

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	GTPγS	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Fluorescence				
GPCRs																	
5-Hydroxytryptamine																	
5-HT _{1A}	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1			X				X	ES-310-A	8-Hydroxy-DPAT, [propyl-2,3-ring-1,2,3- ³ H]-MPPF, [methyl- ³ H]-	5-HT _{1A} agonist 5-HT _{1A} antagonist agonist agonist endogenous agonist antagonist	NET929 NET1109 NET638 NEX199 NET498 NET1071 NET1164		
			AequoZen frozen ready to screen cells	CHO-K1						X	ES-310-AF	Lysergic acid diethylamide, [N-methyl- ³ H]-					
			ValiScreen transfected cell line	CHO-K1		X	X	X				ES-310-C	Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)				
			cAMPZen frozen ready to screen cells	CHO-K1				X				ES-310-CF	5-Carboxamidotryptamine, [1,2- ³ H]-				
			Membrane preparation	CHO-K1		X						6110501400UA	Hydroxytryptamine creatinine sulfate, 5-[1,2- ³ H(N)]-				
			HEK293 EBNA		X							RBHS1AM400UA	5-Carboxamidotryptamine, [1,2- ³ H]-				
	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X	ES-313-A	Ketanserin hydrochloride, [ethylene- ³ H]-, (R41 468)	5-HT _{2A} antagonist 5-HT _{2A} agonist 5-HT _{2A} antagonist agonist agonist agonist endogenous	NET791 NEX255 NET565 NET638 NEX199 NET1071 NET498		
			AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X	X	ES-313-AF	DOI, [¹²⁵ I]-(+)-				
			ValiScreen transfected cell line	CHO-K1		X					X	ES-313-C	Spiperone, [benzene ring- ³ H]-, (Spiroperidol)				
			Membrane preparation	CHO-K1		X						ES-313-M400UA	Lysergic acid diethylamide, [N-methyl- ³ H]-				
													Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)				
													5-Carboxamidotryptamine, [1,2- ³ H]-				
G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X	ES-314-A	Ketanserin hydrochloride, [ethylene- ³ H]-, (R41 468)	5-HT _{2B} antagonist 5-HT _{2B} agonist 5-HT _{2B} antagonist agonist agonist agonist endogenous	NET791 NEX255 NET565 NET638 NEX199 NET1071 NET498			
		AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X	X	ES-314-C	DOI, [¹²⁵ I]-(+)-					
		ValiScreen transfected cell line	CHO-K1		X						ES-314-M400UA	Spiperone, [benzene ring- ³ H]-, (Spiroperidol)					
		Membrane preparation	CHO-K1		X							Lysergic acid diethylamide, [N-methyl- ³ H]-					
												5-Carboxamidotryptamine, [1,2- ³ H]-					
												Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)					
G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X	ES-318-A	Ketanserin hydrochloride, [ethylene- ³ H]-, (R41 468)	5-HT _{2C} antagonist 5-HT _{2C} agonist 5-HT _{2C} antagonist agonist agonist agonist endogenous	NET791 NEX255 NET565 NET638 NEX199 NET1071 NET498 NET1167			
		AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X	X	ES-318-AF	DOI, [¹²⁵ I]-(+)-					
		ValiScreen transfected cell line	1321N1		X			X			ES-318-C	Spiperone, [benzene ring- ³ H]-, (Spiroperidol)					
		Membrane preparation	HEK293		X						6110548400UA	Lysergic acid diethylamide, [N-methyl- ³ H]-					
			1321N1		X						ES-318-M400UA	Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)					
												5-Hydroxytryptamine, [³ H]-					
G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X	ES-315-A	Ketanserin hydrochloride, [ethylene- ³ H]-, (R41 468)	5-HT _{2C} antagonist 5-HT _{2C} agonist 5-HT _{2C} antagonist agonist agonist agonist endogenous	NET791 NEX255 NET565 NET638 NEX199 NET1071 NET498			
		AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X	X	ES-315-AF	DOI, [¹²⁵ I]-(+)-					
		ValiScreen transfected cell line	CHO-K1		X			X		X	ES-315-C	Spiperone, [benzene ring- ³ H]-, (Spiroperidol)					
		Membrane preparation	CHO-K1		X						ES-315-M400UA	Lysergic acid diethylamide, [N-methyl- ³ H]-					
												5-Carboxamidotryptamine, [1,2- ³ H]-					
												Hydroxytryptamine creatinine sulfate, 5-[1,2- ³ H(N)]-					
5-HT _{2A} (ligand-gated ion channel) see Ion Channels																	
G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X	ES-401-A	Lysergic acid diethylamide, [N-methyl- ³ H]-	agonist agonist endogenous agonist agonist	NET638 NEX199 NET498 NET1071			
		AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X		ES-401-AF	Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)					
		ValiScreen transfected cell line	CHO-K1		X						ES-401-C	Hydroxytryptamine creatinine sulfate, 5-[1,2- ³ H(N)]-					
		Membrane preparation	CHO-K1		X						RBHSSAM400UA	5-Carboxamidotryptamine, [1,2- ³ H]-					
			CHO-K1		X							ES-401-M400UA			Lysergic acid diethylamide, [N-methyl- ³ H]-		
												ES-316-A			Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)		
G _s	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X	ES-316-A	Lysergic acid diethylamide, [N-methyl- ³ H]-	agonist agonist endogenous agonist agonist antagonist	NET638 NEX199 NET498 NET1071 NEX424			
		AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X	X	ES-316-AF	Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)					
		ValiScreen transfected cell line	CHO-K1/1321N1		X			X			ES-316-C	Hydroxytryptamine creatinine sulfate, 5-[1,2- ³ H(N)]-					
		cAMPZen frozen ready to screen cells	1321N1		X			X			ES-316-CF	5-Carboxamidotryptamine, [1,2- ³ H]-					
		Membrane preparation	CHO-K1		X						ES-316-M400UA	5-Carboxamidotryptamine, [1,2- ³ H]-					
			HEK293		X						RBHSM400UA	SB258585, [¹²⁵ I]-					
G _s	Rat	Membrane preparation	HEK293		X						RBRS6M400UA						
		Human	Membrane preparation	CHO-K1		X					6110512400UA	Lysergic acid diethylamide, [N-methyl- ³ H]-	agonist agonist agonist endogenous	NET638 NEX199 NET1071 NET498			
			Membrane preparation	HEK293		X					RBRS7M400UA	Lysergic acid diethylamide, 2-[¹²⁵ I]iodo(-)-, ([¹²⁵ I]-LSD)					
	Rat	Membrane preparation	HEK293		X							5-Carboxamidotryptamine, [1,2- ³ H]-					
												Hydroxytryptamine creatinine sulfate, 5-[1,2- ³ H(N)]-					
												5-Carboxamidotryptamine, [1,2- ³ H]-					
Acetylcholine (Muscarinic)																	
M ₁	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X				X	ES-210-A	Scopolamine methyl chloride, [N-methyl- ³ H]-	M ₁ -M ₂ antagonist endogenous M ₁ -M ₂ antagonist M ₁ -M ₂ agonist M ₁ antagonist M ₁ , M ₂ antagonist	NET636 NET113 NET656 NET671 NET780 NET1040			
			AequoZen frozen ready to screen cells	CHO-K1						X	ES-210-AF	Acetylcholine iodide, [acetyl- ³ H]-					
			ValiScreen transfected cell line	CHO-K1		X				X	ES-210-C	Quinuclidinyl benzilate, L-[benzilic-4,4- ³ H]-					
			Membrane preparation	CHO-K1		X						RBHM1M400UA			Oxtremorine-M acetate, [methyl- ³ H]-		
															Pirenzepine, [N-methyl- ³ H]-		
															4-DAMP, [N-methyl- ³ H]-		
M ₂	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X				X	ES-211-A	Scopolamine methyl chloride, [N-methyl- ³ H]-	M ₁ -M ₂ antagonist endogenous M ₁ -M ₂ antagonist M ₁ -M ₂ agonist M ₂ antagonist	NET636 NET113 NET656 NET671 NET1041			
			AequoZen frozen ready to screen cells	CHO-K1						X	ES-211-AF	Acetylcholine iodide, [acetyl- ³ H]-					
			ValiScreen transfected cell line	CHO-K1		X	X					ES-211-C			Quinuclidinyl benzilate, L-[benzilic-4,4- ³ H]-		
			Membrane preparation	CHO-K1		X						RBHM2M400UA			Oxtremorine-M acetate, [methyl- ³ H]-		
															AF-DX 384, [2, 3-dipropylamino- ³ H]-		
															4-DAMP, [N-methyl- ³ H]-		
M ₃	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X				X	ES-212-A	Scopolamine methyl chloride, [N-methyl- ³ H]-	M ₁ -M ₂ antagonist endogenous M ₁ -M ₂ antagonist M ₁ -M ₂ agonist M ₁ , M ₂ antagonist	NET636 NET113 NET656 NET671 NET1040			
			AequoZen frozen ready to screen cells	CHO-K1						X	ES-212-AF	Acetylcholine iodide, [acetyl- ³ H]-					
			ValiScreen transfected cell line	CHO-K1		X				X	ES-212-C	Quinuclidinyl benzilate, L-[benzilic-4,4- ³ H]-					
			Membrane preparation	CHO-K1		X						RBHM3M400UA			Oxtremorine-M acetate, [methyl- ³ H]-		
															4-DAMP, [N-methyl- ³ H]-		
															Scopolamine methyl chloride, [N-methyl- ³ H]-		
M ₄	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	ES-213-A	Scopolamine methyl chloride, [N-methyl- ³ H]-	M ₁ -M ₂ antagonist endogenous M ₁ -M ₂ antagonist M ₁ -M ₂ agonist M ₁ , M ₂ antagonist	NET636 NET113 NET656 NET671			
			AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X	ES-213-AF	Acetylcholine iodide, [acetyl- ³ H]-					
			ValiScreen transfected cell line	CHO-K1		X	X					ES-213-C			Quinuclidinyl benzilate, L-[benzilic-4,4- ³ H]-		
			Membrane preparation	CHO-K1		X						RBHM4M400UA			Oxtremorine-M acetate, [methyl- ³ H]-		
															Scopolamine methyl chloride, [N-methyl- ³ H]-		
															Acetylcholine iodide, [acetyl- ³ H]-		
M ₅	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X				X	ES-214-A	Scopolamine methyl chloride, [N-methyl- ³ H]-	M ₁ -M ₂ antagonist endogenous M ₁ -M ₂ antagonist M ₁ -M ₂ agonist M ₁ , M ₂ antagonist	NET636 NET113 NET656 NET671			
			AequoZen frozen ready to screen cells	CHO-K1						X	ES-214-AF	Acetylcholine iodide, [acetyl- ³ H]-					
			ValiScreen transfected cell line	CHO-K1		X				X	ES-214-C	Quinuclidinyl benzilate, L-[benzilic-4,4- ³ H]-					
			Membrane preparation	CHO-K1		X						RBHM5M400UA			Oxtremorine-M acetate, [methyl- ³ H]-		
															Acetylcholine iodide, [acetyl- ³ H]-		
															Quinuclidinyl benzilate, L-[benzilic-4,4- ³ H]-		

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	qFP70	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Cytosolic Fluorescence				
Adenosine	A ₁	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-010-A	Cyclopentyl-1,3-dipropylxanthine, 8-[dipropyl-2,3- ³ H(N)]-Ethylcarboxamidoadenosine, 5'-N-[adenine-2,8- ³ H]-, (NECA)	A ₁ antagonist A ₁ , A ₂ agonist	NET974 NET811	
				AequoZen frozen ready to screen cells	CHO-K1				X				ES-010-AF				
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-010-C				
				CAMPZen frozen ready to screen cells	CHO-K1				X				ES-010-CF				
				Membrane preparation	CHO-K1		X	X					ES-010-M400UA				
				Membrane preparation	CHO-K1		X						6110511400UA				
	A _{2A}	G _s	Human	AequoScreen aequorin double-transfected cell line	HEK293	Gα ₁₆	X			X	X		ES-011-A	CGS 21680, [carboxyethyl- ³ H(N)]-Ethylcarboxamidoadenosine, 5'-N-[adenine-2,8- ³ H]-, (NECA)	A _{2A} agonist A ₁ , A ₂ agonist	NET1021 NET811	
				AequoZen frozen ready to screen cells	HEK293	Gα ₁₆	X			X			ES-011-AF				
				ValiScreen transfected cell line	HEK293		X		X				ES-011-C				
				CAMPZen frozen ready to screen cells	HEK293		X		X				ES-011-CF				
				Membrane preparation	HEK293		X						RBHA2AM400UA				
				Membrane preparation	HEK293		X						MCL-511				
	A _{2B}	G _s	Human	ValiScreen transfected cell line	CHO-K1		X				X		ES-013-A	CGS 21680, [carboxyethyl- ³ H(N)]-Ethylcarboxamidoadenosine, 5'-N-[adenine-2,8- ³ H]-, (NECA)	A _{2B} agonist A ₁ , A ₂ agonist	NET1021 NET811	
				AequoScreen aequorin double-transfected cell line	HEK293	Gα ₁₆	X			X			ES-013-A				
				ValiScreen transfected cell line	HEK293		X		X				ES-013-C				
				CAMPZen frozen ready to screen cells	HEK293		X		X				ES-013-CF				
				Membrane preparation	HEK293		X						ES-013-M400UA				
				Membrane preparation	HEK293		X						ES-013-M400UA				
A ₃	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₈	X			X	X		ES-012-A	AB-MECA, (4-amino-3-[¹²⁵ I]iodobenzyl-5'-N)-methylcarboxamideadenosine	A ₃ agonist	NEX312		
			AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X			X			ES-012-AF					
			ValiScreen transfected cell line	CHO-K1		X	X	X				ES-012-C					
			CAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-012-CF					
			PhotoScreen Photina double-transfected cell line	CHO-K1	Gα ₁₆	X			X			AX-001-PCF					
			Membrane preparation	CHO-K1		X	X					ES-012-M400UA					
	Rat			ValiScreen transfected cell line	CHO-K1		X						MCL-512				
				Membrane preparation	HEK293 EBNA		X						RBRA3M400UA				
				Membrane preparation	HEK293 EBNA		X						ES-036-A				
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-036-AF				
				AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-036-C				
				Membrane preparation	CHO-K1		X						ES-036-M400UA				
Adrenoceptors	α _{1A}	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X		ES-037-A	Prazosin, [7-methoxy- ³ H]-HEAT, [¹²⁵ I]-Prazosin analog [¹²⁵ I]-iodoazido-, (iodoarylazido)prazosin	α ₁ antagonist α ₁ antagonist α ₁ antagonist	NET823 NEX182 NEX219		
				AequoZen frozen ready to screen cells	CHO-K1		X			X		ES-037-A					
				ValiScreen transfected cell line	CHO-K1		X		X							ES-038-A	
				Membrane preparation	CHO-K1		X									ES-038-A	
				Membrane preparation	HEK293 EBNA		X									ES-030-A	
				Membrane preparation	HEK293 EBNA		X									ES-030-AF	
	α _{1B}	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-030-A	MK-912, [methyl- ³ H]-Iodoclonidine, p-[¹²⁵ I]-2-[(2,5-dichloro-4-[¹²⁵ I] phenyl) imino] imidazolidine	α ₁ antagonist α ₂ agonist α ₂ agonist endogenous α ₂ antagonist endogenous α ₂ antagonist α ₂ agonist α ₂ antagonist	NET1059 NEX253 NET613 NET623 NET659 NET678 NET722 NET853 NET1153	
				AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X			X	X		ES-030-AF				
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-030-C				
				CAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-030-CF				
				Membrane preparation	CHO-K1		X	X	X				ES-030-M400UA				
				Membrane preparation	Sf9		X						6110113400UA				
	α _{2A}	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-031-A	MK-912, [methyl- ³ H]-Iodoclonidine, p-[¹²⁵ I]-2-[(2,5-dichloro-4-[¹²⁵ I] phenyl) imino] imidazolidine	α ₂ antagonist α ₂ agonist α ₂ agonist endogenous α ₂ antagonist endogenous α ₂ antagonist α ₂ agonist α ₂ antagonist	NET1059 NEX253 NET613 NET623 NET659 NET678 NET722 NET853 NET1153	
				AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-031-AF				
				ValiScreen transfected cell line	CHO-K1		X	X					ES-031-C				
				Membrane preparation	CHO-K1		X	X					ES-031-M400UA				
				Membrane preparation	HEK293 EBNA		X						ES-032-A				
				Membrane preparation	HEK293 EBNA		X						ES-032-AF				
α _{2C}	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-032-A	MK-912, [methyl- ³ H]-Iodoclonidine, p-[¹²⁵ I]-2-[(2,6-dichloro-4-[¹²⁵ I]iodophenyl)imino]imidazolidine	α ₂ antagonist α ₂ agonist α ₂ agonist endogenous α ₂ antagonist endogenous α ₂ antagonist α ₂ agonist α ₂ antagonist	NET1059 NEX253 NET613 NET623 NET659 NET678 NET722 NET853 NET1153		
			ValiScreen transfected cell line	CHO-K1		X		X				ES-032-C					
			CAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-032-CF					
			Membrane preparation	Sf9		X						6110114400UA					
			Membrane preparation	CHO-K1		X						ES-032-M400UA					
			Membrane preparation	HEK293 EBNA		X						ES-033-A					
β ₁	G _s	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X		ES-033-A	(-)-CGP 12177, [5,7- ³ H]-(-)-Cyanopindolol, [¹²⁵ I]-(-)-Pindolol, [¹²⁵ I]-Propranolol, L-[4- ³ H]-Dihydroalprenolol hydrochloride, levo [ring, propyl- ³ H]-(+/-)-Cyanopindolol, [¹²⁵ I]-	antagonist antagonist antagonist antagonist antagonist antagonist	NET1061 NEX189 NEX211 NET515 NET720 NEX174		
			AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X			X			ES-033-AF					
			ValiScreen transfected cell line	CHO-K1		X		X				ES-033-C					
			CAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-033-CF					
			Membrane preparation	CHO-K1		X						ES-033-M400UA					
			Membrane preparation	Sf9		X						6110110400UA					

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Biotinylation	GFPG	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Calcium Fluorescence				
Adrenoceptors (continued)	β ₂	G _s	Human	ValiScreen transfected cell line	CHO-K1		X	X					ES-034-C	(-)-CGP 12177, [5,7- ³ H]-	antagonist	NET1061	
				cAMPZen frozen ready to screen cells	CHO-K1		X	X					ES-034-CF	(-)-Cyanopindolol, [1 ²⁵ I]-	antagonist	NEX189	
				Membrane preparation	CHO-K1		X						ES-034-M400UA	(-)-Pindolol, [1 ²⁵ I]-	antagonist	NEX211	
					SI9		X						6110106400UA	Propranolol, L-[4- ³ H]-	antagonist	NET515	
					HEK293 EBNA		X						RBHBE2M400UA	Dihydroalprenolol hydrochloride, levo [ring, propyl- ³ H]-	antagonist	NET720	
							X							(±)-Cyanopindolol, [1 ²⁵ I]-	antagonist	NEX174	
	β ₃	G _s	Human	ValiScreen transfected cell line	CHO-K1		X	X					ES-035-C	(-)-CGP 12177, [5,7- ³ H]-	antagonist	NET1061	
				cAMPZen frozen ready to screen cells	CHO-K1		X	X					ES-035-CF	(-)-Cyanopindolol, [1 ²⁵ I]-	antagonist	NEX189	
				Membrane preparation	CHO-K1		X						ES-035-M400UA	(-)-Pindolol, [1 ²⁵ I]-	antagonist	NEX211	
					SK-N-MC		X						RBHBE3M400UA	Propranolol, L-[4- ³ H]-	antagonist	NET515	
							X							Dihydroalprenolol hydrochloride, levo [ring, propyl- ³ H]-	antagonist	NET720	
							X							(±)-Cyanopindolol, [1 ²⁵ I]-	antagonist	NEX174	
Anaphylatoxin	C3a	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X		ES-730-A	Complement C3a, [1 ²⁵ I]-, Bolton-Hunter labeled	C3a endogenous	NEX356		
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-730-C				
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-730-CF				
				Membrane preparation	CHO-K1		X	X					ES-730-M400UA				
					CHO-K1		X			X			ES-731-A				
					CHO-K1		X				X		ES-731-AF				
	C5a	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X		ES-731-A	Complement C5a, [1 ²⁵ I]-, Bolton-Hunter labeled	C5a endogenous	NEX250		
				AequoZen frozen ready to screen cells	CHO-K1		X				X	ES-731-AF					
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-731-C				
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-731-CF				
				Membrane preparation	CHO-K1		X	X					ES-731-M400UA				
					CHO-K1		X						6110526400UA				
Angiotensin	AT ₁	G _{q/11} , G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X	ES-072-A	Angiotensin II (Sar ¹ , Tyr ⁸ , Ile ⁸), [1 ²⁵ I]-	AT antagonist	NEX248		
				AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X			X			ES-072-AF	Angiotensin I (Tyr ¹), [1 ²⁵ I]-	endogenous	NEX101	
				ValiScreen transfected cell line	CHO-K1		X		X				ES-072-C	Angiotensin IV (Tyr ³), [1 ²⁵ I]-	Non AT ₁ , non AT ₂ agonist	NEX295	
				Membrane preparation	SI9		X						6110121400UA	Angiotensin II (Tyr ¹), [1 ²⁵ I]-	endogenous	NEX105	
					CHO-K1		X						ES-072-M400UA				
					CHO-K1		X						ES-070-C	Angiotensin II (Sar ¹ , Tyr ⁸ , Ile ⁸), [1 ²⁵ I]-	AT antagonist	NEX248	
	AT ₂	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X					ES-070-M400UA	Angiotensin IV (Tyr ³), [1 ²⁵ I]-	Non AT ₁ , non AT ₂ agonist	NEX295		
				Membrane preparation	CHO-K1		X					ES-071-C	Angiotensin II (Tyr ¹), [1 ²⁵ I]-	endogenous	NEX105		
					CHO-K1		X						ES-071-M400UA	CGP42112A, [1 ²⁵ I]-	AT ₂ antagonist	NEX324	
					CHO-K1		X						ES-460-A	Apelin 13 (Glp ⁶⁹ , Nle ⁷⁵ , Tyr ⁷⁷), [1 ²⁵ I]-	APJ agonist	NEX393	
					CHO-K1		X	X	X				ES-460-C				
					CHO-K1		X		X				ES-460-CF				
Beta Alanine	MRGPRD, TGR7	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X	ES-741-A					
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-460-C				
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-460-CF				
				Membrane preparation	CHO-K1		X	X					ES-460-M400UA				
					CHO-K1		X						ES-741-A				
					CHO-K1		X						ES-581-A	Bombesin (Tyr ⁴), [1 ²⁵ I]-	agonist	NEX258	
	Bombesin	BB ₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X	ES-581-A	Bombesin (Tyr ⁴), [1 ²⁵ I]-	agonist	NEX258	
					Membrane preparation	CHO-K1	Gα ₁₆	X					RBHBS1M400UA				
						CHO-K1	Gα ₁₆	X			X	X		ES-582-A	Bombesin (Tyr ⁴), [1 ²⁵ I]-	agonist	NEX258
		BB ₂	G _i /G ₁₁	Human	AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X				X	ES-582-AF				
					ValiScreen transfected cell line	CHO-K1		X			X			ES-582-C	Gastrin Releasing Peptide (Tyr ¹⁵) (Porcine), [125I]-	agonist	NEX421
					Membrane preparation	HEK293		X						RBHBS2M400UA			
Bradykinin	BB ₃	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X		ES-580-A	Bombesin (6-14) (D-Tyr ⁶ , βAla ¹¹ , Phe ¹³ , Nle ¹⁴), [1 ²⁵ I]-	BB ₃ agonist	NEX377		
				Membrane preparation	CHO-K1	Gα ₁₆	X						RBHBS3M400UA	Bombesin (Tyr ⁴), [1 ²⁵ I]-	agonist	NEX258	
					CHO-K1		X			X	X		ES-091-A	Kalidin (Des-Arg ¹⁰), [3,4-prolyl-3,4- ³ H(N)]-	B ₁ agonist	NET1064	
	B ₁	G _i /G ₁₁	Rat	ValiScreen transfected cell line	CHO-K1		X				X	ES-091-C	Kalidin (Des-Arg ¹⁰ , Leu ⁷), [3,4-prolyl-3,4- ³ H(N)]-	B ₁ antagonist	NET1096		
				Membrane preparation	CHO-K1		X					ES-091-M400UA					
					CHO-K1		X						ES-092-A				
B ₂	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X				X	ES-090-A	Bradykinin, [2,3-prolyl-3,4- ³ H]-	endogenous	NET706			
			ValiScreen transfected cell line	CHO-K1		X					ES-090-C						
			Membrane preparation	CHO-K1		X						ES-090-M400UA					
Calcitonin	CGRP ₁	G _s	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{9s}	X			X	X	ES-420-A	Calcitonin Gene Related Peptide, [1 ²⁵ I]-, ([1 ²⁵ I]-hCGRP)	CGPR ₁ , endogenous	NEX354		
				AequoZen frozen ready to screen cells	CHO-K1	Gα _{9s}	X				X	ES-420-AF	Calcitonin Gene Related Peptide (8-37), [1 ²⁵ I]-, ([1 ²⁵ I]-hCGRP)	CGPR ₁ , antagonist	NEX319		
				ValiScreen transfected cell line	CHO-K1		X		X				ES-420-C	Amylin (r) [125I]-	CGPR ₁ , antagonist	NEX448	
				cAMPZen frozen ready to screen cells	CHO-K1		X						ES-420-CF				
				Membrane preparation	CHO-K1		X						ES-420-M400UA				
					SI9	Gα _{βγ2}	X						6110135400UA				
	CT	G _s	Human	Non-recombinant, membrane preparation	SK-N-MC		X						RBHGRPM400UA				
							X							Calcitonin (Tyr ¹³) (human), [1 ²⁵ I]-	endogenous	NEX422	
							X							Calcitonin (salmon), [1 ²⁵ I]-	endogenous	NEX423	
							X							6110136400UA			
							X							Adrenomedullin (13-52), (Tyr), [1 ²⁵ I]-	AM ₁ , AM ₂ endogenous	NEX385	
							X							Adrenomedullin (rat), [1 ²⁵ I]-	AM ₁ , AM ₂ endogenous	NEX427	
AM ₁	G _s	Human	Membrane preparation	SI9	Gα _{βγ2}	X						6110136400UA					
						X							Adrenomedullin (13-52), (Tyr), [1 ²⁵ I]-	AM ₁ , AM ₂ endogenous	NEX385		
						X							Adrenomedullin (rat), [1 ²⁵ I]-	AM ₁ , AM ₂ endogenous	NEX427		
						X							ES-430-A				
						X							ES-430-C				
						X							ES-430-CF				
AM ₂	G _s , G ₁₀	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-430-A	Adrenomedullin (13-52), (Tyr), [1 ²⁵ I]-	AM ₁ , AM ₂ endogenous	NEX385		
			ValiScreen transfected cell line	CHO-K1		X			X			ES-430-C	Adrenomedullin (rat), [1 ²⁵ I]-	AM ₁ , AM ₂ endogenous	NEX427		
			cAMPZen frozen ready to screen cells	CHO-K1		X						ES-430-CF					
			Membrane preparation	CHO-K1		X						ES-430-M400UA					
						X											
						X											

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	GRTPG	cAMP	Inositol Phosphate	Calcium	Luminescence	Fluorescence				
Cannabinoid	CB ₁	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{i,16}				X	X		ES-110-A	CP 55940, [side chain-2,3,4- ³ H(N)]-WIN 55,212-2, [5,7-naphthyl- ³ H]-Anadamide, [arachidonyl-5,6,8,9,11,12,14,15- ³ H]-SR141716A, [³ H]	agonist agonist CB, endogenous CB, antagonist	NET1051 NET1058 NET1073 NET1158	
				AequoZen frozen ready to screen cells	CHO-K1	Gα _{i,16}				X			ES-110-AF				
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-110-C				
				cAMPZen frozen ready to screen cells	CHO-K1				X				ES-110-CF				
				Membrane preparation	CHO-K1		X	X					ES-110-M400UA				
					SH9	Gα _q β _{1,2}	X						6110129400UA				
	CB ₂	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-111-C	CP 55940, [side chain-2,3,4- ³ H(N)]-WIN 55,212-2, [5,7-naphthyl- ³ H]-Anadamide, [arachidonyl-5,6,8,9,11,12,14,15- ³ H]-	agonist agonist CB, endogenous	NET1051 NET1058 NET1073	
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-111-CF				
				Membrane preparation	SH9	Gα _q β _{1,2}	X						6110130400UA				
					HEK293 EBNA		X						RBXC2M400UA				
					CHO-K1		X	X					ES-111-M400UA				
					CHO-K1				X				ES-661-A				
Chemerin	GPR1	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{i,16}					X	ES-132-C					
Chemokine	CCR1	G _i /G _o	Human	AequoZen frozen ready to screen cells	CHO-K1	Gα _{i,16}					X		ES-132-C	MIP-1α, [¹²⁵ I]-, Bolton-Hunter labeled RANTES, [¹²⁵ I]-	CCR1, CCR4, CCR5, endogenous CCR1, CCR3, CCR4, CCR5, endogenous	NEX298 NEX292	
				ValiScreen transfected cell line	CHO-K1		X	X					ES-132-CF				
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-132-CF				
				Membrane preparation	HEK293		X						RBHCC1M400UA				
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-133-A				
				AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-133-AF				
	CCR2b	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-133-C	MCP-1, [¹²⁵ I]-, Bolton-Hunter labeled JE, [¹²⁵ I]-, Bolton-Hunter labeled (murine, recombinant MCP-1)	CCR2, endogenous CCR2 agonist	NEX332 NEX313	
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-133-C				
				Membrane preparation	CHO-K1		X	X					ES-133-M400UA				
					HEK293		X						6110550400UA				
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-138-A				
				AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-138-AF				
	CCR3	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1	Gα _{i,16}						X	ES-138-C	Eotaxin, [¹²⁵ I]-	CCR3, endogenous	NEX314	
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-138-C				
				Membrane preparation	CHO-K1		X	X					ES-138-M400UA				
					HEK293		X										
					CHO-K1		X						ES-139-A				
					CHO-K1		X						ES-139-AF				
	CCR4	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-139-C	TARC, [¹²⁵ I]- MIP-1β (Leu ³ Gly ²), [¹²⁵ I]- MIP-3α, [¹²⁵ I]-	endogenous endogenous CCR6, endogenous	NEX369 NEX299 NEX371	
				cAMPZen frozen ready to screen cells	CHO-K1		X	X	X				ES-139-CF				
				Membrane preparation	CHO-K1		X	X					ES-139-M400UA				
					CHO-K1		X			X	X		ES-140-A				
				AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{i,16}				X	X		ES-140-AF				
				AequoZen frozen ready to screen cells	CHO-K1	Gα _{i,16}				X			ES-140-AF				
CCR7	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-140-C	MIP-3β, [¹²⁵ I]- 6-Ckine, [¹²⁵ I]-, Bolton-Hunter labeled	CCR7, endogenous CCR7, endogenous	NEX370 NEX387		
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-140-CF					
			Membrane preparation	CHO-K1		X	X					ES-140-M400UA					
				CHO-K1		X			X			ES-136-A					
			AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-136-AF					
			AequoZen frozen ready to screen cells	CHO-K1		X	X	X				ES-136-C					
CCR8	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-136-CF	I309, [¹²⁵ I]-, Bolton-Hunter labeled	CCR8, endogenous	NEX364		
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-136-CF					
			Membrane preparation	CHO-K1		X	X					ES-136-M400UA					
				CHO-K1		X			X			ES-146-AF					
			AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-143-A					
			AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-143-AF					
CCR9a	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X					X	ES-143-AF	CTACK, [¹²⁵ I]-	CCR9, endogenous	NEX391		
			cAMPZen frozen ready to screen cells	CHO-K1		X						ES-143-AF					
			Membrane preparation	CHO-K1		X	X					ES-143-AF					
				CHO-K1		X			X	X		ES-137-A					
			AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{i,16}				X	X		ES-137-AF					
			AequoZen frozen ready to screen cells	CHO-K1	Gα _{i,16}				X			ES-137-AF					
CCR10	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-137-C	Fractalkine, [¹²⁵ I]-	CX ₂ CR1, endogenous	NEX368		
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-137-CF					
			Membrane preparation	CHO-K1		X	X					ES-137-M400UA					
				CHO-K1		X			X	X		ES-145-A					
			AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{i,16}				X	X		ES-145-AF					
			AequoZen frozen ready to screen cells	CHO-K1	Gα _{i,16}				X			ES-145-AF					
CX ₂ CR1	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-145-C	Interleukin-8, [¹²⁵ I]-, [¹²⁵ I]-IL-8 GRO Alpha/MGSA, [¹²⁵ I]-	CXCR1, CXCR2, endogenous CXCR2, endogenous	NEX277 NEX321		
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-145-CF					
			Membrane preparation	CHO-K1		X	X					ES-145-M400UA					
				CHO-K1		X			X	X		ES-142-A					
			AequoScreen aequorin double-transfected cell line	CHO-K1	Gα _{i,16}				X			ES-142-AF					
			AequoZen frozen ready to screen cells	CHO-K1	Gα _{i,16}				X			ES-142-AF					
CXCR3	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-142-C	I-TAC, [¹²⁵ I]- IP-10, [¹²⁵ I]-	CXCR3, endogenous CXCR3, endogenous	NEX376 NEX348		
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-142-C					
			Membrane preparation	CHO-K1		X	X					ES-142-M400UA					
				CHO-K1		X			X			ES-720-A					
			AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-720-AF					
			AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-720-AF					
CXCR4	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X					ES-720-C	SDF1α, [¹²⁵ I]- CXCL16, [¹²⁵ I]-	CXCR4, endogenous CXCR6, endogenous	NEX346 NEX398		
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-720-CF					
			Membrane preparation	CHO-K1		X	X					ES-720-M400UA					
				CHO-K1		X			X			ES-720-M400UA					
			AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			RBHCX6M400UA					
			AequoZen frozen ready to screen cells	HEK293		X			X			RBHCX6M400UA					

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	GRTPG	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Fluorescence				
Chemokine (continued)	XCR1	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆				X			ES-148-A				
				AequoZen frozen ready to screen cells	CHO-K1	Gd ₁₆				X				ES-148-AF			
				ValiScreen transfected cell line	CHO-K1		X	X						ES-148-C			
				GAMPZen frozen ready to screen cells	CHO-K1			X						ES-148-CF			
Cholecystokinin	CCK ₁	G _q /G ₁₁	Human	Membrane preparation	CHO-K1		X						ES-148-M400UA				
				AequoScreen aequorin double-transfected cell line	1321N1		X			X	X		ES-530-A	Cholecystokinin octapeptide, [¹²⁵ I]-, Bolton-Hunter labeled, ([¹²⁵ I]-CCK-8)	endogenous	NEX203	
				AequoZen frozen ready to screen cells	1321N1		X			X			ES-530-AF	Gastrin I (Tyr ³), [¹²⁵ I]-	endogenous	NEX176	
	CCK ₂	G _q /G ₁₁	Human	ValiScreen transfected cell line	1321N1		X			X		ES-530-C	Interleukin-8 (IL-8) (h,r) Receptor Grade, [¹²⁵ I]	agonist	NEX443		
				Membrane preparation	1321N1		X			X	X		ES-530-M400UA	CCK-8, [³ H]	agonist	NET1162	
				AequoScreen aequorin double-transfected cell line	1321N1		X			X	X		ES-531-A	Cholecystokinin octapeptide, [¹²⁵ I]-, Bolton-Hunter labeled, ([¹²⁵ I]-CCK-8)	endogenous	NEX203	
Corticotropin-Releasing Factor	CRF ₁	G _o	Human	AequoZen frozen ready to screen cells	1321N1		X			X		ES-531-AF	Gastrin I (Tyr ³), [¹²⁵ I]-	endogenous	NEX176		
				ValiScreen transfected cell line	1321N1		X			X			ES-531-C	Interleukin-8 (IL-8) (h,r) Receptor Grade, [¹²⁵ I]	agonist	NEX443	
				Membrane preparation	1321N1		X			X			ES-531-M400UA	CCK-8, [³ H]	agonist	NET1162	
				AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆	X			X	X		ES-152-A	Sauvagine (Tyr ⁶), [¹²⁵ I]-	endogenous	NEX306	
				ValiScreen transfected cell line	CHO-K1		X	X					ES-152-C	CRF (Tyr ⁶) (ovine), [¹²⁵ I]-	CRF agonist	NEX217	
				Membrane preparation	HEK293		X						RBNCRF1M400UA	CRF (Tyr ⁶) (human,rat), [¹²⁵ I]-	endogenous	NEX216	
Dopamine	D ₁	G _o	Human	ValiScreen transfected cell line	CHO-K1		X		X			ES-172-C	SCH 23390, [N-methyl- ³ H]-	D ₁ , D ₂ antagonist	NET930		
				CAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-172-CF	Dihydroxyphenylethylamine, 3,4-[ring-2,5,6- ³ H]-, (Dopamine)	endogenous	NET673	
				Membrane preparation	L Cells		X						6110513400UA	Dihydroxyphenylethylamine, 3,4-[7- ³ H]-	endogenous	NET131	
	D ₂	G _q /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆				X		ES-171-A	Dihydroxyphenylethylamine [2,5,6,7,8- ³ H] (Dopamine)	endogenous	NET1160		
				AequoZen frozen ready to screen cells	CHO-K1	Gd ₁₆				X			ES-171-AF	Iodosipiperone, 2'-[¹²⁵ I]-	D ₂ antagonist	NEX284	
				Membrane preparation	SI9	Gd ₁₆	X						6110137400UA	R (+) trans-7-hydroxy-PIPAT, [¹²⁵ I]-	D ₂ , D ₃ agonist	NEX307	
														Spiroperone, [benzene ring- ³ H]-, (Spiroperidol)	D ₂ antagonist	NET565	
														Dihydroxyphenylethylamine, 3,4-[ring-2,5,6- ³ H]-, (Dopamine)	endogenous	NET673	
														Iodosulpride, [¹²⁵ I]-	D ₂ , D ₃ antagonist	NEX441	
														(-)Sulpride, [methoxy- ³ H]-	D ₂ antagonist	NET775	
														Methylspiperone, [N-methyl- ³ H]-	D ₂ , D ₃ antagonist	NET856	
														Raclopride, [methoxy- ³ H]-	D ₂ antagonist	NET975	
														YM-09151-2, [N-methyl- ³ H]-	D ₂ antagonist	NET1004	
														Dihydroxyphenylethylamine, 3,4-[7- ³ H]-	endogenous	NET131	
														Amisulpride (DAN-2163), [¹²⁵ I]-	D ₂ , D ₃ antagonist	NEX432	
														Domperidone, [benzene ring- ³ H]-	D ₂ , D ₃ antagonist	NET662	
														Quinpirole, [N-propyl- ³ H]-	D ₂ , D ₃ agonist	NET1085	
				D _{2S}	G _q /G _o	Human	Membrane preparation	CHO-K1		X						RBHD2CM400UA	Iodosipiperone, 2'-[¹²⁵ I]-
													R (+) trans-7-hydroxy-PIPAT, [¹²⁵ I]-	D ₂ , D ₃ agonist	NEX307		
													Spiroperone, [benzene ring- ³ H]-, (Spiroperidol)	D ₂ antagonist	NET565		
													Dihydroxyphenylethylamine, 3,4-[ring-2,5,6- ³ H]-, (Dopamine)	endogenous	NET673		
													Iodosulpride, [¹²⁵ I]-	D ₂ , D ₃ antagonist	NEX441		
													(-)Sulpride, [methoxy- ³ H]-	D ₂ antagonist	NET775		
													Methylspiperone, [N-methyl- ³ H]-	D ₂ , D ₃ antagonist	NET856		
													Raclopride, [methoxy- ³ H]-	D ₂ , D ₃ antagonist	NET975		
													YM-09151-2, [N-methyl- ³ H]-	D ₂ antagonist	NET1004		
													Dihydroxyphenylethylamine, 3,4-[7- ³ H]-	endogenous	NET131		
													Amisulpride (DAN-2163), [¹²⁵ I]-	D ₂ , D ₃ antagonist	NEX432		
													Domperidone, [benzene ring- ³ H]-	D ₂ , D ₃ antagonist	NET662		
D ₃	G _q /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X	X				ES-173-C	Dihydroxyphenylethylamine, 3,4-[ring-2,5,6- ³ H]-, (Dopamine)	endogenous	NET673		
			Membrane preparation	CHO-K1		X	X					ES-173-M400UA	Dihydroxyphenylethylamine, 3,4-[7- ³ H]-	endogenous	NET131		
			Rat	Membrane preparation	SI9		X						6110139400UA	R (+) trans-7-hydroxy-PIPAT, [¹²⁵ I]-	D ₂ , D ₃ agonist	NEX307	
														Iodosulpride, [¹²⁵ I]-	D ₂ , D ₃ antagonist	NEX441	
														Methylspiperone, [N-methyl- ³ H]-	D ₂ , D ₃ antagonist	NET856	
														Amisulpride (DAN-2163), [¹²⁵ I]-	D ₂ , D ₃ antagonist	NEX432	
										Domperidone, [benzene ring- ³ H]-	D ₂ , D ₃ antagonist	NET662					
										Quinpirole, [N-propyl- ³ H]-	D ₂ , D ₃ agonist	NET1085					
D ₄	G _q /G _o	Rat	ValiScreen transfected cell line	CHO-K1		X		X				ES-170-C	Dihydroxyphenylethylamine, 3,4-[ring-2,5,6- ³ H]-, (Dopamine)	endogenous	NET673		
			Membrane preparation	SI9		X						6110112400UA	Dihydroxyphenylethylamine, 3,4-[7- ³ H]-	endogenous	NET131		
D _{4L2}	G _q /G _o	Human	Membrane preparation	CHO-K1		X						RBHD42M400UA	WAY100635 [METHOXY-3H]				
			ValiScreen transfected cell line	CHO-K1		X						MCL 507					
			Membrane preparation	SI9		X						6110122400UA					
D _{4L4}	G _q /G _o	Human	ValiScreen transfected cell line	CHO-K1		X						RBHD44M400UA					
			Membrane preparation	CHO-K1		X						RBHD47M400UA					
D _{4L7}	G _q /G _o	Human	Membrane preparation	CHO-K1		X							Dihydroxyphenylethylamine, 3,4-[ring-2,5,6- ³ H]-, (Dopamine)	endogenous	NET673		
														SCH 23390, [N-methyl- ³ H]-	D ₁ , D ₂ antagonist	NET930	
D ₅	G _q /G _o	Human	Membrane preparation	CHO-K1		X							Dihydroxyphenylethylamine, 3,4-[7- ³ H]-	endogenous	NET131		

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	qTPyB	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Fluorescence				
Endothelin	ET _A	G ₁₂ /G ₁₃	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-320-A	Endothelin-1 (Tyr ¹³), [¹²⁵ I]-	endogenous	NEX259	
				AequoZen frozen ready to screen cells	CHO-K1				X		ES-320-AF						
				ValiScreen transfected cell line	CHO-K1				X		ES-320-C						
		ET _B	G ₁₂ /G ₁₃	Human	Membrane preparation	CHO-K1		X					ES-320-M400UA				
	AequoScreen aequorin double-transfected cell line				CHO-K1		X		X	X	ES-321-A						
	AequoZen frozen ready to screen cells				CHO-K1		X		X		ES-321-AF						
	Formyl Peptide	G ₁₂ /G ₁₃	Human	ValiScreen transfected cell line	CHO-K1		X			X		ES-321-C	Endothelin-1 (Tyr ¹³), [¹²⁵ I]-	endogenous	NEX259		
Membrane preparation				CHO-K1		X				ES-321-M400UA							
AequoScreen aequorin double-transfected cell line				CHO-K1		X		X	X	ES-810-A							
	Free Fatty Acid	G ₁₂ /G ₁₃	Human	AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X			X	X		ES-810-AF	WKYMVM, [¹²⁵ I]-, Bolton-Hunter labeled	FPRL1 agonist	NEX386	
ValiScreen transfected cell line				CHO-K1		X	X	X		X	ES-810-C						
cAMPZen frozen ready to screen cells				CHO-K1		X	X	X		X	ES-810-CF						
Membrane preparation		CHO-K1		X	X	X		X	ES-810-M400UA								
AequoScreen aequorin double-transfected cell line		CHO-K1		X				X	ES-852-A								
AequoZen frozen ready to screen cells		CHO-K1		X				X	ES-852-AF								
	Free Fatty Acid	G ₁₂ /G ₁₃	Human	ValiScreen transfected cell line	1321N1				X			ES-852-C					
AequoScreen aequorin double-transfected cell line				CHO-K1		X			X	ES-800-A							
AequoZen frozen ready to screen cells				CHO-K1		X			X	ES-800-AF							
	GABA _A	GABA _{α1a}	Human	Membrane preparation	CHO-K1		X					6110545400UA	Baclofen, (-)-[butyl-4- ³ H(N)]-aminobutyric acid, γ-[2,3- ³ H(N)]-aminobutyric acid, γ-[2,3- ³ H(N)]-aminobutyric acid, γ-[2,3- ³ H(N)]-aminobutyric acid, tertiary-[¹²⁵ S]-Butyl bicyclophosphorothionate, tertiary-[¹²⁵ S]-	GABA _A agonist endogenous endogenous Cl-channel non competitive antagonist	NET867 NET191 NET191X NEG049		
ValiScreen transfected cell line				HEK293		X				6110560400UA							
Membrane preparation				CHO-K1		X				6110546400UA							
AequoScreen aequorin double-transfected cell line		HEK293		X				6110557400UA									
ValiScreen transfected cell line		CHO-K1	Gα ₁₆	X			X	X	ES-500-A								
cAMPZen frozen ready to screen cells		CHO-K1		X	X	X			ES-500-C								
	GABA _A	GABA _{α1b}	Human	Membrane preparation	CHO-K1		X					6110559400UA	Galanin (porcine), [¹²⁵ I]- Galanin (human), [¹²⁵ I]- Galanin (2-11), [¹²⁵ I]-	agonist endogenous GAL ₂ , GAL ₃ agonist	NEX243 NEX333 NEX416		
ValiScreen transfected cell line				HEK293		X				ES-500-CF							
Membrane preparation				CHO-K1		X	X				ES-500-M400UA						
AequoScreen aequorin double-transfected cell line		HEK293		X				X	6110556400UA								
ValiScreen transfected cell line		CHO-K1	Gα ₁₆	X			X	X	ES-510-A								
cAMPZen frozen ready to screen cells		CHO-K1		X	X	X			ES-510-AF								
	Galanin	GAL ₁	Human	ValiScreen transfected cell line	CHO-K1	Gα ₁₆	X	X	X			ES-510-C	Galanin (porcine), [¹²⁵ I]- Galanin (human), [¹²⁵ I]- Galanin (2-11), [¹²⁵ I]-	agonist endogenous GAL ₂ , GAL ₃ agonist	NEX243 NEX333 NEX416		
Membrane preparation				CHO-K1		X					ES-510-CF						
AequoScreen aequorin double-transfected cell line				HEK293 EBNA		X					6110537400UA						
ValiScreen transfected cell line		CHO-K1	Gα ₁₆	X	X				ES-510-M400UA								
Membrane preparation		CHO-K1		X			X	X	ES-511-A								
AequoZen frozen ready to screen cells		CHO-K1		X			X		ES-511-AF								
	Ghrelin	Ghrelin	Human	ValiScreen transfected cell line	CHO-K1		X	X	X			ES-511-C	Ghrelin (His), [¹²⁵ I]- Ghrelin (Tyr), [¹²⁵ I]-	endogenous agonist	NEX388 NEX417		
Membrane preparation				CHO-K1		X					AX-008-PCF						
AequoScreen aequorin double-transfected cell line				CHO-K1		X			X	X	ES-511-M400UA						
AequoZen frozen ready to screen cells		CHO-K1		X			X	X	ES-410-A								
ValiScreen transfected cell line		CHO-K1		X			X	X	ES-410-AF								
Membrane preparation		HEK293		X					ES-410-C								
	Glucagon	Glucagon	Human	ValiScreen transfected cell line	CHO-K1		X					ES-410-M400UA	Glucagon, [¹²⁵ I]-	endogenous	NEX207		
Membrane preparation				1321N1		X		X			ES-710-C						
AequoScreen aequorin double-transfected cell line				CHO-K1		X			X		ES-710-A						
ValiScreen transfected cell line		HEK293		X					RBHGIPM400UA								
Membrane preparation		1321N1		X					ES-711-C								
AequoScreen aequorin double-transfected cell line		1321N1		X					ES-711-CF								
	Glutamate	mGlu _{1A}	Human	ValiScreen transfected cell line	CHO-K1		X					ES-712-A	Quisqualic Acid, [³ H]	mGlu ₁ , mGlu ₂ agonist	NET1165		
Membrane preparation				CHO-K1		X					ES-712-AF						
AequoScreen aequorin double-transfected cell line				CHO-K1		X		X	X		ES-555-A						
ValiScreen transfected cell line		CHO-K1		X					ES-555-M400UA								
Membrane preparation		CHO-K1		X		X			ES-553-C								
AequoScreen aequorin double-transfected cell line		CHO-K1		X		X			ES-553-CF								
	Glycoprotein Hormone	TSH	Human	ValiScreen transfected cell line	CHO-K1	Gα ₁₆	X			X		ES-790-A	Luteinizing Hormone-Releasing Hormone (D-Trp ⁶), [¹²⁵ I]-, (1 ²⁵ I)-[D-Trp ⁶]-LHRH Luteinizing Hormone Releasing Hormone (Tyr ⁶), [¹²⁵ I]-	GnRH agonist agonist	NEX365 NEX163		
Membrane preparation				CHO-K1		X					ES-600-A						
AequoScreen aequorin double-transfected cell line				CHO-K1		X					ES-600-M400UA						
ValiScreen transfected cell line		CHO-K1		X					ES-744-A								
Membrane preparation		CHO-K1		X					ES-744-AF								
AequoScreen aequorin double-transfected cell line		CHO-K1		X					ES-800-A								
	GPR	SUCNR1 (Succinate, GPR91)	Human	ValiScreen transfected cell line	CHO-K1		X					ES-800-AF					
Membrane preparation				CHO-K1		X					ES-800-A						
AequoScreen aequorin double-transfected cell line				CHO-K1		X					ES-743-A						
ValiScreen transfected cell line		CHO-K1		X					ES-743-AF								
Membrane preparation		CHO-K1		X					ES-743-AF								
AequoScreen aequorin double-transfected cell line		CHO-K1		X					ES-743-AF								
	Growth Hormone-Releasing Factor	H ₁	Human	ValiScreen transfected cell line	CHO-K1		X		X	X		ES-390-A	GRF 1-44 amide, [¹²⁵ I]- Pyrimidine, [pyridinyl 5- ³ H]-, (Mepyrmine) Histamine dihydrochloride, [ring, methylenes- ³ H(N)]-	agonist H ₁ antagonist endogenous	NEX426 NET594 NET732		
Membrane preparation				CHO-K1		X					ES-390-AF						
AequoScreen aequorin double-transfected cell line				CHO-K1		X					ES-390-C						
ValiScreen transfected cell line		CHO-K1		X					ES-390-M400UA								
Membrane preparation		CHO-K1		X					ES-390-M400UA								
AequoScreen aequorin double-transfected cell line		CHO-K1		X					ES-390-M400UA								

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	GTPγS	cAMP	Inositol Phosphate	Calcium	Luminescence	Ca ²⁺ Fluorescence				
Histamine (continued)	H ₂	G _s	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X	ES-391-A	Tiotidine, [methyl- ³ H]-, (ICI 125,211) Histamine dihydrochloride, [ring, methylenes- ³ H(N)]- Aminopotentidine, [¹²⁵ I]-	H ₂ antagonist endogenous H ₂ antagonist	NET688 NET732 NEX431		
				AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆				X	ES-391-AF						
				ValiScreen transfected cell line	CHO-K1		X		X		ES-391-C						
				CAMPZen frozen ready to screen cells	CHO-K1		X		X		ES-391-CF						
					Membrane preparation	HEK293		X				6110565400UA					
						CHO-K1		X				ES-391-M400UA					
	H ₃	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X	ES-392-A	Methylhistamine dihydrochloride, N-α-[methyl- ³ H]- Proxylan, [¹²⁵ I]-	H ₃ agonist H ₃ protean agonist	NET1027 NEX436		
				AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆				X	ES-392-AF						
				ValiScreen transfected cell line	CHO-K1		X	X	X		ES-392-C						
				CAMPZen frozen ready to screen cells	CHO-K1		X	X	X		ES-392-CF						
					Membrane preparation	CHO-K1		X	X			ES-392-M400UA					
	H ₄	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X		ES-393-A	Histamine dihydrochloride, [ring, methylenes- ³ H(N)]-	endogenous	NET732		
AequoZen frozen ready to screen cells				CHO-K1	Gα ₁₆				X	ES-393-AF							
ValiScreen transfected cell line				CHO-K1		X	X	X		ES-393-C							
CAMPZen frozen ready to screen cells				CHO-K1		X	X	X		ES-393-CF							
				Membrane preparation	CHO-K1		X	X			ES-393-M400UA						
KISS1 (Metastin)	KISS1 (GPR54)	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆			X	X	ES-630-A	Metastin 45-54, [¹²⁵ I]-	agonist	NEX395			
				ValiScreen transfected cell line	MDA-MB-435S		X		X		ES-630-C						
				Membrane preparation	MDA-MB-435S		X				ES-630-M400UA						
					CHO-K1		X				RBHMTSM400UA						
Leukotriene	BLT ₁ (LTB4R1)	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X		X	X	ES-340-A	Leukotriene B ₄ , [5,6,8,9,11,12,14,15- ³ H(N)]-	endogenous	NET852			
				ValiScreen transfected cell line	CHO-K1		X	X	X		ES-340-C						
				CAMPZen frozen ready to screen cells	CHO-K1		X	X			ES-340-CF						
				Membrane preparation	CHO-K1		X	X			ES-340-M400UA						
					CysLT ₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X	X	X	ES-470-A	Leukotriene D ₄ , [14,15,19,20- ³ H(N)]-	endogenous endogenous	NET1019 NET1018
	ValiScreen transfected cell line	CHO-K1		X				X		ES-470-C							
	Membrane preparation	CHO-K1		X						ES-470-M400UA							
					OXE (HM74-like)	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆		X		ES-640-A			
	ValiScreen transfected cell line	CHO-K1						X		ES-640-AF							
	Membrane preparation	CHO-K1								ES-640-AF							
	Lysophospholipid	S ₁ P ₂ (EDG5)	G _i	Human	AequoScreen aequorin double-transfected cell line	CHO-K1				X		ES-594-A					
					ValiScreen transfected cell line	CHO-K1					X		ES-594-AF				
Membrane preparation					CHO-K1						ES-594-AF						
					CHO-K1						ES-594-AF						
	S ₁ P ₄ (EDG6)	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₃₁			X	X	ES-592-A						
				ValiScreen transfected cell line	CHO-K1	Gα ₃₁				X	ES-592-AF						
				Membrane preparation	CHO-K1	Gα ₃₁				X	ES-592-AF						
					CHO-K1	Gα ₃₁				X	ES-592-AF						
	S ₁ P ₁ (EDG8)	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₃₁			X	X	ES-593-A						
				ValiScreen transfected cell line	CHO-K1	Gα ₃₁				X	ES-593-AF						
				Membrane preparation	CHO-K1	Gα ₃₁				X	ES-593-AF						
					CHO-K1	Gα ₃₁				X	ES-593-AF						
Mas related	MrgB3	G _i /G _o	Rat	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₂₅			X		ES-745-A						
				ValiScreen transfected cell line	CHO-K1	Gα ₂₅				X	ES-745-AF						
				Membrane preparation	CHO-K1	Gα ₂₅				X	ES-745-AF						
					CHO-K1	Gα ₂₅				X	ES-745-AF						
Melanin-Concentrating Hormone	MCH ₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X		X	X	ES-370-A	S36057 (Tyr), [¹²⁵ I]- Melanin-Concentrating Hormone, [¹²⁵ I]- Melanin-Concentrating Hormone (Phe ³ , Tyr ¹⁹), [¹²⁵ I]- MCH-1R Antagonist, [¹²⁵ I]-	MCH receptor agonist MCH receptor agonist MCH receptor agonist MCH ₁ antagonist	NEX396 NEX373 NEX375 NEX406			
				ValiScreen transfected cell line	CHO-K1	Gα ₁₆	X	X	X		ES-370-AF						
				CAMPZen frozen ready to screen cells	CHO-K1	Gα ₁₆	X	X	X		ES-370-C						
				Membrane preparation	CHO-K1	Gα ₁₆	X	X	X		ES-370-CF						
					Membrane preparation	CHO-K1		X	X			ES-370-M400UA					
	MCH ₂	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X	ES-371-A	S36057 (Tyr), [¹²⁵ I]- Melanin-Concentrating Hormone, [¹²⁵ I]- Melanin-Concentrating Hormone (Phe ³ , Tyr ¹⁹), [¹²⁵ I]-	MCH receptor agonist MCH receptor agonist MCH receptor agonist	NEX396 NEX373 NEX375		
				ValiScreen transfected cell line	CHO-K1	Gα ₁₆	X			X	ES-371-AF						
				CAMPZen frozen ready to screen cells	1321N1		X			X	ES-372-C						
				Membrane preparation	1321N1		X				ES-372-M400UA						
	Melanocortin	MC ₁	G _s	Human	ValiScreen transfected cell line	CHO-K1		X	X			ES-195-C	SHU9119, [¹²⁵ I]- α-MSH (Nle ⁴ , D-Phe ⁷), [¹²⁵ I]-	MC ₁ , MC ₅ agonist MC agonist	NEX362 NEX352		
					CAMPZen frozen ready to screen cells	CHO-K1		X	X			ES-195-CF					
					Membrane preparation	CHO-K1		X	X			ES-195-M400UA					
					CHO-K1		X	X			ES-195-M400UA						
MC ₂		G _s	Human	ValiScreen transfected cell line	CHO-K1							ACTH (1-39) (Tyr ²³), [¹²⁵ I]- ACTH (1-39) (Tyr ²³), [¹²⁵ I]-	MC ₂ agonist MC ₂ agonist	NEX165 NEX342			
				CAMPZen frozen ready to screen cells	CHO-K1												
				Membrane preparation	CHO-K1												
					CHO-K1												
MC ₃		G _s	Human	ValiScreen transfected cell line	CHO-K1		X	X				ES-193-C	α-MSH (Nle ⁴ , D-Phe ⁷), [¹²⁵ I]- SHU9119, [¹²⁵ I]-	MC agonist MC ₃ , MC ₄ antagonist	NEX352 NEX362		
				CAMPZen frozen ready to screen cells	CHO-K1		X	X				ES-193-CF					
				Membrane preparation	HEK293		X	X				RBXMC3M400UA					
					CHO-K1		X	X				ES-193-M400UA					
MC ₄	G _s	Human	ValiScreen transfected cell line	CHO-K1		X	X				ES-190-C	α-MSH (Nle ⁴ , D-Phe ⁷), [¹²⁵ I]- SHU9119, [¹²⁵ I]- AGRP (Ac 87-132), [¹²⁵ I]- MT II (His), [¹²⁵ I]-	MC agonist MC ₃ , MC ₄ antagonist MC ₂ agonist MC ₂ agonist	NEX352 NEX362 NEX372 NEX374			
			CAMPZen frozen ready to screen cells	CHO-K1		X	X				ES-190-CF						
			Membrane preparation	CHO-K1		X	X				ES-190-M400UA						
				CHO-K1		X	X				ES-191-A						

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	qTPP	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Fluorescence				
Melanocortin (continued)	MC ₅	G _s	Human	ValiScreen transfected cell line	CHO-K1		X	X					ES-194-C	α-MSH (Nle ⁶ , D-Phe ⁷), [¹²⁵ I]-SHU9119, [¹²⁵ I]-	MC ₁ , MC ₅ agonist	NEX352 NEX362	
				cAMPZn frozen ready to screen cells	CHO-K1			X				ES-194-CF					
				Membrane preparation	HEK293 EBNA		X					RBXMC5M400UA					
	Mouse	ValiScreen transfected cell line	CHO-K1		X						ES-194-M400UA						
		cAMPZn frozen ready to screen cells	CHO-K1		X		X				ES-192-C						
		Membrane preparation	CHO-K1		X		X				ES-192-CF						
Melatonin	MT ₁	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆	X			X			ES-620-A	Melatonin, 2-[¹²⁵ I]-	agonist	NEX236	
				AequoZn frozen ready to screen cells	CHO-K1	Gd ₁₆			X			ES-620-AF					
				ValiScreen transfected cell line	CHO-K1		X	X				ES-620-C					
	MT ₂	G _i /G _o	Human	Membrane preparation	CHO-K1			X	X				ES-620-M400UA				
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-621-A				
				AequoZn frozen ready to screen cells	CHO-K1		X			X			ES-621-AF				
Motilin	motilin	G _i /G ₁₁	Human	ValiScreen transfected cell line	CHO-K1		X						ES-621-C	Motilin, [¹²⁵ I]-	endogenous	NEX378	
				AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆	X			X	X		ES-380-A				
				ValiScreen transfected cell line	CHO-K1		X			X	X		ES-380-C				
	Membrane preparation	HEK293		X							RBHMOTM400UA						
		CHO-K1		X							ES-380-M400UA						
Muscarinic (see Acetylcholine) Neurokinin (see Tachykinin) Neuromedin U	NMU1	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆	X			X	X		ES-450-A	Neuromedin U-25 (Tyr ¹⁸), [¹²⁵ I]-, ([¹²⁵ I]-NMU-25)	NMU1, NMU2 agonist	NEX383 NEX392	
				ValiScreen transfected cell line	CHO-K1		X		X	X		ES-450-C					
				Membrane preparation	HEK293		X						RBHNU1M400UA				
	NMU2	G _i /G ₁₁	Human	ValiScreen transfected cell line	CHO-K1			X			X		ES-451-A	Neuromedin U-8, [¹²⁵ I]-	NMU1, NMU2 agonist	NEX392 NEX383	
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-451-C				
				Membrane preparation	HEK293		X						RBHNU2M400UA				
Membrane preparation	CHO-K1		X							ES-451-M400UA							
Neuropeptide S	NPS	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₁₆	X			X	X		ES-770-A	Neuropeptide S (Tyr ¹⁵), [¹²⁵ I]-	endogenous	NEX411	
				ValiScreen transfected cell line	CHO-K1		X			X	X		ES-770-C				
Neuropeptide FF	NPFF1	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X		X				ES-491-C	Neuropeptide FF (D-Tyr ¹ , N-MePhe ³), [¹²⁵ I]-	NPFF1, NPFF2 agonist	NEX381	
				cAMPZn frozen ready to screen cells	CHO-K1				X			ES-491-CF					
				Membrane preparation	CHO-DHFR(-)		X						RBHNF1M400UA				
	NPFF2	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1	Gd ₃₁		X			X	X		ES-490-A	Neuropeptide FF (D-Tyr ¹ , N-MePhe ³), [¹²⁵ I]-	NPFF1, NPFF2 agonist	NEX381
				AequoScreen aequorin double-transfected cell line	CHO-K1		X		X	X			ES-490-C				
				Membrane preparation	CHO-K1		X			X			ES-490-CF				
Membrane preparation	HEK293		X							RBHNF2M400UA							
										RBHGPM400UA							
Neuropeptide W/B	NPBW2 (GPR8)	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X						ES-351-A	Neuropeptide W23, [¹²⁵ I]- Neuropeptide B23 (Des-Br), [¹²⁵ I]-	NPBW2, endogenous NPBW2, endogenous	NEX403 NEX405	
				ValiScreen transfected cell line	CHO-K1		X		X				ES-351-C				
	Y ₁	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1			X						ES-351-A	Peptide YY (porcine), [¹²⁵ I]-, ([¹²⁵ I]-PYY) Peptide YY (Leu ³ , Pro ³), [¹²⁵ I]- Pancreatic polypeptide, [¹²⁵ I]- Peptide YY (human), [¹²⁵ I]-, ([¹²⁵ I]-PYY) Neuropeptide Y (porcine), [¹²⁵ I]- Peptide YY (porcine), [¹²⁵ I]-, ([¹²⁵ I]-PYY) Pancreatic polypeptide, [¹²⁵ I]- Peptide YY (human), [¹²⁵ I]-, ([¹²⁵ I]-PYY) Neuropeptide Y (porcine), [¹²⁵ I]-	NPY agonist Y ₁ agonist endogenous endogenous agonist	NEX240 NEX334 NEX315 NEX341 NEX222
				Non-recombinant, membrane preparation	SK-N-MC		X						RBHNP1M400UA				
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-352-A				
				AequoZn frozen ready to screen cells	CHO-K1		X			X			ES-352-AF				
Y ₂	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1			X		X				ES-352-C				
			cAMPZn frozen ready to screen cells	CHO-K1			X				ES-352-CF						
			Non-recombinant, membrane preparation	KAN-TS		X						RBHNP2M400UA					
			Membrane preparation	HEK293		X						ES-690-A					
Neurotensin	NTS ₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-690-A	Neurotensin (Tyr ⁸), [¹²⁵ I]- Neurotensin, [3,11-tyrosyl-3,5- ³ H(N)]-	endogenous endogenous	NEX198 NET605	
				AequoZn frozen ready to screen cells	CHO-K1		X			X		ES-690-AF					
				ValiScreen transfected cell line	CHO-K1		X			X		ES-690-C					
	Membrane preparation	CHO-K1		X							6110518400UA						
		HEK293		X							RBXNT1M400UA						
		CHO-K1		X							ES-690-M400UA						
NTS ₂	G _i /G ₁₁	Human	ValiScreen transfected cell line	1321N1			X					ES-691-C	Neurotensin (Tyr ⁸), [¹²⁵ I]- Neurotensin, [3,11-tyrosyl-3,5- ³ H(N)]-	endogenous endogenous	NEX198 NET605		
			Membrane preparation	HEK293		X						6110566400UA					
			ValiScreen transfected cell line	1321N1		X						ES-691-M400UA					
Nicotinic	GPR109A (HM74A)	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gd ₃₁	X			X	X		ES-760-A				
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-760-C				
				cAMPZn frozen ready to screen cells	CHO-K1				X				ES-760-CF				
	Membrane preparation	CHO-K1		X	X						ES-760-M400UA						
		IMR32		X							RBHNCM400UA						
Obestatin	delta	G _i /G _o	Human	Membrane preparation	CHO-K1		X						RBHODM400UA	Obestatin, [¹²⁵ I]- Naltrindole, [5,7- ³ H]- Met (O ² -E)kephalin, [¹²⁵ I]-Tyr ¹ - E)kephalin (2-D-Alanine-5-D-Leucine), [Tyrosyl-3,5- ³ H(N)]- E)kephalin (Tyrosyl-2,6- ³ H(N)), (2-D-Penicillamine, 5-D-Penicillamine) (DPDPE) Deltorphin II (2-D-Ala), [Tyrosyl-3,5- ³ H]- Diprenorphine [15,16- ³ H]-	endogenous delta antagonist	NEX415 NET1065	
				Membrane preparation	HEK293		X						6110549400UA				
Opioid	delta	G _i /G _o	Mouse	Membrane preparation	Sf9		X						6110115400UA	delta agonist delta agonist delta agonist delta agonist delta agonist antagonist	NEX149 NET648 NET922 NET1087 NET1121		

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	GTPγS	cAMP	Inositol Phosphate	Calcium	Calcineurin	Fluorescence				
Opioid (continued)	kappa	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-541-A	Diprenorphine, [15,16- ³ H]-U-69,593, [phenyl-3,4- ³ H]-	antagonist kappa agonist	NET1121 NET952	
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-541-C				
				Membrane preparation	HEK293		X						6110558400UA				
					CHO-K1		X	X					ES-541-M400UA				
					CHO-K1	Gd ₁₆	X				X	X	ES-542-A				
					CHO-K1	Gd ₁₆	X	X	X				ES-542-AF				
	mu	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1			X			X	X		Diprenorphine, [15,16- ³ H]-Naloxone, [N-allyl-2,3- ³ H]-DAMGO, [tyrosyl-3,5- ³ H(N)]-	antagonist mu antagonist mu agonist	NET1121 NET719 NET902	
				AequoZen frozen ready to screen cells	CHO-K1		X				X		ES-542-C				
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-542-CF				
				cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-542-CF				
				Membrane preparation	CHO-K1		X	X					ES-542-M400UA				
					CHO-K1		X				X	X	ES-542-A				
NOP (ORL1)	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1			X			X	X		Nociceptin (Tyr ¹⁴), [¹²⁵ I]-Nociceptin, [leucyl-3,4,5- ³ H]-	NOP agonist NOP agonist	NEX338 NET1130		
			ValiScreen transfected cell line	CHO-K1		X	X	X				ES-230-C					
			cAMPZen frozen ready to screen cells	CHO-K1		X		X				ES-230-CF					
			Membrane preparation	CHO-K1		X		X				6110540400UA					
				HEK293		X						RBHRLM400UA					
				CHO-K1		X				X	X	ES-330-A					
Orexin	OX ₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		Orexin-A, [¹²⁵ I]-	OX ₁ , OX ₂ agonist	NEX367		
				AequoZen frozen ready to screen cells	CHO-K1		X				X					ES-330-AF	
				ValiScreen transfected cell line	CHO-K1		X				X					ES-330-C	
	OX ₂	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		Gd ₁₆	X			X	X		Orexin-A, [¹²⁵ I]-	OX ₁ , OX ₂ agonist	NEX367	
				AequoZen frozen ready to screen cells	CHO-K1		Gd ₁₆	X			X		ES-331-AF				
				ValiScreen transfected cell line	CHO-K1		X				X		ES-331-C				
Oxytocin												Ornithine Vasotocin Analog, [¹²⁵ I]-Oxytocin (Tyr ⁷), [¹²⁵ I]-Oxytocin, [Tyr ² , 6- ³ H]-	antagonist endogenous endogenous	NEX254 NEX187 NET858			
Parathyroid Hormone												Parathyroid Hormone (1-34) amide (Nle ⁶² , Tyr ⁶⁴) (Rat), [¹²⁵ I]-Parathyroid Hormone (1-34) (Nle ⁶ , Tyr ⁶⁴) (Human), [¹²⁵ I]-	PTH ₁ , PTH ₂ agonist endogenous	NEX353 NEX397			
Peptide P518	QRFP (OX ₂ -like, GPR103)	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X	ES-670-A	QRFP-43, [¹²⁵ I]-	QRFP, endogenous	NEX408		
				ValiScreen transfected cell line	CHO-K1		X				X	ES-670-C					
				Membrane preparation	CHO-K1		X					ES-670-M400UA					
					HEK293		X					RBH103M400UA					
Platelet Activating Factor													Hexadecyl-2-acetyl-sn-glycyl-3-phosphocholine, 1-O-acetyl- ³ H(N)]-	endogenous	NET910		
Prokineticin	PKR ₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		Gd ₁₆	X			X	X	ES-750-A	MIT-1, [¹²⁵ I]-, Bolton-Hunter labeled	PKR ₁ , PKR ₂ antagonist	NEX410	
				AequoZen frozen ready to screen cells	CHO-K1		Gd ₁₆	X			X		ES-750-AF				
				Membrane preparation	CHO-K1		X						ES-750-M400UA				
	PKR ₂	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		MIT-1, [¹²⁵ I]-, Bolton-Hunter labeled	PKR ₁ , PKR ₂ antagonist	NEX410		
				AequoZen frozen ready to screen cells	CHO-K1		X			X		ES-751-A					
				Membrane preparation	CHO-K1		X				X	ES-751-AF					
					CHO-K1		X					ES-751-M400UA					
					HEK293		X					RBHPK2M400UA					
					CHO-K1		X			X		ES-751-C					
				ValiScreen transfected cell line	CHO-K1		X			X		ES-751-CF					
Prolactin Releasing Peptide	PRRP	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		Prolactin-Releasing Peptide (Tyr ²⁰), [¹²⁵ I]-	PRRP agonist	NEX384		
				ValiScreen transfected cell line	CHO-K1		X			X		ES-480-A					
				Membrane preparation	CHO-K1		X			X		ES-480-M400UA					
Prostanoid	DP ₁	G _s	Human	AequoScreen aequorin double-transfected cell line	1321N1		X			X	X		Prostaglandin D ₂ , [5,6,8,9,12,14,15- ³ H(N)]-	endogenous	NET616		
				ValiScreen transfected cell line	1321N1		X			X		ES-560-A					
				cAMPZen frozen ready to screen cells	1321N1		X			X		ES-560-C					
				Membrane preparation	1321N1		X			X		ES-560-CF					
					1321N1		X					ES-560-M400UA					
	DP ₂ (ORTH2)	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		Gd ₁₆	X			X	X	ES-561-A	Prostaglandin D ₂ , [5,6,8,9,12,14,15- ³ H(N)]- Prostaglandin E ₂ , [5,6,8,11,12,14,15- ³ H(N)]-	endogenous endogenous	NET616 NET428	
				ValiScreen transfected cell line	CHO-K1		X	X	X			ES-561-C					
				cAMPZen frozen ready to screen cells	CHO-K1		X			X		ES-561-CF					
				Membrane preparation	HEK293		X					RBHT2M400UA					
					CHO-K1		X	X				ES-561-M400UA					
	EP ₂	G _s	Human	ValiScreen transfected cell line	HEK293		X		X				ES-562-C	Prostaglandin E ₂ , [5,6,8,11,12,14,15- ³ H(N)]-	endogenous	NET428	
				cAMPZen frozen ready to screen cells	HEK293		X		X			ES-562-CF					
				Membrane preparation	HEK293		X					ES-562-M400UA					
	EP ₄	G _s	Human	AequoScreen aequorin double-transfected cell line	HEK293		Gd ₁₆	X			X	X	ES-563-A	Prostaglandin E ₂ , [5,6,8,11,12,14,15- ³ H(N)]-	endogenous	NET428	
				ValiScreen transfected cell line	HEK293		X		X			ES-563-C					
				Membrane preparation	HEK293		X					ES-563-M400UA					
	FP	G _i /G ₁₁	Human	ValiScreen transfected cell line	1321N1		X				X		ES-564-C	Prostaglandin F _{2α} , [5,6,8,9,11,12,14,15- ³ H(N)]-	endogenous	NET433	
				Membrane preparation	1321N1		X					ES-564-M400UA					
	TP	G _i /G ₁₁	Human	ValiScreen transfected cell line	HEK293 EBNA		X						RBHTPC	SQ 29548, [5,6- ³ H]-	TP receptor antagonist	NET936	
				Membrane preparation	HEK293 EBNA		X					RBHTPM400UA					
Protease-activated	PAR2	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-780-A				
	PAR4	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X			ES-782-A				
				AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-782-AF				
Purinergic P2Y	P2Y ₂	G _i /G ₁₁	Human	ValiScreen transfected cell line	1321N1		X			X			ES-290-C				
	P2Y ₆	G _i /G ₁₁	Human	ValiScreen transfected cell line	1321N1		X			X			ES-292-C				
	P2Y ₁₁	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	1321N1		X			X			ES-293-A				
				AequoZen frozen ready to screen cells	1321N1		X			X			ES-293-AF				
	P2Y ₁₂	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X		X				RBHP12C				

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.		
							Binding	GTPγS	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Fluorescence						
Relaxin	RXFP3	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1					X			ES-656-A	INSL-3, [¹²⁵ I]-	RXFP2, RXFP3 agonist	NEX419			
				ValiScreen transfected cell line	CHO-K1		X	X				ES-656-C							
				cAMPZen frozen ready to screen cells	CHO-K1			X				ES-656-CF							
	RXFP4	G _i /G _o	Human	Membrane preparation	CHO-K1		X						ES-656-M400UA	INSL-3, [¹²⁵ I]-	RXFP2, RXFP3 agonist	NEX419			
				ValiScreen transfected cell line	CHO-K1		X	X				ES-659-C							
				Membrane preparation	CHO-K1		X					ES-659-M400UA							
	RXFP1 & RXFP2				AequoScreen aequorin double-transfected cell line	CHO-K1					X		ES-659-A	H2 Relaxin, [¹²⁵ I]-, Bolton-Hunter Labeled H3 Relaxin, [¹²⁵ I]-	RXFP1 (LGR7) agonist RXFP1 (LGR7) agonist	NEX409 NEX418			
	Serotonin (see 5-hydroxytryptamine)																		
	Somatostatin																		
Sst	sst ₁	G _i /G _o	Human	ValiScreen transfected cell line	CHO-K1		X	X					ES-520-C	Somatostatin (Tyr ¹¹), [¹²⁵ I]- Somatostatin-14 (Tyr ¹¹) (human, recombinant), [¹²⁵ I]-	sst ₁ -sst ₆ agonist sst ₁ -sst ₆ agonist	NEX389 NEX446			
				Membrane preparation	CHO-K1		X	X					ES-520-M400UA						
				AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X		ES-521-A						
	sst _{2a}	G _i /G _o	Human	AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆						X		ES-521-AF	Somatostatin (Tyr ¹¹), [¹²⁵ I]- Somatostatin-14 (Tyr ¹¹) (human, recombinant), [¹²⁵ I]-	sst ₁ -sst ₆ agonist	NEX389		
				ValiScreen transfected cell line	CHO-K1	Gα ₁₆	X	X	X				ES-521-C						
				cAMPZen frozen ready to screen cells	CHO-K1		X	X	X				ES-521-CF						
	sst ₂	G _i /G _o	Human	Membrane preparation	CHO-K1		X	X						ES-521-M400UA	Somatostatin (Tyr ¹¹), [¹²⁵ I]- Somatostatin-14 (Tyr ¹¹) (human, recombinant), [¹²⁵ I]-	sst ₁ -sst ₆ agonist	NEX389 NEX446		
				AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X		ES-523-A						
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-523-C						
	sst ₃	G _i /G _o	Human	cAMPZen frozen ready to screen cells	CHO-K1					X				ES-523-CF	Somatostatin (Tyr ¹¹), [¹²⁵ I]- Somatostatin-14 (Tyr ¹¹) (human, recombinant), [¹²⁵ I]-	sst ₁ -sst ₆ agonist	NEX389 NEX446		
				Membrane preparation	CHO-K1		X	X					ES-523-M400UA						
				AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X			X	X		ES-524-A						
	sst ₄	G _i /G _o	Human	AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆		X			X			ES-524-AF	Somatostatin (Tyr ¹¹), [¹²⁵ I]- Somatostatin-14 (Tyr ¹¹) (human, recombinant), [¹²⁵ I]-	sst ₁ -sst ₆ agonist	NEX389 NEX446		
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-524-C						
				cAMPZen frozen ready to screen cells	CHO-K1				X				ES-524-CF						
				Membrane preparation	HEK293 EBNA		X						RBHST4M400UA						
					CHO-K1		X	X					ES-524-M400UA						
					CHO-K1		X	X					ES-524-M400UA						
	sst ₅	G _i /G _o	Human	AequoScreen aequorin double-transfected cell line	CHO-K1	Gα ₁₆	X				X	X		ES-522-A	Somatostatin (Tyr ¹¹), [¹²⁵ I]- Somatostatin-14 (Tyr ¹¹) (human, recombinant), [¹²⁵ I]-	sst ₁ -sst ₆ agonist	NEX389 NEX446		
				AequoZen frozen ready to screen cells	CHO-K1	Gα ₁₆					X		ES-522-AF						
				ValiScreen transfected cell line	CHO-K1		X	X	X				ES-522-C						
				cAMPZen frozen ready to screen cells	CHO-K1				X				ES-522-CF						
				Membrane preparation	HEK293 EBNA		X						RBHST5M400UA						
					CHO-K1		X	X					ES-522-M400UA						
Succinate (see GPR91)																			
Tachykinin																			
NK	NK ₁	G _i /G ₁₁	Human	Non-recombinant, membrane preparation	UC11		X						6110551400UA	Substance P (Lys ³), [¹²⁵ I]-, Bolton-Hunter labeled Substance P (Tyr ⁹), [¹²⁵ I]- Substance-P (Sar ² , 11-Met(O ₂)), [2-prolyl-3,4- ³ H]- Hemokinin-1, [¹²⁵ I]-, Bolton-Hunter labeled Substance P, [leucyl-3,4,5- ³ H(N)]- Substance-P (9-Sar, 11-Met(O ₂)), [2-Prolyl-3, 4-3H]-	NK ₁ agonist NK ₁ agonist NK ₁ agonist NK ₁ agonist NK ₁ agonist NK ₁ agonist	NEX190 NEX152 NET1025 NEX414 NET1115 NET1025			
				NK ₂	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X				X		ES-251-A
							ValiScreen transfected cell line	CHO-K1		X							X		ES-251-C
	Membrane preparation	CHO-K1					X					ES-251-M400UA							
	NK ₃	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-252-A	SR 48968, [benzamide-4- ³ H]- Eledoisin (Lys ⁵), [¹²⁵ I]-, Bolton-Hunter labeled Neurokinin A, [¹²⁵ I]-, [¹²⁵ I]-Substance K, [¹²⁵ I]-, NKA)	NK ₂ antagonist NK ₃ , NK ₃ agonist NK ₃ agonist	NET1077 NEX218 NEX252			
				ValiScreen transfected cell line	CHO-K1		X			X		ES-252-C							
				Membrane preparation	CHO-K1		X			X		ES-252-M400UA							
	NK ₃	G _i /G ₁₁	Human	ValiScreen transfected cell line	CHO-K1		X			X			ES-252-C	Neurokinin B (MePhe ⁷), [¹²⁵ I]-, ([¹²⁵ I]-NKB) Senkide, [phenylalanyl-3,4,5- ³ H]- SR142801, [³ H]-	NK ₃ , NK ₃ agonist NK ₃ agonist NK ₃ agonist	NEX218 NEX285 NET997 NET1183			
				Membrane preparation	CHO-K1		X					ES-252-M400UA							
					CHO-K1		X					ES-252-M400UA							
	Thyrotropin Releasing Hormone																		
	TRH	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X			ES-700-A	Thyrotropin Releasing Hormone (3-Methyl-histidine ²), [L-histidyl-4- ³ H(N)]- Thyrotropin Releasing Hormone (His ⁵), [¹²⁵ I]-	TRH agonist TRH agonist	NET705 NEX153		
AequoZen frozen ready to screen cells				CHO-K1		X				X		ES-700-AF							
ValiScreen transfected cell line				CHO-K1		X				X		ES-700-C							
Membrane preparation				CHO-K1		X						ES-700-M400UA							
Urotensin																			
UT (Uro II, GPR14)	G _i /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X			ES-440-A	Urotensin II (human), [¹²⁵ I]-Tyr ² -	endogenous	NEX379			
			ValiScreen transfected cell line	CHO-K1		X				X		ES-440-C							
			PhotoScreen Phdina double-transfected cell line	CHO-K1		X			X			AX-007-PCF							
			Membrane preparation	CHO-K1		X						ES-440-M400UA							
			Rat	AequoScreen aequorin double-transfected cell line	CHO-K1	X												ES-441-A	
																			AequoZen frozen ready to screen cells
Vasoactive Intestinal Peptide	PAC ₁ (PACAP)	G _s	Human	Membrane preparation	CHO-K1		X						RBMUR2M400UA	PACAP 27, [¹²⁵ I]-	PAC ₁ , endogenous	NEX294			
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-272-A						
				ValiScreen transfected cell line	CHO-K1		X			X	X		ES-273-A						
	VPAC ₁ (VIP)	G _s	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X			ES-273-A	Vasoactive Intestinal Polypeptide, [¹²⁵ I]-	VPAC ₁ , endogenous	NEX192		
				Non-recombinant, membrane preparation	HT29		X						RBHVIPM400UA						
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-274-A						
Vasopressin	VPAC ₂	G _s	Human	Non-recombinant, membrane preparation	HT29		X						ES-274-A	Vasoactive Intestinal Polypeptide, [¹²⁵ I]-	V _{1a} antagonist	NEX310			
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-361-A						
				AequoZen frozen ready to screen cells	CHO-K1		X			X			ES-361-AF						
	V _{1a}	G _i /G ₁₁	Human	ValiScreen transfected cell line	1321N1		X				X		ES-361-C	Vasopressin (linear), V _{1a} antagonist (Phenylacetyl ¹ , 0-Me-D-Tyr ² , [¹²⁵ I]-Arg ⁸)- Vasopressin, 8-arginine, [phenylalanyl-3,4,5- ³ H(N)]-	endogenous	NET800			
				Membrane preparation	HEK293		X						RBHV1AM400UA						
					1321N1		X						ES-361-M400UA						

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	qTPG	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Fluorescence				
Vasopressin (continued)	V _{1b}	G _q /G ₁₁	Human	AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X		ES-362-A	Vasopressin, 8-arginine, [phenylalanyl-3,4,5- ³ H(N)]-	endogenous	NET800	
				AequoZen frozen ready to screen cells	CHO-K1				X		ES-362-AF						
				ValiScreen transfected cell line	1321N1		X			X		ES-362-C					
				Membrane preparation	CHO-K1		X					6110543400UA					
					HEK293		X					RBHV1BM400UA					
	V ₂	G _s	Human	ValiScreen transfected cell line	CHO-K1		X				X		ES-362-M400UA				
				Membrane preparation	CHO-K1		X					X	ES-360-C				
				AequoScreen aequorin double-transfected cell line	CHO-K1		X			X	X	ES-363-A					
				ValiScreen transfected cell line	1321N1		X		X			ES-363-C					
				cAMPZen frozen ready to screen cells	1321N1		X		X			ES-363-CF					
Membrane preparation	CHO-K1		X					6110541400UA									
					1321N1		X				ES-363-M400UA						
Cytokine & Growth Hormone Receptors																	
Growth hormone Please inquire for receptor availability																	
Interleukin Please inquire for receptor availability																	
Interleukin-1β, [¹²⁵ I]-																	
Interleukin-2, [¹²⁵ I]-																	
Interleukin-6, [¹²⁵ I]-																	
Leptin, [¹²⁵ I]-																	
Leptin (murine), [¹²⁵ I]-																	
Prolactin, [¹²⁵ I]-																	
Natriuretic Peptide Receptors																	
Atrial Natriuretic Factor Please inquire for receptor availability																	
ANF (rat), [¹²⁵ I]-																	
ANF (human), [¹²⁵ I]-																	
Tyrosine Kinase Receptors																	
Epidermal Growth Factor																	
EGF Human Non-recombinant, membrane preparation A431 X																	
EGFR1 Human AequoScreen aequorin double-transfected cell line CHO-K1 upon request X																	
Fibroblast Growth Factor																	
FGFR1 Human AequoScreen aequorin double-transfected cell line CHO-K1 upon request X																	
Granulocyte Macrophage-Colony Stimulating-Factor																	
Please inquire for receptor availability																	
Insulin-like Growth Factor																	
Please inquire for receptor availability																	
Insulin-like Growth Factor I, [¹²⁵ I]-																	
Insulin-like Growth Factor II, [¹²⁵ I]-																	
Nerve-Growth factor																	
Please inquire for receptor availability																	
Nerve Growth Factor (human), [¹²⁵ I]-																	
Nerve Growth Factor (human), recombinant, [¹²⁵ I]-																	
Nerve Growth Factor (rat)																	
Platelet-Derived Growth Factor																	
Please inquire for receptor availability																	
PDGF-BB, [¹²⁵ I]-																	
Transforming Growth Factor																	
Please inquire for receptor availability																	
Transforming Growth Factor β1, [¹²⁵ I]-																	
Tumor Necrosis Factor																	
Please inquire for receptor availability																	
Tumor Necrosis Factor-α, [¹²⁵ I]-																	
Vascular Endothelial Growth factor																	
Please inquire for receptor availability																	
Vascular Endothelial Growth factor, [¹²⁵ I]-																	
Insulin																	
Please inquire for receptor availability																	
Insulin (porcine), [¹²⁵ I]-																	
Insulin (human), [¹²⁵ I]-																	
Insulin (Tyr ^{A14}) (porcine), [¹²⁵ I]-																	
Transporters																	
Acetylcholine																	
Dopamine n.a. Human Membrane preparation CHO-K1 X RBHDATM400UA Hemicholinium-3 diacetate salt, [methyl- ³ H]-																	
WIN 35428, [N-methyl- ³ H]-																	
RTI-55, [¹²⁵ I]-																	
RTI-121, [¹²⁵ I]-																	
Cocaine, levo-[benzoyl-3,4- ³ H(N)]-																	
Mazindol, [4- ³ H]-																	
GBR 12935, [propylene-2,3- ³ H]-																	
Methyl-4-phenylpyridinium acetate, [N-methyl- ³ H]-, (MPP(+))																	
Norepinephrine n.a. Human Membrane preparation MDCK X RBHNETM400UA Nisoxetine hydrochloride, [N-methyl- ³ H]-																	
Desmethylimipramine hydrochloride, [benzene ring, 10,11- ³ H]-																	
Serotonin n.a. Human Membrane preparation HEK293 X RBHSTM400UA Imipramine hydrochloride, [benzene ring- ³ H(N)]-																	
Paroxetine, [phenyl-6- ³ H]-																	
Citalopram, [N-methyl-3H]-																	
Transferrin																	
Transferrin, [¹²⁵ I]-																	
Ion Channels																	
5-Hydroxytryptamine																	
5-HT _{2A} AequoScreen aequorin double-transfected cell line HEK293 X ES-402-A																	
AequoZen frozen ready to screen cells HEK293 X ES-402-AF																	
5-HT ₃ Human Membrane preparation HEK293 X RBHSM400UA GR 65630, [N-methyl- ³ H]-																	
BRL-43694, [9-methyl- ³ H]-																	
5-HT ₃ blocker																	
5-HT ₃ blocker																	

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	qTPG	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Calcium Fluorescence				
Aspartate	Please inquire for receptor/channel availability													Aspartic acid, D-[2,3- ³ H]-	endogenous	NET581	
Benzodiazepine (GABA-A)	Please inquire for receptor/channel availability													Diazepam, [methyl- ³ H]- Flunitrazepam, [methyl- ³ H]- Muscimol, 3-hydroxy-5-aminomethylisoxazole, [methylene- ³ H(N)]- SR 95531, [butyryl-2,3- ³ H]- Pregnan-3α-ol-20-one, 5α-[9,11,12- ³ H(N)]- RO 5-4864, [N-methyl- ³ H]- PK 11195, [N-methyl- ³ H]- RO 15-1788, [N-methyl- ³ H]- RO 15-4513, [7,9- ³ H]-	GABA-A agonist GABA-A agonist GABA-A agonist GABA-A agonist GABA-A modulator peripheral antagonist peripheral agonist central antagonist photoaffinity antagonist	NET564 NET567 NET574 NET946 NET1047 NET704 NET885 NET757 NET925	
Calcium Channel	Please inquire for receptor/channel availability													Calcium-45 Radionuclide ω-Conotoxin MVIIC, [125I]- ω-Conotoxin GVIA (Tyr ²⁵), [125I]- Diltiazem, cis-(+)-[N-methyl- ³ H]- Nitrendipine, [5-methyl- ³ H]- PN200-110, (+)-[5-methyl- ³ H]- Verapamil hydrochloride, [N-methyl- ³ H]- Ryanodine, [9,21- ³ H(N)]-	Non-L-type antagonist N-type antagonist L-type antagonist L-type antagonist L-type antagonist L-type antagonist Ryanodine agonist	NEZ013 NEX323 NEX239 NET847 NET741 NET863 NET810 NET950	
Calcium Release Activated	CRAC (Stim1 + Ora1)			PhotoScreen Photina double-transfected cell line validated in e-phys	HEK293					X			AX-013-PCL				
Glutamate	Please inquire for receptor/channel availability													GCP 39653, [propyl-2,3- ³ H]- MK-801, (+)-[3- ³ H]- Ifenprodil, [phenyl- ³ H]- Amino-3-hydroxy-5-methylisoxazole-4-propionic acid, DL-α-[5-methyl- ³ H]- Kainic acid, [vinylidene- ³ H]- TCP, N-(1-(2-Thienyl)cyclohexyl)-3,4-piperidine, [piperidyl-3,4- ³ H(N)]- NAAAG, [glutamate-3,4- ³ H]- Glutamic Acid, L-[3,4- ³ H]-	NMDA agonist non-competitive NMDA agonist NMDA; polyamine site AMPA agonist Kainate agonist Non-competitive glutamate antagonist endogenous endogenous	NET1050 NET972 NET1089 NET833 NET875 NET886 NET1082 NET490	
Glycine	Please inquire for receptor/channel availability													Serine, D-[³ H(G)]- Strychnine, [benzene ring- ³ H]-	endogenous antagonist	NET1092 NET490	
HERG	K ⁺ channel		Human	Membrane preparation	HEK293		X							RBHERGM400UA BeKm-1 (Tyr ¹), [125I]- Astemizole, [O-methyl- ³ H]- Dotelidide, [N-methyl- ³ H]-	ERG blocker ERG blocker ERG blocker	NEX412 NET1140 NET1144	
Potassium Channel	Please inquire for receptor/channel availability													Apamin (His ¹⁸ NH ₂), [125I]- Glibenclamide Gliburide, [cyclohexyl-2,3- ³ H(N)]- α-Dendrotoxin, [125I]- Rubidium-86 Radionuclide	antagonist ATP-dependent antagonist blocker	NET242 NET1024 NEX434 NEZ072	
Nicotinic	n.a.		Human	Non-recombinant, membrane preparation	IMR32		X							RBHNICM400UA Epibatidine, [125I]-, [(³ H)-IPH]- (+/-)-Epibatidine, [5,6-bicycloheptyl- ³ H]- Choline chloride, [methyl- ³ H]- A-85380, [125I]- α-Bungarotoxin (Tyr ²⁶), [125I]- Cytisine hydrochloride, [3,5-3H(N)]- Nicotine, L-(-)-[N-ethyl- ³ H]-	CNS, skeletal muscle and autonomic ganglion activation CNS, skeletal muscle and autonomic ganglion activation Neuronal CNS alpha4 beta2 Neuromuscular blocker CNS activator CNS activator	NEX358 NET1102 NET109 NEX413 NEX126 NET1054 NET827	
Purinergic	P2RX _{2/3} / p2rx ₃		Human/rat	PhotoScreen Photina double-transfected cell line validated in e-phys	CHO-K1						X		AX-014-PCL				
	P2RX ₁		Human	PhotoScreen Photina double-transfected cell line validated in e-phys	CHO-K1							X	AX-009-PCF				
	P2RX ₄		Human	PhotoScreen Photina double-transfected cell line validated in e-phys	HEK293							X	AX-015-PCF				
Sodium Channel	Please inquire for receptor/channel availability													Batrachotoxinin A 20-α-Benzoate, [benzoyl-2,5- ³ H]- Sodium-22 Radionuclide	activator	NET876 NEZ081	
Transient receptor potential	TRPA1		Human	PhotoScreen Photina double-transfected cell line validated in e-phys	HEK293							X	AX-004-PCL				
	TRPC3		Human	VaiScreen transfected cell line validated in e-phys and membrane potential assay	HEK293								AX-011-C				
	TRPC6		Human	VaiScreen transfected cell line validated in e-phys and membrane potential assay	HEK293								AX-012-C				
Vanilloid														Resiniferatoxin, [125I]- Resiniferatoxin, [³ H]-	agonist agonist	NEX394 NEX1132	
Voltage gated potassium channel	K _v 1.3 (KCNA3)		Human	VaiScreen transfected cell line validated in e-phys and membrane potential assay	CHO-DUKX								AX-010-C				
Nuclear Receptors																	
Aldosterone	Please inquire for receptor availability													Aldosterone, D-[1,2,6,7- ³ H(N)]-	MR, endogenous	NET419	
Androgen	Please inquire for receptor availability													Dihydrotestosterone (5α androstan-17β-ol-3-one), [1,2,4,5,6,7- ³ H(N)]- Testosterone, [1, 2, 6, 7, 16, 17- ³ H(N)]- Testosterone, [1,2,6,7- ³ H(N)]- Methyltrienolone (R1881) Methyltrienolone, [17α-methyl- ³ H]- Mibolerone Mibolerone, [17α-methyl- ³ H]-	AR, endogenous AR, endogenous AR, endogenous AR agonist AR agonist AR agonist AR agonist	NET453 NET553 NET370 NLP005 NET590 NLP024 NET919	

Receptor type	Subtype	G-protein coupling	Species	Product Type	Host cell	Promiscuous G-Protein Added	Validated for							Cat. No.	Complementary radioligands	Pharmacological action	Cat. No.
							Binding	GTPγS	cAMP	Inositol Phosphate	Calcium	Calcium Luminescence	Calcium Fluorescence				
Ecdysteroid	Please inquire for receptor availability														Ecdysone, α-[23,24- ³ H(N)]- Ponasterone A, [24,25,26,27- ³ H(N)]-	Ecdysteroid agonist Ecdysteroid agonist	NET621 NET1070
Estrogen	Please inquire for receptor availability														Estradiol, [2,4,6,7,16,17- ³ H(N)]- Estradiol, [2,4,6,7- ³ H(N)]- Estradiol 17β-D-glucuronide, [Estradiol-6,7- ³ H(N)]- Estradiol, [16α-[¹²⁵ I]iodo-3,17β]- Estradiol, [16α-[¹²⁵ I]iodo-3,17β]- Estradiol, [¹²⁵ I]- Dexamethasone, [6,7- ³ H(N)]- Cortisol, [¹²⁵ I]- Corticosterone, [1,2,6,7- ³ H(N)]- Hydrocortisone [1,2,6,7- ³ H(N)]-, (Cortisol, [1,2,6,7- ³ H(N)]-)	ER, endogenous ER, endogenous ER, endogenous ER, endogenous ER, endogenous ER, endogenous GR agonist GR, endogenous GR, endogenous GR, endogenous PR, endogenous photoaffinity agonist photoaffinity agonist	NET517 NET317 NET1106 NEX144 NEX144L NEX438 NET467 NEX435 NET399 NET396 NET381 NET555 NLP004
Glucocorticoid	Please inquire for receptor availability														Progesterone, [1,2,6,7- ³ H(N)]- Promegestone (R5020*), [17α-methyl- ³ H]- Promegestone (R5020*) Solid	GR agonist GR, endogenous GR, endogenous GR, endogenous	NET467 NEX435 NET399 NET396
Progestin	Please inquire for receptor availability														Thyroxine, L-[¹²⁵ I]-, ([¹²⁵ I]-T4) Thyroxine, L-[¹²⁵ I]-, ([¹²⁵ I]-T4) Thyroxine, L-[¹²⁵ I]-, ([¹²⁵ I]-T4) Triiodothyronine, ([¹²⁵ I]-Reverse T3), L-3,3',5'-[¹²⁵ I]- Triiodothyronine, ([¹²⁵ I]-Reverse T3), L-3,5',3'-[¹²⁵ I]- Triiodothyronine, ([¹²⁵ I]-Reverse T3), L-3,5',3'-[¹²⁵ I]- Triiodothyronine, ([¹²⁵ I]-Reverse T3), L-3,5',3'-[¹²⁵ I]-	THR, endogenous THR, endogenous THR, endogenous THR, endogenous THR, endogenous THR, endogenous	NEX1111 NEX111X NEX1109 NEX110 NEX110H NEX110X
Thyroid hormones	Please inquire for receptor availability														Dihydroxyvitamin D ₃ , 1α, 25-[26,27- ³ H]- Hydroxyvitamin D ₃ , 25-[26,27- ³ H]-	VDR, endogenous endogenous	NET626 NET349
Vitamin D₃	Please inquire for receptor availability																
Wild-type membranes																	
1321N1 cells	Wild Type Control Membranes		Human	Membrane preparation	1321N1										RBH1321M010MG		
A9 cells	Wild Type Control Membranes		Mouse	Membrane preparation	A9										RBMA9M010MG		
CHO-K1 cells	Wild Type Control Membranes		Hamster	Membrane preparation	CHO-K1										RBCCOHOM010MG		
HEK293 cells	Wild Type Control Membranes		Human	Membrane preparation	HEK293										RBHHEKM010MG		
K562 cells	Wild Type Control Membranes		Human	Membrane preparation	K562										RBHK56M010MG		
THP-1 cells	Wild Type Control Membranes		Human	Membrane preparation	THP-1										RBHTHPM010MG		
U373 cells	Wild Type Control Membranes		Human	Membrane preparation	U373										RBHU37M010MG		
Signal transduction																	
Inositol triphosphate															Inositol-1,3,4,5-tetrakisphosphate, [inositol-1- ³ H(N)]- Inositol-1,4,5-trisphosphate, D-[inositol-1- ³ H(N)]-	endogenous endogenous	NET941 NET911
Nitric oxide															Arginine monohydrochloride L-[2,3,4- ³ H]-	NOS substrate	NET1123
Phospholipase															Phosphatidylinositol 4,5-bisphosphate (PIP ₂), [inositol-2- ³ H(N)]- Phosphatidylinositol 4,5-bisphosphate, [inositol-1- ³ H]-	PLC, endogenous PLC, endogenous	NET895 NET1135
Sphingomyelinase															Sphingomyelin (egg), [choline methyl- ³ H]-	Sphingomyelinase substrate	NET1134
Fatty acid															Myristic acid (tetradecanoic acid), [9,10- ³ H(N)]-	Endogenous substrate	NET830
GTP analogue															GTPγS	Non-hydrolyzable GTP analogue	NEG030
Parental cell lines																	
				AequoScreen aequorin parental cell line (medium level)	CHO-K1	Gα _{1β}									X	ES-000-A2	
				AequoZen parental Frozen Cells (medium level)	CHO-K1	Gα _{1β}									X	ES-000-A2F	
				AequoScreen aequorin parental cell line	CHO-K1	Gα _{1β}									X	ES-000-A12	
				AequoZen parental Frozen Cells	CHO-K1	Gα _{1β}									X	ES-000-A12F	
				AequoScreen aequorin parental cell line	CHO-K1	Gα _{1β}									X	ES-000-A21	
				AequoZen parental Frozen Cells	CHO-K1	Gα _{1β}									X	ES-000-A21F	
				AequoScreen aequorin parental cell line	CHO-K1	Gα _{1β}									X	ES-000-A24	
				AequoZen parental Frozen Cells	CHO-K1	Gα _{1β}									X	ES-000-A24F	
				AequoScreen aequorin parental cell line	1321N1	Gα _{1β}									X	ES-000-A27	
				AequoZen parental Frozen Cells	1321N1	Gα _{1β}									X	ES-000-A27F	
				AequoScreen aequorin parental cell line	1321N1	Gα _{1β}									X	ES-000-A28	
				AequoZen parental Frozen Cells	1321N1	Gα _{1β}									X	ES-000-A28F	
				AequoScreen aequorin parental cell line	HEK293	Gα _{1β}									X	ES-000-A30	
				AequoZen parental Frozen Cells	HEK293	Gα _{1β}									X	ES-000-A30F	
				AequoScreen aequorin parental cell line	HEK293	Gα _{1β}									X	ES-000-A26	
				AequoZen parental Frozen Cells	HEK293	Gα _{1β}									X	ES-000-A26F	
Plasmids																	
				AequoScreen aequorin plasmid		Gα _{1β}									X	ES-002-AC	
				AequoScreen aequorin plasmid		Gα _{1β}									X	ES-003-AC	
Starter Kits																	
Muscarinic and Histamine	M ₃ and H ₁			AequoZen frozen cells Starter Kit	CHO-K1									X	ES-001-AF		
Muscarinic and Histamine	M ₃ and H ₁			AequoScreen Starter Kit	CHO-K1									X	ES-001-A		
Adenosin	A ₃			PhotoScreen Starter Kit	CHO-K2									X	AX-001-PCF		

©2008 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer Inc. AequoScreen, AequoZen, cAMPZen and NEN are registered trademarks of PerkinElmer Inc. PhotoScreen and ValiScreen are trademarks of PerkinElmer Inc. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at anytime without notice and disclaims liability for editorial, pictorial or typographical errors.