

# LANCE and LANCE *Ultra* TR-FRET Technology Product List

## ORDERING INFORMATION

Please refer to the following for ordering LANCE® TR-FRET and LANCE *Ultra* TR-FRET products. LANCE TR-FRET and LANCE *Ultra* TR-FRET are for research purposes only and are not intended for human diagnostic applications.

### Table of Contents

LANCE <i>Ultra</i> Biomarker Detection Kits .....	2
LANCE and LANCE <i>Ultra</i> Binding Kits.....	4
cAMP, LANCE and Secondary Messenger Assay Kits .....	4
LANCE <i>Ultra</i> Cellular Kinase Assay Kits .....	4
LANCE <i>Ultra</i> KinaSelect Assays .....	4
General LANCE Toolbox Donor Fluorophores .....	5
General LANCE Toolbox Acceptor Fluorophores .....	6
LANCE <i>Ultra</i> Kinase Assay <i>ULight</i> -substrates and Matching Europium Anti-phospho Antibodies .....	6
LANCE <i>Ultra</i> Epigenetic Toolbox Reagents .....	8
LANCE Labeling Reagents .....	8
Buffers and Solutions .....	8

## LANCE *Ultra* Biomarker Detection Kits

Kit name	Target	500 Assay Points	10,000 Assay Points
<i>Animal Health</i>			
Immunoglobulin G (bovine)	IgG (bovine)	TRF1530C	TRF1530M
<i>Angiogenesis</i>			
Angiopoietin 1	ANGPT1	TRF3014C	TRF3014M
Angiopoietin-like 3	ANGPTL3	TRF3016C	TRF3016M
Angiopoietin-like 4	ANGPTL4	TRF3017C	TRF3017M
<i>Biologics/Bioprocess</i>			
CHO Host Cell Proteins	CHO-P	TRF1301C	TRF1301M
CHO Host Cell Proteins (broad reactivity)	CHO-P (broad reactivity)	TRF1210C	TRF1210M
E.coli Host Cell Proteins	E.coli HCP	TRF1261C	TRF1261M
Immunoglobulin E (human)	IgE (human)	TRF1292C	TRF1292M
Immunoglobulin G (rabbit)	IgG (rabbit)	TRF1104C	TRF1104M
Immunoglobulin G (human)	IgG (human)	TRF1205C	TRF1205M
Immunoglobulin G (mouse)	IgG (mouse)	TRF1529C	TRF1529M
Immunoglobulin M (human)	IgM (human)	TRF1263C	TRF1263M
Kappa light chain (human)	κ chain	TRF3023C	TRF3023M
Lambda light chain (human)	λ chain	TRF3024C	TRF3024M
PER.C6 Host Cell Proteins	PER.C6 HCP	TRF1302C	TRF1302M
Residual DNA contaminants	Residual DNA contaminants	TRF1331C	TRF1331M
Residual Protein A	Residual Protein A	TRF1287C	TRF1287M
<i>Cancer</i>			
Alpha-1 type I collagen	COL1A1	TRF1371C	TRF1371M
Bcl-2	Bcl-2	TRF1384C	TRF1384M
COX-2	COX-2	TRF1376C	TRF1376M
E-cadherin	E-cadherin	TRF1370C	TRF1370M
Epidermal Growth Factor	EGF	TRF1339C	TRF1339M
Epidermal Growth Factor Receptor	EGFR	TRF1340C	TRF1340M
Fibronectin	Fibronectin	TRF1351C	TRF1351M
Matrix Metalloproteinase-3	MMP-3	TRF1284C	TRF1284M
Matrix Metalloproteinase-7	MMP-7	TRF1378C	TRF1378M
Matrix Metalloproteinase-8	MMP-8	TRF1374C	TRF1374M
Matrix Metalloproteinase-12	MMP-12	TRF1375C	TRF1375M
Mesencephalic Astrocyte Derived Neurotrophic Factor	MANF	TRF3009C	TRF3009M
Nidogen-1	Nidogen-1	TRF1367C	TRF1367M
Oncostatin M	Oncostatin M	TRF1341C	TRF1341M
Perlecan	Perlecan	TRF1377C	TRF1377M
Phosphatase and tensin homolog	PTEN	TRF1380C	TRF1380M
PIP Collagen	PIP Collagen	TRF1353C	TRF1353M
Programmed cell death protein 1	PD-1	TRF1343C	TRF1343M
Programmed death ligand 1	PD-L1	TRF1355C	TRF1355M
Vascular endothelial growth factor	VEGF	TRF1201C	TRF1201M
Vascular endothelial growth factor (mouse)	VEGF (mouse)	TRF1520C	TRF1520M
Receptor Tyrosine-Protein Kinase ErbB-2 (HER2)	ErbB2 (HER2)	TRF1235C	TRF1235M
Tumor protein p53	p53	TRF1388C	TRF1388M
<i>Cardiovascular</i>			
Fibrinogen	Fibrinogen	TRF1352C	TRF1352M
Tissue plasminogen activator	tPA	TRF1250C	TRF1250M
Transferrin	Transferrin	TRF1311C	TRF1311M
<i>Central Nervous System</i>			
Apolipoprotein E	Apo-E	TRF1395C	TRF1395M
Amyloid β 1-40	Aβ(1-40)	TRF1345C	TRF1345M
Brain Derived Neurotrophic Factor	BDNF	TRF1317C	TRF1317M
Frataxin	Frataxin	TRF1322C	TRF1322M
Glutamic acid decarboxylase 67	GAD67	TRF1397C	TRF1397M
Microtubule-associated protein tau	Tau	TRF1271C	TRF1271M
TAR DNA-binding protein 43	TDP-43	TRF1387C	TRF1387M

## LANCE Ultra Biomarker Detection Kits (Continued)

Kit name	Target	500 Assay Points	10,000 Assay Points
<i>Inflammation and Cytokine</i>			
C-C motif chemokine 2 (Monocyte Chemoattractant Protein-1)	CCL2/MCP-1	TRF1244C	TRF1244M
C-C motif chemokine 2 (Monocyte Chemoattractant Protein-1) (mouse)	CCL2/MCP-1 (mouse)	TRF1509C	TRF1509M
C-C motif chemokine 3 (Macrophage inflammatory protein 1- $\alpha$ )	CCL3/MIP1 $\alpha$	TRF1257C	TRF1257M
C-C motif chemokine 4 (Macrophage inflammatory protein 1- $\beta$ )	CCL4/MIP-1 $\beta$	TRF1258C	TRF1258M
C-C motif chemokine 5 (T-cell-specific protein RANTES)	CCL5/RANTES	TRF1289C	TRF1289M
C-Reactive Protein	CRP	TRF1233C	TRF1233M
C-X-C Motif Chemokine 1/Growth-regulated Oncogene $\alpha$	CXCL1/GRO- $\alpha$	TRF1279C	TRF1279M
C-X-C motif chemokine 9 (Gamma-interferon-induced monokine)	CXCL9/MIG	TRF1280C	TRF1280M
C-X-C motif chemokine 10 (10 kDa IFN $\gamma$ -induced protein)	CXCL10/IP-10	TRF1259C	TRF1259M
Interferon $\alpha$	IFN- $\alpha$	TRF1297C	TRF1297M
Interferon $\lambda$ 2	IFN- $\lambda$ 2	TRF3039C	TRF3039M
Interferon $\gamma$	IFN- $\gamma$	TRF1217C	TRF1217M
Interferon $\gamma$ (mouse)	IFN- $\gamma$ (mouse)	TRF1501C	TRF1501M
Interleukin 1 receptor antagonist	IL-1RA	TRF1358C	TRF1358M
Interleukin 1 $\alpha$	IL-1 $\alpha$	TRF1238C	TRF1238M
Interleukin 1 $\beta$	IL-1 $\beta$	TRF1220C	TRF1220M
Interleukin 1 $\beta$ (mouse)	IL-1 $\beta$ (mouse)	TRF1503C	TRF1503M
Interleukin 2	IL-2	TRF1221C	TRF1221M
Interleukin 4	IL-4	TRF1266C	TRF1266M
Interleukin 5	IL-5	TRF1267C	TRF1267M
Interleukin 6	IL-6	TRF1223C	TRF1223M
Interleukin 6 (mouse)	IL-6 (mouse)	TRF1504C	TRF1504M
Interleukin 8	IL-8	TRF1224C	TRF1224M
Interleukin 11	IL-11	TRF1293C	TRF1293M
Interleukin 12/p40	IL-12/p40	TRF1382C	TRF1382M
Interleukin 17A	IL-17A	TRF1219C	TRF1219M
Interleukin 17A (mouse)	IL-17A (mouse)	TRF1514C	TRF1514M
Interleukin 17F	IL-17F	TRF1391C	TRF1391M
Interleukin 23A	IL-23A	TRF3006C	TRF3006M
Interleukin 28B	IL-28B	TRF1365C	TRF1365M
Interleukin 31	IL-31	TRF3003C	TRF3003M
Tumor Necrosis Factor $\alpha$	TNF- $\alpha$	TRF1208C	TRF1208M
Tumor Necrosis Factor $\alpha$ (mouse)	TNF- $\alpha$ (mouse)	TRF1505C	TRF1505M
<i>Metabolic</i>			
Adiponectin	Adiponectin	TRF1209C	TRF1209M
Albumin	Albumin	TRF1294C	TRF1294M
Apolipoprotein A1	Apo-A1	TRF1389C	TRF1389M
Apolipoprotein C3	Apo-CIII	TRF1396C	TRF1396M
C-peptide (mouse/rat)	C-peptide (mouse/rat)	TRF1510C	TRF1510M
Chorionic Gonadotropin	hCG	TRF3001C	TRF3001M
Cortisol	Cortisol	TRF1314C	TRF1314M
Follicle Stimulating Hormone	FSH	TRF1394C	TRF1394M
Glucagon-Like Peptide 1 (7-36)	GLP-1 (7-36)	TRF1215C	TRF1215M
Glucagon	Glucagon	TRF1312C	TRF1312M
Insulin	Insulin	TRF1204C	TRF1204M
Insulin-like Growth Factor 1	IGF-1	TRF1236C	TRF1236M
Insulin-like Growth Factor 1 (rat)	IGF-1 (rat)	TRF1552C	TRF1552M
Leptin	Leptin	TRF1225C	TRF1225M
Leptin (mouse)	Leptin (mouse)	TRF1521C	TRF1521M
Muscleblind-like protein 1	MBNL1	TRF3008C	TRF3008M
Plasminogen Activator Inhibitor-1	PAI-1	TRF1286C	TRF1286M
Proprotein Convertase Subtilisin/Kexin 9	PCSK9	TRF1270C	TRF1270M
Thyroid Stimulating Hormone	TSH	TRF3005C	TRF3005M
<i>Autophagy</i>			
LC3B	LC3B	TRF1306C	TRF1306M

## LANCE and LANCE Ultra Binding Kits

Kit name	500 Assay Points	10,000 Assay Points
<i>LANCE Ultra Binding Kits</i>		
EGF-EGFR LANCE Ultra Binding Kit	TRF1366C	TRF1366M
<i>LANCE Fc-receptor Binding Kits</i>		
FCGR3A/CD16a (176Phe/F158) LANCE Ultra Binding Kit	TRF1347C	TRF1347M
FCGR3A/CD16a (176Val/V158) LANCE Ultra Binding Kit	TRF1348C	TRF1348M

## cAMP, LANCE Secondary Messenger Assay Kits

Kit name	500 assays	1,000 assays	10,000 assays	50,000 assays
LANCE Ultra cAMP Kit		TRF0262	TRF0263	TRF0264
LANCE cAMP Kit	AD0262	AD0262E	AD0263	AD0264

## LANCE Ultra Cellular Kinase Assay Kits

Kit name	500 assays	10,000 assays
Phospho-AKT1 (Ser473) Cellular Detection Kit	TRF4013C	TRF4013M
Phospho-AKT1 (Thr308) Cellular Detection Kit	TRF4006C	TRF4006M
Phospho-AKT1/2/3 (Ser473) Cellular Detection Kit	TRF4002C	TRF4002M
Phospho-AKT1/2/3 (Thr308) Cellular Detection Kit	TRF4003C	TRF4003M
Total AKT1 Cellular Detection Kit	TRF4007C	TRF4007M
Phospho-AMPK $\alpha$ (Thr172) Cellular Detection Kit	TRF4027C	TRF4027M
Phospho-BTK (Tyr223) Cellular Detection Kit	TRF4021C	TRF4021M
Total BTK Cellular Detection Kit	TRF4022C	TRF4022M
Phospho-eIF2A (Ser51) Cellular Detection Kit	TRF4008C	TRF4008M
Total eIF2 $\alpha$ Cellular Detection Kit	TRF4023C	TRF4023M
Phospho-EGFR (Tyr1068) Cellular Detection Kit	TRF4015C	TRF4015M
Total EGFR Cellular Detection Kit	TRF4016C	TRF4016M
Phospho-ERK1/2 (Thr202/Tyr204) Cellular Detection Kit	TRF4000C	TRF4000M
Total ERK1/2 Cellular Detection Kit	TRF4001C	TRF4001M
Phospho-LCK (Tyr394) Cellular Detection Kit	TRF4026C	TRF4026M
Phospho-NF $\kappa$ B (Ser536) Cellular Detection Kit	TRF4019C	TRF4019M
Total NF $\kappa$ B Cellular Detection Kit	TRF4020C	TRF4020M
Phospho-SLP76 (Ser376) Cellular Detection Kit	TRF4010C	TRF4010M
Total SLP76 Cellular Detection Kit	TRF4017C	TRF4017M
Phospho-SMAD1 (Ser463/465) Cellular Detection Kit	TRF4011C	TRF4011M
Phospho-SMAD2 (Ser465/467) Cellular Detection Kit	TRF4012C	TRF4012M
Phospho-SMAD3 (Ser423/425) Cellular Detection Kit	TRF4014C	TRF4014M
Phospho-Src (Tyr419) Cellular Detection Kit	TRF4018C	TRF4018M
Phospho-STAT1 (Tyr701) Cellular Detection Kit	TRF4028C	TRF4028M
Total STAT1 Cellular Detection Kit	TRF4029C	TRF4029M
Phospho-STAT3 (Tyr705) Cellular Detection Kit	TRF4004C	TRF4004M
Total STAT3 Cellular Detection Kit	TRF4005C	TRF4005M
Phospho-STAT5 (Tyr694/699) Cellular Detection Kit	TRF4009C	TRF4009M
Phospho-TBK1 (Ser172)	TRF4031C	TRF4031M
Phospho-ZAP-70 (Tyr319) Cellular Detection Kit	TRF4024C	TRF4024M
Total ZAP-70 Cellular Detection Kit	TRF4025C	TRF4025M
TREM2 Cellular Detection Kit	TRF4030C	TRF4030M

## LANCE Ultra KinaSelect™ Assays

Kit name	1,000 assays	5 x 100 assays
LANCE Ultra KinaSelect™ Kit for Ser/Thr Kinases		TRF0300-C
LANCE Ultra KinaSelect™ Kit for Tyr Kinases	TRF0301-D	

## LANCE TECHNOLOGY TOOLBOX PRODUCTS

ToolBox reagents, substrates and buffers are available for you to setup your own assay and controls. Our local Field Application Specialists team is available to support you in establishing your own assays, or you can ask our OnPoint Custom Assays Development team to develop the assay you need to your specifications. Alternatively, our custom labeling team is available to label your antibodies, proteins, peptides or small molecules in the technology of your choice.

### General LANCE Toolbox Donor Fluorophores

Product	10 µg	50 µg	1 mg	10 nmoles
LANCE Eu-anti-6X His antibody	AD0205	AD0110	AD0111	
LANCE Eu-anti-c-myc antibody	AD0111	AD0206	AD0114	
LANCE Eu-anti-HA antibody		AD0084	AD0085	
LANCE Eu-anti-GST antibody	AD0252	AD0253	AD0254	
LANCE Eu-anti-FLAG Antibody	AD0273	AD0274	AD0275	
LANCE Eu-anti-FITC Antibody	AD0270	AD0271	AD0272	
LANCE Eu-anti-Protein C tag antibody	AD0279	AD0280	AD0281	
LANCE Eu-anti-V5 Antibody	AD0276	AD0277	AD0278	
LANCE Eu Protein G	AD0211	AD0070	AD0071	
LANCE Eu-anti-human IgG	AD0212	AD0074	AD0075	
LANCE Eu-anti-mouse IgG		AD0076	AD0077	
LANCE Eu-anti-rabbit IgG		AD0082	AD0083	
LANCE Eu-W1024-streptavidin		AD0062	AD0063	
LANCE Eu-W8044-streptavidin		AD0060	AD0061	
LANCE Eu-biotin				CR91-100
LANCE Eu-antiphosphotyrosine PT66		AD0068	AD0069	
LANCE Eu-antiphosphotyrosine P-Tyr-100	AD0203	AD0161	AD0162	
LANCE Eu-antiphosphotyrosine PY20		AD0066	AD0067	
LANCE Eu-antiphosphothreonine	AD0094			
LANCE Eu-antiphospho-threonine-proline	AD0099			
LANCE Eu-anti-phospho-serine/threonine antibody	AD0176			
LANCE Eu-anti-phospho-serine/threonine-phenylalanine antibody	AD0178			
LANCE Eu-anti-phospho-serine/threonine -proline antibody	AD0180			
LANCE Eu-anti-phospho-PKA substrate antibody	AD0182			
LANCE Eu-anti-phospho-(Ser) PKC substrate antibody	AD0188			
LANCE Eu-anti-phospho-(Ser) 14-3-3 binding motif antibody	AD0190			
LANCE Eu-anti-phospho-(Ser) 14-3-3 binding motif 4E2 antibody	AD0192			

## General LANCE Toolbox Acceptor Fluorophores

Product	1 nmole	10 nmoles	100 nmoles
<i>ULight</i> <sup>TM</sup> -streptavidin	TRF0102-D	TRF0102-M	TRF0102-R
<i>ULight</i> anti-His antibody	TRF0105-D	TRF0105-M	TRF0105-R
<i>ULight</i> anti-FLAG antibody	TRF0059D	TRF0059M	TRF0059R
<i>ULight</i> anti-GST antibody	TRF0104-D	TRF0104-M	TRF0104-R
<i>ULight</i> Protein A	TRF0103-D	TRF0103-M	
	0.2 nmoles	2 nmoles	20 nmoles
<i>ULight</i> anti-FITC antibody	TRF503D	TRF503M	TRF503R
<i>ULight</i> anti-c-myc	TRF504D	TRF504M	TRF504R
<i>ULight</i> anti-human IgG	TRF500D	TRF500M	TRF500R
<i>ULight</i> anti-mouse IgG	TRF501D	TRF501M	TRF501R
<i>ULight</i> anti-rabbit IgG	TRF502D	TRF502M	TRF502R
<i>ULight</i> anti-V5 antibody	TRF505D	TRF505M	TRF505R
<i>ULight</i> anti-Protein C tag antibody	TRF506D	TRF506M	TRF506R
	1 mg	30 mg	50 mg
SureLight <sup>TM</sup> APC-streptavidin for kinase	CR130-100		CR130-150
SureLight <sup>TM</sup> APC-streptavidin for binding assays	AD0201	AD0202	
SureLight <sup>TM</sup> APC anti-c-myc	AD0059C		
SureLight <sup>TM</sup> APC anti-FLAG	AD0059F		
SureLight <sup>TM</sup> APC anti-GST	AD0059G		
SureLight <sup>TM</sup> APC anti-6x His	AD0059H		
SureLight <sup>TM</sup> APC anti-mouse IgG	AD0059M		
SureLight <sup>TM</sup> APC anti-rabbit IgG	AD0059R		
SureLight <sup>TM</sup> APC wheat germ agglutinin (WGA)	AD0059W		

## LANCE *Ultra* Kinase Assay *ULight*-substrates and Matching Europium Anti-phospho Antibodies

Substrate	Substrate + Matching Antibody	Size (for 20 $\mu$ L reaction)	Catalog number	Phospho-Motif
<i>ULight</i> -CREBtide	<i>ULight</i> -CREBtide (Ser133)	1,000 assay points	TRF0107-D	RRP $\underline{S}$ YRK
		10,000 assay points	TRF0107-M	
	Europium-anti-phospho-CREB (Ser133)	1,562 assay points	TRF0200-D	
		15,625 assay points	TRF0200-M	
<i>ULight</i> MBP	<i>ULight</i> -Myelin Basic Protein Peptide	1,000 assay points	TRF0109-D	VTPRIPPP
		10,000 assay points	TRF0109-M	
	Europium-anti-phospho-Myelin Basic Protein	1,562 assay points	TRF0201-D	
		15,625 assay points	TRF0201-M	
<i>ULight</i> -Crosstide	<i>ULight</i> -Crosstide (GSK-3 $\alpha$ Ser21)	1,000 assay points	TRF0106-D	PRTSS $\underline{S}$ FAE
		10,000 assay points	TRF0106-M	
	Europium-anti-phospho-Crosstide (GSK-3 $\alpha$ Ser21)	1,562 assay points	TRF0202-D	
		15,625 assay points	TRF0202-M	
<i>ULight</i> -PLK	<i>ULight</i> -PLK (Ser137) Peptide	1,000 assay points	TRF0110-D	RRR $\underline{S}$ LLE
		10,000 assay points	TRF0110-M	
	Europium-anti-phospho-PLK (Ser137)	1,562 assay points	TRF0203-D	
		15,625 assay points	TRF0203-M	
<i>ULight</i> -PKC	<i>ULight</i> -PKC Substrate (Ala25Ser)	1,000 assay points	TRF0108-D	RKG $\underline{S}$ LRQ
		10,000 assay points	TRF0108-M	
	Europium anti-phospho-PKC Peptide (Ala25Ser)	1,562 assay points	TRF0207-D	
		15,625 assay points	TRF0207-M	
<i>ULight</i> -I $\kappa$ B $\alpha$	<i>ULight</i> -I $\kappa$ B $\alpha$ (Ser32/36) Peptide	1,000 assay points	TRF0113-D	RHD $\underline{S}$ GLD $\underline{S}$ M
		10,000 assay points	TRF0113-M	
	Europium-anti-phospho-I $\kappa$ B $\alpha$ (Ser32/36)	1,562 assay points	TRF0206-D	
		15,625 assay points	TRF0206-M	

**LANCE *Ultra* Kinase Assay *ULight*-substrates and Matching Europium Anti-phospho Antibodies (Continued)**

Substrate	Substrate + Matching Antibody	Size (for 20 $\mu$ L reaction)	Catalog number	Phospho-Motif
<i>ULight</i> -4E-BP1	<i>ULight</i> -eIF4E-binding protein 1 (Thr37/46) Peptide	1,000 assay points	TRF0128-D	STIPGGTLFSTIPG
		10,000 assay points	TRF0128-M	
	Europium-anti-phospho-eIF4E-binding protein 1 (Thr37/46)	1,562 assay points	TRF0216-D	
		15,625 assay points	TRF0216-M	
<i>ULight</i> -Acetyl-CoA Carboxylase	<i>ULight</i> -Acetyl-CoA Carboxylase (Ser79) SAMS Peptide	1,000 assay points	TRF0133-D	RSAMS <sub>S</sub> GLHL
		10,000 assay points	TRF0133-M	
	Europium-anti-phospho-Acetyl-CoA Carboxylase (Ser79)	1,562 assay points	TRF0208-D	
		15,625 assay points	TRF0208-M	
<i>ULight</i> -Cdc25C	<i>ULight</i> -Cdc25C (Ser216) Peptide	1,000 assay points	TRF0123-D	YRSP <sub>S</sub> MPEN
		10,000 assay points	TRF0123-M	
	Europium-anti-phospho-(Ser) 14-3-3 binding motif, mAb 4E2	1,562 assay points	AD0192	
<i>ULight</i> -CDK1	<i>ULight</i> -CDK1 (Tyr15) Peptide	1,000 assay points	TRF0122-D	GEGTY <sub>V</sub> VVY
		10,000 assay points	TRF0122-M	
	Europium-anti-phosphotyrosine (PT66)	15,625 assay points	AD0068	
		312,500 assay points	AD0069	
<i>ULight</i> -Histone H3 (Thr3/Ser10)	<i>ULight</i> -Histone H3 (Thr3/Ser10) Peptide	1,000 assay points	TRF0125-D	ARIKQTARK <sub>S</sub> TG
		10,000 assay points	TRF0125-M	
	Europium-anti-phospho-Histone H3 (Ser10)	1,562 assay points	TRF0210-D	
		15,625 assay points	TRF0210-M	
	Europium-anti-phospho-Histone H3 (Thr3)	1,562 assay points	TRF0211-D	
		15,625 assay points	TRF0211-M	
<i>ULight</i> -GS	<i>ULight</i> -Glycogen Synthase (Ser641/pSer657) Peptide	1,000 assay points	TRF0131-D	PAP <sub>S</sub> VPPSPSL
		10,000 assay points	TRF0131-M	
	Europium-anti-phospho-Glycogen Synthase (Ser641) Antibody	1,562 assay points	TRF0220-D	
		15,625 assay points	TRF0220-M	
<i>ULight</i> -IRS-1	<i>ULight</i> -IRS-1 (Tyr983) Peptide	1,000 assay points	TRF0120-D	SRGDY <sub>M</sub> TMQ
		10,000 assay points	TRF0120-M	
	Europium-anti-phosphotyrosine (PT66)	15,625 assay points	AD0068	
		312,500 assay points	AD0069	
<i>ULight</i> -JAK-1	<i>ULight</i> -JAK-1 (Tyr1023) Peptide	1,000 assay points	TRF0121-D	DKEYY <sub>T</sub> VKD
		10,000 assay points	TRF0121-M	
	Europium-anti-phosphotyrosine (PT66)	15,625 assay points	AD0068	
		312,500 assay points	AD0069	
<i>ULight</i> -mTOR	<i>ULight</i> -mTOR (Ser2448) Peptide	1,000 assay points	TRF0119-D	TRTDS <sub>S</sub> YSAG
		10,000 assay points	TRF0119-M	
	Europium-anti phospho-mTOR (Ser2448)	1,562 assay points	TRF0209-D	
		15,625 assay points	TRF0209-M	
<i>ULight</i> -p70S6K	<i>ULight</i> -p70S6K (Thr389) Peptide	1,000 assay points	TRF0126-D	FLGFIYVAP
		10,000 assay points	TRF0126-M	
	Europium-anti-phospho-p70S6K (Thr389)	1,562 assay points	TRF0214-D	
		15,625 assay points	TRF0214-M	
<i>ULight</i> -poly GT	<i>ULight</i> -poly GT	1,000 assay points	TRF0100-D	[EY(4:1)] <sub>n</sub>
		10,000 assay points	TRF0100-M	
	Europium-anti-phosphotyrosine (PT66)	15,625 assay points	AD0068	
		312,500 assay points	AD0069	
<i>ULight</i> -poly GAT	<i>ULight</i> -poly GAT	1,000 assay points	TRF0101-D	[EAY(1:1:1)] <sub>n</sub>
		10,000 assay points	TRF0101-M	
	Europium-anti-phosphotyrosine (PT66)	15,625 assay points	AD0068	
		312,500 assay points	AD0069	

## LANCE *Ultra* Kinase Assay *ULight*-substrates and Matching Europium Anti-phospho Antibodies (Continued)

Substrate	Substrate + Matching Antibody	Size (for 20 µL reaction)	Catalog number	Phospho-Motif
<i>ULight</i> -rpS6	<i>ULight</i> -40S Ribosomal Protein S6 (pSer235/Ser236) Peptide	1,000 assay points	TRF0129-D	RRL(pS)SSLRA
		10,000 assay points	TRF0129-M	
	Europium-anti-phospho-40S Ribosomal Protein S6 (Ser235/236)	1,562 assay points	TRF0217-D	
		15,625 assay points	TRF0217-M	
<i>ULight</i> -Topo IIα	<i>ULight</i> -DNA Topoisomerase 2-α (Thr1342) Peptide	1,000 assay points	TRF0130-D	DEKIDDE
		10,000 assay points	TRF0130-M	
	Europium-anti-phospho-DNA Topoisomerase 2-α (Thr1342)	1,562 assay points	TRF0218-D	
		15,625 assay points	TRF0218-M	
<i>ULight</i> -TK	<i>ULight</i> -TK Peptide	1,000 assay points	TRF0127-D	Y
		10,000 assay points	TRF0127-M	
	Europium-anti-phosphotyrosine (PT66)	15,625 assay points	AD0068	
		312,500 assay points	AD0069	

## LANCE *Ultra* Epigenetic Toolbox Reagents

Product	10 µg	100 µg
LANCE <i>Ultra</i> Eu-anti-acetyl-Histone H3 Lysine 9 (H3K9ac)	TRF0400-D	TRF0400-M
LANCE <i>Ultra</i> Eu-anti-methyl-Histone H3 Lysine 4 (H3K4me1-2)	TRF0402-D	TRF0402-M
LANCE <i>Ultra</i> Eu-anti-methyl-Histone H3 Lysine 9 (H3K9me2)	TRF0403-D	TRF0403-M
LANCE <i>Ultra</i> Eu-anti-unmodified Histone H3 Lysine 4 (H3K4) Antibody	TRF0404-D	TRF0404-M
LANCE <i>Ultra</i> Eu-anti-acetyl-Histone H3 Lysine 27 (H3K27ac) Antibody	TRF0405-D	TRF0405-M
LANCE <i>Ultra</i> Eu-anti-di/mono-methyl-Histone H3 Lysine 27 (H3K27me2-1) Antibody	TRF0406-D	TRF0406-M
LANCE <i>Ultra</i> Eu-anti-tri-methyl-Histone H3 Lysine 27 (H3K27me3) Antibody	TRF0407-D	TRF0407-M
LANCE <i>Ultra</i> Eu-anti-di-methyl-Histone H3 Lysine 36 (H3K36me2) Antibody	TRF0408-D	TRF0408-M
LANCE <i>Ultra</i> Eu-anti-acetyl-p53 Lysine 382 (p53K382ac) Antibody	TRF0409-D	TRF0409-M
LANCE <i>Ultra</i> Eu-anti-unmodified Histone H3 Lysine 9/Lysine 27 (H3K9/K27) Antibody	TRF0411-D	TRF0411-M
LANCE <i>Ultra</i> Eu-anti-acetyl-Histone H4 Lysine (H4Kac pan) Antibody	TRF0412-D	TRF0412-M
LANCE <i>Ultra</i> Eu-anti-O-linked-GlcNAc Antibody	TRF0413-D	TRF0413-M
LANCE <i>Ultra</i> Eu-anti-methyl-Histone H4 Arginine 3 (H4R3me) Antibody	TRF0414-D	TRF0414-M
LANCE <i>Ultra</i> Eu-anti-methyl-Histone H3/H4 Arginine (H3/H4Rme pan) Antibody	TRF0415-D	TRF0415-M

## LANCE Labeling Reagents

Product	100 µg	1 mg
LANCE Eu-W1024-ITC labeling reagent	AD0096	AD0013
LANCE Eu-W1284-iodoacetamide labeling reagent	AD0107	AD0014
LANCE Eu-W8044-DTA labeling reagent		AD0020

## Buffers and Solutions

Product	Catalog number	Size
10X LANCE Detection Buffer (not for use in LANCE cAMP assays)	CR97-100	250 mL
7.5% BSA Solution	CR84-100	50 mL
LANCE Assay Buffer	TRFLAB002	2 mL
	TRFLAB100	100 mL
<i>Ultra</i> HiBlock Buffer 5X	TRF1011C	10 mL
	TRF1011F	100 mL

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

013073\_01

PKI April 2019