

Alpha Technology Product List

ORDERING INFORMATION

Please refer to the following for ordering Alpha Technology products. AlphaLISA®, AlphaScreen®, AlphaLISA® *SureFire*® and AlphaLISA® *SureFire*® *Ultra* are for research purposes or agronomy/veterinary fields diagnostics only and are not intended for human diagnostic applications.

Table of Contents

Biomarkers and Cytokines	2
AlphaLISA kits	2
AlphaPlex kits.....	8
Biologic – AlphaLISA kits.....	9
Binding Assays – AlphaLISA kits.....	10
Epigenetics Reagents	10
Epigenetics Biochemical Assay Reagents (AlphaLISA)	10
Cell-based Epigenetic Assays.....	11
Protein Phosphorylation Analysis	12
Alpha <i>SureFire</i> kits.....	12
Terbium <i>SureFire Ultra</i>	15
Beads for Biochemical kinase assays.....	17
Alpha Technology Products	18
Alpha ToolBox.....	18
Alpha Donor Beads	19
Unconjugated Beads.....	19
Antibody capture solutions	19
Tag-capture solutions.....	19
AlphaLISA Acceptor Beads	20
Unconjugated Beads.....	20
Antibody capture solutions	20
Tag-capture solutions.....	20
AlphaPlex Acceptor Beads	21
AlphaScreen Kits and Kit Components.....	21
Alpha Buffers.....	22
Control beads – TruHits	22
Control beads – Omnibeads	22

Biomarker and Cytokine AlphaLISA Kits

Agriculture and Plant Science

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Cry1F	Cry1F	AL402C	AL402F	AL402S	AL402HV
Cry2A	Cry2A	AL403C	AL403F	AL403S	AL403HV
Cry1Ab/1Ac	Cry1Ab/1Ac	AL401C	AL401F	AL401S	AL401HV

Angiogenesis

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Angiopoietin 1	ANGPT1	AL3014C	AL3014F	AL3014S	AL3014HV
Angiopoietin 2	ANGPT1	AL3015C	AL3015F	AL3015S	AL3015HV
Angiopoietin-like 3	ANGPTL3	AL3016C	AL3016F	AL3016S	AL3016HV
Angiopoietin-like 4	ANGPTL4	AL3017C	AL3017F	AL3017S	AL3017HV
Tumor Necrosis Factor α	TNF- α (human)	AL208C	AL208F	AL208S	
Tumor Necrosis Factor α	TNF- α (mouse)	AL505C	AL505F	AL505S	
Tumor Necrosis Factor α (cynomolgus)	TNF- α (cynomolgus)	AL562C	AL562F	AL562S	AL562HV
Tumor Necrosis Factor α (human) Biotin-Free	TNF- α (human)	AL325C	AL325F		AL325HV
Tumor Necrosis Factor α (mouse) Biotin-Free	TNF- α (mouse)	AL541C	AL541F		AL541HV
Transforming Growth Factor β 1 (human)	TGF- β 1 (human)	AL336C	AL336F	AL336S	AL336HV
Vascular Endothelial Growth Factor	VEGF	AL201C	AL201F	AL201S	

Animal Health

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Immunoglobulin A (bovine)	IgA (bovine)	AL536C	AL536F	AL536S	AL536HV
Immunoglobulin G (bovine)	IgG (bovine)	AL530C	AL530F	AL530S	AL530HV
Immunoglobulin G1 (bovine)	IgG1 (bovine)	AL531C	AL531F	AL531S	AL531HV
Immunoglobulin G2 (bovine)	IgG2 (bovine)	AL532C	AL532F	AL532S	AL532HV
Immunoglobulin M (bovine)	IgM (bovine)	AL533C	AL533F	AL533S	AL533HV
Interferon alpha (porcine)	IFN- α (porcine)	AL576C	AL576F	AL576S	AL576HV
Interferon beta (porcine)	IFN- β (porcine)	AL577C	AL577F	AL577S	AL577HV
Interferon gamma (bovine)	IFN- γ (bovine)	AL535C	AL535F	AL535S	AL535HV
Interferon gamma (porcine)	IFN- γ (porcine)	AL549C	AL549F	AL549S	AL549HV
Interferon gamma (ovine)	IFN- γ (ovine)	AL560C	AL560F	AL560S	AL560HV
Interleukin 1 α (bovine)	IL-1 α (bovine)	AL537C	AL537F	AL537S	AL537HV
Interleukin 1 β (bovine)	IL-1 β (bovine)	AL539C	AL539F	AL539S	AL539HV
Interleukin 1 β (porcine)	IL-1 β (porcine)	AL575C	AL575F	AL575S	AL575HV
Interleukin 2 (bovine)	IL-2 (bovine)	AL543C	AL543F	AL543S	AL543HV
Interleukin 4 (bovine)	IL-4 (bovine)	AL542C	AL542F	AL542S	AL542HV
Interleukin 4 (porcine)	IL-4 (porcine)	AL578C	AL578F	AL578S	AL578HV
Interleukin 6 (bovine)	IL-6 (bovine)	AL538C	AL538F	AL538S	AL538HV
Interleukin 6 (porcine)	IL-6 (porcine)	AL573C	AL573F	AL573S	AL573HV
Interleukin 8 (bovine/ovine)	IL-8 (bovine/ovine)	AL545C	AL545F	AL545S	AL545HV
Interleukin 8 (porcine)	IL-8 (porcine)	AL571C	AL571F	AL571S	AL571HV
Interleukin 10 (bovine/ovine)	IL-10 (bovine/ovine)	AL546C	AL546F	AL546S	AL546HV
Interleukin 10 (porcine)	IL-10 (porcine)	AL574C	AL574F	AL574S	AL574HV
Interleukin 12 (porcine)	IL-12 (porcine)	AL554C	AL554F	AL554S	AL554HV
Interleukin 12/p40 (bovine)	IL-12 p40 (bovine)	AL553C	AL553F	AL553S	AL553HV
Interleukin 13 (bovine)	IL-13 (bovine)	AL557C	AL557F	AL557S	AL557HV
Interleukin 13 (porcine)	IL-13 (porcine)	AL551C	AL551F	AL551S	AL551HV
Interleukin 17A (bovine)	IL-17A (bovine)	AL547C	AL547F	AL547S	AL547HV
Interleukin 17A (porcine)	IL-17A (porcine)	AL572C	AL572F	AL572S	AL572HV
Interleukin 18 (bovine)	IL-18 (bovine)	AL558C	AL558F	AL558S	AL558HV
Interleukin 18 (porcine)	IL-18 (porcine)	AL559C	AL559F	AL559S	AL559HV
TNF α (bovine)	TNF- α (bovine)	AL534C	AL534F	AL534S	AL534HV
TNF α (porcine)	TNF- α (porcine)	AL548C	AL548F	AL548S	AL548HV

Autophagy

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
p62	p62	AL3097C	AL3097F	AL3097S	AL3097HV
LC3B	LC3B	AL306C	AL306F	AL306S	

Cancer

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
4-1BB ligand	4-1BBL	AL3108C	AL3108F	AL3108S	AL3108HV
α -tubulin	α -tubulin	AL383C	AL383F	AL383S	AL383HV
Alpha-1 Anti-Trypsin	A1AT	AL344C	AL344F	AL344S	AL344HV
B- and T-lymphocyte attenuator	BTLA, CD272	AL3062C	AL3062F	AL3062S	AL3062HV
BCL-2	BCL-2	AL384C	AL384F	AL384S	AL384HV
Beta-2 microglobulin	Beta-2 microglobulin	AL3067C	AL3067F	AL3067S	
Beta-secretase 1	BACE-1	AL3033C	AL3033F	AL3033S	AL3033HV
Bone morphogenetic protein 7	BMP-7	AL369C	AL369F	AL369S	AL369HV
Breast cancer type 2 susceptibility protein	BRCA2	AL3036C	AL3036F	AL3036S	AL3036HV
β Nerve Growth Factor	β NGF	AL246C	AL246F	AL246S	
Cluster of differentiation 28	CD28	AL3044C	AL3044F	AL3044S	AL3044HV
Cluster of differentiation 40	CD40	AL3102C	AL3102F	AL3102S	AL3102HV
Cluster of differentiation 70	CD70	AL3103C	AL3103F	AL3103S	AL3103HV
Cluster of differentiation 80	CD80	AL3055C	AL3055F	AL3055S	AL3055HV
Cluster of differentiation 86	CD86	AL3045C	AL3045F	AL3045S	AL3045HV
Cluster of differentiation 276	CD276, B7-H3	AL3060C	AL3060F	AL3060S	AL3060HV
Collagen IV	Collagen IV	AL3034C	AL3034F	AL3034S	AL3034HV
Collagen, type I, alpha 1	COL1A1	AL371C	AL371F	AL371S	AL371HV
COX-2	COX-2	AL376C	AL376F	AL376S	AL376HV
C-X-C motif chemokine 11 (Interferon inducible T-cell chemoattractant)	CXCL11 / I-TAC	AL260C	AL260F	AL260S	
Cytotoxic T-lymphocyte-associated protein 4	CTLA-4	AL3050C	AL3050F	AL3050S	AL3050HV
E-cadherin	E-cadherin	AL370C	AL370F	AL370S	AL370HV
Epidermal growth factor	EGF	AL339C	AL339F	AL339S	AL339HV
Epidermal growth factor receptor	EGFR	AL340C	AL340F	AL340S	AL340HV
Erythropoietin	EPO	AL206C	AL206F	AL206S	
FAS ligand	FASL	AL3104C	AL3104F	AL3104S	AL3104HV
Fibroblast growth factor 23	FGF23	AL3002C	AL3002F	AL3002S	AL3002HV
Fibronectin	Fibronectin	AL3056C	AL3056F	AL3056S	AL3056HV
Galectin-9	GAL9	AL3051C	AL3051F	AL3051S	AL3051HV
Glucocorticoid-induced TNFR-related protein	GITR	AL3101C	AL3101F	AL3101S	AL3101HV
Herpesvirus entry mediator	HVEM	AL3054C	AL3054F	AL3054S	AL3054HV
Human MHC class I antigen C	HLA-C	AL3061C	AL3061F	AL3061S	AL3061HV
Human MHC class I antigen E	HLA-E	AL3057C	AL3057F	AL3057S	AL3057HV
Human Stem Cell Factor	SCF (KIT-ligand)	AL3019C	AL3019F	AL3019S	AL3019HV
Hyaluronic acid	Hyaluronic acid	AL354C	AL354F	AL354S	AL354HV
Interferon β	IFN- β	AL265C	AL265F	AL265S	
Laminin	Laminin	AL368C	AL368F	AL368S	AL368HV
Lymphocyte-activation gene 3	LAG-3	AL3058C	AL3058F	AL3058S	AL3058HV
Mucin-16 (Ovarian Carcinoma Antigen CA125)	CA125	AL256C	AL256F	AL256S	
Mast/stem cell growth factor receptor	CD117 (SCFR)	AL3021C	AL3021F	AL3021S	AL3021HV
Matrix Metalloproteinase-1	MMP-1	AL242C	AL242F	AL242S	
Matrix Metalloproteinase-2	MMP-2	AL269C	AL269F	AL269S	
Matrix Metalloproteinase-3	MMP-3	AL284C	AL284F	AL284S	
Matrix Metalloproteinase-7	MMP-7	AL378C	AL378F	AL378S	AL378HV
Matrix Metalloproteinase-8	MMP-8	AL374C	AL374F	AL374S	AL374HV
Matrix Metalloproteinase-9	MMP-9	AL243C	AL243F	AL243S	
Matrix Metalloproteinase-9 (mouse)	MMP-9 (mouse)	AL519C	AL519F	AL519S	
Matrix Metalloproteinase-12	MMP-12	AL375C	AL375F	AL375S	AL375HV
Matrix Metalloproteinase-13	MMP-13	AL318C	AL318F	AL318S	AL318HV
Matrix Metalloproteinase-23	MMP-23	AL3115C	AL3115F	AL3115S	AL3115HV
Mesothelin	Mesothelin	AL3063C	AL3063F	AL3063S	AL3063HV
N-cadherin	N-cadherin	AL379C	AL379F	AL379S	AL379HV
Neprilysin	NEP	AL337C	AL337F	AL337S	AL337HV
Nidogen-1	Nidogen-1	AL367C	AL367F	AL367S	AL367HV

Cancer (cont'd)

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Oncostatin M	Oncostatin M	AL341C	AL341F	AL341S	AL341HV
OX40 ligand	OX40L	AL3100C	AL3100F	AL3100S	AL3100HV
Perlecan	Perlecan	AL377C	AL377F	AL377S	AL377HV
Phosphatase and tensin homolog	PTEN	AL380C	AL380F	AL380S	AL380HV
PIP collagen	PIP collagen	AL353C	AL353F	AL353S	AL353HV
Programmed cell death protein 1	PD-1	AL343C	AL343F	AL343S	AL343HV
Programmed death-ligand 1	PD-L1	AL355C	AL355F	AL355S	AL355HV
Programmed death 1 ligand 2	PD-L2	AL3093C	AL3093F	AL3093S	AL3093HV
Prostatic Specific Antigen	PSA	AL228C	AL228F	AL228S	
Prostate-specific Membrane Antigen	PSMA	AL3038C	AL3038F	AL3038S	AL3038HV
Receptor Tyrosine-Protein Kinase ErbB-2 (HER2)*	ERBB2 / HER2	AL235C	AL235F	AL235S	
S100 calcium-binding protein A4	S100A4/FSP1	AL3029C	AL3029F	AL3029S	AL3029HV
T-cell immunoglobulin and mucin-domain containing-3	TIM-3	AL3052C	AL3052F	AL3052S	AL3052HV
Tissue inhibitor of metalloproteinases-1	TIMP-1	AL3090C	AL3090F	AL3090S	AL3090HV
Tissue inhibitor of metalloproteinases-2	TIMP-2	AL3091C	AL3091F	AL3091S	AL3091HV
TNF-related apoptosis-inducing ligand	TRAIL	AL3089C	AL3089F	AL3089S	AL3089HV
Transforming growth factor beta 1	TGF-β1	AL336C	AL336F	AL336S	AL336HV
Transforming growth factor beta 1 Biotin-Free	TGF-β1	AL361C	AL361F	AL361S	AL361HV
Trefoil Factor 3	TFF3	AL273C	AL273F	AL273S	
Tumor Necrosis Factor α	TNF-α (human)	AL208C	AL208F	AL208S	
Tumor Necrosis Factor α	TNF-α (mouse)	AL505C	AL505F	AL505S	
Tumor Necrosis Factor α Biotin-Free (human)	TNF-α (human)	AL325C	AL325F		AL325HV
Tumor Necrosis Factor α Biotin-Free (mouse)	TNF-α (mouse)	AL541C	AL541F		AL541HV
V-domain Ig suppressor of T cell activation	VISTA	AL3059C	AL3059F	AL3059S	AL3059HV
Vascular Cell Adhesion Molecule-1	VCAM-1	AL338C	AL338F	AL338S	AL338HV
Vimentin	Vimentin	AL372C	AL372F	AL372S	AL372HV

Cardiovascular

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Cardiac Troponin I	cTnI	AL295C	AL295F	AL295S	
Cardiac Troponin T	cTnT	AL3011C	AL3011F	AL3011S	AL3011HV
Fibrin D-dimer	D-dimer	AL290C	AL290F	AL290S	
Fibrinogen	Fibrinogen	AL352C	AL352F	AL352S	AL352HV
Intercellular Adhesion Molecule 1	ICAM-1	AL282C	AL282F	AL282S	
N-terminal pro b-type natriuretic peptide	NT-proBNP	AL3114C	AL3114F	AL3114S	AL3114S
Myeloperoxidase	MPO	AL245C	AL245F	AL245S	
Myoglobin	MB	AL285C	AL285F	AL285S	
Plasminogen Activator Inhibitor-1	PAI-1	AL286C	AL286F	AL286S	
Proprotein Convertase Subtilisin/Kexin 9	PCSK9	AL270C	AL270F	AL270S	
Tissue Plasminogen Activator	TPA	AL250C	AL250F	AL250S	
Transferrin (human)	TF (human)	AL311C	AL311F	AL311S	AL311HV
Vascular endothelial growth factor receptor 1	FLT1/VEGFR1	AL3018C	AL3018F	AL3018S	AL3018HV

Central Nervous System

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Alpha-synuclein	Alpha-synuclein	AL321C	AL321F	AL321S	AL321HV
Amyloid β 1-40	Aβ 1-40	AL202C	AL202F	AL202S	
Amyloid β 1-42	Aβ 1-42	AL203C	AL203F	AL203S	
Amyloid β 1-42 (high specificity)	Aβ 1-42 (high specificity)	AL276C	AL276F	AL276S	
Amyloid β 1-x	Amyloid β 1-x	AL288C	AL288F		
Apolipoprotein E	ApoE	AL395C	AL395F	AL395S	AL395HV
Beta-Nerve Growth Factor	β-NGF	AL246C	AL246F	AL246S	
Brain Derived Neurotrophic Factor	BDNF	AL317C	AL317F	AL317S	AL317HV
Frataxin	Frataxin	AL322C	AL322F	AL322S	AL322HV
GAD67	GAD67	AL397C	AL397F	AL397S	AL397HV

Central Nervous System (cont'd)

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Glial Derived Neurotrophic Factor	GDNF	AL319C	AL319F	AL319S	AL319HV
Mesencephalic astrocyte-derived neurotrophic factor	MANF	AL3009C	AL3009F	AL3009S	AL3009HV
Microtubule-Associated Protein Tau	Tau	AL271C	AL271F	AL271S	
Phosphorylated neurofilament heavy chain	pNFH	AL3112C	AL3112F	AL3112S	AL3112HV
TAR DNA-binding protein 43	TDP-43	AL387C	AL387F	AL387S	AL387HV
Triggering receptor expressed on myeloid cells 2	TREM2	AL3109C	AL3109F	AL3109S	AL3109HV

Food Testing

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Aflatoxin	Aflatoxin	AL390C	AL390F	AL390S	AL390HV
Beef tropinin	Beef Troponin	AL405C	AL405F	AL405S	AL405HV
Chloramphenicol	Chloramphenicol	AL385C	AL385F		AL385HV
Chloramphenicol (Biotin-Free)	Chloramphenicol	AL393C	AL393F		AL393HV
Soybean Agglutinin	Soybean Agglutinin	AL404C	AL404F	AL404S	AL404HV

Inflammation

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Bruton's tyrosine kinase	BTK	AL3110C	AL3110F	AL3110S	AL3110HV
C-C motif chemokine ligand 1	CCL1	AL3099C	AL3099F	AL3099S	AL3099HV
C-C motif chemokine ligand 2 (Monocyte Chemoattractant Protein-1)	CCL2 / MCP1	AL244C	AL244F	AL244S	
C-C motif chemokine ligand 2 (Monocyte Chemoattractant Protein 1)	CCL2 / MCP-1 (mouse/rat)	AL509C	AL509F	AL509S	
C-C motif chemokine ligand 3 (Macrophage inflammatory protein 1- α)	CCL3 / MIP-1 α	AL257C	AL257F	AL257S	
C-C motif chemokine ligand 4 (Macrophage inflammatory protein 1- β)	CCL4 / MIP-1 β	AL258C	AL258F	AL258S	
C-C motif chemokine ligand 5 (T-cell-specific protein RANTES)	CCL5 / RANTES	AL289C	AL289F	AL289S	
C-C motif chemokine ligand 5 (T-cell-specific protein RANTES)	CCL5 / RANTES (mouse)	AL516C	AL516F	AL516S	
C-C motif chemokine ligand 8	CCL8/MCP2 (mouse)	AL582C	AL582F	AL582S	AL582HV
C-C motif chemokine ligand 18	CCL18	AL3106C	AL3106F	AL3106S	AL3106HV
C-Reactive Protein	CRP	AL233C	AL233F	AL233S	
C-X-C motif chemokine 1, GRO- α	CXCL1/GRO- α	AL349C	AL349F	AL349S	
C-X-C motif chemokine 9 (Gamma-interferon-induced monokine)	CXCL9 / MIG	AL280C	AL280F	AL280S	
C-X-C motif chemokine 10 (10 kD α IFN γ -induced protein)	CXCL10 / IP-10	AL259C	AL259F	AL259S	
C-X-C motif chemokine 10 (10 kD α IFN γ -induced protein) (mouse)	CXCL10 / IP-10 (mouse)	AL556C	AL556F	AL556S	AL556HV
C-X-C motif chemokine ligand 4/PF4	CXCL4/PF4	AL398C	AL398F	AL398S	AL398HV
C-X-C motif chemokine ligand 13	CXCL13	AL3098C	AL3098F	AL3098S	AL3098HV
C-X-C motif chemokine ligand 13 (mouse)	CXCL13 (mouse)	AL581C	AL581F	AL581S	AL581HV
C-X-C motif chemokine 12, SDF-1	CXCL12/SDF-1	AL3028C	AL3028F	AL3028S	AL3028HV
Granulocyte-Colony Stimulating Factor	G-CSF	AL213C	AL213F	AL213S	
Granulocyte-Monocyte Colony-Stimulating Factor	GM-CSF (human)	AL216C	AL216F	AL216S	
Granulocyte-Macrophage Colony-Stimulating Factor	GM-CSF (mouse)	AL517C	AL517F	AL517S	
Interferon α	IFN- α	AL297C	AL297F	AL297S	
Interferon α (mouse)	IFN- α (mouse)	AL550C	AL550F	AL550S	AL550HV
Interferon λ 2	IFN- λ 2	AL3031C	AL3031F	AL3031S	AL3031HV
Interferon γ	IFN- γ	AL217C	AL217F	AL217S	
Interferon γ Biotin-Free (human)	IFN- γ	AL327C	AL327F		AL327HV

Inflammation (cont'd)

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points HV
Interferon γ (cynomolgus)	IFN- γ (cynomolgus)	AL561C	AL561F	AL561S	AL561HV
Interferon γ (mouse)	IFN- γ (mouse)	AL501C	AL501F	AL501S	
Interleukin 1 Receptor Antagonist	IL-1RA	AL358C	AL358F	AL358S	AL358HV
Interleukin 1 α	IL-1 α	AL238C	AL238F	AL238S	
Interleukin 1 β	IL-1 β	AL220C	AL220F	AL220S	
Interleukin 1 β (mouse)	IL-1 β (mouse)	AL503C	AL503F	AL503S	
Interleukin 1 β (rat)	IL-1 β (rat)	AL518C	AL518F	AL518S	
Interleukin 2	IL-2	AL221C	AL221F	AL221S	
Interleukin 2 Biotin-Free	IL-2	AL333C	AL333F		AL333HV
Interleukin 2 (mouse)	IL-2 (mouse)	AL511C	AL511F	AL511S	
Interleukin 4	IL-4	AL266C	AL266F	AL266S	
Interleukin 4 (mouse)	IL-4 (mouse)	AL567C	AL567F	AL567S	AL567HV
Interleukin 4 (rhesus)	IL-4 (rhesus)	AL564C	AL564F	AL564S	AL564HV
Interleukin 5	IL-5	AL267C	AL267F	AL267S	
Interleukin 5 (mouse)	IL-5 (mouse)	AL569C	AL569F	AL569S	AL569HV
Interleukin 6	IL-6	AL223C	AL223F	AL223S	
Interleukin 6 Biotin-Free	IL-6	AL3025C	AL3025F	AL3025S	AL3025HV
Interleukin 6 (mouse)	IL-6 (mouse)	AL504C	AL504F	AL504S	
Interleukin 6 (monkey)	IL-6 (monkey)	AL555C	AL555F	AL555S	AL555HV
Interleukin 7	IL-7	AL268C	AL268F	AL268S	
Interleukin 8	IL-8	AL224C	AL224F	AL224S	
Interleukin 8 Biotin-Free (human)	IL-8	AL328C	AL328F		AL328HV
Interleukin 10	IL-10	AL218C	AL218F	AL218S	
Interleukin 10 (mouse)	IL-10 (mouse)	AL502C	AL502F	AL502S	
Interleukin 11	IL-11	AL293C	AL293F	AL293S	
Interleukin 12 p70	IL-12 p70	AL3116C	AL3116F	AL3116S	AL3116HV
Interleukin 13	IL-13	AL240C	AL240F	AL240S	
Interleukin 13 (mouse)	IL-13 (mouse)	AL565C	AL565F	AL565S	AL565HV
Interleukin 15 Biotin-Free	IL-15	AL360C	AL360F		AL360HV
Interleukin 12/Interleukin 23	IL-12/IL-23	AL382C	AL382F	AL382S	AL382HV
Interleukin 17F	IL-17F	AL391C	AL391F	AL391S	AL391HV
Interleukin 17A	IL-17A	AL219C	AL219F	AL219S	
Interleukin 17A Biotin-Free	IL-17A	AL346C	AL346F		AL346HV
Interleukin 17A (mouse/rat)	IL-17A (mouse/rat)	AL514C	AL514F	AL514S	
Interleukin 17A (mouse)	IL-17 (mouse)	AL540C	AL540F	AL540S	AL540HV
Interleukin 18	IL-18	AL241C	AL241F	AL241S	
Interleukin 18-binding protein	IL-18BP	AL3096C	AL3096F	AL3096S	AL3096HV
Interleukin 21	IL-21	AL3043C	AL3043F	AL3043S	AL3043HV
Interleukin 22 (human/rat)	IL-22	AL320C	AL320F	AL320S	AL320HV
Interleukin 23	IL-23	AL3006C	AL3006F	AL3006S	AL3006HV
Interleukin 25	IL-25	AL3048C	AL3048F	AL3048S	AL3048HV
Interleukin 25 (mouse)	IL-25 (mouse)	AL570C	AL570F	AL570S	AL570HV
Interleukin 28B	IL-28B	AL365C	AL365F	AL365S	AL365HV
Interleukin 31	IL-31	AL3003C	AL3003F	AL3003S	AL3003HV
Interleukin 33	IL-33	AL3042C	AL3042F	AL3042S	AL3042HV
Interleukin 33 (mouse)	IL-33 (mouse)	AL566C	AL566F	AL566S	AL566HV
IP10	IP-10	AL326C	AL326F	AL326S	AL326HV
Periostin	Periostin	AL3105C	AL3105F	AL3105S	AL3105HV
Platelet-derived growth factor	PDGF (BB dimer)	AL3049C	AL3049F	AL3049S	AL3049HV
Thymic stromal lymphopoietin	TSLP	AL3040C	AL3040F	AL3040S	AL3040HV
Tumor Necrosis Factor α	TNF- α	AL208C	AL208F	AL208S	
Tumor Necrosis Factor α (cynomolgus)	TNF- α (cynomolgus)	AL562C	AL562F	AL562S	AL562HV
Tumor Necrosis Factor α (mouse)	TNF- α (mouse)	AL505C	AL505F	AL505S	
Tumor Necrosis Factor α Biotin-Free	TNF- α	AL325C	AL325F		AL325HV
Tumor Necrosis Factor α Biotin-Free (mouse)	TNF- α (mouse)	AL541C	AL541F		AL541HV
Tumor Necrosis Factor receptor 1 (soluble)	sTNFR1	AL3088C	AL3088F	AL3088S	AL3088HV
Tumor Necrosis Factor receptor 1 (soluble, mouse)	sTNFR1 (mouse)	AL579C	AL579F	AL579S	AL579HV

Kidney

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Neutrophil gelatinase-associated lipocalin	NGAL	AL364C	AL364F	AL364S	AL364HV

Metabolic and Hormones

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Adiponectin	Adiponectin	AL209C	AL209F	AL209S	
Apolipoprotein A1	Apo-A1	AL389C	AL389F	AL389S	AL389HV
Apolipoprotein B	Apo-B	AL392C	AL392F	AL392S	AL392HV
Aspartate transaminase	AST	AL3078C	AL3078F	AL3078S	AL3078HV
Serum Albumin (human)	HSA	AL294C	AL294F	AL294S	
Serum Albumin (human) Biotin-Free	HSA	AL363C	AL363F		AL363HV
Serum Albumin (mouse)	Albumin (mouse)	AL508C	AL508F	AL508S	
Bone Morphogenetic Protein 9	BMP-9	AL3013C	AL3013F	AL3013S	AL3013HV
C-Peptide (human)	C-Peptide (human)	AL299C	AL299F	AL299S	
C-Peptide (mouse/rat)	C-Peptide (mouse/rat)	AL510C	AL510F	AL510S	
Chorionic Gonadotropin (human)	hCG	AL3001C	AL3001F	AL3001S	AL3001HV
Corticosterone	Corticosterone	AL3020C	AL3020F	AL3020S	AL3020HV
Cortisol	Cortisol	AL314C	AL314F		AL314HV
Cysteine rich angiogenic inducer 61	CYR61	AL3076C	AL3076F	AL3076S	AL3076HV
Erythropoietin	EPO	AL206C	AL206F	AL206S	
Follicle Stimulating Hormone	FSH	AL394C	AL394F	AL394S	AL394HV
Ghrelin (active)	Ghrelin (active)	AL3007C	AL3007F	AL3007S	AL3007HV
Ghrelin (total)	Ghrelin (total)	AL3012C	AL3012F	AL3012S	AL3012HV
Glucagon	Glucagon	AL312C	AL312F	AL312S	AL312HV
Growth Hormone	GH	AL214C	AL214F	AL214S	
Heme oxygenase 1	HO-1	AL3026C	AL3026F	AL3026S	AL3026HV
Heme oxygenase 2	HO-2	AL3027C	AL3027F	AL3027S	AL3027HV
Human Factor VIII	Factor VIII	AL342C	AL342F	AL342S	AL342HV
Insulin	Insulin	AL204C	AL204F	AL204S	
Insulin (plasma)	Insulin (plasma)	AL350C	AL350F	AL350S	AL350HV
Insulin-like Growth Factor 1	IGF-1	AL236C	AL236F	AL236S	
Insulin-like Growth Factor 1 (rat)	IGF-1 (rat)	AL552C	AL552F	AL552S	AL552HV
Insulin-like Growth Factor 2	IGF-2	AL237C	AL237F	AL237S	
Kidney Injury Molecule 1	KIM-1	AL3075C	AL3075F	AL3075S	AL3075HV
Lactate dehydrogenase subunit A	LDHA	AL3077C	AL3077F	AL3077S	AL3077HV
Leptin	Leptin	AL225C	AL225F	AL225S	
Leptin (mouse)	Leptin (mouse)	AL521C	AL521F	AL521S	
Muscleblind-like protein 1	MBNL1	AL3008C	AL3008F	AL3008S	
Nicotinamide N-Methyltransferase	NNMT	AL3079C	AL3079F	AL3079S	AL3079HV
Paraoxonase 1	PON-1	AL3071C	AL3071F	AL3071S	AL3071HV
Paraoxonase 2	PON-2	AL3072C	AL3072F	AL3072S	AL3072HV
Placental Growth Factor	PIGF	AL3030C	AL3030F	AL3030S	AL3030HV
Progesterone	Progesterone	AL335C	AL335F	AL335S	AL335HV
Osteoprotegerin	Osteoprotegerin	AL3010C	AL3010F	AL3010S	AL3010HV
Src-homology domain-2 containing protein tyrosine phosphatase 2	SHP-2	AL3074C	AL3074F	AL3074S	AL3074HV
Testosterone	Testosterone	AL324C	AL324F	AL324S	AL324HV
Thyroid stimulating hormone	TSH	AL3005C	AL3005F	AL3005S	AL3005HV
Transthyretin	TTR	AL3111C	AL3111F	AL3111S	AL3111HV
Utrophin	Utrophin	AL3004C	AL3004F	AL3004S	AL3004HV

Virology

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Hepatitis B virus e-antigen	HBeAg	AL3082C	AL3082F	AL3082S	AL3082HV
Hepatitis B virus surface antigen	HBsAg	AL3083C	AL3083F	AL3083S	AL3083HV
Hepatitis C virus core antigen	HCV cAg	AL3092C	AL3092F	AL3092S	AL3092HV
HIV p24	HIV p24	AL207C	AL207F	AL207S	
HIV p24 (high sensitivity)	HIV p24 (high sensitivity)	AL291C	AL291F	AL291S	
HIV p24 Biotin-Free	HIV p24	AL332C	AL332F		AL332HV
Influenza A nucleoprotein	Influenza A NP	AL3084C	AL3084F	AL3084S	AL3084HV

AlphaPlex Multiplexing Biomarker Kits

Product Description	Part Number	Part Number	Part Number
	500 data points	5,000 data points	100 data points
AlphaPlex α -tubulin Tb	AP383TB-C	AP383TB-F	AP383TB-HV
AlphaPlex ICAM-1 Sm	AP282SM-C	AP282SM-F	AP282SM-HV
AlphaPlex ICAM-1 Tb	AP282TB-C	AP282TB-F	AP282TB-HV
AlphaPlex IFN- γ Sm	AP217SM-C	AP217SM-F	AP217SM-HV
AlphaPlex IFN- γ Tb	AP217TB-C	AP217TB-F	AP217TB-HV
AlphaPlex IL-1 β Sm	AP220SM-C	AP220SM-F	AP220SM-HV
AlphaPlex IL-1 β Tb	AP220TB-C	AP220TB-F	AP220TB-HV
AlphaPlex IL-2 Sm	AP221SM-C	AP221SM-F	AP221SM-HV
AlphaPlex IL-2 Tb	AP221TB-C	AP221TB-F	AP221TB-HV
AlphaPlex IL-6 Tb	AP223TB-C	AP223TB-F	AP223TB-HV
AlphaPlex IL-17 Sm	AP219SM-C	AP219SM-F	AP219SM-HV
AlphaPlex IL-17 Tb	AP219TB-C	AP219TB-F	AP219TB-HV
AlphaPlex Rat IL-22 Tb	AP320TB-C	AP320TB-F	AP320TB-HV
AlphaPlex Insulin Sm	AP204SM-C	AP204SM-F	AP204SM-HV
AlphaPlex Insulin Tb	AP204TB-C	AP204TB-F	AP204TB-HV
AlphaPlex IP-10 Sm	AP326SM-C	AP326SM-F	AP326SM-HV
AlphaPlex IP-10 Tb	AP326TB-C	AP326TB-F	AP326TB-HV
AlphaPlex TNF- α Sm	AP208SM-C	AP208SM-F	AP208SM-HV
AlphaPlex TNF- α Tb	AP208TB-C	AP208TB-F	AP208TB-HV

AlphaLISA Biologics Kits

Bioprocess

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Serum Albumin (human)	HSA	AL294C	AL294F	AL294S	
Serum Albumin (human) Biotin-Free	HSA	AL363C	AL363F		AL363HV
Serum Albumin (mouse)	Albumin (mouse)	AL508C	AL508F	AL508S	
Anti-Peg IgM Kit	IgM (Anti-PEG)	AL313C	AL313F		AL313HV
CHO Host Cell Proteins	CHO-P	AL210C	AL210F	AL210S	
CHO Host Cell Proteins (broad reactivity)	CHO-P (broad reactivity)	AL301C	AL301F	AL301S	
Host Cell Residual DNA	Residual DNA	AL331C	AL331F	AL331S	AL331HV
PER.C6 Host Cell Proteins	PER.C6 HCP	AL302C	AL302F	AL302S	
Residual Protein A	Residual Protein A	AL287C	AL287F	AL287S	

Ig Detection

Product Description		Part Number	Part Number	Part Number	Part Number
		500 data points	5,000 data points	Standard	100 data points
Immunoglobulin E (mouse)	IgE (mouse)	AL526C	AL526F	AL526S	AL526HV
Immunoglobulin G (human)	IgG (human)	AL205C	AL205F	AL205S	
Immunoglobulin G (mouse)	IgG (mouse)	AL529C	AL529F	AL529S	AL529HV
Immunoglobulin G1 (mouse)	IgG1 (mouse)	AL522C	AL522F	AL522S	AL522HV
Immunoglobulin G2a (mouse)	IgG2a (mouse)	AL523C	AL523F	AL523S	AL523HV
Immunoglobulin G2b (mouse)	IgG2b (mouse)	AL524C	AL524F	AL524S	AL524HV
Immunoglobulin G3 (mouse)	IgG3 (mouse)	AL525C	AL525F	AL525S	AL525HV
Immunoglobulin G4 (human)	IgG4 (human)	AL310C	AL310F	AL310S	AL310HV
Immunoglobulin G4 (pharmacokinetic) (human)	IgG4 (human)	AL304C	AL304F	AL304S	
Immunoglobulin M (human)	IgM (human)	AL263C	AL263F	AL263S	
Immunoglobulin M (mouse)	IgM (mouse)	AL527C	AL527F	AL527S	AL527HV
Immunoglobulin Kappa light chain	Kappa light chain	AL3023C	AL3023F	AL3023S	AL3023HV
Lambda light chain	Lambda light chain	AL3042C	AL3042F	AL3042S	AL3042HV

AlphaLISA Binding Kits

Product Description	Part Number		
	500 data points	5,000 data points	100 data points
4-1BB/4-1BBL	AL3113C	AL3113F	
CTLA4/CD80	AL3046C	AL3046F	
CTLA4/CD86	AL3047C	AL3047F	
EGF/EGFR	AL366C	AL366F	AL366HV
FcGR1 (CD64)	AL3081C	AL3081F	
FcGR2A (CD32a), 167H	AL3086C	AL3086F	
FcGR2A (CD32a), 167R	AL3087C	AL3087F	
FcGR2B (CD32b)	AL3080C	AL3080F	
FcGR3A (CD16a), 176Phe/F158	AL347C	AL347F	AL347HV
FcGR3A (CD16a), 176Val/V158	AL348C	AL348F	AL348HV
FcγR3B (CD16b)	AL3094C	AL3094F	
FcRn	AL3095C	AL3095F	
HLA-DRA/LAG-3	AL3066C	AL3066F	
PD-1/PD-L1	AL356C	AL356F	AL356HV
PD-1/PD-L1 (mouse)	AL580C	AL580F	
PDGF-BB/PDGFRα	AL3064C	AL3064F	
PDGF-BB/PDGFRβ	AL3065C	AL3065F	
PCSK9/LDLR	AL3119C	AL3119F	

AlphaLISA Epigenetic Biochemical Assay Reagents

Product Description	Part Number		
	250 µg	5 mg	25 mg
Anti-acetyl-Histone H3 Lysine 9 (H3K9ac) Acceptor Beads	AL114C	AL114M	AL114R
Anti-methyl-Histone H3 Lysine 4 (H3K4me1-2) Acceptor Beads	AL116C	AL116M	AL116R
Anti-methyl-Histone H3 Lysine 9 (H3K9me2) Acceptor Beads	AL117C	AL117M	AL117R
Biotinylated anti-Histone H3 (C-ter) Antibody	AL118C	AL118M	AL118R
Anti-unmodified-Histone H3 Lysine 4 (H3K4) Acceptor Beads	AL119C	AL119M	AL119R
Anti-acetyl-Histone H3 Lysine 27 (H3K27ac) Acceptor Beads	AL120C	AL120M	AL120R
Anti-di/mono-methyl-Histone H3 Lysine 27 (H3K27me2-1) Acceptor Beads	AL121C	AL121M	AL121R
Anti-tri-methyl-Histone H3 Lysine 27 (H3K27me3) Acceptor Beads	AL122C	AL122M	AL122R
Anti-di-methyl-Histone H3 Lysine 36 (H3K36me2) Acceptor Beads	AL123C	AL123M	AL123R
Anti-unmodified Histone H3 Lysine 9/ Lysine 27 (H3K9/K27) Acceptor Beads	AL138C	AL138M	AL138R
Anti-methyl-Histone H3 Arginine 2 (H3R2me) Acceptor Beads	AL139C	AL139M	AL139R
Anti-Histone H3 (C-ter) Acceptor Beads	AL147C	AL147M	AL147R
Biotinylated anti-di-methyl-Histone H3 Lysine 79 (H3K79me2) Antibody	AL148C	AL148M	AL148R
Anti-acetyl-Histone H4 Lysine pan Acceptor Beads	AL143C	AL143M	AL143R
Anti-mono-methyl-Histone H4 Lysine 20 (H4K20me1) Acceptor Beads	AL145C	AL145M	AL145R
Biotinylated anti-Histone H4 Antibody	AL146C	AL146M	AL146R
Anti-methyl-Histone H4 Arginine 3 (H4R3me) Acceptor Beads	AL150C	AL150M	AL150R
Anti-methyl-Histone H3/H4 Arginine pan Acceptor Beads	AL151C	AL151M	AL151R
Anti-acetyl-p53 Lysine 382 (p53K382ac) Acceptor Beads	AL124C	AL124M	AL124R
Anti-O-GlcNAc Acceptor Beads	AL144C	AL144M	AL144R

AlphaLISA® Epigenetic Cellular Detection Kits

Product Description	Part Number	Part Number	Part Number
	500 data points	5,000 data points	100 data points
Total H2a	AL725C	AL725F	AL725HV
Acetyl-Histone H3 Lysine 9 (H3K9ac)	AL714C	AL714F	AL714HV
Di-methyl-Histone H3 Lysine 4 (H3K4me2)	AL716C	AL716F	AL716HV
Di-methyl-Histone H3 Lysine 9 (H3K9me2)	AL717C	AL717F	AL717HV
Unmodified Histone H3 Lysine 4 (H3K4)*	AL719C	AL719F	AL719HV
Acetyl-Histone H3 Lysine 27 (H3K27ac)	AL720C	AL720F	AL720HV
Di/mono-methyl-Histone H3 Lysine 27 (H3K27me2-1)	AL721C	AL721F	AL721HV
Tri-methyl-Histone H3 Lysine 27 (H3K27me3)	AL722C	AL722F	AL722HV
Di-methyl-Histone H3 Lysine 36 (H3K36me2)	AL723C	AL723F	AL723HV
Di-methyl-Histone H3 Lysine 79 (H3K79me2)	AL748C	AL748F	AL748HV

*Suitable for measuring total H3 content

Alpha SureFire Kits

For AlphaScreen® *SureFire*® products, the corresponding AlphaScreen *SureFire* IgG (Protein A based) Detection Kit needs to be ordered in addition to the AlphaScreen *SureFire* Phospho or total kit. Please note that the sum of the prices of both the AlphaScreen *SureFire* and the IgG detection kits is superior to the price of the corresponding AlphaLISA *SureFire Ultra* kit. If you have not already shifted to the AlphaLISA *SureFire Ultra* we highly recommend to do that shift as, in addition to cost-savings, there is on average a 10x gain in sensitivity and dynamic range for the AlphaLISA *SureFire Ultra* assays compared to the AlphaScreen *SureFire* assays.

For AlphaLISA® *SureFire*® *Ultra*™, the Alpha Donor and Acceptor beads are included in the kits.

The Alpha *SureFire Ultra* Multiplex dual targets kits contain all the reagents necessary to measure two targets from the same well, using the AlphaPlex technology.

The Alpha terbium *SureFire Ultra* kits are intended to be combined with some AlphaLISA *SureFire Ultra* kits in order to multiplex the measurement of two analytes. The terbium *SureFire Ultra* kits include the antibody-containing Reaction Buffers and the beads and cannot be used as standalone kits because the cell lysis buffer, the activation buffer and the dilution buffer are provided only in the AlphaLISA *SureFire Ultra* kits.

If you need a target in the AlphaLISA *SureFire Ultra* or in the terbium *SureFire Ultra* format and do not find it in the listing below, please contact us – we may already have developed it recently. We would also be happy to discuss the possibility of doing this on a custom basis.

All kits are available in three sizes for immunoassays to be performed in 384-well plate format, and the *Ultra* are also available in “High Volume” format: these kits contain 100 data points, in a convenient 96-well plate assay format, for low throughput applications.

Global Overview of Alpha *SureFire* Kits by Pathway

PI3K/Akt/mTOR Pathway	JAK/STAT Pathway	NFκB Pathway	MAPK Pathway	Receptors	TGFβ/BMP/SMAD Pathway	Others
4E-BP1	STAT1	IκBα	ATF-1	ALK	SMAD1	Acetyl CoA carboxylase
Akt	STAT3	IKKα	ATF-2	c-Met	SMAD2	AMPKα 1/2
BAD	STAT4	IKKβ	c-Jun	EGF Receptor	SMAD3	Beta catenin
Caspase-9	STAT5	NFκB p65	ELK1	ErbB2		Chk1
eIF4E	STAT6		ERK 1/2	IGF-1 Receptor β		Chk2
GSK-3α			HSP27	Insulin Receptor β		CREB
GSK-3β			JNK1/2/3	TrkA/B		Cofilin
mTOR			MEK1	VEGF Receptor 2		eIF2α
p70S6K			MKK3/6			GAPDH
PDK1			MKK4			Histone H3
S6RP			P38 MAPK			P53
						SLP-76
						SYK
						Topoisomerase IIα

Alphabetical Listing of Alpha *SureFire* Products

Product Description	Part Number						
	AlphaScreen <i>SureFire</i>			AlphaLISA <i>SureFire Ultra</i>			
	500 data points	10,000 data points	50,000 data points	500 data points	10,000 data points	50,000 data points	100 datapoints
Phospho-4EBP1 (Thr37/46)	TGR4ES500	TGR4ES10K	TGR4ES50K	ALSU-P4EBP-A500	ALSU-P4EBP-A10K	ALSU-P4EBP-A50K	ALSU-P4EBP-A-HV
Phospho-4EBP1 (Thr70)	TGR4E2S500	TGR4E2S10K	TGR4E2S50K				
Phospho-acetyl CoA carboxylase (Ser79)				ALSU-PACC-A500	ALSU-PACC-A10K	ALSU-PACC-A50K	ALSU-PACC-A-HV
Phospho-Akt1 (Thr308)	TGRA2S500	TGRA2S10K	TGRA2S50K				
Phospho-Akt1 (Ser473)	TGRA5S500	TGRA5S10K	TGRA5S50K				
Total Akt1	TGRTA1S500	TGRTA1S10K	TGRTA1S50K	ALSU-TAKT1-A500	ALSU-TAKT1-A10K	ALSU-TAKT1-A50K	ALSU-TAKT1-A-HV
Phospho-Akt 1/2/3 (Thr308)	TGRA3S500	TGRA3S10K	TGRA3S50K	ALSU-PAKT-A500	ALSU-PAKT-A10K	ALSU-PAKT-A50K	ALSU-PAKT-A-HV
Phospho-Akt1/2/3 (Ser473)	TGRAS500	TGRAS10K	TGRAS50K				
Phospho-Akt 1/2/3 (Ser473)	TGRA4S500	TGRA4S10K	TGRA4S50K	ALSU-PAKT-B500	ALSU-PAKT-B10K	ALSU-PAKT-B50K	ALSU-PAKT-B-HV
Total Akt 1/2/3	TGRTAPS500	TGRTAPS10K	TGRTAPS50K	ALSU-TAKT-B500	ALSU-TAKT-B10K	ALSU-TAKT-B50K	ALSU-TAKT-B-HV
Phospho-ALK (Tyr1586)	TGRALS500	TGRALS10K	TGRALS50K				
Phospho-ALK (Tyr1604)	TGRAL2S500	TGRAL2S10K	TGRAL2S50K				
Total ALK	TGRTALS500	TGRTALS10K	TGRTALS50K				

Product Description	Part Number						
	AlphaScreen <i>SureFire</i>			AlphaLISA <i>SureFire Ultra</i>			
	500 data points	10,000 data points	50,000 data points	500 data points	10,000 data points	50,000 data points	100 data points
Phospho-AMPK α 1/2 (Thr172)				ALSU-PAMPK-A500	ALSU-PAMPK-A10K	ALSU-PAMPK-A50K	ALSU-PAMPK-A-HV
Phospho-ATF-1 (Ser63)	TGRAF1S500	TGRAF1S10K	TGRTALS50K				
Phospho-ATF-2 (Thr71)	TGRAF2S500	TGRAF2S10K	TGRAF2S50K				
Phospho-BAD (Ser112)	TGRBS500	TGRBS10K	TGRBS50K				
Phospho-BAD (Ser136)	TGRB2S500	TGRB2S10K	TGRB2S50K				
Phospho-beta-catenin (Ser45)	TGRBCS500	TGRBCS10K	TGRBCS50K				
Phospho-Caspase 9 (Ser196)	TGRC9S500	TGRC9S10K	TGRC9S50K				
Phospho-Chk1 (Ser345)	TGRCHK1S500	TGRCHK1S10K	TGRCHK1S50K				
Phospho-Chk2 (Thr68)	TGRCHK2S500	TGRCHK2S10K	TGRCHK2S50K				
Phospho-c-Jun (Ser63)	TGRCJS500	TGRCJS10K	TGRCJS50K				
Phospho-c-Jun (Ser73)	TGRCJ2S500	TGRCJ2S10K	TGRCJ2S50K				
Phospho-c-Met (Tyr1234/1235)	TGRCMS500	TGRCMS10K	TGRCMS50K	ALSU-PCMET-A500	ALSU-PCMET-A10K	ALSU-PCMET-A50K	ALSU-PCMET-A-HV
Phospho-cofilin (Ser3)	TGRCOFS500	TGRCOFS10K	TGRCOFS50K	ALSU-PCOF-A500	ALSU-PCOF-A10K	ALSU-PCOF-A50K	ALSU-PCOF-A-HV
Total cofilin				ALSU-TCOF-A500	ALSU-TCOF-A10K	ALSU-TCOF-A50K	ALSU-TCOF-A-HV
Phospho-CREB-1 (Ser133)	TGRCBS500	TGRCBS10K	TGRCBS50K	ALSU-PCREB-A500	ALSU-PCREB-A10K	ALSU-PCREB-A50K	ALSU-PCREB-A-HV
Phospho-EGF Receptor (Tyr1068)	TGRERS500	TGRERS10K	TGRERS50K	ALSU-PEGFR-B500	ALSU-PEGFR-B10K	ALSU-PEGFR-B50K	ALSU-PEGFR-B-HV
Total EGF Receptor				ALSU-TEGFR-A500	ALSU-TEGFR-A10K	ALSU-TEGFR-A50K	ALSU-TEGFR-A-HV
Phospho-eIF2 α (Ser51)				ALSU-PEIF2-B500	ALSU-PEIF2-B10K	ALSU-PEIF2-B50K	ALSU-PEIF2-B-HV
Total eIF2 α				ALSU-TEIF2-A500	ALSU-TEIF2-A10K	ALSU-TEIF2-A50K	ALSU-TEIF2-A-HV
Phospho-eIF4E (Ser209)	TGREIF4S500	TGREIF4S10K	TGREIF4S50K	ALSU-PEIF4-A500	ALSU-PEIF4-A10K	ALSU-PEIF4-A50K	ALSU-PEIF4-A-HV
Phospho-Elk-1 (Ser383)	TGRELKS500	TGRELKS10K	TGRELKS50K				
Phospho-ErbB2 (Tyr1221/1222)	TGREB2S500	TGREB2S10K	TGREB2S50K	ALSU-PEB2-A500	ALSU-PEB2-A10K	ALSU-PEB2-A50K	ALSU-PEB2-A-HV
Phospho-ERK1/2 (Thr202/Tyr204)	TRGESB500	TGRESB10K	TGRESB50K	ALSU-PERK-A500	ALSU-PERK-A10K	ALSU-PERK-A50K	ALSU-PERK-A-HV
Total ERK1/2	TGRTEB500	TGRTEB10K	TGRTEB50K	ALSU-TERK-A500	ALSU-TERK-A10K	ALSU-TERK-A50K	ALSU-TERK-A-HV
Total GAPDH (human)	TGRGDS500	TGRGDS10K	TGRGDS50K	ALSU-TGAPD-A500	ALSU-TGAPD-A10K	ALSU-TGAPD-A50K	ALSU-TGAPD-A-HV
Total GAPDH (mouse & human)				ALSU-TGAPD-B500	ALSU-TGAPD-B10K	ALSU-TGAPD-B50K	ALSU-TGAPD-B-HV
Phospho-GSK3 α (Ser 21)	TGRGAS500	TGRGAS10K	TGRGAS50K				
Phospho-GSK3 β (Ser9)	TGRGBS500	TGRGBS10K	TGRGBS50K	ALSU-PGS3B-A500	ALSU-PGS3B-A10K	ALSU-PGS3B-A50K	ALSU-PGS3B-A-HV
Phospho-Histone H3 (Ser10)*	TGRH3S500	TGRH3S10K	TGRH3S50K				
Phospho-HSP27 (Ser82)	TGRHS27S500	TGRHS27S10K	TGRHS27S50K				
Phospho-IGF1 Receptor β (Tyr1135/1136)				ALSU-PIGFR-B500	ALSU-PIGFR-B10K	ALSU-PIGFR-B50K	ALSU-PIGFR-B-HV
Phospho-I κ B α (Ser32/36)	TGRIKS500	TGRIKS10K	TGRIKS50K				
Total I κ B α	TGRTIKS500	TGRTIKS10K	TGRTIKS50K				
Phospho-IKK α (Ser176/180)	TGRKAS500	TGRKAS10K	TGRKAS50K	ALSU-PIKKA-A500	ALSU-PIKKA-A50K	ALSU-PIKKA-A10K	ALSU-PIKKA-A-HV
Phospho-IKK β (Ser177/181)	TGRKBS500	TGRKBS10K	TGRKBS50K				
Phospho-Insulin Receptor β (Tyr1150/1151)	TGRIIRS500	TGRIIRS10K	TGRIIRS50K	ALSU-PINR-A500	ALSU-PINR-A10K	ALSU-PINR-A50K	ALSU-PINR-A-HV
Phospho-JNK1/3 (Thr183/Tyr185)	TGRJS500	TGRJS10K	TGRJS50K				

* Please refer to our Epigenetics Listing for the total H3 (Unmodified H3K4 Cellular Kit) AL719

Product Description				Part Number			
AlphaScreen <i>SureFire</i>				AlphaLISA <i>SureFire Ultra</i>			
	500 data points	10,000 data points	50,000 data points	500 data points	10,000 data points	50,000 data points	100 data points
Phospho-JNK1/2/3 (Thr183/Tyr185)	TGRJPS500	TGRJPS10K	TGRJPS50K	ALSU-PJNK-A500	ALSU-PJNK-A10K	ALSU-PJNK-A50K	ALSU-PJNK-A-HV
Total JNK1/2/3	TGRTJS500	TGRTJS10K	TGRTJS50K				
Phospho-MEK-1 (Ser218/222)				ALSU-PMEK1-A500	ALSU-PMEK1-A10K	ALSU-PMEK1-A50K	ALSU-PMEK1-A-HV
Total MEK-1				ALSU-TMEK1-A500	ALSU-TMEK1-A10K	ALSU-TMEK1-A50K	ALSU-TMEK1-A-HV
Phospho-MKK3/6 (Ser189/Ser207)	TGRMK3S500	TGRMK3S10K	TGRMK3S50K				
Phospho-MKK4 (Ser257)				ALSU-PMMK4-A-HV	ALSU-PMMK4-A500	ALSU-PMMK4-A10K	ALSU-PMMK4-A50K
Phospho-MKK4 (Ser257/Thr261)	TGRMK4S500	TGRMK4S10K	TGRMK4S50K				
Total MKK4				ALSU-TMKK4-A-HV	ALSU-TMKK4-A500	ALSU-TMKK4-A10K	ALSU-TMKK4-A50K
Phospho-mTOR (Ser2448)				ALSU-PMTOR-A500	ALSU-PMTOR-A10K	ALSU-PMTOR-A50K	ALSU-PMTOR-A-HV
Phospho-mTOR (Ser2481)				ALSU-PMTOR-B-HV	ALSU-PMTOR-B500	ALSU-PMTOR-B10K	ALSU-PMTOR-B50K
Total mTOR				ALSU-TMTOR-A500	ALSU-TMTOR-A10K	ALSU-TMTOR-A50K	ALSU-TMTOR-A-HV
Phospho-NFκB p65 (Ser536)	TGRNFS500	TGRNFS10K	TGRNFS50K	ALSU-PNFKB-A500	ALSU-PNFKB-A10K	ALSU-PNFKB-A50K	ALSU-PNFKB-A-HV
Total NFκB p65	TGRTNFS500	TGRTNFS10K	TGRTNFS50K	ALSU-TNFKB-A500	ALSU-TNFKB-A10K	ALSU-TNFKB-A50K	ALSU-TNFKB-A-HV
Phospho-p38 MAPK (Thr180/Tyr182)				ALSU-PP38-B500	ALSU-PP38-B10K	ALSU-PP38-B50K	ALSU-PP38-B-HV
Total p38 MAPK				ALSU-TP38-B500	ALSU-TP38-B10K	ALSU-TP38-B50K	ALSU-TP38-B-HV
Phospho-p53 (Ser15)	TGRP53S500	TGRP53S10K	TGRP53S50K				
Phospho-p53 (Ser392)	TGRP532S500	TGRP532S10K	TGRP532S50K				
Total p53	TGRT53S500	TGRT53S10K	TGRT53S50K				
Phospho-p70S6K (Thr229)	TGR703S500	TGR703S10K	TGR703S50K				
Phospho-p70S6K (Thr389)	TGR70S500	TGR70S10K	TGR70S50K	ALSU-PP70-A500	ALSU-PP70-A10K	ALSU-PP70-A50K	ALSU-PP70-A-HV
Phospho-p70S6K (Thr421/Ser424)	TGR702S500	TGR702S10K	TGR702S50K				
Total p70S6K	TGRT70S500	TGRT70S10K	TGRT70S50K	ALSU-TP90-A500	ALSU-TP90-A10K	ALSU-TP90-A50K	ALSU-TP90-A-HV
Phospho-PDK1 (Ser241)	TGRPS500	TGRPS10K	TGRPS50K				
Phospho-S6 RP (Ser235/236)	TGRS6PS500	TGRS6PS10K	TGRS6PS50K				
Phospho-S6 RP (Ser240/244)	TGRS6P2S500	TGRS6P2S10K	TGRS6P2S50K	ALSU-PS6R-A500	ALSU-PS6R-A10K	ALSU-PS6R-A50K	ALSU-PS6R-A-HV
Phospho-SHP1 (Tyr536)				ALSU-PSHP1-A500	ALSU-PSHP1-A10K	ALSU-PSHP1-A50K	ALSU-PSHP1-A-HV
Phospho-SHP1 (Tyr 564)				ALSU-PSHP1-B500	ALSU-PSHP1-B10K	ALSU-PSHP1-B50K	ALSU-PSHP1-B-HV
Phospho-SHP2 (Tyr542)				ALSU-PSHP2-A500	ALSU-PSHP2-A10K	ALSU-PSHP2-A50K	ALSU-PSHP2-A-HV
Phospho-SLP-76 (Ser376)				ALSU-PSLP-A500	ALSU-PSLP-A10K	ALSU-PSLP-A50K	ALSU-PSLP-A-HV
Total SLP-76				ALSU-TSLP-A500	ALSU-TSLP-A10K	ALSU-TSLP-A50K	ALSU-TSLP-A-HV
Phospho-SMAD1 (Ser463/465)	TGRSM1S500	TGRSM1S10K	TGRSM1S50K	ALSU-PSM1-A500	ALSU-PSM1-A10K	ALSU-PSM1-A50K	ALSU-PSM1-A-HV
Phospho-SMAD2 (Ser465/467)	TGRSM2S500	TGRSM2S10K	TGRSM2S50K	ALSU-PSM2-A500	ALSU-PSM2-A10K	ALSU-PSM2-A50K	ALSU-PSM2-A-HV
Total SMAD2				ALSU-TSM2-A500	ALSU-TSM2-A10K	ALSU-TSM2-A50K	ALSU-TSM2-A-HV

Product Description				Part Number			
AlphaScreen <i>SureFire</i>				AlphaLISA <i>SureFire Ultra</i>			
500 data points		10,000 data points	50,000 data points	500 data points	10,000 data points	50,000 data points	100 data points
Phospho-SMAD3 (Ser423/425)	TGRSM3S500	TGRSM3S10K	TGRSM3S50K	ALSU-TSM2-A10K	ALSU-PSM3-A10K	ALSU-PSM3-A50K	ALSU-PSM3-A-HV
Phospho-Src (Tyr419)				ALSU-TSM2-A50K	ALSU-PSRC-A500	ALSU-PSRC-A10K	ALSU-PSRC-A50K
Phospho-STAT1 (Tyr701)	TGRS1S500	TGRS1S10K	TGRS1S50K	ALSU-TSM2-A-HV	ALSU-PST1-A10K	ALSU-PST1-A50K	ALSU-PST1-A-HV
Phospho-STAT1 (Ser727)				ALSU-PST1-B500	ALSU-PST1-B10K	ALSU-PST1-B50K	ALSU-PST1-B-HV
Phospho-STAT3 (Tyr705)	TGRS3S500	TGRS3S10K	TGRS3S50K	ALSU-PST3-A500	ALSU-PST3-A10K	ALSU-PST3-A50K	ALSU-PST3-A-HV
Total STAT3				ALSU-TST3-A500	ALSU-TST3-A10K	ALSU-TST3-A50K	ALSU-TST3-A-HV
Phospho-STAT4 (Tyr693)				ALSU-PST4-A500	ALSU-PST4-A10K	ALSU-PST4-A50K	ALSU-PST4-A-HV
Phospho-STAT5 (Tyr694/699)	TGRS5S500	TGRS5S10K	TGRS5S50K	ALSU-PST5-B500	ALSU-PST5-B10K	ALSU-PST5-B50K	ALSU-PST5-B-HV
Total STAT5				ALSU-TST5-A500	ALSU-TST5-A10K	ALSU-TST5-A50K	ALSU-TST5-A-HV
Phospho-STAT6 (Tyr641)				ALSU-PST6-A500	ALSU-PST6-A10K	ALSU-PST6-A50K	ALSU-PST6-A-HV
Phospho-Syk (Tyr352)				ALSU-PSYK-B-HV	ALSU-PSYK-B500	ALSU-PSYK-B10K	ALSU-PSYK-B50K
Phospho-SYK (Tyr525/526)				ALSU-PSYK-A500	ALSU-PSYK-A10K	ALSU-PSYK-A50K	ALSU-PSYK-A-HV
Total SYK				ALSU-TSYK-A500	ALSU-TSYK-A10K	ALSU-TSYK-A50K	ALSU-TSYK-A-HV
Phospho-Topoisomerase II α (Ser1106)	TGRT2A2S500	TGRT2A2S10K	TGRT2A2S50K				
Phospho-Topoisomerase II α (Thr1343)	TGRT2AS500	TGRT2AS10K	TGRT2AS50K				
Total Topoisomerase II α	TGRTT2AS500	TGRTT2AS10K	TGRTT2AS50K				
Phospho-TrkA (Tyr674/675) / TrkB (Tyr706/707)				ALSU-PTRKAB-A500	ALSU-PTRKAB-A10K	ALSU-PTRKAB-A50K	ALSU-PTRKAB-A-HV
Phospho-VEGFR2 (Tyr951)				ALSU-PVGFR-B-HV	ALSU-PVGFR-B500	ALSU-PVGFR-B10K	ALSU-PVGFR-B50K
Phospho-VEGFR2 (Tyr1214)				ALSU-PVGFR-C-HV	ALSU-PVGFR-C500	ALSU-PVGFR-C10K	ALSU-PVGFR-C50K
Phospho-VEGFR2 (Tyr1175)	TGRVRS500	TGRVRS10K	TGRVRS50K	ALSU-PVGFR-A500	ALSU-PVGFR-A10K	ALSU-PVGFR-A50K	ALSU-PVGFR-A-HV
Total VEGFR2				ALSU-TVGFR-A500	ALSU-TVGFR-A10K	ALSU-TVGFR-A50K	ALSU-TVGFR-A-HV
Phospho-ZAP-70 (Tyr319)				ALSU-PZAP-A-HV	ALSU-PZAP-A500	ALSU-PZAP-A10K	ALSU-PZAP-A50K
Total ZAP-70				ALSU-TZAP-A-HV	ALSU-TZAP-A500	ALSU-TZAP-A10K	ALSU-TZAP-A50K

Alpha *SureFire Ultra* Multiplex

Product Description		Part Number	Part Number	Part Number	Part Number
		100 data points	500 data points	10,000 data points	50,000 data points
p-ERK (Eu)/Total ERK (Tb)	Dual Target	MPSU-PTERK-K-HV	MPSU-PTERK-K500	MPSU-PTERK-K10K	MPSU-PTERK-K50K
p-Akt 1/2/3 (S473) (Eu)/Total Akt1 (Tb)	Dual Target	MPSU-PTAKT-K-HV	MPSU-PTAKT-K500	MPSU-PTAKT-K10K	MPSU-PTAKT-K50K

Terbium *SureFire Ultra*

The terbium *SureFire Ultra* (TBSU) kits measure a single phosphoprotein target and are intended to be combined with a standard AlphaLISA *SureFire Ultra* kit (ALSU) to assess the second target. The TBSU kits allow for a pick-and-pack combination of any two phosphoprotein targets in the assay lists, providing measurement of two separate targets with ultimate flexibility of selection. The TBSU kits will report the target of interest through the Terbium (Tb) Acceptor beads, while the ALSU kit reports via the Europium (Eu) AlphaLISA Acceptor beads.

Product Description	Part Number	Part Number	Part Number	Part Number
	500 data points	10,000 data points	50,000 data points	100 data points
Phospho-4EBP1 (Thr37/46)	TBSU-P4EBP-A500	TBSU-P4EBP-A10K	TBSU-P4EBP-A50K	TBSU-P4EBP-A-HV
Phospho-Akt1/2/3 (Thr308)	TBSU-PAKT-A500	TBSU-PAKT-A10K	TBSU-PAKT-A50K	TBSU-PAKT-A-HV
Phospho-Akt1/2/3 (Ser473)	TBSU-PAKT-B500	TBSU-PAKT-B10K	TBSU-PAKT-B50K	TBSU-PAKT-B-HV
Total cofilin	TBSU-TCOF-A500	TBSU-TCOF-A10K	TBSU-TCOF-A50K	TBSU-TCOF-A-HV
Phospho-CREB-1 (Ser133)	TBSU-PCREB-A500	TBSU-PCREB-A10K	TBSU-PCREB-A50K	TBSU-PCREB-A-HV
Phospho-EGF Receptor (Tyr1068)	TBSU-PEGFR-A500	TBSU-PEGFR-A10K	TBSU-PEGFR-A50K	TBSU-PEGFR-A-HV
Phospho-eIF2alpha (Ser51)	TBSU-PEIF2-B500	TBSU-PEIF2-B10K	TBSU-PEIF2-B50K	TBSU-PEIF2-B-HV
Phospho-eIF4E (Ser209)	TBSU-PEIF4-A500	TBSU-PEIF4-A10K	TBSU-PEIF4-A50K	TBSU-PEIF4-A-HV
Phospho-ERK1/2 (T202/Y204)	TBSU-PERK-A500	TBSU-PERK-A10K	TBSU-PERK-A50K	TBSU-PERK-A-HV
Phospho-GSK3β (Ser9)	TBSU-PGS3B-A500	TBSU-PGS3B-A10K	TBSU-PGS3B-A50K	TBSU-PGS3B-A-HV
Phospho-IGF1 Receptor β (Tyr1135/1136)	TBSU-PIGFR-B500	TBSU-PIGFR-B10K	TBSU-PIGFR-B50K	TBSU-PIGFR-B-HV
Phospho-IKKα (Ser176/180)	TBSU-PIKKA-A500	TBSU-PIKKA-A10K	TBSU-PIKKA-A50K	TBSU-PIKKA-A-HV
Phospho-Insulin Receptor β (Tyr1150/1151)	TBSU-PINR-A500	TBSU-PINR-A10K	TBSU-PINR-A50K	TBSU-PINR-A-HV
Phospho-JNK1/2/3 (Thr183/Tyr185)	TBSU-PJNK-A500	TBSU-PJNK-A10K	TBSU-PJNK-A50K	TBSU-PJNK-A-HV
Phospho-MEK1 (Ser218/222)	TBSU-PMEK1-A500	TBSU-PMEK1-A10K	TBSU-PMEK1-A50K	TBSU-PMEK1-A-HV
Phospho-mTOR (Ser2448)	TBSU-PMTOR-A500	TBSU-PMTOR-A10K	TBSU-PMTOR-A50K	TBSU-PMTOR-A-HV
Phospho-NFκB p65 (Ser536)	TBSU-PNFKB-A500	TBSU-PNFKB-A10K	TBSU-PNFKB-A50K	TBSU-PNFKB-A-HV
Phospho-p38 MAPK (Thr180/Tyr182)	TBSU-PP38-B500	TBSU-PP38-B10K	TBSU-PP38-B50K	TBSU-PP38-B-HV
Phospho-p70S6K (Thr389)	TBSU-PP70-A500	TBSU-PP70-A10K	TBSU-PP70-A50K	TBSU-PP70-A-HV
Phospho-RPS6 (Ser240/244)	TBSU-PS6R-A500	TBSU-PS6R-A10K	TBSU-PS6R-A50K	TBSU-PS6R-A-HV
Phospho-SLP76 (Ser376)	TBSU-PSLP-A500	TBSU-PSLP-A10K	TBSU-PSLP-A50K	TBSU-PSLP-A-HV
Phospho-SMAD1 (Ser463/465)	TBSU-PSM1-A500	TBSU-PSM1-A10K	TBSU-PSM1-A50K	TBSU-PSM1-A-HV
Phospho-SMAD3 (Ser423/425)	TBSU-PSM3-A500	TBSU-PSM3-A10K	TBSU-PSM3-A50K	TBSU-PSM3-A-HV
Phospho-STAT1 (Tyr701)	TBSU-PST1-A500	TBSU-PST1-A10K	TBSU-PST1-A50K	TBSU-PST1-A-HV
Phospho-STAT1 (Ser727)	TBSU-PST1-B500	TBSU-PST1-B10K	TBSU-PST1-B50K	TBSU-PST1-B-HV
Phospho-STAT3 (Tyr705)	TBSU-PST3-A500	TBSU-PST3-A10K	TBSU-PST3-A50K	TBSU-PST3-A-HV
Phospho-STAT4 (Tyr693)	TBSU-PST4-A500	TBSU-PST4-A10K	TBSU-PST4-A50K	TBSU-PST4-A-HV
Phospho-STAT5 (Tyr694/699)	TBSU-PST5-B500	TBSU-PST5-B10K	TBSU-PST5-B50K	TBSU-PST5-B-HV
Phospho-STAT6 (Tyr641)	TBSU-PST6-A500	TBSU-PST6-A10K	TBSU-PST6-A50K	TBSU-PST6-A-HV
Phospho-SYK (Tyr525/526)	TBSU-PSYK-A500	TBSU-PSYK-A10K	TBSU-PSYK-A50K	TBSU-PSYK-A-HV
Phospho-VEGF Receptor 2 (Tyr1175)	TBSU-PVGFR-A500	TBSU-PVGFR-A10K	TBSU-PVGFR-A50K	TBSU-PVGFR-A-HV

AlphaScreen *SureFire* IgG (Protein A Based) Detection Kits

These kits contain Protein A Acceptor Beads and Streptavidin Donor Beads that are needed when working with AlphaScreen *SureFire* kits.

Product Description	Part Number	Part Number	Part Number	Part Number
	500 data points	10,000 data points	50,000 data points	100 data points
AlphaScreen General IgG (Protein A based) Detection Kit (384-well format)	6760617C	6760617M	6760617R	
AlphaScreen <i>SureFire</i> HV IgG (Protein A) Detection Kit (96-well format)				6760617HV

AlphaScreen *SureFire* Supplementary Buffers

Enough buffer is provided in the kits if following the standard protocol, but you may want to order more if using a different protocol.

Product Description	Part Number	
	10 mL	100 mL
<i>SureFire</i> Lysis Buffer (5X)	TGRLB10ML	TGRLB100ML
<i>SureFire</i> Activation Buffer (50X)	TGRAB10ML	TGRAB100ML
<i>SureFire</i> Activation Buffer B (50X)	TGRABB10ML	TGRABB100ML
<i>SureFire</i> Activation Buffer C (50X)	TGRABC10ML	TGRABC100ML
<i>SureFire</i> Dilution Buffer (1X)	TGRDB10ML	TGRDB100ML

AlphaScreen *SureFire* Supplementary Control Lysates

Enough control lysates are provided in the kits if following the standard protocol, but if you need more for specific applications please contact us to request custom AlphaScreen *SureFire* control lysates.

AlphaLISA *SureFire Ultra* Supplementary Buffers

Enough buffer is provided in the kits if following the standard protocol, but you may want to order more if using a different protocol.

Product Description	Part Number	
	10 mL	100 mL
<i>SureFire Ultra</i> Lysis Buffer (5X)	ALSU-LB-10mL	ALSU-LB-100mL
<i>SureFire Ultra</i> Activation Buffer (50X)	ALSU-AB-10mL	ALSU-AB-100mL
<i>SureFire Ultra</i> Activation Buffer B (50X)	ALSU-ABB-10mL	ALSU-ABB-100mL
<i>SureFire Ultra</i> Activation Buffer C (50X)	ALSU-ABC-10mL	ALSU-ABC-100mL
<i>SureFire Ultra</i> Dilution Buffer (1X)	ALSU-DB-10mL	ALSU-DB-100mL

AlphaLISA *SureFire Ultra* Supplementary Beads

Enough beads are provided in the kits if following the standard protocol, but you may want to order more if using a different protocol.

Product Description	Part Number		
	500 data points (60 µL)	10,000 data points (1.2 mL)	50,000 data points (6 mL)
Alpha Streptavidin Donor Beads (2 mg/mL)	ALSU-ASDB-0.06mL	ALSU-ASDB-1.2mL	ALSU-ASDB-6mL
AlphaLISA <i>SureFire Ultra</i> CaptSure Acceptor Beads	ALSU-ACAB-0.06mL	ALSU-ACAB-1.2mL	ALSU-ACAB-6mL

AlphaLISA *SureFire*® *Ultra* Kits Supplementary Control lysates

Enough control lysates are provided in the kits if following the standard protocol, but if you need more for specific applications, they are available separately. For example, the phospho-ERK control lysate Part# is ALSU-PERK-A-L. Please consult our WEB site to find the references for other kits control lysates.

Alpha *SureFire*® *Ultra* Multiplex Supplementary Beads

Enough beads are provided in the kits if following the standard protocol, but you may want to order more if using a different protocol. These beads can also be substituted to the AlphaLISA (Eu) beads in order to read the target in another wavelength channel.

Product Description	Part Number		
	500 data points (60 µL)	10,000 data points (1.2 mL)	50,000 data points (6 mL)
Alpha 545 (Tb) <i>SureFire Ultra</i> CaptSure Acceptor Beads	MPSU-ACAT-0.06mL	MPSU-ACAT-1.2mL	MPSU-ACAT-6mL

Beads for Biochemical Kinase Assays

Product Description	Part Number			
	500 data points	10K data points	50k data points	1K data points
Antibody-free PhosphoSensor Kit		6760307M	6760307R	6760307D
PY20 Assay Kit	6760601C	6760601M	6760601R	
PT66 Assay Kit	6760602C	6760602M	6760602R	

Alpha Technology Products

ToolBox beads, buffers, and control beads 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. Please contact your PerkinElmer account manager or PerkinElmer sales office for more information.

Alpha ToolBox

	Acceptor Beads				
	Donor Beads 10 ₂	AlphaLISA 615	AlphaPlex (Tb) 545	AlphaPlex (Sm) 645	AlphaScreen 520-620
Unconjugated	6762011/2/3	6772001/2/3	6782001/2/3	6792001/2/3	6762001/2/3
<i>Fusion Tag Detection</i>					
Streptavidin	6760002	AL125	AP125TB	AP125SM	
Strep Tactin®	AS106	AL136			
Anti-DIG	AS108	AL113			6760604
Anti-FITC		AL127			6760605
Anti-DNP	AS111	AL173			
Anti-HRP	AS109	AL171			
Nickel Chelate	AS101	AL108			6760619
Anti-6x His		AL128	AP128TB	AP128SM	
Glutathione (GSH)	6765300/1/2/3	AL109	AP109TB		
Anti-GST		AL110			6760603
Anti-c-myc		AL111			6760611
Anti-Protein C tag		AL172			
Anti-FLAG®	AS103	AL112	AP112TB	AP112SM	6760613
Anti-GFP	AS112	AL133			
Anti-HA		AL170	AP170TB		6760612
LCA (Lens Culinaris Agglutinin)		AL140			
Anti-Maltose-Binding Protein (MBP)		AL134			
Anti-V5		AL129			
<i>Antibody Capture</i>					
Protein A	AS102	AL101	AP101TB	AP101SM	6760137
Protein G		AL102			
Protein L		AL126			
Anti-bovine IgA		AL169			
Anti-chicken IgY		AL131			
Anti-goat IgG		AL107			
Anti-human IgE		AL176			
Anti-human IgG (Fc-specific)	AS114	AL103	AP103TB	AP103SM	
Anti-human IgG1		AL141			
Anti-human IgG2 (isotyping)		AL154			
Anti-human IgG4		AL142			
Anti-mouse IgG	AS104	AL105	AP105TB	AP105SM	6760606
Anti-mouse IgM		AL130			
Anti-rabbit IgG	AS105	AL104	AP104TB		6760607
Anti-rat IgG	AS110	AL106			
Anti-sheep IgG		AL132			
Human IgG (Fc-fragment)	AS113	AL174			
Human IgG (whole)		AL175			

Alpha Donor Beads

Product Description	Part Number	Part Number	Part Number	Part Number
	1 mg	5 mg	25 mg	50 mg
<i>Unconjugated</i>				
Unconjugated Donor Beads	6762013	6762011		6762012
<i>Tag capture</i>				
Streptavidin Donor Beads	6760002S	6760002		6760002B
Strep-Tactin® Donor Beads	AS106D	AS106M	AS106R	
Anti-DIG Donor Beads	AS108D	AS108M	AS108R	
Anti-DNP (dinitrophenol) Donor Beads	AS111D	AS111M	AS111R	
Anti-FITC Donor Beads	AS115D	AS115M	AS115R	
Anti-FLAG Donor Beads	AS103D	AS103M	AS103R	
Anti-GFP Donor Beads	AS112D	AS112M	AS112R	
Anti-HRP Donor Beads	AS109D	AS109M	AS109R	
GSH (glutathione) Donor Beads	6765300	6765301	6765301	
Nickel Chelate Donor Beads	AS101D	AS101M	AS101R	
<i>Antibody capture</i>				
Anti-human IgG Donor beads (Fc-specific)	AS114D	AS114M	AS114R	
Anti-Mouse IgG Donor Beads	AS104D	AS104M	AS104R	
Anti-Rabbit IgG Donor Beads	AS105D	AS105M	AS105R	
Anti-Rat IgG Donor Beads	AS110D	AS110M	AS110R	
Human IgG Fc-fragment	AS113D	AS113M	AS113R	
Protein A Donor Beads	AS102D	AS102M	AS102R	

AlphaScreen Acceptor Beads

Product Description	Part Number	Part Number	Part Number	Part Number
	1 mg	5 mg	25 mg	50 mg
Unconjugated	6762003	6762001		6762002
Protein A		6760137M	6760137R	

AlphaLISA Acceptor Beads

Product Description	Part Number	Part Number	Part Number	Part Number	Part Number
	250 µg	1 mg	5 mg	25 mg	50 mg
<i>Unconjugated</i>					
Unconjugated Acceptor Beads		6772001	6772002		6772003
<i>Tag capture</i>					
Streptavidin Acceptor Beads	AL125C		AL125M		AL125R
Strep-Tactin® Acceptor Beads	AL136C		AL136M		AL136R
Anti-6X His Acceptor Beads	AL128C		AL128M		AL128R
Anti-c-myc Acceptor Beads	AL111C		AL111M		AL111R
Anti-DIG Acceptor Beads	AL113C		AL113M		AL113R
Anti-DNP (dinitrophenol) Acceptor Beads	AL173C		AL173M	AL173R	
Anti-FITC Acceptor Beads	AL127C		AL127M		AL127R
Anti-FLAG Acceptor Beads	AL112C		AL112M		AL112R
Anti-GFP Acceptor Beads	AL133C		AL133M		AL133R
Anti-GST Acceptor Beads	AL110C		AL110M		AL110R
Anti-HA Acceptor Beads	AL170C		AL170M	AL170R	
Anti-HRP Acceptor Beads					
Anti-Maltose-Binding Protein (MBP) Acceptor Beads	AL134C		AL134M		AL134R
Anti-Protein C Tag Acceptor Beads	AL172C		AL172M	AL172R	
Anti-V5 Acceptor Beads	AL129C		AL129M		AL129R
GSH (glutathione) Acceptor Beads	AL109C		AL109M		AL109R
Nickel Chelate Acceptor Beads	AL108C		AL108M		AL108R
LCA Acceptor Beads	AL140C		AL140M		AL140R
<i>Antibody capture</i>					
Anti-Bovine IgA Acceptor Beads	AL169C		AL169M	AL169R	
Anti-Bovine IgG Acceptor Beads	AL165C		AL165M	AL165R	
Anti-Bovine IgG1 Acceptor Beads	AL166C		AL166M	AL166R	
Anti-Bovine IgG2 Acceptor Beads	AL167C		AL167M	AL167R	
Anti-Bovine IgM Acceptor Beads	AL168C		AL168M	AL168R	
Anti-Chicken IgY Acceptor Beads	AL131C		AL131M	AL131R	
Anti-Goat (Fc-specific) Acceptor Beads	AL107C		AL107M	AL107R	
Anti-Human (Fc-specific) Acceptor Beads	AL103C		AL103M	AL103R	
Anti-Human IgE	AL176C		AL176M	AL176R	
Anti-Human IgG1 Acceptor Beads	AL141C		AL141M	AL141R	
Anti-Human IgG2 Acceptor Beads	AL154C		AL154M	AL154R	
Anti-Human IgG4 Acceptor Beads	AL142C		AL142M	AL142R	
Anti-Mouse IgE Acceptor Beads	AL161C		AL161M		AL161R
Anti-Mouse (Fc-specific) IgG Acceptor Beads	AL105C		AL105M	AL105R	
Anti-Mouse IgG Acceptor Beads	AL164C		AL164M	AL164R	
Anti-Mouse IgG1 Acceptor Beads	AL157C		AL157M		AL157R
Anti-Mouse IgG2a Acceptor Beads	AL158C		AL158M		AL158R
Anti-Mouse IgG2b Acceptor Beads	AL159C		AL159M		AL159R
Anti-Mouse IgG3 Acceptor Beads	AL160C		AL160M		AL160R
Anti-Mouse IgM Acceptor Beads	AL130C		AL130M	AL130R	
Anti-Mouse IgM Acceptor Beads	AL162C		AL162M		AL162R
Anti-Rabbit IgG Acceptor Beads	AL104C		AL104M	AL104R	
Anti-Rat IgG Acceptor Beads	AL106C		AL106M	AL106R	
Anti-Sheep IgG Acceptor Beads	AL132C		AL132M	AL132R	
Human IgG Fc-fragment Acceptor Beads	AL174C		AL174M	AL174R	
Human IgG (whole)	AL175C		AL175M	AL175R	
Protein A Acceptor Beads	AL101C		AL101M	AL101R	
Protein G Acceptor Beads	AL102C		AL102M	AL102R	
Protein L Acceptor Beads	AL126C		AL126M	AL126R	

AlphaPlex Acceptor Beads

Product Description	Part Number	Part Number	Part Number	Part Number
	250 µg	1 mg	5 mg	25 mg
<i>Unconjugated</i>				
Unconjugated AlphaPlex 545 (Tb) Acceptor Beads		6782001	6782002	6782003
Unconjugated AlphaPlex 645 (Sm) Acceptor Beads		6792001	6792002	6792003
<i>Tag capture</i>				
Streptavidin AlphaPlex 545 (Tb) Acceptor Beads	AP125TB-C		AP125TB-M	AP125TB-R
Streptavidin AlphaPlex 645 (Sm) Acceptor Beads	AP125SM-C		AP125SM-M	AP125SM-R
Anti-6X His AlphaPlex 545 (Tb) Acceptor Beads	AP128TB-C		AP128TB-M	AP128TB-R
Anti-6X His AlphaPlex 645 (Sm) Acceptor Beads	AP128SM-C		AP128SM-M	AP128SM-R
Anti-FLAG AlphaPlex 545 (Tb) Acceptor Beads	AP112TB-C		AP112TB-M	AP112TB-R
Anti-FLAG AlphaPlex 645 (Sm) Acceptor Beads	AP112SM-C		AP112SM-M	AP112SM-R
Anti-HA AlphaPlex 545 (Tb) Acceptor Beads	AP170TB-C		AP170TB-M	AP170TB-R
GSH (glutathione) AlphaPlex 545 (Tb) Acceptor Beads	AP109TB-C		AP109TB-M	AP109TB-R
<i>Antibody capture</i>				
Anti-Human IgG AlphaPlex 545 (Tb) Acceptor Beads	AP103TB-C		AP103TB-M	AP103TB-R
Anti-Human IgG AlphaPlex 645 (Sm) Acceptor Beads	AP103SM-C		AP103SM-M	AP103SM-R
Anti-Mouse IgG AlphaPlex 545 (Tb) Acceptor Beads	AP105TB-C		AP105TB-M	AP105TB-R
Anti-Mouse IgG AlphaPlex 645 (Sm) Acceptor Beads	AP105SM-C		AP105SM-M	AP105SM-R
Anti-Rabbit IgG AlphaPlex 545 (Tb) Acceptor Beads	AP104TB-C		AP104TB-M	AP104TB-R
Protein A AlphaPlex 545 (Tb) Acceptor Beads	AP101TB-C		AP101TB-M	AP101TB-R
Protein A AlphaPlex 645 (Sm) Acceptor Beads	AP101SM-C		AP101SM-M	AP101SM-R

AlphaScreen Kits and Kit Components

Product Description	Part Number	Part Number	Part Number
	500 data points	10,000 data points	50,000 data points
<i>Unconjugated</i>			
AlphaScreen Conjugation Kit	676000K		
<i>Tag capture</i>			
Anti-c-myc Assay Kit	6760611C	6760611M	6760611R
Anti-DIG Assay Kit	6760604C	6760604M	6760604R
Anti-FITC Assay Kit	6760605C	6760605M	6760605R
Anti-FLAG Assay Kit	6760613C	6760613M	6760613R
Anti-GST Assay Kit	6760603C	6760603M	6760603R
Anti-HA Assay Kit	6760612C	6760612M	6760612R
Nickel Chelate Assay Kit	6760619C	6760619M	6760619R
Biotinylated HIS Tag		6760303M	
Biotinylated GST		6760305M	
<i>Antibody capture</i>			
Anti-mouse IgG Assay Kit	6760606C	6760606M	6760606R
Anti-rabbit IgG Assay Kit	6760607C	6760607M	6760607R
Protein A Assay Kit	6760617C	6760617M	6760617R

Alpha Buffers

Product Description	Part Number	Part Number
	10 mL	100 mL
Bovine Immunoassay Buffer	AL010C	AL010F
Casein Buffer, 5X	AL014C	AL014F
Dissociation Buffer, 3X	AL006C	AL006F
Epigenetics Buffer, 5X	AL008C	AL008F
HiBlock Buffer, 10X	AL004C	AL004F
High Salt (NaCl) Buffer, 5X	AL007C	AL007F
Immunoassay Buffer, 10X	AL000C	AL000F
Lysis Buffer, 5X	AL003C	AL003F
MES Buffer, 5X	AL016C	AL016F
PPI Buffer, 5X	AL015C	AL015F
Universal Assay Buffer, 5X	AL001C	AL001F

Control Beads: TruHits

Product Description	Part Number	Part Number	Part Number	Part Number
	500 data points	1,000 data points	5,000 data points	10,000 data points
AlphaScreen TruHits		6760627D		6760627M
AlphaLISA TruHits		AL900D		AL900M
AlphaPlex TruHits Sm	AP900SM-D		AP900SM-M	
AlphaPlex TruHits Tb	AP900TB-D		AP900TB-M	

Alpha Control Kits

Product Description	Part Number	Part Number	Part Number
	1,000 data points	10,000 data points	50,000 data points
AlphaScreen Omnibeads	6760626D	6760626M	6760626R

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



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2011-2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

H (009841G_01) PKI April 2019