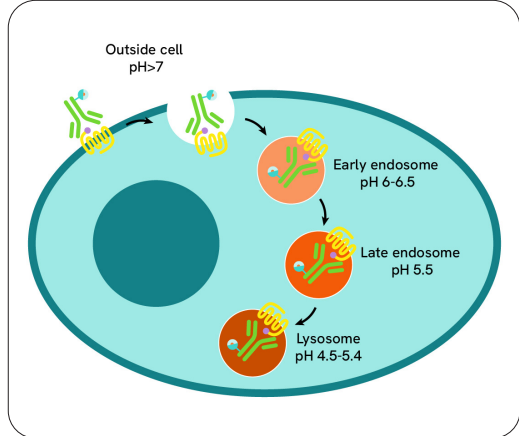


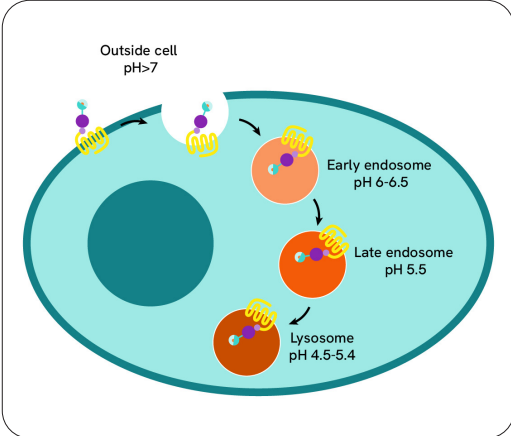


Obesity, Diabetes and Metabolic Disease Reagents

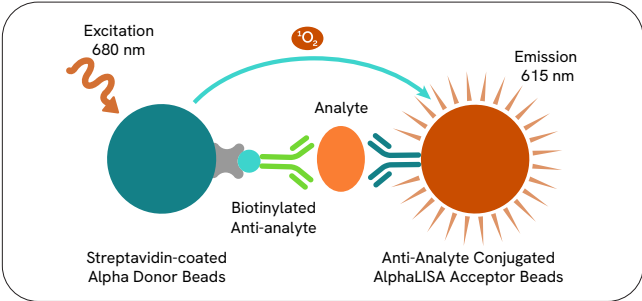
Revvity has supported the research and development of therapies for metabolic disorders with 360° research solutions. We offer an extensive portfolio of GPCRs, cell signaling, cytokines and biomarkers research reagents and assays for obesity, diabetes, MASH, and MASLD, which can be used to investigate the mechanisms at play in these critical diseases.



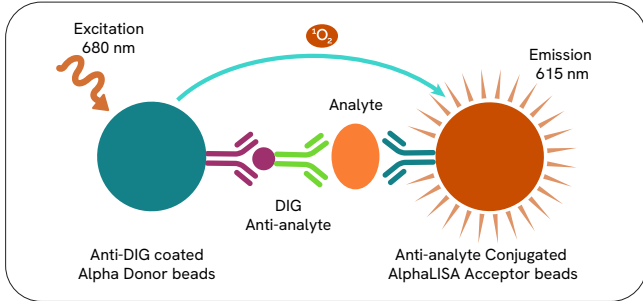
| pHSense™ Eu Anti-FLAG



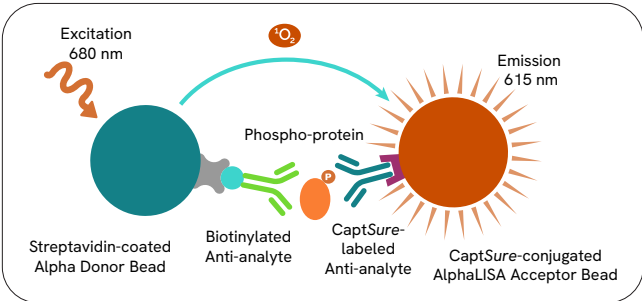
| pHSense™ Eu SNAP



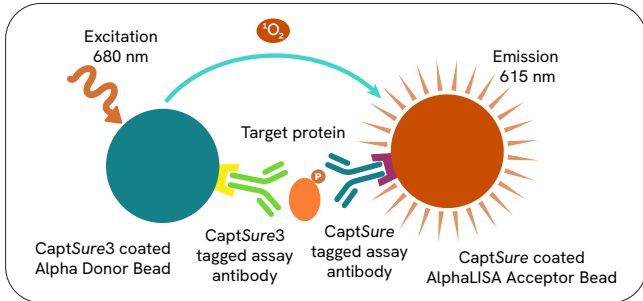
| AlphaLISA™



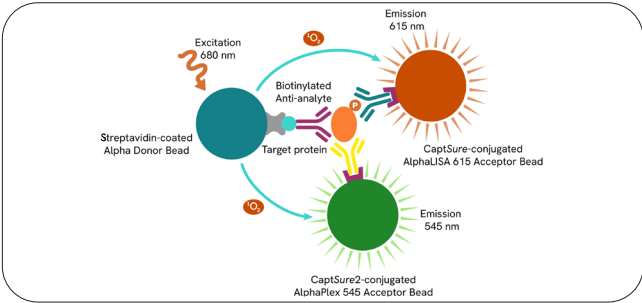
| AlphaLISA™ Biotin-Free



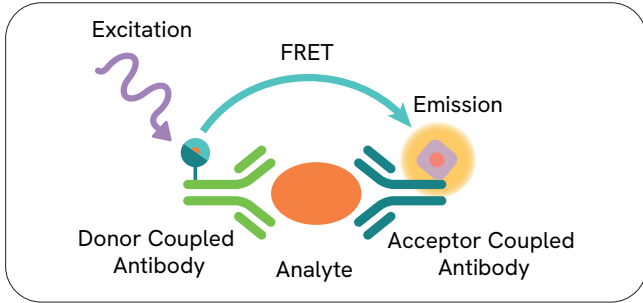
| AlphaLISA™ SureFire® Ultra™



| AlphaLISA SureFire® Biotin-Free



| Multiplex SureFire® Ultra™



| HTRF™

Obesity

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
β-Arrestin 1	Total	64BAR1TPEB						
β-Arrestin 2	Recruitment	62BDBAR2PEB						
	Total	64BAR2TPEB						
β-Catenin	Phospho-T41/S33/37	64CATPEG	Phospho-S37	ALSU-PBCAT-A500				
			Phospho-S675	ALSU-PBCAT-C500				
	Total	64NCAPEG	Phospho-T41/S45	ALSU-PBCAT-B500				
			Total	ALSU-TBCAT-A500				
AKT1/2/3	Phospho-S473	64AKSPEG	Phospho-S473	ALSU-PAKT-B500	Phospho-S473 & AKT1 Total	MPSU-PTAKT-M500		
				ASBF-PAKT-A500*				
	Phospho-T308	64AKTPEG	Phospho-T308	ALSU-PAKT-A500	Phospho-T308 & Total	MPSU-PTAKT-O500		
				Phospho-T450				
	Total	64NKTPEG	Total	ALSU-TAKT-B500	Phospho-T450 & Total	MPSU-PTAKT-N500		
				ASBF-TAKT-A500*				
AKT3	Phospho-S473	63ADK082PEG						
	Total	63ADK083PEG						
AMPK	Phospho-T172	64MPKPEG						
	Total	63ADK060PEG						
cAMP	Gi	62AM9PEB						
	Gs Dynamic	62AM4PEB						
	Gs HiRange	62AM6PEB						
c-RAF	Phospho-S338	63ADK018PEG						
	Phospho-S43	64CRAFS43PEG						
		64CRAFS43V2PEG						
	Total	64CRAFTPEG						
FOXO1	Phospho-S256	64FOXPEG						
	Total	64NFOPEG						

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
GSK3α	Phospho-S21	64GPAPEG						
	Total	64GTAPEG						
GSK3β	Phospho-S9	64GPBPEG	Phospho-S9	ALSU-PGS3B-A500				
	Total	64GTBPEG	Total	ALSU-TGS3B-A500				
GTP / Gi protein		62GTPPEG						
Insulin	High range	62IN1PEG					Human and Mouse	AL3184C
	Ultra sensitive	62IN2PEG						
	Mouse Serum	62IN3PEB						
IP-One	Gq	62IPAPEB						AL3145D
IR-β (Insulin Receptor Beta)	Phospho-Y1150/1151	63ADK016PEG	Phospho-Y1150/1151	ALSU-PINR-B500	Phospho-T1150/1151 & Total	MPSU-PTINR-K500		
	Total	63ADK019PEG	Total	ALSU-TINR-A500				
IRS1	Phospho-S307/S312	64IRSPEG	Phospho-S612	ALSU-PIRS1-A500				
	Total	64NRSPEG	Total	ALSU-TIRS1-A500				
JNK	Phospho-T183/Y185	64JNKPEG						
mTOR	Phospho-S2448	64TORPEG	Phospho-S2448	ALSU-PMTOR-C500				
			Total	ALSU-TMTOR-B500				

Diabetes

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
β-Arrestin 1	Total	64BAR1TPEB						
β-Arrestin 2	Recruitment	62BDBAR2PEB						
	Total	64BAR2TPEB						
4EBP1	Phospho-T37/46	64BP1PEG	Phospho-T37/46	ALSU-P4EBP-A500				
			Total	ALSU-T4EBP-A500				
AKT1	Phospho-S473	63ADK078PEG	Phospho-S129	ALSU-PAKT1-D500				
			Phospho-S473	ALSU-PAKT1-C500				
	Total	63ADK079PEG	Total	ALSU-TAKT1-B500				

* Biotin Free kit

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
AKT2	Phospho-S473	63ADK080PEG						
	Total	63ADK081PEG						
AKT3	Phospho-S473	63ADK082PEG						
	Total	63ADK083PEG						
BAD	Phospho-S112	64BADPEG	Phospho-S112	ALSU-PBAD-A500				
			Phospho-S136	ALSU-PBAD-B500				
			Total	ALSU-TBAD-A500				
BCL-2			Phospho-S70	ALSU-PBCL2-A500			Human	AL384C
			Total	ALSU-TBCL2-A500				
BCL-XL			Total	ALSU-TBCLXL-A500				
cAMP	Gi	62AM9PEB						
	Gs Dynamic	62AM4PEB						
	Gs HiRange	62AM6PEB						
CCL2	Human	62HCCL2PEG					Human	AL244C
	Mouse	62MCCL2PEG					Mouse & Rat	AL509C
CCL3	Human	62HCCL3PEG					Human	AL257C
CREB	Phospho-S133	64CREPEG	Phospho-S133	ALSU-PCREB-A500				
	Total	63ADK052PEG	Total	ALSU-TCREB-A500				
c-Src	Phospho-Y419	64SRCPEG						
	Total	64NSRPEG						
DNA PKcs			Phospho-S2056	ALSU-PDNA-A500				
			Total	ALSU-TDNA-A500				
EGFR	Phospho-Pan-Y	64HR1PEG	Phospho-Y1068	ALSU-PEGFR-B500	Phospho-Y1068 & Total	MPSU-PTEGFR-K500	Total	AL340C
	Phospho-Y1068	64EG1PEG						
	Total	64NG1PEG	Total	ALSU-TEGFR-A500				
EGFR DEL19	Total	64EGFRD19PEG						
EGFR L858R	Total	64EGFRL8PEG	Total	ALSU-TEGFR-B500				
eIF2α	Phospho-S52	64EF2PEG	Phospho-S51	ALSU-PEIF2-B500	Phospho-S51 & Total	MPSU-PTEIF2-K500		
	Total	64NEFPEG	Total	ALSU-TEIF2-A500				

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
eIF4E	Phospho-S209	64EF4PEG	Phospho-S209	ALSU-PEIF4-A500				
			Total	ALSU-TEIF4-A500				
ERK	Phospho-T202/TY204	64ERKPEG						
	Phospho-T202/Y204 Advanced	64AERPEG						
	Total	64NRKPEG						
ERK1/2			Phospho-T202/Y204	ALSU-PERK-A500	Phospho-T202/Y204 & Total	MPSU-PTERK-M500		
				ASBF-PERK-A500*				
			Total	ALSU-TERK-A500				
				ASBF-TERK-A500*				
FOXO1	Phospho-S256	64FOXPEG						
	Total	64NFOPEG						
GLP1	Active	62GLPPEG					Human 7-36 amide	AL359C
Glucagon		62CGLPEG					Human	AL3140C
		62SGLPEB						
GSK3α	Phospho-S21	64GPAPEG						
	Total	64GTAPEG						
GSK3β	Phospho-S9	64GPBPEG	Phospho-S9	ALSU-PGS3B-A500				
	Total	64GTBPEG	Total	ALSU-TGS3B-A500				
IκKβ	Phospho-S177/181	64KKBPEG	Phospho-S177/181	ALSU-PIKKB-A500				
	Total	63ADK097PEG	Total	ALSU-TIKKB-A500				

* Biotin-free

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
IFN-γ	Human	62HIFNGPEG					Bovine	AL535C
							Cynomolgus	AL561C
							Human	AL217C
								AL327C*
							Human HP	AL3153C
	AL3154C*							
	Mouse	62MIFNGPEG					Mouse	AL501C
							Mouse HP	AL593C
							Ovine	AL560C
							Porcine	AL549C
Rat			AL3195C					
IGF1 Receptor β			Phospho-Y1135/1136	ALSU-PIGFR-B500	Phospho-Y1135/1136 & Total	MPSU-PTIGFR-K500		
			Total	ALSU-TIGFR-A500				
IL-1β	Human	62HIL1BPEG					Bovine	AL539C
							Human	AL220C
	Human HP	62HIL1B2PEG					Human HP	AL3160C
							Mouse & Rat HP	AL592C
	Mouse	62MIL1BPEG					Mouse	AL503C
							Porcine	AL575C
							Rat	AL518C
IL-6	Human	62HIL06PEG					Bovine	AL538C
							Cynomolgus	AL555C
	Human HP	62HIL6V2PEG					Human	AL223C
								AL3025C*
							Human HP	AL3202C
	Mouse	62MIL06PEG						AL3203C*
							Mouse	AL504C
							Porcine	AL573C

* Biotin-free
HP = High-performance

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
IL-8	Human	62HIL08PEG					Bovine & Ovine	AL545C
							Human	AL224C
								AL328C*
							Human HP	AL3164C
								AL3165C*
							Porcine	AL571C
Insulin	High range	62IN1PEG					Human and Mouse	AL3184C
	Ultra sensitive	62IN2PEG						
	Mouse Serum	62IN3PEB						
IP-One	Gq	62IPAPEB						AL3145D
IR-β (Insulin Receptor Beta)	Phospho-Y1150/1151	63ADK016PEG	Phospho-Y1150/1151	ALSU-PINR-B500	Phospho-T1150/1151 & Total	MPSU-PTINR-K500		
	Total	63ADK019PEG	Total	ALSU-TINR-A500				
IRS1	Phospho-mouse S307/ human S312	64IRSPEG	Phospho-S612	ALSU-PIRS1-A500				
	Total	64NRSPEG	Total	ALSU-TIRS1-A500				
IRS2	Total	64IRTPEG						
JNK	Phospho-T183/Y185	64JNKPEG						
JNK1/2			Total	ALSU-TJNK-A500				
JNK1/3			Phospho-T183/Y185	ALSU-PJNK-A500				
MEK1	Phospho-S218/222	64ME1PEG	Phospho-S218/222	ALSU-PMEK1-A500	Human and Mouse Phospho-S218/222 & Total	MPSU-PTMEK1-K500		
	Phospho-S298	64MEK1S98PEG						
	Total	64NE1PEG	Total	ALSU-TMEK1-A500				
MEK1/2	Phospho-S218/222	64ME2PEG						
MEK2			Phospho-S217/221	ALSU-PMEK2-A500				
			Total	ALSU-TMEK2-A500				
mTOR	Phospho-S2448	64TORPEG	Phospho-S2448	ALSU-PMTOR-C500				
			Total	ALSU-TMTOR-B500				
NF-kB	Phospho-S536	64NFBPEG			Phospho-S536 & Total	MPSU-PTNFKB-K500		
	Total	64NFTPEG						
NF-kB p65			Phospho-S536	ALSU-PNFKB-A500				
			Total	ALSU-TNFKB-A500				

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
NF-kB2			Total	ALSU-TNFKB2-A500				
P38 MAPK	Phospho-T180/Y182	64P38PEG	Phospho-T180/Y182	ALSU-PP38-B500				
			Total	ALSU-TP38-B500				
p70 S6K	Phospho-T389	64S6KPEG	Phospho-T389	ALSU-PP70-A500	Phospho-T389 & total	MPSU-PTP70-K500		
			Total	ALSU-TP70-A500				
PERK			Phospho-T980	ALSU-PPERK-A500				
			Phospho-T982	ALSU-PPERK-B500				
			Total	ALSU-TPERK-A500				
PI3K p85			Phospho-Y458	ALSU-PP13K-A500				
			Phospho-Y467	ALSU-PPI3K-B500				
			Human Total	ALSU-TPI3K-B500				
			Human and Mouse Total	ALSU-TP13K-A500				
PKC-θ			Phospho-T538	ALSU-PPKC-A500				
PLC-γ	Phospho-Y783	64PLCG1Y3PEG	Phospho-S1248	ALSU-PPLCG1-B500				
			Phospho-Y1253	ALSU-PPLCG1-C500				
	Total	64PLCG1TPEG	Phospho-Y783	ALSU-PPLCG1-A500				
			Total	ALSU-TPLCG1-A500				
PLC-γ2	Phospho-Y1217	64PLCG2Y7PEG	Phospho-Y1197	ALSU-PPLCG2-C500				
			Phospho-Y1217	ALSU-PPLCG2-D500				
	Total	64PLCG2TPEG	Phospho-Y753	ALSU-PPLCG2-A500				
			Phospho-Y759	ALSU-PPLCG2-B500				
			Total	ALSU-TPLCG2-A500				
Raf1			Phospho-S259	ALSU-PRAF-B500				
			Phospho-S338	ALSU-PRAF-A500				
			Total	ALSU-TRAF-A500				
S6RP	Phospho-S235/236	64RP6PEG						
SRC			Phospho-Y419	ALSU-PSRC-A500				

* Biotin-free
HP = High-performance

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
STAT1	Phospho-Y701	63ADK026PEG	Phospho-S727	ALSU-PST1-B500	Phospho-S727 & Total	MPSU-PTST1-M500		
			Phospho-Y701 HP	ALSU-PST1-C500				
			Phospho-Tyr701 Biotin-free	ASBF-PST1-A500				
	Total	63ADK096PEG	Phospho-Y701	ALSU-PST1-A500	Phospho-Y701 & Total	MPSU-PTST1-K500		
			Total	ALSU-TST1-A500				
				ASBF-TST1-A500*				
TAK1			Phospho-T184/187	ALSU-PTAK-A500				
			Total	ALSU-TTAK-A500				
TNF-α	Cynomolgus	62CTNFAPEG					Bovine	AL534C
							Cynomolgus	AL562C
	Human HP	62HTNFAV2PEG					Human	AL208C
								AL325C*
	Human	62HTNFAPEG					Human HP	AL3157C
								AL3158C*
	Mouse	62MTNFAPEG					Mouse	AL505C
			AL541C*					
			Porcine	AL548C				
TNF-α/ TNFR1		64BDTNFPEG						
TNF-α/ TNFR2							Human	AL3121C
VEGF Receptor 2 (VEGFR2)	Phospho-Y1175	63ADK041PEG	Phospho-Y1175	ALSU-PVGFR-A500				
			Phospho-Y1214	ALSU-PVGFR-C500				
			Phospho-Y951	ALSU-PVGFR-B500				
			Total	ALSU-TVGFR-A500				

* Biotin-free
HP = High-performance

MASH/MASLD

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
A-SMA	Human	62ASMAPEG						
ACC	Phospho-S79	64ACCPEG	Phospho-S79	ALSU-PACC-A500				
	Total	64NACPEG	Total	ALSU-TACC-A500				
AKT1/2/3	Phospho-S473	64AKSPEG	Phospho-S473	ALSU-PAKT-B500 ASBF-PAKT-A500*	Phospho-S473 & AKT1 & Total	MPSU-PTAKT-M500		
	Phospho-T308	64AKTPEG	Phospho-T308 Phospho-T450	ALSU-PAKT-A500 ALSU-PAKT-C500		MPSU-PTAKT-O500		
	Total	64NKTPEG	Total	ALSU-TAKT-B500 ASBF-TAKT-A500*	Phospho-T450 & Total	MPSU-PTAKT-N500		
AKT1	Phospho-S473	63ADK078PEG	Phospho-S129 Phospho-S473	ALSU-PAKT1-D500 ALSU-PAKT1-C500				
	Total	63ADK079PEG	Total	ALSU-TAKT1-B500				
AKT2	Phospho-S473	63ADK080PEG						
	Total	63ADK081PEG						
AKT3	Phospho-S473	63ADK082PEG						
	Total	63ADK083PEG						
AMPK	Phospho-T172	64MPKPEG						
	Total	63ADK060PEG						
AMPKα			Phospho-T172	ALSU-PAMPK-A500				
AMPKα1/2			Total	ALSU-TAMPK-A500	Phospho-T172 & Total	MPSU-PTAMPK-K500		
ApoB	Human	64APOBPEG						
ATF2	Phospho-T71	63ADK015PEG	Phospho-T71	ALSU-PATF2-A500				
			Total	ALSU-TATF2-A500				
ATF4	Total	64ATF4TPEG	Total	ALSU-TATF4-A500				
BAD	Phospho-S112	64BADPEG	Human Phospho-S112	ALSU-PBAD-A500				
			Human Phospho-S136	ALSU-PBAD-B500				
			Human Total	ALSU-TBAD-A500				

* Biotin-free

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
cAMP	Gi	62AM9PEB						
	Gs Dynamic	62AM4PEB						
	Gs HiRange	62AM6PEB						
CCL2	Human	62HCCL2PEG					Human	AL244C
	Mouse	62MCCL2PEG					Mouse & Rat	AL509C
CCL5	Human	62HCCL5PEG					Human	AL289C
							Mouse	AL589C
c-Jun	Phospho-S63	64JUNPEG	Phospho-S63	ALSU-PCJUN-A500				
			Phospho-S73	ALSU-PCJUN-B500				
			Total	ALSU-TCJUN-A500				
Cleaved PARP1			Total	ALSU-CPARP1-A500				
eIF2α	Phospho-S52	64EF2PEG	Phospho-S51	ALSU-PEIF2-B500	Phospho-S51 & Total	MPSU-PTEIF2-K500		
	Total	64NEFPEG	Total	ALSU-TEIF2-A500				
ERK	Phospho-T202/Y204	64ERKPEG						
	Phospho-T202/y204 Advanced	64AERPEG						
	Total	64NRKPEG						
ERK1			Phospho-T202/Y204	ALSU-PERK-A500				
				ASBF-PERK-A500*				
			Total	ALSU-TERK-A500 ASBF-TERK-A500*				
ERK1/2			Phospho-T202/Y204	ALSU-PERK-A500 ASBF-PERK-A500*	Phospho-T202/Y204 & Total	MPSU-PTERK-M500		
				ALSU-TERK-A500 ASBF-TERK-A500*				
			Total					
FOXO1	Phospho-S256	64FOXPEG						
	Total	64NFOPEG						
GSK3β	Phospho-S9	64GPBPEG	Phospho-S9	ALSU-PGS3B-A500				
	Total	64GTBPEG	Total	ALSU-TGS3B-A500				
Hyaluronic acid		63ADK089PEG					Human	AL354C

* Biotin-free

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™		
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number	
IFN-α	Human	62HIFNAPEG					Human	AL297C	
							Mouse	AL550C	
							Porcine	AL576C	
IFN-β	Human	62HIFNBPEG					Human HP	AL3191C	
	Mouse	62MIFNBPEG					Mouse	AL586C	
								Porcine	AL577C
IL-1β	Human	62HIL1BPEG					Bovine	AL539C	
	Human HP	62HIL1B2PEG					Human	AL220C	
								Human HP	AL3160C
								Mouse & Rat HP	AL592C
	Mouse	62MIL1BPEG					Mouse	AL503C	
								Porcine	AL575C
								Rat	AL518C
IL-12	Human	62HIL12PEG							
IL-18	Human	62HIL18PEG					Bovine	AL558C	
							Human	AL3137C	
							Mouse	AL3189C	
							Porcine	AL559C	
IL-6	Human	62HIL06PEG					Bovine	AL538C	
	Human HP	62HIL6V2PEG					Cynomolgus	AL555C	
								Human HP	AL3202C
									AL3203C*
	Mouse	62MIL06PEG					Human	AL223C	
									AL3025C*
								Mouse	AL504C
						Porcine	AL573C		

* Biotin-free
HP = High-performance

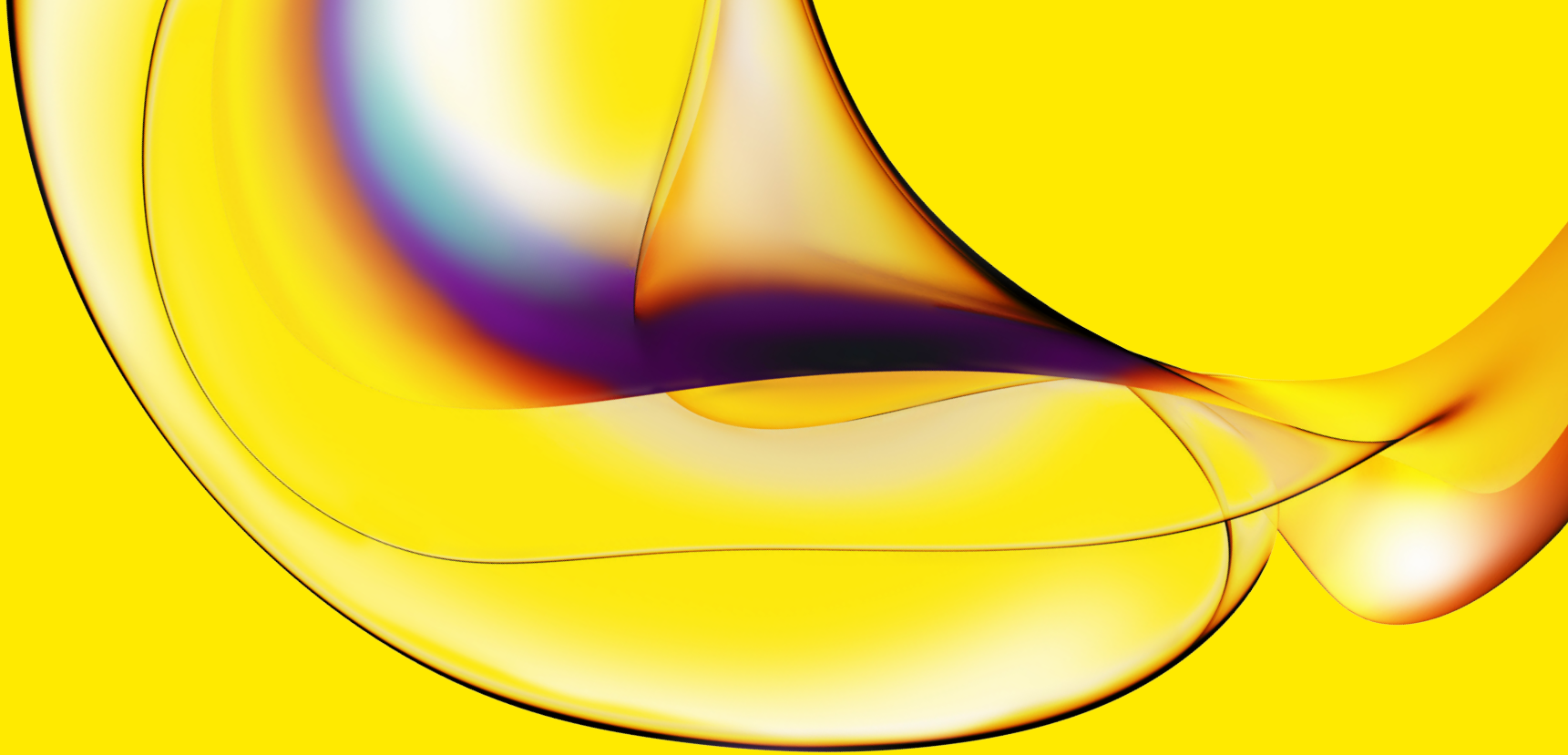
Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
IL-8	Human	62HIL08PEG					Bovine & Ovine	AL545C
							Human	AL224C
								AL328C*
							Human HP & Biotin-free	AL3165C
Insulin	High range	62IN1PEG					Human and Mouse	AL3184C
	Ultra sensitive	62IN2PEG						
	Mouse Serum	62IN3PEB						
IP-One	Gq	62IPAPEB						AL3145D
IR-β (Insulin Receptor Beta)	Phospho-Y1150/1151	63ADK016PEG	Phospho-Y1150/1151	ALSU-PINR-B500	Phospho-T1150/1151 & Total	MPSU-PTINR-K500		
	Total	63ADK019PEG	Total	ALSU-TINR-A500				
IRF3	Phospho-S385/386	6FRF3PEG	Phospho-S386	ALSU-PIRF3-A500				
			Total	ALSU-TIRF3-A500				
IRS1	Phospho-mouse S307/ human S312	64IRSPEG	Phospho-S612	ALSU-PIRS1-A500				
	Total	64NRSPEG	Total	ALSU-TIRS1-A500				
IRS2	Total	64IRTPEG						
JNK	Phospho-T183/Y185	64JNKPEG						
JNK1/2			Total	ALSU-TJNK-A500				
JNK1/3			Phospho-T183/Y185	ALSU-PJNK-A500				
MEK1	Phospho-S218/222	64ME1PEG	Phospho-S218/222	ALSU-PMEK1-A500	Phospho-S218/222 & Total	MPSU-PTMEK1-K500		
	Phospho-S298	64MEK1S98PEG						
	Total	64NE1PEG	Total	ALSU-TMEK1-A500				
MEK1/2	Phospho-S218/222	64ME2PEG						
MEK2			Phospho-S217/221	ALSU-PMEK2-A500				
			Total	ALSU-TMEK2-A500				
MMP2	Human	62MM2PEG						
MMP9	Human	62MM9PEG					Human	AL3138C
							Mouse	AL519C

* Biotin-free
HP = High-performance

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
mTOR	Phospho-S2448	64TORPEG	Phospho-S2448	ALSU-PMTOR-C500				
			Total	ALSU-TMTOR-B500				
Myd88	Total	64MYD88PEG	Total	ALSU-TMYD-A500				
NF-κB	Phospho-S536	64NFBPEG			Phospho-S536 & Total	MPSU-PTNFKB-K500		
	Total	64NFTPEG						
NF-κB p65			Phospho-S536	ALSU-PNFKB-A500				
			Total	ALSU-TNFKB-A500				
NF-κB2			Total	ALSU-TNFKB2-A500				
P38 MAPK	Human and Mouse Phospho-T180/Y182	64P38PEG	Phospho-T180/Y182	ALSU-PP38-B500				
			Total	ALSU-TP38-B500				
PARP	Human and Mouse cleaved-D214	64PARPEG						
PDGF Receptor (PDGFR)			Total	ALSU-TPDGF-A500				
PDGF-BB							Human	AL3049C
PDGF-BB / PDGFR-α							Human	AL3064C
PDGF-BB / PDGFR-β							Human	AL3065C
PDGFR-β (PDGF Receptor Beta)	Phospho-Y751	64PDGPEG	Phospho-Y751	ALSU-PPDGF-A500	Phospho-Y751 & Total	MPSU-PTPDGF-K500		
PERK			Mouse Phospho-T980	ALSU-PPERK-A500				
			Human Phospho-T982	ALSU-PPERK-B500				
			Human and Mouse Total	ALSU-TPERK-A500				
PI3K p85			Mouse Phospho-Y458	ALSU-PP13K-A500				
			Human Phospho-Y467	ALSU-PPI3K-B500				
			Human Total	ALSU-TPI3K-B500				
			Human and Mouse Total	ALSU-TP13K-A500				
PICP (Pro-Collagen 1 C-Terminal Pro-Peptide)							Human	AL3208C
PKC-θ			Phospho-T538	ALSU-PPKC-A500				
Pro-Collagen Type 1	Human	63ADK014PEG						

Analyte	HTRF™		AlphaLISA™ Surefire® Ultra™		AlphaLISA™ Surefire® Ultra™ Multiplex		AlphaLISA™	
	Variant	Part Number	Variant	Part Number	Variant	Part Number	Variant	Part Number
Raf1			Phospho-S259	ALSU-PRAF-B500				
			Phospho-S338	ALSU-PRAF-A500				
			Total	ALSU-TRAF-A500				
SMAD2	Phospho-S465/467	64SMAD2S5PEG	Phospho-S465/467	ALSU-PSM2-A500	Phospho-S465/467	MPSU-PTSM2-K500		
	Total	64SMAD2TPEG	Total	ALSU-TSM2-A500	Total			
SMAD3	Phospho-S423/425	63ADK025PEG	Phospho-S423/425	ALSU-PSM3-A500				
	Total	64ND3PEG	Total	ALSU-TSM3-A500				
SMAD4	Phospho-T277	64SMAD4T7PEG	Total	ALSU-TSM4-A500				
	Total	64SMAD4TPEG						
TAK1			Phospho-T184/187	ALSU-PTAK-A500				
			Total	ALSU-TTAK-A500				
TBK1	Phospho-S172	64TBKPEG	Phospho-S172	ALSU-PTBK1-A500				
	Total	64NTBPEG	Total	ALSU-TTBK1-A500				
TGF-β1	Human and Mouse	62HTGFBPEG					Human	AL336C
								AL361C*
TIMP1	Human	64TIMPEG						AL272C
TNF-α	Cynomolgus	62CTNFAPEG					Bovine	AL534C
							Cynomolgus	AL562C
	Human	62HTNFAPEG					Human	AL208C
								AL325C*
	Mouse	62MTNFAPEG					Human HP	AL3157C
							Mouse	AL541C*
		Porcine	AL548C					
YAP	Phospho-S127	64YAPPEG	Phospho-S127	ALSU-PYAP-B500				
	Total	64YATPEG	Total	ALSU-TYAP-A500				

* Biotin-free
HP = High-performance



www.revivity.com

revvity

Copyright ©2025, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

1906451

For research use only. Not for use in diagnostic procedures.