

SPECIES COMPATIBILITY

JULY 2022

Cell Signaling: Phospho- & Total-protein assays

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

	HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™ 4EBP1, Phospho-T37/46	●	●	●	○	○	○	○	○
HTRF® 4EBP1, Phospho-T37/46	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™ 4EBP1, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ Alpha-Synuclein Aggregate Specific	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ Alpha-Synuclein, Phospho-S129	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ Alpha-Synuclein, Total	●	○	○	○	○	○	○	○
HTRF® Alpha-SMA	●	●	●	○	○	○	○	○
HTRF® Alpha-Synuclein-S129	●	●	○	○	○	○	○	○
HTRF® Alpha-Synuclein, Total	●	✗	○	○	○	○	○	○
HTRF® Alpha-Tubulin Housekeeping	●	●	●	●	●	○	●	●
AlphaLISA® SureFire® Ultra™ ACC, Phospho-S79	●	✗	✗	○	○	○	○	○
HTRF® ACC, Phospho-S79	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ ACC, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ AKT1/2/3, Phospho-S473	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	AKT1/2/3, Phospho-S473	●	●	●	●	●	●	○	○
AlphaLISA® SureFire® Ultra™	AKT1/2/3, Phospho-T308	●	●	●	○	○	○	○	○
HTRF®	AKT1/2/3, Phospho-T308	●	●	●	●	●	●	○	○
AlphaLISA® SureFire® Ultra™	AKT 1/2/3, Phospho-T450	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	AKT1/2/3, Total	●	●	○	○	○	○	○	○
HTRF®	AKT 1/2/3total	●	●	●	○	●	○	○	○
Alpha SureFire® Ultra™ Multiplex	AKT 1/2/3, Phospho-T308 + AKT 1/2/3, Total	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	AKT 1/2/3, Phospho-T450 + AKT 1/2/3, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	AKT1, Phospho-S473	●	●	●	○	○	○	○	○
HTRF®	AKT1, Phospho-S473	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	AKT1, Total	●	●	●	○	○	○	○	○
HTRF®	AKT1, Total	●	●	●	●	○	○	○	○
HTRF®	AKT2, Phospho-S473	●	●	●	●	○	○	○	○
HTRF®	AKT2, Total	●	●	●	●	○	○	○	○
HTRF®	AKT3, Phospho-S473	●	●	●	○	○	○	○	○
HTRF®	AKT3, Total	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	ALK, Phospho-Y1604 + ALK, Total	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ALK, Phospho-Y1586	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ALK, Phospho-Y1604	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ALK, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	AMPKα1/2, Phospho-T172	●	●	●	○	○	○	○	○
HTRF®	AMPK, Phospho-T172	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	AMPKα1/2, Total	●	●	●	○	○	○	○	○
HTRF®	AMPK, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Androgen Receptor, Total NEW	●	●	○	○	○	○	○	○
HTRF®	Androgen receptor, Total	●	✗	✗	✗	✗	✗	✗	✗

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	AP2, Total	●	●	●	○	●	○	○	○
HTRF®	ATF2, Phospho-T71	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATG16L1, Phospho-S278	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATG16L1, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATM, Phospho-S1981	●	○	○	○	○	○	○	○
HTRF®	ATM, Phospho-S1981 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATM, Total	●	●	●	○	○	○	○	○
HTRF®	ATM, Total NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Aurora A, Phospho-T288 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Aurora A, Total NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Aurora B, Phospho-T232 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Aurora B, Total NEW	●	X	○	○	○	○	○	○
HTRF®	BAD, Phospho-S112	●	●	●	●	○	○	○	○
HTRF®	B-arrestin 1, Total	●	●	●	○	●	○	○	○
HTRF®	B-arrestin 2, Total	●	●	●	○	●	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin, Phospho-S37	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin, Phospho-T41/S45	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin, Phospho-S675	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Beclin-1, Total	●	●	●	○	○	○	○	○
HTRF®	Beta Catenin, Phospho-T41/S33/37	●	●	●	●	○	○	●	○
HTRF®	Beta Catenin, Total	●	●	●	●	○	○	●	○
HTRF®	BRD2, Total	●	○	○	○	○	○	○	○
HTRF®	BRD3, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	BRD4, Total	●	X	X	○	○	○	○	○
HTRF®	BRD4, Total	●	●	○	○	○	○	○	○
HTRF®	BRD7, Total	●	●	○	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	BRD9, Total	●	●	●	○	○	○	○	○
HTRF®	BRD9, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	BTK, Phospho-Y223	●	○	○	○	○	○	○	○
HTRF®	BTK, Phospho-Y223	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	BTK, Phospho-Y551	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	BTK, Total	●	○	○	○	○	○	○	○
HTRF®	BTK, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CaMKII, Phospho-T286	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CaMKII, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CD3 Zeta, Phospho-Y142	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CD3 Zeta, Phospho-Y83	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CD3 Zeta, Total	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK1, Phospho-T14	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK1, Phospho-T161	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK1, Phospho-Y15	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK1, Total	●	X	X	○	○	○	○	○
HTRF®	CDK1, Total	●	X	X	●	X	X	X	X
Alpha SureFire® Ultra™ Multiplex	CDK1, Phospho-T14 + Total CDK1 NEW	●	X	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	CDK1, Phospho-Y15 + Total CDK1 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK2, Phospho-T160 NEW	●	●	○	○	○	○	○	○
HTRF®	CDK2, Phospho-Y15	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK2, Total	●	○	○	○	○	○	○	○
HTRF®	CDK2, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CDK4, Total	●	○	○	○	○	○	○	○
HTRF®	CDK4, Total	●	●	○	○	○	○	○	○
HTRF®	CDK6, Total	●	X	X	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	CDK7, Total	●	✗	✗	○	○	○	○	○
HTRF®	CDK7, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	CDK9, Total	●	●	●	○	○	○	○	○
HTRF®	CDK9, Total	●	●	●	✗	✗	✗	✗	✗
HTRF®	CDK12, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	Cereblon, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK1, Phospho-S345	●	●	○	○	○	○	○	○
HTRF®	CHK1, Phospho-S345	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK1, Total	●	○	○	○	○	○	○	○
HTRF®	CHK1, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK2, Phospho-T68	●	○	○	○	○	○	○	○
HTRF®	CHK2, Phospho-T68	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK2, Total	●	○	○	○	○	○	○	○
HTRF®	CHK2, Total	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Jun, Phospho-S63	●	●	●	○	○	○	○	○
HTRF®	cJun, Phospho-S63	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Jun, Phospho-S73	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Jun, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Met, Phospho-Y1234/1235	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Myb, Phospho-S11	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Myb, Total	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Myc, Phospho-S62	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Myc, Phospho-T58	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Myc, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Cofilin, Phospho-S3	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Cofilin, Total	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	cRAF, Phospho-S338	●	●	●	○	○	○	○	○
HTRF®	cRaf, Phospho S43	●	○	○	○	○	○	○	○
HTRF®	cRaf, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CREB, Phospho-S133	●	●	●	○	○	○	○	○
HTRF®	CREB, Phospho-S133	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	CREB, Total	●	●	●	○	○	○	○	○
HTRF®	CREB, Total	●	●	●	●	○	●	○	○
AlphaLISA® SureFire® Ultra™	Cyclin D1, Phospho-T286	●	X	X	○	○	○	○	○
HTRF®	Cyclin D1, Phospho-T286	●	X	○	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	Cyclin D1, Total	●	X	X	○	○	○	○	○
HTRF®	Cyclin D1, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p-DAP12 (panY)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	DAP12, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	DNA PKcs, Phospho-S2056	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	DNA PKcs, Total	●	X	X	○	○	○	○	○
HTRF®	EGF Receptor, Phospho-Y1068	●	X	X	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	EGF Receptor, Phospho-Y1068	●	●	○	○	○	○	○	○
HTRF®	EGF Receptor, Total	●	●	X	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	EGF Receptor, Total	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	EGFR, Phospho-Y1068 + EGFR Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	eIF2α, Phospho-S51	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	eIF2α, Phospho-S51 - OLD	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	eIF2α, Total	●	●	○	○	○	○	○	○
HTRF®	eIF2α, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	eIF4E, Phospho-S209	●	●	○	○	○	○	○	○
HTRF®	eIF4E, Phospho-S209	●	●	●	●	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	eIF4E, Total	●	○	○	○	○	○	○	○
HTRF®	eNOS, Phospho-S1177	●	●	●	○	○	○	○	○
HTRF®	eNOS, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ErbB2, Phospho-Y1221/1222	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ErbB2, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ErbB3, Phospho-Y1289	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ErbB3, Total	●	✗	✗	○	○	○	○	○
HTRF®	Advanced ERK, Phospho-T202/Y204	●	●	●	●	●	●	●	●
HTRF®	ERK1/2, Phospho-T202/Y204	●	●	●	●	●	●	●	●
HTRF®	ERK1/2, Total	●	●	●	●	●	●	●	●
AlphaLISA® SureFire® Ultra™	ERK1/2, Phospho-T202/Y204	●	●	●	○	●	○	○	○
AlphaLISA® SureFire® Ultra™	ERK1/2, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Estrogen Receptor α, Phospho-S118	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Estrogen Receptor α, Phospho-S167	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Estrogen Receptor α, Total	●	✗	✗	○	○	○	○	○
HTRF®	Estrogen Receptor α, Total	●	○	○	○	○	○	○	○
HTRF®	FAK, Phospho-Y397	●	●	○	○	○	○	○	○
HTRF®	FAK, Total	●	●	○	○	○	○	○	○
HTRF®	FGFR1, Phospho Y653/654	●	●	○	○	○	○	○	○
HTRF®	FGFR1, Total	●	✗	○	○	○	○	○	○
HTRF®	FGFR2, Phospho Y653/654	●	✗	○	○	○	○	○	○
HTRF®	FGFR2, Total	●	✗	○	○	○	○	○	○
HTRF®	FGFR3, Phospho Y653/654	●	○	○	○	○	○	○	○
HTRF®	FGFR3, Total	●	✗	○	○	○	○	○	○
HTRF®	FGFR4, Phospho Y642	●	✗	○	○	○	○	○	○
HTRF®	FGFR4, Total	●	✗	○	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

	HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™ FGF Receptor 1, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FGF Receptor 2, Phospho-Y653/654	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FGF Receptor 2, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FGF Receptor 3, Phospho-Y724	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FGF Receptor 3, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FGF Receptor 4, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FLT3, Phospho-Y591	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FLT3, Phospho-Y842	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ FLT3, Total	●	✗	✗	○	○	○	○	○
HTRF® FoxO1, Phospho-S256	●	●	●	●	○	○	○	○
HTRF® FoxO1, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™ GAPDH, Total	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ GAPDH, Total (Mouse / Human)	●	●	○	○	○	○	○	○
HTRF® GAPDH Housekeeping	●	●	●	●	●	○	○	✗
AlphaLISA® SureFire® Ultra™ GCN2, Phospho-T899	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ GCN2, Total	●	○	○	○	○	○	○	○
HTRF® GSK3 alpha, Phospho-S21	●	●	●	●	○	○	○	○
HTRF® GSK3 alpha, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™ GSK3 beta, Phospho-S9	●	●	●	○	○	○	○	○
HTRF® GSK3 beta, Phospho-S9	●	●	●	●	○	○	○	○
HTRF® GSK3 beta, Total	●	●	●	●	○	○	○	○
HTRF® H2AX, Phospho-S139	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™ p-HDAC4 (Ser246)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™ HDAC4, Total	●	●	○	○	○	○	○	○
HTRF® HER2, Phospho-Y1221/1222	●	✗	✗	✗	✗	✗	✗	✗
HTRF® HER2, Total	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	HER3, Phospho-Y1289	●	●	●	○	○	○	○	○
HTRF®	HER3, Total	●	●	●	○	○	○	○	○
HTRF®	HIF-1α	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p-Histone H2AX (Ser139)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Histone H2AX, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Histone H3, Phospho-S10/T11	●	●	○	○	○	○	○	○
HTRF®	Histone H3, Phospho-T3	●	●	●	○	○	○	○	○
HTRF®	Histone H3, Phospho-S10	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	Histone H3, Total	●	●	○	○	○	○	○	○
HTRF®	Histone H3, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	HSP27, Phospho-S15 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	HSP27, Phospho-S82 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	HSP27, Phospho-S78 NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	HSP27, Total NEW	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IGF-1 Rec.β, Phospho-Y1135/1136	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IGF-1 Rec.β, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IκBα, Phospho-S32/36	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IκBα, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IKKα, Phospho-S176/180	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IKKα, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IKKβ, Phospho-S177/181	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IKKβ, Total	●	X	X	○	○	○	○	○
HTRF®	IKKB, Phospho-S177/181	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	Insulin Rec. β, Phospho-Y1150/1151	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Insulin Rec. β, Total	●	●	●	○	○	○	○	○
HTRF®	IRAK3, Total	●	X	○	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	IRAK4, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IRAK4, Total	●	●	○	○	○	○	○	○
HTRF®	IRβ, Phospho Y1150/1151	●	●	●	○	○	○	○	○
HTRF®	IRβ, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IRF3, Phospho-S386	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IRF3, Total	●	○	○	○	○	○	○	○
HTRF®	IRF3, Phospho-S386	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	IRF5	●	✗	✗	✗	○	○	○	○
HTRF®	IRS1, Phospho-S307 mouse/S312 human	●	●	○	○	○	○	○	○
HTRF®	IRS1, Total	●	●	○	○	○	○	○	○
HTRF®	IRS2, Total	●	●	●	○	○	○	○	○
HTRF®	JAK1, Phospho-Y1034/35	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	JAK1, Total	●	✗	✗	○	○	○	○	○
HTRF®	JAK1, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	JAK2, Phospho-Y1007/08	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	JAK2, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	JAK2, Total	●	●	●	○	○	○	○	○
HTRF®	JAK3, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	JNK1/2/3, Phospho-T183/Y185	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	JNK 1/2total	●	○	○	○	○	○	○	○
HTRF®	JNK, Phospho-T183/Y185	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	KAP1, Phospho-S824	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	KAP1, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Lck, Phospho-Y505	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Lck, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	LRRK2, Phospho-S935	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	LRRK2, Phospho-S935	●	●	●	○	○	○	○	○
HTRF®	LRRK2, Total	●	●	●	○	○	○	○	○
HTRF®	MDA5, Total	●	○	○	○	○	○	○	○
HTRF®	MDA5 aggregation	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MEK1, Phospho-S218/222	●	●	○	○	○	○	○	○
HTRF®	MEK1, Phospho-S298	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MEK1, Total	●	●	○	○	○	○	○	○
HTRF®	MEK1, Total	●	●	●	●	●	○	○	○
HTRF®	MEK1/2, Phospho-S218/222	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	MKK4, Phospho-S257	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MKK4, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p-MLKL(Ser358)	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MLKL, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MLKL, Phospho-Ser345 (Mouse)	X	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MLKL, Total (Mouse)	●	●	○	○	○	○	○	○
HTRF®	MLKL, Phospho S35	●	○	○	○	○	○	○	○
HTRF®	MLKL, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	mTOR, Phospho-S2448	●	●	○	○	○	○	○	○
HTRF®	mTOR, Phospho-S2448	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	mTOR, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	NF-kB p65, Phospho-S536	●	●	●	○	○	○	○	○
HTRF®	NF-kB, Phospho-S536	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	NF-kB p65, Total	●	●	●	○	○	○	○	○
HTRF®	NF-kB, Total	●	●	●	○	●	○	○	●
HTRF®	PDGFRβ, Phospho-Y751	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PDHA1, Phospho-S293	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PDHA1, Total	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	PERK, Phospho-T980 (Mouse)	✗	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PERK, Total	●	●	●	○	○	○	○	○
HTRF®	Phospholamban, Phospho-S16	●	●	●	○	○	○	○	○
HTRF®	Phospholamban, Phospho-T17	●	●	●	○	○	○	○	○
HTRF®	Phospholamban, Total	●	●	●	○	○	○	○	○
HTRF®	P21	●	✗	✗	✗	○	○	○	○
AlphaLISA® SureFire® Ultra™	p38 MAPK, Phospho-T180/Y182	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p38 MAPK, Total	●	●	●	○	○	○	○	○
HTRF®	p38 MAPK, Phospho-T180/Y182	●	●	●	●	●	○	●	○
AlphaLISA® SureFire® Ultra™	p53 Ac-K382	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p53, Phospho-S15	●	○	○	○	○	○	○	○
HTRF®	p53, Phospho-S15	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	p53, Total	●	○	○	○	○	○	○	○
HTRF®	p53, Total	●	●	○	○	○	○	○	○
HTRF®	p62/SQSTM1, Phospho-S403	●	●	●	○	○	○	○	○
HTRF®	p62/SQSTM1, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p70 S6K, Phospho-T389	●	●	●	○	○	○	○	○
HTRF®	p70S6K, Phospho-T389	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	p70 S6K, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PD-1, Phospho-panY	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PD-1, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PDGFRβ, Phospho-Y751	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PDGFRβ, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p-PI3 Kinase p85 (Y458) (Mouse Specific)	○	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PI3 Kinase p85, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PKCθ, Phospho-T538	●	✗	✗	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	PKR, Phospho-T451	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PKR, Total	●	○	○	○	○	○	○	○
HTRF®	PLCg1, Phospho-Y783	●	●	●	○	○	○	○	○
HTRF®	PLCg1, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ1, Phospho-Y783	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ1, Phospho-S1248	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ1, Phospho-Y1253	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ1, Total	●	●	●	○	○	○	○	○
HTRF®	PLCg2, Phospho-Y1217	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ2, Phospho-Y753	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ2, Phospho-Y759	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ2, Phospho-Y1197	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ2, Phospho-Y1217	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCγ2, Total	●	○	○	○	○	○	○	○
HTRF®	PLCg2, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLK1, Phospho-T210	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLK1, Total	●	●	●	○	○	○	○	○
HTRF®	Progesterone Receptor, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RAB10, Phospho-T73	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RAB10, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Raf-1, Phospho-S338	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Raf-1, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Ras, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RB, Phospho-S780	●	●	○	○	○	○	○	○
HTRF®	RB, Phospho-S780	●	●	○	○	○	○	○	○
HTRF®	RB, Phospho-S807/811	●	●	○	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	RB, Total	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Rb, Phospho-S807/811	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Rb, Phospho-T821/826	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Rb, Total	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Rb, Phospho-S780 (Eu) +, Total Rb (Tb)	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Rb, Phospho-S807/811 (Eu) +, Total Rb (Tb)	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Rb, Phospho-T821/826 (Eu) +, Total Rb (Tb)	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Ret, Phospho-panY + Ret, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p-Ret (panY)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Ret, Total	●	●	○	○	○	○	○	○
HTRF®	RET, Phospho pan Y	●	●	●	○	○	○	○	○
HTRF®	RET, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIP3, Phospho-S227	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIP3, Total (Human)	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIP3, Phospho-T231/S232 (Mouse)	✗	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIP3, Total (Mouse)	✗	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIP3, Phospho-S166	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIPK1, Total	●	○	○	○	○	○	○	○
HTRF®	RIPK1, Phospho-S166	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	RIPK1, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	RPS6, Phospho-S235/236	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RPS6, Total	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	p-RPS6 (Ser235/236) (Eu) +, Total RPS6 (Tb)	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	p-RPS6 (SerS240/244) (Eu) +, Total RPS6 (Tb)	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	RSK1, Phospho-S380	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1, Phospho-T359	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1, Phospho-T573	●	○	○	○	○	○	○	○
HTRF®	RSK1, Phospho S380	●	●	●	○	○	○	○	○
HTRF®	RSK1, Phospho T359	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1, Total	●	●	●	○	○	○	○	○
HTRF®	RSK1, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	S6RP, Phospho-S240/244	●	●	●	○	○	○	○	○
HTRF®	S6RP, Phospho-S235/236	●	●	●	●	●	○	○	○
AlphaLISA® SureFire® Ultra™	SHP-1, Phospho-Y536	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SHP-1, Phospho-Y564	●	✗	✗	○	○	○	○	○
HTRF®	SHP-1, Phospho-Y564	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SHP-1, Total	●	✗	✗	○	○	○	○	○
HTRF®	SHP-1, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SHP-2, Phospho-Y542	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SHP-2, Phospho-Y580	●	●	●	○	○	○	○	○
HTRF®	SHP-2, Phospho-Y542	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SHP-2, Total	●	●	●	○	○	○	○	○
HTRF®	SHP-2, Total	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	SLP-76, Phospho-S376 + SLP-76, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76, Phospho-S376	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76, Phospho-S376 (Mouse, Human)	●	●	○	○	○	○	○	○
HTRF®	SLP-76, Phospho-S376	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	SLP-76, Phospho-S376 (mouse)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76, Phospho-Y113	●	○	○	○	○	○	○	○

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	SLP-76, Phospho-Y145	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76, Phospho-Y145 (Mouse) NEW	●	●	○	○	○	○	○	○
HTRF®	SLP-76, Phospho-Y145	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	SLP-76, Total	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76, Total (Mouse, Human)	●	●	○	○	○	○	○	○
HTRF®	SLP-76, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	SLP-76, Total (mouse)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD1, Phospho-S463/465	●	●	●	○	○	○	○	○
HTRF®	SMAD1, Phospho-S463/465	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD1, Total	●	●	●	○	○	○	○	○
HTRF®	SMAD1, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD2, Phospho-S465/467	●	●	●	○	○	○	○	○
HTRF®	SMAD2, Phospho-S465/467	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD2, Total	●	●	●	○	○	○	○	○
HTRF®	SMAD2, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD3, Phospho-S423/425	●	●	●	○	○	○	○	○
HTRF®	SMAD3, Phospho-S423/425	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD3, Total	●	●	●	○	○	○	○	○
HTRF®	SMAD3, Total	●	●	○	○	○	○	○	○
HTRF®	SMAD4, Phospho-T277	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMAD4, Total	●	●	●	○	○	○	○	○
HTRF®	SMAD4, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	SMARCA4, Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SQSTM1 p62, Phospho-S351	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SQSTM1 p62, Total	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Src, Phospho-Y419	●	○	○	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	c-Src, Phospho-Y419	●	●	●	●	○	○	○	○
HTRF®	c-Src, Total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT1, Phospho-S727	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT1, Phospho-Y701	●	●	○	○	○	○	○	○
HTRF®	STAT1, Phospho-Y701	●	●	X	X	X	X	X	X
AlphaLISA® SureFire® Ultra™	STAT1, Total	●	○	○	○	○	○	○	○
HTRF®	STAT1, Total	●	●	X	X	X	X	X	X
Alpha SureFire® Ultra™ Multiplex	p-STAT1 (Y701) (Eu) +, Total STAT1 (Tb)	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	p-STAT1 (Ser727) (Eu) +, Total STAT1 (Tb)	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT3, Phospho-Y705	●	●	●	○	○	○	○	○
HTRF®	STAT3, Phospho-Y705	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT3, Total	●	●	●	○	○	○	○	○
HTRF®	STAT3, Total	●	●	●	●	●	○	○	○
AlphaLISA® SureFire® Ultra™	STAT4, Phospho-Y693	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT4, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT5, Phospho-Y694/699	●	○	○	○	○	○	○	○
HTRF®	STAT5, Phospho-Y694	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT5, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT6, Phospho-Y641	●	●	○	○	○	○	○	○
HTRF®	STAT6, Phospho-Y641	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT6, Total	●	○	○	○	○	○	○	○
HTRF®	STAT6, Total NEW	●	●	X	X	X	X	X	X
HTRF®	STING, Phospho-S366	●	X	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STING, Total	●	○	○	○	○	○	○	○
HTRF®	STING (m), Phospho-S365	X	●	X	X	X	X	X	X
HTRF®	STING (m), Total	●	●	X	X	X	X	X	X

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	STING, Total	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SYK, Phospho-Y352	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SYK, Phospho-Y525/526	●	●	●	○	○	○	○	○
HTRF®	SYK, Phospho-Y525	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	SYK, Total	●	✗	✗	○	○	○	○	○
HTRF®	SYK, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	Tau, Phospho-S396 NEW	●	●	○	○	○	○	○	○
HTRF®	Tau, Phospho-S202/T205	●	●	●	○	○	○	○	○
HTRF®	Tau, Phospho-S356	●	●	●	○	○	○	○	○
HTRF®	Tau, Phospho-S404	●	●	●	○	○	○	○	○
HTRF®	Tau, Phospho-S422	●	●	●	○	○	○	○	○
HTRF®	Tau, Phospho-T181	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Tau, Total NEW	●	●	○	○	○	○	○	○
HTRF®	Tau, Total	●	✗	○	○	○	○	○	○
HTRF®	TDP-43 aggregated	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	TDP43, Phospho-S409/410	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	TDP43, Total	●	✗	✗	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	TBK1, Phospho-S172	●	●	●	○	○	○	○	○
HTRF®	TBK1, Phospho-S172	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	TBK1, Total	●	●	●	○	○	○	○	○
HTRF®	TBK1, Total	●	●	●	○	○	○	○	○
HTRF®	Tie2, Phospho-Y992	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	TrkA, Phospho-Y674/675 / TrkB, Phospho-Y706/707	●	●	●	○	○	○	○	○
HTRF®	TYK2, Phospho-Y1054/55	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	TYK2, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	Ubiquitin, Phospho-S65	●	●	●	○	○	○	○	○

● Compatible / ● Predicted / X Not compatible / ○ Unknown

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	Ubiquitin, Total	●	●	○	○	○	○	○	○
HTRF®	Ubiquitin, Phospho-S65	●	●	●	●	●	○	●	●
AlphaLISA® SureFire® Ultra™	ULK1, Phospho-S556	●	X	X	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ULK1, Phospho-S758	●	X	X	○	○	○	○	○
HTRF®	ULK1 phosho-S757 mouse/S758 human	●	●	○	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	ULK1, Total	●	X	X	○	○	○	○	○
HTRF®	VASP, Phospho-S157	●	●	○	○	○	○	○	○
HTRF®	VASP, Phospho-S239 kit	●	●	○	○	○	○	○	○
HTRF®	VASP, Total	●	●	○	○	○	○	○	○
HTRF®	VEGFR2, Phospho-Y1175	●	●	○	○	○	○	○	○
HTRF®	VEGFR2, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	VEGFR2, Phospho-Y1175	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	VEGFR2, Phospho-Y1214	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	VEGFR2, Total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	YAP, Phospho-S127	●	●	●	○	○	○	○	○
HTRF®	YAP, Phospho-S127	●	●	●	○	●	○	○	○
AlphaLISA® SureFire® Ultra™	YAP, Total	●	X	X	○	○	○	○	○
HTRF®	YAP, Total	●	●	○	●	●	○	○	○
AlphaLISA® SureFire® Ultra™	ZAP-70, Phospho-Y319	●	●	○	○	○	○	○	○
HTRF®	ZAP-70, Phospho-Y319	●	X	X	X	X	X	X	X
AlphaLISA® SureFire® Ultra™	ZAP-70, Total	●	●	○	○	○	○	○	○
HTRF®	ZAP-70, Total	●	X	X	X	X	X	X	X

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 ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.