



Species compatibility.

October 2024

Cell signaling: phospho- and total-protein assays

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

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	4E-BP1, Phospho-T37/46	●	●	○	○	○	○	○	○
HTRF™	4E-BP1, Phospho-T37/46	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	4E-BP1, Total	●	●	○	○	○	○	○	○
HTRF™	α-SMA	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	α-Synuclein Aggregate Specific	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	α-Synuclein, Phospho-S129	●	●	○	○	○	○	○	○
HTRF™	α-Synuclein-S129 Aggregation	●	●	○	○	○	○	○	○
HTRF™	α-Synuclein, Phospho-S129	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	α-Synuclein, Total	●	●	○	○	○	○	○	○
HTRF™	α-Synuclein, Total	●	X	○	○	○	○	○	○
HTRF™	α-Tubulin Housekeeping	●	●	●	●	●	○	●	●
AlphaLISA™ SureFire® Ultra™	ACC, Phospho-S79	●	X	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	ACC, Phospho-S79	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ACC, Total	●	●	○	○	○	○	○	○
HTRF™	ADAR1, Total	●	○	○	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Aiolos, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1/2/3, Phospho-S473	●	○	○	○	○	○	○	○
HTRF™	AKT 1/2/3, Phospho-S473	●	●	●	●	●	●	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1/2/3, Phospho-T308	●	○	○	○	○	○	○	○
HTRF™	AKT 1/2/3, Phospho-T308	●	●	●	●	●	●	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1/2/3, Phospho-T450	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1/2/3, Total	●	○	○	○	○	○	○	○
HTRF™	AKT 1/2/3 Total	●	●	●	○	●	○	○	○
Alpha SureFire® Ultra™ Multiplex	AKT 1/2/3, Phospho-S473 + AKT 1, Total	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	AKT 1/2/3, Phospho-T308 + AKT, Total	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	AKT 1/2/3, Phospho-T450 + AKT, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1, Phospho-S129	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1, Phospho-S473	●	○	○	○	○	○	○	○
HTRF™	AKT 1, Phospho-S473	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AKT 1, Total	●	○	○	○	○	○	○	○
HTRF™	AKT 1, Total	●	●	●	●	○	○	○	○
HTRF™	AKT 2, Phospho-S473	●	●	●	●	○	○	○	○
HTRF™	AKT 2, Total	●	●	●	●	○	○	○	○
HTRF™	AKT 3, Phospho-S473	●	●	●	○	○	○	○	○
HTRF™	AKT 3, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ALK, Phospho-Y1586	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ALK, Phospho-Y1604	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	ALK, Phospho-Y1604 + ALK, Total	●	✗	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	ALK, Phospho-Y1604	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	ALK, Total	●	○	○	○	○	○	○	○
HTRF™	ALK, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AMPKα1/2, Phospho-T172	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	AMPKα1/2, Phospho-T172 + AMPKα1/2, Total	●	○	○	○	○	○	○	○
HTRF™	AMPK, Phospho-T172	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	AMPKα1/2, Total	●	✗	○	○	○	○	○	○
HTRF™	AMPK, Total	●	●	●	○	○	○	○	○
HTRF™	Androgen Receptor, Full-length	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Androgen Receptor, Total	●	●	○	○	○	○	○	○
HTRF™	Androgen Receptor, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	Androgen Receptor, Variant 7	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	AP2, Total	●	●	●	○	●	○	○	○
AlphaLISA™ SureFire® Ultra™	ATF-2, Phospho-T71	●	●	○	○	○	○	○	○
HTRF™	ATF2, Phospho-T71	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATF-2, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATF-4, Total	●	●	○	○	○	○	○	○
HTRF™	ATF4, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATG14, Phospho-S29	●	●	○	○	○	○	○	○
HTRF™	ATG14, Phospho-S29	●	●	●	○	○	○	○	○
HTRF™	ATG14, Total	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATG16L1, Phospho-S278	●	●	○	○	○	○	○	○
HTRF™	ATG16L1, Phospho-S278	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATG16L1, Total	●	●	○	○	○	○	○	○
HTRF™	ATG16L1, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATM, Phospho-S1981	●	✗	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	ATM, Phospho-S1981	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ATM, Total	●	●	○	○	○	○	○	○
HTRF™	ATM, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Aurora A, Phospho-T288	●	✗	○	○	○	○	○	○
HTRF™	Aurora A, Phospho-T288	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Aurora A, Total	●	✗	○	○	○	○	○	○
HTRF™	Aurora A, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Aurora B, Phospho-T232	●	○	○	○	○	○	○	○
HTRF™	Aurora B, Phospho-T232	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Aurora B, Total	●	○	○	○	○	○	○	○
HTRF™	Aurora B, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BAD, Phospho-S112 NEW	●	●	○	○	○	○	○	○
HTRF™	BAD, Phospho-S112	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BAD, Phospho-S136 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BAD, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BCL-2, Phospho-S70	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BCL-2, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BCL-XL, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Beclin-1, Total	●	●	○	○	○	○	○	○
HTRF™	β-arrestin 1, Total	●	●	●	○	●	○	○	○
HTRF™	β-arrestin 2, Total	●	●	●	○	●	○	○	○
AlphaLISA™ SureFire® Ultra™	β-Catenin, Phospho-S37	●	●	○	○	○	○	○	○
HTRF™	β-Catenin, Phospho-T41/S33/37	●	●	●	●	○	○	●	○
AlphaLISA™ SureFire® Ultra™	β-Catenin, Phospho-T41/S45	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	β-Catenin, Phospho-S675	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	β-Catenin, Total	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	β-Catenin, Total	●	●	●	●	○	○	●	○
AlphaLISA™ SureFire® Ultra™	BRD2, Phospho-S37	●	●	○	○	○	○	○	○
HTRF™	BRD2, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BRD2, Total	●	●	○	○	○	○	○	○
HTRF™	BRD3, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BRD3, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BRD4, Total	●	✗	○	○	○	○	○	○
HTRF™	BRD4, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BRD7, Total	●	●	○	○	○	○	○	○
HTRF™	BRD7, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BRD9, Total	●	●	○	○	○	○	○	○
HTRF™	BRD9, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	BTK, Phospho-Y223	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	BTK, Phospho-Y223 + BTK, Total	●	○	○	○	○	○	○	○
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-Y83	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CD3 Zeta, Phospho-Y142	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CD3 Zeta, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDC6, Phospho-S54	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDC6, Phospho-S106	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDC6, Total	●	✗	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	CDC25C, Phospho-S216	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDC25C, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK1, Phospho-T14	●	✗	○	○	○	○	○	○
HTRF™	CDK1, Phospho-T14	●	●	●	✗	✗	✗	✗	✗
Alpha SureFire® Ultra™ Multiplex	CDK1, Phospho-T14 + CDK1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK1, Phospho-T161	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK1, Phospho-Y15	●	●	○	○	○	○	○	○
HTRF™	CDK1, Phospho-Y15	●	●	●	✗	✗	✗	✗	✗
Alpha SureFire® Ultra™ Multiplex	CDK1, Phospho-Y15 + CDK1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK1, Total	●	✗	✗	○	○	○	○	○
HTRF™	CDK1, Total	●	✗	○	●	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	CDK2, Phospho-T160	●	●	○	○	○	○	○	○
HTRF™	CDK2, Phospho-Y15	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK2, Total	●	●	○	○	○	○	○	○
HTRF™	CDK2, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK4, Total	●	✗	○	○	○	○	○	○
HTRF™	CDK4, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK5, Total	●	●	○	○	○	○	○	○
HTRF™	CDK5, Total	●	●	●	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	CDK6, Total	●	✗	○	○	○	○	○	○
HTRF™	CDK6, Total	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	CDK7, Total	●	✗	○	○	○	○	○	○
HTRF™	CDK7, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	CDK8, Total	●	●	○	○	○	○	○	○
HTRF™	CDK8, Total	●	●	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	CDK9, Total	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	CDK9, Total	●	●	●	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	CDK12, Total	●	●	○	○	○	○	○	○
HTRF™	CDK12, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	Cereblon, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-FOS, Phospho-S32	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-FOS, Total	●	✗	○	○	○	○	○	○
HTRF™	cGAS, 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™	c-Jun, Phospho-S63	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-Jun, Phospho-S73	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-Jun, Total	●	●	○	○	○	○	○	○
HTRF™	CK1α, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-Met, Phospho-Y1234/1235	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-Met, Total	●	✗	○	○	○	○	○	○
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	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	c-Myc, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cofilin, Phospho-S3	●	●	○	○	○	○	○	○
HTRF™	Cofilin, Phospho-S3	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Cofilin, Phospho-S3 + Cofilin, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cofilin, Total	●	●	○	○	○	○	○	○
HTRF™	Cofilin, Total	●	●	●	○	○	○	○	○
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™	CREBBP, Total	●	●	○	○	○	○	○	○
HTRF™	CREBBP, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin A2, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin B1, Phospho-S126	●	✗	○	○	○	○	○	○
HTRF™	Cyclin B1, Phospho-S126	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin B1, Total	●	✗	○	○	○	○	○	○
HTRF™	Cyclin B1, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin D1, Phospho-T286	●	●	○	○	○	○	○	○
HTRF™	Cyclin D1, Phospho-T286	●	✗	○	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin D1, Total	●	●	○	○	○	○	○	○
HTRF™	Cyclin D1, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin D2, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin D3, Total	●	✗	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	Cyclin E1, Phospho-T77	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Cyclin E1, Total	●	✗	○	○	○	○	○	○
HTRF™	Cyclin E1, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	DAP12, Aggregate	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DAP12, Phospho-panY	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DAP12, Phospho-Y91	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	DAP12, Phospho-Y91 + DAP12, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DAP12, Phospho-Y92 (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DAP12, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DAP12, Total (mouse)	✗	●	○	○	○	○	○	○
HTRF™	DDR1, Phospho-PanY	●	✗	○	○	○	○	○	○
HTRF™	DDR1, Phospho-Y513	●	✗	○	○	○	○	○	○
HTRF™	DDR1, Phospho-Y796	●	✗	○	○	○	○	○	○
HTRF™	DDR1, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DJ1, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DNA PKcs, Phospho-S2056	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	DNA PKcs, Total	●	✗	○	○	○	○	○	○
HTRF™	EGF Receptor DEL19, Total	●	●	●	●	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	EGF Receptor L858R, Total	●	✗	○	○	○	○	○	○
HTRF™	EGF Receptor L858R, Total	●	●	●	●	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	EGF Receptor, Phospho-Y1068	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	EGF Receptor Phospho-Y1068 + EGF Receptor, Total	●	●	○	○	○	○	○	○
HTRF™	EGF Receptor, Phospho-Y1068	●	✗	✗	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	EGF Receptor, Total	●	✗	○	○	○	○	○	○
HTRF™	EGF Receptor, Total	●	●	✗	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eIF2α, Phospho-S51	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
Alpha SureFire® Ultra™ Multiplex	eIF2α, Phospho-S51 + eIF2α, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eIF2α, Total	●	●	○	○	○	○	○	○
HTRF™	eIF2α, Total	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eIF4E, Phospho-S209	●	●	○	○	○	○	○	○
HTRF™	eIF4E, Phospho-S209	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eIF4E, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eNOS, Phospho-S1177 NEW	●	✗	○	○	○	○	○	○
HTRF™	eNOS, Phospho-S1177	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eNOS, Phospho-T495 NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	eNOS, Total NEW	●	✗	○	○	○	○	○	○
HTRF™	eNOS, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ErbB2, Phospho-Y1221/1222	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	ErbB2 Phospho-Y1221/1222 + ErbB2, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ErbB2, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ErbB3, Phospho-Y1289	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Erb3, Phospho-T1289 + Erb3, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ErbB3, Total	●	✗	○	○	○	○	○	○
HTRF™	ERK advanced, Phospho-T202/Y204	●	●	●	●	●	●	●	●
AlphaLISA™ SureFire® Ultra™	ERK1/2, Phospho-T202/Y204	●	●	○	○	●	○	○	○
HTRF™	ERK1/2, Phospho-T202/Y204	●	●	●	●	●	●	●	●
Alpha SureFire® Ultra™ Multiplex	ERK1/2 Phospho-T202/Y204 + ERK1/2, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ERK1/2, Total	●	●	○	○	○	○	○	○
HTRF™	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	●	○	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	FAK, Phospho-Y397	●	●	○	○	○	○	○	○
HTRF™	FAK, Phospho-Y397	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FAK, Phospho-Y576/577	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FAK, Total	●	●	○	○	○	○	○	○
HTRF™	FAK, Total	●	●	○	○	○	○	○	○
HTRF™	FGFR1, Phospho-Y653/654	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 1, Total	●	○	○	○	○	○	○	○
HTRF™	FGFR1, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 2, Phospho-Y653/654	●	○	○	○	○	○	○	○
HTRF™	FGFR2, Phospho Y653/654	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 2, Total	●	○	○	○	○	○	○	○
HTRF™	FGFR2, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 3, pan-Tyr	●	✗	○	○	○	○	○	○
HTRF™	FGFR3, Phospho-Y653/654	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 3, Phospho-Y724	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 3, Total	●	○	○	○	○	○	○	○
HTRF™	FGFR3, Total	●	✗	○	○	○	○	○	○
HTRF™	FGFR4, Phospho-Y642	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FGF Receptor 4, Total	●	○	○	○	○	○	○	○
HTRF™	FGFR4, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FLT3, Phospho-Y591	●	✗	○	○	○	○	○	○
HTRF™	FLT3, Phospho-Y589/591	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FLT3, Phospho-Y842	●	✗	○	○	○	○	○	○
HTRF™	FLT3, Phospho-Y842	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	FLT3, Total	●	✗	○	○	○	○	○	○
HTRF™	FLT3, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	FOXO1, Phospho-T600	●	✗	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	FOXM1, Total	●	✗	○	○	○	○	○	○
HTRF™	FoxO1, Phospho-S256	●	●	●	●	○	○	○	○
HTRF™	FoxO1, Total	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	GAPDH, Total (human)	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	GAPDH, Total	●	●	○	○	○	○	○	○
HTRF™	GAPDH Housekeeping	●	●	●	●	●	○	○	✗
AlphaLISA™ SureFire® Ultra™	GCN2, Phospho-T899	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	GCN2, Total	●	○	○	○	○	○	○	○
HTRF™	GSK3α, Phospho-S21	●	●	●	●	○	○	○	○
HTRF™	GSK3α, Total	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	GSK3β, Phospho-S9	●	●	○	○	○	○	○	○
HTRF™	GSK3β, Phospho-S9	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	GSK3β, Total	●	●	○	○	○	○	○	○
HTRF™	GSK3β, Total	●	●	●	●	○	○	○	○
HTRF™	GSPT1, Total	●	○	○	○	○	○	○	○
HTRF™	H2AX, Phospho-S139	●	●	●	●	○	○	○	○
HTRF™	H2AX, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HDAC4, Phospho-S246	●	●	○	○	○	○	○	○
HTRF™	HDAC4, Phospho-S246	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HDAC4, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HDAC5, Phospho-S259	●	●	○	○	○	○	○	○
HTRF™	HDAC5, phospho-S259	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HDAC5, Total	●	●	○	○	○	○	○	○
HTRF™	HDAC5, Total	●	●	○	○	○	○	○	○
HTRF™	HER2, Phospho-Y1221/1222	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	HER2, Total	●	●	●	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	HER3, Phospho-Y1289	●	●	●	○	○	○	○	○
HTRF™	HER3, Total	●	●	●	○	○	○	○	○
HTRF™	HIF-1α	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Histone H2AX, Phospho-S139	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Histone H2AX, Phospho-S139 + Histone H2AX, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Histone H2AX, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Histone H3, Phospho-S10/T11	●	●	○	○	○	○	○	○
HTRF™	Histone H3, Phospho-S10	●	●	●	●	○	○	○	○
HTRF™	Histone H3, Phospho-T3	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Histone H3, Total	●	●	○	○	○	○	○	○
HTRF™	Histone H3, Total	●	●	●	●	○	○	○	○
HTRF™	HPK1, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HSP27, Phospho-S15	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HSP27, Phospho-S78	●	○	○	○	○	○	○	○
HTRF™	HSP27, Phospho-S78	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HSP27, Phospho-S82	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	HSP27, Total	●	○	○	○	○	○	○	○
HTRF™	HSP27, Total	●	✗	○	○	○	○	○	○
HTRF™	IDO1, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IGF-1 Rec.β, Phospho-Y1135/1136	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	IGF-1 Rec.β, Phospho-Y1135/1136 + IGF-1 Rec.β, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IGF-1 Rec.β, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IKAROS, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IKKα, Phospho-S176/180	●	●	○	○	○	○	○	○
HTRF™	IKKα, Phospho-S176/180	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IKKα, Total	●	○	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	IKKα, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IκBα, Phospho-S32/36	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IκBα, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IKKβ, Phospho-S177/181	●	✗	○	○	○	○	○	○
HTRF™	IKKβ, Phospho-S177/181	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IKKβ, Total	●	✗	○	○	○	○	○	○
HTRF™	IKZF1, Total	●	●	○	○	○	○	○	○
HTRF™	IKZF3, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Insulin Rec. β, Phospho-Y1150/1151	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Insulin Rec. β, Phospho-Y1150/1151 + Insulin Rec.β, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Insulin Rec. β, Total	●	●	○	○	○	○	○	○
HTRF™	IRAK3, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IRAK4, Phospho-T345/S346	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IRAK4, Total	●	●	○	○	○	○	○	○
HTRF™	IRAK4, Total	●	●	○	○	○	○	○	○
HTRF™	IRβ, Phospho-Y1150/1151	●	●	●	○	○	○	○	○
HTRF™	IRβ, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IRF3, Phospho-S386	●	✗	○	○	○	○	○	○
HTRF™	IRF3, Phospho-S386	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	IRF3, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IRF5, Phospho-S446	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	IRF5, Total	●	●	○	○	○	○	○	○
HTRF™	IRF5, Total	●	○	○	○	○	○	○	○
HTRF™	IRS1, Phospho-S307 mouse/S312	●	●	○	○	○	○	○	○
HTRF™	IRS1, Total	●	●	○	○	○	○	○	○
HTRF™	IRS2, Total	●	●	●	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	JAK1, Phospho-Y1034/35	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	JAK1, Total	●	✗	○	○	○	○	○	○
HTRF™	JAK1, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	JAK2, Phospho-Y1007/08	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	JAK2, Total	●	●	○	○	○	○	○	○
HTRF™	JAK2, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	JAK3, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	JNK1/2/3, Phospho-T183/Y185	●	●	○	○	○	○	○	○
HTRF™	JNK, Phospho-T183/Y185	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	JNK 1/2, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	KAP1, Phospho-S824	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	KAP1, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	KRAS, Total	●	✗	○	○	○	○	○	○
HTRF™	KRAS, Total	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Lamin A/C, Phospho-S22	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Lamin A/C, Total	●	●	○	○	○	○	○	○
HTRF™	LC3B-II, Total	●	●	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Lck, Phospho-Y505	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Lck, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	LRRK2, Phospho-S910	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	LRRK2, Phospho-S935	●	●	○	○	○	○	○	○
HTRF™	LRRK2, Phospho-S935	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	LRRK2, Phospho-S955	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	LRRK2, Total	●	●	○	○	○	○	○	○
HTRF™	LRRK2, Total	●	●	●	○	○	○	○	○
HTRF™	MDA5 aggregation	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	MDA5, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MDM2, Phospho-S166	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MDM2, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MEK1, Phospho-S218/222	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	MEK1, Phospho-S218/222 + MEK1, Total	●	●	○	○	○	○	○	○
HTRF™	MEK1, Phospho-S298	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MEK1, Total	●	●	○	○	○	○	○	○
HTRF™	MEK1, Total	●	●	●	●	●	○	○	○
HTRF™	MEK1/2, Phospho-S218/222	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MEK2, Phospho-S217/221 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MEK2, Total NEW	●	●	○	○	○	○	○	○
HTRF™	MerTK, Phospho-Y749	●	✗	○	○	○	○	○	○
HTRF™	MerTK, Total	●	✗	○	○	○	○	○	○
HTRF™	MET, Phospho-Y1234/1235	●	✗	○	○	○	○	○	○
HTRF™	MET, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MKK4, Phospho-S257	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MKK4, Total	●	○	○	○	○	○	○	○
HTRF™	MLKL, Phospho-S35	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MLKL, Phospho-S358	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MLKL, Phospho-S345 (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MLKL, Total (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MLKL, Total	●	✗	○	○	○	○	○	○
HTRF™	MLKL, Total	●	○	○	○	○	○	○	○
HTRF™	MSH3, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	mTOR, Phospho-S2448	●	●	○	○	○	○	○	○
HTRF™	mTOR, Phospho-S2448	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	mTOR, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	MYD88, Total	●	✗	○	○	○	○	○	○
HTRF™	MYD88, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	NF-κB p65, Phospho-S536	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	NF-κB p65, Phospho-S536 + NF-κB p65, Total	●	●	○	○	○	○	○	○
HTRF™	NF-κB, Phospho-S536	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	NF-κB p65, Total	●	●	○	○	○	○	○	○
HTRF™	NF-κB, Total	●	●	●	○	●	○	○	●
AlphaLISA™ SureFire® Ultra™	NF-κB2 p100, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	NF-κB2 p100/p52, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	NPM, Phospho-T199	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	NPM, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Nrf2, Phospho-S40	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Nrf2, Total	●	✗	○	○	○	○	○	○
HTRF™	Nrf2, Total	●	●	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	p16, INK4A, Total NEW	●	✗	○	○	○	○	○	○
HTRF™	p21	●	✗	✗	✗	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p21 CiP1, Total NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p27 KIP1, Phospho-S10	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p27 KIP1, Total	●	●	○	○	○	○	○	○
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	●	○	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Acc p53 Phospho-K382 + p53, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p53, Phospho-S15	●	○	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	p53, Phospho-S15	●	✗	✗	✗	✗	✗	✗	✗
Alpha SureFire® Ultra™ Multiplex	p53, Phospho-S15 + p53, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p53, Total	●	○	○	○	○	○	○	○
HTRF™	p53, Total	●	●	○	○	○	○	○	○
HTRF™	p62/SQSTM1, Phospho-S403	●	●	●	○	○	○	○	○
HTRF™	p62/SQSTM1, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p70 S6K, Phospho-T389	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	p70 S6K, Phospho-T389 + p70S6K, Total	●	●	○	○	○	○	○	○
HTRF™	p70S6K, Phospho-T389	●	●	●	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p70S6K, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	p300, Total	●	✗	○	○	○	○	○	○
HTRF™	p300, Total	●	●	○	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Parkin, Phospho-S65	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Parkin, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PARP1 Cleaved (Asp 214) NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PD-1, Phospho-panY	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PD-1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PDGF Rec β, Phospho-Y751	●	●	○	○	○	○	○	○
HTRF™	PDGF Rec β, Phospho-Y751	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	PDGF Rec β, Phospho-Y751 + PDGF Rec β, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PDGF Rec β, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PDHA1, Phospho-S293	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PDHA1, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PERK, Phospho-T980 (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PERK, Phospho-T982 NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PERK, Total	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	Phospholamban, Phospho-S16	●	●	●	○	○	○	○	○
HTRF™	Phospholamban, Phospho-T17	●	●	●	○	○	○	○	○
HTRF™	Phospholamban, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PI3 Kinase p85, Phospho-Y458 (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PI3 Kinase p85, Phospho-Y467 NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PI3 Kinase p85, Total (human) NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PI3 Kinase p85, Total (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PINK1, Total	●	✗	○	○	○	○	○	○
HTRF™	PINK1, Total	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PKCθ, Phospho-T538	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PKR, Phospho-T451	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PKR, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ1, Phospho-Y783	●	●	○	○	○	○	○	○
HTRF™	PLCγ1, Phospho-Y783	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ1, Phospho-S1248	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ1, Phospho-Y1253	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ1, Total	●	●	○	○	○	○	○	○
HTRF™	PLCγ1, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ2, Phospho-Y753	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ2, Phospho-Y759	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ2, Phospho-Y1197	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ2, Phospho-Y1217	●	✗	○	○	○	○	○	○
HTRF™	PLCγ2, Phospho-Y1217	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLCγ2, Total	●	✗	○	○	○	○	○	○
HTRF™	PLCγ2, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLK1, Phospho-T210	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	PLK1, Phospho-T210	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PLK1, Total	●	●	○	○	○	○	○	○
HTRF™	PLK1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PRAS40, Phospho-T246 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PRAS40, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Progesterone Receptor, Phospho-S190 NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Progesterone Receptor, Total NEW	●	✗	○	○	○	○	○	○
HTRF™	Progesterone Receptor, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PTEN, Phospho-T366	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	PTEN, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB8A, Phospho-S111	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB8A, Phospho-T72	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB8A, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB10, Phospho-T73	●	●	○	○	○	○	○	○
HTRF™	RAB10, Phospho-T73	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB10, Total	●	●	○	○	○	○	○	○
HTRF™	RAB10, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB12, Phospho-S106	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB29, Phospho-S72	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAB29, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Raf-1, Phospho-S259	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Raf-1, Phospho-S338	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Raf-1, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RAPTOR, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Ras, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Rb, Phospho-S780	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	Rb, Phospho-S780	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Rb, Phospho-S780 + Rb, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Rb, Phospho-S807/811	●	●	○	○	○	○	○	○
HTRF™	Rb, Phospho-S807/811	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Rb, Phospho-S807/811 + Rb, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Rb, Phospho-T821 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Rb, Phospho-T821/826	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	Rb, Phospho-S821/826 + Rb, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Rb, Phospho-T826 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Rb, Total	●	✗	○	○	○	○	○	○
HTRF™	Rb, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RBM39, Total	●	●	○	○	○	○	○	○
HTRF™	RBM39, Total	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	RET, Phospho-panY + Ret, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RET, Phospho-panY	●	✗	○	○	○	○	○	○
HTRF™	RET, Phospho-panY	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RET, Total	●	✗	○	○	○	○	○	○
HTRF™	RET, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIP3, Phospho-S166	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIP3, Phospho-S227	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIP3, Phospho-T231/S232 (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIP3, Total (human)	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIP3, Total (mouse)	✗	●	○	○	○	○	○	○
HTRF™	RIPK1, Phospho-S166	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	RIPK1, Phospho-S320 NEW	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIPK1, Phospho-S321 NEW	✗	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	RIPK1, Total (human)	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RIPK1, Total (mouse) NEW	✗	●	○	○	○	○	○	○
HTRF™	RIPK1, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	RNA POL II, Phospho-S2	●	●	○	○	○	○	○	○
HTRF™	RNA POL II, Phospho-S2 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RNA POL II, Phospho-S5	●	●	○	○	○	○	○	○
HTRF™	RNA POL II, Phospho-S5 NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RNA POL II, Total	●	●	○	○	○	○	○	○
HTRF™	RNA POL II, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RPS6, Phospho-S240/244	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RPS6, Phospho-S235/236	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	RPS6, Phospho-S235/236 + RPS6, Total	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	RPS6, Phospho-S240/244 + RPS6, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RPS6, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RSK1, Phospho-S380	●	●	○	○	○	○	○	○
HTRF™	RSK1, Phospho-S380	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RSK1, Phospho-T359	●	●	○	○	○	○	○	○
HTRF™	RSK1, Phospho-T359	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RSK1, Phospho-T573	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	RSK1, Total	●	●	○	○	○	○	○	○
HTRF™	RSK1, Total	●	●	●	○	○	○	○	○
HTRF™	S6RP, Phospho-S235/236	●	●	●	●	●	○	○	○
HTRF™	SALL4, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SHP-1, Phospho-Y536	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SHP-1, Phospho-Y564	●	●	○	○	○	○	○	○
HTRF™	SHP-1, Phospho-Y564	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	SHP-1, Total	●	✗	○	○	○	○	○	○
HTRF™	SHP-1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SHP-2, Phospho-Y542	●	●	○	○	○	○	○	○
HTRF™	SHP-2, Phospho-Y542	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SHP-2, Phospho-Y580	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SHP-2, Total	●	●	○	○	○	○	○	○
HTRF™	SHP-2, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SLP-76, Phospho-S376 (human)	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SLP-76, Phospho-S376	●	●	○	○	○	○	○	○
HTRF™	SLP-76, Phospho-S376	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	SLP-76, Phospho-S376 (mouse)	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	SLP-76, Phospho-S376 + SLP-76, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SLP-76, Phospho-Y113	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SLP-76, Phospho-Y145	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SLP-76, Phospho-Y145	●	●	○	○	○	○	○	○
HTRF™	SLP-76, Phospho-Y145	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	SLP-76, Total (human)	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SLP-76, Total	●	●	○	○	○	○	○	○
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	●	●	●	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
Alpha SureFire® Ultra™ Multiplex	SMAD2, Phospho-S465/467 + SMAD2, Total	●	●	○	○	○	○	○	○
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™	SMARCA2, Total	●	✗	○	○	○	○	○	○
HTRF™	SMARCA2, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SMARCA4, Total	●	●	○	○	○	○	○	○
HTRF™	SMARCA4, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SQSTM1 p62, Phospho-S351	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SQSTM1 p62, Phospho-S403	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SQSTM1 p62, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	c-Src, Phospho-Y419	●	○	○	○	○	○	○	○
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	●	●	✗	✗	✗	✗	✗	✗
Alpha SureFire® Ultra™ Multiplex	STAT1, Phospho-Y701 + STAT1, Total	●	●	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	STAT1, Phospho-S727 + STAT1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	STAT1, Total	●	●	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
HTRF™	STAT1, Total	●	●	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	STAT3, Phospho-Y705	●	●	○	○	○	○	○	○
HTRF™	STAT3, Phospho-Y705	●	●	●	●	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	STAT3, Phospho-Y705 + STAT3, Total	●	○	○	○	○	○	○	○
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	●	●	●	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	STAT5, Phospho-Y694/699 + STAT5, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	STAT5, Total	●	○	○	○	○	○	○	○
HTRF™	STAT5, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	STAT6, Phospho-Y641	●	○	○	○	○	○	○	○
HTRF™	STAT6, Phospho-Y641	●	✗	○	○	○	○	○	○
Alpha SureFire® Ultra™ Multiplex	STAT6, Phospho-Y641 + STAT6, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	STAT6, Total	●	●	○	○	○	○	○	○
HTRF™	STAT6, Total	●	●	✗	✗	✗	✗	✗	✗
HTRF™	STING, Phospho-S365 (mouse)	✗	●	✗	✗	✗	✗	✗	✗
HTRF™	STING, Phospho-S366	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	STING, Total	●	○	○	○	○	○	○	○
HTRF™	STING, Total (mouse)	●	●	✗	✗	✗	✗	✗	✗
HTRF™	STING, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SYK, Phospho-Y352	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SYK, Phospho-Y525/526	●	●	○	○	○	○	○	○
HTRF™	SYK, Phospho-Y525	●	✗	✗	✗	✗	✗	✗	✗

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
Alpha SureFire® Ultra™ Multiplex	SYK, Phospho-Y525/526 + SYK, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SYK, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	SYK, Total (mouse)	✗	●	○	○	○	○	○	○
HTRF™	SYK, Total	●	✗	✗	✗	✗	✗	✗	✗
HTRF™	Tau, Phospho-S202/T205	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Phospho-S214	●	○	○	○	○	○	○	○
HTRF™	Tau, Phospho-S356	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Phospho-S396	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Phospho-S404	●	○	○	○	○	○	○	○
HTRF™	Tau, Phospho-S404	●	●	●	○	○	○	○	○
HTRF™	Tau, Phospho-S422	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Phospho-T181	●	○	○	○	○	○	○	○
HTRF™	Tau, Phospho-T181	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Phospho-T205	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Phospho-T217	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Tau, Total	●	○	○	○	○	○	○	○
HTRF™	Tau, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TAK1, Phospho-T184/187	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TAK1, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TBK1, Phospho-S172	●	●	○	○	○	○	○	○
HTRF™	TBK1, Phospho-S172	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TBK1, Total	●	●	○	○	○	○	○	○
HTRF™	TBK1, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TDP-43 aggregated	●	●	○	○	○	○	○	○
HTRF™	TDP-43 aggregated	●	●	●	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	TDP-43, Phospho-S409/410 (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TDP-43, Phospho-S409/410	●	✗	○	○	○	○	○	○
HTRF™	TDP-43, Phospho-S409/410	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TDP-43, Total (mouse)	✗	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TDP-43, Total	●	✗	○	○	○	○	○	○
HTRF™	TDP-43, Total	●	●	●	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFE3, Phospho-S321	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFE3, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFEB, Phospho-S122	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFEB, Phospho-S122 (mouse)	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFEB, Phospho-S211	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFEB, Phospho-S211 (mouse)	●	●	○	○	○	○	○	○
HTRF™	TFEB, Phospho-S211	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFEB, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TFEB, Total (mouse)	●	●	○	○	○	○	○	○
HTRF™	TFEB, Total	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TIE2, Phospho-Y992	●	○	○	○	○	○	○	○
HTRF™	TIE2, Phospho-Y992	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TIE2, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TRAF3, Total NEW	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TREM2, Aggregate	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TREM2/DAP12 Complex	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TREM2, Total	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TrkA, Phospho-Y674/675 / TrkB, Phospho-Y706/707	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	TYK2, Phospho-Y1054/1055	●	●	○	○	○	○	○	○
HTRF™	TYK2, Phospho-Y1054/1055	●	✗	✗	✗	✗	✗	✗	✗

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	TYK2, Total	●	✗	○	○	○	○	○	○
HTRF™	TYK2, Total	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	Ubiquitin, Phospho-S65	●	●	○	○	○	○	○	○
HTRF™	Ubiquitin, Phospho-S65	●	●	●	●	●	○	●	●
AlphaLISA™ SureFire® Ultra™	Ubiquitin, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ULK1, Phospho-S556	●	✗	○	○	○	○	○	○
HTRF™	ULK1, Phospho-S556	●	●	○	○	○	○	○	○
HTRF™	ULK1, Phospho-S757 mouse/S758 human	●	●	○	●	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ULK1, Phospho-S758	●	✗	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	ULK1, Total	●	✗	○	○	○	○	○	○
HTRF™	ULK1, Total	●	●	○	○	○	○	○	○
HTRF™	VASP, Phospho-S157	●	●	○	○	○	○	○	○
HTRF™	VASP, Phospho-S239	●	●	○	○	○	○	○	○
HTRF™	VASP, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	VAV1, Phospho-Y174	●	✗	○	○	○	○	○	○
HTRF™	VAV1, Phospho-Y174	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	VAV1, Total	●	✗	○	○	○	○	○	○
HTRF™	VAV1, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	VEGF Rec 2, Phospho-Y951	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	VEGF Rec 2, Phospho-Y1175	●	○	○	○	○	○	○	○
HTRF™	VEGFR2, Phospho-Y1175	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	VEGF Rec 2, Phospho-Y1214	●	○	○	○	○	○	○	○

		Human	Mouse	Rat	Monkey	Hamster	Drosophila	Pig	Dog
AlphaLISA™ SureFire® Ultra™	VEGF Rec 2, Total	●	○	○	○	○	○	○	○
HTRF™	VEGFR2, Total	●	○	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	WDR5, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	Wee1, Total	●	●	●	○	○	○	○	○
HTRF™	Wee1, Total	●	●	○	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	XBP1s, Total	●	✗	○	○	○	○	○	○
HTRF™	XBP1s, Total	●	✗	✗	○	○	○	○	○
HTRF™	XBP1, Total	●	✗	✗	○	○	○	○	○
AlphaLISA™ SureFire® Ultra™	YAP, Phospho-S127	●	●	○	○	○	○	○	○
HTRF™	YAP, Phospho-S127	●	●	●	○	●	○	○	○
AlphaLISA™ SureFire® Ultra™	YAP, Total	●	●	○	○	○	○	○	○
HTRF®	YAP, Total	●	●	○	●	●	○	○	○
AlphaLISA™ SureFire® Ultra™	ZAP-70, Phospho-Y319	●	○	○	○	○	○	○	○
HTRF®	ZAP-70, Phospho-Y319	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA™ SureFire® Ultra™	ZAP-70, Total	●	●	○	○	○	○	○	○
HTRF®	ZAP-70, Total	●	✗	✗	✗	✗	✗	✗	✗

