

Species compatibility

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	●	●	●	○	○	○	○	○
HTRF®	AKT1/2/3 phospho-S473	●	●	●	●	●	●	○	○
AlphaLISA® SureFire® Ultra™	AKT1/2/3 phospho-T308	●	●	●	○	○	○	○	○
HTRF®	AKT1/2/3 phospho-T308	●	●	●	●	●	●	○	○
AlphaLISA® SureFire® Ultra™	AKT1/2/3 total	●	●	○	○	○	○	○	○
HTRF®	AKT 1/2/3total	●	●	●	○	●	○	○	○
AlphaLISA® SureFire® Ultra™	AKT1 phospho-S473	●	●	●	○	○	○	○	○
HTRF®	AKT1 phospho-S473	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	AKT1 total	●	●	●	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	AKT1 total	●	●	●	●	○	○	○	○
HTRF®	AKT2 phospho-S473	●	●	●	●	○	○	○	○
HTRF®	AKT2 total	●	●	●	●	○	○	○	○
HTRF®	AKT3 phospho-S473	●	●	●	○	○	○	○	○
HTRF®	AKT3 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	●	●	●	○	○	○	○	○
HTRF®	AP2 total	●	●	●	○	●	○	○	○
HTRF®	ATF2 phospho-T71	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATG16L1 phospho-S278 NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATG16L1 total NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATM phospho-S198	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	ATM Total	●	●	●	○	○	○	○	○
HTRF®	BAD phospho-S112	●	●	●	●	○	○	○	○
HTRF®	B-arrestin 1 total	●	●	●	○	●	○	○	○
HTRF®	B-arrestin 2 total	●	●	●	○	●	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin phospho-S37 NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin phospho-T41/S45 NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin phospho-S675 NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	β-Catenin total NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Beclin-1 total	●	●	●	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	Beta Catenin phospho-T41/S33/37	●	●	●	●	○	○	●	○
HTRF®	Beta Catenin 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	●	●	○	○	○	○	○	○
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	○	○	○	○	○
HTRF®	Cereblon total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK1 phospho-S345	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK1 total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK2 phospho-T68	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	CHK2 total	●	○	○	○	○	○	○	○
HTRF®	cJun phospho-S63	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	c-Met phospho-Y1234/1235	●	○	○	○	○	○	○	○
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	●	●	●	○	○	○	○	○
HTRF®	cRAF phospho-S338	●	●	●	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	ERK1/2 phospho-T202/Y204	●	●	●	○	●	○	○	○
AlphaLISA® SureFire® Ultra™	ERK1/2 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	●	✗	○	○	○	○	○	○
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	●	○	○	○	○	○	○	○
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	●	●	●	●	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
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	●	●	●	○	○	○	○	○
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 NEW	●	●	○	○	○	○	○	○
HTRF®	Histone H3 phospho-T3	●	●	●	○	○	○	○	○
HTRF®	Histone H3 phospho-S10	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	Histone H3 total NEW	●	●	○	○	○	○	○	○
HTRF®	Histone H3 total	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	IGF-1 Rec.β phospho-Y1135/1136	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IGF-1 Rec.β Total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IKKα phospho-S176/180	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	IKKα Total	●	○	○	○	○	○	○	○
HTRF®	IKKB phospho-S177/181	●	●	●	●	○	○	○	○
AlphaLISA® SureFire® Ultra™	Insulin Rec. β phospho-Y1150/1151	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Insulin Rec. β Total	●	●	●	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	IRAK3 total	●	✗	○	○	○	○	○	○
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®	IRS1 phospho-S307 mouse/S312 human	●	●	○	○	○	○	○	○
HTRF®	IRS1 total	●	●	○	○	○	○	○	○
HTRF®	IRS2 total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	JAK2 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	●	●	●	○	○	○	○	○
HTRF®	LRRK2 phospho-S935	●	●	●	○	○	○	○	○
HTRF®	LRRK2 total	●	●	●	○	○	○	○	○
HTRF®	MDA5 total	●	○	○	○	○	○	○	○
HTRF®	MDA5 aggregation	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MEK1 phospho-S218/222	●	●	○	○	○	○	○	○
HTRF®	MEK1 phospho-S218/222	●	●	●	●	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
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™	p-MLKL (Ser345) (Mouse)	X	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	MLKL Total (Mouse)	●	●	○	○	○	○	○	○
HTRF®	MLKL phospho S358 NEW	●	○	○	○	○	○	○	○
HTRF®	MLKL total NEW	●	○	○	○	○	○	○	○
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	●	●	●	○	○	○	○	○
HTRF®	Phospholamban phospho-S16	●	●	●	○	○	○	○	○
HTRF®	Phospholamban phospho-T17	●	●	●	○	○	○	○	○
HTRF®	Phospholamban 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	●	○	○	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
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	●	●	●	○	○	○	○	○
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™	PLCy2 phospho-Y753	●	○	○	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	PLCy2 phospho-Y759	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCy2 phospho-Y1197	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCy2 phospho-Y1217	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	PLCy2 Total	●	○	○	○	○	○	○	○
HTRF®	PLCg2 total	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RAB10 phospho-T73 NEW	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RAB10 total NEW	●	●	●	○	○	○	○	○
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	●	●	○	○	○	○	○	○
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)	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	p-Ret (panY)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Ret Total	●	●	○	○	○	○	○	○
HTRF®	RET phospho pan Y NEW	●	●	●	○	○	○	○	○
HTRF®	RET total NEW	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RIP3 phospho-S227	●	✗	✗	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
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 (Ser240/244) (Eu) + Total RPS6 (Tb)	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1 phospho-S380	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1 phospho-T359/S363	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1 phospho-T573	●	○	○	○	○	○	○	○
HTRF®	RSK1 phospho S380 NEW	●	●	●	○	○	○	○	○
HTRF®	RSK1 phospho T359/S363 NEW	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	RSK1 total	●	●	●	○	○	○	○	○
HTRF®	RSK1 total NEW	●	●	●	○	○	○	○	○
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	●	○	○	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
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	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76 phospho-S376	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76 phospho-S376 (Mouse, Human) NEW	●	●	○	○	○	○	○	○
HTRF®	SLP-76 phospho-S376	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	SLP-76 phospho-S376 (mouse)	●	●	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76 phospho-Y113	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76 phospho-Y145	●	○	○	○	○	○	○	○
HTRF®	SLP-76 phospho-Y145	●	✗	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	SLP-76 total	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	SLP-76 total (Mouse, Human) NEW	●	●	○	○	○	○	○	○
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	●	●	○	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	SMAD3 total	●	●	●	○	○	○	○	○
HTRF®	SMAD3 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	●	●	✗	✗	✗	✗	✗	✗
AlphaLISA® SureFire® Ultra™	STAT1 total	●	○	○	○	○	○	○	○
HTRF®	STAT1 total	●	●	✗	✗	✗	✗	✗	✗
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	●	✗	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	STAT6 total	●	○	○	○	○	○	○	○
HTRF®	STING phospho-S366	●	✗	○	○	○	○	○	○

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
AlphaLISA® SureFire® Ultra™	STING total	●	○	○	○	○	○	○	○
HTRF®	STING (m) phospho-S365	✗	●	✗	✗	✗	✗	✗	✗
HTRF®	STING (m) Total	●	●	✗	✗	✗	✗	✗	✗
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	●	✗	✗	✗	✗	✗	✗	✗
HTRF®	Tau phospho-S202/T205	●	●	●	○	○	○	○	○
HTRF®	Tau phospho-S356	●	●	●	○	○	○	○	○
HTRF®	Tau phospho-S404	●	●	●	○	○	○	○	○
HTRF®	Tau phospho-S422	●	●	●	○	○	○	○	○
HTRF®	Tau phospho-T181	●	●	●	○	○	○	○	○
HTRF®	Tau total	●	✗	○	○	○	○	○	○
HTRF®	TDP-43 aggregated	●	●	●	○	○	○	○	○
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	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Ubiquitin phospho-S65	●	●	●	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	Ubiquitin Total	●	●	○	○	○	○	○	○
HTRF®	Ubiquitin phospho-S65	●	●	●	●	●	○	●	●

Species compatibility

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

		HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
HTRF®	ULK1 phosho-S757 mouse/S758 human	●	●	○	●	○	○	○	○
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 phospho-Y951	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	VEGFR2 total	●	○	○	○	○	○	○	○
AlphaLISA® SureFire® Ultra™	YAP phospho-S127	●	●	●	○	○	○	○	○
HTRF®	YAP phospho-S127	●	●	●	○	●	○	○	○
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	●	✗	✗	✗	✗	✗	✗	✗

FOR MORE INFORMATION

Europe and other countries +33(0)466-796-705 U.S. and Canada 1-888-963-4567 China +86-21-5018-9880 Japan +81-(0)43-306-8712

Visit www.cisbio.com to find a list of our regional distributors