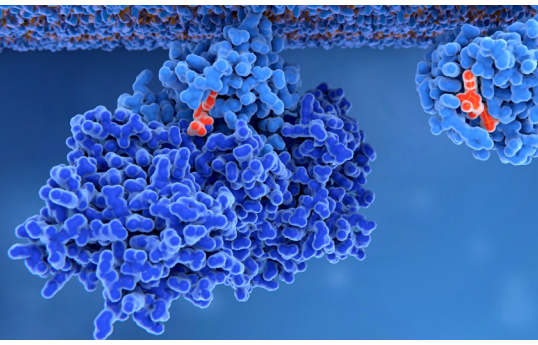


# Alpha SureFire Technology

Assays for phosphorylated targets



## The Ultimate Tools for Phosphoprotein Research

Alpha SureFire® assays are no-wash cellular kinase assays. While they exist in different formats (AlphaScreen® SureFire, AlphaLISA® SureFire Ultra, and Alpha SureFire Ultra Multiplex, which include the Terbium SureFire Ultra kits), all of them use PerkinElmer's exclusive bead-based technology and sandwich immunoassays for the detection of phosphorylated proteins in cells. Alpha SureFire assays are a quantitative alternative to Western Blotting, automation-friendly, easy to miniaturize, and detect both endogenous and recombinant proteins.

### AlphaScreen SureFire

The original format available since 2005, AlphaScreen SureFire has a rich track record of small molecule screening successes

- Uses protein A to capture one of the kit's antibodies
- 4-hour incubation protocol
- Need to order the AlphaScreen protein A kit in addition to the AlphaScreen SureFire kit

### AlphaLISA SureFire Ultra

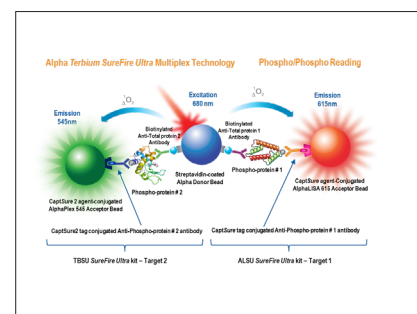
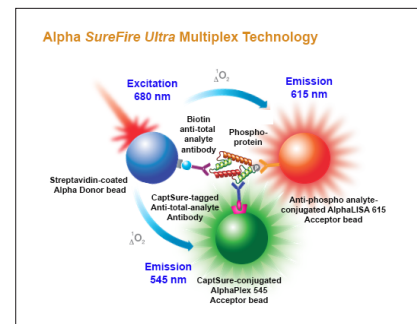
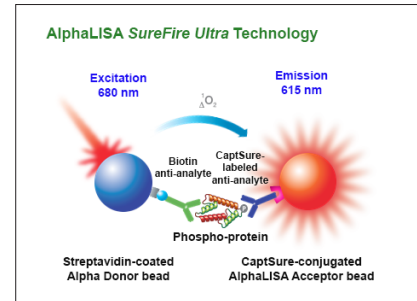
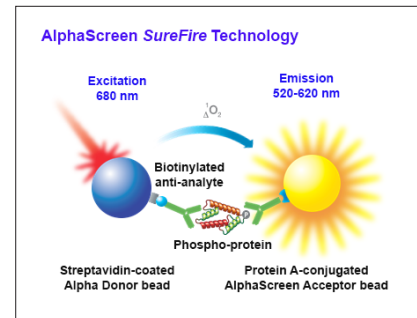
A significant improvement over the AlphaScreen SureFire format; one of the antibodies is labeled with CaptSure™, a small peptide tag, which is recognized by CaptSure-antibody-coated beads

- Lower background signal, improved sensitivity, and wider dynamic range
- No interference from extraneous antibodies; ideal for therapeutic antibody screening
- AlphaLISA beads are optimal for hemoglobin-containing samples
- 2-hour incubation protocol
- Beads are included in the kit, eliminating the need to purchase any additional components separately
- Cost-effective alternative to AlphaScreen SureFire

### Alpha SureFire Ultra Multiplex

This format utilizes AlphaPlex™ technology to read multiple signals from the same well

- Offers the option of either improved data quality with internal controls (total protein) or the ability to gather more data from a single experiment (multiplex phosphorylation events), both options results in significant time and cost savings
- Same features as the AlphaLISA SureFire Ultra, beads are included in the kit
- Dual-target kits include all the reagents needed to measure both total and phosphorylated target
- Terbium SureFire Ultra kits include reagents needed to be combined with those of an AlphaLISA SureFire Ultra kit to measure two phosphoprotein targets or one phosphoprotein target and total cofilin (as a control housekeeping protein)



# ORDERING INFORMATION

July 2018

Analyte	AlphaScreen SureFire Kits			AlphaLISA SureFire Ultra Kits				
	500 data points	10,000 data points	50,000 data points	500 data points	10,000 data points	50,000 data points	HV (100 data points)	Additional Control Lysate
Phospho-4EBP1 (Thr37/46)	TGR4E5500	TGR4E510K	TGR4E550K	ALSU-P4EBP-A500	ALSU-P4EBP-A10K	ALSU-P4EBP-A50K	ALSU-P4EBP-A-HV	ALSU-P4EBP-A-L
Phospho-4EBP1 (Thr70)	TGR4E25500	TGR4E2510K	TGR4E2550K					
<b>Phospho-ACC (Ser79)</b>				<b>ALSU-PACC-A500</b>	<b>ALSU-PACC-A10K</b>	<b>ALSU-PACC-A50K</b>	<b>ALSU-PACC-A-HV</b>	<b>ALSU-PACC-A-L</b>
Phospho-AKT1 (Thr308)	TGRA25500	TGRA2510K	TGRA2550K					
Phospho-AKT1 (Ser473)	TGRA55500	TGRA5510K	TGRA5550K					
Total AKT1	TGRTA15500	TGRTA1510K	TGRTA1550K	ALSU-TAKT1-A500	ALSU-TAKT1-A10K	ALSU-TAKT1-A50K	ALSU-TAKT1-A-HV	ALSU-TAKT1-A-L
Phospho-AKT1/2/3 (Thr308)	TGRA35500	TGRA3510K	TGRA3550K	ALSU-PAKT-A500	ALSU-PAKT-A10K	ALSU-PAKT-A50K	ALSU-PAKT-A-HV	ALSU-PAKT-A-L
Phospho-AKT1/2/3 (Ser473)	TGRA5500	TGRA510K	TGRA550K					
Phospho-AKT1/2/3 (Ser473)	TGRA45500	TGRA4510K	TGRA4550K	ALSU-PAKT-B500	ALSU-PAKT-B10K	ALSU-PAKT-B50K	ALSU-PAKT-B-HV	ALSU-PAKT-B-L
Total AKT1/2/3	TGRTAPS500	TGRTAPS10K	TGRTAPS50K					
Phospho-ALK (Tyr1586)	TGRALS500	TGRALS10K	TGRALS50K					
Phospho-ALK (Tyr1604)	TGRAL25500	TGRAL2510K	TGRAL2550K					
Total ALK	TGRTALS500	TGRTALS10K	TGRTALS50K					
<b>Phospho-AMPK<math>\alpha</math> 1/2 (Thr172)</b>				<b>ALSU-PAMPK-A500</b>	<b>ALSU-PAMPK-A10K</b>	<b>ALSU-PAMPK-A50K</b>	<b>ALSU-PAMPK-A-HV</b>	<b>ALSU-PAMPK-A-L</b>
Phospho-ATF-1 (Ser63)	TGRAF15500	TGRAF1510K	TGRAF1550K					
Phospho-ATF-2 (Thr71)	TGRAF25500	TGRAF2510K	TGRAF2550K					
Phospho-BAD (Ser112)	TGRB5500	TGRB510K	TGRB550K					
Phospho-BAD (Ser136)	TGRB25500	TGRB2510K	TGRB2550K					
Phospho- $\beta$ -Catenin (Ser45)	TGRBCS500	TGRBCS10K	TGRBCS50K					
Phospho-Caspase 9 (Ser196)	TGRC95500	TGRC9510K	TGRC9550K					
Phospho-Chk1 (Ser345)	TGRCHK15500	TGRCHK1510K	TGRCHK1550K					
Phospho-Chk2 (Thr68)	TGRCHK25500	TGRCHK2510K	TGRCHK2550K					
Phospho-c-Jun (Ser63)	TGRCJ5500	TGRCJ510K	TGRCJ550K					
Phospho-c-Jun (Ser73)	TGRCJ25500	TGRCJ2510K	TGRCJ2550K					
Phospho-c-Met (Tyr1234/1235)	TGRCMS500	TGRCMS10K	TGRCMS50K	<b>ALSU-PCMET-A500</b>	<b>ALSU-PCMET-A10K</b>	<b>ALSU-PCMET-A50K</b>	<b>ALSU-PCMET-A-HV</b>	<b>ALSU-PCMET-A-L</b>
Phospho-Cofilin (Ser3)	TGRCOFS500	TGRCOFS10K	TGRCOFS50K	<b>ALSU-PCOF-A500</b>	<b>ALSU-PCOF-A10K</b>	<b>ALSU-PCOF-A50K</b>	<b>ALSU-PCOF-A-HV</b>	<b>ALSU-PCOF-A-L</b>
<b>Total Cofilin</b>				<b>ALSU-TCOF-A500</b>	<b>ALSU-TCOF-A10K</b>	<b>ALSU-TCOF-A50K</b>	<b>ALSU-TCOF-A-HV</b>	<b>ALSU-TCOF-A-L</b>
Phospho-CREB-1 (Ser133)	TGRCBS500	TGRCBS10K	TGRCBS50K	ALSU-PCREB-A500	ALSU-PCREB-A10K	ALSU-PCREB-A50K	ALSU-PCREB-A-HV	ALSU-PCREB-A-L
Phospho-EGF Receptor (Tyr1068)	TGRERS500	TGRERS10K	TGRERS50K	ALSU-PEGFR-A500	ALSU-PEGFR-A10K	ALSU-PEGFR-A50K	ALSU-PEGFR-A-HV	ALSU-PEGFR-A-L
Phospho-eIF2 $\alpha$ (Ser51)				ALSU-PEIF2-B500	ALSU-PEIF2-B10K	ALSU-PEIF2-B50K	ALSU-PEIF2-B-HV	ALSU-PEIF2-B-L
Phospho-eIF4E (Ser209)	TGREIF45500	TGREIF4510K	TGREIF4550K	ALSU-PEIF4-A500	ALSU-PEIF4-A10K	ALSU-PEIF4-A50K	ALSU-PEIF4-A-HV	ALSU-PEIF4-A-L
Phospho-Elk-1 (Ser383)	TGRELKS500	TGRELKS10K	TGRELKS50K					
Phospho-ErbB2 (Tyr1221/1222)	TGREB25500	TGREB2510K	TGREB2550K	<b>ALSU-PEB2-A500</b>	<b>ALSU-PEB2-A10K</b>	<b>ALSU-PEB2-A50K</b>	<b>ALSU-PEB2-A-HV</b>	<b>ALSU-PEB2-A-L</b>
Phospho-ERK1/2 (T202/Y204)	TGRESB500	TGRESB10K	TGRESB50K	ALSU-PERK-A500	ALSU-PERK-A10K	ALSU-PERK-A50K	ALSU-PERK-A-HV	ALSU-PERK-A-L
Total ERK1/2	TGRTESB500	TGRTESB10K	TGRTESB50K	ALSU-TERK-A500	ALSU-TERK-A10K	ALSU-TERK-A50K	ALSU-TERK-A-HV	ALSU-TERK-A-L
Total GAPDH (human)	TGRGDS500	TGRGDS10K	TGRGDS50K	ALSU-TGAPD-A500	ALSU-TGAPD-A10K	ALSU-TGAPD-A50K	ALSU-TGAPD-A-HV	ALSU-TGAPD-A-L
Total GAPDH (human, mouse, rat)				ALSU-TGAPD-B500	ALSU-TGAPD-B10K	ALSU-TGAPD-B50K	ALSU-TGAPD-B-HV	ALSU-TGAPD-B-L
Phospho-GSK3 $\alpha$ (Ser21)	TGRGAS500	TGRGAS10K	TGRGAS50K					
Phospho-GSK3 $\beta$ (Ser9)	TGRGBS500	TGRGBS10K	TGRGBS50K	ALSU-PGS3B-A500	ALSU-PGS3B-A10K	ALSU-PGS3B-A50K	ALSU-PGS3B-A-HV	ALSU-PGS3B-A-L
Phospho-Histone H3 (Ser10)	TGRH35500	TGRH3510K	TGRH3550K					
Phospho-HSP27 (Ser82)	TGRHS275500	TGRHS27510K	TGRHS27550K					
Phospho-IGF1 Receptor $\beta$ (Tyr1135/1136)				ALSU-PIGFR-B500	ALSU-PIGFR-B10K	ALSU-PIGFR-B50K	ALSU-PIGFR-B-HV	ALSU-PIGFR-B-L
Phospho-I $\kappa$ B $\alpha$ (Ser32/36)	TGRIKS500	TGRIKS10K	TGRIKS50K					
Total I $\kappa$ B $\alpha$	TGRTIKS500	TGRTIKS10K	TGRTIKS50K					
Phospho-I $\kappa$ K $\alpha$ (Ser176/180)	TGRKAS500	TGRKAS10K	TGRKAS50K	ALSU-PIKKA-A500	ALSU-PIKKA-A10K	ALSU-PIKKA-A50K	ALSU-PIKKA-A-HV	ALSU-PIKKA-A-L
Phospho-I $\kappa$ K $\beta$ (Ser177/181)	TGRKBS500	TGRKBS10K	TGRKBS50K					
Phospho-Insulin Receptor $\beta$ (Tyr1150/1151)	TGRIRS500	TGRIRS10K	TGRIRS50K	ALSU-PINR-A500	ALSU-PINR-A10K	ALSU-PINR-A50K	ALSU-PINR-A-HV	ALSU-PINR-A-L
Phospho-JNK1/3 (Thr183/Tyr185)	TGRJ5500	TGRJ510K	TGRJ550K					
Phospho-JNK1/2/3 (Thr183/Tyr185)	TGRJP5500	TGRJP510K	TGRJP550K	ALSU-PJNK-A500	ALSU-PJNK-A10K	ALSU-PJNK-A50K	ALSU-PJNK-A-HV	ALSU-PJNK-A-L
Total JNK1/2/3	TGRTJ5500	TGRTJ510K	TGRTJ550K					
Phospho-MEK1 (Ser218/222)				ALSU-PMEK1-A500	ALSU-PMEK1-A10K	ALSU-PMEK1-A50K	ALSU-PMEK1-A-HV	ALSU-PMEK1-A-L
Total MEK1				ALSU-TMEK1-A500	ALSU-TMEK1-A10K	ALSU-TMEK1-A50K	ALSU-TMEK1-A-HV	ALSU-TMEK1-A-L
Phospho-MKK3/6 (Ser189/Ser207)	TGRMK35500	TGRMK3510K	TGRMK3550K					
Phospho-MKK4 (Ser257/Thr261)	TGRMK45500	TGRMK4510K	TGRMK4550K					
Phospho-mTOR (Ser2448)				ALSU-PMTOR-A500	ALSU-PMTOR-A10K	ALSU-PMTOR-A50K	ALSU-PMTOR-A-HV	ALSU-PMTOR-A-L
Total mTOR				ALSU-TMTOR-A500	ALSU-TMTOR-A10K	ALSU-TMTOR-A50K	ALSU-TMTOR-A-HV	ALSU-TMTOR-A-L
Phospho-NF $\kappa$ B p65 (Ser536)	TGRNFS500	TGRNFS10K	TGRNFS50K	ALSU-PNFKB-A500	ALSU-PNFKB-A10K	ALSU-PNFKB-A50K	ALSU-PNFKB-A-HV	ALSU-PNFKB-A-L
Total NF $\kappa$ B p65	TGRTNFS500	TGRTNFS10K	TGRTNFS50K					
Phospho-p38 MAPK (Thr180/Tyr182)				ALSU-PP38-B500	ALSU-PP38-B10K	ALSU-PP38-B50K	ALSU-PP38-B-HV	ALSU-PP38-B-L
Total p38 MAPK				ALSU-TP38-B500	ALSU-TP38-B10K	ALSU-TP38-B50K	ALSU-TP38-B-HV	ALSU-TP38-B-L
Phospho-p53 (Ser15)	TGRP535500	TGRP53510K	TGRP53550K					
Phospho-p53 (Ser392)	TGRP5325500	TGRP532510K	TGRP532550K					
Total p53	TGRT535500	TGRT53510K	TGRT53550K					
Phospho-p70S6K (Thr229)	TGR7035500	TGR703510K	TGR703550K					
Phospho-p70S6K (Thr389)	TGR705500	TGR70510K	TGR70550K	ALSU-PP70-A500	ALSU-PP70-A10K	ALSU-PP70-A50K	ALSU-PP70-A-HV	ALSU-PP70-A-L
Phospho-p70S6K (Thr421/Ser424)	TGR7025500	TGR702510K	TGR702550K					
Total p70S6K	TGRT705500	TGRT70510K	TGRT70550K					

Kits launched recently are highlighted in **bold and red**

Analyte	AlphaScreen SureFire Kits			AlphaLISA SureFire Ultra Kits				
	500 data points	10,000 data points	50,000 data points	500 data points	10,000 data points	50,000 data points	HV (100 data points)	Additional Control Lysate
Phospho-PDK1 (Ser241)	TGRPS500	TGRPS10K	TGRPS50K					
Phospho-S6RP (Ser235/236)	TGRS6PS500	TGRS6PS10K	TGRS6PS50K					
Phospho-RPS6 (Ser240/244)	TGRS6P2S500	TGRS6P2S10K	TGRS6P2S50K	ALSU-PS6R-A500	ALSU-PS6R-A10K	ALSU-PS6R-A50K	ALSU-PS6R-A-HV	ALSU-PS6R-A-L
Phospho-SLP76 (Ser376)				ALSU-PSLP-A500	ALSU-PSLP-A10K	ALSU-PSLP-A50K	ALSU-PSLP-A-HV	ALSU-PSLP-A-L
Total-SLP76				ALSU-TSLP-A500	ALSU-TSLP-A10K	ALSU-TSLP-A50K	ALSU-TSLP-A-HV	ALSU-TSLP-A-L
Phospho-SMAD1 (Ser463/465)	TGRSM1S500	TGRSM1S10K	TGRSM1S50K	ALSU-PSM1-A500	ALSU-PSM1-A10K	ALSU-PSM1-A50K	ALSU-PSM1-A-HV	ALSU-PSM1-A-L
Phospho-SMAD2 (Ser465/467)	TGRSM2S500	TGRSM2S10K	TGRSM2S50K					
Phospho-SMAD3 (Ser423/425)	TGRSM3S500	TGRSM3S10K	TGRSM3S50K	ALSU-PSM3-A500	ALSU-PSM3-A10K	ALSU-PSM3-A50K	ALSU-PSM3-A-HV	ALSU-PSM3-A-L
Phospho-STAT1 (Tyr701)	TGRS1S500	TGRS1S10K	TGRS1S50K	ALSU-PST1-A500	ALSU-PST1-A10K	ALSU-PST1-A50K	ALSU-PST1-A-HV	ALSU-PST1-A-L
Phospho-STAT1 (Ser727)				ALSU-PST1-B500	ALSU-PST1-B10K	ALSU-PST1-B50K	ALSU-PST1-B-HV	ALSU-PST1-B-L
Phospho-STAT3 (Tyr705)	TGRS3S500	TGRS3S10K	TGRS3S50K	ALSU-PST3-A500	ALSU-PST3-A10K	ALSU-PST3-A50K	ALSU-PST3-A-HV	ALSU-PST3-A-L
Total STAT3				ALSU-TST3-A500	ALSU-TST3-A10K	ALSU-TST3-A50K	ALSU-TST3-A-HV	ALSU-TST3-A-L
Phospho-STAT4 (Tyr693)				ALSU-PST4-A500	ALSU-PST4-A10K	ALSU-PST4-A50K	ALSU-PST4-A-HV	ALSU-PST4-A-L
Phospho-STAT5 (Tyr694/699)	TGRS5S500	TGRS5S10K	TGRS5S50K	ALSU-PST5-B500	ALSU-PST5-B10K	ALSU-PST5-B50K	ALSU-PST5-B-HV	ALSU-PST5-B-L
Total STAT5				ALSU-TST5-A500	ALSU-TST5-A10K	ALSU-TST5-A50K	ALSU-TST5-A-HV	ALSU-TST5-A-L
Phospho-STAT6 (Tyr641)				ALSU-PST6-A500	ALSU-PST6-A10K	ALSU-PST6-A50K	ALSU-PST6-A-HV	ALSU-PST6-A-L
Phospho-Syk (Tyr525/526)				ALSU-PSYK-A500	ALSU-PSYK-A10K	ALSU-PSYK-A50K	ALSU-PSYK-A-HV	ALSU-PSYK-A-L
Total Syk				ALSU-TSYK-A500	ALSU-TSYK-A10K	ALSU-TSYK-A50K	ALSU-TSYK-A-HV	ALSU-TSYK-A-L
Phospho-Topoisomerase IIa (Ser1106)	TGRT2A2S500	TGRT2A2S10K	TGRT2A2S50K					
Phospho-Topoisomerase IIa (Thr1343)	TGRT2AS500	TGRT2AS10K	TGRT2AS50K					
Total Topoisomerase IIa	TGRT2AS500	TGRT2AS10K	TGRT2AS50K					
<b>Phospho-TrkA (Tyr674/675) / B (Tyr 706/707)</b>				<b>ALSU-PTRKAB-A500</b>	<b>ALSU-PTRKAB-A10K</b>	<b>ALSU-PTRKAB-A50K</b>	<b>ALSU-PTRKAB-A-HV</b>	<b>ALSU-PTRKAB-A-L</b>
Phospho-VEGF Receptor 2 (Tyr1175)	TGRVRS500	TGRVRS10K	TGRVRS50K	ALSU-PVGFR-A500	ALSU-PVGFR-A10K	ALSU-PVGFR-A50K	ALSU-PVGFR-A-HV	ALSU-PVGFR-A-L
<b>Total VEGF Receptor 2</b>				<b>ALSU-TVGFR-A500</b>	<b>ALSU-TVGFR-A10K</b>	<b>ALSU-TVGFR-A50K</b>	<b>ALSU-TVGFR-A-HV</b>	<b>ALSU-TVGFR-A-L</b>

Kits launched recently are highlighted in **bold and red**

Analyte	Alpha SureFire Ultra Multiplex Kits Total/Phospho Dual Kits			
	500 data points	10,000 data points	50,000 data points	HV (100 data points)
Phospho-AKT 1/2/3 (Ser473) (Eu) + Total AKT1 (Tb)	MPSU-PTAKT-K500	MPSU-PTAKT-K10K	MPSU-PTAKT-K50K	MPSU-PTAKT-K-HV
Phospho-ERK 1/2 (Eu) + Total ERK (Tb)	MPSU-PTERK-K500	MPSU-PTERK-K10K	MPSU-PTERK-K50K	MPSU-PTERK-K-HV

Analyte	Terbium SureFire Ultra Kits			
	500 data points	10,000 data points	50,000 data points	HV (100 data points)
Phospho-4EBP1 (Thr37/46) (Tb)	TBSU-P4EBP-A500	TBSU-P4EBP-A10K	TBSU-P4EBP-A50K	TBSU-P4EBP-A-HV
Phospho-Akt (Thr308) (Tb)	TBSU-PAKT-A500	TBSU-PAKT-A10K	TBSU-PAKT-A50K	TBSU-PAKT-A-HV
Phospho-Akt (Ser473) (Tb)	TBSU-PAKT-B500	TBSU-PAKT-B10K	TBSU-PAKT-B50K	TBSU-PAKT-B-HV
Phospho-CREB-1 (Ser133) (Tb)	TBSU-PCREB-A500	TBSU-PCREB-A10K	TBSU-PCREB-A50K	TBSU-PCREB-A-HV
Phospho-EGF Receptor (Tyr1068) (Tb)	TBSU-PEGFR-A500	TBSU-PEGFR-A10K	TBSU-PEGFR-A50K	TBSU-PEGFR-A-HV
Phospho-eIF2a (Ser51) (Tb)	TBSU-PEIF2-B500	TBSU-PEIF2-B10K	TBSU-PEIF2-B50K	TBSU-PEIF2-B-HV
Phospho-eIF4E (Ser209) (Tb)	TBSU-PEIF4-A500	TBSU-PEIF4-A10K	TBSU-PEIF4-A50K	TBSU-PEIF4-A-HV
Phospho-ERK1/2 (Thr202/Tyr204) (Tb)	TBSU-PERK-A500	TBSU-PERK-A10K	TBSU-PERK-A50K	TBSU-PERK-A-HV
Phospho-GSK3β (Ser9) (Tb)	TBSU-PGS3B-A500	TBSU-PGS3B-A10K	TBSU-PGS3B-A50K	TBSU-PGS3B-A-HV
Phospho-IGF1 Receptor β (Tyr1135/1136) (Tb)	TBSU-PIGFR-B500	TBSU-PIGFR-B10K	TBSU-PIGFR-B50K	TBSU-PIGFR-B-HV
Phospho-IKKa (Ser176/180) (Tb)	TBSU-PIKKA-A500	TBSU-PIKKA-A10K	TBSU-PIKKA-A50K	TBSU-PIKKA-A-HV
Phospho-Insulin Receptor β (Tyr1150/1151) (Tb)	TBSU-PINR-A500	TBSU-PINR-A10K	TBSU-PINR-A50K	TBSU-PINR-A-HV
Phospho-JNK1/2/3 (Thr183/Tyr185) (Tb)	TBSU-PJNK-A500	TBSU-PJNK-A10K	TBSU-PJNK-A50K	TBSU-PJNK-A-HV
Phospho-MEK-1 (Ser218/222) (Tb)	TBSU-PMEK1-A500	TBSU-PMEK1-A10K	TBSU-PMEK1-A50K	TBSU-PMEK1-A-HV
Phospho-mTOR (Ser2448) (Tb)	TBSU-PMTOR-A500	TBSU-PMTOR-A10K	TBSU-PMTOR-A50K	TBSU-PMTOR-A-HV
Phospho-NFκB p65 (Ser536) (Tb)	TBSU-PNFKB-A500	TBSU-PNFKB-A10K	TBSU-PNFKB-A50K	TBSU-PNFKB-A-HV
Phospho-p38 MAPK (Thr180/Tyr182) (Tb)	TBSU-PP38-B500	TBSU-PP38-B10K	TBSU-PP38-B50K	TBSU-PP38-B-HV
Phospho-p70S6K (Thr389) (Tb)	TBSU-PP70-A500	TBSU-PP70-A10K	TBSU-PP70-A50K	TBSU-PP70-A-HV
Phospho-S6 RP (Ser240/244) (Tb)	TBSU-PS6R-A500	TBSU-PS6R-A10K	TBSU-PS6R-A50K	TBSU-PS6R-A-HV
Phospho-SLP-76 (Ser376) (Tb)	TBSU-PSLP-A500	TBSU-PSLP-A10K	TBSU-PSLP-A50K	TBSU-PSLP-A-HV
Phospho-SMAD1 (Ser463/465) (Tb)	TBSU-PSM1-A500	TBSU-PSM1-A10K	TBSU-PSM1-A50K	TBSU-PSM1-A-HV
Phospho-SMAD3 (Ser423/425) (Tb)	TBSU-PSM3-A500	TBSU-PSM3-A10K	TBSU-PSM3-A50K	TBSU-PSM3-A-HV
Phospho-STAT1 (Tyr701) (Tb)	TBSU-PST1-A500	TBSU-PST1-A10K	TBSU-PST1-A50K	TBSU-PST1-A-HV
Phospho-STAT1 (Ser727) (Tb)	TBSU-PST1-B500	TBSU-PST1-B10K	TBSU-PST1-B50K	TBSU-PST1-B-HV
Phospho-STAT3 (Tyr705) (Tb)	TBSU-PST3-A500	TBSU-PST3-A10K	TBSU-PST3-A50K	TBSU-PST3-A-HV
Phospho-STAT4 (Tyr693) (Tb)	TBSU-PST4-A500	TBSU-PST4-A10K	TBSU-PST4-A50K	TBSU-PST4-A-HV
Phospho-STAT5 (Tyr694/699) (Tb)	TBSU-PST5-B500	TBSU-PST5-B10K	TBSU-PST5-B50K	TBSU-PST5-B-HV
Phospho-STAT6 (Tyr641) (Tb)	TBSU-PST6-A500	TBSU-PST6-A10K	TBSU-PST6-A50K	TBSU-PST6-A-HV
Phospho-SYK (Tyr525/526) (Tb)	TBSU-PSYK-A500	TBSU-PSYK-A10K	TBSU-PSYK-A50K	TBSU-PSYK-A-HV
Phospho-VEGF Receptor 2 (Tyr1175) (Tb)	TBSU-PVGFR-A500	TBSU-PVGFR-A10K	TBSU-PVGFR-A50K	TBSU-PVGFR-A-HV
Total Cofilin (Tb)	TBSU-TCOF-A500	TBSU-TCOF-A10K	TBSU-TCOF-A50K	TBSU-TCOF-A-HV

# ORDERING INFORMATION

## Alpha SureFire Ultra Multiplex Supplementary Beads

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

Beads	500 data points (60 µL)	10,000 data points (1.2 mL)	50,000 data points (6 mL)
Alpha 545 (Tb) SureFire Ultra CaptSure Acceptor Beads	MPSU-ACAT-0.06mL	MPSU-ACAT-1.2mL	MPSU-ACAT-6mL

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

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

Beads	500 data points	10,000 data points	50,000 data points	HV (100 data points)
AlphaScreen General IgG (Protein A based) Detection Kit (384-well format)	6760617C	6760617M	6760617R	
AlphaScreen SureFire HV IgG (Protein A) Detection Kit (96-well format)				6760617HV

## AlphaScreen SureFire Supplementary Buffers

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

Buffers	10 mL	100 mL
SureFire Lysis Buffer (5X)	TGRLB10ML	TGRLB100ML
SureFire Activation Buffer (50X)	TGRAB10ML	TGRAB100ML
SureFire Activation Buffer B (50X)	TGRABB10ML	TGRABB100ML
SureFire Activation Buffer C (50X)	TGRABC10ML	TGRABC100ML
SureFire Dilution Buffer (1X)	TGRDB10ML	TGRDB100ML

## AlphaScreen SureFire Supplementary Control Lysates

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

## AlphaLISA SureFire Ultra Supplementary Beads

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

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

## AlphaLISA SureFire Ultra Supplementary Buffers

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

Buffers	10 mL	100 mL
SureFire Ultra Lysis Buffer (5X)	ALSU-LB-10mL	ALSU-LB-100mL
SureFire Ultra Activation Buffer (50X)	ALSU-AB-10mL	ALSU-AB-100mL
SureFire Ultra Activation Buffer B (50X)	ALSU-ABB-10mL	ALSU-ABB-100mL
SureFire Ultra Activation Buffer C (50X)	ALSU-ABC-10mL	ALSU-ABC-100mL
SureFire Ultra Dilution Buffer (1X)	ALSU-DB-10mL	ALSU-DB-100mL

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2016-2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.