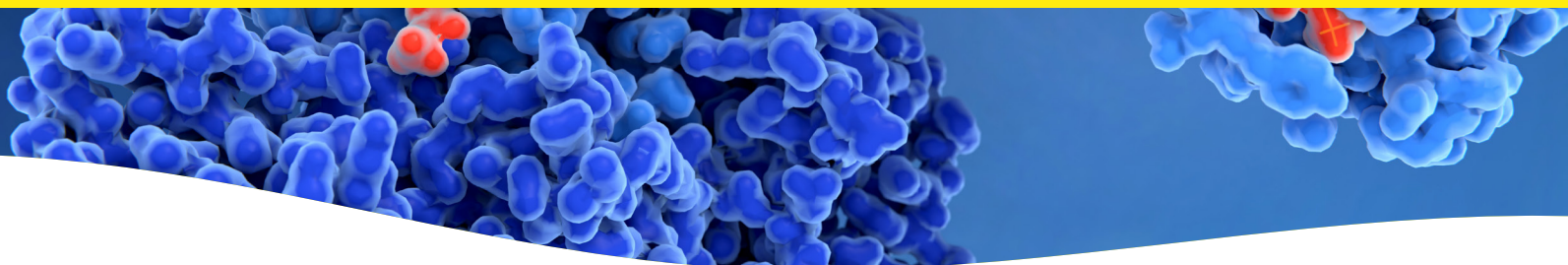




## Alpha SureFire technology assays for phosphorylated targets.



Exclusive Supplier of SureFire™ Technology to Revvity.

### The ultimate tools for phosphoprotein research

Alpha SureFire™ assays are no-wash cellular kinase assays that use Revvity's exclusive bead-based technology and sandwich immunoassays for the detection of phosphorylated proteins in cells. A quantitative alternative to Western Blotting, Alpha SureFire assays are automation-friendly, easy to miniaturize, and detect both endogenous and recombinant proteins.

#### AlphaLISA SureFire Ultra™ (ALSU)

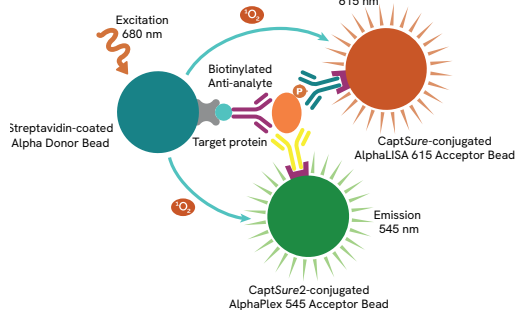
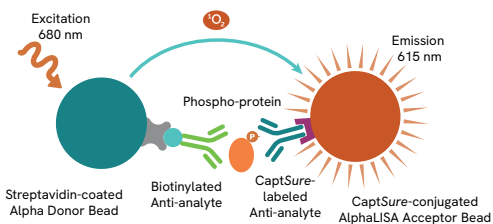
For single target analysis.

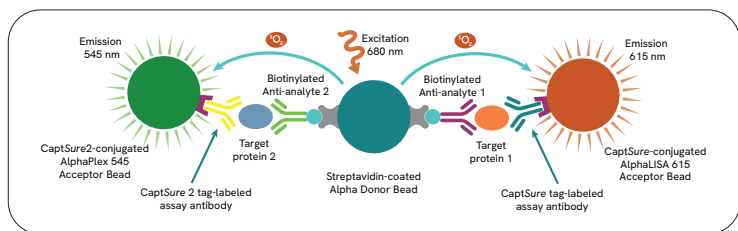
- Lower background signal, improved sensitivity, and wider dynamic range than traditional phospho-protein detection
- No interference from extraneous antibodies; ideal for therapeutic antibody screening
- Kits include all reagents

#### Alpha SureFire Ultra Multiplex (MPSU)

For phospho/total target analysis.

- 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 result in significant time and cost savings
- Uses same assay plate well from a single sample
- Dual-target kits include all the reagents needed to measure both total and phosphorylated target





## AlphaLISA SureFire Ultra Terbium Multiplex (TBSU)

For dual phospho/phospho target analysis.

- Kits include reagents needed to be combined with those of an AlphaLISA SureFire Ultra kit to measure two different phosphoproteins
- Total kit available for normalization of AlphaLISA SureFire Ultra phosphorylated targets

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
4E-BP1, Phospho-T37/46	ALSU-P4EBP-A500	ALSU-P4EBP-A10K	ALSU-P4EBP-A50K	ALSU-P4EBP-A-HV
4EBP1, Total	ALSU-T4EBP-A500	ALSU-T4EBP-A10K	ALSU-T4EBP-A50K	ALSU-T4EBP-A-HV
α-Synuclein, Aggregate Specific	ALSU-AASYN-A500	ALSU-AASYN-A10K	ALSU-AASYN-A50K	ALSU-AASYN-A-HV
α-Synuclein, Phospho-S129	ALSU-PASYN-B500	ALSU-PASYN-B10K	ALSU-PASYN-B50K	ALSU-PASYN-B-HV
α-Synuclein, Total	ALSU-TASYN-B500	ALSU-TASYN-B10K	ALSU-TASYN-B50K	ALSU-TASYN-B-HV
β-Catenin, Phospho-S37	ALSU-PBCAT-A500	ALSU-PBCAT-A10K	ALSU-PBCAT-A50K	ALSU-PBCAT-A-HV
β-Catenin, Phospho-T41/S45	ALSU-PBCAT-B500	ALSU-PBCAT-B10K	ALSU-PBCAT-B50K	ALSU-PBCAT-B-HV
β-Catenin, Phospho-S675	ALSU-PBCAT-C500	ALSU-PBCAT-C10K	ALSU-PBCAT-C50K	ALSU-PBCAT-C-HV
β-Catenin, Total	ALSU-TBCAT-A500	ALSU-TBCAT-A10K	ALSU-TBCAT-A50K	ALSU-TBCAT-A-HV
ACC, Phospho-S79	ALSU-PACC-A500	ALSU-PACC-A10K	ALSU-PACC-A50K	ALSU-PACC-A-HV
ACC, Total	ALSU-TACC-A500	ALSU-TACC-A10K	ALSU-TACC-A50K	ALSU-TACC-A-HV
<b>AIOLOS, Total</b>	<b>ALSU-TAIOL-A500</b>	<b>ALSU-TAIOL-A10K</b>	<b>ALSU-TAIOL-A50K</b>	<b>ALSU-TAIOL-A-HV</b>
AKT1, Phospho-S129	ALSU-PAKT1-D500	ALSU-PAKT1-D10K	ALSU-PAKT1-D50K	ALSU-PAKT1-D-HV
AKT1, Phospho-S473	ALSU-PAKT1-C500	ALSU-PAKT1-C10K	ALSU-PAKT1-C50K	ALSU-PAKT1-C-HV
AKT1, Total	ALSU-TAKT1-B500	ALSU-TAKT1-B10K	ALSU-TAKT1-B50K	ALSU-TAKT1-B-HV
AKT1/2/3, Phospho-S473	ALSU-PAKT-B500	ALSU-PAKT-B10K	ALSU-PAKT-B50K	ALSU-PAKT-B-HV
AKT 1/2/3, Phospho-T450	ALSU-PAKT-C500	ALSU-PAKT-C10K	ALSU-PAKT-C50K	ALSU-PAKT-C-HV
AKT1/2/3, Phospho-T308	ALSU-PAKT-A500	ALSU-PAKT-A10K	ALSU-PAKT-A50K	ALSU-PAKT-A-HV
AKT1/2/3, Total	ALSU-TAKT-B500	ALSU-TAKT-B10K	ALSU-TAKT-B50K	ALSU-TAKT-B-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
ALK, Phospho-Y1586	ALSU-PALK-A500	ALSU-PALK-A10K	ALSU-PALK-A50K	ALSU-PALK-A-HV
ALK, Phospho-Y1604	ALSU-PALK-C500	ALSU-PALK-C10K	ALSU-PALK-C50K	ALSU-PALK-C-HV
ALK, Total	ALSU-TALK-A500	ALSU-TALK-A10K	ALSU-TALK-A50K	ALSU-TALK-A-HV
AMPK $\alpha$ 1/2, Phospho-T172	ALSU-PAMPK-A500	ALSU-PAMPK-A10K	ALSU-PAMPK-A50K	ALSU-PAMPK-A-HV
AMPK $\alpha$ 1/2, Total	ALSU-TAMPK-A500	ALSU-TAMPK-A10K	ALSU-TAMPK-A50K	ALSU-TAMPK-A-HV
Androgen Receptor, Total	ALSU-TANDR-A500	ALSU-TANDR-A10K	ALSU-TANDR-A50K	ALSU-TANDR-A-HV
<b>ATF-2, Phospho-T71</b>	<b>ALSU-PATF2-A500</b>	<b>ALSU-PATF2-A10K</b>	<b>ALSU-PATF2-A50K</b>	<b>ALSU-PATF2-A-HV</b>
<b>ATF-2, Total</b>	<b>ALSU-TATF2-A500</b>	<b>ALSU-TATF2-A10K</b>	<b>ALSU-TATF2-A50K</b>	<b>ALSU-TATF2-A-HV</b>
ATF-4, Total	ALSU-TATF4-A500	ALSU-TATF4-A10K	ALSU-TATF4-A50K	ALSU-TATF4-A-HV
<b>ATG14, Phospho S29</b>	<b>ALSU-PATG14-A500</b>	<b>ALSU-PATG14-A10K</b>	<b>ALSU-PATG14-A50K</b>	<b>ALSU-PATG14-A-HV</b>
ATG16L1, Phospho-S278	ALSU-PATG-A500	ALSU-PATG-A10K	ALSU-PATG-A50K	ALSU-PATG-A-HV
ATG16L1, Total	ALSU-TATG-A500	ALSU-TATG-A10K	ALSU-TATG-A50K	ALSU-TATG-A-HV
ATM, Phospho-S1981	ALSU-PATM-A500	ALSU-PATM-A10K	ALSU-PATM-A50K	ALSU-PATM-A-HV
ATM, Total	ALSU-TATM-A500	ALSU-TATM-A10K	ALSU-TATM-A50K	ALSU-TATM-A-HV
Aurora A, Phospho-T288	ALSU-PAURA-A500	ALSU-PAURA-A10K	ALSU-PAURA-A50K	ALSU-PAURA-A-HV
Aurora A, Total	ALSU-TAURA-A500	ALSU-TAURA-A10K	ALSU-TAURA-A50K	ALSU-TAURA-A-HV
Aurora B, Phospho-T232	ALSU-PAURB-A500	ALSU-PAURB-A10K	ALSU-PAURB-A50K	ALSU-PAURB-A-HV
Aurora B, Total	ALSU-TAURB-A500	ALSU-TAURB-A10K	ALSU-TAURB-A50K	ALSU-TAURB-A-HV
BCL-2, Phospho-S70	ALSU-PBCL2-A500	ALSU-PBCL2-A10K	ALSU-PBCL2-A50K	ALSU-PBCL2-A-HV
BCL-2, Total	ALSU-TBCL2-A500	ALSU-TBCL2-A10K	ALSU-TBCL2-A50K	ALSU-TBCL2-A-HV
BCL-XL	ALSU-TBCLXL-A500	ALSU-TBCLXL-A10K	ALSU-TBCLXL-A50K	ALSU-TBCLXL-A-HV
Beclin-1, Total	ALSU-TBEC-A500	ALSU-TBEC-A10K	ALSU-TBEC-A50K	ALSU-TBEC-A-HV
BRD2, Phospho-S37	ALSU-PBRD2-A500	ALSU-PBRD2-A10K	ALSU-PBRD2-A50K	ALSU-PBRD2-A-HV
BRD2, Total	ALSU-TBRD2-A500	ALSU-TBRD2-A10K	ALSU-TBRD2-A50K	ALSU-TBRD2-A-HV
BRD4, Total	ALSU-TBRD4-A500	ALSU-TBRD4-A10K	ALSU-TBRD4-A50K	ALSU-TBRD4-A-HV
BRD3, Total	ALSU-TBRD3-A500	ALSU-TBRD3-A10K	ALSU-TBRD3-A50K	ALSU-TBRD3-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
BRD7, Total	ALSU-TBRD7-A500	ALSU-TBRD7-A10K	ALSU-TBRD7-A50K	ALSU-TBRD7-A-HV
BRD9, Total	ALSU-TBRD9-A500	ALSU-TBRD9-A10K	ALSU-TBRD9-A50K	ALSU-TBRD9-A-HV
Btk, Phospho-Y223	ALSU-PBTK-A500	ALSU-PBTK-A10K	ALSU-PBTK-A50K	ALSU-PBTK-A-HV
Btk, Phospho-Y551	ALSU-PBTK-B500	ALSU-PBTK-B10K	ALSU-PBTK-B50K	ALSU-PBTK-B-HV
Btk, Total	ALSU-TBTK-A500	ALSU-TBTK-A10K	ALSU-TBTK-A50K	ALSU-TBTK-A-HV
CaMKII, Phospho-T286	ALSU-PCAM-A500	ALSU-PCAM-A10K	ALSU-PCAM-A50K	ALSU-PCAM-A-HV
CaMKII, Total	ALSU-TCAM-A500	ALSU-TCAM-A10K	ALSU-TCAM-A50K	ALSU-TCAM-A-HV
CD3 Zeta, Phospho-Y142	ALSU-PCD3-B500	ALSU-PCD3-B10K	ALSU-PCD3-B50K	ALSU-PCD3-B-HV
CD3 Zeta, Phospho-Y83	ALSU-PCD3-A500	ALSU-PCD3-A10K	ALSU-PCD3-A50K	ALSU-PCD3-A-HV
CD3 Zeta, Total	ALSU-TCD3-A500	ALSU-TCD3-A10K	ALSU-TCD3-A50K	ALSU-TCD3-A-HV
CDC25C, Phospho-S216	ALSU-PCDC25-A500	ALSU-PCDC25-A10K	ALSU-PCDC25-A50K	ALSU-PCDC25-A-HV
CDC25C, Total	ALSU-TCDC25-A500	ALSU-TCDC25-A10K	ALSU-TCDC25-A50K	ALSU-TCDC25-A-HV
CDC6, Phospho-S54	ALSU-PCDC6-A500	ALSU-PCDC6-A10K	ALSU-PCDC6-A50K	ALSU-PCDC6-A-HV
CDC6, Phospho-S106	ALSU-PCDC6-B500	ALSU-PCDC6-B10K	ALSU-PCDC6-B50K	ALSU-PCDC6-B-HV
CDC6, Total	ALSU-TCDC6-A500	ALSU-TCDC6-A10K	ALSU-TCDC6-A50K	ALSU-TCDC6-A-HV
CDK1, Phospho-T14	ALSU-PCDK1-A500	ALSU-PCDK1-A10K	ALSU-PCDK1-A50K	ALSU-PCDK1-A-HV
CDK1, Phospho-Y15	ALSU-PCDK1-B500	ALSU-PCDK1-B10K	ALSU-PCDK1-B50K	ALSU-PCDK1-B-HV
CDK1, Phospho-T161	ALSU-PCDK1-C500	ALSU-PCDK1-C10K	ALSU-PCDK1-C50K	ALSU-PCDK1-C-HV
CDK1, Total	ALSU-TCDK1-A500	ALSU-TCDK1-A10K	ALSU-TCDK1-A50K	ALSU-TCDK1-A-HV
CDK2, Phospho-T160	ALSU-PCDK2-A500	ALSU-PCDK2-A10K	ALSU-PCDK2-A50K	ALSU-PCDK2-A-HV
CDK2, Total	ALSU-TCDK2-A500	ALSU-TCDK2-A10K	ALSU-TCDK2-A50K	ALSU-TCDK2-A-HV
CDK4, Total	ALSU-TCDK4-A500	ALSU-TCDK4-A10K	ALSU-TCDK4-A50K	ALSU-TCDK4-A-HV
CDK5, Total	ALSU-TCDK5-A500	ALSU-TCDK5-A10K	ALSU-TCDK5-A50K	ALSU-TCDK5-A-HV
CDK6, Total	ALSU-TCDK6-A500	ALSU-TCDK6-A10K	ALSU-TCDK6-A50K	ALSU-TCDK6-A-HV
CDK7, Total	ALSU-TCDK7-A500	ALSU-TCDK7-A10K	ALSU-TCDK7-A50K	ALSU-TCDK7-A-HV
<b>CDK8, Total</b>	<b>ALSU-TCDK8-A500</b>	<b>ALSU-TCDK8-A10K</b>	<b>ALSU-TCDK8-A50K</b>	<b>ALSU-TCDK8-A-HV</b>

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
CDK9, Total	ALSU-TCDK9-A500	ALSU-TCDK9-A10K	ALSU-TCDK9-A50K	ALSU-TCDK9-A-HV
CDK12, Total	ALSU-TCDK12-A500	ALSU-TCDK12-A10K	ALSU-TCDK12-A50K	ALSU-TCDK12-A-HV
CHK1, Phospho-S345	ALSU-PCHK1-A500	ALSU-PCHK1-A10K	ALSU-PCHK1-A50K	ALSU-PCHK1-A-HV
CHK1, Total	ALSU-TCHK1-A500	ALSU-TCHK1-A10K	ALSU-TCHK1-A50K	ALSU-TCHK1-A-HV
CHK2, Phospho-T68	ALSU-PCHK2-A500	ALSU-PCHK2-A10K	ALSU-PCHK2-A50K	ALSU-PCHK2-A-HV
CHK2, Total	ALSU-TCHK2-A500	ALSU-TCHK2-A10K	ALSU-TCHK2-A50K	ALSU-TCHK2-A-HV
<b>c-FOS, Phospho-S32</b>	<b>ALSU-PCFOS-A500</b>	<b>ALSU-PCFOS-A10K</b>	<b>ALSU-PCFOS-A50K</b>	<b>ALSU-PCFOS-A-HV</b>
<b>c-FOS, Total</b>	<b>ALSU-TCFOS-A500</b>	<b>ALSU-TCFOS-A10K</b>	<b>ALSU-TCFOS-A50K</b>	<b>ALSU-TCFOS-A-HV</b>
c-Jun, Phospho-S63	ALSU-PCJUN-A500	ALSU-PCJUN-A10K	ALSU-PCJUN-A50K	ALSU-PCJUN-A-HV
c-Jun, Phospho-S73	ALSU-PCJUN-B500	ALSU-PCJUN-B10K	ALSU-PCJUN-B50K	ALSU-PCJUN-B-HV
c-Jun, Total	ALSU-TCJUN-A500	ALSU-TCJUN-A10K	ALSU-TCJUN-A50K	ALSU-TCJUN-A-HV
c-Met, Phospho-Y1234/1235	ALSU-PCMET-A500	ALSU-PCMET-A10K	ALSU-PCMET-A50K	ALSU-PCMET-A-HV
c-Met, Total	ALSU-TCMET-A500	ALSU-TCMET-A10K	ALSU-TCMET-A50K	ALSU-TCMET-A-HV
c-Myb, Phospho-S11	ALSU-PCMYB-A500	ALSU-PCMYB-A10K	ALSU-PCMYB-A50K	ALSU-PCMYB-A-HV
c-Myb, Total	ALSU-TCMYB-A500	ALSU-TCMYB-A10K	ALSU-TCMYB-A50K	ALSU-TCMYB-A-HV
c-Myc, Phospho-S62	ALSU-PCMYC-A500	ALSU-PCMYC-A10K	ALSU-PCMYC-A50K	ALSU-PCMYC-A-HV
c-Myc, Phospho-T58	ALSU-PCMYC-B500	ALSU-PCMYC-B10K	ALSU-PCMYC-B50K	ALSU-PCMYC-B-HV
c-Myc, Total	ALSU-TCMYC-A500	ALSU-TCMYC-A10K	ALSU-TCMYC-A50K	ALSU-TCMYC-A-HV
Cofilin, Phospho-S3	ALSU-PCOF-A500	ALSU-PCOF-A10K	ALSU-PCOF-A50K	ALSU-PCOF-A-HV
Cofilin, Total	ALSU-TCOF-A500	ALSU-TCOF-A10K	ALSU-TCOF-A50K	ALSU-TCOF-A-HV
CREB, Phospho-S133	ALSU-PCREB-A500	ALSU-PCREB-A10K	ALSU-PCREB-A50K	ALSU-PCREB-A-HV
CREB, Total	ALSU-TCREB-A500	ALSU-TCREB-A10K	ALSU-TCREB-A50K	ALSU-TCREB-A-HV
CREBBP, Total	ALSU-TCRBP-A500	ALSU-TCRBP-A10K	ALSU-TCRBP-A50K	ALSU-TCRBP-A-HV
Cyclin A2, Total	ALSU-TCYCA2-A500	ALSU-TCYCA2-A10K	ALSU-TCYCA2-A50K	ALSU-TCYCA2-A-HV
Cyclin B1, Phospho-S126	ALSU-PCYCB1-A500	ALSU-PCYCB1-A10K	ALSU-PCYCB1-A50K	ALSU-PCYCB1-A-HV
Cyclin B1, Total	ALSU-TCYCB1-A500	ALSU-TCYCB1-A10K	ALSU-TCYCB1-A50K	ALSU-TCYCB1-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
Cyclin D1, Phospho-T286	ALSU-PCYCD1-A500	ALSU-PCYCD1-A10K	ALSU-PCYCD1-A50K	ALSU-PCYCD1-A-HV
Cyclin D1, Total	ALSU-TCYCD1-A500	ALSU-TCYCD1-A10K	ALSU-TCYCD1-A50K	ALSU-TCYCD1-A-HV
<b>Cyclin D2, Total</b>	<b>ALSU-TCYCD2-A500</b>	<b>ALSU-TCYCD2-A10K</b>	<b>ALSU-TCYCD2-A50K</b>	<b>ALSU-TCYCD2-A-HV</b>
<b>Cyclin D3, Total</b>	<b>ALSU-TCYCD3-A500</b>	<b>ALSU-TCYCD3-A10K</b>	<b>ALSU-TCYCD3-A50K</b>	<b>ALSU-TCYCD3-A-HV</b>
Cyclin E1, Phospho-T77	ALSU-PCYCE1-A500	ALSU-PCYCE1-A10K	ALSU-PCYCE1-A50K	ALSU-PCYCE1-A-HV
Cyclin E1, Total	ALSU-TCYCE1-A500	ALSU-TCYCE1-A10K	ALSU-TCYCE1-A50K	ALSU-TCYCE1-A-HV
DAP12, Phospho-Y91	ALSU-PDAP12-B500	ALSU-PDAP12-B10K	ALSU-PDAP12-B50K	ALSU-PDAP12-B-HV
DAP12, Phospho-Y92 (Mouse)	ALSU-PDAP12-C500	ALSU-PDAP12-C10K	ALSU-PDAP12-C50K	ALSU-PDAP12-C-HV
DAP12, Phospho-parY	ALSU-PDAP12-A500	ALSU-PDAP12-A10K	ALSU-PDAP12-A50K	ALSU-PDAP12-A-HV
DAP12 Aggregate	ALSU-ADAP12-A500	ALSU-ADAP12-A10K	ALSU-ADAP12-A50K	ALSU-ADAP12-A-HV
DAP12, Total	ALSU-TDAP12-A500	ALSU-TDAP12-A10K	ALSU-TDAP12-A50K	ALSU-TDAP12-A-HV
DAP12, Total (Mouse)	ALSU-TDAP12-B500	ALSU-TDAP12-B10K	ALSU-TDAP12-B50K	ALSU-TDAP12-B-HV
<b>DJ1, Total</b>	<b>ALSU-TDJ1-A500</b>	<b>ALSU-TDJ1-A10K</b>	<b>ALSU-TDJ1-A50K</b>	<b>ALSU-TDJ1-A-HV</b>
DNA PKcs, Phospho-S2056	ALSU-PDNA-A500	ALSU-PDNA-A10K	ALSU-PDNA-A50K	ALSU-PDNA-A-HV
DNA PKcs, Total	ALSU-TDNA-A500	ALSU-TDNA-A10K	ALSU-TDNA-A50K	ALSU-TDNA-A-HV
EGF Receptor, Phospho-Y1068	ALSU-PEGFR-B500	ALSU-PEGFR-B10K	ALSU-PEGFR-B50K	ALSU-PEGFR-B-HV
EGF Receptor, Total	ALSU-TEGFR-A500	ALSU-TEGFR-A10K	ALSU-TEGFR-A50K	ALSU-TEGFR-A-HV
EGF Receptor L858R, Total	ALSU-TEGFR-B500	ALSU-TEGFR-B10K	ALSU-TEGFR-B50K	ALSU-TEGFR-B-HV
eIF2 $\alpha$ , Phospho-S51	ALSU-PEIF2-B500	ALSU-PEIF2-B10K	ALSU-PEIF2-B50K	ALSU-PEIF2-B-HV
eIF2 $\alpha$ , Total	ALSU-TEIF2-A500	ALSU-TEIF2-A10K	ALSU-TEIF2-A50K	ALSU-TEIF2-A-HV
eIF4E, Phospho-S209	ALSU-PEIF4-A500	ALSU-PEIF4-A10K	ALSU-PEIF4-A50K	ALSU-PEIF4-A-HV
eIF4E, Total	ALSU-TEIF4-A500	ALSU-TEIF4-A10K	ALSU-TEIF4-A50K	ALSU-TEIF4-A-HV
ER $\alpha$ , Phospho-S118	ALSU-PERA-A500	ALSU-PERA-A10K	ALSU-PERA-A50K	ALSU-PERA-A-HV
ER $\alpha$ , Phospho-S167	ALSU-PERA-B500	ALSU-PERA-B10K	ALSU-PERA-B50K	ALSU-PERA-B-HV
ER $\alpha$ , Total	ALSU-TERA-A500	ALSU-TERA-A10K	ALSU-TERA-A50K	ALSU-TERA-A-HV
ErbB2, Phospho-Y1221/1222	ALSU-PEB2-A500	ALSU-PEB2-A10K	ALSU-PEB2-A50K	ALSU-PEB2-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
ErbB2, Total	ALSU-TEB2-A500	ALSU-TEB2-A10K	ALSU-TEB2-A50K	ALSU-TEB2-A-HV
ErbB3, Phospho-Y1289	ALSU-PERB3-A500	ALSU-PERB3-A10K	ALSU-PERB3-A50K	ALSU-PERB3-A-HV
ErbB3, Total	ALSU-TERB3-A500	ALSU-TERB3-A10K	ALSU-TERB3-A50K	ALSU-TERB3-A-HV
ERK1/2, Phospho-T202/Y204	ALSU-PERK-A500	ALSU-PERK-A10K	ALSU-PERK-A50K	ALSU-PERK-A-HV
ERK1/2, Total	ALSU-TERK-A500	ALSU-TERK-A10K	ALSU-TERK-A50K	ALSU-TERK-A-HV
FAK, Phospho-Y397	ALSU-PFAK-A500	ALSU-PFAK-A10K	ALSU-PFAK-A50K	ALSU-PFAK-A-HV
FAK, Phospho-Y576/577	ALSU-PFAK-B500	ALSU-PFAK-B10K	ALSU-PFAK-B50K	ALSU-PFAK-B-HV
FAK, Total	ALSU-TFAK-A500	ALSU-TFAK-A10K	ALSU-TFAK-A50K	ALSU-TFAK-A-HV
FGFR1, Total	ALSU-TFGFR1-A500	ALSU-TFGFR1-A10K	ALSU-TFGFR1-A50K	ALSU-TFGFR1-A-HV
FGFR2, Phospho-Y653/654	ALSU-PFGFR2-A500	ALSU-PFGFR2-A10K	ALSU-PFGFR2-A50K	ALSU-PFGFR2-A-HV
FGFR2, Total	ALSU-TFGFR2-A500	ALSU-TFGFR2-A10K	ALSU-TFGFR2-A50K	ALSU-TFGFR2-A-HV
FGFR3, Phospho-Y724	ALSU-PFGFR3-A500	ALSU-PFGFR3-A10K	ALSU-PFGFR3-A50K	ALSU-PFGFR3-A-HV
FGFR3, Phospho-panY	ALSU-PFGFR3-B500	ALSU-PFGFR3-B10K	ALSU-PFGFR3-B50K	ALSU-PFGFR3-B-HV
FGFR3, Total	ALSU-TFGFR3-A500	ALSU-TFGFR3-A10K	ALSU-TFGFR3-A50K	ALSU-TFGFR3-A-HV
FGFR4, Total	ALSU-TFGFR4-A500	ALSU-TFGFR4-A10K	ALSU-TFGFR4-A50K	ALSU-TFGFR4-A-HV
FLT3, Phospho-Y591	ALSU-PFLT3-A500	ALSU-PFLT3-A10K	ALSU-PFLT3-A50K	ALSU-PFLT3-A-HV
FLT3, Phospho-Y842	ALSU-PFLT3-B500	ALSU-PFLT3-B10K	ALSU-PFLT3-B50K	ALSU-PFLT3-B-HV
FLT3, Total	ALSU-TFLT3-A500	ALSU-TFLT3-A10K	ALSU-TFLT3-A50K	ALSU-TFLT3-A-HV
FOXO1, Phospho-T600	ALSU-PFOXO1-A500	ALSU-PFOXO1-A10K	ALSU-PFOXO1-A50K	ALSU-PFOXO1-A-HV
FOXO1, Total	ALSU-TFOXO1-A500	ALSU-TFOXO1-A10K	ALSU-TFOXO1-A50K	ALSU-TFOXO1-A-HV
GAPDH, Total	ALSU-TGAPD-A500	ALSU-TGAPD-A10K	ALSU-TGAPD-A50K	ALSU-TGAPD-A-HV
GAPDH, Total (Mouse/Human)	ALSU-TGAPD-B500	ALSU-TGAPD-B10K	ALSU-TGAPD-B50K	ALSU-TGAPD-B-HV
GCN2, Phospho-T899	ALSU-PGCN2-A500	ALSU-PGCN2-A10K	ALSU-PGCN2-A50K	ALSU-PGCN2-A-HV
GCN2, Total	ALSU-TGCN2-A500	ALSU-TGCN2-A10K	ALSU-TGCN2-A50K	ALSU-TGCN2-A-HV
GSK3 $\beta$ , Phospho-S9	ALSU-PGS3B-A500	ALSU-PGS3B-A10K	ALSU-PGS3B-A50K	ALSU-PGS3B-A-HV
GSK3 $\beta$ , Total	ALSU-TGS3B-A500	ALSU-TGS3B-A10K	ALSU-TGS3B-A50K	ALSU-TGS3B-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
Histone H2A.X, Phospho S139	ALSU-PH2X-A500	ALSU-PH2X-A10K	ALSU-PH2X-A50K	ALSU-PH2X-A-HV
Histone H3, Phospho-S10/T11	ALSU-PHISH3-A500	ALSU-PHISH3-A10K	ALSU-PHISH3-A50K	ALSU-PHISH3-A-HV
Histone H3, Total	ALSU-THISH3-A500	ALSU-THISH3-A10K	ALSU-THISH3-A50K	ALSU-THISH3-A-HV
Histone H2A.X, Total	ALSU-TH2X-A500	ALSU-TH2X-A10K	ALSU-TH2X-A50K	ALSU-TH2X-A-HV
HDAC4, Phospho-S246	ALSU-PHDAC-A500	ALSU-PHDAC-A10K	ALSU-PHDAC-A50K	ALSU-PHDAC-A-HV
HDAC4, Total	ALSU-THDAC-A500	ALSU-THDAC-A10K	ALSU-THDAC-A50K	ALSU-THDAC-A-HV
HDAC5, Phospho-S259	ALSU-PHDAC5-A500	ALSU-PHDAC5-A10K	ALSU-PHDAC5-A50K	ALSU-PHDAC5-A-HV
HDAC5, Total	ALSU-THDAC5-A500	ALSU-THDAC5-A10K	ALSU-THDAC5-A50K	ALSU-THDAC5-A-HV
HSP27, Phospho-S15	ALSU-PHSP-A500	ALSU-PHSP-A10K	ALSU-PHSP-A50K	ALSU-PHSP-A-HV
HSP27, Phospho-S78	ALSU-PHSP-C500	ALSU-PHSP-C10K	ALSU-PHSP-C50K	ALSU-PHSP-C-HV
HSP27, Phospho-S82	ALSU-PHSP-B500	ALSU-PHSP-B10K	ALSU-PHSP-B50K	ALSU-PHSP-B-HV
HSP27, Total	ALSU-THSP-A500	ALSU-THSP-A10K	ALSU-THSP-A50K	ALSU-THSP-A-HV
IGF-1 Rec. $\beta$ , Phospho-Y1135/1136	ALSU-PIGFR-B500	ALSU-PIGFR-B10K	ALSU-PIGFR-B50K	ALSU-PIGFR-B-HV
IGF-1 Rec. $\beta$ Total	ALSU-TIGFR-A500	ALSU-TIGFR-A10K	ALSU-TIGFR-A50K	ALSU-TIGFR-A-HV
IKAROS, Total	ALSU-TIKARO-A500	ALSU-TIKARO-A10K	ALSU-TIKARO-A50K	ALSU-TIKARO-A-HV
I $\kappa$ B $\alpha$ , Phospho-S32/36	ALSU-PIKBA-A500	ALSU-PIKBA-A10K	ALSU-PIKBA-A50K	ALSU-PIKBA-A-HV
I $\kappa$ B $\alpha$ , Total	ALSU-TIKBA-A500	ALSU-TIKBA-A10K	ALSU-TIKBA-A50K	ALSU-TIKBA-A-HV
IKK $\alpha$ , Phospho-S176/180	ALSU-PIKKA-A500	ALSU-PIKKA-A10K	ALSU-PIKKA-A50K	ALSU-PIKKA-A-HV
IKK $\alpha$ , Total	ALSU-TIKKA-A500	ALSU-TIKKA-A10K	ALSU-TIKKA-A50K	ALSU-TIKKA-A-HV
IKK $\beta$ , Phospho-S177/181	ALSU-PIKKB-A500	ALSU-PIKKB-A10K	ALSU-PIKKB-A50K	ALSU-PIKKB-A-HV
IKK $\beta$ , Total	ALSU-TIKKB-A500	ALSU-TIKKB-A10K	ALSU-TIKKB-A50K	ALSU-TIKKB-A-HV
Insulin Rec. $\beta$ , Phospho-Y1150/1151	ALSU-PINR-B500	ALSU-PINR-B10K	ALSU-PINR-B50K	ALSU-PINR-B-HV
Insulin Rec. $\beta$ , Total	ALSU-TINR-A500	ALSU-TINR-A10K	ALSU-TINR-A50K	ALSU-TINR-A-HV
IRAK4, Phospho-T345/S346	ALSU-PIRAK-A500	ALSU-PIRAK-A10K	ALSU-PIRAK-A50K	ALSU-PIRAK-A-HV
IRAK4, Total	ALSU-TIRAK-A500	ALSU-TIRAK-A10K	ALSU-TIRAK-A50K	ALSU-TIRAK-A-HV
IRF3, Phospho-S386	ALSU-PIRF3-A500	ALSU-PIRF3-A10K	ALSU-PIRF3-A50K	ALSU-PIRF3-A-HV

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



## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
IRF3, Total	ALSU-TIRF3-A500	ALSU-TIRF3-A10K	ALSU-TIRF3-A50K	ALSU-TIRF3-A-HV
<b>IRF5, Phospho-S446</b>	<b>ALSU-PIRF5-A500</b>	<b>ALSU-PIRF5-A10K</b>	<b>ALSU-PIRF5-A50K</b>	<b>ALSU-PIRF5-A-HV</b>
<b>IRF5, Total</b>	<b>ALSU-TIRF5-A500</b>	<b>ALSU-TIRF5-A10K</b>	<b>ALSU-TIRF5-A50K</b>	<b>ALSU-TIRF5-A-HV</b>
JAK1, Total	ALSU-TJAK1-A500	ALSU-TJAK1-A10K	ALSU-TJAK1-A50K	ALSU-TJAK1-A-HV
JAK2, Total	ALSU-TJAK2-A500	ALSU-TJAK2-A10K	ALSU-TJAK2-A50K	ALSU-TJAK2-A-HV
JNK1/2, Total	ALSU-TJNK-A500	ALSU-TJNK-A10K	ALSU-TJNK-A50K	ALSU-TJNK-A-HV
JNK1/2/3, Phospho-T183/Y185	ALSU-PJNK-A500	ALSU-PJNK-A10K	ALSU-PJNK-A50K	ALSU-PJNK-A-HV
KAP1, Phospho-S824	ALSU-PKAP1-A500	ALSU-PKAP1-A10K	ALSU-PKAP1-A50K	ALSU-PKAP1-A-HV
KAP1, Total	ALSU-TKAP1-A500	ALSU-TKAP1-A10K	ALSU-TKAP1-A50K	ALSU-TKAP1-A-HV
<b>KRAS, Total</b>	<b>ALSU-TKRAS-A500</b>	<b>ALSU-TKRAS-A10K</b>	<b>ALSU-TKRAS-A50K</b>	<b>ALSU-TKRAS-A-HV</b>
Lamin A/C, Phospho-S22	ALSU-PLAM-A500	ALSU-PLAM-A10K	ALSU-PLAM-A50K	ALSU-PLAM-A-HV
Lamin A/C, Total	ALSU-TLAM-A500	ALSU-TLAM-A10K	ALSU-TLAM-A50K	ALSU-TLAM-A-HV
Lck, Phospho-Y505	ALSU-PLCK-A500	ALSU-PLCK-A10K	ALSU-PLCK-A50K	ALSU-PLCK-A-HV
Lck, Total	ALSU-TLCK-A500	ALSU-TLCK-A10K	ALSU-TLCK-A50K	ALSU-TLCK-A-HV
LRRK2, Phospho-S910	ALSU-PLRRK2-C500	ALSU-PLRRK2-C10K	ALSU-PLRRK2-C50K	ALSU-PLRRK2-C-HV
LRRK2, Phospho-S935	ALSU-PLRRK2-A500	ALSU-PLRRK2-A10K	ALSU-PLRRK2-A50K	ALSU-PLRRK2-A-HV
LRRK2, Phospho-S955	ALSU-PLRRK2-B500	ALSU-PLRRK2-B10K	ALSU-PLRRK2-B50K	ALSU-PLRRK2-B-HV
LRRK2, Total	ALSU-TLRRK2-A500	ALSU-TLRRK2-A10K	ALSU-TLRRK2-A50K	ALSU-TLRRK2-A-HV
MDM2, Phospho-S166	ALSU-PMDM2-A500	ALSU-PMDM2-A10K	ALSU-PMDM2-A50K	ALSU-PMDM2-A-HV
MDM2, Total	ALSU-TMDM2-A500	ALSU-TMDM2-A10K	ALSU-TMDM2-A50K	ALSU-TMDM2-A-HV
MEK1, Phospho-S218/222	ALSU-PMEK1-A500	ALSU-PMEK1-A10K	ALSU-PMEK1-A50K	ALSU-PMEK1-A-HV
MEK1, Total	ALSU-TMEK1-A500	ALSU-TMEK1-A10K	ALSU-TMEK1-A50K	ALSU-TMEK1-A-HV
MKK4, Phospho-S257	ALSU-PMKK4-A500	ALSU-PMKK4-A10K	ALSU-PMKK4-A50K	ALSU-PMKK4-A-HV
MKK4, Total	ALSU-TMKK4-A500	ALSU-TMKK4-A10K	ALSU-TMKK4-A50K	ALSU-TMKK4-A-HV
MLKL, Phospho-S345 (Mouse)	ALSU-PMLKL-A500	ALSU-PMLKL-A10K	ALSU-PMLKL-A50K	ALSU-PMLKL-A-HV
MLKL, Total (Mouse)	ALSU-TMLKL-B500	ALSU-TMLKL-B10K	ALSU-TMLKL-B50K	ALSU-TMLKL-B-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
MLKL, Phospho S358	ALSU-PMLKL-B500	ALSU-PMLKL-B10K	ALSU-PMLKL-B50K	ALSU-PMLKL-B-HV
MLKL, Total	ALSU-TMLKL-A500	ALSU-TMLKL-A10K	ALSU-TMLKL-A50K	ALSU-TMLKL-A-HV
mTOR, Phospho-S2448	ALSU-PMTOR-C500	ALSU-PMTOR-C10K	ALSU-PMTOR-C50K	ALSU-PMTOR-C-HV
mTOR, Total	ALSU-TMTOR-B500	ALSU-TMTOR-B10K	ALSU-TMTOR-B50K	ALSU-TMTOR-B-HV
NFκB p65, Phospho-S536	ALSU-PNFKB-A500	ALSU-PNFKB-A10K	ALSU-PNFKB-A50K	ALSU-PNFKB-A-HV
NFκB p65, Total	ALSU-TNFKB-A500	ALSU-TNFKB-A10K	ALSU-TNFKB-A50K	ALSU-TNFKB-A-HV
NPM, Phospho-T199	ALSU-PNPM-A500	ALSU-PNPM-A10K	ALSU-PNPM-A50K	ALSU-PNPM-A-HV
NPM, Total	ALSU-TNPM-A500	ALSU-TNPM-A10K	ALSU-TNPM-A50K	ALSU-TNPM-A-HV
Nrf2, Phospho-S40	ALSU-PNRF2-A500	ALSU-PNRF2-A10K	ALSU-PNRF2-A50K	ALSU-PNRF2-A-HV
Nrf2, Total	ALSU-TNRF2-A500	ALSU-TNRF2-A10K	ALSU-TNRF2-A50K	ALSU-TNRF2-A-HV
p27 KIP1, Phospho-S10	ALSU-PP27K-A500	ALSU-PP27K-A10K	ALSU-PP27K-A50K	ALSU-PP27K-A-HV
p27 KIP1, Total	ALSU-TP27K-A500	ALSU-TP27K-A10K	ALSU-TP27K-A50K	ALSU-TP27K-A-HV
p38 MAPK, Phospho-T180/Y182	ALSU-PP38-B500	ALSU-PP38-B10K	ALSU-PP38-B50K	ALSU-PP38-B-HV
p38 MAPK, Total	ALSU-TP38-B500	ALSU-TP38-B10K	ALSU-TP38-B50K	ALSU-TP38-B-HV
p53, Ac-K382	ALSU-ACP53-A500	ALSU-ACP53-A10K	ALSU-ACP53-A50K	ALSU-ACP53-A-HV
p53, Phospho-S15	ALSU-PP53-A500	ALSU-PP53-A10K	ALSU-PP53-A50K	ALSU-PP53-A-HV
p53, Total	ALSU-TP53-A500	ALSU-TP53-A10K	ALSU-TP53-A50K	ALSU-TP53-A-HV
p70 S6K, Phospho-T389	ALSU-PP70-A500	ALSU-PP70-A10K	ALSU-PP70-A50K	ALSU-PP70-A-HV
p70 S6K, Total	ALSU-TP70-A500	ALSU-TP70-A10K	ALSU-TP70-A50K	ALSU-TP70-A-HV
p300, Total	ALSU-TP300-A500	ALSU-TP300-A10K	ALSU-TP300-A50K	ALSU-TP300-A-HV
Parkin, Phospho-S65	ALSU-PPARK-A500	ALSU-PPARK-A10K	ALSU-PPARK-A50K	ALSU-PPARK-A-HV
Parkin, Total	ALSU-TPARK-A500	ALSU-TPARK-A10K	ALSU-TPARK-A50K	ALSU-TPARK-A-HV
PD-1, Phospho-panY	ALSU-PPD1-A500	ALSU-PPD1-A10K	ALSU-PPD1-A50K	ALSU-PPD1-A-HV
PD-1, Total	ALSU-TPD1-A500	ALSU-TPD1-A10K	ALSU-TPD1-A50K	ALSU-TPD1-A-HV
PDGFRβ, Phospho-Y751	ALSU-PPDGF-A500	ALSU-PPDGF-A10K	ALSU-PPDGF-A50K	ALSU-PPDGF-A-HV
PDGFRβ, Total	ALSU-TPDGF-A500	ALSU-TPDGF-A10K	ALSU-TPDGF-A50K	ALSU-TPDGF-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
PDHA1, Phospho-S293	ALSU-PPDHA1-A500	ALSU-PPDHA1-A10K	ALSU-PPDHA1-A50K	ALSU-PPDHA1-A-HV
PDHA1, Total	ALSU-TPDHA1-A500	ALSU-TPDHA1-A10K	ALSU-TPDHA1-A50K	ALSU-TPDHA1-A-HV
PERK, Phospho-T980 (Mouse)	ALSU-PPERK-A500	ALSU-PPERK-A10K	ALSU-PPERK-A50K	ALSU-PPERK-A-HV
PERK, Total	ALSU-TPERK-A500	ALSU-TPERK-A10K	ALSU-TPERK-A50K	ALSU-TPERK-A-HV
PI3 Kinase p85, Phospho-Y730 (Mouse)	ALSU-PP13K-A500	ALSU-PP13K-A10K	ALSU-PP13K-A50K	ALSU-PP13K-A-HV
P13 Kinase p85, Total	ALSU-TP13K-A500	ALSU-TP13K-A10K	ALSU-TP13K-A50K	ALSU-TP13K-A-HV
PINK1, Total	ALSU-TPINK-A500	ALSU-TPINK-A10K	ALSU-TPINK-A50K	ALSU-TPINK-A-HV
PKCθ, Phospho-T538	ALSU-PPKC-A500	ALSU-PPKC-A10K	ALSU-PPKC-A50K	ALSU-PPKC-A-HV
PKR, Phospho-T451	ALSU-PPKR-B500	ALSU-PPKR-B10K	ALSU-PPKR-B50K	ALSU-PPKR-B-HV
PKR, Total	ALSU-TPKR-A500	ALSU-TPKR-A10K	ALSU-TPKR-A50K	ALSU-TPKR-A-HV
PLC G1, Phospho-Y783	ALSU-PPLCG1-A500	ALSU-PPLCG1-A10K	ALSU-PPLCG1-A50K	ALSU-PPLCG1-A-HV
PLC G1, Phospho-S1248	ALSU-PPLCG1-B500	ALSU-PPLCG1-B10K	ALSU-PPLCG1-B50K	ALSU-PPLCG1-B-HV
PLC G1, Phospho-Y1253	ALSU-PPLCG1-C500	ALSU-PPLCG1-C10K	ALSU-PPLCG1-C50K	ALSU-PPLCG1-C-HV
PLC G1, Total	ALSU-TPLCG1-A500	ALSU-TPLCG1-A10K	ALSU-TPLCG1-A50K	ALSU-TPLCG1-A-HV
PLC G2, Phospho-Y753	ALSU-PPLCG2-A500	ALSU-PPLCG2-A10K	ALSU-PPLCG2-A50K	ALSU-PPLCG2-A-HV
PLC G2, Phospho-Y759	ALSU-PPLCG2-B500	ALSU-PPLCG2-B10K	ALSU-PPLCG2-B50K	ALSU-PPLCG2-B-HV
PLC G2, Phospho-Y1197	ALSU-PPLCG2-C500	ALSU-PPLCG2-C10K	ALSU-PPLCG2-C50K	ALSU-PPLCG2-C-HV
PLC G2, Phospho-Y1217	ALSU-PPLCG2-D500	ALSU-PPLCG2-D10K	ALSU-PPLCG2-D50K	ALSU-PPLCG2-D-HV
PLC G2, Total	ALSU-TPLCG2-A500	ALSU-TPLCG2-A10K	ALSU-TPLCG2-A50K	ALSU-TPLCG2-A-HV
PLK1, Phospho-T210	ALSU-PPLK1-A500	ALSU-PPLK1-A10K	ALSU-PPLK1-A50K	ALSU-PPLK1-A-HV
PLK1, Total	ALSU-TPLK1-A500	ALSU-TPLK1-A10K	ALSU-TPLK1-A50K	ALSU-TPLK1-A-HV
<b>PTEN, Phospho-T366</b>	<b>ALSU-PPTEN-A500</b>	<b>ALSU-PPTEN-A10K</b>	<b>ALSU-PPTEN-A50K</b>	<b>ALSU-PPTEN-A-HV</b>
<b>PTEN, Total</b>	<b>ALSU-TPTEN-A500</b>	<b>ALSU-TPTEN-A10K</b>	<b>ALSU-TPTEN-A50K</b>	<b>ALSU-TPTEN-A-HV</b>
RAB 8A, Phospho-S111	ALSU-PRAB8-A500	ALSU-PRAB8-A10K	ALSU-PRAB8-A50K	ALSU-PRAB8-A-HV
RAB 8A, Phospho-T72	ALSU-PRAB8-B500	ALSU-PRAB8-B10K	ALSU-PRAB8-B50K	ALSU-PRAB8-B-HV
RAB 8A, Total	ALSU-TRAB8-A500	ALSU-TRAB8-A10K	ALSU-TRAB8-A50K	ALSU-TRAB8-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
RAB10, Phospho-T73	ALSU-PRAB-A500	ALSU-PRAB-A10K	ALSU-PRAB-A50K	ALSU-PRAB-A-HV
RAB10, Total	ALSU-TRAB-A500	ALSU-TRAB-A10K	ALSU-TRAB-A50K	ALSU-TRAB-A-HV
RAB12, Phospho-S106	ALSU-PRAB12-A500	ALSU-PRAB12-A10K	ALSU-PRAB12-A50K	ALSU-PRAB12-A-HV
RAB29, Phospho-S72	ALSU-PRAB29-A500	ALSU-PRAB29-A10K	ALSU-PRAB29-A50K	ALSU-PRAB29-A-HV
RAB29, Total	ALSU-TRAB29-A500	ALSU-TRAB29-A10K	ALSU-TRAB29-A50K	ALSU-TRAB29-A-HV
Raf-1, Phospho-S338	ALSU-PRAF-A500	ALSU-PRAF-A10K	ALSU-PRAF-A50K	ALSU-PRAF-A-HV
Raf-1, Total	ALSU-TRAF-A500	ALSU-TRAF-A10K	ALSU-TRAF-A50K	ALSU-TRAF-A-HV
RAPTOR, Total	ALSU-TRAPT-A500	ALSU-TRAPT-A10K	ALSU-TRAPT-A50K	ALSU-TRAPT-A-HV
Ras, Total	ALSU-TRAS-A500	ALSU-TRAS-A10K	ALSU-TRAS-A50K	ALSU-TRAS-A-HV
Rb, Phospho-S780	ALSU-PRB-C500	ALSU-PRB-C10K	ALSU-PRB-C50K	ALSU-PRB-C-HV
Rb, Phospho-S807/811	ALSU-PRB-A500	ALSU-PRB-A10K	ALSU-PRB-A50K	ALSU-PRB-A-HV
Rb, Phospho-T821/826	ALSU-PRB-B500	ALSU-PRB-B10K	ALSU-PRB-B50K	ALSU-PRB-B-HV
Rb, Total	ALSU-TRB-A500	ALSU-TRB-A10K	ALSU-TRB-A50K	ALSU-TRB-A-HV
RBM39, Total	ALSU-TRBM39-A500	ALSU-TRBM39-A10K	ALSU-TRBM39-A50K	ALSU-TRBM39-A-HV
RET, Phospho panY	ALSU-PRET-A500	ALSU-PRET-A10K	ALSU-PRET-A50K	ALSU-PRET-A-HV
RET, Total	ALSU-TRET-A500	ALSU-TRET-A10K	ALSU-TRET-A50K	ALSU-TRET-A-HV
Ribosomal Protein S6, Phospho-S235/236	ALSU-PS6R-B500	ALSU-PS6R-B10K	ALSU-PS6R-B50K	ALSU-PS6R-B-HV
Ribosomal Protein S6, Phospho-S240/244	ALSU-PS6R-A500	ALSU-PS6R-A10K	ALSU-PS6R-A50K	ALSU-PS6R-A-HV
Ribosomal Protein S6, Total	ALSU-TS6R-A500	ALSU-TS6R-A10K	ALSU-TS6R-A50K	ALSU-TS6R-A-HV
RIP3, Phospho-S227 (Human)	ALSU-PRIP3-A500	ALSU-PRIP3-A10K	ALSU-PRIP3-A50K	ALSU-PRIP3-A-HV
RIP3, Total (Human)	ALSU-TRIP3-A500	ALSU-TRIP3-A10K	ALSU-TRIP3-A50K	ALSU-TRIP3-A-HV
RIP3, Phospho-T231/S232 (Mouse)	ALSU-PRIP3-B500	ALSU-PRIP3-B10K	ALSU-PRIP3-B50K	ALSU-PRIP3-B-HV
RIP3, Total (Mouse)	ALSU-TRIP3-B500	ALSU-TRIP3-B10K	ALSU-TRIP3-B50K	ALSU-TRIP3-B-HV
RIPK1, Phospho-S166	ALSU-PRIPK1-A500	ALSU-PRIPK1-A10K	ALSU-PRIPK1-A50K	ALSU-PRIPK1-A-HV
RIPK1, Total	ALSU-TRIPK1-A500	ALSU-TRIPK1-A10K	ALSU-TRIPK1-A50K	ALSU-TRIPK1-A-HV
RNA POL II, Phospho-S2	ALSU-PRNAP-A500	ALSU-PRNAP-A10K	ALSU-PRNAP-A50K	ALSU-PRNAP-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
RNA POL II, Phospho-S5	ALSU-PRNAP-B500	ALSU-PRNAP-B10K	ALSU-PRNAP-B50K	ALSU-PRNAP-B-HV
RNA POL II, Total	ALSU-TRNAP-A500	ALSU-TRNAP-A10K	ALSU-TRNAP-A50K	ALSU-TRNAP-A-HV
RSK1, Phospho-T359	ALSU-PRSK1-B500	ALSU-PRSK1-B10K	ALSU-PRSK1-B50K	ALSU-PRSK1-B-HV
RSK1, Phospho-S380	ALSU-PRSK1-C500	ALSU-PRSK1-C10K	ALSU-PRSK1-C50K	ALSU-PRSK1-C-HV
RSK1, Phospho-T573	ALSU-PRSK1-A500	ALSU-PRSK1-A10K	ALSU-PRSK1-A50K	ALSU-PRSK1-A-HV
RSK1, Total	ALSU-TRSK1-A500	ALSU-TRSK1-A10K	ALSU-TRSK1-A50K	ALSU-TRSK1-A-HV
SHP-1, Phospho-Y536	ALSU-PSHP1-A500	ALSU-PSHP1-A10K	ALSU-PSHP1-A50K	ALSU-PSHP1-A-HV
SHP-1, Phospho-Y564	ALSU-PSHP1-B500	ALSU-PSHP1-B10K	ALSU-PSHP1-B50K	ALSU-PSHP1-B-HV
SHP-1, Total	ALSU-TSHP1-A500	ALSU-TSHP1-A10K	ALSU-TSHP1-A50K	ALSU-TSHP1-A-HV
SHP-2, Phospho-Y542	ALSU-PSHP2-A500	ALSU-PSHP2-A10K	ALSU-PSHP2-A50K	ALSU-PSHP2-A-HV
SHP-2, Phospho-Y580	ALSU-PSHP2-B500	ALSU-PSHP2-B10K	ALSU-PSHP2-B50K	ALSU-PSHP2-B-HV
SHP-2, Total	ALSU-TSHP2-A500	ALSU-TSHP2-A10K	ALSU-TSHP2-A50K	ALSU-TSHP2-A-HV
SLP-76, Phospho-S376	ALSU-PSLP-A500	ALSU-PSLP-A10K	ALSU-PSLP-A50K	ALSU-PSLP-A-HV
SLP76, Phospho-S376 (Mouse)	ALSU-PSLP-D500	ALSU-PSLP-D10K	ALSU-PSLP-D50K	ALSU-PSLP-D-HV
SLP76, Phospho-Y113	ALSU-PSLP-B500	ALSU-PSLP-B10K	ALSU-PSLP-B50K	ALSU-PSLP-B-HV
SLP76, Phospho-Y145	ALSU-PSLP-C500	ALSU-PSLP-C10K	ALSU-PSLP-C50K	ALSU-PSLP-C-HV
SLP76, Phospho-Y145 (Mouse)	ALSU-PSLP-E500	ALSU-PSLP-E10K	ALSU-PSLP-E50K	ALSU-PSLP-E-HV
SLP-76, Total	ALSU-TSLP-A500	ALSU-TSLP-A10K	ALSU-TSLP-A50K	ALSU-TSLP-A-HV
SLP76, Total (Mouse)	ALSU-TSLP-B500	ALSU-TSLP-B10K	ALSU-TSLP-B50K	ALSU-TSLP-B-HV
SMAD1, Phospho-S463/465	ALSU-PSM1-A500	ALSU-PSM1-A10K	ALSU-PSM1-A50K	ALSU-PSM1-A-HV
SMAD1, Total	ALSU-TSM1-A500	ALSU-TSM1-A10K	ALSU-TSM1-A50K	ALSU-TSM1-A-HV
SMAD2, Phospho-S465/467	ALSU-PSM2-A500	ALSU-PSM2-A10K	ALSU-PSM2-A50K	ALSU-PSM2-A-HV
SMAD2, Total	ALSU-TSM2-A500	ALSU-TSM2-A10K	ALSU-TSM2-A50K	ALSU-TSM2-A-HV
SMAD3, Phospho-S423/425	ALSU-PSM3-A500	ALSU-PSM3-A10K	ALSU-PSM3-A50K	ALSU-PSM3-A-HV
SMAD3, Total	ALSU-TSM3-A500	ALSU-TSM3-A10K	ALSU-TSM3-A50K	ALSU-TSM3-A-HV
SMAD4, Total	ALSU-TSM4-A500	ALSU-TSM4-A10K	ALSU-TSM4-A50K	ALSU-TSM4-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
SMARCA4, Total	ALSU-TBRG1-A500	ALSU-TBRG1-A10K	ALSU-TBRG1-A50K	ALSU-TBRG1-A-HV
SQSTM1 p62, Phospho-S351	ALSU-PSQSTM-A500	ALSU-PSQSTM-A10K	ALSU-PSQSTM-A50K	ALSU-PSQSTM-A-HV
SQSTM1 p62, Phospho-S403	ALSU-PSQSTM-B500	ALSU-PSQSTM-B10K	ALSU-PSQSTM-B50K	ALSU-PSQSTM-B-HV
SQSTM1 p62, Total	ALSU-TSQSTM-A500	ALSU-TSQSTM-A10K	ALSU-TSQSTM-A50K	ALSU-TSQSTM-A-HV
Src, Phospho-Y419	ALSU-PSRC-A500	ALSU-PSRC-A10K	ALSU-PSRC-A50K	ALSU-PSRC-A-HV
STAT1, Phospho-S727	ALSU-PST1-B500	ALSU-PST1-B10K	ALSU-PST1-B50K	ALSU-PST1-B-HV
STAT1, Phospho-Y701	ALSU-PST1-A500	ALSU-PST1-A10K	ALSU-PST1-A50K	ALSU-PST1-A-HV
STAT1, Total	ALSU-TST1-A500	ALSU-TST1-A10K	ALSU-TST1-A50K	ALSU-TST1-A-HV
STAT3, Phospho-Y705	ALSU-PST3-A500	ALSU-PST3-A10K	ALSU-PST3-A50K	ALSU-PST3-A-HV
STAT3, Total	ALSU-TST3-A500	ALSU-TST3-A10K	ALSU-TST3-A50K	ALSU-TST3-A-HV
STAT4, Phospho-Y693	ALSU-PST4-A500	ALSU-PST4-A10K	ALSU-PST4-A50K	ALSU-PST4-A-HV
STAT4, Total	ALSU-TST4-A500	ALSU-TST4-A10K	ALSU-TST4-A50K	ALSU-TST4-A-HV
STAT5, Phospho-Y694/699	ALSU-PST5-B500	ALSU-PST5-B10K	ALSU-PST5-B50K	ALSU-PST5-B-HV
STAT5, Total	ALSU-TST5-A500	ALSU-TST5-A10K	ALSU-TST5-A50K	ALSU-TST5-A-HV
STAT6, Phospho-Y641	ALSU-PST6-A500	ALSU-PST6-A10K	ALSU-PST6-A50K	ALSU-PST6-A-HV
STAT6, Total	ALSU-TST6-A500	ALSU-TST6-A10K	ALSU-TST6-A50K	ALSU-TST6-A-HV
STING, Total	ALSU-TSTNG-A500	ALSU-TSTNG-A10K	ALSU-TSTNG-A50K	ALSU-TSTNG-A-HV
SYK, Phospho-Y352	ALSU-PSYK-B500	ALSU-PSYK-B10K	ALSU-PSYK-B50K	ALSU-PSYK-B-HV
SYK, Phospho-Y525/526	ALSU-PSYK-A500	ALSU-PSYK-A10K	ALSU-PSYK-A50K	ALSU-PSYK-A-HV
SYK, Total	ALSU-TSYK-A500	ALSU-TSYK-A10K	ALSU-TSYK-A50K	ALSU-TSYK-A-HV
SYK, Total (Mouse)	ALSU-TSYK-B500	ALSU-TSYK-B10K	ALSU-TSYK-B50K	ALSU-TSYK-B-HV
TAK1, Phospho-T184/187	ALSU-PTAK-A500	ALSU-PTAK-A10K	ALSU-PTAK-A50K	ALSU-PTAK-A-HV
TAK1, Total	ALSU-TTAK-A500	ALSU-TTAK-A10K	ALSU-TTAK-A50K	ALSU-TTAK-A-HV
Tau, Phospho-S214	ALSU-PTAU-A500	ALSU-PTAU-A10K	ALSU-PTAU-A50K	ALSU-PTAU-A-HV
Tau, Phospho-S404	ALSU-PTAU-C500	ALSU-PTAU-C10K	ALSU-PTAU-C50K	ALSU-PTAU-C-HV
Tau, Phospho-S396	ALSU-PTAU-B500	ALSU-PTAU-B10K	ALSU-PTAU-B50K	ALSU-PTAU-B-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
Tau, Phospho-T181	ALSU-PTAU-D500	ALSU-PTAU-D10K	ALSU-PTAU-D50K	ALSU-PTAU-D-HV
Tau, Phospho-T205	ALSU-PTAU-E500	ALSU-PTAU-E10K	ALSU-PTAU-E50K	ALSU-PTAU-E-HV
Tau, Total	ALSU-TTAU-A500	ALSU-TTAU-A10K	ALSU-TTAU-A50K	ALSU-TTAU-A-HV
TBK1, Phospho-S172	ALSU-PTBK1-A500	ALSU-PTBK1-A10K	ALSU-PTBK1-A50K	ALSU-PTBK1-A-HV
TBK1, Total	ALSU-TTBK1-A500	ALSU-TTBK1-A10K	ALSU-TTBK1-A50K	ALSU-TTBK1-A-HV
TDP43, Phospho-S409/410	ALSU-PTDP43-A500	ALSU-PTDP43-A10K	ALSU-PTDP43-A50K	ALSU-PTDP43-A-HV
TDP43, Phospho-S409/410 (Mouse)	ALSU-PTDP43-B500	ALSU-PTDP43-B10K	ALSU-PTDP43-B50K	ALSU-PTDP43-B-HV
TDP43 Aggregate	ALSU-ATDP43-A500	ALSU-ATDP43-A10K	ALSU-ATDP43-A50K	ALSU-ATDP43-A-HV
TDP43, Total (Mouse)	ALSU-TTDP43-B500	ALSU-TTDP43-B10K	ALSU-TTDP43-B50K	ALSU-TTDP43-B-HV
TDP43, Total	ALSU-TTDP43-A500	ALSU-TTDP43-A10K	ALSU-TTDP43-A50K	ALSU-TTDP43-A-HV
<b>TFE3, Phospho-S321</b>	<b>ALSU-PTFE3-A500</b>	<b>ALSU-PTFE3-A10K</b>	<b>ALSU-PTFE3-A50K</b>	<b>ALSU-PTFE3-A-HV</b>
<b>TFE3, Total</b>	<b>ALSU-TTFE3-A500</b>	<b>ALSU-TTFE3-A10K</b>	<b>ALSU-TTFE3-A50K</b>	<b>ALSU-TTFE3-A-HV</b>
TFEB, Phospho-S122	ALSU-PTFEB-A500	ALSU-PTFEB-A10K	ALSU-PTFEB-A50K	ALSU-PTFEB-A-HV
TFEB, Phospho-S122 (Mouse)	ALSU-PTFEB-C500	ALSU-PTFEB-C10K	ALSU-PTFEB-C50K	ALSU-PTFEB-C-HV
TFEB, Phospho-S211	ALSU-PTFEB-B500	ALSU-PTFEB-B10K	ALSU-PTFEB-B50K	ALSU-PTFEB-B-HV
TFEB, Phospho-S211 (Mouse)	ALSU-PTFEB-D500	ALSU-PTFEB-D10K	ALSU-PTFEB-D50K	ALSU-PTFEB-D-HV
TFEB, Total	ALSU-TTFEB-A500	ALSU-TTFEB-A10K	ALSU-TTFEB-A50K	ALSU-TTFEB-A-HV
TFEB, Total (Mouse)	ALSU-TTFEB-B500	ALSU-TTFEB-B10K	ALSU-TTFEB-B50K	ALSU-TTFEB-B-HV
TREM2 Aggregate	ALSU-ATREM2-A500	ALSU-ATREM2-A10K	ALSU-ATREM2-A50K	ALSU-ATREM2-A-HV
TREM2, Total	ALSU-TTREM2-A500	ALSU-TTREM2-A10K	ALSU-TTREM2-A50K	ALSU-TTREM2-A-HV
TREM2/DAP12, Total	ALSU-TTRMDP-A500	ALSU-TTRMDP-A10K	ALSU-TTRMDP-A50K	ALSU-TTRMDP-A-HV
TrkA, Phospho-Y674/675 / TrkB, Phospho-Y706/707	ALSU-PTRKAB-A500	ALSU-PTRKAB-A10K	ALSU-PTRKAB-A50K	ALSU-PTRKAB-A-HV
<b>TYK2, Phospho-Y1054/1055</b>	<b>ALSU-PTYK2-A500</b>	<b>ALSU-PTYK2-A10K</b>	<b>ALSU-PTYK2-A50K</b>	<b>ALSU-PTYK2-A-HV</b>
<b>TYK2, Total</b>	<b>ALSU-TTYK2-A500</b>	<b>ALSU-TTYK2-A10K</b>	<b>ALSU-TTYK2-A50K</b>	<b>ALSU-TTYK2-A-HV</b>
Ubiquitin, Phospho-S65	ALSU-PUBI-A500	ALSU-PUBI-A10K	ALSU-PUBI-A50K	ALSU-PUBI-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra kits			
	500 assay points	10,000 assay points	50,000 assay points	100 assay points
Ubiquitin, Total	ALSU-TUBI-A500	ALSU-TUBI-A10K	ALSU-TUBI-A50K	ALSU-TUBI-A-HV
ULK1, Phospho-S556	ALSU-PULK1-A500	ALSU-PULK1-A10K	ALSU-PULK1-A50K	ALSU-PULK1-A-HV
ULK1, Phospho-S758	ALSU-PULK1-B500	ALSU-PULK1-B10K	ALSU-PULK1-B50K	ALSU-PULK1-B-HV
ULK1, Total	ALSU-TULK1-A500	ALSU-TULK1-A10K	ALSU-TULK1-A50K	ALSU-TULK1-A-HV
VAV1, Phospho-Y174	ALSU-PVAV1-A500	ALSU-PVAV1-A10K	ALSU-PVAV1-A50K	ALSU-PVAV1-A-HV
VAV1, Total	ALSU-TVAV1-A500	ALSU-TVAV1-A10K	ALSU-TVAV1-A50K	ALSU-TVAV1-A-HV
VEGF Receptor 2, Phospho-Y951	ALSU-PVGFR-B500	ALSU-PVGFR-B10K	ALSU-PVGFR-B50K	ALSU-PVGFR-B-HV
VEGF Receptor 2, Phospho-Y1175	ALSU-PVGFR-A500	ALSU-PVGFR-A10K	ALSU-PVGFR-A50K	ALSU-PVGFR-A-HV
VEGF Receptor 2, Phospho-Y1214	ALSU-PVGFR-C500	ALSU-PVGFR-C10K	ALSU-PVGFR-C50K	ALSU-PVGFR-C-HV
VEGF Receptor 2, Total	ALSU-TVGFR-A500	ALSU-TVGFR-A10K	ALSU-TVGFR-A50K	ALSU-TVGFR-A-HV
WDR5, Total	ALSU-TWDR-A500	ALSU-TWDR-A10K	ALSU-TWDR-A50K	ALSU-TWDR-A-HV
Wee1, Total	ALSU-TWEE-A500	ALSU-TWEE-A10K	ALSU-TWEE-A50K	ALSU-TWEE-A-HV
<b>XBP1s, Total</b>	<b>ALSU-TXBP1-A500</b>	<b>ALSU-TXBP1-A10K</b>	<b>ALSU-TXBP1-A50K</b>	<b>ALSU-TXBP1-A-HV</b>
YAP, Phospho-S127	ALSU-PYAP-B500	ALSU-PYAP-B10K	ALSU-PYAP-B50K	ALSU-PYAP-B-HV
YAP, Total	ALSU-TYAP-A500	ALSU-TYAP-A10K	ALSU-TYAP-A50K	ALSU-TYAP-A-HV
ZAP-70, Phospho-Y319	ALSU-PZAP-A500	ALSU-PZAP-A10K	ALSU-PZAP-A50K	ALSU-PZAP-A-HV
ZAP-70, Total	ALSU-TZAP-A500	ALSU-TZAP-A10K	ALSU-TZAP-A50K	ALSU-TZAP-A-HV

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



## Ordering information

Analyte	Alpha SureFire Ultra Multiplex			
	500 Assay Points	10,000 Assay Points	50,000 Assay Points	100 Assay Points
AKT 1/2/3, Phospho-T308 + AKT 1/2/3, Total	MPSU-PTAKT-O500	MPSU-PTAKT-O10K	MPSU-PTAKT-O50K	MPSU-PTAKT-O-HV
AKT 1/2/3, Phospho-T450 + AKT 1/2/3, Total	MPSU-PTAKT-N500	MPSU-PTAKT-N10K	MPSU-PTAKT-N50K	MPSU-PTAKT-N-HV
AKT1/2/3, Phospho-S473 + AKT1, Total	MPSU-PTAKT-M500	MPSU-PTAKT-M10K	MPSU-PTAKT-M50K	MPSU-PTAKT-M-HV
ALK, Phospho-Y1604 + ALK, Total	MPSU-PTALK-K500	MPSU-PTALK-K10K	MPSU-PTALK-K50K	MPSU-PTALK-K-HV
AMPK $\alpha$ 1/2, Phospho-T172 + AMPK $\alpha$ 1/2, Total	MPSU-PTAMPK-K500	MPSU-PTAMPK-K10K	MPSU-PTAMPK-K50K	MPSU-PTAMPK-K-HV
Btk, Phospho-Y223 + Btk, Total	MPSU-PTBTK-K500	MPSU-PTBTK-K10K	MPSU-PTBTK-K50K	MPSU-PTBTK-K-HV
CDK1, Phospho-T14 + CDK1, Total	MPSU-PTCDK1-K500	MPSU-PTCDK1-K10K	MPSU-PTCDK1-K50K	MPSU-PTCDK1-K-HV
CDK1, Phospho-Y15 + CDK1, Total	MPSU-PTCDK1-M500	MPSU-PTCDK1-M10K	MPSU-PTCDK1-M50K	MPSU-PTCDK1-M-HV
Cofilin, Phospho-S3 + Cofilin, Total	MPSU-PTCOF-K500	MPSU-PTCOF-K10K	MPSU-PTCOF-K50K	MPSU-PTCOF-K-HV
DAP12, Phospho-Y91 + DAP12, Total	MPSU-PTDAP12-K500	MPSU-PTDAP12-K10K	MPSU-PTDAP12-K50K	MPSU-PTDAP12-K-HV
EGFR, Phospho-Y1068 + EGFR, Total	MPSU-PTEGFR-K500	MPSU-PTEGFR-K10K	MPSU-PTEGFR-K50K	MPSU-PTEGFR-K-HV
eIF2a, Phospho-S51 + eIF2a, Total	MPSU-PTEIF2-K500	MPSU-PTEIF2-K10K	MPSU-PTEIF2-K50K	MPSU-PTEIF2-K-HV
ErbB2, Phospho-Y1221/1222 + ErbB2, Total	MPSU-PTEB2-K500	MPSU-PTEB2-K10K	MPSU-PTEB2-K50K	MPSU-PTEB2-K-HV
ErbB3, Phospho-Y1289 + ErbB3, Total	MPSU-PTERB3-K500	MPSU-PTERB3-K10K	MPSU-PTERB3-K50K	MPSU-PTERB3-K-HV
ERK1/2, Phospho-T202/Y204 + ERK1/2, Total	MPSU-PTERK-M500	MPSU-PTERK-M10K	MPSU-PTERK-M50K	MPSU-PTERK-M-HV
Histone H2A.X, Phospho p-S139 + Histone H2A.X, Total	MPSU-PTH2X-K500	MPSU-PTH2X-K10K	MPSU-PTH2X-K50K	MPSU-PTH2X-K-HV
IGF-1 Rec. $\beta$ , Phospho-Y1135/1136 + IGF-1 Rec. $\beta$ , Total	MPSU-PTIGFR-K500	MPSU-PTIGFR-K10K	MPSU-PTIGFR-K50K	MPSU-PTIGFR-K-HV
Insulin Rec. $\beta$ , Phospho-Y1150/1151 + Insulin Rec. $\beta$ , Total	MPSU-PTINR-K500	MPSU-PTINR-K10K	MPSU-PTINR-K50K	MPSU-PTINR-K-HV
MEK,1 Phospho-S218/222 + MEK1, Total	MPSU-PTMEK1-K500	MPSU-PTMEK1-K10K	MPSU-PTMEK1-K50K	MPSU-PTMEK1-K-HV
NFkB p65, Phospho-S536 + NFkB p65, Total	MPSU-PTNFKB-K500	MPSU-PTNFKB-K10K	MPSU-PTNFKB-K50K	MPSU-PTNFKB-K-HV
p53, Acetyl-K382 + p53, Total	MPSU-ATP53-K500	MPSU-ATP53-K10K	MPSU-ATP53-K50K	MPSU-ATP53-K-HV
p53, Phospho-S15 + p53, Total	MPSU-PTP53-K500	MPSU-PTP53-K10K	MPSU-PTP53-K50K	MPSU-PTP53-K-HV
p70S6K, Phospho-T389 + p70S6K, Total	MPSU-PTP70-K500	MPSU-PTP70-K10K	MPSU-PTP70-K50K	MPSU-PTP70-K-HV
PDGFR $\beta$ , Phospho-Y751 + PDGFR $\beta$ , Total	MPSU-PTPDGF-K500	MPSU-PTPDGF-K10K	MPSU-PTPDGF-K50K	MPSU-PTPDGF-K-HV

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

## Ordering information

Analyte	Alpha SureFire Ultra Multiplex			
	500 Assay Points	10,000 Assay Points	50,000 Assay Points	100 Assay Points
Rb, Phospho-S780 + Rb, Total	MPSU-PTRB-K500	MPSU-PTRB-K10K	MPSU-PTRB-K50K	MPSU-PTRB-K5-HV
Rb, Phospho-S807/811 + Rb, Total	MPSU-PTRB-M500	MPSU-PTRB-M10K	MPSU-PTRB-M50K	MPSU-PTRB-M-HV
Rb, Phospho-T821/826 + Rb, Total	MPSU-PTRB-N500	MPSU-PTRB-N10K	MPSU-PTRB-N50K	MPSU-PTRB-N-HV
Rb, Phospho-T821/826 + Rb, Total	MPSU-PTRPS6-B500	MPSU-PTRPS6-B10K	MPSU-PTRPS6-B50K	MPSU-PTRPS6-B-HV
Ret, Phospho-panY + Ret, Total	MPSU-PTRET-K500	MPSU-PTRET-K10K	MPSU-PTRET-K50K	MPSU-PTRET-K-HV
Ribosomal Protein S6, Phospho-S235/236 + Ribosomal Protein S6, Total	MPSU-PTRPS6-A500	MPSU-PTRPS6-A10K	MPSU-PTRPS6-A50K	MPSU-PTRPS6-A-HV
SLP-76, Phospho-S376 + SLP-76, Total	MPSU-PTSPLP-K500	MPSU-PTSPLP-K10K	MPSU-PTSPLP-K50K	MPSU-PTSPLP-K-HV
SMAD2, Phospho-S465/467 + SMAD2, Total	MPSU-PTSM2-K500	MPSU-PTSM2-K10K	MPSU-PTSM2-K50K	MPSU-PTSM2-K-HV
STAT1, Phospho-Y701 + STAT1, Total	MPSU-PTST1-K500	MPSU-PTST1-K10K	MPSU-PTST1-K50K	MPSU-PTST1-K-HV
STAT1, Phospho-S727 + STAT1, Total	MPSU-PTST1-M500	MPSU-PTST1-M10K	MPSU-PTST1-M50K	MPSU-PTST1-M-HV
STAT3, Phospho-Y705 + STAT3, Total	MPSU-PTST3-K500	MPSU-PTST3-K10K	MPSU-PTST3-K50K	MPSU-PTST3-K-HV
STAT5, Phospho-Y694/699 + STAT5, Total	MPSU-PTST5-K500	MPSU-PTST5-K10K	MPSU-PTST5-K50K	MPSU-PTST5-K-HV
<b>STAT6, Phospho-Y641 + STAT6, Total</b>	<b>MPSU-PTST6-K500</b>	<b>MPSU-PTST6-K10K</b>	<b>MPSU-PTST6-K50K</b>	<b>MPSU-PTST6-K-HV</b>
SYK, Phospho-Y525/526 + SYK, Total	MPSU-PTSYK-K500	MPSU-PTSYK-K10K	MPSU-PTSYK-K50K	MPSU-PTSYK-K-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra Terbium Multiplex			
	500 Assay Points	10,000 Assay Points	50,000 Assay Points	100 Assay Points
4E-BP1, Phospho-T37/46	TBSU-P4EBP-A500	TBSU-P4EBP-A10K	TBSU-P4EBP-A50K	TBSU-P4EBP-A-HV
AKT1/2/3, Phospho-T308	TBSU-PAKT-A500	TBSU-PAKT-A10K	TBSU-PAKT-A50K	TBSU-PAKT-A-HV
AKT1/2/3, Phospho-S473	TBSU-PAKT-B500	TBSU-PAKT-B10K	TBSU-PAKT-B50K	TBSU-PAKT-B-HV
Cofilin, Total	TBSU-TCOF-A500	TBSU-TCOF-A10K	TBSU-TCOF-A50K	TBSU-TCOF-A-HV
CREB, Phospho-S133	TBSU-PCREB-A500	TBSU-PCREB-A10K	TBSU-PCREB-A50K	TBSU-PCREB-A-HV
EGF Receptor, Phospho-Y1068	TBSU-PEGFR-A500	TBSU-PEGFR-A10K	TBSU-PEGFR-A50K	TBSU-PEGFR-A-HV
eIF2 $\alpha$ , Phospho-S51	TBSU-PEIF2-B500	TBSU-PEIF2-B10K	TBSU-PEIF2-B50K	TBSU-PEIF2-B-HV
eIF4E, Phospho-S209	TBSU-PEIF4-A500	TBSU-PEIF4-A10K	TBSU-PEIF4-A50K	TBSU-PEIF4-A-HV
ERK1/2, Phospho-T202/Y204	TBSU-PERK-A500	TBSU-PERK-A10K	TBSU-PERK-A50K	TBSU-PERK-A-HV
GSK3 $\beta$ , Phospho-S9	TBSU-PGS3B-A500	TBSU-PGS3B-A10K	TBSU-PGS3B-A50K	TBSU-PGS3B-A-HV
IGF-1 Receptor $\beta$ , Phospho-Y1135/1136	TBSU-PIGFR-B500	TBSU-PIGFR-B10K	TBSU-PIGFR-B50K	TBSU-PIGFR-B-HV
IKK $\alpha$ , Phospho-S176/180	TBSU-PIKKA-A500	TBSU-PIKKA-A10K	TBSU-PIKKA-A50K	TBSU-PIKKA-A-HV
Insulin Receptor $\beta$ , Phospho-Y1150/1151	TBSU-PINR-A500	TBSU-PINR-A10K	TBSU-PINR-A50K	TBSU-PINR-A-HV
JNK1/2/3, Phospho-T183/Y185	TBSU-PJNK-A500	TBSU-PJNK-A10K	TBSU-PJNK-A50K	TBSU-PJNK-A-HV
MEK1, Phospho-S218/222	TBSU-PMEK1-A500	TBSU-PMEK1-A10K	TBSU-PMEK1-A50K	TBSU-PMEK1-A-HV
mTOR, Phospho-S2448	TBSU-PMTOR-A500	TBSU-PMTOR-A10K	TBSU-PMTOR-A50K	TBSU-PMTOR-A-HV
NF- $\kappa$ B p65, Phospho-S536	TBSU-PNFKB-A500	TBSU-PNFKB-A10K	TBSU-PNFKB-A50K	TBSU-PNFKB-A-HV
p38 MAPK, Phospho-T180/Y182	TBSU-PP38-B500	TBSU-PP38-B10K	TBSU-PP38-B50K	TBSU-PP38-B-HV
p70 S6K, Phospho-T389	TBSU-PP70-A500	TBSU-PP70-A10K	TBSU-PP70-A50K	TBSU-PP70-A-HV
Ribosomal Protein S6, Phospho-S240/244	TBSU-PS6R-A500	TBSU-PS6R-A10K	TBSU-PS6R-A50K	TBSU-PS6R-A-HV
SLP-76, Phospho-S376	TBSU-PSLP-A500	TBSU-PSLP-A10K	TBSU-PSLP-A50K	TBSU-PSLP-A-HV
SMAD1, Phospho-S463/465	TBSU-PSM1-A500	TBSU-PSM1-A10K	TBSU-PSM1-A50K	TBSU-PSM1-A-HV
SMAD3, Phospho-S423/425	TBSU-PSM3-A500	TBSU-PSM3-A10K	TBSU-PSM3-A50K	TBSU-PSM3-A-HV
STAT1, Phospho-S727	TBSU-PST1-B500	TBSU-PST1-B10K	TBSU-PST1-B50K	TBSU-PST1-B-HV
STAT1, Phospho-Y701	TBSU-PST1-A500	TBSU-PST1-A10K	TBSU-PST1-A50K	TBSU-PST1-A-HV
STAT3, Phospho-Y705	TBSU-PST3-A500	TBSU-PST3-A10K	TBSU-PST3-A50K	TBSU-PST3-A-HV

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

## Ordering information

Analyte	AlphaLISA SureFire Ultra Terbium Multiplex			
	500 Assay Points	10,000 Assay Points	50,000 Assay Points	100 Assay Points
STAT4, Phospho-Y693	TBSU-PST4-A500	TBSU-PST4-A10K	TBSU-PST4-A50K	TBSU-PST4-A-HV
STAT5, Phospho-Y694/699	TBSU-PST5-B500	TBSU-PST5-B10K	TBSU-PST5-B50K	TBSU-PST5-B-HV
STAT6, Phospho-Y641	TBSU-PST6-A500	TBSU-PST6-A10K	TBSU-PST6-A50K	TBSU-PST6-A-HV
SYK, Phospho-Y525/526	TBSU-PSYK-A500	TBSU-PSYK-A10K	TBSU-PSYK-A50K	TBSU-PSYK-A-HV
VEGF Receptor 2, Phospho-Y1175	TBSU-PVGFR-A500	TBSU-PVGFR-A10K	TBSU-PVGFR-A50K	TBSU-PVGFR-A-HV

### 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-0.06 mL	ALSU-ASDB-1.2 mL	ALSU-ASDB-6 mL
AlphaLISA SureFire Ultra CaptSure™ Acceptor Beads (2 mg/mL)	ALSU-ACAB-0.06 mL	ALSU-ACAB-1.2mL	ALSU-ACAB-6 mL

### 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
AlphaLISA SureFire Ultra Lysis Buffer (5X)	ALSU-LB-10 mL	ALSU-LB-100 mL
AlphaLISA SureFire Ultra Lysis Buffer B (5X)	ALSU-LBB-10ML	ALSU-LBB-100ML
AlphaLISA SureFire Ultra Lysis Buffer C (5X)	ALSU-LBC-10ML	ALSU-LBC-100ML
AlphaLISA SureFire Ultra Activation Buffer	ALSU-AB-10 mL	ALSU-AB-100 mL
AlphaLISA SureFire Ultra Activation Buffer B	ALSU-ABB-10 mL	ALSU-ABB-100 mL
AlphaLISA SureFire Ultra Activation Buffer C	ALSU-ABC-10 mL	ALSU-ABC-100 mL
AlphaLISA SureFire Ultra Dilution Buffer	ALSU-DB-10 mL	ALSU-DB-100 mL

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

## Ordering information

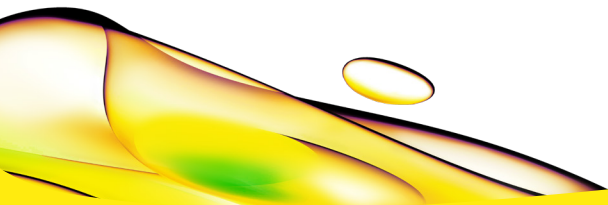
### Alpha SureFire Ultra Multiplex supplementary beads

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 $\mu$ L)	10,000 Data Points (1.2 mL)	50,000 Data Points (6 mL)
<b>SureFire Ultra MultiPlex Tb (545) CaptSure Acceptor Beads (2 mg/mL)</b>	MPSU-ACAT-0.06 mL	MPSU-ACAT-1.2 mL	MPSU-ACAT-6 mL
SureFire Ultra MultiPlex Tb (545) CaptSure 2 Acceptor Beads (2 mg/mL)	MPSU-CS2B-0.06 mL	MPSU-CS2B-1.2 mL	MPSU-CS2B-6 mL

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

SureFire™ is a registered trademark of TGR BioSciences Pty Ltd, Australia. SureFire® Ultra™ and CaptSure™ are trademarks of TGR BioSciences Pty Ltd, Australia.



revvity