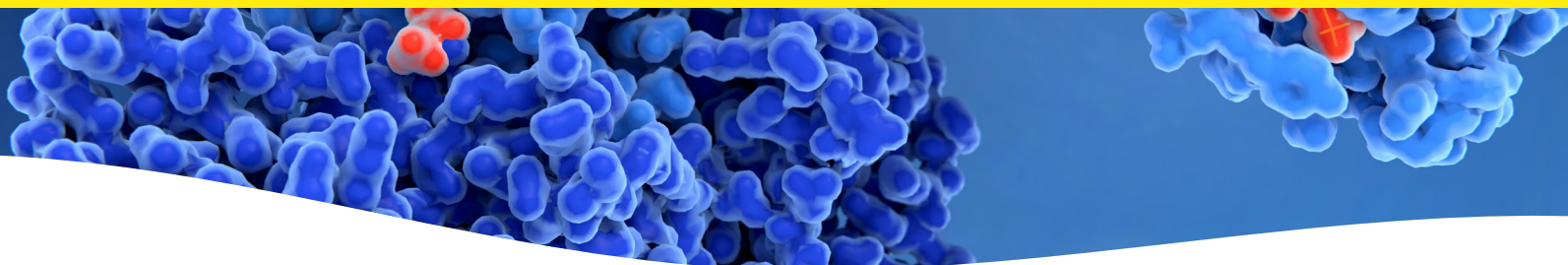




## Alpha SureFire technology assays for phosphorylated targets.



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)

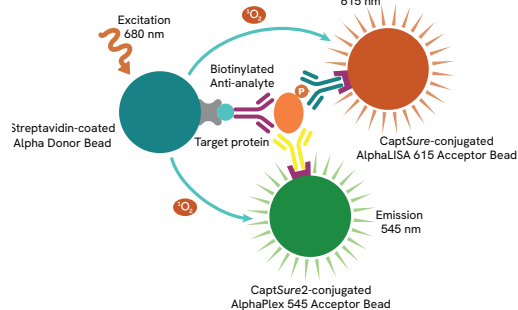
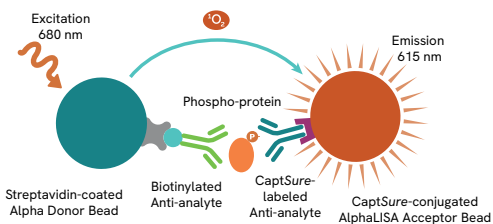
For single target analysis.

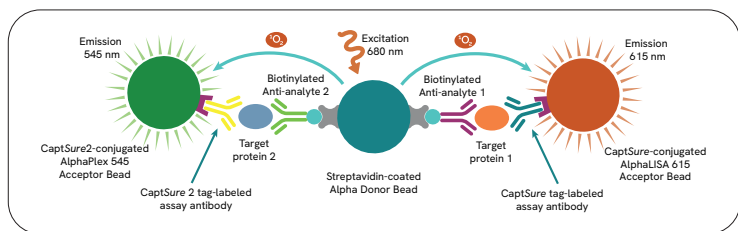
- Lower background signal, improved sensitivity, and wider dynamic range than traditional phospho-protein detection
- No interference from extraneous antibodies; ideal for therapeutic antibody screening
- Kits include all reagents

#### 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

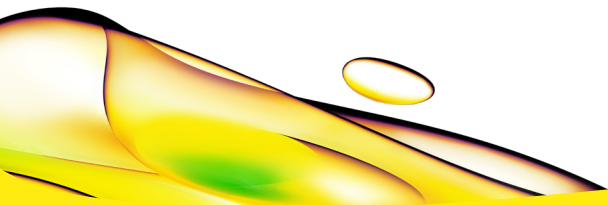
### Alpha SureFire Ultra Multiplex supplementary beads

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 µL)	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