



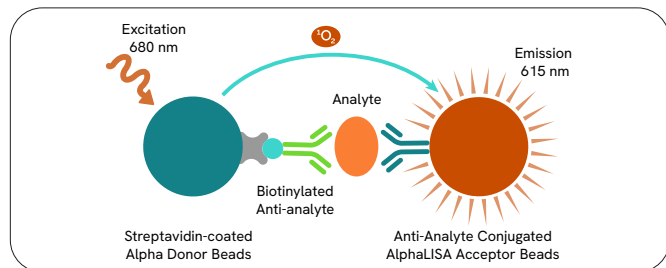
New products to meet your research and drug discovery needs.

AlphaLISA and HTRF as your science evolves, our product offering expands!

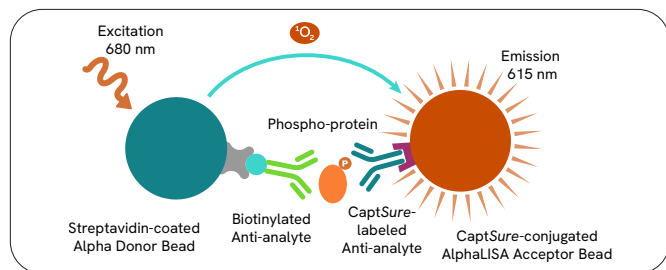
Our R&D scientists have been working tirelessly to bring you fully validated and ready-to-use immunoassay Kits for the new targets you need most.

AlphaLISA and HTRF no-wash technologies provide a larger dynamic range, improved sensitivity, and higher signal-to-noise ratios than traditional wash-based ELISAs.

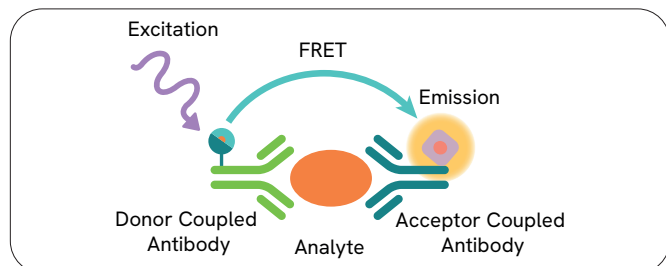
We are pleased to announce our newest AlphaLISA and HTRF assays. Benefit from Revvity's broad offering of reagent technologies and ditch your ELISA wash steps for good!



| AlphaLISA™ technology



| AlphaLISA™ SureFire® Ultra™ technology



| HTRF™ technology

Ordering information – July 2024

| Cellular phospho-protein assays | |
|---|------------------|
| AlphaLISA™ SureFire® Ultra™ products | Part – 500 tests |
| MYD88, Total Kit | ALSU-TMYD-A500 |
| Raf-1, phospho-S259 Kit | ALSU-PRAF-B500 |
| SMARCA2, Total Kit | ALSU-TSMCA2-A500 |
| Tau, phospho-T217 Kit | ALSU-PTAU-G500 |
| TIE2, phospho-Y992 Kit | ALSU-PTIE2-A500 |
| TIE2, Total Kit | ALSU-TTIE2-A500 |
| HTRF™ products | Part – 500 tests |
| ATF4, Total Kit | 64ATF4TPEG |
| c-Raf, phospho-S43 Kit | 64CRAFS43V2PEG |
| HPK1, Total Kit | 64HPK1TPEG |
| HSP27 phospho-S78 Kit | 64HSPS78PEG |
| HSP27, Total Kit | 64HSPTPEG |
| KRAS, Total Kit | 64KRASTPEG |
| MerTK, phospho-Y749 Kit | 64MERTK49PEG |
| MerTK, Total Kit | 64MERTKTPEG |
| STAT5, Total Kit | 64AT5TPEG |

| Protein-protein interaction / toolbox assays | |
|--|--------|
| AlphaLISA™ products | Part |
| Mab anti-DIG Donor Beads - 1 mg | AS118D |

| Cytokine & Biomarkers assays | |
|------------------------------|-------------|
| HTRF™ products | Part |
| dsRNA (IVT) Kit | 64DSRNAPEG |
| Diluent 14 (4X) | 62DL14DDE |
| HP Human IL-6 Kit | 62HIL6V2PEG |

HTRF™ products also available under 10000 tests & bulk size
AlphaLISA™ products also available under 100 – 5000 – 10000 – 50000 tests

Ordering information – April 2024

| Cellular phospho-protein assays | |
|---|------------------|
| AlphaLISA™ SureFire® Ultra™ products | Part - 500 tests |
| Aiolos, Total Kit | ALSU-TAIOL-A500 |
| ATG14, phospho-S29 Kit | ALSU-PATG14-A500 |
| ATF2, phospho-T71 Kit | ALSU-PATF2-A500 |
| ATF2, Total Kit | ALSU-TATF2-A500 |
| CDK8, Total Kit | ALSU-TCDK8-A500 |
| c-FOS, phospho-S32 Kit | ALSU-PCFOS-A500 |
| c-FOS, Total Kit | ALSU-TCFOS-A500 |
| Cyclin D2, Total Kit | ALSU-TCYCD2-A500 |
| Cyclin D3, Total Kit | ALSU-TCYCD3-A500 |
| DJ1, Total Kit | ALSU-TDJ1-A500 |
| IRF5, phospho-S446 Kit | ALSU-PIRF5-A500 |
| IRF5, Total Kit | ALSU-TIRF5-A500 |
| KRAS, Total Kit | ALSU-TKRAS-A500 |
| PTEN, phospho-T366 Kit | ALSU-PPTEN-A500 |
| PTEN, Total Kit | ALSU-TPTEN-A500 |
| TFE3, phospho-S321 Kit | ALSU-PTFE3-A500 |
| TFE3, Total Kit | ALSU-TTFE3-A500 |
| TYK2, phospho-Y1054/1055 Kit | ALSU-PTYK2-A500 |
| TYK2, Total Kit | ALSU-TTYK2-A500 |
| XBP1s, Total Kit | ALSU-TXBP1-A500 |
| Alpha SureFire® Ultra™ Multiplex products | Part - 500 tests |
| STAT6, phospho-Y641 + STAT6, Total Kit | MPSU-PTST6-K500 |
| HTRF™ products | Part - 500 tests |
| ALK, Phospho-Y1604 Kit | 64ALKY04PEG |
| ALK, Total Kit | 64ALKTPEG |
| Myd88, Total Kit | 64MYD88PEG |
| DDR1, Phospho-Pan-Y Kit | 64DDR1YPEG |
| DDR1, Phospho-Y513 Kit | 64DDR1Y13PEG |
| DDR1, Phospho-Y796 Kit | 64DDR1Y96PEG |
| DDR1, Total Kit | 64DDR1TPEG |

| Protein-protein interaction assays | |
|--|---------------|
| AlphaLISA™ products | Part |
| Alpha TE Lysis Buffer 1, 100 mL | TE-BUF1-100ML |
| Alpha TE Lysis Buffer 2, 100 mL | TE-BUF2-100ML |
| Alpha TE Lysis Buffer 3, 100 mL | TE-BUF3-100ML |
| Alpha TE Lysis Buffer 4, 100 mL | TE-BUF4-100ML |
| Alpha TE Lysis Buffer 5, 100 mL | TE-BUF5-100ML |
| Alpha TE Lysis Buffer Selection Kit | TE-BUF-KIT |
| HTRF™ products | Part |
| Mab Anti Flag-Eu | 61FGBKLA |
| PPI Europium Low Ionic Strength detection buffer | 61DB11RDF |

| Cytokine & Biomarkers assays | |
|------------------------------|-----------|
| HTRF™ products | Part |
| MMP1 Kit | 62MMP1PEG |
| Apolipoprotein B Kit | 64APOBPEG |

HTRF™ products also available under 10000 tests & bulk size
AlphaLISA™ products also available under 100 – 5000 – 10000 – 50000 tests

Ordering information – January 2024

| Cellular phospho-proteins assays | |
|---|-------------------|
| AlphaLISA™ SureFire® Ultra™ products | Part - 500 tests |
| AKT1, phospho-S129 Kit | ALSU-PAKT1-D500 |
| CDC25C, phospho-S216 Kit | ALSU-PCDC25-A500 |
| CDC25C, Total Kit | ALSU-TCDC25-A500 |
| Cyclin A2, Total Kit | ALSU-TCYCA2-A500 |
| Mouse DAP12, phospho-Y92 Kit | ALSU-PDAP12-C500 |
| Mouse DAP12, Total Kit | ALSU-TDAP12-B500 |
| RAB29, phospho-S72 Kit | ALSU-PRAB29-A500 |
| RAB29, Total Kit | ALSU-TRAB29-A500 |
| Mouse TFEB, phospho-S122 Kit | ALSU-PTFEB-C500 |
| Mouse TFEB, phospho-S211 Kit | ALSU-PTFEB-D500 |
| Mouse TFEB, Total Kit | ALSU-TTFEB-B500 |
| Multiplex DAP12 phospho-Y91/ Total DAP12 Kit | MPSU-PTDAP12-K500 |
| HTRF™ products | Part - 500 tests |
| Androgen Receptor, Full-length Kit | 64ARFLPEG |
| Androgen Receptor, Variant 7 Kit | 64ARV7PEG |
| Cyclin B1, phospho-S126 Kit | 64CYCB1S1PEG |
| Cyclin B1, Total Kit | 64CYCB1TPEG |
| IDO1, Total Kit | 64IDO1TPEG |
| IKZF1, Total Kit | 64IKZF1TPEG |
| IKZF3, Total Kit | 64IKZF3TPEG |
| LC3B II, Kit | 64LC3B2PEG |
| MSH3, Total Kit | 64MSH3TPEG |
| VAV1, phospho-Tyr174 Kit | 64VAV1Y174PEG |
| VAV1, Total Kit | 64VAV1TPEG |

| Protein-protein interaction assays | |
|------------------------------------|-----------|
| HTRF™ products | Part |
| Anti-VHH-d2 Kit | 610VHHDLA |
| Anti-VHH XL665 Kit | 610VHHXLA |

| Cytokine and biomarkers assays | |
|--------------------------------|--------------|
| AlphaLISA™ products | Part |
| GFAP Kit | AL3185C |
| Mutant Ataxin-2 Kit | AL3197C |
| HEK293 Host Cell Protein Kit | AL3198C |
| Rat IFN γ Kit | AL3195C |
| IL-15 Kit | AL3196C |
| HTRF™ products | Part |
| Mutant Human Ataxin 2 Kit | 64ATA2MPEG |
| IL-2 High Performance Kit | 62HIL02V2PEG |
| IL-15 Kit | 62HIL15PEG |

HTRF™ products also available under 10000 tests and bulk size
AlphaLISA™ products also available under 100 - 5000 - 10000 - 50000 tests.

Ordering information – October 2023

| Cellular phospho-proteins assays | |
|---|------------------|
| AlphaLISA™ SureFire® Ultra™ products | Part - 500 tests |
| BCL-2, phospho-S70 Kit | ALSU-PBCL2-A500 |
| BCL-2, Total Kit | ALSU-TBCL2-A500 |
| BCL-XL, Total Kit | ALSU-TBCLXL-A500 |
| DAP12, Aggregate Kit | ALSU-ADAP12-A500 |
| DAP12, phospho-Y91 Kit | ALSU-PDAP12-B500 |
| IKAROS, Total Kit | ALSU-TIKARO-A500 |
| Lamin A/C, phospho-S22 Kit | ALSU-PLAM-A500 |
| Lamin A/C, Total Kit | ALSU-TLAM-A500 |
| Parkin, phospho-S65 Kit | ALSU-PPARK-A500 |
| Parkin, Total Kit | ALSU-TPARK-A500 |
| RAB12, phospho-S106 Kit | ALSU-PRAB12-A500 |
| RAPTOR, Total Kit | ALSU-TRAPT-A500 |
| TAK1, phospho-T184/187 Kit | ALSU-PTAK-A500 |
| TAK1, Total Kit | ALSU-TTAK-A500 |
| TREM2, Aggregate Kit | ALSU-ATREM2-A500 |
| TREM2, Total Kit | ALSU-TTREM2-A500 |
| TREM2/DAP12 Complex Kit | ALSU-TTRMDP-A500 |
| WDR5, Total Kit | ALSU-TWDR-A500 |
| Wee1, Total Kit | ALSU-TWEE-A500 |
| HTRF™ products | Part - 500 tests |
| Cofilin, phospho-S3 Kit | 64COFS3PEG |
| Cofilin, Total Kit | 64COFTPEG |
| CREBBP, Total Kit | 64CREBBPTPEG |
| GSPT1, Total Kit | 64GSPT1TPEG |
| HDAC5, phospho-S259 Kit | 64HDAC5S9PEG |
| HDAC5, Total Kit | 64HDAC5TPEG |
| p300, Total Kit | 64P300TPEG |
| PLK1, phospho-T210 Kit | 64PLKT2PEG |
| PLK1, Total Kit | 64PLKTPEG |

| Protein-protein interaction assays | |
|------------------------------------|--------------|
| AlphaLISA™ products | Part |
| Anti-VHH Donor Beads – 1mg | AS117D |
| Anti-VHH Acceptor Beads – 250 µg | AL180C |
| KEAP1 Binding Kit – 500 tests | AL3194C |
| HTRF™ products | Part |
| KEAP1 Binding Kit – 500 tests | 64BDKEAP1PEG |
| KEAP1 Red Ligand – 1 unit | 64BDKEAP1RED |

| Cytokine assays | |
|-------------------------------|------------------|
| HTRF™ products | Part - 500 tests |
| IL1-beta High Performance Kit | 62HIL1B2PEG* |

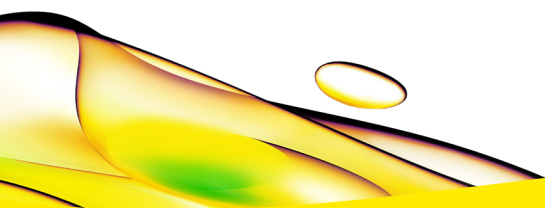
HTRF™ products also available under 10000 tests & bulk size
AlphaLISA™ products also available under 100 - 5000 - 10000 - 50000 tests

Ordering information – July 2023

| Cellular phospho-proteins assays | |
|---|-------------------------|
| AlphaLISA™ SureFire® Ultra™ products | Part - 500 tests |
| CREBBP, Total Kit | ALSU-TCRBP-A |
| IRAK4, Phospho- T345/S346 Kit | ALSU-PIRAK-A |
| MDM2, Phospho-S166 Kit | ALSU-PMDM2-A |
| MDM2, Total Kit | ALSU-TMDM2-A |
| NRF2, Phospho-S40 Kit | ALSU-PNRF2-A |
| NRF2, Total Kit | ALSU-TNRF2-A |
| p27 KIP1, Phospho-S10 Kit | ALSU-PP27K-A |
| p27 KIP1, Total Kit | ALSU-TP27K-A |
| p300, Total Kit | ALSU-TP300-A |
| PINK1, Total Kit | ALSU-TPINK-A |
| SYK, Total Kit (Mouse) | ALSU-TSYK-B |
| HTRF™ products | Part - 500 tests |
| Alpha-Synuclein, Phospho-S129 Aggregation Kit | 64SYN29APEG |
| CDK5, Total Kit | 64CDK5TPEG |
| CDK8, Total Kit | 64CDK8TPEG |
| ULK1, Phospho-S556 Kit | 64ULK556PEG |
| ULK1, Total Kit | 64ULK1TPEG |
| Wee1, Total Kit | 64WEE1TPEG |

| Protein-protein interaction assays | |
|------------------------------------|--------------------------|
| HTRF™ products | Part - 5000 tests |
| Anti-VHH Eu-cryptate | 610VHHKLA |
| Anti-VHH Tb-cryptate | 610VHHTLA |

HTRF™ products also available under 10000 tests & bulk size
AlphaLISA™ products also available under 100 – 5000 – 10000 – 50000 tests



revvity