



Physiologically relevant results for fast flowing research.



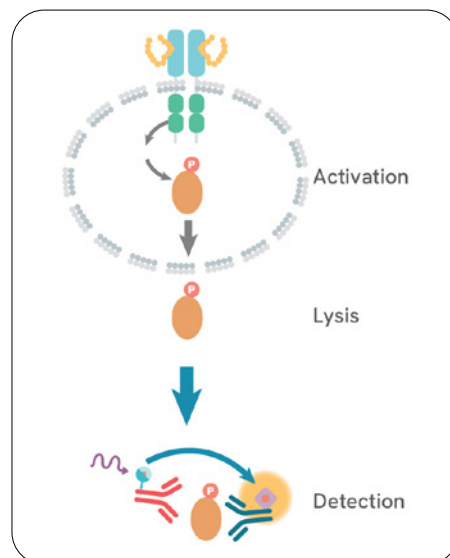
Time may be money, but relevant results are golden

Revvity's smart pathway readout tools simplify and accelerate your research projects, from basic biological investigations to high-throughput drug screening. They are ideal for both small molecule and biotherapeutic pharmacology.

Based on HTRF®, Revvity's unique Homogeneous Time- Resolved Fluorescence detection platform, this no-wash cell-based system offers add-and-read assays for physiologically relevant results, ideal for studying signaling pathways in the cell.

Progress faster, with more relevant cell signaling results using cell-based HTRF phospho-assays.

- Greater specificity and high Z' scores for robust, reproducible results
- Fewer false negatives and repeat runs
- Adapts easily to all throughput requirements
- Simple wash free add-and-read protocol
- High species cross-reactivity



Ordering information

| ANALYTE | HTRF |
|------------------------------------------|--------------|
| 4EBP1 phospho T37/46 kit | 64BP1PEG |
| ACC phospho S79 kit | 64ACCPEG |
| ACC total kit | 64NACPEG |
| ADAR1 total kit | 64ADAR1TPEG |
| AKT phospho S473 kit | 64AKSPEG |
| AKT phospho T308 kit | 64AKTPEG |
| AKT total kit | 64NKTPEG |
| AKT1 phospho S473 kit | 63ADK078PEG |
| AKT1 total kit | 63ADK079PEG |
| AKT2 phospho S473 kit | 63ADK080PEG |
| AKT2 total kit | 63ADK081PEG |
| AKT3 phospho S473 kit | 63ADK082PEG |
| AKT3 total kit | 63ADK083PEG |
| AMPK phospho T172 kit | 64MPKPEG |
| AMPK total kit | 63ADK060PEG |
| Androgen Receptor total kit | 64ANDRPEG |
| a-synuclein phospho S129 aggregation kit | 64SYN29AGPEG |
| a-synuclein phospho S129 kit | 6FSYNPEG |
| a-synuclein total kit | 6FNSYPEG |
| ATF2 phospho T71 kit | 63ADK015PEG |
| ATG14 phospho S29 kit | 64ATG14S9PEG |
| ATG14 total kit | 64ATG14TPEG |
| ATG16L1 phospho S278 kit | 64ATG16S8PEG |
| ATG16L1 total kit | 64ATG16TPEG |
| ATM phospho S1981 kit | 64ATMS1PEG |
| ATM total kit | 64ATMTPEG |
| Aurora A phospho T288 kit | 64AURAT2PEG |
| Aurora A total kit | 64AURATPEG |
| Aurora B phospho T232 kit | 64AURBT2PEG |
| Aurora B total kit | 64AURBTPEG |
| BAD phospho S112 kit | 64BADPEG |
| b-Catenin phospho T41/ S33/37 kit | 64CATPEG |
| b-Catenin total kit | 64NCAPEG |
| BRD2 total kit | 64BRD2TPEG |
| BRD3 total kit | 64BRD3TPEG |
| BRD4 total kit | 64BRD4TPEG |

| ANALYTE | HTRF |
|----------------------------|--------------|
| BRD7 total kit | 64BRD7TPEG |
| BRD9 total kit | 64BRD9TPEG |
| BTK phospho Y223 kit | 63ADK017PEG |
| BTK total kit | 63ADK064PEG |
| CDK1 phospho T14 kit | 64CDK1T4PEG |
| CDK1 phospho T15 kit | 64CDK1Y5PEG |
| CDK1 total kit | 64CDK1TPEG |
| CDK12 total kit | 64CDK12TPEG |
| CDK2 phospho Y15 kit | 64CDK2Y15PEG |
| CDK2 total kit | 64CDK2TPEG |
| CDK4 total kit | 64CDK4TPEG |
| CDK5 total kit | 64CDK5TPEG |
| CDK6 total kit | 64CDK6TPEG |
| CDK7 total kit | 64CDK7TPEG |
| CDK8 total kit | 64CDK8TPEG |
| CDK9 total kit | 64CDK9TPEG |
| Cereblon total kit | 64CRBNPEG |
| cGAS total kit | 64CGASTPEG |
| CHK1 phospho S345 kit | 64CHK1S5PEG |
| CHK1 total kit | 64CHK1TPEG |
| CHK2 phospho T68 kit | 64CHK2T8PEG |
| CHK2 total kit | 64CHK2TPEG |
| c-Jun phospho S63 kit | 64JUNPEG |
| COFILIN phospho S3 kit | 64COFS3PEG |
| COFILIN total kit | 64COFTPEG |
| c-Raf phospho S338 kit | 63ADK018PEG |
| c-RAF phospho S43 kit | 64CRAFS43PEG |
| c-RAF total kit | 64CRAFTPEG |
| CREB phospho S133 kit | 64CREPEG |
| CREB total kit | 63ADK052PEG |
| CREBBP total kit | 64CREBBPTPEG |
| c-Src phospho Y419 kit | 64SRCPEG |
| c-Src total kit | 64NSRPEG |
| Cyclin D1 phospho T286 kit | 64CYCD1T6PEG |
| Cyclin D1 total kit | 64CYCD1TPEG |
| Cyclin E1 total kit | 64CYCE1TPEG |

Ordering information (continued)

| ANALYTE | HTRF |
|-------------------------------------|--------------|
| EGFR DEL19 total kit | 64EGFRD19PEG |
| EGFR L858R total kit | 64EGFRL8PEG |
| EGFR phospho pan kit | 64HR1PEG |
| EGFR phospho Y1068 kit | 64EG1PEG |
| EGFR total kit | 64NG1PEG |
| EIF2a phospho S52 kit | 64EF2PEG |
| EIF2a total kit | 64NEFPEG |
| EIF4E phospho S209 kit | 64EF4PEG |
| eNOS phospho S1177 kit | 64ENOSS7PEG |
| eNOS total kit | 64ENOSTPEG |
| ErbB2 (HER2) phospho Y1221/1222 kit | 64HR2PEG |
| ErbB2 (HER2) total kit | 64NR2PEG |
| ERK phospho T202/Y204 kit | 64AERPEG |
| ERK total kit | 64NRKPEG |
| FAK phospho Y397 kit | 64FAKPEG |
| FAK total kit | 64NAKPEG |
| FGFR1 phospho Y653/654 kit | 64FGFR1Y6PEG |
| FGFR1 total kit | 64FGFR1TPEG |
| FGFR2 phospho Y653/654 kit | 63ADK054PEG |
| FGFR2 total kit | 63ADK057PEG |
| FGFR3 phospho Y653/654 kit | 64FGFR3Y6PEG |
| FGFR3 total kit | 64FGFR3TPEG |
| FGFR4 phospho Y642 kit | 64FGFR4Y6PEG |
| FGFR4 total kit | 64FGFR4TPEG |
| FLT3 phospho Y589/591 kit | 64FLT3Y1PEG |
| FLT3 phospho Y842 kit | 64FLT3Y2PEG |
| FLT3 total kit | 64FLT3TPEG |
| FoxO1 phospho S256 kit | 64FOXPEG |
| FoxO1 total kit | 64NFOPEG |
| GAPDH housekeeping kit | 64GAPDHPEG |
| GSK3a phospho S21 kit | 64GPAPEG |
| GSK3a total kit | 64GTAPEG |
| GSK3b phospho S9 kit | 64GPBPEG |
| GSK3b total kit | 64GTBPEG |
| GSPT1 total kit | 64GSPT1TPEG |
| H2AX phospho S139 kit | 64H2XPEG |

| ANALYTE | HTRF |
|---------------------------------|--------------|
| H2AX total kit | 64H2AXTPEG |
| HDAC4 phospho S246 kit | 64HDAC4S6PEG |
| HDAC4 total kit | 64HDAC4TPEG |
| HDAC5 phospho S259 kit | 64HDAC5S9PEG |
| HDAC5 total kit | 64HDAC5TPEG |
| HER3 phospho Y1289 kit | 64HR3PEG |
| HER3 total kit | 64NR3PEG |
| Histone H3 phospho S10 kit | 64HH3PEG |
| Histone H3 phospho T3 kit | 63ADK061PEG |
| Histone H3 total kit | 64NH3PEG |
| HTT total kit | 64HTTTPEG |
| IKKA phospho S176/180 kit | 64KKAS6PEG |
| IKKA total kit | 64KKATPEG |
| IKKB phospho S177/181 kit | 64KKBPEG |
| IKKB total kit | 63ADK097PEG |
| IR beta phospho Y1150 /1151 kit | 63ADK016PEG |
| IRAK3 total kit | 63ADK101PEG |
| IRAK4 total kit | 63ADK108PEG |
| IRb total kit | 63ADK019PEG |
| IRF3 phospho S385/386 kit | 6FRF3PEG |
| IRF5 total kit | 64HIRF5PEG |
| IRS1 phospho Ser312 kit | 64IRSPEG |
| IRS1 total kit | 64NRSPEG |
| IRS2 total kit | 64IRTPEG |
| JAK1 phospho Y1034/1035 kit | 64JAK1Y10PEG |
| JAK1 total kit | 64JAK1TPEG |
| JAK2 phospho Y1007/1008 kit | 64JAK2Y10PEG |
| JAK2 total kit | 64JAK2TPEG |
| JAK3 total kit | 64JAK3TPEG |
| JNK phospho T183/Y185 kit | 64JNKPEG |
| KRAS total kit | 64KRASTPEG |
| LRRK2 phospho S935 kit | 6FLRKPEG |
| LRRK2 total kit | 6FNRKPEG |
| MEK1 phospho S218/222 kit | 64ME1PEG |
| MEK1 phospho S298 kit | 63ADK021PEG |
| MEK1 total kit | 64NE1PEG |

Ordering information (continued)

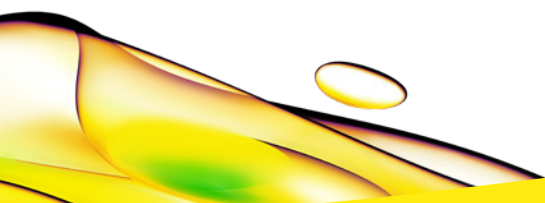
| ANALYTE | HTRF |
|-------------------------------------------------|--------------|
| MEK1/2 phospho S218/222 kit | 64ME2PEG |
| MET phospho T1234/1235 kit | 64METY34PEG |
| MET total kit | 64METTPEG |
| MKK4 phospho S257 kit | 64MK4PEG |
| MKK4 total kit | 64NK4PEG |
| MLKL phospho S358 kit | 64MLKLS8PEG |
| MLKL total kit | 64MLKLTPEG |
| MNK phospho T197/202 kit | 63ADK022PEG |
| mTOR phospho S2448 kit | 64TORPEG |
| NDRG1 phospho S330 kit | 63ADK093PEG |
| NDRG1 phospho T346 kit | 63ADK092PEG |
| NDRG1 total kit | 63ADK094PEG |
| NFκB phospho S536 kit | 64NFBPEG |
| NFκB total kit | 64NFTPEG |
| NRF2 total kit | 64NRF2TPEG |
| p21 total kit | 64HP21PEG |
| p300 total kit | 64P300TPEG |
| P38 MAPK phospho T180/Y182 kit | 64P38PEG |
| p53 phospho S15 kit | 64P53PEG |
| P53 total kit | 64T53PEG |
| p62 phospho S403 kit | 64P62S4PEG |
| p62 total kit | 64N62PEG |
| p70 S6K phospho T389 kit | 64S6KPEG |
| PDGF receptor beta (PDGFRβ) phospho Y751 kit | 63ADK040PEG |
| Phospholamban phospho S16 kit | 64PLN16PEG |
| Phospholamban phospho T17 kit | 63ADK074PEG |
| Phospholamban total kit | 63ADK075PEG |
| PINK1 total kit | 64PINKTPEG |
| PLCg1 phospho Y783 kit | 64PLCG1Y3PEG |
| PLCg1 total kit | 64PLCG1TPEG |
| PLCg2 phospho Y1217 kit | 64PLCG2Y7PEG |
| PLCg2 total kit | 64PLCG2TPEG |
| PLK1 phospho T210 kit | 64PLKT2PEG |
| PLK1 total kit | 64PLKTPEG |
| PR total kit | 64HPRTPPEG |
| pRB Ser780 phospho S780 kit | 63ADK104PEG |

| ANALYTE | HTRF |
|----------------------------|--------------|
| RAB10 phospho T73 kit | 64RABT73PEG |
| RAB10 total kit | 64RAB10TPEG |
| Rb phospho S780 kit | 64RBS780PEG |
| Rb phospho S807/811 kit | 63ADK105PEG |
| RB total kit | 63ADK106PEG |
| RBM39 total kit | 64RBM39TPEG |
| RET phospho pan kit | 64RETYPEG |
| RET total kit | 64RETTPEG |
| RIPK1 phospho S166 kit | 64RIPK1S1PEG |
| RIPK1 total kit | 64RIPK1TPEG |
| RSK1 phospho S380 kit | 64RSK1S0PEG |
| RSK1 phospho T359/S363 kit | 64RSK1T9PEG |
| RSK1 total kit | 64RSK1TPEG |
| S6RP phospho S235/236 kit | 64RP6PEG |
| SHP1 phospho Y564 kit | 64SH1PEG |
| SHP1 total kit | 64NH1PEG |
| SHP2 phospho Y542 kit | 64SH2PEG |
| SHP2 total kit | 64NH2PEG |
| SLP76 phospho S376 kit | 63ADK076PEG |
| SLP76 phospho Y145 kit | 64SLPY5PEG |
| SLP76 total kit | 63ADK077PEG |
| SMAD1 phospho S463/465 kit | 63ADK062PEG |
| SMAD1 total kit | 63ADK063PEG |
| SMAD2 phospho S465/467 kit | 63ADK102PEG |
| SMAD2 total kit | 63ADK103PEG |
| SMAD3 phospho S423/425 kit | 63ADK025PEG |
| SMAD3 total kit | 64ND3PEG |
| SMAD4 phospho T277 kit | 64SMAD4T7PEG |
| SMAD4 total kit | 64SMAD4TPEG |
| SMARCA-2 total kit | 64SMARC2TPEG |
| SMARCA-4 total kit | 64SMARC4TPEG |
| STAT1 phospho Y701 kit | 63ADK026PEG |
| STAT1 total kit | 63ADK096PEG |
| STAT3 phospho Y705 kit | 62AT3PEG |
| STAT3 total kit | 64NT3PEG |
| STAT5 phospho Y694 kit | 64AT5PEG |

Ordering information (continued)

| ANALYTE | HTRF |
|---------------------------|--------------|
| STAT6 phospho Y641 kit | 64AT6PEG |
| STAT6 total kit | 64STAT6TPEG |
| STING phospho S365 kit | 64MSTGPEG |
| STING phospho S366 kit | 64STGPEG |
| STING total kit | 64NTGPEG |
| SYK phospho Y525/526 kit | 64SYKY525PEG |
| SYK total kit | 64SYKTPEG |
| Tau phospho S181 kit | 64TT181PEG |
| Tau phospho S202/T205 kit | 64TS2PEG |
| Tau phospho S356 kit | 64TS356PEG |
| Tau phospho S404 kit | 64TS4PEG |
| Tau phospho S422 kit | 64TS422PEG |
| Tau total kit | 64NTAUPEG |
| TBK1 phospho S172 kit | 64TBKPEG |
| TBK1 total kit | 64NTBPEG |
| TDP-43 phospho S409 kit | 64TDPS4PEG |
| TDP-43 total kit | 64TDPTPEG |
| TFEB phospho S211 kit | 64TFEBS2PEG |
| TFEB total kit | 64TFEBTPEG |
| Tie2 phospho Y992 kit | 64TIE2Y2PEG |

| ANALYTE | HTRF |
|--------------------------------------------|--------------|
| TYK2 phospho Y1054/1055 kit | 64TYK2Y10PEG |
| TYK2 total kit | 64TYK2TPEG |
| Ubiquitin phospho S65 kit | 64UBIS65PEG |
| ULK1 phospho S556 kit | 64ULKS5PEG |
| ULK1 phospho S758 kit | 64ULKPEG |
| ULK1 total kit | 64ULKTPPEG |
| VASP phospho S157 kit | 63ADK066PEG |
| VASP phospho S239 kit | 63ADK065PEG |
| VASP total kit | 63ADK067PEG |
| VEGF receptor 2 (VEGFR2) phospho Y1175 kit | 63ADK041PEG |
| VEGFR2 total kit | 64VEGFR2TPEG |
| Wee1 total kit | 64WEE1TPEG |
| XBP1 total kit | 64XBPPTPEG |
| XBP1s total kit | 64XBPSTPEG |
| YAP phospho S127 kit | 64YAPPEG |
| YAP total kit | 64YATPEG |
| ZAP70 phospho Y319 kit | 64ZAPPEG |
| ZAP70 total kit | 64ZATPEG |



revvity