

Species compatibility

Cell Signaling: Biomarkers, Phospho- & total-protein assays

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILA | PIG | DOG |
|---------------------------------|-------|-------|-----|--------|---------|------------|-----|-----|
| Alpha-SMA | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Alpha-Synuclein-S129 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| Alpha-Synuclein total | ● | ✗ | ○ | ○ | ○ | ○ | ○ | ○ |
| Alpha-Tubulin Housekeeping | ● | ● | ● | ● | ● | ○ | ● | ● |
| 4EBP1 phospho-T37/46 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| ACC phospho-S79 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| AKT phospho-S473 | ● | ● | ● | ● | ● | ● | ○ | ○ |
| AKT phospho-T308 | ● | ● | ● | ● | ● | ● | ○ | ○ |
| AKT total | ● | ● | ● | ○ | ● | ○ | ○ | ○ |
| AKT1 phospho-S473 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| AKT1 total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| AKT2 phospho-S473 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| AKT2 total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| AKT3 phospho-S473 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| AKT3 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| AMPK phospho-T172 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| AMPK total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| ATF2 phospho-T71 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| BAD phospho-S112 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| Beta Catenin phospho-T41/S33/37 | ● | ● | ● | ● | ○ | ○ | ● | ○ |
| Beta Catenin total | ● | ● | ● | ● | ○ | ○ | ● | ○ |
| BTK phospho-Y223 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| BTK total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |

Species compatibility

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILA | PIG | DOG |
|--------------------------------|-------|-------|-----|--------|---------|------------|-----|-----|
| cJun phospho-S63 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| cRAF phospho-S338 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| CREB phospho-S133 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| CREB total | ● | ● | ● | ● | ○ | ● | ○ | ○ |
| EGFR phospho-Y1068 | ● | ✗ | ✗ | ● | ○ | ○ | ○ | ○ |
| EGFR total | ● | ● | ✗ | ● | ○ | ○ | ○ | ○ |
| EIF2a Phospho-S52 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| EIF2a total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| EIF4E phospho-S209 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| Advanced ERK phospho-T202/Y204 | ● | ● | ● | ● | ● | ● | ● | ● |
| ERK phospho-T202/Y204 | ● | ● | ● | ● | ● | ● | ● | ● |
| ERK total | ● | ● | ● | ● | ● | ● | ● | ● |
| FAK phospho-Y397 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| FAK total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| FoxO1 phospho-S256 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| FoxO1 total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| GSK3 alpha phospho-S21 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| GSK3 alpha total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| GSK3 beta phospho-S9 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| GSK3 beta total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| H2AX phospho-S139 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| HER2 phospho-Y1221/1222 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| HER2 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| HER3 phospho-Y1289 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| HER3 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| HIF-1 α | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |

Species compatibility

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILA | PIG | DOG |
|------------------------------------|-------|-------|-----|--------|---------|------------|-----|-----|
| Histone H3 phospho-T3 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Histone H3 phospho-S10 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| Histone H3 total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| IKKB phospho-S177/181 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| IRAK3 total | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |
| IRAK4 total NEW | ● | ✗ | ○ | ○ | ○ | ○ | ○ | ○ |
| IRF3 phospho-S386 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| IRS1 phospho-S307 mouse/S312 human | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| IRS1 total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| IRS2 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| JNK phospho-T183/Y185 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| LRRK2 phospho-S935 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| LRRK2 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| MEK1 phospho-S218/222 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| MEK1 total | ● | ● | ● | ● | ● | ○ | ○ | ○ |
| MEK1/2 phospho-S218/222 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| mTOR phospho-S2448 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| NFκB phospho-S536 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| NFκB total | ● | ● | ● | ○ | ● | ○ | ○ | ● |
| PDGFRβ phospho-Y751 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Phospholamban phospho-T17 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Phospholamban total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| p38 phospho-T180/Y182 | ● | ● | ● | ● | ● | ○ | ● | ○ |
| p53 phospho-S15 | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |
| p53 total | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |

Species compatibility

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILA | PIG | DOG |
|------------------------------------|-------|-------|-----|--------|---------|------------|-----|-----|
| p62/SQSTM1 phospho-S403 NEW | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| p62/SQSTM1 total NEW | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| p70S6K phospho-T389 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| RB phospho-S780 | ● | ● | ○ | ● | ○ | ○ | ○ | ○ |
| RB phospho-S807/811 | ● | ● | ○ | ● | ○ | ○ | ○ | ○ |
| RB total | ● | ○ | ○ | ● | ○ | ○ | ○ | ○ |
| S6RP phospho-S235/236 | ● | ● | ● | ● | ● | ○ | ○ | ○ |
| SHP1 phospho-Y564 | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |
| SHP1 total | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |
| SHP2 phospho-Y542 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| SHP2 total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| SLP-76 phospho-S376 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| SLP-76 total | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| SMAD1 phospho-S463/465 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| SMAD1 total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| SMAD2 phospho-S465/467 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| SMAD2 total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| SMAD3 phospho-S423/425 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| SMAD3 total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| c-Src phospho-Y419 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| c-Src total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| STAT1 phospho-Y701 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| STAT1 total | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| STAT3 phospho-Y705 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| STAT3 total | ● | ● | ● | ● | ● | ○ | ○ | ○ |
| STAT5 phospho-Y694 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |

Species compatibility

● Compatible / ● Predicted / ✗ Not compatible / ○ Unknown

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILA | PIG | DOG |
|------------------------------------|-------|-------|-----|--------|---------|------------|-----|-----|
| STAT6 phospho-Y641 | ● | ✗ | ○ | ○ | ○ | ○ | ○ | ○ |
| STING phospho-S366 | ● | ✗ | ○ | ○ | ○ | ○ | ○ | ○ |
| STING total | ● | ✗ | ○ | ○ | ○ | ○ | ○ | ○ |
| Tau phospho-S202/T205 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Tau phospho-S404 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| TBK1 phospho-S172 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| TBK1 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| ULK1 phospho-S757 mouse/S758 human | ● | ● | ○ | ● | ○ | ○ | ○ | ○ |
| VASP Phospho-S157 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| VASP Phospho-S239 kit | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| VASP total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| YAP phospho-S127 | ● | ● | ● | ○ | ● | ○ | ○ | ○ |
| VEGFR2 phospho-Y1175 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| YAP total | ● | ● | ● | ○ | ● | ○ | ○ | ○ |
| ZAP-70 phospho-Y319 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| ZAP-70 total | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 21 5018 9880 Japan +81-(0)43-306-8712

Visit www.cisbio.com to find a list of our regional distributors