

Species compatibility

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

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILIA | PIG | DOG |
|------------------------------------|-------|-------|-----|--------|---------|-------------|-----|-----|
| 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 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Histone H3 phospho-T3 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| Histone H3 phospho-S10 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| Histone H3 total | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| IKKB phospho-S177/181 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| IRF3 phospho-S386 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| IRS1 phospho-S307 mouse/S312 human | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| IRS1 total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| IRS2 total | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| JNK phospho-T183/Y185 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| MEK1 phospho-S218/222 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| MEK1 total | ● | ● | ● | ● | ● | ○ | ○ | ○ |
| MEK1/2 phospho-S218/222 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| mTOR phospho-S2448 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| NFkB phospho-S536 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| NFkB total | ● | ● | ● | ○ | ● | ○ | ○ | ● |

Species compatibility

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

| | HUMAN | MOUSE | RAT | MONKEY | HAMSTER | DROSOPHILIA | PIG | DOG |
|------------------------------------|-------|-------|-----|--------|---------|-------------|-----|-----|
| p38 phospho-T180/Y182 | ● | ● | ● | ● | ● | ○ | ● | ○ |
| P53 phospho-S15 | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |
| p53 total | ● | ○ | ○ | ○ | ○ | ○ | ○ | ○ |
| p70S6K phospho-T389 | ● | ● | ● | ● | ○ | ○ | ○ | ○ |
| S6RP phospho-S235/236 | ● | ● | ● | ● | ● | ○ | ○ | ○ |
| SLP-76 phospho-S376 | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| SLP-76 total | ● | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ | ✗ |
| SMAD1 phospho-S463/465 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| SMAD1 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 | ● | ● | ● | ○ | ○ | ○ | ○ | ○ |
| ULK1 phospho-S757 mouse/S758 human | ● | ● | ○ | ● | ○ | ○ | ○ | ○ |
| VASP Phospho-S157 | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| VASP Phospho-S239 kit | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| VASP total | ● | ● | ○ | ○ | ○ | ○ | ○ | ○ |
| YAP phospho-S127 | ● | ● | ● | ○ | ● | ○ | ○ | ○ |
| 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