

# 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-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	●	●	○	○	○	○	○	○
BAD phospho-S112	●	●	●	●	○	○	○	○
Beta Catenin phospho-T41/S33/37	●	●	●	●	○	○	●	○
Beta Catenin total	●	●	●	●	○	○	●	○
BTK phospho-Y223	●	●	●	○	○	○	○	○
BTK total	●	●	○	○	○	○	○	○
cJun phospho-S63	●	●	●	●	○	○	○	○
CREB phospho-S133	●	●	●	●	○	○	○	○
CREB total	●	●	●	●	○	●	○	○

## Species compatibility

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

	HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
EGFR phospho-Y1068	●	✗	✗	●	○	○	○	○
EGFR total	●	●	✗	●	○	○	○	○
EIF2a Phospho-S52	●	●	●	●	○	○	○	○
EIF2a total	●	●	●	●	○	○	○	○
EIF4E phospho-S209	●	●	●	●	○	○	○	○
Advanced ERK phospho-T202/Y204	●	●	●	●	●	●	●	●
ERK phospho-T202/Y204	●	●	●	●	●	●	●	●
ERK 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 $\alpha$	●	●	●	○	○	○	○	○
Histone H3 phospho-T3	●	●	●	○	○	○	○	○
Histone H3 phospho-S10	●	●	●	●	○	○	○	○
Histone H3 total	●	●	●	●	○	○	○	○
IKKB phospho-S177/181	●	●	●	●	○	○	○	○
IRF3 phospho-S386	●	✗	✗	✗	✗	✗	✗	✗

## Species compatibility

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

	HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
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	●	●	●	○	○	○	○	○
NFκB phospho-S536	●	●	●	○	○	○	○	○
NFκB total	●	●	●	○	●	○	○	●
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	●	●	●	●	○	○	○	○

## Species compatibility

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

	HUMAN	MOUSE	RAT	MONKEY	HAMSTER	DROSOPHILA	PIG	DOG
STAT1 phospho-Y701	●	✗	✗	✗	✗	✗	✗	✗
STAT1 total	●	✗	✗	✗	✗	✗	✗	✗
STAT3 phospho-Y705	●	●	●	●	○	○	○	○
STAT3 total	●	●	●	●	●	○	○	○
STAT5 phospho-Y694	●	●	●	○	○	○	○	○
ULK1 phosho-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](http://www.cisbio.com) to find a list of our regional distributors