

Lysis buffer compatibility

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Alpha-SMA	X	○	LB	X	○
Alpha-Synuclein phospho-S129	●	●	●	●	LB
Alpha-Synuclein total	●	●	●	●	LB
Alpha-Tubulin Housekeeping	●	●	●	LB	○
4EBP1 phospho-T37/46	LB	○	○	○	○
ACC phospho-S79	●	○	X	LB	○
ACC total	X	X	●	LB	○
AKT1 phospho-S473	LB	●	○	○	○
AKT1 total	LB	●	○	○	○
AKT2 phospho-S473	LB	●	○	○	○
AKT2 total	LB	●	○	○	○
AKT3 phospho-S473	LB	●	○	○	○
AKT3 total	LB	●	○	○	○
AKT phospho-S473 & AKT phospho-T308	LB	○	●	●	○
AKT total	LB	○	○	●	○
AMPK phospho-T172	LB	○	○	●	○
AMPK total	LB	○	○	●	○
Androgen receptor Total NEW	●	●	LB	●	○
AP2 total	●	●	●	LB	○

-  Included in the kit
-  High performance
-  Good performance
-  Low performance
-  Not suitable
-  Not determined



Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
ATF2 phospho-T71	LB	●	X	X	○
BAD phospho-S112	●	○	LB	○	○
B-arrestin 1 total	●	●	●	LB	○
B-arrestin 2 total	●	●	●	LB	○
Beta Catenin phospho-T41/S33/37	○	○	LB	●	○
Beta Catenin total	○	○	LB	●	○
BTK phospho-Y223	LB	○	○	●	○
BTK total	LB	○	○	●	○
Cereblon total	●	●	●	LB	○
CDK2 phospho-Y15	●	LB	●	●	○
CDK2 total	X	LB	●	●	○
CDK4 total	LB	X	X	●	○
CDK6 total	LB	X	X	●	○
c-Jun phospho-S63	LB	○	X	●	○
c-RAF phospho-S338		LB	○	●	○
cRaf phospho-S43	●	●	○	○	○
cRaf total	●	●	○	○	○
CREB phospho-S133	LB	○	●	X	○
CREB total	LB	○	○	●	○
EGFR phospho-pan	○	○	●	LB	○
EGFR phospho-Y1068	X	○	●	LB	○

- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
EGFR total	○	○	●	LB	○
EIF2α phospho-S52	●	○	○	LB	○
EIF2α total	●	○	○	LB	○
EIF4E phospho-S209	LB	○	X	○	○
eNOS phospho-S1177	●	●	LB	●	○
eNOS total	●	●	LB	●	○
Advanced ERK phospho-T202/Y204	LB	○	X	●	○
ERK phospho-T202/Y204	LB	○	X	●	○
ERK total	LB	○	X	●	○
FAK phospho-Y397	X	●	LB	X	○
FAK total	●	●	LB	●	○
FGFR1 phospho Y653/654	X	X	X	LB	○
FGFR1 total	●	●	●	LB	○
FGFR2 phospho-Y653/654	X	●	●	LB	○
FGFR2 total	●	●	●	LB	○
FGFR3 phospho Y653/654	X	●	X	LB	○
FGFR3 total	●	●	X	LB	○
FGFR4 phospho Y642	X	●	●	LB	○
FGFR4 total	X	●	●	LB	○
FoxO1 phospho-S256	LB	○	X	●	○
FoxO1 total	LB	○	X	●	○

-  Included in the kit
-  High performance
-  Good performance
-  Low performance
-  Not suitable
-  Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
GAPDH Housekeeping	LB	●	●	●	○
GSK3 alpha phospho-S21	●	○	●	LB	○
GSK3 alpha total	●	○	●	LB	○
GSK3 beta phospho-S9	●	○	●	LB	○
GSK3 beta total	●	○	●	LB	○
H2AX phospho-S139	●	○	○	LB	○
HER2 phospho-Y1221/1222	X	○	●	LB	○
HER2 total	○	○	●	LB	○
HER3 phospho-Y1289	●	○	○	LB	○
HER3 total	○	○	○	LB	○
HIF-1 alpha	●	○	●	LB	○
Histone H3 phospho-T4	LB	○	○	○	○
Histone H3 phospho-S10	LB	○	○	○	○
Histone H3 total	LB	○	X	○	○
Insulin receptor phospho-Y1150/1151	LB	●	○	○	○
Insulin receptor total	LB	●	○	○	○
IRAK3 total	LB	●	●	●	●
IRAK4 total	LB	●	●	●	●
IRF3 phospho-S386	X	LB	●	X	○
IRβ phospho Y1150/1151	LB	●	●	●	○
IRβ total	LB	●	●	●	○

- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
IRS1 phospho-S307 mouse / S312 human	LB	○	○	●	○
IRS1 total	LB	○	○	●	○
IRS2 total	LB	○	○	●	○
IKKB phospho-S177/181	LB	○	X	●	○
IKKB total	LB	●	●	●	○
JAK1 Phospho Y1034/35 NEW	X	●	●	LB	○
JAK1 total NEW	●	●	●	LB	○
JAK2 Phospho Y1007/08 NEW	LB	●	●	●	○
JAK2 total NEW	LB	●	●	●	○
JAK3 total NEW	●	●	●	LB	○
JNK phospho-T183/Y185	X	○	LB	○	○
Ki-67	●	●	X	LB	○
LRRK2 phospho-S9358	X	X	X	LB	○
LRRK2 total	X	X	X	LB	○
MDA5 total	●	●	LB	●	○
MDA5 aggregation	X	X	LB	X	○
MEK1 phospho-S298	LB	○	○	○	○
MEK1 phospho-S218/222	LB	○	○	○	○
MEK1 total	LB	○	○	○	○
MEK1/2 phospho-S218/222	LB	○	●	○	○
MKK4 phospho-S257	LB	●	●	●	○

-  Included in the kit
-  High performance
-  Good performance
-  Low performance
-  Not suitable
-  Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
MKK4 total	LB	●	●	●	○
MLKL phospho-S358	●	●	●	LB	○
MLKL total	●	●	●	LB	○
MNK phospho-T197/202	LB	X	○	○	○
mTOR phospho-S2448	LB	○	●	●	○
NDRG1 phospho-T346	○	○	○	LB	○
NDRG1 phospho-S330	○	○	○	LB	○
NDRG1 total	○	○	○	LB	○
NFκB phospho-S536	○	○	●	LB	○
NFκB total	○	○	●	LB	○
P38 phospho-T180/Y182	LB	○	X	X	○
P53 phospho-S15	LB	○	X	X	○
P53 total	LB	○	X	X	○
P62/SQSTM1-S403	●	●	●	LB	○
P62/SQSTM1 total	●	●	●	LB	○
p70S6K phospho-T389	LB	○	●	○	○
PDGFRβ phospho-Y751	●	●	●	LB	○
PLCγ1 phospho-Y1217	X	X	LB	X	○
PLCγ1 total	●	●	LB	●	○
PLCγ2 phospho-Y1217	●	●	●	LB	○
PLCγ2 total	●	X	X	LB	○

- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Phospholamban phospho-S16	●	●	LB ●	●	○
Phospholamban total	●	●	LB ●	●	○
RB phospho-S780	●	LB ●	●	●	○
RB phospho-S807/811	●	LB ●	●	●	○
RB total	●	LB ●	●	●	○
RET phospho-pan Y	●	●	●	LB ●	○
RET total	●	●	●	LB ●	○
RIPK1 phospho-S166	LB ●	●	●	●	○
RIPK1 total	LB ●	●	●	●	○
RSK1 phospho-S380	X	LB ●	○	X	●
RSK1 phospho-T359/S363	●	LB ●	○	○	○
RSK1 total	X	LB ●	○	○	○
S6RP phospho-S235/236	LB ●	○	●	●	○
SARS-CoV-2 capsid	●	●	●	●	LB ●
SHP1 phospho-Y564	X	X	●	LB ●	○
SHP1 total	X	●	●	LB ●	○
SHP2 phospho-Y542	X	LB ●	●	X	○
SHP2 total	X	LB ●	●	X	○
SLP76 phospho-S376	LB ●	○	○	●	○
SLP76 phospho-Y145	X	●	●	LB ●	○
SLP76 total	LB ●	●	●	●	○

- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
SLP76 phospho-S376 mouse	LB	●	○	○	○
SLP76 total mouse	LB	●	○	○	○
SMAD1 phospho-S463/465	●	○	○	LB	○
SMAD1 total	●	○	○	LB	○
SMAD2 phospho-S465/467	●	X	●	LB	○
SMAD2 total	●	●	●	LB	○
SMAD3 phospho-S463/465	●	○	●	LB	○
SMAD3 total	●	○	●	LB	○
Spike S1 protein	●	○	○	●	LB
c-Src phospho-Y419	X	○	LB	●	○
c-Src total	●	○	LB	●	○
STAT1 phospho-Y701	●	LB	●	●	○
STAT1 total	●	LB	●	●	○
STAT3 phospho-Y705	LB	○	●	●	○
STAT3 total	LB	○	○	○	○
STAT5 phospho-Y694	●	○	X	LB	○
STAT6 phospho-Y641	X	X	●	LB	○
STING phospho-S366	●	●	●	LB	○
STING total	●	●	●	LB	○
STING (m) phospho-S365	X	X	●	LB	○
STING (m) Total	●	●	●	LB	○



- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
SYK phospho-Y525	●	●	LB	●	○
SYK total	●	●	LB	●	○
TAU phospho-S202/T205	LB	●	●	●	○
TAU phospho-S356	LB	●	●	●	○
TAU phospho-S404	LB	X	●	●	○
TAU phospho-S422	LB	X	●	●	○
TAU phospho-T181	LB	●	●	●	○
TAU total	LB	●	●	●	○
TBK1 phospho-S172	●	X	X	LB	○
TBK1 total	●	X	X	LB	○
TDP43 aggregated	LB	●	●	●	○
Tie2 phospho-Y992	LB	X	X	●	●
TOM20	●	●	X	LB	○
Ubiquitin phospho-S65	LB	X	●	●	○
ULK1 phospho-S757 mouse / S758 human	LB	○	●	●	○
VASP phospho-S157	●	LB	●	●	○
VASP phospho-S239	●	LB	●	●	○
VASP total	●	LB	●	●	○
VEGFR2 phospho-Y1175	●	LB	●	●	○
VEGFR2 total	●	LB	X	●	○
YAP phospho-S127	●	○	X	LB	○

- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
ZAP-70 phospho-Y319	●	○	●		○
ZAP-70 total	●	○	●		○

-  Included in the kit
- High performance
- Good performance
- Low performance
- ✗ Not suitable
- Not determined

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