







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	○	○	●	○
ATF2 phospho-T71	LB	●	X	X	○
BAD phospho-S112	●	○	LB	○	○

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







Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Beta Catenin phospho-T41/S33/37	○	○	LB	●	○
Beta Catenin total	○	○	LB	●	○
BTK phospho-Y223	LB	○	○	●	○
BTK total	LB	○	○	●	○
Cereblon total NEW	●	●	●	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	○	●	○
CREB phospho-S133	LB	○	●	X	○
CREB total	LB	○	○	●	○
EGFR phospho-pan	○	○	●	LB	○
EGFR phospho-Y1068	X	○	●	LB	○
EGFR total	○	○	●	LB	○
EIF2 α phospho-S52	●	○	○	LB	○
EIF2 α total	●	○	○	LB	○
EIF4E phospho-S209	LB	○	X	○	○
eNOS phospho-S1177 NEW	●	●	LB	●	○
eNOS total NEW	●	●	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
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	●	○
FGFR2 phospho-Y653/654	X	LB	○	○	○
FGFR2 total	X	LB	○	○	○
FoxO1 phospho-S256	LB	○	X	●	○
FoxO1 total	LB	○	X	●	○
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	○	○	○	○

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







Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
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	○
IRS1 phospho-S307 mouse / S312 human	LB	○	○	●	○
IRS1 total	LB	○	○	●	○
IRS2 total	LB	○	○	●	○
IKKB phospho-S177/181	LB	○	X	●	○
IKKB total	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	○	●	○	○







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

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
MKK4 phospho-S257	LB	●	●	●	○
MKK4 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	○
PLCg1 phospho-Y1217 NEW	X	X	LB	X	○
PLCg1 total NEW	●	●	LB	●	○
PLCg2 phospho-Y1217	●	●	●	LB	○
PLCg2 total	●	X	X	LB	○







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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Phospholamban phospho-S16	●	●	LB ●	●	○
Phospholamban total	●	●	LB ●	●	○
RB phospho-S780	LB	○	○	○	○
RB phospho-S807/811	LB	○	○	○	○
RB total	LB	○	○	○	○
S6RP phospho-S235/236	LB	○	●	●	○
SHP1 phospho-Y564	X	X	●	LB	○
SHP1 total	X	●	●	LB	○
SHP2 phospho-Y542	X	LB	●	X	○
SHP2 total	X	LB	●	X	○
SLP-76 phospho-S376	LB	○	○	●	○
SLP-76 total	LB	○	○	●	○
SMAD1 phospho-S463/465	●	○	○	LB	○
SMAD1 total	●	○	○	LB	○
SMAD2 phospho-S465/467	X	●	LB	○	○
SMAD2 total	X	●	LB	○	○
SMAD3 phospho-S463/465	●	○	●	LB	○
SMAD3 total	●	○	●	LB	○

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

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
c-Src phospho-Y419	X	○	LB	●	○
c-Src total	●	○	LB	●	○
SLP76 phospho-S376 mouse	LB	●	○	○	○
SLP76 total mouse	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	○
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	●	●	○

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

Lysis buffer compatibility

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
TAU phospho-T181					
TAU total					
TBK1 phospho-S172					
TBK1 total					
TDP-43 aggregated					
ULK1 phospho-S757 mouse / S758 human					
VASP phospho-S157					
VASP phospho-S239					
VASP total					
VEGFR2 phospho-Y1175					
YAP phospho-S127					
YAP total					
ZAP-70 phospho-Y319					
ZAP-70 total					

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

July 2020 - Copyright Cisbio. All rights reserved. All trademarks are property of Cisbio Bioassays.

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



cisbio