

# Cisbio lysis buffer compatibility

## Cell Signaling: Biomarkers, Phospho- & total-protein Assays

CELL SIGNALING

CISBIO LYSIS BUFFER COMPATIBILITY	64K11FDF	64K12FDF	64K13FDF	64K14FDF
Alpha-SMA	-	ND	LB+++	-
Alpha-Tubulin Housekeeping	++	+++	+++	LB+++
4EBP1 phospho-T37/46	LB+++	ND	ND	ND
ACC phospho-S79	+++	ND	-	LB+++
AKT1 phospho-S473	LB+++	+	ND	ND
AKT1 total	LB+++	+	ND	ND
AKT2 phospho-S473	LB+++	+	ND	ND
AKT2 total	LB+++	+	ND	ND
AKT3 phospho-S473	LB+++	+	ND	ND
AKT3 total	LB+++	+	ND	ND
AKT phospho-S473 & AKT phospho-T308	LB+++	ND	+	+++
AKT total	LB+++	ND	ND	+++
AMPK phospho-T172	LB+++	ND	ND	+
AMPK total	LB+++	ND	ND	+++
ATF2 phospho-T711	LB+++	ND	ND	ND
BAD phospho-S112	+	ND	LB+++	ND
Beta Catenin phospho-T41/S33/37	ND	ND	LB+++	++
Beta Catenin total	ND	ND	LB+++	++
BTK phospho-Y223	LB+++	ND	ND	++
BTK total	LB+++	ND	ND	++
c-Jun phospho-S63	LB+++	ND	-	++
c-RAF phospho-S338	+	LB+++	ND	+
CREB phospho-S133	LB+++	ND	++	-
CREB total	LB+++	ND	ND	++
EGFR phospho-pan	ND	ND	+++	LB+++
EGFR phospho-Y1068	-	ND	++	LB+++
EGFR total	ND	ND	+++	LB+++
EIF2 $\alpha$ phospho-S52	+++	ND	ND	LB+++
EIF2 $\alpha$ total	+++	ND	ND	LB+++
EIF4E phospho-S209	LB+++	ND	-	ND

+++ High performance    ++ Good performance    + Low performance    - Not suitable  
 ND Not determined    LB Lysis buffer included in the kit

CISBIO LYSIS BUFFER COMPATIBILITY	64K11FDF	64K12FDF	64K13FDF	64K14FDF
Advanced ERK phospho-T202/Y204	LB+++	ND	-	+++
ERK phospho-T202/Y204	LB+++	ND	-	+++
ERK total	LB+++	ND	-	+++
FGFR2 phospho-Y653/654	-	LB+++	ND	ND
FGFR2 total	-	LB+++	ND	ND
FoxO1 phospho-S256	LB+++	ND	-	++
FoxO1 total	LB+++	ND	-	++
GSK3 alpha phospho-S21	+++	ND	+	LB+++
GSK3 alpha total	+++	ND	+	LB+++
GSK3 beta phospho-S9	+++	ND	+	LB+++
GSK3 beta total	+++	ND	+	LB+++
H2AX phospho-S139	+++	ND	ND	LB+++
HER2 phospho-Y1221/1222	-	ND	+	LB+++
HER2 total	ND	ND	++	LB+++
HER3 phospho-Y1289	+	ND	ND++	LB+++
HER3 total	ND	ND	ND	LB+++
HIF-1 alpha	++	ND	++	LB+++
Histone H3 phospho-T4	LB+++	ND	ND	ND
Histone H3 phospho-S10	LB+++	ND	ND	ND
Histone H3 total	LB+++	ND	-	ND
Insulin receptor phospho-Y1150/1151	LB+++	+	ND	ND
Insulin receptor total	LB+++	+	ND	ND
IRF3 phospho-S386	-	LB+++	++	-
IRS1 phospho-S307 mouse / S312 human	LB+++	ND	ND	+
IRS1 total	LB+++	ND	ND	+
IRS2 total	LB+++	ND	ND	+++
IKKB phospho-S177/181	LB+++	ND	-	+
IKKB total <b>NEW</b>	LB+++	+	+	+++
JNK phospho-T183/Y185	-	ND	LB+++	ND
LRRK2 phospho-S9358	-	-	-	LB+++
LRRK2 total	-	-	-	LB+++
MEK1 phospho-S298	LB+++	ND	ND	ND
MEK1 phospho-S218/222	LB+++	ND	ND	ND
MEK1 total	LB+++	ND	ND	ND
MEK1/2 phospho-S218/222	LB+++	ND	++	ND
MNK phospho-T197/202	LB+++	-	ND	ND
mTOR phospho-S2448	LB+++	ND	++	+++
NDRG1 phospho-T346	ND	ND	ND	LB+++
NDRG1 phospho-S330	ND	ND	ND	LB+++
NDRG1 total	ND	ND	ND	LB+++
NFkB phospho-S536	ND	ND	+++	LB+++
NFkB total	ND	ND	+++	LB+++
P38 phospho-T180/Y182	LB+++	ND	-	-
P53 phospho-S15	LB+++	ND	-	-
P53 total	LB+++	ND	-	-

+++ High performance    ++ Good performance    + Low performance    - Not suitable  
ND Not determined    LB Lysis buffer included in the kit  
July 2018

CISBIO LYSIS BUFFER COMPATIBILITY	64K11FDF	64K12FDF	64K13FDF	64K14FDF
p70S6K phospho-T389	LB+++	ND	++	ND
PDGFRβ phospho-Y751 <b>NEW</b>	++	++	++	LB+++
Phospholamban total	LB+++	-	ND	ND
S6RP phospho-S235/236	LB+++	ND	+	+++
SLP-76 phospho-S376	LB+++	ND	ND	++
SLP-76 total	LB+++	ND	ND	++
SMAD1 phospho-S463/465	+++	ND	ND	LB+++
SMAD1 total	+++	ND	ND	LB+++
SMAD3 phospho-S463/465	+++	ND	+++	LB+++
SMAD3 total	+++	ND	+++	LB+++
c-Src phospho-Y419	-	ND	LB+++	++
c-Src total	+++	ND	LB+++	+++
STAT1 phospho-Y701	++	LB+++	+++	+++
STAT1 total	+++	LB+++	++	+++
STAT3 phospho-Y705	LB+++	ND	++	+++
STAT3 total	LB+++	ND	ND	ND
STAT5 phospho-Y694	+	ND	-	LB+++
TAU phospho-S202/T205	LB+++	ND	ND	ND
TBK1 phospho-S172 <b>NEW</b>	+++	-	-	LB+++
TBK1 total <b>NEW</b>	+++	-	-	LB+++
ULK1 phospho-S757 mouse / S758 human	LB+++	ND	+++	+++
VASP phospho-S157	+++	LB++	+++	++
VASP phospho-S239	++	LB+++	+	++
VASP total	+	LB+++	++	++
VEGFR2 phospho-Y1175	ND	LB+++	ND	ND
YAP phospho-S127	+++	ND	-	LB+++
YAP total	+++	ND	-	LB+++
ZAP-70 phospho-Y319	+	ND	++	LB+++
ZAP-70 total	+	ND	++	LB+++