







# LYSIS BUFFER COMPATIBILITY HTRF ASSAYS

JANUARY 2023

## 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	○
ADAR1 total <b>NEW</b>	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	●	○	○	○


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


	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
AKT2 total		●	○	○	○
AKT3 phospho-S473		●	○	○	○
AKT3 total		●	○	○	○
AKT phospho-S473 & AKT phospho-T308		○	●	●	○
AKT total		○	○	●	○
AMPK phospho-T172		○	○	●	○
AMPK total		○	○	●	○
Androgen receptor total	●	●		●	○
AP2 total	●	●	●		○
ATF2 phospho-T71		●	X	X	○
ATM phospho-S1981		●	○	●	○
ATM total		●	○	●	○
ATG14 phospho-S29	●	●	●		○
ATG14 total	●	X	●		○
ATG16L1 phospho-S278 <b>NEW</b>	●	●		●	○
ATG16L1 total <b>NEW</b>	●	●		●	○
Aurora A phospho-T288		X	X	●	○
Aurora A total		X	X	●	○
Aurora B, phospho-T232		X	X	●	○
Aurora B total		X	X	●	○
BAD phospho-S112	●	○		○	○
B-arrestin 1 total	●	●	●		○

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







	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
B-arrestin 2 total	●	●	●	LB	○
Beta Catenin phospho-T41/S33/37	○	○	LB	●	○
Beta Catenin total	○	○	LB	●	○
BRD2 total	●	●	●	LB	○
BRD3 total	●	●	●	LB	○
BRD4 total	●	●	●	LB	○
BRD7 total	●	X	X	LB	○
BRD9 total	●	●	●	LB	○
BTK phospho-Y223	LB	○	○	●	○
BTK total	LB	○	○	●	○
Cereblon total	●	●	●	LB	○
CDK1 phospho-T14 <b>NEW</b>	LB	●	●	●	○
CDK1 phospho-Y15 <b>NEW</b>	LB	●	●	●	○
CDK1 total	LB	●	●	●	○
CDK2 phospho-Y15	●	LB	●	●	○
CDK2 total	X	LB	●	●	○
CDK4 total	LB	X	X	●	○
CDK6 total	LB	X	X	●	○
CDK7 total	LB	●	●	●	○
CDK9 total	LB	●	X	●	○
CDK12 total	●	LB	●	●	○
cGAS total <b>NEW</b>	LB	●	○	○	○

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


	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
CHK1 phospho-S345	 ●	●	●	●	○
CHK1 total	 ●	●	●	●	○
CHK2 phospho-T68	 ●	●	●	●	○
CHK2 total	 ●	●	●	●	○
c-Jun phospho-S63	 ●	○	✗	●	○
c-RAF phospho-S338		 ●	○	●	○
cRaf phospho-S43	●	●	○	○	○
cRaf total	●	●	○	○	○
CREB phospho-S133	 ●	○	●	✗	○
CREB total	 ●	○	○	●	○
Cyclin D1 phospho-T286	 ●	✗	✗	●	○
Cyclin D1 total	 ●	●	●	●	○
EGFR DEL19 total	●	●	●	 ●	○
EGFR L858R total	●	●	●	 ●	○
EGFR phospho-pan	○	○	●	 ●	○
EGFR phospho-Y1068	✗	○	●	 ●	○
EGFR total	○	○	●	 ●	○
EIF2 $\alpha$ phospho-S52	●	○	○	 ●	○
EIF2 $\alpha$ total	●	○	○	 ●	○
EIF4E phospho-S209	 ●	○	✗	○	○
eNOS phospho-S1177	●	●	 ●	●	○
eNOS total	●	●	 ●	●	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
ERa 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	○
Fibronectin	●	●	LB	●	○
Fibronectin EDA	●	●	LB	●	○
FLT3 total	●	●	X	LB	○
FoxO1 phospho-S256	LB	○	X	●	○
FoxO1 total	LB	○	X	●	○
GAPDH Housekeeping	LB	●	●	●	○
GSK3 alpha phospho-S21	●	○	●	LB	○
GSK3 alpha total	●	○	●	LB	○









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


	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
GSK3 beta phospho-S9	●	○	●	LB	○
GSK3 beta total	●	○	●	LB	○
H2AX phospho-S139	●	○	○	LB	○
H2AX total	●	●	●	LB	○
HDAC4 phospho-S246	●	●	X	LB	○
HDAC4 total	●	●	●	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	○
IRF5	●	LB	○	○	○
IRβ phospho Y1150/1151	LB	●	●	●	○
IRβ total	LB	●	●	●	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
IRS1 phospho-S307 mouse / S312 human		○	○	●	○
IRS1 total		○	○	●	○
IRS2 total		○	○	●	○
IKKB phospho-S177/181		○	X	●	○
IKKB total		●	●	●	○
JAK1 Phospho Y1034/35	X	●	●		○
JAK1 total	●	●	●		○
JAK2 Phospho Y1007/08		●	●	●	○
JAK2 total		●	●	●	○
JAK3 total	●	●	●		○
JNK phospho-T183/Y185	X	○		○	○
Ki-67	●	●	X		○
LRRK2 phospho-S9358	X	X	X		○
LRRK2 total	X	X	X		○
MDA5 total	●	●		●	○
MDA5 aggregation	X	X		X	○
MEK1 phospho-S298	●		●	●	●
MEK1 phospho-S218/222		○	○	○	○
MEK1 total		○	○	○	○
MEK1/2 phospho-S218/222		○	●	○	○
MET phospho-Y1234/1235 <b>NEW</b>	●		○	○	○
MET total <b>NEW</b>	●		○	○	○







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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
MKK4 phospho-S257		●	●	●	○
MKK4 total		●	●	●	○
MLKL phospho-S358	●	●	●		○
MLKL total	●	●	●		○
MNK phospho-T197/202		X	○	○	○
mTOR phospho-S2448		○	●	●	○
NDRG1 phospho-T346	○	○	○		○
NDRG1 phospho-S330	○	○	○		○
NDRG1 total	○	○	○		○
NFκB phospho-S536	○	○	●		○
NFκB total	○	○	●		○
P21	●		○	○	○
P38 phospho-T180/Y182		○	X	X	○
P53 phospho-S15		○	X	X	○
P53 total		○	X	X	○
P62/SQSTM1-S403	●	●	●		○
P62/SQSTM1 total	●	●	●		○
p70S6K phospho-T389		○	●	○	○
PDGFRβ phospho-Y751	●	●	●		○
PLCg1 phospho-Y1217	X	X		X	○
PLCg1 total	●	●		●	○
PLCg2 phospho-Y1217	●	●	●		○







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



	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
PLCg2 total	●	✗	✗	LB	○
Phospholamban phospho-S16	●	●	LB	●	○
Phospholamban total	●	●	LB	●	○
PR total	●	●	●	LB	○
RAB10 phospho-T73	✗	●	LB	✗	○
RAB10 total	●	●	LB	●	○
RBM39 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	✗	LB	○	✗	●
RSK1 phospho-T359/S363	●	LB	○	○	○
RSK1 total	✗	LB	○	○	○
S6RP phospho-S235/236	LB	○	●	●	○
SARS-CoV-2 capsid	●	●	●	●	LB
SHP1 phospho-Y564	✗	✗	●	LB	○
SHP1 total	✗	●	●	LB	○
SHP2 phospho-Y542	✗	LB	●	✗	○

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







	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
SHP2 total	X	LB	●	X	○
SLP76 phospho-S376	LB	○	○	●	○
SLP76 phospho-Y145	X	●	●	LB	○
SLP76 total	LB	●	●	●	○
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	○
SMAD4 phospho-T277	●	●	●	LB	●
SMAD4 total	●	●	●	LB	●
SMARCA2 total	LB	●	●	●	○
SMARCA4 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	○	●	●	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
STAT3 total		○	○	○	○
STAT5 phospho-Y694	●	○	X		○
STAT6 phospho-Y641	X	X	●		○
STAT6 total	●	●	●		○
STING phospho-S366	●	●	●		○
STING total	●	●	●		○
STING (m) phospho-S365	X	X	●		○
STING (m) Total	●	●	●		○
SYK phospho-Y525	●	●		●	○
SYK total	●	●		●	○
TAU phospho-S202/T205		●	●	●	○
TAU phospho-S356		●	●	●	○
TAU phospho-S404		X	●	●	○
TAU phospho-S422		X	●	●	○
TAU phospho-T181		●	●	●	○
TAU total		●	●	●	○
TBK1 phospho-S172	●	X	X		○
TBK1 total	●	X	X		○
TDP43 aggregated		●	●	●	○
Tie2 phospho-Y992		X	X	●	●
TOM20	●	●	X		○
TYK2 phospho-Y1054/55	X	X	●		○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
TYK2 total	●	●	●	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	○
ZAP-70 phospho-Y319	●	○	●	LB	○
ZAP-70 total	●	○	●	LB	○

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