



# Lysis buffer compatibility HTRF<sup>®</sup> assays

Cell signaling: Phospho- and total-protein assays: April 2023

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Alpha-SMA	X	○	LB	X	○
Alpha-Synuclein phospho-S129	●	●	●	●	LB
Alpha-Synuclein phospho-S129 Aggregation <b>NEW</b>	X	X	X	X	LB
Alpha-Synuclein total	●	●	●	●	LB
Alpha-Tubulin Housekeeping	●	●	●	LB	○
ADAR1 total	LB	●	●	○	○
4EBP1 phospho-T37/46	LB	○	○	○	○
ACC phospho-S79	●	○	X	LB	○
ACC total	X	X	●	LB	○
AKT1 phospho-S473	LB	●	○	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
AKT1 total					
AKT2 phospho-S473					
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					
ATM phospho-S1981					
ATM total					
ATG14 phospho-S29					
ATG14 total					
ATG16L1 phospho-S278					
ATG16L1 total					
Aurora A phospho-T288					
Aurora A total					
Aurora B, phospho-T232					

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







	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Aurora B total		X	X	●	○
BAD phospho-S112	●	○		○	○
B-arrestin 1 total	●	●	●		○
B-arrestin 2 total	●	●	●		○
Beta Catenin phospho-T41/S33/37	○	○		●	○
Beta Catenin total	○	○		●	○
BRD2 total	●	●	●		○
BRD3 total	●	●	●		○
BRD4 total	●	●	●		○
BRD7 total	●	X	X		○
BRD9 total	●	●	●		○
BTK phospho-Y223		○	○	●	○
BTK total		○	○	●	○
Cereblon total	●	●	●		○
CDK1 phospho-T14		●	●	●	○
CDK1 phospho-Y15		●	●	●	○
CDK1 total		●	●	●	○
CDK2 phospho-Y15	●		●	●	○
CDK2 total	X		●	●	○
CDK4 total		X	X	●	○
CDK5 total <b>NEW</b>		●	●	●	○
CDK6 total		X	X	●	○

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







	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
CDK7 total		●	●	●	○
CDK8 total <b>NEW</b>		●	●	●	○
CDK9 total		●	X	●	○
CDK12 total	●		●	●	○
cGAS total		●	○	○	○
CHK1 phospho-S345		●	●	●	○
CHK1 total		●	●	●	○
CHK2 phospho-T68		●	●	●	○
CHK2 total		●	●	●	○
c-Jun phospho-S63		○	X	●	○
c-RAF phospho-S338			○	●	○
cRaf phospho-S43	●	●	○	○	○
cRaf total	●	●	○	○	○
CREB phospho-S133		○	●	X	○
CREB total		○	○	●	○
Cyclin D1 phospho-T286		X	X	●	○
Cyclin D1 total		●	●	●	○
Cyclin E1 total	●		●	●	○
EGFR DEL19 total	●	●	●		○
EGFR L858R total	●	●	●		○
EGFR phospho-pan	○	○	●		○
EGFR phospho-Y1068	X	○	●		○

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

	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	●	○
ERα 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	●	○

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







	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Fibronectin EDA	●	●	LB	●	○
FLT3 phospho Y842 <i>New</i>	X	●	●	LB	○
FLT3 phospho Y589/591 <i>New</i>	X	●	●	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	○
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	○	○	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
Histone H3 phospho-S10		○	○	○	○
Histone H3 total		○	X	○	○
Insulin receptor phospho-Y1150/1151		●	○	○	○
Insulin receptor total		●	○	○	○
IRAK3 total		●	●	●	●
IRAK4 total		●	●	●	●
IRF3 phospho-S386	X		●	X	○
IRF5	●		○	○	○
IRβ phospho Y1150/1151		●	●	●	○
IRβ total		●	●	●	○
IRS1 phospho-S307 mouse / S312 human		○	○	●	○
IRS1 total		○	○	●	○
IRS2 total		○	○	●	○
IKKβ phospho-S177/181		○	X	●	○
IKKα phospho-S176/180		X	●	●	○
IKKα total		X	X	●	○
IKKβ total		●	●	●	○
JAK1 Phospho Y1034/35	X	●	●		○
JAK1 total	●	●	●		○
JAK2 Phospho Y1007/08		●	●	●	○
JAK2 total		●	●	●	○







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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
JAK3 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	○	●	○	○
MET phospho-Y1234/1235	●	LB	○	○	○
MET total	●	LB	○	○	○
MKK4 phospho-S257	LB	●	●	●	○
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	○





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



	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
NDRG1 total	○	○	○	LB	○
NFκB phospho-S536	○	○	●	LB	○
NFκB total	○	○	●	LB	○
NRF2 total	LB	●	●	●	○
P21	●	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	○
PINK1 total	LB	X	X	●	○
PLCg1 phospho-Y1217	X	X	LB	X	○
PLCg1 total	●	●	LB	●	○
PLCg2 phospho-Y1217	●	●	●	LB	○
PLCg2 total	●	X	X	LB	○
Phospholamban phospho-S16	●	●	LB	●	○
Phospholamban total	●	●	LB	●	○
PR total	●	●	●	LB	○
RAB10 phospho-T73	X	●	LB	X	○

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







	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
RAB10 total	●	●	LB	●	○
RBM39 total	●	X	X	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	●	●	●	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
SLP76 phospho-S376 mouse					
SLP76 total mouse					
SMAD1 phospho-S463/465					
SMAD1 total					
SMAD2 phospho-S465/467					
SMAD2 total					
SMAD3 phospho-S463/465					
SMAD3 total					
SMAD4 phospho-T277					
SMAD4 total					
SMARCA2 total					
SMARCA4 total					
Spike S1 protein					
c-Src phospho-Y419					
c-Src total					
STAT1 phospho-Y701					
STAT1 total					
STAT3 phospho-Y705					
STAT3 total					
STAT5 phospho-Y694					
STAT6 phospho-Y641					

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
STAT6 total	●	●	●	LB	○
STING phospho-S366	●	●	●	LB	○
STING total	●	●	●	LB	○
STING (m) phospho-S365	X	X	●	LB	○
STING (m) 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	●	●	○
TAU phospho-T181	LB	●	●	●	○
TAU total	LB	●	●	●	○
TBK1 phospho-S172	●	X	X	LB	○
TBK1 total	●	X	X	LB	○
TDP43 aggregated	LB	●	●	●	○
TDP43 phospho-S409/410	LB	●	●	●	○
TDP43 total	LB	●	●	●	○
TFEB phospho-S211	X	●	●	LB	○
TFEB total	●	●	●	LB	○
Tie2 phospho-Y992	LB	X	X	●	●

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF
TOM20	●	●	X	LB	○
TYK2 phospho-Y1054/55	X	X	●	LB	○
TYK2 total	●	●	●	LB	○
Ubiquitin phospho-S65	LB	X	●	●	○
ULK1 phospho-S556 <b>NEW</b>	●	●	●	LB	○
ULK1 phospho-S757 mouse / S758 human	LB	○	●	●	○
ULK1 total <b>NEW</b>	●	●	●	LB	○
VASP phospho-S157	●	LB	●	●	○
VASP phospho-S239	●	LB	●	●	○
VASP total	●	LB	●	●	○
VEGFR2 phospho-Y1175	●	LB	●	●	○
VEGFR2 total	●	LB	X	●	○
WEE1 total <b>NEW</b>	LB	X	X	X	X
XBP1s	LB	●	●	●	○
XBP1 total	LB	●	●	●	○
YAP phospho-S127	●	○	X	LB	○
ZAP-70 phospho-Y319	●	○	●	LB	○
ZAP-70 total	●	○	●	LB	○

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

