



# Lysis buffer compatibility HTRF assays.

October 2024

## Cell signaling: phospho- and total-protein assays

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
Alpha-SMA	X	○	LB	X	○	○
Alpha-Synuclein phospho-S129	●	●	●	●	LB	○
Alpha-Synuclein phospho-S129 Aggregation	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	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
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	○	○	●	○	○
ALK phospho-Y1604	X	●	X	LB	○	○
ALK total	●	●	●	LB	○	○
AMPK phospho-T172	LB	○	○	●	○	○
AMPK total	LB	○	○	●	○	○
Androgen receptor full-length	●	●	LB	●	○	○
Androgen receptor total	●	●	LB	●	○	○
Androgen receptor variant 7	●	●	LB	●	○	○
AP2 total	●	●	●	LB	○	○
ATF2 phospho-T71	LB	●	X	X	○	○
ATF4 total	●	●	●	LB	○	○
ATG14 phospho-S29	●	●	●	LB	○	○
ATG14 total	●	X	●	LB	○	○
ATG16L1 phospho-S278	●	●	LB	●	○	○
ATG16L1 total	●	●	LB	●	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
ATM phospho-S1981	LB	●	○	●	○	○
ATM total	LB	●	○	●	○	○
Aurora A phospho-T288	LB	X	X	●	○	○
Aurora A total	LB	X	X	●	○	○
Aurora B, phospho-T232	LB	X	X	●	○	○
Aurora B total	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	●	○	○
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	LB	●	●	●	○	○
CDK1 phospho-Y15	LB	●	●	●	○	○
CDK1 total	LB	●	●	●	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
CDK2 phospho-Y15	●	LB	●	●	○	○
CDK2 total	X	LB	●	●	○	○
CDK4 total	LB	X	X	●	○	○
CDK5 total	LB	●	●	●	○	○
CDK6 total	LB	X	X	●	○	○
CDK7 total	LB	●	●	●	○	○
CDK8 total	LB	●	●	●	○	○
CDK9 total	LB	●	X	●	○	○
CDK12 total	●	LB	●	●	○	○
cGAS total	LB	●	○	○	○	○
CHK1 phospho-S345	LB	●	●	●	○	○
CHK1 total	LB	●	●	●	○	○
CHK2 phospho-T68	LB	●	●	●	○	○
CHK2 total	LB	●	●	●	○	○
c-Jun phospho-S63	LB	○	X	●	○	○
CK1α total <b>NEW</b>	●	●	X	LB	○	○
Cofilin phospho-S3	X	X	●	LB	●	○
Cofilin total	●	●	●	LB	●	○
c-RAF phospho-S338	X	LB	●	X	○	○
c-RAF phospho-S43	●	LB	●	●	●	○
c-RAF Total	●	LB	●	●	○	○
CREB phospho-S133	LB	○	●	X	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
CREB total	LB	○	○	●	○	○
CREBBP total	LB	●	●	●	○	○
Cyclin B1 phospho-S126	LB	X	X	●	○	○
Cyclin B1 total	LB	●	X	●	○	○
Cyclin D1 phospho-T286	LB	X	X	●	○	○
Cyclin D1 total	LB	●	●	●	○	○
Cyclin E1 total	●	LB	●	●	○	○
DDR1 phospho-panY	●	●	●	LB	○	○
DDR1 phospho-Y513	X	X	●	LB	○	○
DDR1 phospho-Y796	●	●	●	LB	○	○
DDR1 total	●	●	X	LB	○	○
EGFR DEL19 total	●	●	●	LB	○	○
EGFR L858R total	●	●	●	LB	○	○
EGFR phospho-pan	○	○	●	LB	○	○
EGFR phospho-Y1068	X	○	●	LB	○	○
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	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
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 phospho Y842	X	●	●	LB	○	○
FLT3 phospho Y589/591	X	●	●	LB	○	○
FLT3 total	●	●	X	LB	○	○
FoxO1 phospho-S256	LB	○	X	●	○	○
FoxO1 total	LB	○	X	●	○	○
GAPDH Housekeeping	LB	●	●	●	○	○
GSK3 alpha phospho-S21	●	○	●	LB	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
GSK3 alpha total	●	○	●	LB	○	○
GSK3 beta phospho-S9	●	○	●	LB	○	○
GSK3 beta total	●	○	●	LB	○	○
GSPT1 total	●	●	●	LB	○	○
H2AX phospho-S139	●	○	○	LB	○	○
H2AX total	●	●	●	LB	○	○
HDAC4 phospho-S246	●	●	X	LB	○	○
HDAC4 total	●	●	●	LB	○	○
HDAC5 phospho-S259	X	X	●	LB	○	○
HDAC5 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	○	○	○
HPK1 total	●	●	●	LB	○	○
HSP27 phospho-S78	LB	●	●	●	○	○
HSP27 total	LB	●	●	●	○	○
IDO1 total	LB	●	●	●	○	○







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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
IKZF1 total	●	●	●	LB	○	○
IKZF3 total	●	●	●	LB	○	○
Insulin receptor phospho-Y1150/1151	LB	●	○	○	○	○
Insulin receptor total	LB	●	○	○	○	○
IRAK3 total	LB	●	●	●	●	○
IRAK4 total	LB	●	●	●	●	○
IRF3 phospho-S386	X	LB	●	X	○	○
IRF5 total <small>UPDATE</small>	●	LB	●	●	○	○
IRβ phospho Y1150/1151	LB	●	●	●	○	○
IRβ total	LB	●	●	●	○	○
IRS1 phospho-S307 mouse / S312 human	LB	○	○	●	○	○
IRS1 total	LB	○	○	●	○	○
IRS2 total	LB	○	○	●	○	○
IKKβ phospho-S177/181	LB	○	X	●	○	○
IKKα phospho-S176/180	LB	X	●	●	○	○
IKKα total	LB	X	X	●	○	○
IKKβ total	LB	●	●	●	○	○
JAK1 Phospho Y1034/35	X	●	●	LB	○	○
JAK1 total	●	●	●	LB	○	○
JAK2 Phospho Y1007/08	LB	●	●	●	○	○
JAK2 total	LB	●	●	●	○	○
JAK3 total	●	●	●	LB	○	○

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



	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
JNK phospho-T183/Y185	X	○	LB	○	○	○
Ki-67	●	●	X	LB	○	○
KRAS total	X	X	X	X	X	LB
LC3B-II total	X	X	LB	X	X	○
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	○	●	○	○	○
MerTK phospho-Y749	●	●	●	LB	○	○
MerTK total	●	X	X	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	○	○	○	○
MSH3 total	●	LB	○	○	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
mTOR phospho-S2448	LB	○	●	●	○	○
Myd88 total	LB	●	●	●	○	○
NDRG1 phospho-T346	○	○	○	LB	○	○
NDRG1 phospho-S330	○	○	○	LB	○	○
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	○	○
P300 total	LB	●	●	●	○	○
p70S6K phospho-T389	LB	○	●	○	○	○
PDGFRb 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	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
PLK1 phospho-T210	LB	●	●	●	○	○
PLK1 total	LB	●	●	●	○	○
Phospholamban phospho-S16	●	●	LB	●	○	○
Phospholamban total	●	●	LB	●	○	○
PR total	●	●	●	LB	○	○
RAB10 phospho-T73	X	●	LB	X	○	○
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	●	●	●	○	○
RNA pol II phospho-S2 <b>NEW</b>	LB	●	●	●	○	○
RNA pol II phospho-S5 <b>NEW</b>	LB	●	●	●	○	○
RNA pol II total <b>NEW</b>	LB	●	●	●	○	○
RSK1 phospho-S380	X	LB	○	X	●	○
RSK1 phospho-T359/S363	●	LB	○	○	○	○
RSK1 total	X	LB	○	○	○	○
S6RP phospho-S235/236	LB	○	●	●	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
SALL4 total <b>NEW</b>	●	●	X	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	●	●	●	○	○
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	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
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	○	○
STAT5 total	●	●	●	LB	○	○
STAT6 phospho-Y641	X	X	●	LB	○	○
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	●	●	●	○	○

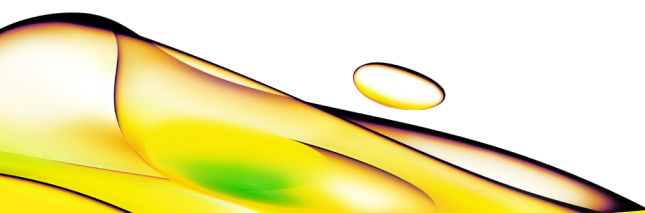
- LB Included in the kit
- High performance
- Good performance
- Low performance
- X Not suitable
- Not determined

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
TBK1 phospho-S172	●	✗	✗	LB ●	○	○
TBK1 total	●	✗	✗	LB ●	○	○
TDP43 aggregated	LB ●	●	●	●	○	○
TDP43 phospho-S409/410	LB ●	●	●	●	○	○
TDP43 total	LB ●	●	●	●	○	○
TFEB phospho-S211	✗	●	●	LB ●	○	○
TFEB total	●	●	●	LB ●	○	○
Tie2 phospho-Y992	LB ●	✗	✗	●	●	○
TOM20	●	●	✗	LB ●	○	○
TYK2 phospho-Y1054/55	✗	✗	●	LB ●	○	○
TYK2 total	●	●	●	LB ●	○	○
Ubiquitin phospho-S65	LB ●	✗	●	●	○	○
ULK1 phospho-S556	●	●	●	LB ●	○	○
ULK1 phospho-S757 mouse / S758 human	LB ●	○	●	●	○	○
ULK1 total	●	●	●	LB ●	○	○
VASP phospho-S157	●	LB ●	●	●	○	○
VASP phospho-S239	●	LB ●	●	●	○	○
VASP total	●	LB ●	●	●	○	○
VAV1 phospho-Y174	LB ●	✗	○	○	○	○
VAV1 total	LB ●	●	○	○	○	○
VEGFR2 phospho-Y1175	●	LB ●	●	●	○	○
VEGFR2 total	●	LB ●	✗	●	○	○

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

	64KL1FDF	64KL2FDF	64KL3FDF	64KL4FDF	64KL5FDF	64KL6FDF
WEE1 total	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



revvity