

Lysates Provided With Bio-Plex™ Assays

Refer to the table below to identify the lysates to be used with each Bio-Plex phosphoprotein and total target assay. Then follow the instructions in the Bio-Plex phosphoprotein detection instruction manual.

NOTE: These lysates are used to determine the assay performance and should not be considered as references. Activation signals and ratios should be based on experimental control lysates. For determining total protein concentrations, consider Bio-Rad's *DC*™ protein assay kit (Bio-Rad catalog #500-0112).

Lysate	Phospho-Specific Targets
EGF-Treated HEK293	Akt, ERK1, ERK2, ERK1/2, GSK-3 α/β , MEK1
EGF-Treated HeLa	EGFR, HSP27, S6 Ribosomal Protein
IFN- α -Treated HeLa	Stat2, Stat3
IL-4-Treated HeLa	Stat6
M Phase-Treated HeLa	Histone H3
NGF β -Treated PC12	IRS-1, p70 S6 kinase, p90RSK, TrkA
PDGF-Treated NIH3T3	PDGF Receptor β
TNF- α -Treated HeLa	I κ B- α , NF- κ B p65
UV-Treated HEK293	ATF-2, c-Jun, CREB, JNK, p38 MAPK, p53
Untreated HeLa	Background control for all phosphoprotein assays. Positive control for all total target assays.