

Acute Phase Response
Cancer
Cardiovascular Disease
Diabetes
Cytokines, Chemokines,
Growth Factors
Immunology/Inflammation
Immunoglobulin Isotyping
Cell Signaling
Toxicology

Bio-Plex Pro™ Cell Signaling Assays

Akt, ATF-2, β -Actin, BAD, Btk, c-Abl, c-Jun, CREB, EGFR, Erk1/2, GSK-3 α/β , HER-2, HSP27, GAPDH, IGF-1R, I κ B- α , IR- β , IRS-1, JNK, Lyn, MEK1, mTOR, NF- κ B p65, p38 MAPK, p53, P70 S6 Kinase, p90 RSK, PDGFR- α , PDGFR- β , PI3K p85, PTEN, S6 Ribosomal Protein, Smad2, Src, Stat1, Stat3, Syk, VEGFR-2, ZAP-70

MAGNETIC SEPARATION ENABLED

- Magnetic or vacuum assay separation
- Convenient kit format
- Flexible ordering options
- Minimal sample volume
- Antibodies from CST



Cell Signaling Assays with Exceptional Sensitivity

The Bio-Plex Pro cell signaling assays are magnetic bead-based immunoassays designed to meet the sensitivity needs of the most discerning scientists. The multiplex format enables robust, reproducible, and simultaneous measurement of proteins involved in key intracellular signaling pathways. Design your own assay choosing from a broad selection of phosphoprotein and total protein targets to investigate pathways associated with:

- Cancer
- Inflammation
- Diabetes
- Neurological disorders
- Cardiovascular disorders
- Drug mechanism of action
- Toxicology

These assays incorporate several features to enhance both quality and ease of use:

- Panel includes house keeping proteins GAPDH and β -Actin
- Assay quick guide to get you started right away
- Assay protocol optimized for exceptional sensitivity and broad dynamic range
- Flexible ordering options — order singleplex sets or visit www.bio-rad.com/assaybuilder to configure a custom premixed kit

Benefits of Magnetic Bead-Based Assays

Magnetic bead-based assays enable automation of wash steps with a Bio-Plex Pro series or similar wash station. This innovation simplifies assay processing and eliminates the need for a vacuum manifold. After adopting the magnetic assay workflow, many users experience improved assay precision, in particular with viscous samples.

Rigorous Assay Validation

All Bio-Plex Pro cell signaling assays undergo a rigorous evaluation that includes the following parameters:

- Specificity and cross-reactivity
- Inter- and intra-assay precision
- Sensitivity (limit of detection, LOD)
- Dynamic range
- Performance characteristics in real samples

Assay Performance Definitions

Precision — the coefficient of variation (%CV) at concentrations within the assay working range

Sensitivity (LOD) — the amount of cell lysate protein for which the fluorescence intensity signal of the specific analyte is two standard deviations above the background signal

BIO-RAD

Available Assays

Table 1 shows a list of available Bio-Plex Pro cell signaling assays now offered for the semiquantitative detection of total target proteins and phosphoproteins in cell and tissue lysates.

Table 1. Available assays.

Total Target Assays	Bead Regions	Total Target Assays	Bead Regions	Phosphoprotein Assays	Bead Regions	Phosphoprotein Assays	Bead Regions
Akt*	75	JNK	34	Akt (Ser ⁴⁷³)	75	mTOR (Ser ²⁴⁴⁸)*	46
Btk *	39	MEK1	27	Akt (Thr ³⁰⁸)	75	NF-κB p65 (Ser ⁵³⁶)	37
c-Jun	56	mTOR*	46	ATF-2 (Thr ⁷¹)*	20	p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	36
CREB*	19	p38 MAPK	36	BAD (Ser ¹³⁶)*	26	p53 (Ser ¹⁵)*	53
Erk1/2	38	P70 S6 Kinase	55	Btk (Tyr ²²³)*	39	p70 S6 Kinase (Thr ³⁸⁹)	55
GSK-3β	18	PTEN*	22	c-Abl (Tyr ²⁴⁵)*	45	p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)	55
HER-2	30	Smad2	14	c-Jun (Ser ⁶³)	56	p90 RSK (Ser ³⁸⁰)*	35
IGF-1R	43	Src*	42	CREB (Ser ¹³³)*	19	PDGFR-α (Tyr ⁷⁵⁴)	28
IκB-α	67	ZAP-70*	64	EGFR (Tyr ¹⁰⁶⁸)	44	PDGFR-β (Tyr ⁷⁵¹)	57
				EGFR (Tyr ¹¹⁷³)	44	PI3K p85 (Tyr ⁴⁵⁸)*	54
Housekeeping Proteins	Bead Regions	Housekeeping Proteins	Bead Regions	Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	38	PTEN (Ser ³⁸⁰)*	22
Human GAPDH	21	β-Actin	47	GSK-3α/β (Ser ²¹ /Ser ⁹)	18	S6 Ribosomal Protein* (Ser ²³⁵ /Ser ²³⁶)	74
				HER-2 (Tyr ¹²⁴⁸)	30	Smad2 (Ser ⁴⁶⁵ /Ser ⁴⁶⁷)	14
				HSP27 (Ser ⁷⁸)*	51	Src (Tyr ⁴¹⁶)*	42
				IGF-1R (Tyr ¹¹³¹)	43	Stat1 (Tyr ⁷⁰¹)	61
				IR-β (Tyr ¹¹⁴⁶)*	43	Stat3 (Ser ⁷²⁷)	52
				IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)*	76	Stat3 (Tyr ⁷⁰⁵)	52
				IκB-α (Ser ³² /Ser ³⁶)	67	Syk (Tyr ³⁵²)*	65
				JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	34	VEGFR-2 (Tyr ¹¹⁷⁵)	29
				Lyn (Tyr ⁵⁰⁷)*	33	ZAP-70 (Tyr ³¹⁹)*	64
				MEK1 (Ser ²¹⁷ /Ser ²²¹)	27		

* Latest additions to the Bio-Plex[®] Pro cell signaling panel.

Assay Performance Characteristics

The Bio-Plex Pro cell signaling assays were evaluated for sensitivity (P/B ratio and LOD), intra- and inter-assay %CV, and species specificity (Table 2).

Table 2. Assay performance characteristics.

Total Target Assays	Intra-assay %CV	Inter-assay %CV	Sensitivity (LOD), µg/well	Reaction Species*
β-Actin	3.4	13.0	0.15	H, M, R, NHP
Akt	3.8	8.4	0.15	H, M, R
Btk	7.6	10.0	0.15	H, M
c-Jun	4.0	4.7	0.15	H
CREB	0.7	10.4	0.15	H, M, R, NHP
Erk1/2	3.4	9.2	0.15	H, M
GAPDH	1.9	4.8	0.15	H
GSK-3β	2.9	4.1	0.8	H, M, R
HER-2	5.6	4.2	0.04	H, M, R
IGF-1R	4.7	4.4	0.15	H
IκB-α	1.9	3.1	0.15	H
JNK	1.9	4.5	0.3	H
MEK 1	1.8	3.9	0.15	H, M, R
mTOR	4.1	3.3	0.15	H, M, R, NHP
p38 MAPK	4.3	9.6	0.3	H, M, R
p70 S6 Kinase	2.0	3.6	0.15	H, M, R
PTEN	5.0	3.2	0.15	H, M, R
Smad2	3.9	16.9	0.15	H
Src	11.3	5.4	0.15	H, M, R
ZAP-70	3.1	11.2	0.15	H

* H, human; M, mouse; R, rat; NHP, nonhuman primate.

Table 2. Assay performance characteristics (cont).

Phosphoprotein Assays	P/B Ratio*	Intra-assay %CV	Inter-assay %CV	Sensitivity (LOD), µg/well	Reaction Species**
Akt (Ser ⁴⁷³)	65	2.9	4.5	0.15	H, M, R
Akt (Thr ³⁰⁸)	34	3.3	6.1	0.15	H, M, R
ATF-2 (Thr ⁷¹)	68	4.0	6.3	0.15	H, M, R, NHP
BAD (Ser ¹³⁶)	12	9.7	9.4	0.15	H, M, R, NHP
Btk (Tyr ²²³)	99	3.2	1.3	0.15	H
c-Abl (Tyr ²⁴⁵)	210	1.3	10.0	0.15	H, M
c-Jun (Ser ⁶³)	295	5.8	8.4	0.15	H, M, R
CREB (Ser ¹³³)	50	4.0	8.9	0.15	H, M, R
EGFR (Tyr ¹⁰⁶⁸)	98	14.1	17.7	0.15	H
EGFR (Tyr ¹¹⁷³)	30	3.9	6.3	0.15	H
Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	48	4.3	4.9	0.15	H, M
GSK-3α/β (Ser ²¹ /Ser ⁹)	19	2.9	5.4	0.15	H, M, R
HER-2 (Tyr ¹²⁴⁹)	49	10.1	9.1	0.15	H, M, R
HSP27 (Ser ⁷⁸)	58	2.7	4.4	0.15	H, NHP
IGF-1R (Tyr ¹¹³¹)	68	4.8	6.5	0.15	H
IR-β (Tyr ¹¹⁴⁶)	74	1.5	3.2	0.15	H, M, R
IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	52	5.5	9.3	0.3	H, M, R
IκB-α (Ser ³² /Ser ³⁶)	103	6.8	8.9	0.15	H
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	73	4.7	4.1	0.15	H
Lyn (Tyr ⁵⁰⁷)	125	4.2	4.4	0.15	H, M, R
MEK1 (Ser ²¹⁷ /Ser ²²¹)	44	1.9	5.5	0.15	H, M, R
mTOR (Ser ²⁴⁴⁸)	40	3.3	9.2	0.3	H, M, R, NHP
NF-κB p65 (Ser ⁵³⁶)	20	10.7	13.5	0.3	H, M, R
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	62	3.1	6.4	0.15	H, M, R
p53 (Ser ¹⁵)	59	8.0	12.2	0.15	H
p70 S6 Kinase (Thr ³⁸⁹)	196	8.4	10.3	0.15	H, M, R
p70 S6 Kinase (Thr ⁴²¹ / Ser ⁴²⁴)	11	2.9	20.1	0.15	H, M, R
p90 RSK (Ser ³⁸⁰)	80	1.7	4.2	0.15	H, M, R, NHP
PDGFR-α (Tyr ⁷⁵⁴)	40	5.0	6.2	0.04	H, M
PDGFR-β (Tyr ⁷⁵¹)	142	8.1	10.3	0.04	H, M, R
PI3K p85 (Tyr ⁴⁵⁸)	42	9.3	11.4	1.6	H, M, R
PTEN (Ser ³⁸⁰)	39	9.0	6.3	0.3	H, M, R
S6 Ribosomal Protein (Ser ²³⁵ /Ser ²³⁶)	26	4.3	5.2	0.15	H, M, R
Smad2 (Ser ⁴⁶⁵ /Ser ⁴⁶⁷)	14	4.4	8.8	0.3	H, M, R
Src (Tyr ⁴¹⁶)	178	5.7	9.1	0.15	H, M, R
Stat1 (Tyr ⁷⁰¹)	18	3.6	7.8	0.15	H
Stat3 (Ser ⁷²⁷)	89	7.0	10.1	0.15	H, M, R
Stat3 (Tyr ⁷⁰⁵)	18	2.5	6.8	0.15	H, M, R
Syk (Tyr ³⁵²)	232	1.4	2.0	0.15	H, M, R
VEGFR-2 (Tyr ¹¹⁷⁵)	614	8.3	11.7	0.15	H, M
ZAP-70 (Tyr ³¹⁹)	138	2.6	5.1	0.15	H, M

* Signal ratio (positive/background) of stimulated cell lysate/background cell lysate.

** H, human; M, mouse; R, rat; NHP, nonhuman primate.

Correlation and Sensitivity Comparison

Western blotting is a traditional technique for detecting phosphorylated proteins within a variety of cell culture and tissue sample lysates. Comparison of the Bio-Plex Pro cell signaling assays to that of western blotting shows a correlation in the expression pattern of phosphorylated proteins (Figure 1) as well as a greater sensitivity in the detection of these phosphorylated proteins (Figure 2).

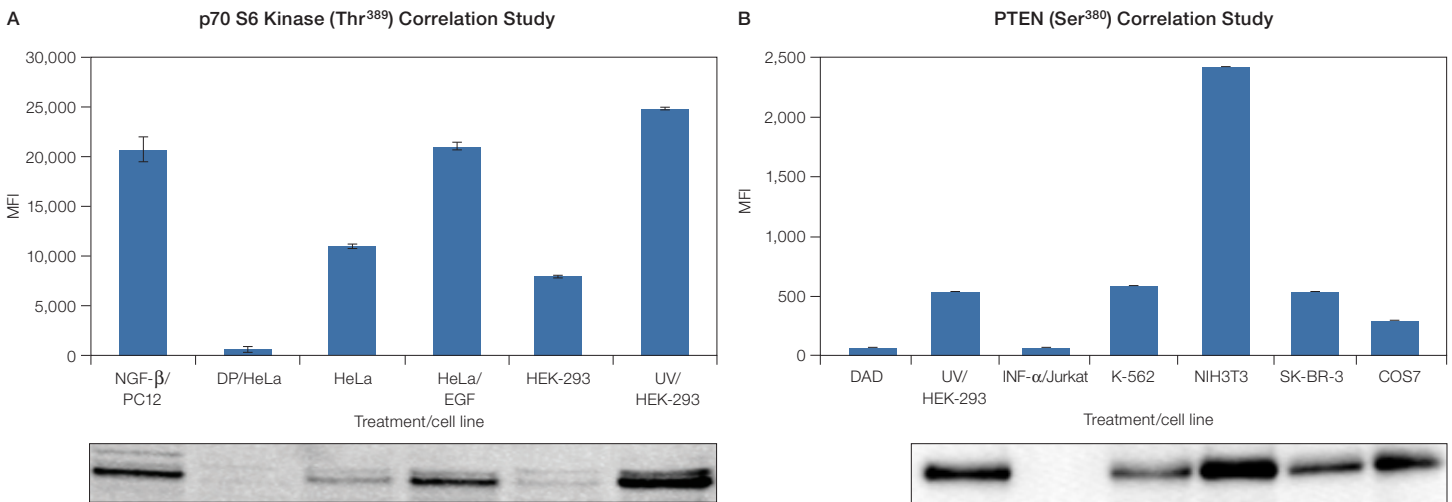


Fig. 1. Performance of Bio-Plex Pro cell signaling assays compared to western blotting. Measurements of p70 S6 Kinase (Thr³⁸⁹) (A) and PTEN (Ser³⁸⁰) (B) in a variety of cell lines. DAD indicates detection antibody diluent alone. In panel A, DP/HeLa indicates phosphatase-treated HeLa cells. This condition serves as a negative control.

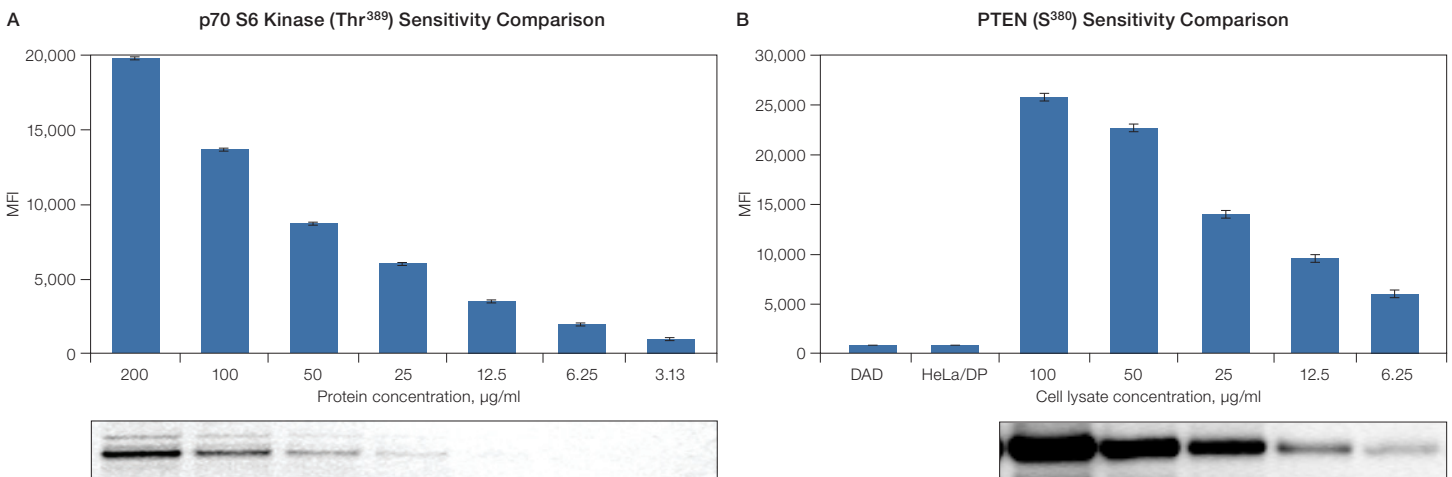


Fig. 2. Sensitivity of Bio-Plex Pro cell signaling assays. Phosphorylated p70 S6 Kinase (Thr³⁸⁹) levels were measured in PC12 cells stimulated with NGF- β (A). The Bio-Plex assay was able to detect the analyte in as little as 3 μ g/ml (150 ng/well) of total lysate protein. Phosphorylated PTEN (Ser³⁸⁰) levels were measured in HeLa cells stimulated with TNF- α (B). The Bio-Plex assay was able to detect the analyte in as little as 6 μ g/ml (300 ng/well) of total lysate protein.

Flexible Ordering Options

Premixed Panels

One kit with everything you need to run an experiment.

Individual Components

Select compatible singleplex sets and lysates. Use Table 3 to check for assay cross-reactivity. Refer to Table 4 to select the appropriate lysate controls for your assays. Complete your order with the Bio-Plex Pro cell signaling reagent kit for a mix-it-yourself multiplex solution.

x-Plex™ Custom Assay Service

A great choice for premium custom-mixed assays. Simply select your analytes of interest using the online Bio-Plex assay builder at www.bio-rad.com/assaybuilder. Assays are mixed for you at Bio-Rad and delivered as an all-in-one kit.

Ordering Information

Catalog # Description

Premixed, All-in-One Kits

LQ0-0000S6KL81S **Bio-Plex Pro Cell Signaling MAPK Panel**, 9-plex, 1 x 96 includes coupled magnetic beads and detection antibodies, cell lysis buffer, cell lysis factor QG, cell wash buffer, bead resuspension buffer, detection antibody diluent, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions for detecting ATF-2 (Thr⁷¹), ERK1/2 (Thr²⁰²/Tyr²⁰⁴, Thr¹⁸⁵/Tyr¹⁸⁷), HSP27 (Ser⁷⁹), JNK (Thr¹⁸³/Tyr¹⁸⁵), MEK1 (Ser²¹⁷/Ser²²¹), p38 MAPK (Thr¹⁸⁰/Tyr¹⁸²), p53 (Ser¹⁵), p90 RSK (Ser³⁸⁰), and Stat3 (Ser⁷²⁷)

LQ0-0006JK0K0RR **Bio-Plex Pro Cell Signaling Akt Panel**, 8-plex, 1 x 96 includes coupled magnetic beads and detection antibodies, cell lysis buffer, cell lysis factor QG, cell wash buffer, bead resuspension buffer, detection antibody diluent, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions for detecting Akt (Ser⁴⁷³), BAD (Ser¹³⁶), GSK-3 α / β (Ser²¹/Ser⁹), IRS-1 (Ser⁶³⁶/Ser⁶³⁹), mTOR (Ser²⁴⁴⁸), p70 S6 Kinase (Thr³⁸⁹), PTEN (Ser³⁸⁰), and S6 Ribosomal Protein (Ser²³⁵/Ser²³⁶)

Bio-Plex Pro Cell Signaling Reagents

171-304006M **Bio-Plex Pro Cell Signaling Reagent Kit**, 1 x 96-well, includes cell lysis buffer, cell lysis factor QG, cell wash buffer, bead resuspension buffer, detection antibody diluent, wash buffer, streptavidin-PE, 96-well flat bottom plate, sealing tape, and instructions

171-304515 **Bio-Plex Pro Cell Signaling Wash Buffer**, 330 ml, for 1 x 96-well assay, for use with Bio-Plex Pro cell signaling assays only, compatible with both magnetic and vacuum separation methods

Wash Stations

300-34376 **Bio-Plex Pro Wash Station**, includes magnetic plate carrier, waste bottle, 2 buffer bottles

Accessories

171-020100 **Bio-Plex Handheld Magnetic Washer**, includes magnetic washer and adjustment hex tools for use in manual wash steps for all Bio-Plex magnetic assays

171-304502 **Filter Plate**, pkg of 1, 96-well plate, with clear plastic lid and tray, for Bio-Plex assays using the vacuum wash method, sealing tape not included

Software

171-001510 **Bio-Plex Data Pro™ Software with Bio-Plex Manager™ Software**, Bio-Plex Data Pro software (5 seats), for multi-experiment analysis and advanced data visualization, and Bio-Plex Manager software (5 seats), for instrument data evaluation and optimization. CDs and security HASP key included

171-001513 **Bio-Plex Data Pro Software**, (5 seats), for multi-experiment analysis and advanced data visualization

171-STND01 **Bio-Plex Manager Software**, includes 1 user desktop license, to analyze Bio-Plex data and generate protocols, does not operate the instrument

Catalog # Description

Phosphoprotein Singleplex Sets*

1 x 96-well, includes coupled magnetic beads and detection antibodies for detecting phosphorylated protein, requires cell signaling reagent kit and optional lysate controls

171-V50001M Akt (Ser⁴⁷³)**
 171-V50002M Akt (Thr³⁰⁸)**
 171-V50024M ATF-2 (Thr⁷¹)**
 171-V50025M BAD (Ser¹³⁶)
 171-V50026M Btk (Tyr²²³)**
 171-V50027M c-Abl (Tyr²⁴⁵)**
 171-V50003M c-Jun (Ser⁶³)**
 171-V50028M CREB (Ser¹³³)
 171-V50004M EGFR (Tyr¹⁰⁶⁸)**
 171-V50005M EGFR (Tyr¹¹⁷³)**
 171-V50006M ERK1/2 (Thr²⁰²/Tyr²⁰⁴, Thr¹⁸⁵/Tyr¹⁸⁷)**
 171-V50007M GSK-3 α / β (Ser²¹/Ser⁹)
 171-V50008M HER-2 (Tyr¹²⁴⁸)**
 171-V50029M HSP27 (Ser⁷⁹)
 171-V50009M IGF-1R (Tyr¹¹³¹)**
 171-V50031M IR- β (Tyr¹¹⁴⁶)**
 171-V50030M IRS-1 (Ser⁶³⁶/Ser⁶³⁹)**
 171-V50010M I κ B- α (Ser³²/Ser³⁶)**
 171-V50011M JNK (Thr¹⁸³/Tyr¹⁸⁵)
 171-V50032M Lyn (Tyr⁵⁰⁷)**
 171-V50012M MEK1 (Ser²¹⁷/Ser²²¹)**
 171-V50033M mTOR (Ser²⁴⁴⁸)
 171-V50013M NF- κ B p65 (Ser⁵³⁶)**
 171-V50014M p38 MAPK (Thr¹⁸⁰/Tyr¹⁸²)**
 171-V50034M p53 (Ser¹⁵)
 171-V50016M p70 S6 Kinase (Thr³⁸⁹)**
 171-V50015M p70 S6 Kinase (Thr⁴²¹/Ser⁴²⁴)**
 171-V50035M p90 RSK (Ser³⁸⁰)
 171-V50017M PDGFR- α (Tyr⁷⁵⁴)**
 171-V50018M PDGFR- β (Tyr⁷⁵¹)**
 171-V50036M PI3K p85 (Tyr⁴⁵⁸)**
 171-V50037M PTEN (Ser³⁸⁰)
 171-V50038M S6 Ribosomal Protein (Ser²³⁵/Ser²³⁶)
 171-V50019M Smad2 (Ser⁴⁶⁵/Ser⁴⁶⁷)
 171-V50039M Src (Tyr⁴¹⁶)**
 171-V50020M Stat1 (Tyr⁷⁰¹)**
 171-V50021M Stat3 (Ser⁷²⁷)**
 171-V50022M Stat3 (Tyr⁷⁰⁵)**
 171-V50040M Syk (Tyr³⁵²)**
 171-V50023M VEGFR-2 (Tyr¹¹⁷⁵)
 171-V50041M ZAP-70 (Tyr³¹⁹)**

Total Target Singleplex Sets*

171-V60001M Akt**
 171-V60012M Btk
 171-V60002M c-Jun
 171-V60013M CREB
 171-V60003M Erk1/2
 171-V60004M GSK-3 β
 171-V60005M Her2
 171-V60014M IGF-1R
 171-V60006M I κ B- α
 171-V60007M JNK
 171-V60008M MEK1
 171-V60015M mTOR
 171-V60009M p38 MAPK**
 171-V60010M p70 S6 Kinase
 171-V60016M PTEN
 171-V60011M Smad2
 171-V60017M Src
 171-V60018M ZAP-70

Housekeeping Protein Sets*

171-V60020M β -Actin**
 171-V60019M Human GAPDH

* Refer to Table 4 to select lysate controls for multiplexing.

** Analytes with known cross-reactivity. Refer to Table 3 when selecting analytes for multiplexing.

Assay Multiplexability

Cross-reactivity between antibodies should be considered when multiplexing immunoassays to obtain the best results. While these assays have been extensively optimized to minimize nonspecific binding, homologies among proteins in the same family, especially in the phosphorylated domains, make the multiplexing of some assays impractical. It is not recommended to multiplex assays with shared bead regions and 10% or greater cross-reactivity. Refer to Table 3 for assay multiplexability.

Table 3. Assay multiplexability. Only assays with known cross-reactivity are shown.

Assays	Cross-Reactivity >10%
β-Actin	Akt (Ser ⁴⁷³), Akt (Thr ³⁰⁸), Stat3 (Tyr ⁷⁰⁵)
Akt (Total)	Smad2
Akt (Ser ⁴⁷³)	β-Actin, Akt (Thr ³⁰⁸)
Akt (Thr ³⁰⁸)	β-Actin, ATF-2 (Thr ⁷¹), Stat3 (Tyr ⁷⁰⁵), Akt (Ser ⁴⁷³)
ATF-2 (Thr ⁷¹)	Akt (Thr ³⁰⁸), c-Jun (Ser ⁶³)
Btk (Tyr ²²³)	HER-2 (Tyr ¹²⁴⁸), Lyn (Tyr ⁵⁰⁷), PI3K p85 (Tyr ⁴⁵⁸), Src (Tyr ⁴¹⁶)
c-Abl (Tyr ²⁴⁵)	EGFR (Tyr ¹¹⁷³), IR-β (Tyr ¹¹⁴⁶), PI3K p85 (Tyr ⁴⁵⁸)
c-Jun (Ser ⁶³)	ATF-2 (Thr ⁷¹), P70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)
EGFR (Tyr ¹⁰⁶⁸)	HER-2 (Tyr ¹²⁴⁸), IGF-1R (Tyr ¹¹³¹), Stat1 (Tyr ⁷⁰¹), Stat3 (Tyr ⁷⁰⁵), EGFR (Tyr ¹¹⁷³)
EGFR (Tyr ¹¹⁷³)	c-Abl (Tyr ²⁴⁵), HER-2 (Tyr ¹²⁴⁸), IGF-1R (Tyr ¹¹³¹), Stat1 (Tyr ⁷⁰¹), EGFR (Tyr ¹⁰⁶⁸)
Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	HER-2 (Tyr ¹²⁴⁸)
HER-2 (Tyr ¹²⁴⁸)	Btk (Tyr ²²³), EGFR (Tyr ¹⁰⁶⁸), EGFR (Tyr ¹¹⁷³), Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷), IGF-1R (Tyr ¹¹³¹), Lyn (Tyr ⁵⁰⁷), PDGFR-α (Tyr ⁷⁵⁴), PDGFR-β (Tyr ⁷⁵¹), PI3K p85 (Tyr ⁴⁵⁸), Stat1 (Tyr ⁷⁰¹)
IGF-1R (Tyr ¹¹³¹)	EGFR (Tyr ¹⁰⁶⁸), EGFR (Tyr ¹¹⁷³), HER-2 (Tyr ¹²⁴⁸), IR-β (Tyr ¹¹⁴⁶), PDGFR-α (Tyr ⁷⁵⁴), PDGFR-β (Tyr ⁷⁵¹)
IR-β (Tyr ¹¹⁴⁶)	c-Abl (Tyr ²⁴⁵), IGF-1R (Tyr ¹¹³¹), Lyn (Tyr ⁵⁰⁷), PI3K p85 (Tyr ⁴⁵⁸), Src (Tyr ⁴¹⁶), ZAP-70 (Tyr ³¹⁹)
IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	PI3K p85 (Tyr ⁴⁵⁸)
IκB-α (Ser ³² /Ser ³⁶)	NF-κB p65 (Ser ⁵³⁶)
Lyn (Tyr ⁵⁰⁷)	Btk (Tyr ²²³), HER-2 (Tyr ¹²⁴⁸), IR-β (Tyr ¹¹⁴⁶), PI3K p85 (Tyr ⁴⁵⁸)
MEK1 (Ser ²¹⁷ /Ser ²²¹)	P70 S6 Kinase (Thr ³⁸⁹), P70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)
NF-κB p65 (Ser ⁵³⁶)	IκB-α (Ser ³² /Ser ³⁶), Stat3 (Tyr ⁷⁰⁵)
p38 MAPK (Total)	c-Jun
p70 S6 Kinase (Thr ³⁸⁹)	MEK1 (Ser ²¹⁷ /Ser ²²¹), p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)
p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)	c-Jun (Ser ⁶³), MEK1 (Ser ²¹⁷ /Ser ²²¹), p70 S6 Kinase (Thr ³⁸⁹)
PDGFR-α (Tyr ⁷⁵⁴)	HER-2 (Tyr ¹²⁴⁸), IGF-1R (Tyr ¹¹³¹), PDGFR-β (Tyr ⁷⁵¹)
PDGFR-β (Tyr ⁷⁵¹)	HER-2 (Tyr ¹²⁴⁸), IGF-1R (Tyr ¹¹³¹), PDGFR-α (Tyr ⁷⁵⁴)
PI3K p85 (Tyr ⁴⁵⁸)	Btk (Tyr ²²³), c-Abl (Tyr ²⁴⁵), HER-2 (Tyr ¹²⁴⁸), IR-β (Tyr ¹¹⁴⁶), IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹), Lyn (Tyr ⁵⁰⁷), Syk (Tyr ³⁵²), ZAP-70 (Y ³¹⁹)
Src (Tyr ⁴¹⁶)	Btk (Tyr ²²³), IR-β (Tyr ¹¹⁴⁶)
Stat1 (Tyr ⁷⁰¹)	EGFR (Tyr ¹⁰⁶⁸), EGFR (Tyr ¹¹⁷³), HER-2 (Tyr ¹²⁴⁸), Stat3 (Ser ⁷²⁷), Stat3 (Tyr ⁷⁰⁵)
Stat3 (Ser ⁷²⁷)	Stat1 (Tyr ⁷⁰¹), Stat3 (Tyr ⁷⁰⁵)
Stat3 (Tyr ⁷⁰⁵)	β-Actin, Akt (Thr ³⁰⁸), EGFR (Tyr ¹⁰⁶⁸), NF-κB p65 (Ser ⁵³⁶), Stat1 (Tyr ⁷⁰¹), Stat3 (Ser ⁷²⁷)
Syk (Tyr ³⁵²)	PI3K p85 (Tyr ⁴⁵⁸)
ZAP-70 (Tyr ³¹⁹)	IR-β (Tyr ¹¹⁴⁶), PI3K p85 (Tyr ⁴⁵⁸)

Cell Lysate Controls

To identify the appropriate controls for your Bio-Plex Pro cell signaling phosphoprotein, total target, or housekeeping protein assay of interest, refer to Table 4.

Table 4. Bio-Rad lysate controls for phosphoprotein, total target, and housekeeping protein assays.

Phosphoprotein	Lysate Control	Catalog #	Total Target	Lysate Control	Catalog #
Akt (Ser ⁴⁷³)			Total Akt		
Akt (Thr ³⁰⁸)			Total Erk1/2		
Erk1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	EGF-treated HEK-293	171-YZ0001	Total GSK-3β		
GSK-3α/β (Ser ²¹ /Ser ⁹)			Total IκB-α		
MEK1 (Ser ²¹⁷ /Ser ²²¹)			Total JNK		
ATF-2 (Thr ⁷¹)			Total MEK1	Untreated HeLa	171-YZT002
c-Jun (Ser ⁶³)			Total mTOR		
CREB (Ser ¹³³)	UV-treated HEK-293	171-YZ0009	Total p38 MAPK		
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)			Total p70 S6 Kinase		
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)			Total PTEN		
p53 (Ser ¹⁵)			Total Smad2		
BAD (Ser ¹³⁶)			Total IGF-1R		
IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)			Total Btk	H ₂ O ₂ -treated Ramos	171-YZ0011
mTOR (Ser ²⁴⁴⁸)	PDGF-treated NIH3T3	171-YZ0007	Total ZAP-70	H ₂ O ₂ -treated Jurkat	171-YZ0012
PDGFR-α (Tyr ⁷⁵⁴)			Total c-Jun	Untreated HEK-293	171-YZT001
PDGFR-β (Tyr ⁷⁵¹)			Total CREB		
PTEN (Ser ³⁸⁰)			Total HER-2	EGF-treated SK-BR-3	171-YZ0003
Btk (Tyr ²²³)			Total Src	Src-transfected NIH3T3	171-YZ0013
Lyn (Tyr ⁵⁰⁷)	H ₂ O ₂ -treated Ramos	171-YZ0011	Negative control for all total target assays	Detection antibody diluent*	
PI3K p85 (Tyr ⁴⁵⁸)					
Syk (Tyr ³⁵²)					
c-Abl (Tyr ²⁴⁵)	Untreated K-562	171-YZT003			
EGFR (Tyr ¹⁰⁶⁸)					
EGFR (Tyr ¹¹⁷³)	EGF-treated HeLa	171-YZ0002			
HER-2 (Tyr ¹²⁴⁸)					
HSP27 (Ser ⁷⁸)					
p90 RSK (Ser ³⁸⁰)	EGF-treated SK-BR-3	171-YZ0003			
S6 Ribosomal Protein (Ser ²³⁵ /Ser ²³⁶)					
IGF-1R (Tyr ¹¹³¹)	IGF-1-treated HEK-293	171-YZ0005			
IR-β (Tyr ¹¹⁴⁶)					
IκB-α (Ser ³² /Ser ³⁶)					
NF-κB p65 (Ser ⁵³⁶)	TNF-α-treated HeLa	171-YZ0008			
Smad2 (Ser ⁴⁶⁵ /Ser ⁴⁶⁷)					
p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴)	NGFβ-treated PC12	171-YZ0006			
p70 S6 Kinase (Thr ³⁸⁹)					
Src (Tyr ⁴¹⁶)	Src-transfected NIH3T3	171-YZ0013			
Stat1 (Tyr ⁷⁰¹)					
Stat3 (Ser ⁷²⁷)	IFN-α-treated HeLa	171-YZ0004			
Stat3 (Tyr ⁷⁰⁵)					
VEGFR-2 (Tyr ¹¹⁷⁵)	VEGF-treated HUVEC	171-YZ0010			
ZAP-70 (Tyr ³¹⁹)	H ₂ O ₂ -treated Jurkat	171-YZ0012			
Negative control for all phosphoprotein assays	Phosphatase-treated HeLa	171-YZB001			

Housekeeping Protein	Lysate Control	Catalog #
β-Actin	Untreated HeLa	171-YZT002
Human GAPDH	Untreated HeLa	171-YZT002
Negative control for all housekeeping protein assays	Detection antibody diluent*	

* Detection antibody diluent is included with the Bio-Plex Pro cell signaling reagent kit (171-304006M).

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.



CST antibodies developed and validated for Bio-Plex cell signaling, phosphoprotein, and total target assays.

HASP is a trademark of Aladdin Knowledge Systems, Ltd.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000