


Ordering Information

Bio-Plex™ Phosphoprotein Assays and Bio-Plex Total Target Assays

What to Order

		
<p>Bio-Plex phosphoprotein assays and Bio-Plex total target assays — includes coupled beads, detection antibodies, and lysates to test for specific proteins in lysate samples.</p>	<p>Bio-Plex cell lysis kit — includes cell lysis buffer, cell wash buffer, factor 1, and factor 2; essential for optimal lysis of your culture or tissue samples.</p>	<p>Bio-Plex phosphoprotein detection reagent kit — includes assay buffer, wash buffer, streptavidin-PE, 96-well filter plate, sealing tape, and instructions required to prepare assays and acquire data.</p>

Note: Phosphoprotein and total target assays use the same reagent kit and cell lysis kit.

Available As Individual or x-Plex™ Assays

Bio-Plex phosphoprotein and total target assays are available individually or as x-Plex assays. Individual assays can be used independently or combined to test for a specific set of proteins in a single sample (catalog numbers listed below).

x-Plex assays are custom-premixed multiplex assays that are quality-tested at Bio-Rad and are a convenient format for repeat sample testing; catalog numbers are available online at www.bio-rad.com/bio-plex/x-plex/

Catalog #	Description
Reagent Kits	
171-304004	Bio-Plex Phosphoprotein Detection Reagent Kit, 1 x 96-well
171-304005	Bio-Plex Phosphoprotein Detection Reagent Kit, 10 x 96-well
Sample Preparation Kits	
171-304011	Bio-Plex Cell Lysis Kit, 1 x 96-well
171-304012	Bio-Plex Cell Lysis Kit, 10 x 96-well
Phosphoprotein Assays	
171-V21075	Bio-Plex Phospho-Akt (Ser ⁴⁷³) Assay, 1 x 96-well
171-V21620*	Bio-Plex Phospho-ATF-2 (Thr ⁷¹) Assay, 1 x 96-well
171-V24356*	Bio-Plex Phospho-c-Jun (Ser ⁶³) Assay, 1 x 96-well
171-V26119	Bio-Plex Phospho-CREB (Ser ¹³³) Assay, 1 x 96-well
171-V23120**	Bio-Plex Phospho-EGFR (Tyr) Assay, 1 x 96-well
171-V21938	Bio-Plex Phospho-ERK1 (Thr ²⁰² /Tyr ²⁰⁴) Assay, 1 x 96-well
171-V20438	Bio-Plex Phospho-ERK2 (Thr ¹⁸⁵ /Tyr ¹⁸⁷) Assay, 1 x 96-well
171-V22238	Bio-Plex Phospho-ERK1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷) Assay, 1 x 96-well
171-V23318	Bio-Plex Phospho-GSK-3 α/β (Ser ²¹ /Ser ⁹) Assay, 1 x 96-well
171-V24777*	Bio-Plex Phospho-Histone H3 (Ser ¹⁰) Assay, 1 x 96-well
171-V24551	Bio-Plex Phospho-HSP27 (Ser ⁷⁸) Assay, 1 x 96-well
171-V20758*	Bio-Plex Phospho-I κ B- α (Ser ³² /Ser ³⁶) Assay, 1 x 96-well
171-V25576	Bio-Plex Phospho-IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹) Assay, 1 x 96-well
171-V21034	Bio-Plex Phospho-JNK (Thr ¹⁸³ /Tyr ¹⁸⁵) Assay, 1 x 96-well
171-V25340	Bio-Plex Phospho-MEK1 (Ser ²¹⁷ /Ser ²²¹) Assay, 1 x 96-well
171-V24937*	Bio-Plex Phospho-NF- κ B p65 (Ser ⁵³⁶) Assay, 1 x 96-well
171-V21336	Bio-Plex Phospho-p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²) Assay, 1 x 96-well
171-V25153	Bio-Plex Phospho-p53 (Ser ¹⁵) Assay, 1 x 96-well
171-V24155	Bio-Plex Phospho-p70 S6 Kinase (Thr ⁴²¹ /Ser ⁴²⁴) Assay, 1 x 96-well

Catalog #	Description
Phosphoprotein Assays (cont.)	
171-V23535	Bio-Plex Phospho-p90RSK (Thr ³⁵⁹ /Ser ³⁶³) Assay, 1 x 96-well
171-V25957*	Bio-Plex Phospho-PDGFR (Tyr ⁷⁵¹) Assay, 1 x 96-well
171-V25374**	Bio-Plex Phospho-S6 Ribosomal Protein (Ser ²³⁵ /Ser ²³⁶) Assay, 1 x 96-well
171-V23732*	Bio-Plex Phospho-STAT2 (Tyr ⁶⁸⁹) Assay, 1 x 96-well
171-V22552*	Bio-Plex Phospho-STAT3 (Tyr ⁷⁰⁵) Assay, 1 x 96-well
171-V25772*	Bio-Plex Phospho-STAT6 (Tyr ⁶⁴¹) Assay, 1 x 96-well
171-V23973	Bio-Plex Phospho-TrkA (Tyr ⁴⁹⁰) Assay, 1 x 96-well
Total Target Assays	
171-V31075	Bio-Plex Total Akt Assay, 1 x 96-well
171-V31620*	Bio-Plex Total ATF-2 Assay, 1 x 96-well
171-V34356*	Bio-Plex Total c-Jun Assay, 1 x 96-well
171-V36119	Bio-Plex Total CREB Assay, 1 x 96-well
171-V30438	Bio-Plex Total ERK2 Assay, 1 x 96-well
171-V32238	Bio-Plex Total ERK1/2 Assay, 1 x 96-well
171-V34551	Bio-Plex Total HSP27 Assay, 1 x 96-well
171-V30758	Bio-Plex Total I κ B- α Assay, 1 x 96-well
171-V31034	Bio-Plex Total JNK Assay, 1 x 96-well
171-V35340	Bio-Plex Total MEK1 Assay, 1 x 96-well
171-V31336	Bio-Plex Total p38 MAPK Assay, 1 x 96-well
171-V33535	Bio-Plex Total p90RSK Assay, 1 x 96-well

* Refer to multiplexing matrix for multiplexing capabilities of specific assay combinations.

** This assay cannot be multiplexed.

Bio-Plex Phosphoprotein Assays

Multiplexable Assay Matrix

Phosphoprotein Assay	Tested Antibody Species*	Akt	ATF-2	c-Jun	CREB	ERK1	ERK2	ERK1/2	GSK-3 α/β	Histone H3	HSP27	I κ B- α	IRS-1	JNK	MEK1	NF- κ B p65	p38 MAPK	p53	p70 S6 Kinase	p90RSK	PDGF Receptor- β	STAT2	STAT3	STAT6	TrkA	
Akt	H, R																									
ATF-2	H, M																									
c-Jun	H, M, R																									
CREB	H, M, R																									
ERK1	H, M, R																									
ERK2	H, M, R																									
ERK1/2	H, M																									
GSK-3 α/β	H, M, R																									
Histone H3	H, M, R, K																									
HSP27	H, K																									
I κ B- α	H																									
IRS-1	H, M, R																									
JNK	H, M																									
MEK1	H, M, R, K																									
NF- κ B p65	H																									
p38 MAPK	H, M, R																									
p53	H																									
p70 S6 kinase	H, M, R																									
p90RSK	H, M, R																									
PDGF receptor- β	H, M, R																									
STAT2	H																									
STAT3	H, M, R																									
STAT6	H, M, B																									
TrkA	H, M, R																									

Singleplex Assays

Phosphoprotein Assay	Tested Antibody Species*
EGFR	H
S6 ribosomal protein	H, M, R

* B, bovine; H, human; K, monkey; M, mouse; R, rat.

Bio-Plex Total Target Assays

Multiplexable Assay Matrix

Total Target Assay	Tested Antibody Species*	Akt	ATF-2	c-Jun	CREB	ERK2	ERK1/2	HSP27	I κ B- α	JNK	MEK1	p38 MAPK	p90RSK
Akt	H, M, R												
ATF-2	H												
c-Jun	H, M, R												
CREB	H, M, R												
ERK2	H, M, R												
ERK1/2	H, M												
HSP27	H, K												
I κ B- α	H												
JNK	H												
MEK1	H, M, R												
p38 MAPK	H, M												
p90RSK	H, M, R												

* H, human; K, monkey; M, mouse; R, rat.



** Representative data based on lysates provided with the assay (HEK293, HeLa, PC12, and NIH 3T3 cell lines).



Bio-Rad Laboratories, Inc.

Life Science Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000