

**Catalog #** \_\_\_\_\_

<input checked="" type="checkbox"/>	Assay	Bead Region
<input type="checkbox"/>	Phosphoprotein	
<input type="checkbox"/>	Total Target	
<input type="checkbox"/>	Akt (Ser473)	75
<input type="checkbox"/>	ATF-2 (Thr71)	20
<input type="checkbox"/>	c-Jun (Ser63)	56
<input type="checkbox"/>	CREB (Ser133)	19
<input type="checkbox"/>	EGFR (Tyr)	20
<input type="checkbox"/>	ERK1 (Thr202/Tyr204)	38
<input type="checkbox"/>	ERK2 (Thr185/Tyr187)	38
<input type="checkbox"/>	ERK1/2 (Thr202/Tyr204, Thr185/Tyr187)	38
<input type="checkbox"/>	GSK-3 $\alpha/\beta$ (Ser21/Ser9)	18
<input type="checkbox"/>	Histone H3 (Ser10)	77
<input type="checkbox"/>	HSP27 (Ser78)	51

<input checked="" type="checkbox"/>	Assay	Bead Region
<input type="checkbox"/>	I $\kappa$ B- $\alpha$ (Ser32/Ser36)	58
<input type="checkbox"/>	IRS-1 (Ser636/Ser639)	76
<input type="checkbox"/>	JNK (Thr183/Tyr185)	34
<input type="checkbox"/>	MEK1 (Ser217/Ser221)	40
<input type="checkbox"/>	NF- $\kappa$ B p65 (Ser536)	37
<input type="checkbox"/>	PDGFR (Tyr751)	57
<input type="checkbox"/>	p38 MAPK (Thr180/Tyr182)	36
<input type="checkbox"/>	p53 (Ser15)	53
<input type="checkbox"/>	p70 S6 kinase (Thr421/Ser424)	55
<input type="checkbox"/>	p90RSK (Thr359/Ser363)	35
<input type="checkbox"/>	S6 Ribosomal Protein (Ser235/Ser236)	74
<input type="checkbox"/>	Stat2 (Tyr689)	32
<input type="checkbox"/>	Stat3 (Tyr705)	52
<input type="checkbox"/>	Stat6 (Tyr641)	72
<input type="checkbox"/>	TrkA (Tyr490)	73

For technical support, call your local Bio-Rad office, or in the US, call 1-800-4BIORAD (1-800-424-6723)

## Bio-Plex™ Phosphoprotein Detection Product Insert

For use with:  
 Bio-Plex phosphoprotein detection reagent kits  
 Catalog #171-304004 (one 96-well assay)  
 Catalog #171-304005 (ten 96-well assays)  
 and Bio-Plex cell lysis kits  
 Catalog #171-304011 (one 96-well assay)  
 Catalog #171-304012 (ten 96-well assays)

For research use only.  
 Not for diagnostic procedures.



## Bio-Plex™ Phosphoprotein Assays

Multiplexable Assay Matrix

Phosphoprotein Assay	Tested Antibody Species*	Akt	ATF-2	c-Jun	CREB	ERK1	ERK2	ERK1/2	GSK-3 $\alpha/\beta$	Histone H3	HSP27	I $\kappa$ B- $\alpha$	IRS-1	JNK	MEK1	NF- $\kappa$ B p65	p38 MAPK	p53	p70 S6 Kinase	p90RSK	PDGFR	STAT2	STAT3	STAT6	TrkA
Akt	H, M, R																								
ATF-2	H, M, R																								
c-Jun	H, M, R, K																								
CREB	H, M, R																								
ERK1	H, M, R																								
ERK2	H, M, R																								
ERK1/2	H, M																								
GSK-3 $\alpha/\beta$	H, M, R																								
Histone H3	H, M, R, K																								
HSP27	H, K																								
I $\kappa$ B- $\alpha$	H, M																								
IRS-1	H, M, R																								
JNK	H, M																								
MEK1	H, M, R, K																								
NF- $\kappa$ B p65	H, M, R																								
p38 MAPK	H, M, R																								
p53	H, M, R																								
p70 S6 kinase	H, M, R																								
p90RSK	H, M, R																								
PDGFR	H, M, R																								
STAT2	H, M, R																								
STAT3	H																								
STAT6	H, M																								
TrkA	H, M, R																								

### Singleplex Assays

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

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

## 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 $\kappa$ B- $\alpha$	JNK	MEK1	p38 MAPK	p90RSK
Akt	H, M, R												
ATF-2	H, M, R												
c-Jun	H												
CREB	H, M, R												
ERK2	H, M, R												
ERK1/2	H, M												
HSP27	H, K												
I $\kappa$ B- $\alpha$	H												
JNK	H												
MEK1	H, M, R												
p38 MAPK	H, M												
p90RSK	H, M, R												

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

### Multiplexing Key\*\*

Multiplexable  Not multiplexable

\*\* Representative data based on lysates provided with the assay.

For expanded views of these matrices, request bulletin 3071 or visit us on the Web at [www.bio-rad.com/products/phosphoproteins](http://www.bio-rad.com/products/phosphoproteins)

## Instructions

See Bio-Plex phosphoprotein detection manual.

## Specifications†

Reactive species

Human, mouse, rat

Matrices

Cell culture lysate, tissue lysate

Sample volume

25 $\mu$ l

Limit of detection

<10  $\mu$ g/ml of total protein

Precision

Intra-assay CV <15%  
Intra-assay CV <25%

Multiplexing capability

Yes

Storage conditions

2–8°C for coupled beads and detection antibodies

-20°C for lysates (shipped separately)

Stability††

12 months

† Values vary by assay.

†† For x-Plex assays, check label for expiration date.