

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 α/β (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 κ B- α (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- κ 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-RAD

4110019 Rev F 0506

