

Bio-Plex® Phosphoprotein Assays

BIO-RAD

Multiplexable Assay Matrix

Phosphoprotein Assay	Tested Antibody Species*	Akt (Ser ⁴⁷³)	ATF-2 (Thr ⁷¹)	Bcr-Abl (Tyr ²⁴⁹)	c-Abl (Tyr ²⁴⁹)	c-Abl (Tyr ¹¹²)	c-Jun (Ser ⁶³)	CREB (Ser ¹³³)	ERK1 (Thr ²⁰² /Tyr ²⁰⁴)	ERK2 (Thr ¹⁸⁵ /Tyr ¹⁸⁷)	ERK1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	GSK-3α/β (Ser ²¹ /Ser ⁹)	Histone H3 (Ser ¹⁰)	HSP27 (Ser ⁷⁰)	IGF-IR (Tyr ¹¹³¹)	IκB-α (Ser ³² /Ser ³⁶)	IR-β (Tyr ¹¹⁴⁶)	IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	MEK1 (Ser ²¹⁷ /Ser ²²¹)	NF-κB p65 (Ser ⁵³⁶)	p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	p53 (Ser ¹⁵)	p53 (Ser ³⁷)	p53 (Ser ³⁹)	p70 S6 kinase (Thr ⁴²¹ /Ser ⁴²⁴)	p90RSK (Thr ³⁵⁹ /Ser ³⁶³)	PDGF receptor-β (Tyr ⁷⁵¹)	Src (Tyr ⁴¹⁶)	STAT2 (Tyr ⁶⁸⁹)	STAT3 (Ser ²⁷)	STAT3 (Tyr ⁷⁰⁵)	STAT6 (Tyr ⁶⁴¹)	TrkA (Tyr ⁴⁹⁰)	Tyk2 (Tyr ¹⁰⁵⁴ /Tyr ¹⁰⁵⁵)
Akt (Ser ⁴⁷³)	H, R																																		
ATF-2 (Thr ⁷¹)	H, M																																		
Bcr-Abl (Tyr ²⁴⁵)	H, M																																		
c-Abl (Tyr ²⁴⁵)	H, M																																		
c-Abl (Tyr ¹¹²)	H																																		
c-Jun (Ser ⁶³)	H, M, R																																		
CREB (Ser ¹³³)	H, M, R																																		
ERK1 (Thr ²⁰² /Tyr ²⁰⁴)	H, M, R																																		
ERK2 (Thr ¹⁸⁵ /Tyr ¹⁸⁷)	H, M, R																																		
ERK1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	H, M																																		
GSK-3α/β (Ser ²¹ /Ser ⁹)	H, M, R																																		
Histone H3 (Ser ¹⁰)	H, M, R, K																																		
HSP27 (Ser ⁷⁸)	H, K																																		
IGF-IR (Tyr ¹¹³¹)	H																																		
IκB-α (Ser ³² /Ser ³⁶)	H																																		
IR-β (Tyr ¹¹⁴⁶)	H, M, R																																		
IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	H, M, R																																		
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	H, M																																		
MEK1 (Ser ²¹⁷ /Ser ²²¹)	H, M, R, K																																		
NF-κB p65 (Ser ⁵³⁶)	H																																		
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	H, M, R																																		
p53 (Ser ¹⁵)	H																																		
p53 (Ser ³⁷)	H, K																																		
p53 (Ser ³⁹)	H, K																																		
p70 S6 kinase (Thr ⁴²¹ /Ser ⁴²⁴)	H, M, R																																		
p90RSK (Thr ³⁵⁹ /Ser ³⁶³)	H, M, R																																		
PDGF receptor-β (Tyr ⁷⁵¹)	H, M, R																																		
Src (Tyr ⁴¹⁶)	H, M, R																																		
STAT2 (Tyr ⁶⁸⁹)	H																																		
STAT3 (Ser ²⁷)	H, M, R																																		
STAT3 (Tyr ⁷⁰⁵)	H, M, R																																		
STAT6 (Tyr ⁶⁴¹)	H, M																																		
TrkA (Tyr ⁴⁹⁰)	H, M, R																																		
Tyk2 (Tyr ¹⁰⁵⁴ /Tyr ¹⁰⁵⁵)	H																																		

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

Phosphoprotein Singleplex Assay

Tested Antibody Species*	EGFR (Tyr)	S6 ribosomal protein (Ser ²³⁵ /Ser ²³⁶)
	H	H, M, R

Multiplexable Assay Matrix

Total Target Assay	Tested Antibody Species*													
		Akt	ATF-2	Bcr-Abl	c-Jun	CREB	ERK2	ERK1/2	HSP27	I _K B- α	JNK	MEK1	p38 MAPK	p53
Akt	H, M, R													
ATF-2	H													
Bcr-Abl	H, M, R													
c-Jun	H, M, R													
CREB	H, M, R													
ERK2	H, M, R													
ERK1/2	H, M													
HSP27	H, K													
I _K B- α	H													
JNK	H													
MEK1	H, M, R													
p38 MAPK	H, M													
p53	H													
p90RSK	H, M, R													

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

Multiplexing Key**

Multiplexable				Not multiplexable

** Representative data are based on lysates provided with the assay (in-house cell lines).