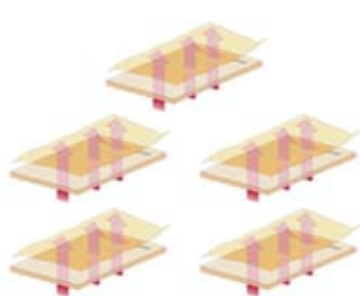
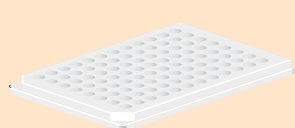

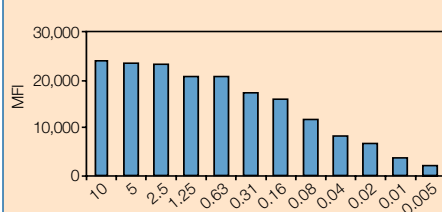


Bio-Plex[®] Phosphoprotein Detection

How would you like to have higher sensitivity than western blots?

The Power of xMAP Technology

Side-by-Side Comparison: Analyzing 5 Phosphoproteins in 12 Samples

	Western Blot	xMAP Technology
Number of phosphoproteins	5	5
Number of samples	12	12
Total data points	60	60
Number of gels or 96-well plates	5 	1 
Data points per gel or plate	12	60
Total time required	~8 hr + one overnight incubation	3 hr + one overnight incubation
Sample volume	125 µl	25 µl
Dynamic range	Limited	50–900 µg/ml total protein
Sensitivity*	 Total protein, µg Western blot of total HSP27 (COS-7 cell lysate)	 MFI Total protein, µg Bio-Plex total HSP27 assay (COS-7 cell lysate)

* For more details, see bulletin 5483.

Bio-Plex Phosphoprotein and Total Target Assays

Available Assays*

Assays	Phosphoprotein	Total
Akt Signaling		
Akt (Ser ⁴⁷³)	●	●
GSK-3 α / β (Ser ²¹ /Ser ⁹)	●	
Cell Cycle/Checkpoint Control		
p53 (Ser ¹⁵)	●	● new
p53 (Ser ³⁷)	● new	● new
p53 (Ser ⁴⁶)	● new	● new
Chromatin Regulation/Acetylation		
Histone H3 (Ser ¹⁰)	●	
Immunology/Inflammation		
I κ B- α (Ser ³² /Ser ³⁶)	●	●
NF- κ B p65 (Ser ⁵³⁶)	●	
STAT2 (Tyr ⁶⁸⁹)	●	
STAT3 (Ser ⁷²⁷)	● new	
STAT3 (Tyr ⁷⁰⁵)	●	
STAT6 (Tyr ⁶⁴¹)	●	
Tyk2 (Tyr ¹⁰⁵⁴ /Tyr ¹⁰⁵⁵)	● new	
Glucose/Energy Metabolism		
IGF-IR (Tyr ¹¹³¹)	● new	
IR- β (Tyr ¹¹⁴⁶)	● new	
IRS-1 (Ser ⁶³⁶ /Ser ⁶³⁹)	●	
MAP Kinase Signaling		
ATF-2 (Thr ⁷¹)	●	●
c-Jun (Ser ⁶³)	●	●
ERK1 (Thr ²⁰² /Tyr ²⁰⁴)	●	
ERK2 (Thr ¹⁸⁵ /Tyr ¹⁸⁷)	●	●
ERK1/2 (Thr ²⁰² /Tyr ²⁰⁴ , Thr ¹⁸⁵ /Tyr ¹⁸⁷)	●	●
HSP27 (Ser ⁷⁸)	●	●
JNK (Thr ¹⁸³ /Tyr ¹⁸⁵)	●	●
MEK1 (Ser ²¹⁷ /Ser ²²¹)	●	●
p38 MAPK (Thr ¹⁸⁰ /Tyr ¹⁸²)	●	●
p90RSK (Thr ³⁵⁹ /Ser ³⁶³)	●	●
Neuroscience		
CREB (Ser ¹³³)	●	●
TrkA (Tyr ⁴⁹⁰)	●	
Translational Control		
p70 S6 kinase (Thr ⁴²¹ /Ser ⁴²⁴)	●	
S6 ribosomal protein (Ser ²³⁵ /Ser ²³⁶)	●	
Tyrosine Kinases		
Bcr-Abl (Tyr ²⁴⁵)	● new	● new
c-Abl (Tyr ²⁴⁵)	● new	
c-Abl (Tyr ⁴¹²)	● new	
EGFR (Tyr)	●	
PDGF receptor- β (Tyr ⁷⁵¹)	●	
Src (Tyr ⁴¹⁶)	● new	

* You can design an x-Plex assay for combinations of assays. For more information, go to www.bio-rad.com/bio-plex/x-plex/ or contact your local Bio-Rad sales representative.

For more information, visit us on the Web at www.bio-rad.com/products/phosphoproteins/
Call toll free at 1-800-4BIORAD (1-800-424-6723);
outside the U.S., contact your local sales office.

Cell Signaling Technology and the Cell Signaling Technology logo are trademarks of Cell Signaling Technology, Inc. xMAP is a trademark of Luminex Corporation. The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.



CST antibodies exclusively developed and validated for Bio-Plex phosphoprotein and total target assays.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 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 Israel 03 963 6050
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
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 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000