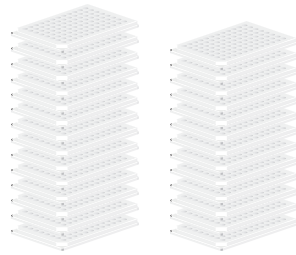



Bio-Plex® Cytokine Assays

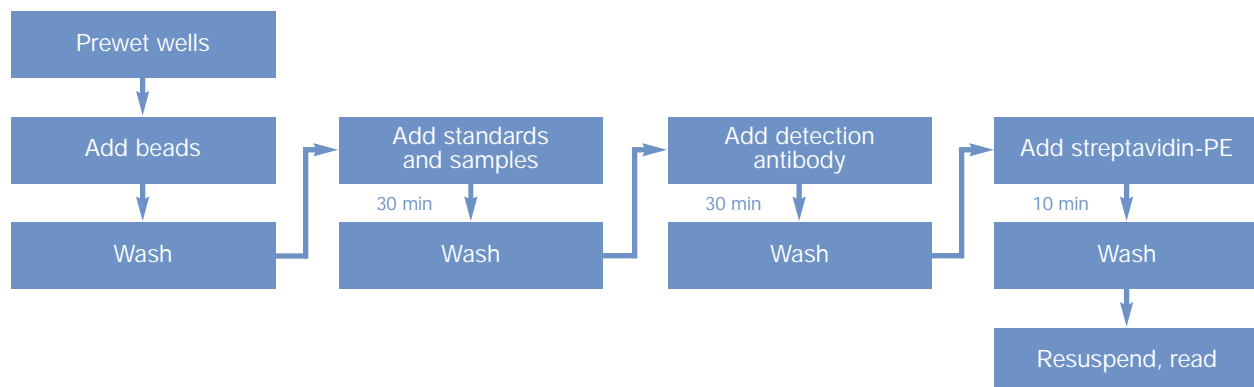
How would you like to have 2,160 data points in one 96-well plate?

The Power of the Bio-Plex System

Side-by-Side Comparison: Analyzing 27 Cytokines in 80 Samples

	ELISA	Bio-Plex
Number of cytokines	27	27
Number of samples	80	80
Total data points	2,160	2,160
Number of 96-well plates	27 	1 
Data points per plate	80	2,160
Total time required	> 60 hr	3 hr
Sample volume	Serum or plasma, >1 ml* Cell culture supernatant, >1 ml*	Serum or plasma, 12.5 µl Cell culture supernatant, 50 µl
Assay range	Serum or plasma, 2–3,000 pg/ml Cell culture supernatant, 2–3,000 pg/ml	Serum or plasma, ~0.2–3,200 pg/ml Cell culture supernatant, ~2–32,000 pg/ml

* Based on 50 µl/well of sample.



Bio-Plex cytokine assay methodology.



Cytokines Available in Bio-Plex Assays

Singleplex assays, preconfigured multiplex panels, and custom-mixed multiplex panels are available.

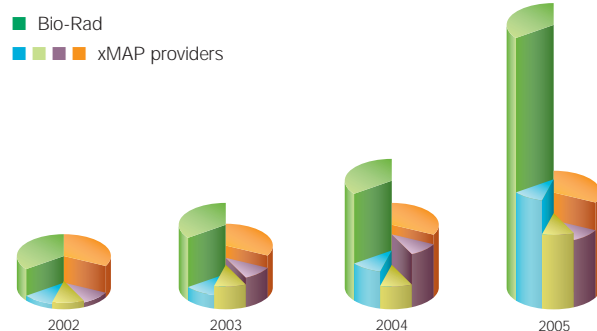
Assays	Human	Mouse	Rat
IL-1 α	●	●	●
IL-1 β	●	●	●
IL-1ra	●		
IL-2	●	●	●
IL-2R α	●		
IL-3	●	●	
IL-4	●	●	●
IL-5	●	●	
IL-6	●	●	●
IL-7	●		
IL-8	●		
IL-9	●	●	
IL-10	●	●	●
IL-12 (p40)	●	●	
IL-12 (p70)	●	●	
IL-13	●	●	
IL-15	●	●	
IL-16	●		
IL-17	●	●	
IL-18*	●	●	
Basic FGF	●	●	
CTACK	●	●	
Eotaxin	●	●	
G-CSF	●	●	
GM-CSF	●	●	●
GRO- α	●		

Assays	Human	Mouse	Rat
HGF	●		
ICAM-1	●		
IFN- α 2	●		
IFN- γ	●	●	●
IP-10	●		
KC		●	
LIF	●	●	
MCP-1 (MCAF)	●	●	
MCP-3	●		
M-CSF	●	●	
MIF	●		
MIG	●	●	
MIP-1 α	●	●	
MIP-1 β	●	●	
MIP-2		●	
β -NGF	●		
PDGF-BB	●	●	
RANTES	●	●	
SCF	●		
SCGF- β	●		
SDF-1 α	●		
TNF- α	●	●	●
TNF- β	●		
TRAIL	●		
VCAM-1	●		
VEGF	●	●	

* IL-18 is not available as a singleplex assay, only as part of an assay panel.

Visit us on the Web at www.bio-rad.com/bio-plex/.
Call toll free at 1-800-4BIORAD (1-800-424-6723);
outside the US, contact your local sales office.

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.



xMAP cytokine assay citations published from 2002 through 2005.



**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 83282000