



## Bio-Plex® Assays — Washing Steps Made Easy

The Bio-Plex Pro Wash Station eliminates manual wash steps from Bio-Plex Assays, making Bio-Plex Assays as easy as enzyme-linked immunosorbent assays (ELISAs). Benefits of the Bio-Plex Pro Wash Station include:

- Improved laboratory productivity
- Reliable and reproducible results
- Optimized onboard protocols

Visit [bio-rad.com/web/Bio-PlexProWash](http://bio-rad.com/web/Bio-PlexProWash) for more information.



## Performance

The Bio-Plex Pro Wash Station helps decrease variability and increase consistency in results between experiments. The wash station is specifically designed to perform Bio-Plex Assay wash steps, but is compatible with any standard xMAP Assay.

### Representative results using the Bio-Plex® Precision Pro™

**Human Cytokine Assay.** Experiments determining intra- and inter-assay variability were carried out using the magnetic plate carrier on the Bio-Plex Pro Wash Station. Intra-assay variability was calculated from triplicate experiments. Inter-assay variability was determined from three separate experiments performed on different days.

Human Target	Mean Observed Concentration, pg/ml	Standard Deviation	%CV
<b>Intra-Assay Variability</b>			
IL-1 $\beta$	6.39	0.17	2.63
IL-4	13.22	0.59	4.50
IL-6	465.09	13.63	2.93
IL-12 (p70)	10.12	1.14	11.29
IFN- $\gamma$	25.03	0.86	3.45
TNF- $\alpha$	11.84	0.78	6.63
<b>Inter-Assay Variability</b>			
IL-1 $\beta$	6.23	0.40	6.45
IL-4	16.68	0.21	1.28
IL-6	467.62	17.12	3.66
IL-12 (p70)	16.34	1.51	9.26
IFN- $\gamma$	32.40	2.30	7.11
TNF- $\alpha$	13.91	0.22	1.56

CV, coefficient of variation.

## Specifications

Magnetic field	4,500 $\pm$ 200 gauss
Number of dispensing channels	8
Residual volume magnetic wash	<4 $\mu$ l
Dispensing accuracy	$\leq$ 2%
Dispensing uniformity	CV $\leq$ 4% over the plate
Dimensions (W $\times$ D $\times$ H)	28 $\times$ 37 $\times$ 18 cm
Weight	6.6 kg

Note: Additional features include plate-shaking capability and PC-based software for custom protocol development.

## Ordering Information

Catalog #	Description
<b>Bio-Plex Pro Wash Station</b> 30034376	<b>Bio-Plex Pro Wash Station</b> , includes magnetic plate carrier, waste bottle, 2 buffer bottles
<b>Bio-Plex Accessories</b> 171020100	<b>Bio-Plex Handheld Magnetic Washer</b> , includes magnetic washer and hex tools for use in manual wash steps for all Bio-Plex Magnetic Assays
171025001*	<b>Bio-Plex Pro Flat Bottom Plates</b> , 40 $\times$ 96-well plates
<b>Bulk Buffer</b> 171304500	<b>Bio-Plex Wash Buffer</b> , 1.5 L

\* Required for washing Bio-Plex Pro Assays or other magnetic bead-based assays using the magnetic plate carrier.

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.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](http://bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550  
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00  
France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 972 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 (0)318 540 666  
New Zealand 64 9 415 2280 Norway 47 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 (0) 861 246 723 Spain 34 91 590 5200 Sweden 46 08 555 12700 Switzerland 41 026 674 55 05  
Taiwan 886 2 2578 7189 Thailand 66 662 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

