

# Bio-Plex Pro™ Human Cytokine Screening Panel

## Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Human Cytokine Screening Panel	10000092045

New users can download the complete manual, which includes detailed instructions and a list of kit components, at [bio-rad.com/bio-plex](http://bio-rad.com/bio-plex).

## Initial Preparation

1. Plan the plate layout.
2. Start up/warm up the Bio-Plex® System (30 min).
  - Bring diluents, including wash buffer, assay buffer, standard diluent HB, detection antibody diluent HB, and sample diluent HB, to room temperature (RT). Keep the other items on ice until needed
    - Mix by inversion to ensure all salts are in solution
    - Prepare 1x wash buffer: dilute **1 part** 10x wash buffer (60 ml) with **9 parts** dH<sub>2</sub>O (540 ml)
  - Begin to thaw frozen samples
3. Prepare sample dilution according to the guidelines provided in the table below. It is important to centrifuge serum or plasma samples at **1,000 x g** for **15 min** at **4°C** to remove particulates from all samples prior to use.

Sample Type	Recommended Sample Dilution	Diluent
Serum and plasma	(1:4)	Sample diluent
Culture media and fluids	User defined	Diluent + 0.5% BSA (w/v)

**Note:** ICAM-1 and VCAM-1 require higher dilution for serum and plasma (recommended 100-fold).

4. Calibrate the Bio-Plex System within Bio-Plex Manager™ Software.
5. Reconstitute the standards and control by adding 250 µl of standard diluent HB to each. Vortex at medium speed for 5 sec and incubate all vials on ice for precisely 30 min.

## Bio-Plex Pro Human Cytokine Screening Panel Quick Guide

- Prepare a fourfold standard dilution series and blank as shown below. Vortex at medium speed for 5 sec between liquid transfers.

**Note:** Standards are at S1 concentration after reconstitution and the controls are ready to use after reconstitution. Controls are included with the fixed panel only.

- Vortex** coupled beads at medium speed for **30 sec** and **dilute to 1x** in Bio-Plex Assay Buffer as shown in the following table. Protect from light.

### Premixed Panels

# of Wells	10x Beads, $\mu\text{l}$	Assay Buffer, $\mu\text{l}$	Total Volume, $\mu\text{l}$
96	570	5,130	5,700

### Singleplex Assays

# of Wells	Singleplex #1 and #2 20x Beads, $\mu\text{l}$	Assay Buffer, $\mu\text{l}$	Total Volume, $\mu\text{l}$
96	285	5,130	5,700

**Note:** 20x singleplex beads allow multiplexing up to 20 analytes.

Transfer Volume, $\mu\text{l}$	250	50	50	50	50	50	50	50	50	
Reconstituted Standard										
Standard Diluent, $\mu\text{l}$	0	150	150	150	150	150	150	150	150	150

## Running the Assay

- Vortex** the diluted (1x) beads. **Add 50  $\mu\text{l}$**  to each well of the assay plate.
- Wash the plate two times** with **100  $\mu\text{l}$**  Bio-Plex Wash Buffer.
- Vortex** samples, standards, blank, and control. **Add 50  $\mu\text{l}$**  to each well.
- Cover plate with sealing tape. Incubate on shaker at **850  $\pm$  50 rpm** at RT for **30 min**.
- With 10 min left in the incubation, **vortex** detection antibodies for **5 sec** and quick-spin to collect liquid. **Dilute to 1x** as shown below.

## Bio-Plex Pro Human Cytokine Screening Panel Quick Guide

### Premixed Panels

# of Wells	10x Detection Ab, $\mu\text{l}$	10x Detection Ab Diluent HB, $\mu\text{l}$	Total Volume, $\mu\text{l}$
96	300	2,700	3,000

### Singleplex Assays

# of Wells	Singleplex #1 and #2 20x Detection Ab, $\mu\text{l}$	Detection Ab Diluent HB, $\mu\text{l}$	Total Volume, $\mu\text{l}$
96	150	2,700	3,000

**Note:** 20x singleplex beads allow multiplexing up to 20 analytes.

6. Wash the plate three times with 100  $\mu\text{l}$  wash buffer.
7. Vortex the diluted (1x) detection antibodies. Add 25  $\mu\text{l}$  to each well.
8. Cover plate with sealing tape and incubate at  $850 \pm 50$  rpm for 30 min at RT. Meanwhile, prepare Bio-Plex Manager Software protocol; enter standard S1 values and units provided in the assay kit.
9. With 10 min left in the incubation, vortex 100x streptavidin-PE (SA-PE) for 5 sec and quick-spin to collect liquid. Dilute to 1x as shown below and protect from light.

# of Wells	100x SA-PE, $\mu\text{l}$	Assay Buffer, $\mu\text{l}$	Total Volume, $\mu\text{l}$
96	60	5,940	6,000

10. Wash the plate three times with 100  $\mu\text{l}$  wash buffer.
11. Vortex the diluted (1x) SA-PE. Add 50  $\mu\text{l}$  to each well.
12. Cover plate with sealing tape and incubate at  $850 \pm 50$  rpm for 10 min at RT.
13. Wash the plate three times with 100  $\mu\text{l}$  wash buffer.
14. Resuspend beads in 125  $\mu\text{l}$  assay buffer. Cover and shake at  $850 \pm 50$  rpm for 30 sec.
15. Remove the sealing tape and read plate using the settings below.

Instrument	RP1 (PMT)	DD Gates	Bead Events
Bio-Plex 3D*	Standard	Select MagPlex Beads	50
Bio-Plex 100, 200*	Low	5,000 (low); 25,000 (high)	50
Bio-Plex® MAGPIX™	N/A, use default instrument settings		

\* Or similar Luminex-based system.

The observed concentration ranges of the control apply only when standards and controls are prepared using the provided Bio-Plex Standard Diluent HB.

# Bio-Plex Pro Human Cytokine Screening Panel Quick Guide

## Assay Workflow

Add 50  $\mu$ l 1x beads to wells

Wash buffer: 2 x 200  $\mu$ l

Add 50  $\mu$ l standards, samples, controls; incubate on shaker at 850 rpm for 30 min at RT

Wash buffer: 3 x 100  $\mu$ l

Add 25  $\mu$ l 1x detection antibody; incubate on shaker at 850 rpm for 30 min at RT

Wash buffer: 3 x 100  $\mu$ l

Add 50  $\mu$ l 1x streptavidin-PE; incubate on shaker at 850 rpm for 10 min at RT

Wash buffer: 3 x 100  $\mu$ l

Resuspend in 125  $\mu$ l assay buffer; shake at 850 rpm for 30 sec

Acquire data on Bio-Plex® System

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation. Luminex, MAGPIX, and MagPlex are trademarks of Luminex Corporation.

**BIO-RAD**

**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 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192  
**Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190  
**India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000  
**Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 310 318 540 666  
**New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191  
**Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188  
**South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00  
**Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311  
**United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

