

Bio-Plex Pro Human Cytokine Assays

Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Cytokine, Chemokine, and Growth Factor Assays	10000111560

This guide can be used to prepare and run a full 1 x 96-well assay plate. New users can download the complete manual, which includes detailed instructions and a list of kit components, at bio-rad.com/bio-plex.

Initial Preparation

1. Plan the plate layout.
2. Start up/warm up the Bio-Plex Multiplex Immunoassay 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** distilled water (540 ml)
 - Begin to thaw frozen samples
3. Prepare the sample dilution according to the guidelines provided in the following table. 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% bovine serum albumin (BSA) (w/v)

Note: ICAM-1 and VCAM-1 require higher dilution for serum and plasma (recommended 100-fold). Refer to the Bio-Plex Pro Cytokine, Chemokine, and Growth Factor Assays Instruction Manual (#10000111560) for detailed sample preparation recommendations.

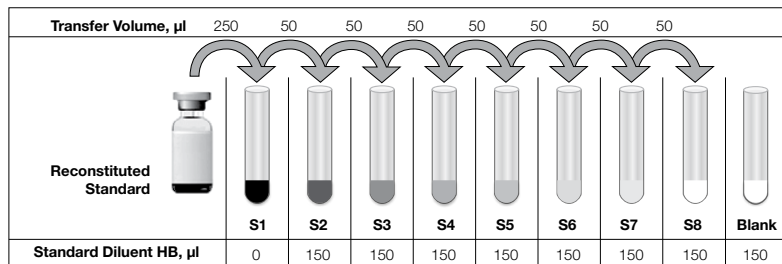
4. Calibrate the Bio-Plex System within Bio-Plex Manager Software.
5. Reconstitute the standards and control by adding 250 µl of standard diluent HB to each. **Vortex** at medium speed for **5 sec** and incubate all vials on ice for precisely **30 min**.

Bio-Plex Pro Human Cytokine Assays Quick Guide

6. Prepare a fourfold standard dilution series and blank as shown.

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.



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

Premixed Panels

Number of Wells	10x Beads, μl	Assay Buffer, μl	Total Volume, μl
96	570	5,130	5,700

Singleplex Assays

Number of Wells	Singleplex #1	Singleplex #2	Assay Buffer, μl	Total Volume, μl
	20x Beads, μl	20x Beads, μl		
96	285	285	5,130	5,700

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

Running the Assay

- Vortex** the diluted (1x) beads. **Add 50 μl** to each well of the assay plate.
- Wash the plate two times** with **100 μl** Bio-Plex Wash Buffer.
- Vortex** the samples, standards, blank, and control. **Add 50 μl** to each well.
- Cover the 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** the detection antibodies for **5 sec** and quick-spin to collect liquid. **Dilute to 1x** as shown.

Bio-Plex Pro Human Cytokine Assays Quick Guide

Premixed Panels

Number of Wells	10x Detection Antibodies, μl	10x Detection Antibody Diluent HB, μl	Total Volume, μl
96	300	2,700	3,000

Singleplex Assays

Number of Wells	Singleplex #1 20x Detection Antibodies, μl	Singleplex #2 20x Detection Antibodies, μl	Detection Antibody Diluent HB, μl	Total Volume, μl
96	150	150	2,700	3,000

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

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

Number of Wells	100x SA-PE, μl	Assay Buffer, μl	Total Volume, μl
96	60	5,940	6,000

- Wash the plate three times with 100 μl wash buffer.
- Vortex the diluted (1x) SA-PE. Add 50 μl to each well.
- Cover the plate with sealing tape and incubate at 850 ± 50 rpm for 10 min at RT.
- Wash the plate three times with 100 μl wash buffer.
- Resuspend the beads in 125 μl assay buffer. Cover and shake at 850 ± 50 rpm for 30 sec.
- Remove the sealing tape and read plate using the following settings:

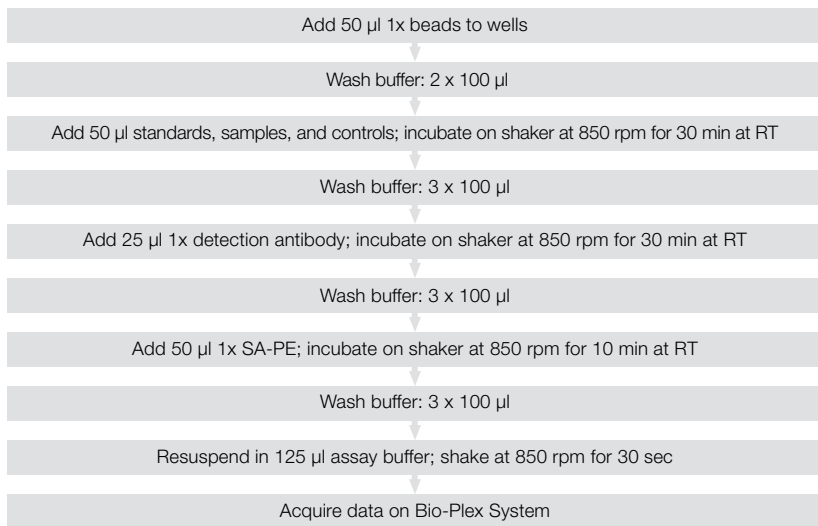
Instrument	RP1 (PMT)	DD Gates	Bead Events
Bio-Plex 3D*	Standard	Select MagPlex Beads	50
Bio-Plex 100, 200*	Low	5,000 (low); 25,000 (high)	50
Bio-Plex MAGPIX	N/A, use default instrument settings		

* Or similar Luminex System.

Bio-Plex Pro Human Cytokine Assays Quick Guide

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

Assay Workflow



BIO-RAD and BIO-PLEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation. Luminex is a trademark of Luminex Corporation. All trademarks used herein are the property of their respective owner.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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** 31 0 318 540 666
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

