

Bio-Plex Pro Human Inflammation Panel I Assays

Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Human Inflammation and Treg Cytokine Assays	10044282

This guide can be used to prepare and run a full 1 x 96-well assay plate. For more information on a given step, refer to the complete instruction manual. New users can go to bio-rad.com/bio-plex and download the manual, which includes detailed instructions and a list of kit components.

IMPORTANT! Pay close attention to **vortexing**, **shaking**, and **incubation** instructions. Deviation from the protocol may result in low assay signal and assay variability.

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
 - Begin to thaw frozen samples
3. Prime the wash station for a flat bottom plate or set a vacuum manifold to -1 to -3" Hg for filter plate.
4. Calibrate the Bio-Plex System by following the prompts in Bio-Plex Manager Software. This can be done now or during an assay incubation step.
5. Prepare 1x wash buffer. Mix 10x stock by inversion to ensure all salts are in solution. Then dilute **1 part** 10x wash buffer with **9 parts** distilled water.
6. Reconstitute the vial of standards in **781 µl** standard diluent HB (or a diluent similar to your sample matrix). Reconstitute the vial of control in **250 µl** of standard diluent HB, as shown. **Vortex** at medium speed for **5 sec** and incubate all vials **on ice** for precisely **30 min**.

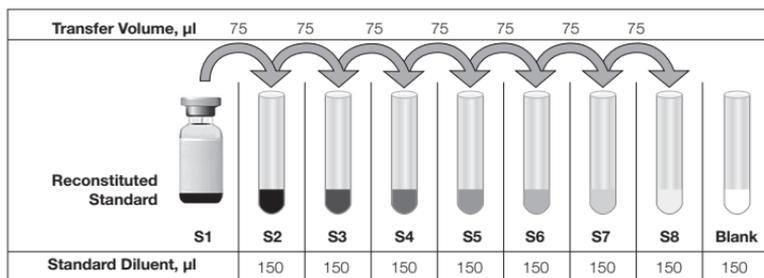
Bio-Plex Pro Human Inflammation Panel I Assays Quick Guide

Sample Type	Diluent for Standard and Controls*	Add Bovine Serum Albumin (BSA)
Serum and plasma	Standard diluent HB	None
Culture media, with serum	Culture media	None
Culture media, serum-free	Culture media	To 0.5% final
Lavage, lysate, other fluids	Sample diluent HB	To 0.5% final

* If using diluents other than the standard diluent HB provided, users must establish their own control ranges.

7. Prepare a threefold standard dilution series and blank as shown. **Vortex** at medium speed for **5 sec** between liquid transfers.

Note: The controls are ready to use after reconstitution. No dilution is needed. Controls are included with the fixed panel only.



8. After thawing samples, prepare them according to the following guidelines.

Sample Type	Dilution Factor	Diluent
Serum and plasma	1:4	Sample diluent HB
Fluids	User defined	Diluent + 0.5% BSA w/v

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

Number of Wells	10x Beads, µl	Assay Buffer, µl	Total Volume, µl
96	575	5,175	5,750

Running the Assay

Note: Make sure all assay components are at RT before pipetting. **Vortex** at medium speed.

1. Prewet the filter plate with **100 µl** Bio-Plex Assay Buffer (skip for flat bottom plate).
2. **Vortex** the diluted (1x) beads. Add **50 µl** to each well of the assay plate.
3. **Wash the plate two times** with **100 µl** Bio-Plex Wash Buffer.
4. **Vortex** samples, standards, blank, and control. Add **50 µl** to each well.
5. Cover the plate with sealing tape and protect from light with aluminum foil. Incubate on shaker at **850 ± 50 rpm** for **1 hr** at RT.
6. With 10 min left in the incubation, **vortex** detection antibodies for **15 sec** and quick-spin to collect liquid. Dilute to 1x as shown.

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

7. **Wash the plate three times** with **100 µl** wash buffer.
8. **Vortex** the diluted (1x) detection antibodies. Add **25 µl** to each well.
9. Cover with aluminum foil and incubate at **850 ± 50 rpm** in the dark for **30 min** at RT. Meanwhile, prepare the Bio-Plex Manager Software protocol; enter standard S1 values and units provided in the assay kit.
10. 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

11. **Wash the plate three times** with **100 µl** wash buffer.
12. **Vortex** the diluted (1x) SA-PE. Add **50 µl** to each well.
13. Cover with aluminum foil and incubate at **850 ± 50 rpm** in the dark for **10 min** at RT.

Bio-Plex Pro Human Inflammation Panel I Assays Quick Guide

14. Wash the plate three times with 100 µl wash buffer.
15. Resuspend the beads in 125 µl assay buffer. Cover and shake at 850 ± 50 rpm for 30 sec.
16. Remove the sealing tape and read plate using the following settings.

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

* Or similar Luminex System.

17. Control is included with the fixed panel only. If the control was run, then compare the observed concentration against the ranges provided in the assay kit. Ranges apply only when standards and controls are prepared in Bio-Plex Standard Diluent HB.

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

