

Bio-Plex Pro Human Chemokine Assays

Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Human Chemokine Assays	10031990

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 the 10x wash buffer, assay buffer, and diluents to room temperature (RT). Keep the other items on ice until needed
 - Begin to thaw frozen samples
 - Prepare 1x wash buffer. Mix 10x stock by inversion to ensure all salts are in solution. Then dilute **1 part** 10x wash buffer (60 ml) with **9 parts** distilled water (540 ml)
3. Prime the wash station for a flat bottom 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. After thawing samples, prepare them according to the following guidelines.

Bio-Plex Pro Human Chemokine Assays Quick Guide

- 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	20x Beads, μl	Assay Buffer, μl	Total Volume, μl
96	288	5,472	5,760

Running the Assay

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

- 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** samples, standards, blank, and control. Add **50 μl** to each well.
- Cover the plate with sealing tape and protect from light with aluminum foil. Incubate on shaker at **850 \pm 50 rpm** for **1 hr** at RT.
- 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	20x Detection Antibodies, μl	Detection Antibody Diluent HB, μl	Total Volume, μl
96	150	2,850	3,000

- Wash the plate three times** with **100 μl** wash buffer.
- Vortex** the diluted (1x) detection antibodies. Add **25 μl** to each well.
- Cover and incubate at **850 \pm 50 rpm**, as in step 4, 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.

Bio-Plex Pro Human Chemokine Assays Quick Guide

- Cover and incubate at 850 ± 50 rpm, as in step 4, for 10 min at RT.
- Wash the plate three times with 100 μ l wash buffer.
- Resuspend the beads in 125 μ l 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 200*	Low	5,000 (low); 25,000 (high)	50
Bio-Plex MAGPIX	N/A, use default instrument settings	Select MagPlex Beads	N/A

* Or similar Luminex System.

- Quality controls are included with the fixed panel only. If they were run, then compare the observed concentrations against the ranges provided in the assay kit. Ranges apply only when standard 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

