

Bio-Plex Pro™ Assays

Chemokine 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 download the manual, which includes detailed instructions and a list of kit components, at www.bio-rad.com/bio-plex.

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® system (30 min).
 - Bring the 10x wash buffer, assay buffer, and diluents to room temperature (RT). Keep other items on ice until needed
 - Begin to thaw frozen samples
3. Prime wash station for flat bottom plate or set vacuum manifold to -1 to -3" Hg for filter plate.
4. Calibrate the Bio-Plex system by following the prompts within the 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 dH₂O.

Bio-Plex Pro Human Chemokine Quick Guide

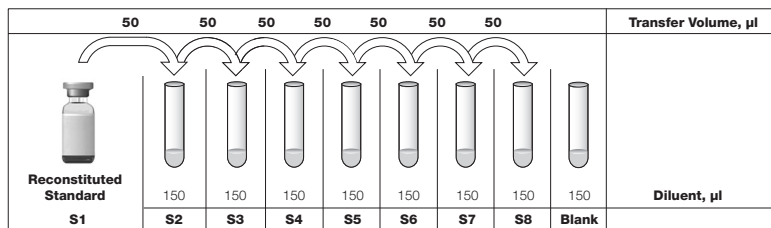
6. Reconstitute the vial of standards in **781 μ l** of a diluent similar to your final sample type or matrix. Reconstitute the vial of quality controls in **250 μ l** of the same diluent, as shown below. **Vortex** at medium speed for **5 sec** and incubate all vials at once **on ice** for **30 min**.

Sample Type	Diluent for Standards and Controls*	Add 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, then users must establish their own control ranges.

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

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



8. After thawing samples, prepare according to the guidelines shown below.

Sample Type	Diluent	Add BSA	Recommended Sample Dilution
Serum and plasma	Sample diluent HB	None	Fourfold (1:4)
Culture media, with serum	Culture media	None	Neat to 1:10
Culture media, serum-free	Culture media	To 0.5% final	Neat to 1:10
Lavage, other fluids	Sample diluent HB	To 0.5% final	User optimized
Lysate	Sample diluent HB	To 0.5% final	User optimized (at least 1:2 for 50 to 500 μ g/ml final protein)

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

# of Wells	20x Beads, μ l	Assay Buffer, μ l	Total Volume, μ l
96	288	5,472	5,760

Running the Assay

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

1. Prewet filter plate with **100 µl** Bio-Plex assay buffer (skip for flat bottom).
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 controls. **Add 50 µl** to each well.
5. Cover plate with sealing tape and protect from light with aluminum foil. Incubate on shaker at **850 ± 50 rpm** at RT for **1 hr**.
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 below.

# of Wells	20x Detection Ab, µl	Detection Ab Diluent HB, µl	Total Volume, µl
96	150	2,850	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 and incubate** at **850 ± 50 rpm**, as described above, in the dark for **30 min** at RT. Meanwhile, prepare 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-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, 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 and incubate** at **850 ± 50 rpm**, as described above, in the dark for **10 min** at RT.

Bio-Plex Pro Human Chemokine Quick Guide

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

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

* A similar Luminex-based system may be used.

17. 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.

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation. Luminex and MagPlex are trademarks of Luminex Corporation.



P10031991

BIO-RAD

**Bio-Rad
Laboratories, Inc.**



Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800
Austria 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500
Czech Republic 420 241 430 532 **Denmark** 44 52 10 00
Finland 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0
Greece 30 210 9532 220 **Hong Kong** 852 2789 3300
Hungary 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460
Mexico 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 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 27 861 246 723 **Spain** 34 91 590 5200
Sweden 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000