

Bio-Plex Pro Mouse Cytokine Assays

Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Mouse Cytokine, Chemokine, and Growth Factor Assays	10000142118

This guide can be used to prepare and run a full 1 x 96-well assay plate. Refer to the complete instruction manual for more information on a given step. 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 assay buffer, wash buffer, and sample diluent to room temperature (RT) and keep other items on ice until needed. Begin to thaw frozen samples.
3. After thawing samples, prepare them according to the following guidelines.

Assay	Serum and Plasma		Culture Supernatant and Other Fluids		Cell and Tissue Lysate	
	Dilution	Diluent	Dilution	Diluent	Dilution	Diluent
Mouse cytokine	1:4	Bio-Plex sample diluent	User optimized	Cell culture medium or buffer similar to sample*	User optimized (1:2 lysates at 200–900 µg/ml protein)	Bio-Plex sample diluent

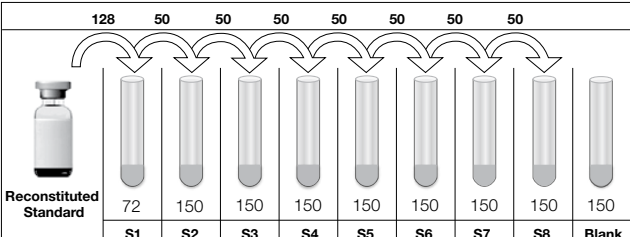
* If samples are serum-free, add bovine serum albumin (BSA) to 0.5% final w/v.

Bio-Plex Pro Mouse Cytokine Assays Quick Guide

- Prime the wash station for a flat bottom plate. 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).
- 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.
- Reconstitute a single vial of standards in **500 µl** of a diluent similar to the final sample type or matrix. **Vortex** for **5 sec** and incubate **on ice** for **30 min**.

Sample Type	Diluent for Standards	Add BSA
Serum and plasma	Standard diluent	None
Culture media with serum	Culture media	None
Culture media, serum-free	Culture media	To 0.5% final (w/v)

- Prepare a fourfold standard dilution series and blank as shown. **Vortex** for **5 sec** between liquid transfers.

	128	50	50	50	50	50	50	50	50	Transfer Volume, µl
	72	150	150	150	150	150	150	150	150	
	S1	S2	S3	S4	S5	S6	S7	S8	Blank	Diluent, µl

Note: Change tips between each dilution.

- Vortex** the 10x coupled beads for **30 sec** and dilute to 1x in Bio-Plex Assay Buffer as shown. Protect from light.

Mouse Cytokine Assays

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

Running the Assay

1. **Vortex** the diluted (1x) beads for **10–20 seconds**. Add **50 µl** to each well of the assay plate.
2. **Wash the plate two times** with **100 µl** Bio-Plex Wash Buffer.
3. **Vortex** samples, standards, and blank. Add **50 µl** to each well.
4. Cover the plate with sealing tape and incubate on shaker at **850 ± 50 rpm** at RT. See the following table for incubation times for sample, detection antibody, and streptavidin-phycoerythrin (SA-PE).

Assay	Incubation Time		
	Sample	Detection Antibody	SA-PE
Bio-Plex Pro Mouse Cytokine	30 min	30 min	10 min

5. With 10 min left in the incubation, **vortex** the 10x detection antibody for **5 sec** and quick-spin to collect liquid. Dilute to 1x as shown.

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

6. **Wash the plate three times** with **100 µl** wash buffer.
7. **Vortex** the diluted (1x) detection antibody. Add **25 µl** to each well.
8. Repeat step 4. See table for incubation time. Meanwhile, prepare the Bio-Plex Manager Software protocol; enter standard S1 values provided in the assay kit.
9. With 10 min left in the incubation, **vortex** the 100x 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

10. **Wash the plate three times** with **100 µl** wash buffer.
11. **Vortex** the diluted (1x) SA-PE. Add **50 µl** to each well.
12. Repeat step 4. See table for incubation time.

Bio-Plex Pro Mouse Cytokine Assays Quick Guide

13. Wash the plate three times with 100 μ l wash buffer.
14. Resuspend the beads in 125 μ l assay buffer. Cover the plate as in step 4 and shake it at 850 \pm 50 rpm for 30 sec.
15. Remove the sealing tape and **read plate** using the following settings for optimal sensitivity. Use low photomultiplier tube (PMT) RP1 setting for the Bio-Plex 100, 200, 3D, or similar system. Use default instrument settings for the Bio-Plex MAGPIX Multiplex Reader.

Assay	Low PMT, RP1	High PMT, RP1
Bio-Plex Pro Mouse Cytokine	•	–

BIO-RAD and BIO-PLEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Lumindex Corporation.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800
Austria 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500
Czech Republic 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23
Finland 00 800 00 24 67 23 **France** 00 800 00 24 67 23
Germany 00 800 00 24 67 23 **Hong Kong** 852 2789 3300
Hungary 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460
Luxembourg 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280
Norway 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23
Portugal 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23
Singapore 65 6415 3188 **South Africa** 00 800 00 24 67 23
Spain 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189
Thailand 66 2 651 8311 **United Arab Emirates** 36 1 459 6150
United Kingdom 00 800 00 24 67 23

