

Bio-Plex Pro™ Assays

Quick Guide

For Use with	Instruction Manual #
Mouse and Rat Cytokine Assays	10014905

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 download the manual, which includes detailed instructions and a list of kit components, at bio-rad.com/web/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 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 as shown below.

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

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

4. Prime wash station for flat bottom plate. Prepare 1x wash buffer.

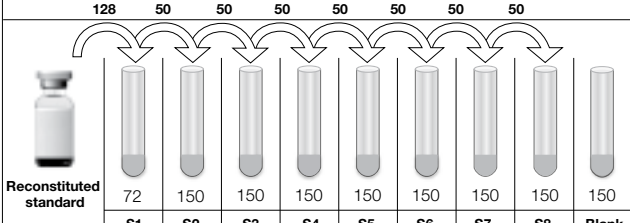
Bio-Plex Pro Assays Quick Guide

Mix 10x stock by inversion to ensure all salts are in solution. Then dilute 1 part 10x wash buffer (60 ml) with 9 parts dH₂O (540 ml).

- Calibrate the Bio-Plex System by following the prompts within 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 below. **Vortex for 5 sec** between liquid transfers. If mixing diabetes assays with cytokine assays, refer to the diabetes instruction manual.

	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 or 20x coupled beads for **30 sec** and dilute to 1x in Bio-Plex Assay Buffer as shown. Protect from light.

Mouse Cytokine Group I and II Assays

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

Mouse Cytokine Group III and Rat Cytokine Group I Assays

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

Running the Assay

- Vortex** the diluted (1x) beads for **10–20 sec**. 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. Add **50 µl** to each well.
- Cover plate with sealing tape and incubate on shaker at **850 ± 50 rpm** at RT. See table for incubation time.

Incubation times for sample, detection Ab, and SA-PE

Assay	Sample	Detection Ab	SA-PE
Bio-Plex Pro Mouse Cytokine (Groups I and II)	30 min	30 min	10 min
Bio-Plex Pro Mouse Cytokine (Group III)	1 hr	30 min	10 min
Bio-Plex Pro Rat Cytokine (Group I)	1 hr	30 min	10 min

- With 10 min left in the incubation, **vortex** the 10x or 20x detection antibodies for **5 sec** and quick-spin to collect liquid. Dilute to 1x as shown below.

# of Wells	Detection Ab, µl		Detection Ab Diluent HB, µl	Total Volume, µl
	10x	20x		
96	300	—	2,700	3,000
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.
- Repeat Step 4. See table for incubation time. Meanwhile, prepare Bio-Plex Manager Software protocol; enter standard S1 values from the assay kit.
- 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.

# 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 Assays Quick Guide

- Repeat Step 4. See table for incubation time.
- Wash the plate three times with **100 μ l** wash buffer.
- Resuspend beads in **125 μ l** assay buffer. Cover plate as in Step 4 and shake the plate at **850 \pm 50 rpm** for **30 sec.**
- Remove sealing tape and **read the plate** using the settings below. Use low PMT, RP1 setting for the Bio-Plex 100, 200, and 3D Systems. Use default instrument settings for the Bio-Plex[®] MAGPIX[™] System.

Settings for optimal sensitivity on the Bio-Plex 100, 200, 3D or similar system.

Assay	Low PMT, RP1	High PMT, RP1
Bio-Plex Pro Mouse Cytokine (Groups I, II, and III)	•	–
Bio-Plex Pro Rat Cytokine (Group I)	–	•

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

MAGPIX is a trademark of Luminex Corporation.

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 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500
Czech Republic 420 241 430 532 **Denmark** 45 44 52 10 00
Finland 358 09 804 22 00 **France** 33 01 47 95 69 65 **Germany** 49 89 31 884 0
Hong Kong 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **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 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 (0) 861 246 723 **Spain** 34 91 590 5200
Sweden 46 08 555 12700 **Switzerland** 41 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 2 651 8311
United Arab Emirates 971 4 8187300 **United Kingdom** 44 020 8328 2000

