



Bio-Plex Pro Mouse Th17 Cytokine Assays

Quick Guide

| For Use with | Instruction Manual # |
|--|----------------------|
| Bio-Plex Pro Cytokine, Chemokine, and Growth Factor Assays | 10000111560 |

This guide can be used to prepare and run a full 1 x 96-well assay plate. New users should download the complete manual, which includes detailed instructions and a list of kit components, at bio-rad.com/bio-plex.

Initial Preparation

- Plan the plate layout.
- Start up/warm up the Bio-Plex Multiplex Immunoassay System (**30 min**).
 - Bring diluents, including wash buffer, assay buffer, standard diluent, detection antibody diluent HP, and sample diluent, to room temperature (RT). Keep the other items on ice until needed
 - Mix by inversion to ensure all salts are in solution
 - Prepare 1x wash buffer: dilute **1 part** 10x wash buffer (60 ml) with **9 parts** distilled water (540 ml)
 - Begin to thaw the frozen samples
- Prepare the sample dilution according to the guidelines provided in the following table. It is important to centrifuge serum or plasma samples at **1,000 x g** for **15 min** at **4°C** to remove particulates from all samples prior to use.

| 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 | 1:100 | Bio-Plex Standard and Sample Diluents | User optimized | | | |

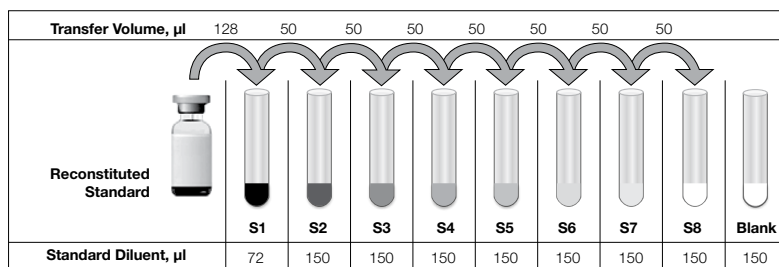
* If samples are serum-free, add bovine serum albumin (BSA) to 0.5% final w/v.

For example:

- For serum or plasma cytokine assays, dilute samples 1:4 by adding 40 µl sample + 120 µl Bio-Plex Sample Diluent
 - For serum or plasma ICAM-1 assays, dilute samples 1:100
 - First dilution (1:4): 10 µl sample + 30 µl Bio-Plex **Sample Diluent**
 - Second dilution (1:25): 5 µl from the first dilution + 120 µl Bio-Plex **Standard Diluent**
- Calibrate the Bio-Plex System in Bio-Plex Manager Software.
 - Reconstitute the standards and control by adding **500 µl** of standard diluent to each. **Vortex** at medium speed for **5 sec** and incubate all vials on ice for precisely **30 min**.
 - Prepare a fourfold standard dilution series and blank as shown. **Vortex** at medium speed for **5 sec** between liquid transfers.

Note: The control is ready to use after reconstitution. Controls are included with the fixed panel only.

Bio-Plex Pro Mouse Th17 Cytokine Assays Quick Guide



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

Premixed Panels

| Number of Wells | 10x Beads, µl | Assay Buffer, µl | Total Volume, µl |
|-----------------|---------------|------------------|------------------|
| 96 | 570 | 5,130 | 5,700 |

Singleplex Assays

| Number of Wells | Singleplex #1 | | Singleplex #2 | | Assay Buffer, µl | Total Volume, µl |
|-----------------|---------------|---------------|---------------|---------------|------------------|------------------|
| | 10x Beads, µl | 10x Beads, µl | 10x Beads, µl | 10x Beads, µl | | |
| 96 | 570 | 570 | 570 | 570 | 4,560 | 5,700 |

Note: 10x singleplex beads allow multiplexing up to ten analytes.

Running the Assay

- 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** the samples, standards, blank, and control. **Add 50 µl** to each well.
- Cover the plate with sealing tape. Incubate on shaker at **850 ± 50 rpm** at RT for **30 min**.
- With 10 min left in the incubation, **vortex** the detection antibodies for **5 sec** and quick-spin to collect liquid. **Dilute to 1x** as shown.

Premixed Panels

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

Singleplex Assays

| Number of Wells | Singleplex #1 | | Singleplex #2 | | Detection Antibody Diluent HP, µl | Total Volume, µl |
|-----------------|------------------------------|------------------------------|------------------------------|------------------------------|-----------------------------------|------------------|
| | 10x Detection Antibodies, µl | 10x Detection Antibodies, µl | 10x Detection Antibodies, µl | 10x Detection Antibodies, µl | | |
| 96 | 300 | 300 | 300 | 300 | 2,400 | 3,000 |

Note: 10x singleplex beads allow multiplexing up to ten analytes.

- Wash the plate three times** with **100 µl** wash buffer.
- Vortex** the diluted (1x) detection antibodies. **Add 25 µl** to each well.
- Cover the plate with sealing tape and incubate at **850 ± 50 rpm** 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.

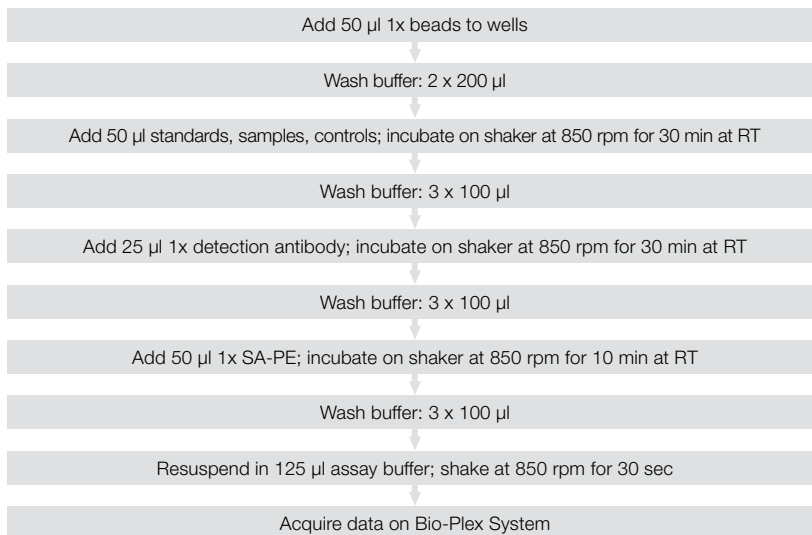
| Number of Wells | 100x SA-PE, µl | Assay Buffer, µl | Total Volume, µl |
|-----------------|----------------|------------------|------------------|
| 96 | 60 | 5,940 | 6,000 |

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. Cover the plate with sealing tape and incubate at 850 ± 50 rpm for 10 min at RT.
13. Wash the plate three times with 100 µl wash buffer.
14. Resuspend the beads in 125 µl assay buffer. Cover and shake at 850 ± 50 rpm for 30 sec.
15. 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 100, 200* | Low | 5,000 (low); 25,000 (high) | 50 |
| Luminex MAGPIX | N/A, use default instrument settings | | |

* Or similar Luminex System.

Assay Workflow



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
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 310 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 470 233 841 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

