

Bio-Plex Pro™ Assays

Quick Guide 1

For Use with*	Instruction Manual #
Bio-Plex Pro Human Cancer Biomarker Panel 1 and Panel 2	10022186
Bio-Plex Pro Human Th17 Cytokine Panel	10023381

* Includes premixed multiplex panels, singleplex sets, custom x-Plex™ assays, and Express assays.

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 corresponding section of 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. (**Section 1** of the complete instruction manual)
2. Start up/warm up the Bio-Plex® system (30 min). (**Section 2**)
 - Meanwhile, bring buffers and diluents to room temperature (RT)
 - Begin to thaw frozen samples
3. Prime wash station or set vacuum manifold to -1 to -3" Hg. (**Section 3**)
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. (**Section 2**)

Bio-Plex Pro Assay Quick Guide 1

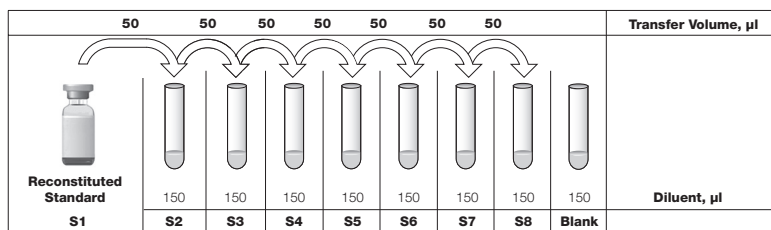
5. Reconstitute a single vial of standard in **781 μ l** of a diluent similar to your final sample type or matrix. Reconstitute the quality controls in **250 μ l** of the same diluent, as shown below. **Vortex** for **5 sec** and incubate all vials at once **on ice** for **30 min.** (Section 4)

Note: The quality controls are ready to use after reconstitution. No dilution is needed. Only premixed panels and x-Plex kits include controls.

Sample Type	Diluent for Standard 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 standard diluent HB, users must establish their own control ranges.

6. Prepare a fourfold standard dilution series and blank as shown below. **Vortex** for **5 sec** between liquid transfers. (Section 4)



7. After thawing samples, prepare as shown below. (Section 5)

Sample Type	Diluent	Add BSA	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, lysate, other fluids	Sample diluent HB	To 0.5% final	User optimized

8. **Vortex** coupled beads for **30 sec** and dilute to 1x in Bio-Plex assay buffer as shown below. Protect from light. (Section 6)

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

Running the Assay (all Section 7 in the manual unless otherwise noted)

Note: Make sure all assay components are at RT before proceeding.

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 and incubate** in the dark for **1 hr** at RT with vigorous shaking.
- Note:** Shake at **850 ± 50 rpm**. Ramp up to speed slowly to avoid splashing.
6. With 10 min left in the incubation, **vortex** detection antibodies for **5 sec** and quick-spin to collect liquid. Dilute to 1x as shown below.

# of Wells	20x Detection Ab, µl	Detection Ab Diluent, µl	Total Volume, µl
96	145	2,755	2,900

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** in the dark for **30 min** at RT. Shake at **850 ± 50 rpm**.
Meanwhile, prepare Bio-Plex Manager software protocol; enter standard S1 values provided in the assay kit. (**Section 8**)
10. With 10 min left in the incubation, **vortex** 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** in the dark for **10 min** at RT. Shake at **850 ± 50 rpm**.

Bio-Plex Pro Assay Quick Guide 1

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. (Section 8)

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

* Or similar Luminex-based system.

16. If quality controls were run, 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.

MAGPIX and MagPlex are trademarks of Luminex Corp.



P10022581



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** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000