

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)

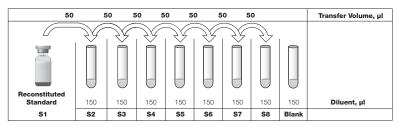
5. Reconstitute a single vial of standard in 781 μl of a diluent similar to your final sample type or matrix. Reconstitute the quality controls in 250 μl of the same diluent, as shown below. Vortex for 5 sec and incubate all vials at once on ice for 30 min. (Section 4)

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.

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, μl	Assay Buffer, µl	Total Volume, µl
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 \muI** 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, μΙ
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 µI 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, µI	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.

- 14. Wash the plate three times with 100 µl wash buffer.
- 15. Resuspend beads in 125 µl assay buffer. Cover and shake at $850 \pm 50 \text{ rpm for } 30 \text{ sec. (Section 8)}$

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.







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

10022581 Rev C Sia 1211