

Bio-Plex Pro™ Assays

Quick Guide 5

For use with	Instruction Manual #
TGF- β Assays	10024984

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 complete instruction manual. New users can download the manual 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

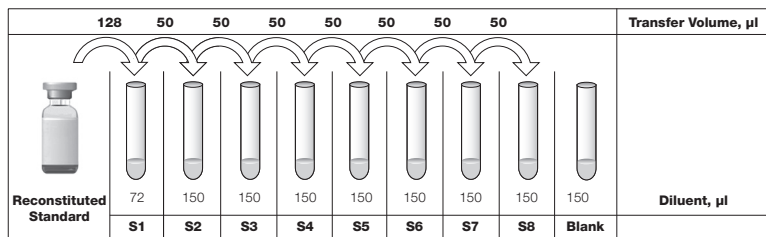
1. Plan the plate layout.
2. Start up/warm up the Bio-Plex® system (**30 min**).
 - Bring assay buffer, wash buffer, and sample diluent to room temperature (RT). Keep other items on ice until needed
 - Begin to thaw frozen samples
3. Prime wash station for flat bottom plate or set vacuum manifold to -1 to -3" Hg for filter plate.
4. Follow the prompts in Bio-Plex Manager™ software to calibrate the system. This can be done now or during an incubation step.
5. Mix 1 volume of Bio-Plex standard diluent with 3 volumes of Bio-Plex sample diluent (each supplied in the kit). The resulting solution is used for reconstitution and subsequent dilution of standards. This results in a serum-matrix based diluent that mimics the matrix in 1:16 diluted serum and plasma samples. For samples in serum-free media and other biological fluids, use a diluent that most closely matches the sample matrix. Add carrier protein such as BSA at a final concentration of 0.5% (w/v).

Bio-Plex Pro Assay Quick Guide 5

6. Reconstitute a single vial of standards in **500 µl** of a diluent similar to the final sample type or matrix as shown below. **Vortex** for **5 sec** and incubate **on ice** for **30 min**.

Sample Type	Diluent for Standard	Add BSA
Serum and plasma	Standard/sample diluent mix (1:3)	None
Culture media, with serum	Culture media	None
Culture media, serum-free	Culture media	To 0.5% final (w/v)
Lavage, lysate, other fluids	Sample diluent	To 0.5% final (w/v)

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



8. After thawing samples, activate by adding 1 volume of 1 N HCl to 5 volumes of sample. **Vortex**, incubate at RT for **10 min**. Neutralize by adding the same volume of base (1.2 N NaOH/0.5 M HEPES). After treatment, dilute samples as shown below.

Sample Type	Diluent	Add BSA	Sample Dilution
Serum and plasma	Sample diluent	None	1:16 final*
Culture media, with serum	Culture media	None	User optimized
Culture media, serum-free	Culture media	To 0.5% final	User optimized
Lavage, lysate, other fluids	Sample diluent	To 0.5% final	User optimized

* For example, activate 25 µl sample, neutralize, and bring to a final volume of 400 µl.

9. **Vortex** the 20x coupled beads for **30 sec** and dilute to 1x in Bio-Plex assay buffer as shown below. Protect from light.

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

Running the Assay

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 two times with **100 µl** Bio-Plex wash buffer.
4. **Vortex** samples, standards, blank. Add **50 µl** to each well.
5. Cover plate with sealing tape and protect from light with aluminum foil. Incubate on shaker at **850 ± 50 rpm** for **2 hr** at RT.

Note: **850 rpm** provides equivalent performance to recommended shaker settings in previous manuals (**1,100 rpm** for **30 sec**, **300 rpm** for incubation).

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

# of Wells	20x Detection Ab, µl	Detection Ab Diluent, µl	Total Volume, µl
96	150	2,850	3,000

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 at **850 ± 50 rpm**, as in Step 5, for **1 hr** at RT. Meanwhile, prepare Bio-Plex Manager software protocol; enter standard S1 values provided in the assay kit.
10. With **10 min** left in the incubation, **vortex** the 100x SA-PE for **5 sec**, quick-spin to collect liquid, and dilute to 1x as shown below. 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.

Bio-Plex Pro Assay Quick Guide 5

- Cover and incubate at **850 ± 50 rpm**, as in Step 5, for **30 min** at RT.
- Wash the plate three times with **100 µl** wash buffer.
- Resuspend beads in **125 µl** assay buffer. Cover and shake at **850 ± 50 rpm**, as in Step 5, for **30 sec**.
- Remove the sealing tape and **read plate** using the settings below.

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
Bio-Plex® MAGPIX™*	N/A, use default instrument settings		Default

* Or similar Luminex-based system.

MagPlex is a trademark of Luminex Corporation.

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



P10024986

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** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000