

Bio-Plex Pro™ Assays

Quick Guide 3

For Use With:

Product Family	Instruction Manual #
Bio-Plex Pro cell signaling assays	10024929

This guide describes how to prepare and run a full 1 x 96-well flat bottom plate using a magnetic assay workflow. 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. **Do Not** mix phosphoprotein assays with corresponding total target assays (e.g., phospho-Akt and total Akt). Use only the reagents supplied with cell signaling assay kits.

Sample Preparation

1. Prepare lysates from cell culture or tissue samples. To remove insoluble debris from lysates (≤ 2 ml), centrifuge at **15,000 x g** for **10 min** at 4°C in a benchtop microfuge. (**Section 1** of the complete instruction manual)
2. Measure protein concentration using the Bio-Rad® DC™ protein assay.
3. Adjust protein concentration to **20–200 µg/ml** in Bio-Plex® cell lysis buffer containing cell lysis factor QG and PMSF. For samples with unknown target expression or phosphorylation levels, start with **200 µg/ml** protein.
4. If necessary, store lysates at –70°C.
5. Thaw frozen lysates and keep on ice. Centrifuge again as above and bring to room temperature (RT) immediately before use.

BIO-RAD

Bio-Plex Pro Assay Quick Guide 3

Assay – Day 1

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

1. Plan the plate layout. (**Section 2** of the complete instruction manual)
2. Prime wash station or prepare handheld magnetic washer. (**Section 3**)
3. Reconstitute Bio-Rad cell lysates in **250 μ l** of diH₂O per vial. **Vortex** for **5 sec** and incubate at RT for **20 min**. Transfer the cell lysate to a centrifuge tube and centrifuge at **15,000 x g** for **10 min** at 4°C. Protein concentration is now **200 μ g/ml**. (**Section 4**)
4. **Vortex** coupled beads for **30 sec** and dilute to 1x in cell signaling wash buffer as shown below. Protect from light. (**Section 4**)

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

5. **Vortex** the diluted (1x) beads. Add **50 μ l** to each well of the assay plate.
6. Wash the plate two times with **200 μ l** wash buffer.
7. Gather samples, Bio-Rad cell lysates, and blank (detection antibody diluent alone). Add **50 μ l** to each well. Recommended total amount of protein to add is **1–10 μ g** per well.
8. Cover and incubate in the dark overnight (**15–18 hrs**) at RT with shaking.

Note: To resuspend the beads, shake vigorously at **900–1,100 rpm** for **30 sec**. Slowly ramp up to speed to avoid splashing. Then turn down to **300–450 rpm** for the specified incubation time.

Assay – Day 2 (Section 4 unless otherwise noted)

1. Start up/warm up the Bio-Plex system (may take up to **30 min**). (**Section 5**)
Meanwhile, bring buffers and diluents to RT.

2. Prime wash station or prepare handheld magnetic washer. (**Section 3**)
3. Calibrate the Bio-Plex system by following the prompts within Bio-Plex Manager™ software. (**Section 5**)
4. With **10 min** left in the overnight incubation, **vortex** 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

5. Wash the plate three times with **200 μl** wash buffer.
6. **Vortex** the diluted (1x) detection antibodies. Add **25 μl** to each well.
7. Cover and incubate in the dark for **30 min** at RT. Shake at **900–1,100 rpm** for **30 sec** and then turn down to **300–450 rpm** for the specified incubation time.
Meanwhile, prepare the Bio-Plex Manager software protocol. (**Section 5**)
8. With **10 min** left, **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	Detection Ab Diluent, μl	Total Volume, μl
96	60	5,940	6,000

9. Wash the plate three times with **200 μl** wash buffer.
10. **Vortex** the diluted (1x) SA-PE. Add **50 μl** to each well.
11. Cover and incubate in the dark for **10 min** at RT. Shake at **900–1,100 rpm** for **30 sec** and then turn down to **300–450 rpm** for the specified incubation time.
12. Wash the plate three times with **200 μl** wash buffer.

Bio-Plex Pro Assay Quick Guide 3

- Resuspend beads in **125 µl** bead resuspension buffer. Cover and shake at **1,100 rpm** for **30 sec**.
- Read the plate using the appropriate settings as shown below. (**Section 5**)

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

* Or similar Luminex-based system.

MAGPIX and MagPlex are trademarks 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.



CST antibodies exclusively developed and validated for Bio-Plex phosphoprotein and total target assays.



P10024930

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** 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000