

# Bio-Plex Pro™ RBM Metabolic and Hormone Assays

## Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro™ RBM Metabolic and Hormone Assays	10041818

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](http://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.

### A. Reagent Preparation

1. Reconstitute the following lyophilized reagents in dH<sub>2</sub>O before use, according to the table below.

Reagent	Volume, $\mu$ l
Standards Mix	150
Control 1	100
Control 2	100

Reagent	Volume, ml
Blocking Buffer	1.5
Standard Diluent	1.0
Detection antibiotics	4.8

- a. Allow vial to sit at room temperature for a minimum of **5 min**, not to exceed **30 min**.
  - b. Mix by **vortexing** at a medium setting.
2. Bring the 10x assay buffer to ambient/room temperature (RT).
    - a. Mix by inversion to ensure all salts are in solution.
    - b. Prepare 1x assay buffer — dilute **1 part** 10x assay buffer (60 ml) with **9 parts** of dH<sub>2</sub>O (540 ml).

### B. Dilution of Standard (1:3 Serial Dilution)

1. Label 9 polypropylene tubes **S1** through **S8** and **Blank**.
2. Transfer the reconstituted standard into the tube labeled **S1**.

## Bio-Plex Pro RBM Metabolic and Hormone Assays Quick Guide

3. Add the appropriate amount of the standard diluent into the labeled tubes according to the table below (this will be sufficient for duplicate standard curves and blanks).

Standard	Volume of Standard Diluent, $\mu$ l	Volume of Standard, $\mu$ l
S1	—	150 from reconstituted vial
S2	100	50 of S1
S3	100	50 of S2
S4	100	50 of S3
S5	100	50 of S4
S6	100	50 of S5
S7	100	50 of S6
S8	100	50 of S7
Blank	100	—

4. Prepare working standards (**S2–S8**) by serial dilution. Transfer the appropriate volume of standard into each of the labeled tubes with standard diluent, as outlined above.
5. **Vortex** each standard at a medium setting before proceeding with the next serial dilution. Change pipet tip at each dilution step.
6. The **Blank** tube consists of standard diluent alone.

### C. Sample Preparation

1. 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.
2. Prepare sample dilutions in **0.5 ml** or **1.0 ml** polypropylene tubes, as required for the assay.

## Bio-Plex Pro RBM Metabolic and Hormone Assays Quick Guide

3. Dilution scenarios provided below are sufficient to run each sample in duplicate.

Panel	Sample Dilution	Volume of Sample, $\mu\text{l}$		Volume of Sample Buffer, $\mu\text{l}$
Metabolic panel 1	1:5	20		80
Metabolic panel 2	1:5	20		80
Metabolic panel 3	1:500,000	(a) Prepare 1:50	5	245
		(b) Prepare 1:100	5 of (a)	495
		Prepare 1:100	5 of (b)	495
Metabolic panel 4	1:500	(c) Prepare 1:10	10	90
		Prepare 1:50	10 of (c)	490
Hormone panel 1	1:5	20		80

**Note:** Controls are ready to use after reconstitution. No further dilution is needed.

### D. Dispensing of Reagents

1. Add **10  $\mu\text{l}$**  of blocker to all wells of the plate.
2. Add **30  $\mu\text{l}$**  of the standard, control, sample, or blank to the appropriate well of the plate.
3. **Vortex** the capture beads at medium speed for **10–20 sec**. Add **10  $\mu\text{l}$**  of the beads to all wells of the plate.
4. Cover plate with plate seal and protect from light with aluminum foil. Incubate on shaker at **850  $\pm$  50 rpm** for **1 hr** at RT.
5. Wash the plate three times with **100  $\mu\text{l}$**  1x assay buffer.
6. **Vortex** the reconstituted detection antibodies at medium speed for **10–20 sec**. Add **40  $\mu\text{l}$**  to each well.
7. Cover and incubate at **850  $\pm$  50 rpm**, as in step 4, for **1 hr** at RT. **Do not aspirate after incubation.**
8. Prepare the required dilution of streptavidin-PE (SA-PE), as outlined in the following table.

**Note:** Volumes in the table are for an entire 96-well plate. Smaller volumes can be prepared, provided that the dilution ratios are maintained.

## Bio-Plex Pro RBM Metabolic and Hormone Assays Quick Guide

9. Add **20 µl** of diluted SA-PE to the required plate wells.

SA-PE Dilution	Volume of SA-PE, µl	Volume of 1x Assay Buffer, µl	Total Volume, µl
1:10	225	2,025	2,250

10. Cover and incubate at **850 ± 50 rpm**, as in step 4, for **30 min** at RT.

11. Wash the plate three times with **100 µl** 1x assay buffer.

12. After the final wash, resuspend the beads in **100 µl** 1x assay buffer. Cover plate, as in step 4, and shake the plate at **850 ± 50 rpm** for **30 sec**.

13. Remove the plate seal and **read plate** at low PMT (Bio-Plex® 200), standard PMT (Bio-Plex 3D), or default settings (Bio-Plex® MAGPIX™).

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by Luminex Corporation. MAGPIX is a trademark of Luminex Corporation.

Bio-Plex Pro RBM kits are manufactured by Myriad RBM. Myriad RBM is a trademark of Myriad RBM, Inc.



P10041817



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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 3065 7550 **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** 81 3 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** 1800 88 22 88  
**United Kingdom** 020 8328 2000