

# Bio-Plex Pro™ RBM IGF and IGFBP Assays

## Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro™ RBM IGF and IGFBP Assays	10042015

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, µl	Reagent	Volume, ml
Standards Mix	150	Blocking Buffer	1.5
Control 1	100	Standard Diluent	1.0
Control 2	100	Detection antibodies	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**.

## Bio-Plex Pro RBM IGF and IGFBP Assays Quick Guide

2. Transfer the reconstituted standard into the tube labeled **S1**.
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

**Note:** Most of the circulating IGF-1 and IGF-2 are in complex with IGFBPs. A sample extraction step is required prior to measuring IGFs in serum and plasma samples. Refer to **Section 5** of the instruction manual for guidance on extraction protocol. The IGF-1 and IGF-2 assays have been optimized to measure total IGF. For users measuring free IGF, please skip the sample extraction step and proceed to dilute the sample 1:5 with the provided sample diluent. The IGFBP assays do not require sample extraction.

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. Refer to instruction manual #10042015 for a detailed sample extraction protocol.
2. Prepare sample dilutions in **0.5 ml** or **1.0 ml** polypropylene tubes, as required for the assay.

## Bio-Plex Pro RBM IGF and IGFBP Assays Quick Guide

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

Panel	Sample Dilution	Volume of Sample	Volume of Extraction Buffer	Volume of Neutralization Buffer	Volume of Sample Dilution Buffer 2
IGF	1:30	20 $\mu$ l	180 $\mu$ l	Mix 50 $\mu$ l of extracted sample with 50 $\mu$ l of neutralization buffer	Mix the neutralized sample with 50 $\mu$ l of sample dilution buffer 2
IGFBP	1:20	10 $\mu$ l	N/A	N/A	190 $\mu$ l

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

### D. Dispensing of Reagents

**Note:** This protocol is optimized with a single wash step after the SA-PE incubation. Do not wash the plate after the sample incubation and the detection antibody incubation steps.

1. Add **10  $\mu$ l** of blocker to all wells of the plate.
2. Add **30  $\mu$ 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  $\mu$ 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. **Do not aspirate after incubation.**
5. **Vortex** the reconstituted detection antibodies at medium speed for **10–20 sec**. Add **40  $\mu$ l** to each well.
6. Cover and incubate at **850  $\pm$  50 rpm**, as in step 4, for **1 hr** at RT. **Do not aspirate after incubation.**
7. 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 IGF and IGFBP Assays Quick Guide

8. 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

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

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

11. 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**.

12. 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.



**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