

Bio-Plex Pro[™] RBM Metabolic and Hormone Assays

Instruction Manual

For technical support, call your local Bio-Rad office, or in the U.S., call 1-800-424-6723.
For research use only. Not for diagnostic procedures.



BIO-RAD

Table of Contents

Introduction	1
Principle	2
Kit Contents and Storage	4
Recommended Materials	5
Assay Workflow	6
Important Considerations	7
Detailed Instructions	7
1. Plan Plate Layout	8
2. Prepare Instrument	9
3. Prepare Wash Method	10
4. Prepare Reagents	11
5. Prepare Samples	13
6. Run the Assay	16
7. Read Plate	18
Troubleshooting Guide	24
Plate Layout Template	29
Safety Considerations	30
Legal Notices	30
Ordering Information	31

Introduction

Approximately 65% of individuals with diabetes die from heart disease and stroke. Individuals with type 2 diabetes also have an increased rate of high blood pressure, lipid problems, and obesity, which contribute to their high rates of cardiovascular disease (CVD). Given the complexity of these conditions, it is not surprising that assessment of a single biomarker alone has not yielded more than an incremental advantage over traditional measures of assessing risk. In diabetes, further complexity is added by the common coincidence of metabolic syndrome, which includes insulin resistance, dyslipidemia, hypertension, and systemic inflammation, all of which contribute to the pathogenesis of CVD (Iqbal et al. 2012). As such, researchers studying CVD and/or diabetes need to monitor multiple biomarkers for cardiovascular risk factors along with biomarkers directly linked to the metabolic disorders mentioned above. These markers include gut and pituitary hormones, adipokines, growth factors, and cardiovascular and acute phase-related proteins. Since these hormones and proteins frequently act in concert, it is advantageous to have the capacity to measure a number of different biomarkers simultaneously in a single blood sample.

Bio-Plex Pro™ RBM Metabolic and Hormone Assays

The Bio-Plex Pro RBM metabolic and hormone assays include a wide range of proteins implicated in the pathophysiology of diabetes, obesity, and metabolic syndrome, as well as biomarkers for associated conditions such as cardiovascular disease and inflammation.

For researchers working with limited sample volume, the capacity to multiplex provides an effective option over the traditional ELISA method. The use of magnetic (MagPlex) beads allows researchers to automate wash steps on a Bio-Plex Pro (or similar) wash station.

References

Iqbal N et al. (2012). Cardiac biomarkers: New tools for heart failure management. *Cardiovas Diagn Ther* 2, 147–164.

Principle

Technology

The Bio-Plex® multiplex system is built upon the three core elements of xMAP technology:

- Fluorescently dyed magnetic microspheres (also called beads), each with a distinct color code or spectral address to permit discrimination of individual tests within a multiplex suspension. This allows simultaneous detection of up to 500 different molecules in a single well of a 96-well microplate on the Bio-Plex® 3D system, up to 100 different molecules on the Bio-Plex® 200 system, and up to 50 different molecules on the Bio-Plex® MAGPIX™ system
- A dedicated plate reader. The Bio-Plex 200 and Bio-Plex 3D systems are flow cytometry–based instruments with two lasers and associated optics to measure the different molecules bound to the surface of the beads. In the Bio-Plex MAGPIX system, the sample is injected into a chamber where the beads are imaged using LED and CCD technology
- A high-speed digital signal processor that efficiently manages the fluorescence data

Assay Format

The Bio-Plex Pro™ RBM metabolic and hormone assays are essentially immunoassays formatted on magnetic beads. The assay principle is similar to that of a sandwich ELISA (Figure 1). Capture antibodies directed against the desired biomarker are covalently coupled to the beads. Coupled beads react with the sample containing the biomarker of interest. After a series of washes to remove unbound protein, a biotinylated detection antibody is added to create a sandwich complex. The final detection complex is formed with the addition of streptavidin-phycoerythrin (SA-PE) conjugate. Phycoerythrin serves as a fluorescent indicator, or reporter. The use of magnetic (MagPlex®) beads allows researchers to automate wash steps on a Bio-Plex Pro (or similar) wash station. Magnetic separation offers greater convenience, productivity, and reproducibility compared to vacuum filtration.

Note: The cortisol assay in Metabolic Panel 1 is a competitive immunoassay (schematic not shown).

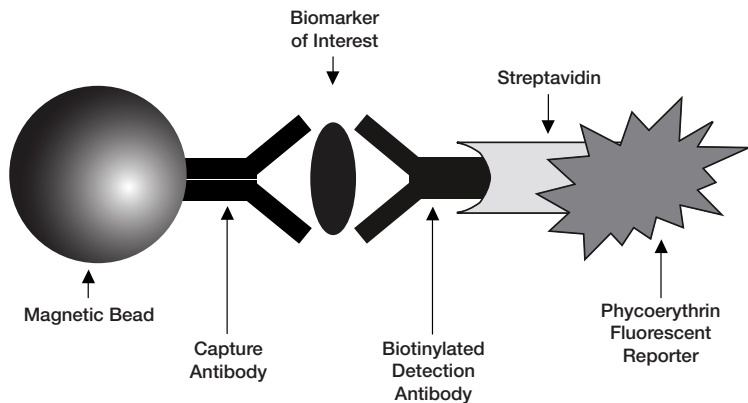


Fig. 1. Bio-Plex sandwich immunoassay.

Data Acquisition and Analysis

Data from the reactions are acquired using a Bio-Plex system or similar Luminex-based reader. When a multiplex assay suspension is drawn into the Bio-Plex 200 reader, for example, a red (635 nm) laser illuminates the fluorescent dyes within each bead to provide bead classification and thus assay identification. At the same time, a green (532 nm) laser excites PE to generate a reporter signal, which is detected by a photomultiplier tube (PMT). A high-speed digital processor manages data output and Bio-Plex Manager™ software presents data as median fluorescence intensity (MFI) as well as concentration. The concentration of analyte bound to each bead is proportional to the MFI of reporter signal.

Kit Contents and Storage

The Bio-Plex Pro™ RBM metabolic and hormone assays are available in a convenient kit format that includes assay, reagent, and diluent components in a single box (Table 1). All other recommended materials are listed in Table 2.

Table 1. Contents of 1 x 96-well kits.

Component	Quantity	Volume	Volume after Reconstitution or Dilution
Capture beads (1x)	1 tube	1.4 ml	
Detection antibodies	1 vial	Lyophilized	4.8 ml
Standards mix	1 vial	Lyophilized	150 µl
Control 1 (high)	1 vial	Lyophilized	100 µl
Control 2 (low)	1 vial	Lyophilized	100 µl
Blocking buffer	1 vial	Lyophilized	1.5 ml
Standard diluent	1 vial	Lyophilized	1.0 ml
Sample dilution buffer	1 bottle	100 ml	
Assay buffer (10x)	1 bottle	60 ml	600 ml
Streptavidin-PE (10x)	1 tube	250 µl	2.5 ml
Assay plate (96-well flat bottom)	1 plate		
Plate seals	1 pack of 3		
Assay quick guide	1 sheet		
Product data sheet	1 sheet		

Storage and Stability

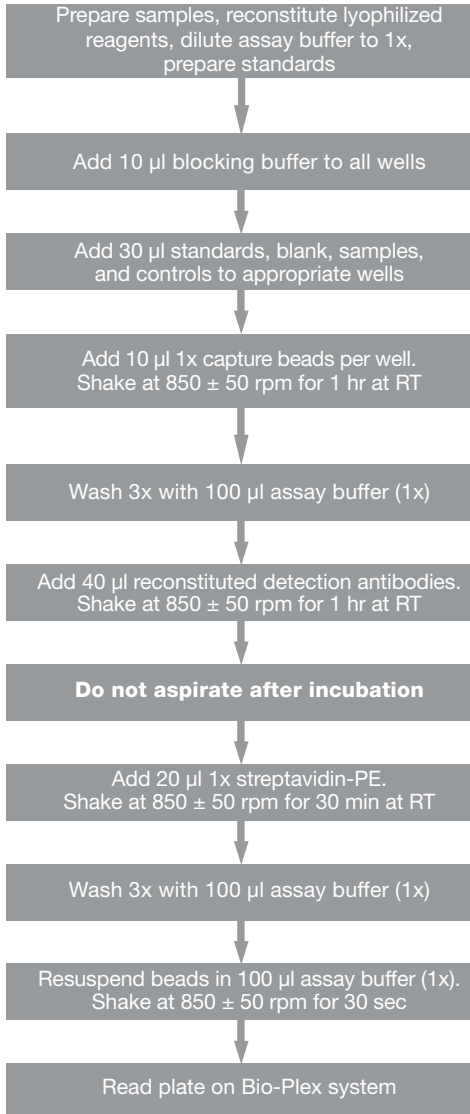
Kit contents should be stored at 2–8°C and never frozen. Coupled magnetic capture beads and streptavidin-PE should be stored in the dark. All components are guaranteed for a minimum of six months from the date of purchase when stored as specified.

Table 2. Recommended materials.

Item	Ordering Information
Bio-Plex® 200 system or Luminex system with HTF	Bio-Rad catalog #171-000205
Bio-Plex validation kit Note: Run the validation kit monthly to ensure optimal performance of fluidics and optics systems	Bio-Rad catalog #171-203001
Bio-Plex calibration kit Note: Run the calibration kit daily to standardize fluorescence signal	Bio-Rad catalog #171-203060
Bio-Plex Pro wash station For use with magnetic bead-based assays only	Bio-Rad catalog #300-34376
Bio-Plex Pro II wash station For use with both polystyrene (nonmagnetic) and magnetic bead-based assays	Bio-Rad catalog #300-34377
Bio-Plex handheld magnetic washer For use with magnetic bead-based assays only	Bio-Rad catalog #170-20100
Bio-Plex Pro flat bottom plates (40 x 96-well) For magnetic separation on the Bio-Plex Pro wash station	Bio-Rad catalog #171-025001
Titertube® micro test tubes For preparing replicate standards, samples, and controls prior to loading the plate	Bio-Rad catalog #223-9390
Microtiter plate shaker IKA MTS 2/4 shaker for 2 or 4 microplates or Barnstead/Lab-Line Model 4625 plate shaker (or equivalent capable of 300–1,100 rpm)	IKA catalog #320-8000 WWR catalog #57019-600
Aurum™ vacuum manifold For vacuum filtration	Bio-Rad catalog #732-6470
BR-2000 vortexer	Bio-Rad catalog #166-0610
Reagent reservoirs, 25 ml For capture beads and detection antibodies	VistaLab catalog #3054-1002 or VistaLab catalog #3054-1004
Reagent reservoir, 50 ml (for reagents and buffers)	VistaLab catalog #3054-1006
Pall Life Science Acrodisc: 25 mm PF syringe filter (0.8/0.2 µm Supor membrane)	Pall Life Sciences catalog #4187
Filter plate, 1 x 96 with clear plastic lid and tray	Bio-Rad catalog #171-304502

Other: 15 ml polypropylene tubes for reagent dilutions, calibrated pipets, pipet tips, sterile distilled water, aluminum foil, absorbent paper towels, 1.5 or 2 ml microcentrifuge tubes, and standard flat bottom microplate (for calibrating vacuum manifold).

Assay Workflow



Important Considerations

Instruments and Software

The assays described in this manual are compatible with all currently available Luminex-based life science research instruments. Assays can be read and analyzed with either Bio-Plex Manager™ software or Luminex xPONENT software.

Assay Procedures

Pay close attention to vortexing, shaking, and incubation times and to Bio-Plex® reader PMT (RP1) setting, as these have been optimized specifically for each assay panel.

Assay Quick Guide

Each assay kit comes complete with a printed Bio-Plex Pro RBM Metabolic and Hormone Assays Quick Guide (bulletin #10041817), which can be used to set up and prepare a full 1 x 96-well assay plate. Users can also download a copy at www.bio-rad.com/bio-plex.

Bead Regions and Multiplexing Compatibility

- Bead regions for all analytes are listed in the Read Plate section
- Do not mix analytes between different panels, or with other Bio-Plex assay panels or reagents

Detailed Instructions

The following pages provide detailed instructions for each step of the assay procedure, including preparation, running the assay, and reading the plate with Bio-Plex Manager™ and Luminex xPONENT software.

1. Plan Plate Layout

Determine the total number of wells in the experiment using the Plate Layout Template on page 29 or the Plate Formatting tab in Bio-Plex Manager™ software. A suggested plate layout is shown in Figure 2, with all conditions in duplicate.

1. Assign standards to columns 1 and 2, with the highest concentration in row A and the lowest concentration in row H.
2. Assign the blank to wells A3 and A4. The blank should consist of standard diluent alone and be processed in the same manner as sample and standard wells. Bio-Plex Manager software automatically subtracts the assay blank (B) MFI value from all other assay wells.
3. User-specific controls, as well as the quality controls supplied in the kits, are assigned to wells in columns 3 and 4.
4. The remainder of the plate is available for samples.

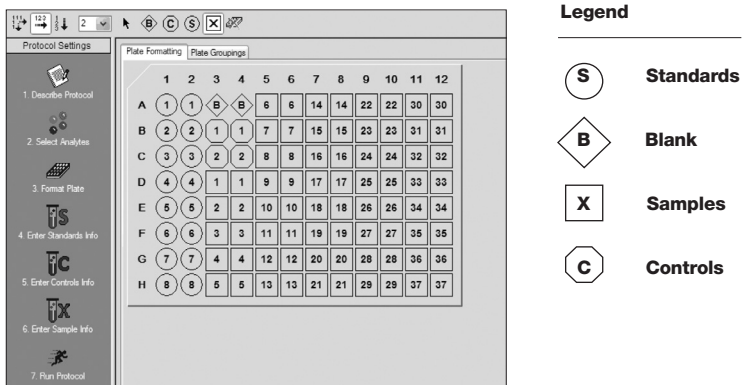


Fig. 2. Suggested plate layout. For detailed instructions on plate formatting in Bio-Plex Manager, see section Read Plate.

2. Prepare Instrument



These directions are specific for the Bio-Plex® 100/200 reader. To prepare either a Bio-Plex 3D or Bio-Plex® MAGPIX™ reader, consult their respective user manuals.

Start up and calibrate the Bio-Plex system with Bio-Plex Manager™ software prior to setting up the assay. The calibration kit should be run daily or before each use of the instrument to standardize the fluorescent signal. For instructions on using other xMAP system software packages, contact Bio-Rad Technical Support.


Note: While the instrument is warming up, bring the 10x assay buffer and sample dilution buffer to room temperature. Keep other items on ice until needed. Also, begin to thaw frozen samples.

The validation kit should be run monthly to ensure performance of fluidics and optics systems. Refer to either the software manual or online Help for directions on how to conduct validation.


Start Up System (Bio-Plex 100, 200, or Similar)

1. Empty the waste bottle and fill the sheath fluid bottle before starting if high throughput fluidics (HTF) are not present. This will prevent fluidic system backup and potential data loss.
2. Turn on the reader, XY platform, and HTF (if included). Allow the system to warm up for 30 min (if not already done).
3. Select Start up  and follow the instructions. If the system is idle for 4 hr without acquiring data, the lasers will automatically turn off. To reset the 4-hr countdown, select Warm up  and wait for the lasers/optics to reach operational temperature.

Calibrate System

1. Select Calibrate  and confirm that the default values for CAL1 and CAL2 are the same as the values printed on the bottle of Bio-Plex calibration beads. Use the Bio-Plex system low RP1 target value.

2. Select OK and follow the software prompts for step-by-step instructions for CAL1 and CAL2 calibration.

Note: In Bio-Plex Manager version 6.1 and higher, startup, warm up, and calibration can be performed together by selecting the “Start up and calibrate” icon. 

3. Prepare Wash Method

Bio-Plex Pro™ assays are compatible with both magnetic separation and vacuum filtration methods. However, for best results, we recommend performing the assays in a flat bottom plate with magnetic separation.

Table 3. Summary of compatible wash stations and plate types.

Wash Method	Wash Station	Assay Plate
Magnetic separation	Bio-Plex Pro Bio-Plex Pro II (use MAG programs) Bio-Plex® handheld magnetic washer	Flat bottom plate
Vacuum filtration	Bio-Plex Pro II (use VAC programs) Vacuum manifold (manual)	Filter plate

Setting Up the Bio-Plex Pro or Bio-Plex Pro II Wash Station

The wash station should be primed before use. For more information, refer to the Bio-Plex Pro Wash Stations Quick Guide (bulletin #5826).

1. Install the appropriate plate carrier on the wash station.
2. Use the prime procedure to prime channel 1 with 1x assay buffer

Setting Up the Bio-Plex Handheld Magnetic Washer

Place an empty flat bottom plate on the magnetic washer by sliding it under the retaining clips. Push the clips inward to secure the plate. Make sure the plate is held securely. If needed, the clips can be adjusted for height and tension. For detailed instructions, refer to the user guide (bulletin #10023087).

Setting Up a Vacuum Manifold

Calibrate the vacuum manifold by placing a standard 96-well flat bottom plate on the unit and adjusting the pressure to -1 to -3" Hg. In general, 100 μ l liquid should take 3-4 sec to clear the well. For more detailed instructions, refer to bulletin #10005042.

4. Prepare Reagents

1. Reconstitute the following lyophilized reagents in dH₂O before use according to the table below.

Table 4. Reagent volume.

Reagent	dH₂O Volume
Standards mix	150 μ l
Control 1	100 μ l
Control 2	100 μ l
Blocking buffer	1.5 ml
Standard diluent	1.0 ml
Detection antibodies	4.8 ml

- 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 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 dH₂O** (540 ml).

Dilution of Standard (1:3 Serial Dilution)

This preparation provides sufficient volume to run duplicate standard dilution curves. Ensure that each new standard is mixed well by vortexing before proceeding to the next dilution. Change tips between each dilution.

Note: The product data sheet in each kit lists the most concentrated point on the standard curve (S1). Enter the values and units into Bio-Plex Manager™ software as instructed in the Read Plate section.

1. Label 9 polypropylene tubes **S1** through **S8** and **Blank**.
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 Figure 3 (this will be sufficient for duplicate standard curves and blanks).
4. Prepare working standards (**S2–S8**) by 1:3 (threefold) serial dilution. Transfer the appropriate volume of standard into each of the labeled tubes containing standard diluent as outlined above.
5. **Vortex** each standard at a medium setting for **5 sec** before proceeding with the next serial dilution. Change pipet tip at each dilution step.
6. The **Blank** tube consists of standard diluent alone.

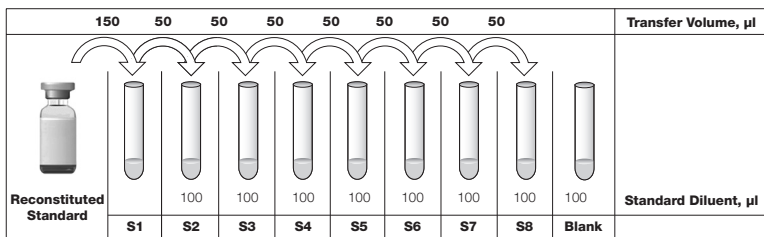


Fig. 3. Preparing a threefold dilution series with a single reconstituted standard.

5. Prepare Samples

The kit has sufficient reagents to run 37 samples in duplicate for a full plate analysis. General guidelines for preparing samples are provided below. For more information, contact Bio-Rad Technical Support.

Serum and Plasma

Serum, EDTA plasma, or citrate plasma are the preferred sample types for these assays. Heparin-treated plasma, while compatible with these assays, may absorb certain soluble proteins of interest.

- Avoid using hemolyzed samples, as this may lead to false results.
- Freezing samples at -70°C immediately after preparation and keeping samples frozen until use should provide adequate protection from degradation. Avoid multiple freeze and thaw cycles.
- Once thawed, keep samples on ice.
- Do not freeze diluted samples

Protease Inhibitors

In general, these biomarkers are detectable in serum or plasma without using protease inhibitors. Users may choose to add protease inhibitors as a precautionary measure if samples will not be frozen immediately after collection. Alternatively, blood sample collection tubes containing inhibitors may be obtained from Becton-Dickinson, www.bd.com.

Note: Protease inhibitors are not recommended for use with serum samples. Compatibility of a given inhibitor cocktail with the assays must be validated by the end user.

1. Draw whole blood into collection tubes containing anticoagulant. Invert tubes several times to mix.
2. For serum, allow blood to clot at room temperature for 30–45 min. For plasma, proceed directly to the centrifugation steps.
3. Perform centrifugation at 1,000 x g for 15 min at 4°C and transfer the serum or plasma to a clean polypropylene tube.
4. To completely remove platelets and precipitates, centrifuge again at 10,000 x g for 10 min at 4°C.
5. Prepare sample dilutions in 0.5 or 1.0 ml polypropylene tubes as required for the assay.
6. Dilution scenarios provided below are sufficient to run each sample in duplicate.

Cell culture media

1. Collect cell culture supernatants and centrifuge at 1,000 x g for 15 min at 4°C. For cell lines cultured in serum-free media, collect samples and add BSA as a carrier protein to a final concentration of at least 0.5%. This is done to stabilize protein analytes and to prevent adsorption to labware.
2. Transfer to a clean polypropylene tube. If cellular debris or precipitates are present, centrifuge again at 10,000 x g for 10 min at 4°C.
3. If high levels of analyte are expected, samples can be diluted in culture media. Optimal dilution factor must be determined by the end user. Supplement serum-free media with at least 0.5% BSA final.
4. Assays samples immediately or aliquot and store at -70°C.

Note: Compatibility of a given media type with the assays must be validated by the end user.

Table 5. Serum and plasma sample dilution guidelines.

Panel	Sample Dilution	Volume of Sample, μ l	Volume of Sample Buffer, μ 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 (b) Prepare 1:100 5 of (a) Prepare 1:100 5 of (b)	245 495 495
Metabolic panel 4	1:500	(c) Prepare 1:10 10 Prepare 1:50 10 of (c)	90 490
Hormone panel 1	1:5	20	80

Note: Sample dilution factor for culture media samples must be optimized by the end user.

6. Run the Assay

Considerations

- Bring all assay components and samples to room temperature before use
- Use calibrated pipets and pipet carefully, avoiding bubbles. Use new pipet tips for every volume transfer
- Pay close attention to vortexing, shaking, and incubation instructions. Deviation from protocol may result in low assay signal and assay variability
- Assay incubations are carried out in the dark on a shaker at **850 ± 50 rpm**. Cover the plate with a plate seal and protect from light with aluminum foil

Table 6. Summary of wash options and protocols. After each assay step, select the appropriate Bio-Plex Pro wash station program or perform the appropriate manual wash step as summarized below.

Assay Step	Bio-Plex Pro or Pro II Wash Station	Bio-Plex Pro II Wash Station	Handheld Magnet or Vacuum Manifold
	Magnetic Program	Vacuum Program	Manual Wash Steps
Sample incubation	MAG x3	VAC x3	3 x 100 µl
SA-PE incubation			

Considerations When Using a Vacuum Manifold

- After each incubation, place the filter plate on a calibrated vacuum apparatus and remove the liquid by vacuum filtration
- To wash, add 100 µl wash buffer to each well and remove the liquid as before. Ensure that all wells are exposed to the vacuum
- Thoroughly blot the bottom of the filter plate with a clean paper towel between each vacuum step to prevent cross contamination
- Place the assay plate on the plastic plate holder/tray as needed
- Before each incubation, gently cover the plate with a new plate seal. Avoid pressing down on the wells to prevent leaking from the bottom

Assay Protocol: Dispensing of Reagents

1. Add **10 μ l** blocker to all wells of the plate.
2. Add **30 μ l** the standard, control, or sample to the appropriate well of the plate.
3. **Vortex** the capture beads at medium speed for **10–20 sec**. Add **10 μ 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 μ l** 1x assay buffer.
6. **Vortex** the reconstituted detection antibodies at medium speed for **10–20 sec**. Add **40 μ 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 Table 7.

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

Table 7. SA-PE dilution.

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

9. Add **20 μ l** diluted SA-PE to the required plate wells.
10. Cover and incubate at **850 \pm 50 rpm**, as in Step 4, for **30 min** at RT.
11. Wash the plate three times with **100 μ l** 1x assay buffer.
12. After the final wash, resuspend the beads in **100 μ l** assay buffer. Cover plate as in Step 4 and shake the plate at **850 \pm 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[™]).

7. Read Plate


Bio-Plex Manager software is recommended for all Bio-Plex Pro assay data acquisition and analysis. Instructions for Luminex xPONENT software are also included. For instructions on using other xMAP system software packages, contact Bio-Rad Technical Support or your regional Bio-Rad field applications specialist.

Prepare Protocol in Bio-Plex Manager Software Version 6.0 and Higher

The protocol should be prepared in advance so that the plate is read as soon as the experiment is complete.

A protocol file specifies the analytes used in the reading, the plate wells to be read, sample information, the values of standards and controls, and instrument settings. Protocols may be obtained from within Bio-Plex Manager software version 6.1 or created from the File menu.

To create a new protocol, select **File**, then **New** from the main menu. Locate and follow the steps under **Protocol Settings**.

1. Click **Describe Protocol** and enter information about the assay (optional).
2. Click **Select Analytes** and create a new panel.
 - a. Click **Add Panel**  in the Select Analytes toolbar. Enter a new panel name. Select **Bio-Plex Pro Assay Magnetic** from the assay dropdown menu. If using Bio-Plex Manager version 5.0 or lower, select **MagPlex** from the assay dropdown menu.
 - b. Click **Add**. Enter the bead region number and name for the first analyte. Click **Add Continue** to repeat for each analyte in the assay. Refer to the bead regions in parentheses () listed on the peel-off label provided with the standards.

For reference, bead regions for the individual assays are listed in Table 8.

- c. Click **Add** when the last analyte has been added and click **OK** to save the new panel.

- d. Highlight analytes from the **Available** list (left) and move to the **Selected** list (right) using the **Add** button. To move all analytes at once, simply click **Add All**.
- e. If some of the analytes need to be removed from the Selected list, highlight them and select **Remove**. If desired, it is possible to rename the panel by clicking on **Rename Panel** and entering a new panel name.

Table 8. Bead regions for Bio-Plex Pro RBM metabolic and hormone assays.

Metabolic Panel 1	Bead Region	Metabolic Panel 2	Bead Region
C-Peptide	45	FGF-21	63
Cortisol	12	FGF-23	47
Pancreatic polypeptide	52	Galectin-3	46
Proinsulin	67	sLeptin R	27
Peptide YY	55	Omentin-1	20
		Pentraxin-3	44
		PON-1	53
		Vaspin	34
Metabolic Panel 3	Bead Region	Metabolic Panel 4	Bead Region
AAT	26	ANGPTL3	66
AGP-1	57	Chemerin	42
Hemopexin	35	DPP4	25
RBP-4	43	Protein S	62
Transferrin	61	SEPP	36
Transthyretin	13		
VDBP	28		

Hormone Panel 1	Bead Region
FSH	18
GH	38
LH	15
Prolactin	19
TSH	22

Note: The cortisol assay is a competitive immunoassay.





3. Click **Format Plate** and format the plate according to the plate layout template (located on page 29) created for the assay. To modify the plate layout, follow the steps below.
 - a. Select the **Plate Formatting** tab.
 - b. Select the standards icon  and drag the cursor over all the wells that contain standards. Repeat this process for blanks , controls , and samples .
4. Click **Enter Standards Info** in the Protocol Settings bar.
 - a. Enter the highest concentration of each analyte in the top row (labeled S1) of the table. S1 concentration information is listed in the product data sheet.
 - b. Enter a dilution factor of 3 and click **Calculate**. The concentrations for each standard point will be populated for all analytes in the table.
 - c. Optional: enter the lot number of the vial of standards into the **Standard Lot** box and click **Save**.
5. Click **Enter Controls Info**.
 - a. For user-specified controls, select an analyte from the dropdown menu, then enter a description and concentration. Repeat for each additional analyte in the assay.
 - b. For the kit controls supplied, format the appropriate wells as controls and enter descriptions, but leave the concentrations blank. Alternatively, the controls can be formatted as samples with clear descriptions such as “quality control.” In any case, the expected control ranges provided on the product data sheet are not entered into Bio-Plex Manager software version 6.1 and earlier.
6. Click **Enter Sample Info** — enter sample information and the appropriate dilution factor.
7. Click **Run Protocol** — confirm that the assay settings follow Table 9.

Table 9. Read the plate using the appropriate instrument settings.



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		

* A similar Luminex-based system may be used.

- a. Confirm that data acquisition is set to **50 beads per region**.
- b. In Bio-Plex Manager software prior to 6.1, go to Advanced Settings, confirm that the bead map is set to **100 region**, the sample size is set to **50 µl**, and the DD gates are set to **5,000 (Low)** and **25,000 (High)**. In Bio-Plex Manager software versions 4.0, 4.1, 4.1.1, and 5.0, check **Override Gates** and set the DD gate values as indicated.

Select **Start**, name and save the .rbx file, and begin data acquisition. The Run Protocol pop-up screen will appear. Click **Eject/Retract** to eject the plate carrier.

Acquire Data

1. Shake the assay plate at **850 ± 50 rpm** for **30 sec**, and visually inspect the plate to ensure that the assay wells are filled with buffer. Slowly remove the sealing tape and any plate cover before placing the plate on the plate carrier.
2. Click **Run Protocol** — on the pop-up screen, select **Load Plate** and click **OK** to start acquiring data.
3. Use the **Wash Between Plates**  command after every plate run to reduce the possibility of clogging the instrument.
4. If acquiring data from more than one plate, empty the waste bottle and refill the sheath bottle after each plate (if HTF are not present). Select **Wash Between Plates** and follow the instructions. Then repeat the Prepare Protocol and Acquire Data instructions.
5. When data acquisition is complete, select **Shut Down**  and follow the instructions.

Data Analysis

Quality Controls

If the quality controls were run in the assay plate, open the results (.rbx) file, click **Report Table**, and locate the control wells. Compare the observed concentrations against the lot-specific control ranges in the product data sheet.

Note: Expected control ranges are provided for reference and should be used as general guidelines. Actual results may vary for some operators. If the controls do not fall within the expected ranges, please refer to the troubleshooting section for possible causes and solutions.

Removing Outliers

Outliers are identified as standard data points that do not meet accuracy or precision requirements and should be considered invalid when performing curve fitting. As such, they should be removed to generate a more realistic and accurate standard curve. This may result in an extended assay working range and allow quantitation of samples that might otherwise be considered out of range (OOR).

In Bio-Plex Manager software version 6.0 and higher, outliers can be automatically removed by selecting the **Optimize** button in the Standard Curve window. In earlier versions of the software, outliers can also be manually selected in the Report Table. Visit online Help to learn more about the standard curve optimizer feature and how outliers are determined.

Previous Versions of Bio-Plex Manager Software

For instructions on using previous versions of Bio-Plex Manager software, please contact Bio-Rad Technical Support.

Luminex xPONENT Software

Luminex xPONENT software may be used to analyze Bio-Plex assays. Although guidelines are provided here, consult the xPONENT software manual for more details. Perform a system initialization with Luminex's calibration and performance verification kit, as directed by Luminex. Select **Batches** to set up the protocol and follow the information under Settings.

Note: The instrument settings described below apply to Luminex 100/200 and FLEXMAP 3D or Bio-Plex 3D instruments. For the Bio-Plex MAGPIX reader, use the default instrument settings.

1. Select MagPlex as the bead type for magnetic beads, which automatically sets the DD gates.
2. Volume = 50 μ l.
3. Low PMT (Standard PMT).
4. Plate name: 96-well plate.
5. Analysis type: Quantitative; 5PL Curve Fit.
6. Number of standards: 8.

Select **Analytes** to set up the panel.

1. Enter "ng/ml" in the Units field.
2. Enter 50 in the Count field.
3. Select the bead region and enter the analyte name.
4. Click **Apply all** for Units and Count.

Select **Stds and Ctrl.**

1. Enter standard concentrations, lot number, dilution factor, and other information, as applicable.

After the assay is complete, select **Results**, then select **Saved Batches**.

Troubleshooting Guide

This troubleshooting guide addresses problems that may be encountered with Bio-Plex Pro™ RBM assays. If you experience any of the problems listed below, review the possible causes and solutions provided. Poor assay performance may also be due to the Bio-Plex® suspension array reader. To eliminate this possibility, use the validation kit to validate all the key functions of the array reader and to assist in determining whether or not the array reader is functioning properly.

Possible Causes

High Inter-Assay CV

Standards were not reconstituted consistently between assays

Reconstituted standards and diluted samples were not stored properly

Bottom of filter plate not dry

Possible Solutions

Incubate the reconstituted standards for for the recommended time period. Always be consistent with the incubation time and temperature.

Diluted samples should be prepared on ice as instructed. Prior to plating, the reconstituted standards and diluted samples should be equilibrated to room temperature.

Dry the bottom of the filter plate with absorbent paper towel (preferably lint-free) to prevent cross-well contamination.

Possible Causes

High Intra-Assay CV

Improper pipetting technique

Reagents and assay components not equilibrated to room temperature prior to pipetting

Contamination with buffer during wash steps

Slow pipetting of samples and reagents across the plate

Low Bead Count

Beads clumped in multiplex bead stock tube

Possible Solutions

Pipet carefully when adding standards, samples, detection antibodies, and streptavidin-PE, especially when using a multichannel pipet. Use a calibrated pipet. Change pipet tip after every volume transfer.

All reagents and assay components should be equilibrated to room temperature prior to pipetting.

During the wash steps, be careful not to splash buffer from one well to another. Be sure that the wells are filtered completely and that no residual volume remains. Ensure that the microplate shaker setting is not too high. Reduce the microplate shaker speed to minimize splashing.

Sample pipetting across the entire plate should take less than 4 min. Reagent pipetting across the entire plate should take less than 1 min.

Vortex for 30 sec at medium speed before aliquoting beads.

Possible Causes

Low Bead Count

Vacuum on for too long when aspirating buffer from wells

Reader is clogged

Low Signal or Poor Sensitivity

Standards reconstituted incorrectly

Detection antibody or streptavidin-PE prepared incorrectly

High Background Signal

Incorrect buffer was used (for example, assay buffer used to dilute standards)

Accidentally spiked blank wells

Detection antibodies or streptavidin-PE incubated too long

Possible Solutions

Do not apply vacuum to the filter plate for longer than 10 sec after the buffer is completely drained from each well.

Refer to the troubleshooting guide in the Bio-Plex system hardware instruction manual (bulletin #10005042).

Follow the standard preparation instructions carefully.

Check your calculations and be careful to add the correct volumes.

Use standard diluent to dilute standards.

Do not add any antigens to the blank wells.

Follow the procedure incubation time precisely.

Possible Causes

Poor Recovery

Expired Bio-Plex reagents were used

Incorrect amounts of components were added

Microplate shaker set to an incorrect speed

High end saturation of the standard curve

Controls do not fall within expected ranges

Improper pipetting technique

Possible Solutions

Check that reagents have not expired. Use new or nonexpired components.

Check your calculations and be careful to add the correct volumes.

Check the microplate shaker speed and use the recommended setting. Setting the speed too high may cause splashing and contamination. Use the recommended plate shaker.

Make sure that correct shaker speed and incubation times are used. Remove S1 for data analysis if needed.

Make sure that the vial of controls is reconstituted at the same time as standards and in the correct diluent. Incubate for times indicated.

Pipet carefully when adding standards, samples, detection antibodies, and streptavidin-PE, especially when using a multichannel pipet. Use a calibrated pipet. Change pipet tip after every volume transfer.

Possible Causes

Impact of Sample Matrix

Negative MFI values in samples

Possible Solutions

If samples contain little or no analyte, negative values observed may be due to statistical variation. If assay drift is suspected, retest the samples by positioning them next to the standards. If contamination of standards is suspected, check the standard replicate value and be careful when adding samples to the wells. Matrix effects could also produce negative sample values.

Bio-Plex Manager™ software automatically subtracts the blank (B) MFI value from all other assay wells. While this has no impact on observed concentrations of samples within the assay working range, it may result in a negative MFI value if the blank's MFI value is greater than either the standard or the sample value. If this is undesirable, then reformat the blank wells as sample (X) or control (C) in the protocol or results file.

Plate Layout Template

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H

Safety Considerations

Eye protection and gloves are recommended when using these products. Consult the MSDS for additional information. The Bio-Plex Pro™ assays contain components of animal origin. This material should be handled as if capable of transmitting infectious agents. Use universal precautions. These components should be handled at Biosafety Level 2 containment (U.S. government publication: Biosafety in Microbiological and Biomedical Laboratories (CDC 1999)).

Legal Notices

Bio-Plex Pro RBM kits are manufactured by Myriad RBM.

Acrodisc and Supor are trademarks of Pall Corporation. MAGPIX, MagPlex, xMAP, xPONENT, FLEXMAP 3D, and Luminex are trademarks of Luminex Corporation. Myriad RBM is a trademark of Myriad RBM, Inc.

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

Ordering Information

Detailed ordering information can be found at www.bio-rad.com/bio-plex.

Catalog #	Premixed 1 x 96-Well All-In-One Multiplex Kit
	Includes premixed magnetic capture beads, premixed detection antibodies, standards mix, 2-level controls, blocking buffer, standard diluent, sample dilution buffer, 10x assay buffer, 10x streptavidin-PE, 96-well flat bottom plate, plate seals, and instructions.
171-AMR1CK	Bio-Plex Pro RBM Human Metabolic Panel 1 , 1 x 96
171-AMR2CK	Bio-Plex Pro RBM Human Metabolic Panel 2 , 1 x 96
171-AMR3CK	Bio-Plex Pro RBM Human Metabolic Panel 3 , 1 x 96
171-AMR4CK	Bio-Plex Pro RBM Human Metabolic Panel 4 , 1 x 96
171-AHR1CK	Bio-Plex Pro RBM Human Hormone Panel 1 , 1 x 96
Catalog #	Wash Stations and Accessories
300-34376	Bio-Plex Pro Wash Station , includes magnetic plate carrier, waste bottle, 2 buffer bottles
300-34377	Bio-Plex Pro II Wash Station , includes magnetic plate carrier, vacuum manifold plate carrier, waste bottle, 2 buffer bottles
171-020100	Bio-Plex Handheld Magnetic Washer , includes magnetic washer and adjustment hex tools for use in manual wash steps for all Bio-Plex magnetic assays
171-025001	Bio-Plex Pro Flat Bottom Plates , 40 x 96-well plates
171-304500	Bio-Plex Wash Buffer , 1.5 L
171-304502	Filter plate , pkg of 1, 96-well plate with clear plastic lid and tray, for Bio-Plex assays using the vacuum wash method, sealing tape not included
Catalog #	Software
171-001510	Bio-Plex Data Pro™ with Bio-Plex Manager Software (5 seats) , for multi-experiment analysis and advanced data visualization
171-001513	Bio-Plex Data Pro Software (5 seats) , for multi-experiment analysis and advanced data visualization
171-001523	Bio-Plex Data Pro Plus Software , contains all the features of Bio-Plex Data Pro software with added visualization, sharing, and analysis functionality
171-STND01	Bio-Plex Manager Software (1 seat) , for instrument data evaluation and optimization



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 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** 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88
United Kingdom 020 8328 2000