

ProcartaPlex[®]

Multiplex Immunoassay

Using Magnetic Beads
For Serum, Plasma, and Cell Culture Supernatant Samples


Instructions for Mouse High Sensitivity Assays

North America

Technical Support:
Research Products:
888-810-6168
858-642-2058
tech@eBioscience.com
Clinical Products:
877-726-8559
858-642-2058
tech@eBioscience.com
Customer Service:
888-999-1371
858-642-2058
info@eBioscience.com
Fax: 858-642-2046

Europe/International*

Technical Support:
+43 1796 40 40 120
tech@eBioscience.com
Customer Service:
+43 1 796 40 40 305
europe@eBioscience.com
Fax: +43 1 796 40 40 400

 Bender MedSystems GmbH
Campus Vienna Biocenter 2
1030 Vienna, Austria
www.eBioscience.com

* Customers in countries where direct sales are not available may contact their eBioscience distributor listed at www.eBioscience.com/distributors *

**For research use only.
Not for use in diagnostic procedures.**

Limited License

Subject to the eBioscience, an Affymetrix Company, terms and conditions that govern your use of eBioscience products, eBioscience grants you a non-exclusive, non-transferable, non-sub licensable license to use this eBioscience product only in accordance with the manual and written instructions provided by eBioscience. You understand and agree that, except as expressly set forth in the eBioscience terms and conditions, no right or license to any patent or other intellectual property owned or licensable by eBioscience is conveyed or implied by this eBioscience product. In particular, no right or license is conveyed or implied to use this eBioscience product in combination with a product not provided, licensed, or specifically recommended by eBioscience for such use.

Citing ProcartaPlex® Immunoassay in Publications

When describing a procedure for publication using this product, please refer to it as the ProcartaPlex® Multiplex Immunoassay from eBioscience.

Disclaimer

eBioscience, Inc. reserves the right to change its products and services at any time to incorporate technological developments. This manual is subject to change without notice.

Although this manual has been prepared with every precaution to ensure accuracy, eBioscience, Inc. assumes no liability for any errors or omissions, not for any damages resulting from the application or use of this information.

Trademarks

©2014 Affymetrix Inc. All rights reserved. Affymetrix®, eBioscience® and ProcartaPlex® are trademarks or registered trademarks of Affymetrix, Inc. Luminex®, xMAP®, MAGPIX®, Luminex® 100/200™, and FLEXMAP 3D® are registered trademarks of the Luminex Corporation. All other trademarks are the property of their respective owners.

Contents

Intended Use	4
How it Works	4
Materials Provided and Storage Conditions	4
Precautions and Technical Hints	5
Required Equipment and Materials Not Supplied	5
Sample Preparation	5
Assay Protocol Overview	6
Preparation of Reagents	7
Assay Protocol	10
Setup of the Luminex Instruments	12
Analyzing Results	12
Troubleshooting	13
Recommended and Blank Plate Layout	14

Intended Use

This user manual is for a ProcartaPlex® High Sensitivity Immunoassay Kit from the eBioscience division of Affymetrix to perform quantitative, multiplexed protein measurements from serum, plasma, and cell culture supernatant samples using magnetic beads technology from Luminex®. Other biological samples might be suitable for use in the assay.

NOTE: For the most current version of user documentation, go to our website at www.ebioscience.com

How it Works

ProcartaPlex Immunoassays incorporate magnetic microsphere technology licensed from the Luminex Corporation to enable the simultaneous detection and quantitation of multiple protein targets in diverse matrices. The platform allows the simultaneous detection from a single sample of up to 100 protein targets on the Luminex 200/100 and FLEXMAP 3D® platforms and 50 protein targets on the MAGPIX® platform.

Materials Provided and Storage Conditions

ProcartaPlex Immunoassay Kits contain the components listed below. Refer to the Certificate of Analysis for quantities and details of components supplied. Expiration date is stated on the kit when stored between 2-8°C. Do not use past kit expiration date.

Components Supplied	Pre-mixed Panels	Simplex Kit	Basic Kit	Mix&Match Panels
Antigen Standards, premixed	√	√		√
Detection Antibody, (50X) ¹	√	√		
Detection Antibody, premixed (1X) ¹				√
Magnetic Beads, Simplexes (50X) ¹		√		
Magnetic Beads, premixed (1X) ¹	√			√
Streptavidin-PE (SA-PE) (1X) ¹	√		√	√
Amplification Reagent 1	√		√	√
Amplification Reagent 2	√		√	√
Wash Buffer Concentrate (10X) ¹	√		√	√
Detection Antibody Diluent ¹	√		√	
Universal Assay Buffer (1X) ¹	√		√	√
Reading Buffer ¹	√		√	√
PCR 8-Tube Strip	√		√	√
96-Well Flat Bottom Plate	√		√	√
Black Microplate Lid	√		√	√
Plate Seals	√		√	√

¹ Contains sodium azide. See WARNING.

WARNING: All chemicals should be considered potentially hazardous. We recommend that this product and its components be handled by those trained in laboratory techniques and be used according to the principles of good laboratory practice. This kit contains small quantities of sodium azide. Sodium azide is highly toxic and reactive in the pure form. At this product's concentration, though not classified as hazardous, build up of sodium azide may react with lead and copper plumbing to form highly reactive explosive metal azide. Dispose of the product in accordance with all State and local regulations.

Precautions and Technical Hints

- Thoroughly read this user manual and Certificate of Analysis that is included with the assay kit. The product insert may contain specific instructions for proper use of your kit.
- For Luminex 100/200 and FLEXMAP 3D instruments initiate the startup protocol to warm up the lasers for at least 30 minutes. Ensure that the Luminex machine is calibrated according to the manufacturer's instructions. MAGPIX instrument doesn't require additional warm up.
- When working with samples and standards, change the pipette tips after every transfer and avoid creating bubbles when pipetting.
- During the incubation steps, cover the 96-Well Flat Bottom Plate with the Black Microplate Lid provided in the kit to minimize exposure of the beads to light.
- Be careful not to invert the 96-Well Flat Bottom Plate during the assay or allow contents from one well to mix with another well.
- Use a multi-channel pipette and reagent reservoirs whenever possible to achieve optimal assay precision.
- Store the reconstituted standards on ice before adding to the 96-Well Flat Bottom Plate

Required Equipment and Materials Not Supplied

- MAGPIX, Luminex® 100/200, FLEXMAP 3D, or Luminex-based Instrument.
- Glass-distilled or deionized water.
- Adjustable single and multi channel pipettes with disposable tips.
- Multichannel pipette reservoir.
- Beakers, flasks, cylinders necessary for preparation of reagents.
- Hand-Held Magnetic Plate Washer, Vortex mixer and Microtiter plate shaker.

IMPORTANT: Use only single use trays for all reagents and buffers.

Sample Preparation

- For frozen samples, thaw samples on ice and mix well by vortexing followed by centrifugation to remove particulates. Avoid multiple freeze/thaw cycles.
- If there is a high lipid content in the sample, centrifuge at 10,000 x g for 10 min at 2-8 °C and transfer contents to a new tube.

For preparation instruction of tissue homogenates and lysates from cultured cells refer to www.ebioscience.com/resources/best-protocols/multiplexing.htm

Plasma Sample Preparation

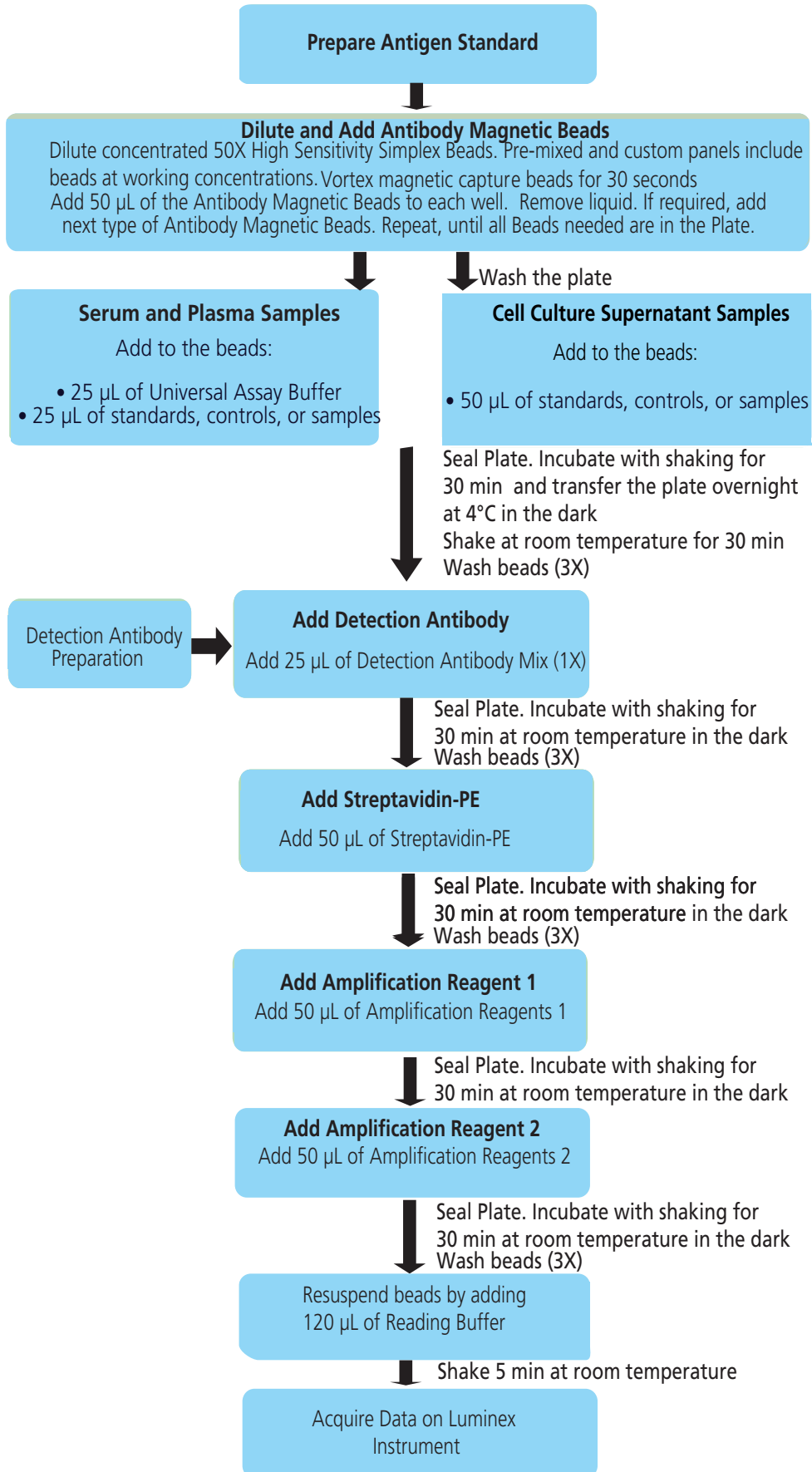
Action
A. Collect samples in sodium citrate or EDTA tubes. When using heparin as an anticoagulant, no more than 10 IU of heparin per mL of blood collected should be used since an excess of heparin may give falsely high values of some of the analytes.
B. Centrifuge samples at 1,000 x g at 4 °C for 10 min within 30 min of blood collection.
C. Collect the plasma fraction.
D. Use immediately or aliquot and store at -80 °C.

Serum Sample Preparation

We recommend to spin down serum samples at 1000 x g for 10 min at 20-25 °C before running the assay.

Action
A. Allow blood to clot for 20-30 min at 20-25 °C.
B. Centrifuge at 1,000 x g for 10 min at 20-25 °C.
C. Use immediately or aliquot and store at -80 °C.

Assay Protocol Overview



Preparation of Reagents

Preparing 1X Wash Buffer

Bring the Wash Buffer Concentrate (10X) to room temperature and vortex for 15 seconds. Mix 20 mL of the Wash Buffer Concentrate (10X) with 180 mL ddH₂O. Wash Buffer (1X) can be stored at 2-8 °C for up to 6 months.

NOTE: Wash Buffer Concentrate volume might not be sufficient if using automated plate washer. For bulk orders use Cat. No. EPX-66666-001.

Preparing 1X High Sensitivity Simplex Beads

High Sensitivity Simplex kits are provided with concentrated 50X beads. Therefore, dilution of concentrated High Sensitivity Simplex beads is required.

- A. Vortex each High Sensitivity Simplex bead vial (50X) for 30 sec and add 100 µL of each High Sensitivity Simplex bead vial (50X) to a mixing bottle if using a whole plate (otherwise adjust the volume accordingly).
- B. Fill up with Wash Buffer (1X) to a final volume of 5 mL. To combine 2 or more different bead vials follow the table below (using a whole plate):

# of different Simplex Bead vials to be mixed	Total Volume of mixed Bead Solution	Volume of Wash Buffer (1X) to add
1	100 µL	4900 µL
2	200 µL	4800 µL
3	300 µL	4700 µL
4	400 µL	4600 µL
5	500 µL	4500 µL
6	600 µL	4400 µL

Preparing 1X Detection Antibody Mixture

For High Sensitivity Simplex and Pre-mixed Panels (see label: 50X), detection antibody is provided at 50X concentration. Add 60 µL of each detection antibody concentrate to the mixing bottle and bring volume to a total of 3 mL using detection antibody diluent if using a whole plate (otherwise adjust the volume accordingly).

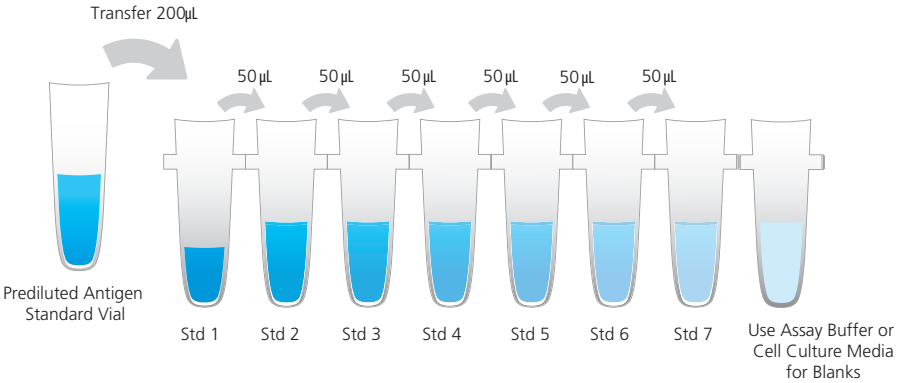
# of Vials of Detection Antibody	Total Volume of Detection Antibody	Volume of Diluent to add
1	60 µL	2940 µL
2	120 µL	2880 µL
3	180 µL	2820 µL
4	240 µL	2760 µL
5	300 µL	2700 µL
6	360 µL	2640 µL

IMPORTANT: High Sensitivity Mix&Match Panels contain 1X detection antibody and no further dilution is required.

Preparing Antigen Standard

The majority of kits is supplied with lyophilized multi-standards containing a mix of multiple standard proteins. Some kits contain multiple sets of standards, each with a unique lot number, that require pooling prior to use. Each kit is shipped with two identical vials of each premixed antigen standard set from the same lot to permit the user to run the assay twice if running a partial plate. When preparing Antigen Standards, the final volume after reconstitution and pooling should be 250 μ L. When combining multiple kits, ensure that the Antigen Standards of your analytes of interest are only present in one of the used standard vials.

Step	Action																														
Step 1. Reconstitution and pooling of Standards	A. Centrifuge each different antigen standard vial(s) at 2000 x g for 10 sec.																														
	B. Add 50 μ L of sample type specific buffer into each standard vial. If you want to measure serum or plasma samples, use Universal Assay Buffer (1X) to reconstitute the standard, if you want to analyse cell culture supernatant samples, use the cell culture medium that was used to culture the cells to dissolve the standard.																														
	C. Gently vortex the vial(s) for 10 seconds and centrifuge at 2000 x g for 10 seconds to collect contents at the bottom of the vial(s).																														
	D. Incubate on ice for 10 min to ensure complete reconstitution.																														
	E. Pool entire contents of each vial into one of the vials and fill up with sample type specific buffer to a total volume of 250 μ L.																														
	F. Gently vortex the vial for 10 seconds and centrifuge at 2000 x g for 10 seconds to collect contents at the bottom of the vial.																														
	<table border="1"> <thead> <tr> <th># of Standard Sets</th> <th>Reconstitution Volume per vial</th> <th>Pooled Volume</th> <th>Buffer to add</th> <th>Total Volume</th> </tr> </thead> <tbody> <tr> <td>1</td> <td>50 μL</td> <td>50 μL</td> <td>200 μL</td> <td>250 μL</td> </tr> <tr> <td>2</td> <td>50 μL</td> <td>100 μL</td> <td>150 μL</td> <td>250 μL</td> </tr> <tr> <td>3</td> <td>50 μL</td> <td>150 μL</td> <td>100 μL</td> <td>250 μL</td> </tr> <tr> <td>4</td> <td>50 μL</td> <td>200 μL</td> <td>50 μL</td> <td>250 μL</td> </tr> <tr> <td>5</td> <td>50 μL</td> <td>250 μL</td> <td>0 μL</td> <td>250 μL</td> </tr> </tbody> </table>	# of Standard Sets	Reconstitution Volume per vial	Pooled Volume	Buffer to add	Total Volume	1	50 μ L	50 μ L	200 μ L	250 μ L	2	50 μ L	100 μ L	150 μ L	250 μ L	3	50 μ L	150 μ L	100 μ L	250 μ L	4	50 μ L	200 μ L	50 μ L	250 μ L	5	50 μ L	250 μ L	0 μ L	250 μ L
# of Standard Sets	Reconstitution Volume per vial	Pooled Volume	Buffer to add	Total Volume																											
1	50 μ L	50 μ L	200 μ L	250 μ L																											
2	50 μ L	100 μ L	150 μ L	250 μ L																											
3	50 μ L	150 μ L	100 μ L	250 μ L																											
4	50 μ L	200 μ L	50 μ L	250 μ L																											
5	50 μ L	250 μ L	0 μ L	250 μ L																											
	G. Predilute the reconstituted antigen standard 1:10 (25 μ L of reconstituted antigen standard with 225 μ L diluent. For serum or plasma samples, use Universal Assay Buffer and for cell culture supernatant samples use the cell culture media that was used to culture the cells.																														

Step	Action
Step 2. Prepare 4-Fold Serial Dilution	<p>A. Prepare a 4-fold serial dilution of the reconstituted and prediluted standard using the PCR 8-tube strip provided. Label tubes Std1, Std2, Std3, Std4, Std5, Std6 and Std7.</p> <p>B. Add 200 μL of the reconstituted and prediluted antigen standards into the first tube of the strip tube and label as Standard 1 (Std1).</p> <p>C. Add 150 μL of sample type specific standard buffer into Std tubes 2-7. Use Universal Assay Buffer for serum or plasma samples and cell culture media for culture supernatant samples.</p> <p>D. Transfer 50 μL of the reconstituted and prediluted antigen standards from Tube 1 into Tube 2.</p> <p>E. Mix by pipetting up and down for a total of 10 times.</p> <p>F. Change the pipette tip and transfer 50 μL of the mixed standards from Tube 2 into Tube 3.</p> <p>G. Mix by pipette up and down 10 times.</p> <p>H. Repeat steps E-H for Std tubes 4-7.</p> <p>I. Add 200 μL of Universal Assay Buffer or cell culture media into tube 8 which serves as a blank. Keep on ice until ready to use.</p>
	

Assay Protocol

IMPORTANT: Use only single use trays for all reagents and buffers.

Step	Action
Step 1. Define the plate map	Mark the standard, sample and blank wells using the plate map at the end of this manual.
Step 2. Addition of Magnetic Beads to the plate	<p>IMPORTANT: If working with High Sensitivity Simplex kits only, add 50 μL per well of diluted 1X High Sensitivity Simplex beads prepared according to instructions provided under "Preparation of Reagents".</p> <p>Note that High Sensitivity Pre-mixed and Mix&Match Panels contain only one bead vial at working concentration 1X.</p> <ol style="list-style-type: none"> Vortex the Magnetic Bead vial for 30 sec. Add 50 μL of the Magnetic Bead solution to each well of the plate. Use a multichannel pipette for this step as well as for the steps below.
Step 3. Wash Magnetic Beads	<ol style="list-style-type: none"> Securely insert the 96-Well Flat Bottom Plate into the Hand-Held Magnetic Plate Washer and ensure that the plate is held in place by the tabs and wait 2 min to allow the Magnetic Beads to accumulate on the bottom of each well. Remove the liquid in the wells by quickly inverting the Hand-Held Magnetic Plate Washer and 96-Well Flat Bottom Plate assembly over a sink or waste container. Do not remove the 96-Well Flat Bottom Plate from the Hand-Held Magnetic Plate Washer. Blot the inverted assembly onto several layers of paper towels or absorbent surface to remove any residual solution. Add 150 μL of Wash Buffer (1X) into each well and wait 30 seconds to allow the beads to accumulate on the bottom of each well. Remove the Wash Buffer in the wells by quickly inverting the plate and then blotting onto an absorbent towel to remove any residual solution. Remove the 96-Well Flat Bottom Plate from the Hand Held Magnetic Plate Washer and proceed to the next step.
Step 4. Add sample type-specific buffer, samples, standards and blanks and incubate	<ol style="list-style-type: none"> Add 25 μL of Universal Assay Buffer (1X) to each well for serum and plasma followed by 25 μL of prepared standards or samples into dedicated wells. For cell culture supernatant samples, add 50 μL prepared standards or samples into dedicated wells. For wells designated as blanks, add an additional 25 μL of Universal Assay Buffer for serum or plasma samples. For cell culture supernatant samples, add 50 μL of cell culture medium. Incubate the 96 well plate overnight. Seal the plate with the provided Plate Seal. Cover the plate with the Black Microplate Lid and shake for 30 min at RT at 500 rpm, then transfer the plate to 4°C and store on a level surface. After overnight incubation, shake the plate for an additional 30 min at RT at 500 rpm.
Step 5. Wash the 96-Well Plate	Wash the plate for a total of three washes using "Step 3. Wash Magnetic Beads".

Step	Action
Step 6. Add Detection Antibody Mixture and incubate	<ul style="list-style-type: none"> A. Add 25 μL of Detection Antibody Mixture (1X) to each well. B. Seal the plate with a new Plate Seal, cover plate with the Black Microplate Lid and incubate 30 min on a plate shaker at RT at 500 rpm.
Step 7. Wash the 96-Well Plate	Wash plate for a total of three washes using "Step 3. Wash Magnetic Beads".
Step 8. Add SAPE and incubate	<ul style="list-style-type: none"> A. Add 50 μL of SAPE solution to each well. B. Seal the plate with a new Plate Seal, cover plate with the Black Microplate Lid and incubate 30 min on a plate shaker at RT at 500 rpm.
Step 9. Wash the 96-Well Plate	Wash plate for a total of three washes using "Step 3. Wash Magnetic Beads".
Step 10. Add Amplification Reagent 1	<ul style="list-style-type: none"> A. Add 50 μL of Amplification Reagent 1 into each well. B. Seal the plate with a new Plate Seal. C. Remove the plate from Hand-Held Magnetic Plate Washer and cover plate with the Black Microplate Lid D. Incubate 30 min on a plate shaker at RT at 500 rpm. After incubation immediately proceed to next step. (DO NOT WASH)
Step 11. Add Amplification Reagent 2	<ul style="list-style-type: none"> A. Add 50 μL of Amplification Reagent 2 into each well. B. Seal the plate with a new Plate Seal and cover plate with the Black Microplate Lid. C. Incubate 30 min on a plate shaker at RT at 500 rpm.
Step 12. Wash the 96-Well Plate	Wash plate for a total of three washes using "Step 3. Wash Magnetic Beads".
Step 13. Prepare the 96-Well Plate for Analysis on a Luminex Instrument	<ul style="list-style-type: none"> A. Add 120 μL of Reading Buffer into each well. B. Seal the plate with a new Plate Seal, cover plate with the Black Microplate Lid and incubate 5 min on a plate shaker at RT at 500 rpm. C. Remove Plate Seal and run the plate on a Luminex Instrument.

Setup of the Luminex Instruments

Instrument	Sample Size	DD Gate	Timeout	Bead Event/Bead Region
Luminex 100/200 FLEXMAP 3D	50 μ L	5,000 - 25,000	60 sec	50-100
MAGPIX	50 μ L	N/A	N/A	50-100

Prior to running the assay, ensure that the probe height has been calibrated with 96-Well Flat Bottom Plate supplied with the kit. Failure to adjust the probe height can cause damage to the instrument or low bead count. The Luminex system allows for calibration of Low and High RP1 target values. We recommend RP1 Low target value settings for ProcartaPlex Immunoassays. When entering the information into the Luminex Acquisition Software please refer to the Certificate of Analysis provided with the kit for bead region and S1 values for each analyte of the current lot.

NOTE: If there is a malfunction of the Luminex Instrument or software during the run, the 96-Well Flat Bottom Plate can be re-read. Remove the 96-Well Flat Bottom Plate from the instrument, insert the 96-Well Flat Bottom Plate into the Hand-Held Magnetic Plate Washer, wait 2 min, then remove the buffer in the wells by quickly inverting the 96-Well Flat Bottom Plate over a sink or waste container. Blot the assembly onto several layers of paper towels to remove any residual solution. Resuspend the beads in 120 μ L of Reading Buffer, remove from the Hand-Held Magnetic Plate Washer, seal the 96-Well Flat Bottom Plate with a new Plate Seal and Lid and shake at 500 rpm for 5 min at room temperature. The assayed samples may take longer to read since there will be less beads in the well.

Analyzing Results

The concentration of the samples can be calculated by plotting the expected concentration of the standards against the MFI generated by each standard. A 4PL or 5PL algorithm is recommended for the best curve fit. Analyze the assayed samples according to the operation manual for the Luminex Instrument (e.g MAGPIX[®], Luminex[®] 100/200[™], FLEXMAP 3D[®]). We offer a free and robust analysis software package for data analysis. ProcartaPlex Analyst 1.0 can be download at:

www.ebioscience.com/resources/procartaplex-analyst-1.0-software.htm

Troubleshooting

Observation	Probable Cause	Recommend Solution
Low Flow Rate	Samples/beads are stuck in flow cell	Remove the 96-Well Plate and perform a wash and rinse cycle.
High CVs	Samples and antigen standards not stored on ice	Prepare the samples and standards on ice before setting up the assay.
	Contamination from re-using the Plate Seal	Use a new Plate Seal for each incubation step.
	Incomplete washing	After adding the standards and samples, it is very important that any excess standards are removed during the wash step.
	Contamination from contents from adjacent wells	Avoid splashing the Wash Buffer during wash steps into adjacent wells.
	Poor pipetting techniques	Use a multichannel pipettor and careful pipette techniques. Avoid touching pipette tips to sides of the wells when adding Wash Buffer.
Limited dynamic range for BioPlex software users	Instrument calibrated at high PMT settings	Calibrate the instrument using the CAL2 Low RP1 target value.
Low bead count	Volume of bead solution is too low	Add 120 μ L Reading Buffer into each well and shake at 500 rpm for 5 min at room temperature to resuspend beads prior to reading on the Luminex Instrument.
	High bead aggregation	Vortex the bead suspension well before using in the assay and ensure that the beads are properly mixed during the incubation steps.
	Dyes contained in the beads are photo-bleached from overexposure to light	Store bead solution and the 96-well plate in the dark.
	Samples causing the instrument to clog	Remove the 96-Well Flat Bottom Plate and perform a wash and rinse to the instrument. Rerun the assay with further dilution of samples
	Probe height is incorrect	Refer to the Luminex Manual for proper adjustment of the needle height.
	Instrument needle is partially clogged	Replace or clean needle according to the manufacturer's recommendations.
	Beads stuck to the bottom of the plate	Confirm that the plate shaker is set to 500 rpm and shaking for at least 5 min before reading.
	Air bubble in the sample loop	Refer to the Luminex manual for proper removal of the air bubble.
Low signal or sensitivity	Standards not reconstituted and diluted correctly	Prepare fresh antigen standards following the instructions in the Preparing Antigen Standards section.
Poor recovery	Did not use appropriate cell culture media to prepare the standards	Use the same cell culture media that is used to culture the cells.
	Samples and antigen standards were not stored on ice	Prepare the samples and standards on ice before setting up the assay.

Recommended and Blank Plate Layout

Standards		Samples									
Standard 1	Standard 1	1	1	9	9	17	17	25	25	33	33
Standard 2	Standard 2	2	2	10	10	18	18	26	26	34	34
Standard 3	Standard 3	3	3	11	11	19	19	27	27	35	35
Standard 4	Standard 4	4	4	12	12	20	20	28	28	36	36
Standard 5	Standard 5	5	5	13	13	21	21	29	29	37	37
Standard 6	Standard 6	6	6	14	14	22	22	30	30	38	38
Standard 7	Standard 7	7	7	15	15	23	23	31	31	39	39
Blank	Blank	8	8	16	16	24	24	32	32	40	40

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