

ProcartaPlex™ Multiplex Immunoassay USER GUIDE

Instructions for Mouse Assays (384 tests)

Using Magnetic Beads for Cell Culture Supernatant Samples

for use with FLEXMAP 3D™

Publication Number MAN0017963

Revision A.0 (30)



The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0017962

Revision	Date	Description
A.0	22 June 2019	New document

Limited Use Label Licence No. 660: Label License/Sticker for Luminex Assay Product: Notice to Purchaser: By opening the packaging containing this Assay Product (which contains fluorescently labeled microsphere beads authorized by Luminex Corporation) or using this Assay Product in any manner, you are consenting and agreeing to be bound by the following terms and conditions. You are also agreeing that the following terms and conditions constitute a legally valid and binding contract that is enforceable against you. If you do not agree to all of the terms and conditions set forth below, you must promptly return this Assay Product for a full refund prior to using it in any manner.

You, the customer, acquire the right under Luminex Corporation's patent rights, if any, to use this Assay Product or any portion of this Assay Product, including without limitation the microsphere beads contained herein, only with Luminex Corporation's fluorescent analytical test instrumentation marketed under the name Luminex Instrument.

Limited Use Label Licence No. 661: Standard Terms And Conditions For Use of Luminex Assay Product: Note to Purchaser: By opening the packaging containing this assay product or cassette ("Product") or by using such Product in any manner, you are consenting and agreeing to be bound by the following terms and conditions. You are also agreeing that the following terms and conditions constitute a legally valid and binding contract that is enforceable against you. If you do not agree to all of the terms and conditions set forth below, you must promptly return the Product for a full refund prior to using them in any manner.

- Acceptance - ALL SALES ARE SUBJECT TO AND EXPRESSLY CONDITIONED UPON THE TERMS AND CONDITIONS CONTAINED HEREIN, AND UPON BUYER'S ASSENT THERETO. NO VARIATION OF THESE TERMS AND CONDITIONS SHALL BE BINDING UPON LUMINEX CORPORATION ("LUMINEX") UNLESS AGREED TO IN WRITING AND SIGNED BY AN AUTHORIZED REPRESENTATIVE OF LUMINEX.

For purposes of this agreement, "Seller" shall mean either Luminex, if the Product is purchased directly from Luminex, or a Luminex authorized reseller. Buyer, by accepting the Product shall be deemed to have assented to the terms and conditions set forth herein, notwithstanding any terms contained in any prior or later communications from Buyer and whether or not Seller shall specifically or expressly object to any such terms.
- Warranties - Notwithstanding Buyer's acceptance thereof, if Product is purchased directly from Luminex, Luminex warrants that the Product shall conform to the quantity and content stated on the label and perform in all material respects consistent with Product specifications accompanying the Product until the expiration date set forth on the Product label. If Product is purchased from a Luminex authorized reseller, any warranty obligations shall be provided in writing directly by such Luminex authorized reseller to Buyer. THIS WARRANTY IS EXCLUSIVE AND LUMINEX MAKES NO OTHER WARRANTY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, ANY IMPLIED WARRANTY OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. Seller's warranties made in connection with this sale shall not be effective if Seller has determined, in its sole discretion, that Buyer has misused the Product in any manner, has failed to use the Product in accordance with industry standards or practices, or has failed to use the Product in accordance with instructions, if any, furnished by Seller.

BUYER'S EXCLUSIVE REMEDY WITH RESPECT TO PRODUCT PROVED TO SELLER'S SATISFACTION TO BE DEFECTIVE OR NONCONFORMING SHALL BE REPLACEMENT OF SUCH PRODUCTS WITHOUT CHARGE OR REFUND OF THE PURCHASE PRICE, IN SELLER'S SOLE DISCRETION, UPON THE RETURN OF SUCH PRODUCTS IN ACCORDANCE WITH SELLER'S INSTRUCTIONS. NEITHER SELLER NOR LUMINEX NOR ITS AFFILIATES SHALL IN ANY EVENT BE LIABLE FOR INCIDENTAL, CONSEQUENTIAL OR SPECIAL DAMAGES OF ANY KIND RESULTING FROM ANY USE OR FAILURE OF THE PRODUCT, EVEN IF SELLER OR LUMINEX OR ITS AFFILIATE HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES, INCLUDING, WITHOUT LIMITATION, LIABILITY FOR LOSS OF WORK IN PROGRESS, DOWN TIME, LOSS OF REVENUE OR PROFITS, FAILURE TO REALIZE SAVINGS, LOSS OF PRODUCTS OF BUYER OR OTHER USE OR ANY LIABILITY OF BUYER TO A THIRD PARTY ON ACCOUNT OF SUCH LOSS, OR FOR ANY LABOR OR ANY OTHER EXPENSE, DAMAGE OR LOSS OCCASIONED BY SUCH PRODUCT INCLUDING PERSONAL INJURY OR PROPERTY DAMAGE UNLESS SUCH PERSONAL INJURY OR PROPERTY DAMAGE IS CAUSED BY SELLER'S GROSS NEGLIGENCE.
- Buyer's Use of Product - Buyer agrees that no rights or licenses under Luminex's patents shall be implied from the sale of the Product, except as expressly provided herein or as specifically agreed to in writing by Luminex, and Buyer does not receive any right under Luminex's patent rights hereunder. Buyer acknowledges and agrees that the Product is sold and licensed only for use with Luminex's instrumentation. In order to maintain the quality of the Product, Buyer may use this Product only once on a single use basis and shall not reuse this Product under any circumstances. Buyer further acknowledges that the Product has not received clearance from the United States Food and Drug Administration or other federal, state or local regulatory agencies and has not been tested by Seller or Luminex for safety or efficacy in food, drug, medical device, cosmetic, commercial or any other use, unless otherwise stated on the Product label or in Seller's technical specifications or material data sheets furnished to Buyer. Buyer expressly represents and warrants to Seller that Buyer will use the Product in accordance with the Product label, if applicable, and will properly test and use any Product in accordance with the practices of a reasonable person who is an expert in the field and in strict compliance with the United States Food and Drug Administration and all applicable domestic and international laws and regulations, now and hereinafter enacted.

BUYER HEREBY GRANTS TO LUMINEX A NONEXCLUSIVE, WORLDWIDE, UNRESTRICTED, ROYALTY-FREE, FULLY PAID-UP LICENSE, WITH THE RIGHT TO GRANT AND AUTHORIZE SUBLICENSES, UNDER ANY AND ALL PATENT RIGHTS IN INVENTIONS COMPRISING MODIFICATIONS, EXTENSIONS, OR ENHANCEMENTS MADE BY BUYER TO THE PRODUCT OR TO THE MANUFACTURE OR USE OF THE PRODUCT ("IMPROVEMENT PATENTS"), TO MAKE, HAVE MADE, USE, IMPORT, OFFER FOR SALE OR SELL ANY AND ALL OF THE PRODUCT; EXPLOIT ANY AND ALL METHODS OR PROCESSES; AND OTHERWISE EXPLOIT IMPROVEMENT PATENTS FOR ALL PURPOSES. NOTWITHSTANDING THE FOREGOING, "IMPROVEMENT PATENTS" SPECIFICALLY EXCLUDES PATENT CLAIMS CONCEIVED AND REDUCED TO PRACTICE BY BUYER

CONSISTING OF METHODS OF SAMPLE PREPARATION, THE COMPOSITION OF MATTER OF THE SPECIFIC CHEMISTRIES OF THE ASSAYS DEVELOPED BY BUYER AND METHODS OF PERFORMING THE ASSAYS (I.E., THE PROTOCOL FOR THE ASSAY).

Buyer has the responsibility and hereby expressly assumes the risk to verify the hazards and to conduct any further research necessary to learn the hazards involved in using the Product. Buyer also has the duty to warn Buyer's customers, employees, agents, assigns, officers, successors and any auxiliary or third party personnel (such as freight handlers, etc.) of any and all risks involved in using or handling the Product. Buyer agrees to comply with instructions, if any, furnished by Seller or Luminex relating to the use of the Product and to not misuse the Product in any manner. Buyer shall not reverse engineer, decompile, disassemble or modify the Product. Buyer acknowledges that Luminex retains ownership of all patents, trademarks, trade secrets and other proprietary rights relating to or residing in the Product and Buyer receives no rights to such intellectual property rights by virtue of its purchase of Product other than as expressly set forth herein. Buyer shall have no right to use any trademarks owned or licensed to Luminex without the express written permission of Luminex.

4. Buyer's Representations, Release and Indemnity - Buyer represents and warrants that it shall use the Product in accordance with Paragraph 3, "Buyer's Use of Product," and that any such use of the Product will not violate any law, regulation, judicial order or injunction. Buyer agrees to release, discharge, disclaim and renounce any and all claims, demands, actions, causes of action and/or suits in law or equity, now existing or hereafter arising, whether known or unknown, against Seller and Luminex, and their respective officers, directors, employees, agents, successors and assigns (collectively the "Released Parties"), with respect to the use of the Product. Buyer agrees to indemnify and hold harmless the Released Parties from and against any suits, losses, claims, demands, liabilities, costs and expenses (including attorney, accounting, expert witness, and consulting fees) that any of the Released Parties may sustain or incur as a result of any claim against such Released Party based upon negligence, breach of warranty, strict liability in tort, contract or any other theory of law or equity arising out of, directly or indirectly, the use of the Product or by reason of Buyer's failure to perform its obligations contained herein. Buyer shall fully cooperate with the Released Parties in the investigation and determination of the cause of any accident involving the Product which results in personal injury or property damage and shall make available to the Released Parties all statements, reports, recordings and tests made by Buyer or made available to Buyer by others.
5. Patent Disclaimer - Neither Seller nor Luminex warrants that the use or sale of the Product will not infringe the claims of any United States or other patents covering the Product itself or the use thereof in combination with other products or in the operation of any process.

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2019 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■ ProcartaPlex™ multiplex immunoassay	6
Product use	6
How it works	6
Materials provided and storage conditions	6
Precautions and technical hints	7
Required equipment and materials not supplied	8
Sample preparation	8
Diluting samples with high concentration of target analytes	8
Assay protocol overview	9
Preparation of reagents	10
Antigen standard	10
Step 1. Reconstitution and pooling of standards	10
Step 2. Prepare 4-Fold serial dilution	10
Assay protocol	11
Step 1. Prepare 1X wash buffer	11
Step 2. Define the plate map	11
Step 3a. Add the antibody magnetic beads (only 1 bead set)	12
Step 3b. Add the antibody magnetic beads (2 or more beads sets)	12
Step 4. Wash antibody magnetic beads	12
Step 5. Add sample type-specific buffer, samples, standards and blanks and incubate	13
Step 6. Wash the 384-Well plate	13
Step 7. Prepare 1X detection antibody mixture	13
Step 8. Add detection antibody mixture and incubate	13
Step 9. Wash the 384-Well plate	14
Step 10. Add SAPE and incubate	14
Step 11. Wash the 384-Well plate	14
Step 12. Prepare the 384-Well plate for analysis on a Luminex™ instrument	14
Setup of the Luminex™ instruments	14
Analyzing results	15
Troubleshooting	15
Blank plate layout	16

■ APPENDIX A Safety	17
Chemical safety	18
Biological hazard safety	19
Documentation and support	20
Customer and technical support	20
Limited product warranty	20



ProcartaPlex™ multiplex immunoassay

Product use

This user manual is for a ProcartaPlex™ Immunoassay Kit from Thermo Fisher Scientific 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.thermofisher.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	Custom Panels
Antigen Standards, premixed	√	√		√
Detection Antibody, premixed (50X) ^[1]	√	√		
Detection Antibody, premixed (1X) ^[1]				√
Antibody Magnetic Beads, premixed ^[1]	√	√		√



Components Supplied	Pre-mixed Panels	Simplex Kit	Basic Kit	Custom Panels
Streptavidin-PE (SA-PE) (1X) ^[1]	✓		✓	✓
Wash Buffer Concentrate (10X) ^[1]	✓		✓	✓
Detection Antibody Diluent ^[1]	✓		✓	
Reading Buffer ^[1]	✓		✓	✓
PCR 8-Tube Strip	✓		✓	✓
Filter Plate ^[2]				
Black Microplate Lid	✓		✓	✓
Plate Seals	✓		✓	✓

^[1] Contains sodium azide. See WARNING on page 7.

^[2] 384-Well Filter Plate (Cat. No. EPX38488-000) is not included into this kit and can be ordered separately.



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.

Note: ProcartaPlex™ Supplemental Products can be also purchased as stand alone items. For more information contact our technical support.

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 FLEXMAP 3D™ instrument 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.
- 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 384-Well Filter Plate with the Black Microplate Lid provided in the kit to minimize exposure of the beads to light.
- Be careful not to invert the 384-Well Filter 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 (including standard diluent sets) on ice before adding to the 384-Well Filter Plate.

Required equipment and materials not supplied

- FLEXMAP 3D™.
- 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.
- Vacuum Filtration Manifold.

Sample preparation

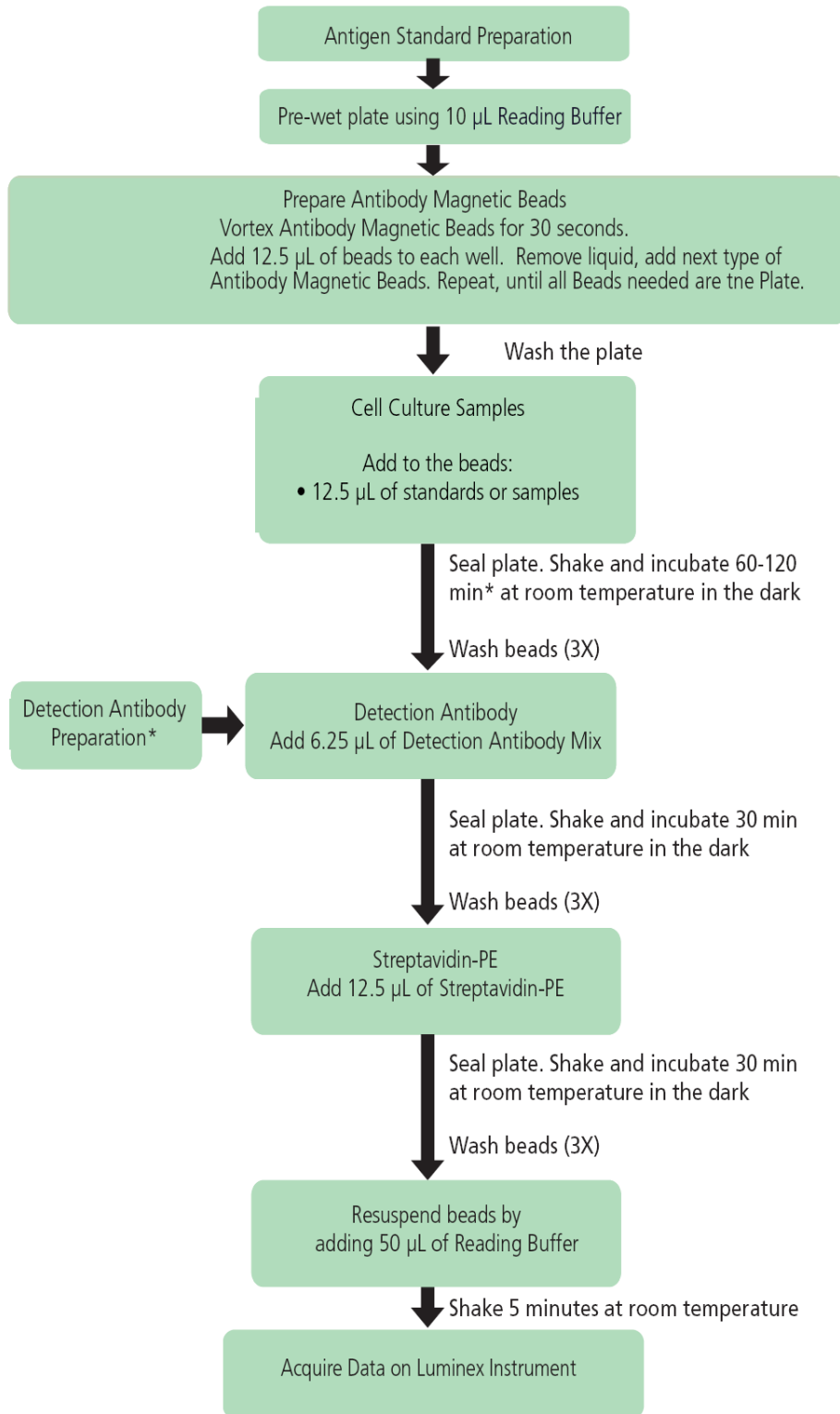
- For frozen samples, thaw samples on ice and mix well by vortexing followed by centrifugation to remove particulates. Avoid multiple freeze/thaw cycles.
- If there is a high lipid content in the sample, centrifuge at 10,000 x g for 10 min at 2-8 °C and transfer contents to a new tube.

Diluting samples with high concentration of target analytes

You may need to further dilute your samples if the analyte concentration is above the assay upper limit of quantitation (ULOQ). For cell culture supernatant samples, we recommend using the medium that was used to culture the cells.



Assay protocol overview



*For assays that require higher sensitivity, 120 min or overnight incubation is recommended.



Preparation of reagents

Antigen standard

Carefully review the Certificate of Analysis for kit specific Antigen Standard preparation instructions. 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 µL. When combining multiple kits, ensure that the Antigen Standards of your analytes of interest are only present in one of the used standard vials. For instructions for combining more than 5 antigen standard sets contact our technical support.

Step 1. Reconstitution and pooling of standards

1. Centrifuge one of the antigen standard set vial(s) at 2000 x g for 10 sec.
2. Add 50 µL of cell culture media that was used to culture the cells.
3. Gently vortex the vial(s) for 30 seconds and centrifuge at 2000 x g for 10 seconds to collect contents at the bottom of the vial(s).
4. Incubate on ice for 10 min to ensure complete reconstitution.
5. Pool entire contents of each vial into one of the vials and add sample type specific buffer to quantity sufficient (q.s.) to 250 µL. (See table below for example).
6. Gently vortex the vial for 30 seconds and centrifuge at 2000 x g for 10 seconds to collect contents at the bottom of the vial.

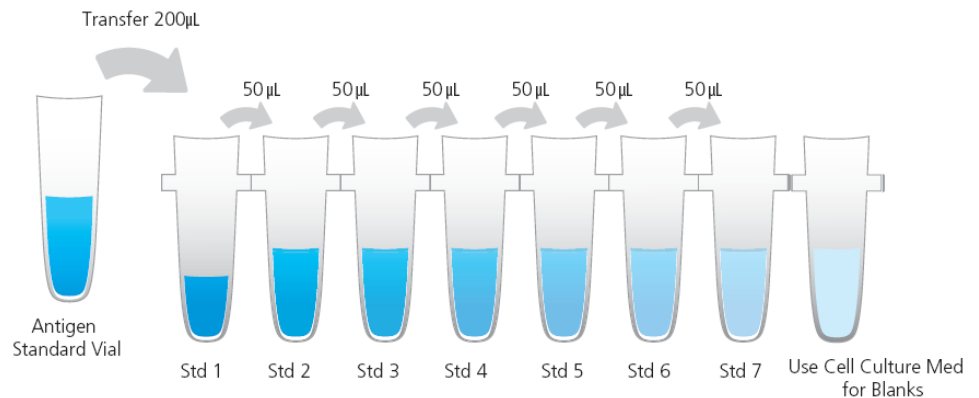
# of Standard Sets	Reconstitution Volume per vial	Pooled Volume	Buffer to q.s.	Total Volume
1	50 µL	50 µL	200 µL	250 µL
2	50 µL	100 µL	150 µL	250 µL
3	50 µL	150 µL	100 µL	250 µL
4	50 µL	200 µL	50 µL	250 µL
5	50 µL	250 µL	0 µL	250 µL

Step 2. Prepare 4-Fold serial dilution

1. Refer to Certificate of Analysis for the value of each premixed standard with assigning S1 values for each analyte for the current lot.
2. Prepare a 4-fold serial dilution of the reconstituted standard(s) using the PCR 8-tube strip provided. Label tubes Std1, Std2, Std3, Std4, Std5, Std6 and Std7.
3. Add 200 µL of the reconstituted antigen standards into the first tube of the strip tube and label as Standard 1 (Std1).
4. Add 150 µL of cell culture media into Std tubes 2-7.



5. Transfer 50 μL of the reconstituted antigen standards from Tube 1 into Tube 2.
6. Mix by pipetting up and down for a total of 10 times.
7. Change the pipette tip and transfer 50 μL of the mixed standards from Tube 2 into Tube 3.
8. Mix by pipette up and down 10 times.
9. Repeat step 5-step 8 for Std tubes 4-7.
10. Add 200 μL of cell culture media into tube 8 which serves as a blank. Keep on ice until ready to use.



Assay protocol

The Antibody Magnetic Beads can be used with a Vacuum Filtration Manifold and the 384-Well Filter Plate. The Vacuum Manifold can be properly calibrated by placing a standard 384-Well Filter Plate on top of the manifold and then turning on the vacuum. Press down on all 4 corners of the standard 384-Well Filter Plate to form a tight seal and adjust the pressure so that it takes 4-6 seconds to evacuate 50 μL of Wash Buffer from the wells. If the vacuum is too high, beads can get trapped or pulled through the filter. Turn off vacuum as soon as the solution filters through the wells and remove the plate from the manifold.

Step 1. Prepare 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.

Step 2. Define the plate map

1. Mark the standard, sample and blank wells using the plate map at the end of this manual.
2. Add 10 μL Reading Buffer (1X) to the 384-Well Filter Plate to pre-wet the wells. Aspirate using the Vacuum Manifold. Blot the bottom of the plate after filtration.



Step 3a. Add the antibody magnetic beads (only 1 bead set)

1. Vortex the Antibody Magnetic Beads for 30 sec.
2. Add 12.5 µL of the Antibody Magnetic Beads to each well. Use a multichannel pipette for this step as well as the steps below.

Step 3b. Add the antibody magnetic beads (2 or more beads sets)

IMPORTANT! ProcartaPlex™ pre-mixed panel and simplex kits can be mixed together for enhanced flexibility. Ensure that the bead regions from your ProcartaPlex™ panels and simplex kits do not overlap. Some analytes use the same bead region and cannot be combined together in one multiplex assay. Please check the compatibility of our analytes using our online panel configurator or contact our technical support.

Both simplex and pre-mixed kits are supplied with Antibody Magnetic Beads at working concentration that require 12.5 µL beads per well or 5 mLs per 384-wells. For running a partial plate adjust the volume accordingly.

1. Vortex each of the Antibody Magnetic Bead sets for 30 sec and add 5 mL of each bead set for 384-wells to an appropriate sized tube. After all of the bead sets are added, vortex the tube for another 30 sec.
2. Add the bead mix to a disposable reservoir and add the appropriate volume of Antibody Magnetic Beads to each well of the 384-Well Filter Plate using the table below:

# of different Magnetic Bead sets to be mixed	Amount added to each well in µL
2	25 µL
3	37.5 µL
4	50 µL
5	62.5 µL
6	75 µL

If more than 6 bead sets are to be mixed proceed to “Step 4. Wash antibody magnetic beads” on page 12 and repeat “Step 3b. Add the antibody magnetic beads (2 or more beads sets)” on page 12 until all Antibody Magnetic Beads have been added and washed.

Step 4. Wash antibody magnetic beads

1. Carefully remove the Plate Seal to avoid splashing the plate contents.
2. Remove the liquid in the wells by aspiration using Vacuum Filtration Manifold. Blot the bottom of the plate after filtration.
3. Add 50 µL of Wash Buffer (1X) into each well.
4. Remove the liquid in the wells by aspiration using Vacuum Filtration Manifold. Blot the bottom of the plate after filtration.



Step 5. Add sample type-specific buffer, samples, standards and blanks and incubate

1. For cell culture supernatant samples add 12.5 µL standards or samples into dedicated wells.
2. For wells designated as blanks add 12.5 µL of cell culture medium.
3. Seal the plate with the provided Plate Seal. Cover the plate with the Black Microplate Lid and shake at 500 rpm for 60 to 120 min at room temperature (RT).
4. Alternatively, the 384-Well Plate can be incubated overnight. Shake the 384-Well Plate 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 6. Wash the 384-Well plate

Wash plate for a total of three times using “Step 4. Wash antibody magnetic beads” on page 12.

Step 7. Prepare 1X detection antibody mixture

1. Prepare fresh prior to use detection antibodies. If intending overnight incubation do not prepare detection antibodies at this point.
2. Detection antibodies for Custom Panels are provided at a 1X concentration and do not require dilution.
3. For simplex and pre-mixed panels, detection antibody is provided at 50X concentration. Add 60 µL of each detection antibody concentrate to the mixing bottle and bring volume of the mixing bottle to a total of 3 mL using detection antibody diluent. Tables below are an example for 384 and 192-wells.

Table 1 Example for using 384-wells

# of Vials of Detection Antibody	Total Volume of Mixed 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

Table 2 Example for using 192-wells

# of Vials of Detection Antibody	Total Volume of Mixed Detection Antibody	Volume of Diluent to Add
1	30 µL	1470 µL
2	60 µL	1440 µL
3	90 µL	1410 µL
4	120 µL	1380 µL

Step 8. Add detection antibody mixture and incubate

1. Add 6.25 µL of Detection Antibody Mixture (1X) to each well.
2. Seal the plate with a new Plate Seal, cover plate with Black Microplate Lid and incubate 30 min on a plate shaker at RT at 500 rpm.



- Step 9. Wash the 384-Well plate** Wash plate for a total of three times using “Step 4. Wash antibody magnetic beads” on page 12.
- Step 10. Add SAPE and incubate**
1. Add 12.5 µL of SAPE solution to each well.
 2. Seal the plate with a new Plate Seal, cover plate with Black Microplate Lid and incubate 30 min on a plate shaker at RT at 500 rpm.
- Step 11. Wash the 384-Well plate** Wash plate for a total of three times using “Step 4. Wash antibody magnetic beads” on page 12.
- Step 12. Prepare the 384-Well plate for analysis on a Luminex™ instrument**
1. Add 50 µL of Reading Buffer into each well.
 2. Seal the plate with a new Plate Seal, cover plate with Black Microplate Lid and incubate 5 min on a plate shaker at RT at 500 rpm.
 3. Remove Plate Seal and run the plate on a Luminex™ Instrument.

Setup of the Luminex™ instruments

Sample Size	DD Gate	Timeout	Bead Event/Bead Region
50 µL	5,000 - 25,000	60 sec	50-100

Verify the probe height for each plate before reading. Failure to adjust the probe height can cause damage to the instrument. The Luminex™ system allows for calibration of Low and High RP1 target values. We recommend RP1 Low target value settings for ProcortaPlex™ Immunoassays.

Please refer to the Certificate of Analysis provided with the kit for bead region and analyte associations when entering the information into the Luminex™ Acquisition Software.

Note: If there is a malfunction of the Luminex™ Instrument or software during the run, the 384-Well Plate can be re-read. Remove the 384-Well Plate from the instrument and remove the liquid in the wells by aspiration using the Vacuum Filtration Manifold. Blot the bottom of the plate after filtration. Resuspend the beads in 50 µL of Reading Buffer, seal the 384-Well Plate with a new Plate Seal 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 wells.



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. For download information visit our website under Life Sciences and Support Centers → Luminex™ Assays or contact our technical support.

Troubleshooting

Observation	Probable Cause	Recommend Solution
Low Flow Rate	Samples/beads are stuck in flow cell	Remove the 384-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 µL 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 384-well plate in the dark.



Observation	Probable Cause	Recommend Solution
Low bead count	Samples causing the instrument to clog	Remove the 384-Well Filter 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.

Blank plate layout

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A																								
B																								
C																								
D																								
E																								
F																								
G																								
H																								
I																								
J																								
K																								
L																								
M																								
N																								
O																								
P																								



Safety



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
 - Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, see the “Documentation and Support” section in this document.
-

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if needed) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



WARNING! HAZARDOUS WASTE (from instruments). Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



WARNING! 4L Reagent and Waste Bottle Safety. Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position.

Biological hazard safety



WARNING! Potential Biohazard. Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:
<https://www.cdc.gov/labs/pdf/CDC-BiosafetymicrobiologicalBiomedicalLaboratories-2009-P.pdf>
 - World Health Organization, *Laboratory Biosafety Manual*, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:
www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf
-

Documentation and support

Customer and technical support

Visit [thermofisher.com/support](http://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

