



## **Human Cytokine / Chemokine Panel II**

### **96 Well Plate Assay**

**Cat. # HCYP2MAG-62K,  
HCP2MAG-62K-PX23,  
HCP2MAG62KPX23BK**

# MILLIPLEX® MAP

## HUMAN CYTOKINE / CHEMOKINE PANEL II 96 Well Plate Assay

# HCYP2MAG-62K,  
# HCP2MAG-62K-PX23,  
# HCP2MAG62KPX23BK

<b><u>TABLE OF CONTENTS</u></b>	<b><u>PAGE</u></b>
Introduction	2
Principle	3
Storage Conditions Upon Receipt	4
Reagents Supplied	4
Materials Required But Not Provided	6
Safety Precautions	6
Technical Guidelines	8
Sample Collection And Storage	9
Preparation of Reagents for Immunoassay	10
Immunoassay Procedure	12
Plate Washing	15
Equipment Settings	15
Quality Controls	16
Assay Characteristics	17
Troubleshooting Guide	20
Replacement Reagents	23
Ordering Information	24
Well Map	25

**For Research Use Only. Not for Use in Diagnostic Procedures.**

By purchasing this product, which contains fluorescently labeled microsphere beads authorized by Luminex® Corporation (“Luminex®”), you, the customer, acquire the right under Luminex®’s patent rights, if any, to use this product or any portion of this product, including without limitation the microsphere beads contained herein, only with Luminex®’s laser based fluorescent analytical test instrumentation marketed under the name of Luminex® 100™ IS, 200™, HTS, FLEXMAP 3D®, MAGPIX®..

# Human Cytokine / Chemokine Panel II

## INTRODUCTION

“Cytokine” is a general term used for a diverse group of soluble proteins and peptides which act as regulators under both normal and pathological conditions to modulate the functional activities of individual cells and tissues. These proteins also mediate direct interactions between cells and regulate processes taking place in the extracellular environment. Cytokines differ from hormones in that they act on a wider spectrum of target cells. Also, unlike hormones, they are not produced by specialized cells which are organized in specialized glands. The cytokine group of proteins includes lymphokines, interferons, colony stimulating factors and chemokines. Cytokine and chemokine research plays a significant role in achieving a deeper understanding of the immune system and its multi-faceted response to most antigens, as well as disease states such as inflammatory disease, allergic reactions, IBD, sepsis, and cancer.

MILLIPLEX<sup>®</sup> MAP offers the broadest selection of analytes across a wide range of disease states and species. Once the analytes of interest have been identified, you can rely on the quality that we build into each kit to produce results you can trust. In addition to the assay characteristics listed in the protocol, other performance criteria evaluated during the validation process include: cross-reactivity, dilution linearity, kit stability, and sample behavior (e.g. detectability and stability).

Each MILLIPLEX<sup>®</sup> MAP panel and kit includes:

- Quality controls (QCs) provided to qualify assay performance
- Comparison of standard (calibrator) and QC lots to a reference lot to ensure lot-to-lot consistency
- Optimized serum matrix to mimic native analyte environment
- Detection antibody cocktails designed to yield consistent analyte profiles within panel

In addition each panel and kit meets stringent manufacturing criteria to ensure batch-to-batch reproducibility. The MILLIPLEX<sup>®</sup> MAP Human Cytokine / Chemokine Panel II Magnetic Bead Panel thus enables you to focus on the therapeutic potential of cytokines as well as the modulation of cytokine expression. Coupled with the Luminex<sup>®</sup> xMAP<sup>®</sup> platform in a magnetic bead format, you receive the advantage of ideal speed and sensitivity, allowing quantitative multiplex detection of dozens of analytes simultaneously, which can dramatically improve productivity.

EMD Millipore's MILLIPLEX<sup>®</sup> MAP Human Cytokine / Chemokine Panel II Magnetic Bead Panel is part of the most versatile system available for cytokine and chemokine research. From our single to multiplex biomarker solutions, we partner with you to design, develop, analytically validate and build the most comprehensive library available for protein detection and quantitation.

- MILLIPLEX<sup>®</sup> MAP offers you:
  - The ability to select a 23-plex or premixed option
  - The ability to choose any combination of analytes from our panel of 23 analytes to design a custom kit that better meets your needs.
  - A convenient “all-in-one” box format that gives you the assurance that you will have all the necessary reagents you need to run your assay.

EMD Millipore's MILLIPLEX<sup>®</sup> MAP Human Cytokine / Chemokine Panel II is to be used for the simultaneous quantification of any or all of the following analytes in serum, plasma, tissue/cell lysate and culture supernatant samples: Eotaxin-2/CCL24, MCP-2, BCA-1/CXCL13, MCP-4, I-309/CCL1, IL-16, TARC/CCL17, 6Ckine/CCL21/Exodus-2, Eotaxin-3/CCL26, LIF, TPO, SCF, TSLP, IL-33/NF-HEV (mature), IL-20, IL-21, IL-23, TRAIL/TNFSF10, CTACK/CCL27, SDF-1 $\alpha$ + $\beta$ /CXCL12, ENA-78/CXCL5, MIP-1 $\delta$ /MIP-5/CCL15, IL-28A/IFN $\lambda$ 2.

***For Research Use Only. Not for Use in Diagnostic Procedures.***

***Please read entire protocol before use.***

***It is important to use same assay incubation conditions throughout your study.***

## **PRINCIPLE**

MILLIPLEX<sup>®</sup> MAP is based on the Luminex<sup>®</sup> xMAP<sup>®</sup> technology — one of the fastest growing and most respected multiplex technologies offering applications throughout the life-sciences and capable of performing a variety of bioassays including immunoassays on the surface of fluorescent-coded magnetic bead (MagPlex<sup>®</sup>-C and non-magnetic bead (MicroPlex<sup>®</sup>) microspheres..

- Luminex<sup>®</sup> uses proprietary techniques to internally color-code microspheres with multiple fluorescent dyes. Through precise concentrations of these dyes, distinctly colored bead sets of 500 5.6  $\mu$ m non-magnetic or 80 6.45  $\mu$ m magnetic polystyrene microspheres can be created, each of which is coated with a specific capture antibody.
- After an analyte from a test sample is captured by the bead, a biotinylated detection antibody is introduced.
- The reaction mixture is then incubated with Streptavidin-PE conjugate, the reporter molecule, to complete the reaction on the surface of each microsphere.
- EMD Millipore provides three Luminex<sup>®</sup> instruments to acquire and analyze data using two detection methods:
  - The Luminex<sup>®</sup> analyzers Luminex<sup>®</sup> 200<sup>™</sup> and FLEXMAP 3D<sup>®</sup>, flow cytometry-based instruments that integrate key xMAP<sup>®</sup> detection components, such as lasers, optics, advanced fluidics and high-speed digital signal processors.
  - The Luminex<sup>®</sup> analyzer (MAGPIX<sup>®</sup>), a CCD-based instrument that integrates key xMAP<sup>®</sup> capture and detection components with the speed and efficiency of magnetic beads.
- Each individual microsphere is identified and the result of its bioassay is quantified based on fluorescent reporter signals. EMD Millipore combines the streamlined data acquisition power of Luminex<sup>®</sup> xPONENT<sup>®</sup> acquisition software with sophisticated analysis capabilities of the new MILLIPLEX<sup>®</sup> Analyst 5.1, integrating data acquisition and analysis seamlessly with all Luminex<sup>®</sup> instruments.

The capability of adding multiple conjugated beads to each sample results in the ability to obtain multiple results from each sample. Open-architecture xMAP<sup>®</sup> technology enables multiplexing of many types of bioassays reducing time, labor and costs over traditional methods.

## STORAGE CONDITIONS UPON RECEIPT

- Recommended storage for kit components is 2 - 8°C.
- For long-term storage, freeze reconstituted standards and controls at ≤ -20°C. Avoid multiple (>2) freeze thaw cycles.
- **DO NOT FREEZE Antibody-Immobilized Beads, Detection Antibody, and Streptavidin-Phycoerythrin.**

## REAGENTS SUPPLIED

**Note: Store all reagents at 2 – 8°C**

REAGENTS SUPPLIED	CATALOG NUMBER	VOLUME	QUANTITY
Human Cytokine / Chemokine Panel II Standard	MXH8062	Lyophilized	1 vial
Human Cytokine Panel II Quality Controls 1 and 2	MXH6062	Lyophilized	2 vials
Serum Matrix Note: Contains 0.08% Sodium Azide	MXHSM-2	Lyophilized	1 vial
Set of one 96-Well Plate with 2 Sealers	-----		1 plate 2 sealers
Assay Buffer	L-AB	30 mL	1 bottle
Bead Diluent	LBD	3,5mL	1bottle
10X Wash Buffer Note: Contains 0.05% Proclin	L-WB	60 mL	1 bottle
Human Cytokine Panel II Detection Antibodies	MXH1062	3.2 mL	1 bottle
Streptavidin-Phycoerythrin	L-SAPE4	3.2 mL	1 bottle
Mixing Bottle (not provided with premixed panel)	-----	-----	1 bottle

## Human Cytokine / Chemokine Panel II Antibody-Immobilized Premixed Magnetic Beads:

Premixed 23-plex Beads	HP2PMX23-MAG	3.5 mL	1 bottle
------------------------	--------------	--------	----------

## REAGENTS SUPPLIED (continued)

### Human Cytokine / Chemokine Panel II Antibody-Immobilized Beads:

Bead/Analyte Name	Luminex® Bead Region	Customizable 23 Analytes (50X concentration, 90µL)		23-Plex Premixed Beads
		Available	Cat. #	
Anti-Human Eotaxin-2 Bead	12	✓	HETXN2-MAG	✓
Anti-Human MCP-2 Bead	13	✓	HMCP2-MAG	✓
Anti-Human BCA-1 Bead	15	✓	HBCA1-MAG	✓
Anti-Human MCP-4 Bead	18	✓	HMCP4-MAG	✓
Anti-Human I-309 Bead	19	✓	HI309-MAG	✓
Anti-Human IL-16 Bead	21	✓	HIL16-MAG	✓
Anti-Human TARC bead	26	✓	HTARC-MAG	✓
Anti-Human 6CKine Bead	28	✓	H6CKINE-MAG	✓
Anti-Human Eotaxin-3 Bead	30	✓	HETXN3-MAG	✓
Anti-Human LIF Bead	34	✓	HLIF-MAG	✓
Anti-Human TPO Bead	36	✓	HTPO-MAG	✓
Anti-Human SCF Bead	38	✓	HCYSCF-MAG	✓
Anti-Human TSLP bead	43	✓	HTSLP-MAG	✓
Anti-Human IL-33 bead	45	✓	HIL33-MAG	✓
Anti-Human IL-20 bead	51	✓	HIL20--MAG	✓
Anti-Human IL-21 bead	52	✓	HIL21-MAG	✓
Anti-Human IL-23 bead	54	✓	HIL23-MAG	✓
Anti-Human TRAIL bead	56	✓	HCYTRAIL-MAG	✓
Anti-Human CTACK bead	62	✓	HCTACK-MAG	✓
Anti-Human SDF-1α+β bead	64	✓	HSDF1AB-MAG	✓
Anti-Human ENA-78 bead	66	✓	HENA78-MAG	✓
Anti-Human MIP-1d bead	76	✓	HMIP1D-MAG	✓
Anti-Human IL-28A bead	77	✓	HIL28A-MAG	✓

## **MATERIALS REQUIRED BUT NOT PROVIDED**

### Reagents

1. Luminex® Sheath Fluid (EMD Millipore Catalog # SHEATHFLUID) or Luminex® Drive Fluid (EMD Millipore Catalog # MPXDF-4PK)

### Instrumentation / Materials

1. Adjustable Pipettes with Tips capable of delivering 25 µL to 1000 µL
2. Multichannel Pipettes capable of delivering 5 µL to 50 µL or 25 µL to 200 µL
3. Reagent Reservoirs
4. Polypropylene Microfuge Tubes
5. Rubber Bands
6. Aluminum Foil
7. Absorbent Pads
8. Laboratory Vortex Mixer
9. Sonicator (Branson Ultrasonic Cleaner Model #B200 or equivalent)
10. Titer Plate Shaker (VWR® Microplate Shaker Cat # 12620-926 or equivalent)
11. Luminex® 200™, HTS, FLEXMAP 3D®, or MAGPIX® with xPONENT® software by Luminex® Corporation
12. Automatic Plate Washer for magnetic beads (BioTek® 405 LS and 405 TS, EMD Millipore Catalog # 40-094, # 40-095, # 40-096, # 40-097 or equivalent) or Handheld Magnetic Separation Block (EMD Millipore Catalog # 40-285 or equivalent).








Note: If a plate washer or handheld magnetic separation block for magnetic beads is not available, one can use a microtiter filter plate (EMD Millipore Catalog # MX-PLATE) to run the assay using a Vacuum Filtration Unit (EMD Millipore Vacuum Manifold Catalog # MSVMHTS00 or equivalent with EMD Millipore Vacuum Pump Catalog # WP6111560 or equivalent).

## **SAFETY PRECAUTIONS**

- All blood components and biological materials should be handled as potentially hazardous. Follow universal precautions as established by the Centers for Disease Control and Prevention and by the Occupational Safety and Health Administration when handling and disposing of infectious agents.
- Sodium Azide or Proclin has been added to some reagents as a preservative. Although the concentrations are low, Sodium Azide and Proclin may react with lead and copper plumbing to form highly explosive metal azides. Dispose of unused contents and waste in accordance with international, federal, state, and local regulations.

**Note: See Full Labels of Hazardous components on next page.**

Full Labels of Hazardous components:

Ingredient, Cat #		Full Label	
Streptavidin-Phycoerythrin	L-SAPE4		<b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
10X Wash Buffer	L-WB		<b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.
Human Cytokine/Chemokine Panel II Detection Antibody	MXH1062		<b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
Human Cytokine/Chemokine Panel II Quality Controls 1 & 2	MXH6062	 	<b>Danger.</b> Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/attention.
Human Cytokine/Chemokine Panel II Standard	MXH8062	 	<b>Danger.</b> Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/attention.
Serum Matrix	MXHSM-2	No Symbol Required	<b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.



## TECHNICAL GUIDELINES

To obtain reliable and reproducible results, the operator should carefully read this entire manual and fully understand all aspects of each assay step before running the assay. The following notes should be reviewed and understood before the assay is set-up.

- FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.
- Do not use beyond the expiration date on the label.
- Do not mix or substitute reagents with those from other lots or sources.
- The Antibody-Immobilized Beads are light sensitive and must be protected from light at all times. Cover the assay plate containing beads with opaque plate lid or aluminum foil during all incubation steps.
- It is important to allow all reagents to warm to room temperature (20-25°C) before use in the assay.
- Incomplete washing can adversely affect the assay outcome. All washing must be performed with the Wash Buffer provided.
- The standards prepared by serial dilution must be used within 1 hour of preparation. Discard any unused standards except the standard stock which may be stored at  $\leq -20^{\circ}\text{C}$  for 1 month and at  $\leq -80^{\circ}\text{C}$  for greater than one month.
- If samples fall outside the dynamic range of the assay, further dilute the samples with the appropriate diluent and repeat the assay.
- Any unused mixed Antibody-Immobilized Beads may be stored in the Bead Mix bottle at 2-8°C for up to one month.
- During the preparation of the standard curve, make certain to mix the higher concentration well before making the next dilution. Use a new tip with each dilution.
- The plate should be read immediately after the assay is finished. If, however, the plate cannot be read immediately, seal the plate, cover with aluminum foil or an opaque lid, and store the plate at 2-8°C for up to 24 hours. Prior to reading, agitate the plate on the plate shaker at room temperature for 10 minutes. Delay in reading a plate may result in decreased sensitivity for some analytes.
- The titer plate shaker should be set at a speed to provide maximum orbital mixing without splashing of liquid outside the wells. For the recommended plate shaker, this would be a setting of 5-7, which is approximately 500-800 rpm.
- Ensure that the needle probe is clean. This may be achieved by sonication and/or alcohol flushes.
- When reading the assay on Luminex® 200™, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended EMD Millipore filter plates using 3 alignment discs. When reading the assay on MAGPIX®, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended EMD Millipore filter plates using 2 alignment discs. When reading the assay on FLEXMAP 3D®, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 1 alignment disc.

For FLEXMAP 3D® when using the solid plate in the kit, the final resuspension should be with 150  $\mu\text{L}$  Sheath Fluid in each well and 75  $\mu\text{L}$  should be aspirated.

## TECHNICAL GUIDELINES (continued)

- For cell culture supernatants or tissue extraction, use the culture or extraction medium as the matrix solution in blank, standard curve and controls. If samples are diluted in assay buffer, use the assay buffer as matrix.
- For serum/plasma sample that require a dilution instead of “Neat”, use the Serum Matrix provided in the kit as the diluent.
- For cell/tissue homogenate, the final cell or tissue homogenate should be prepared in a buffer that has a neutral pH, contains minimal detergents or strong denaturing detergents, and has an ionic strength close to physiological concentration. Avoid debris, lipids, and cell/tissue chunks. Centrifuge samples before use.
- Vortex all reagents well before adding to plate.

## SAMPLE COLLECTION AND STORAGE

### A. Preparation of Serum Samples:

- Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000xg. Remove serum and assay immediately or aliquot and store samples at  $\leq -20^{\circ}\text{C}$ .
- Avoid multiple (>2) freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.
- Neat serum samples are used. If further dilution is required, use Serum Matrix as the diluent. Note: Serum Matrix volume provided in the kit may be insufficient. Additional matrix may be obtained from EMD Millipore.

### B. Preparation of Plasma Samples:

- Plasma collection using EDTA as an anti-coagulant is recommended. Centrifuge for 10 minutes at 1000xg within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at  $\leq -20^{\circ}\text{C}$ .
- Avoid multiple (> 2) freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.
- Neat Plasma samples are used. If further dilution is required, use Serum Matrix as the diluent. Note: Serum Matrix volume provided in the kit may be insufficient. Additional matrix may be obtained from EMD Millipore.

### C. Preparation of Tissue Culture Supernatant:

- Centrifuge the sample to remove debris and assay immediately or aliquot and store samples at  $\leq -20^{\circ}\text{C}$ .
- Avoid multiple (> 2) freeze/thaw cycles.
- Tissue culture supernatant may require a dilution with an appropriate control medium prior to assay.

## SAMPLE COLLECTION AND STORAGE (continued)

### NOTE:

- A maximum of 25  $\mu$ L per well of neat or diluted serum or plasma can be used. Tissue culture or other media may also be used.
- All samples must be stored in polypropylene tubes. **DO NOT STORE SAMPLES IN GLASS.**
- Avoid debris, lipids and cells when using samples with gross hemolysis or lipemia.
- Care must be taken when using heparin as an anticoagulant since an excess of heparin will provide falsely high values. Use no more than 10 IU heparin per mL of blood collected.

## PREPARATION OF REAGENTS FOR IMMUNOASSAY

### A. Preparation of Antibody-Immobilized Beads

If premixed beads are used, sonicate the premixed bead bottle 30 seconds and then vortex for 1 minute before use.

For individual vials of beads, sonicate each antibody-bead vial for 30 seconds; vortex for 1 minute. Add 60  $\mu$ L from each antibody bead vial to the Mixing Bottle and bring final volume to 3.0 mL with Bead Diluent. Vortex the mixed beads well. Unused portion may be stored at 2-8°C for up to one month. (Note: Due to the composition of magnetic beads, you may notice a slight color in the bead solution. This does not affect the performance of the beads or the kit.)

Example 1: When using 5 cytokine antibody-immobilized beads, add 60  $\mu$ L from each of the 5 bead sets to the Mixing Bottle. Then add 2.7mL Bead Diluent.

Example 2: When using 9 cytokine antibody-immobilized beads, add 60  $\mu$ L from each of the 9 bead sets to the Mixing Bottle. Then add 2.46 mL Bead Diluent.

### B. Preparation of Quality Controls

Before use, reconstitute Quality Control 1 and Quality Control 2 with 250  $\mu$ L deionized water. Invert the vial several times to mix and vortex. Allow the vial to sit for 5-10 minutes. Unused portion may be stored at  $\leq$  -20°C for up to one month.

### C. Preparation of Wash Buffer

Bring the 10X Wash Buffer to room temperature and mix to bring all salts into solution. Dilute 60 mL of 10X Wash Buffer with 540 mL deionized water. Store unused portion at 2-8°C for up to one month.

## PREPARATION OF REAGENTS FOR IMMUNOASSAY (continued)

### D. Preparation of Serum Matrix

**This step is required for serum or plasma samples only.**

Add 1.0 mL Assay Buffer to the bottle containing lyophilized Serum Matrix. Mix well. Allow at least 10 minutes for complete reconstitution. Leftover reconstituted Serum Matrix should be stored at  $\leq -20^{\circ}\text{C}$  for up to one month.

### E. Preparation of Human Cytokine Panel II Standard

1.) Prior to use, reconstitute the Human Cytokine Panel II Standard with 250  $\mu\text{L}$  deionized water. Invert the vial several times to mix. Vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes. This will be used as the Std 6; the unused portion may be stored at  $\leq -20^{\circ}\text{C}$  for up to one month.

#### 2). Preparation of Working Standards

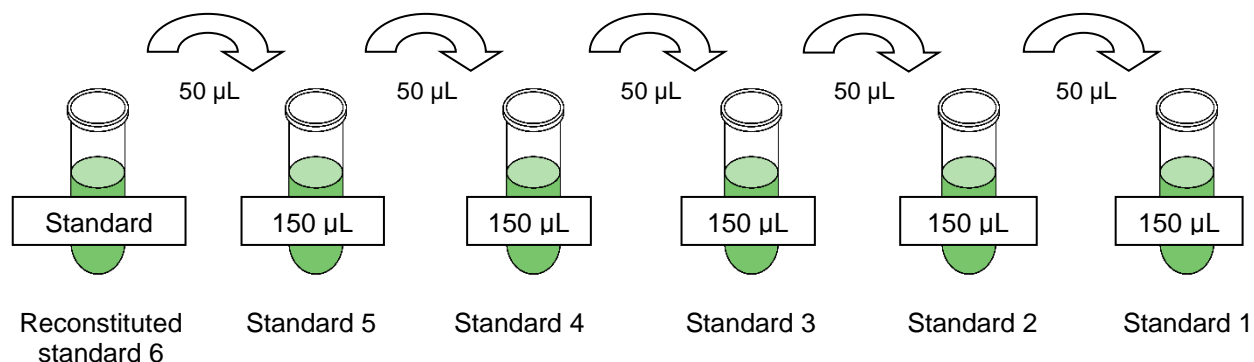
Label five polypropylene microfuge tubes Std 5, Std 4, Std 3, Std 2 and Std 1. Add 150  $\mu\text{L}$  of Assay Buffer to each of the five tubes. Prepare serial dilutions by adding 50  $\mu\text{L}$  of the reconstituted Standard 6 to the Std 5 tube, mix well and transfer 50  $\mu\text{L}$  of the Standard 5 to the Std 4 tube, mix well and transfer 50  $\mu\text{L}$  of the Standard 4 to the Std 3 tube, mix well and transfer 50  $\mu\text{L}$  of the Standard 3 to Std 2 tube, mix well and transfer 50  $\mu\text{L}$  of the Standard 2 to the Std 1 tube and mix well. The Standard 0 (Background) will be Assay Buffer.

Standard	Volume of Deionized Water to Add	Volume of Standard to Add
Original (Std 6)	250 $\mu\text{L}$	0

Standard	Volume of Assay Buffer to Add	Volume of Standard to Add
Standard 5	150 $\mu\text{L}$	50 $\mu\text{L}$ of Standard 6
Standard 4	150 $\mu\text{L}$	50 $\mu\text{L}$ of Standard 5
Standard 3	150 $\mu\text{L}$	50 $\mu\text{L}$ of Standard 4
Standard 2	150 $\mu\text{L}$	50 $\mu\text{L}$ of Standard 3
Standard 1	150 $\mu\text{L}$	50 $\mu\text{L}$ of Standard 2

## PREPARATION OF REAGENTS FOR IMMUNOASSAY (continued)

### Preparation of Standards



After serial dilutions, the tubes should have the following concentrations for constructing standard curves.

Standard Tube #	BCA-1, TARC (pg/mL)	I-309 (pg/mL)	MCP-2, CTACK (pg/mL)	MCP-4, Eotaxin-2, TRAIL, SCF, TSLP, IL-28A, IL-16 (pg/mL)	ENA-78, 6Ckine, LIF, IL-21, IL-33 (pg/mL)	MIP-1 $\delta$ , Eotaxin-3, IL-23, TPO, IL-20 (pg/mL)	SDF-1 $\alpha$ + $\beta$ , (pg/mL)
1	1.0	2.0	4.9	9.8	19.5	48.8	97.7
2	3.9	7.8	19.5	39.1	78.1	195.3	390.6
3	15.6	31.3	78.1	156.3	312.5	781.3	1562.5
4	62.5	125	312.5	625	1250	3125	6250
5	250	500	1250	2500	5000	12500	25,000
6	1000	2000	5000	10,000	20,000	50,000	100,000

### IMMUNOASSAY PROCEDURE

- Prior to beginning this assay, it is imperative to read this protocol completely and to thoroughly understand the Technical Guidelines.
- Allow all reagents to warm to room temperature (20-25°C) before use in the assay.
- Diagram the placement of Standards [0 (Background), Std1, Std2, Std3, Std4, Std5, and Std6], Controls 1 and 2, and Samples on Well Map Worksheet in a vertical configuration. (Note: Most instruments will only read the 96-well plate vertically by default.) It is recommended to run the assay in duplicate.
- If using a filter plate, set the filter plate on a plate holder at all times during reagent dispensing and incubation steps so that the bottom of the plate does not touch any surface.

## IMMUNOASSAY PROCEDURE (continued)

1. Pre-wet the plate by pipetting 200  $\mu\text{L}$  of Wash Buffer into each well of the Plate. Seal and shake on a plate shaker for 10 minutes at room temperature (20-25°C).
2. Decant Wash Buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times.
3. Add 25  $\mu\text{L}$  of each Standard or Control into the appropriate wells. Assay Buffer should be used for 0 pg/mL standard (Background).
4. Add 25  $\mu\text{L}$  of Assay Buffer to the sample wells.
5. Add 25  $\mu\text{L}$  of appropriate matrix solution to the background, standards, and control wells. When assaying serum or plasma, use the Serum Matrix provided in the kit. When assaying tissue culture or other supernatant, use proper control culture medium as the matrix solution.
6. Add 25  $\mu\text{L}$  of Sample (neat) into the appropriate wells.
7. Vortex Mixing Bottle and add 25  $\mu\text{L}$  of the Mixed or Premixed Beads to each well. (Note: During addition of Beads, shake bead bottle intermittently to avoid settling.)
8. Seal the plate with a plate sealer. Wrap the plate with foil and incubate with agitation on a plate shaker overnight at 4°C or 2 hours at room temperature (20-25°C). *An overnight incubation (16-18 hr) may improve assay sensitivity for some analytes.*
9. Gently remove well contents and wash plate 2 times following instructions listed in the **PLATE WASHING** section.
10. Add 25  $\mu\text{L}$  of Detection Antibodies into each well. (Note: Allow the Detection Antibodies to warm to room temperature prior to addition.)

Add 200  $\mu\text{L}$  Wash Buffer per well



Shake 10 min, RT

Decant

- Add 25  $\mu\text{L}$  Standard or Control to appropriate wells
- Add 25  $\mu\text{L}$  Assay Buffer to background and sample wells



- Add 25  $\mu\text{L}$  appropriate matrix to background, standards, and control wells
- Add 25  $\mu\text{L}$  Samples to sample wells
- Add 25  $\mu\text{L}$  Beads to each well



Incubate overnight (16-18 hr) at 4°C or 2 hours at RT with shaking



Remove well contents and wash 2X with 200  $\mu\text{L}$  Wash Buffer

Add 25  $\mu\text{L}$  Detection Antibodies per well

11. Seal, cover with foil and incubate with agitation on a plate shaker for 1 hour at room temperature (20-25°C). **DO NOT ASPIRATE AFTER INCUBATION.**

12. Add 25  $\mu$ L Streptavidin-Phycoerythrin to each well containing the 25  $\mu$ L of Detection Antibodies.

13. Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25°C).

14. Gently remove well contents and wash plate 2 times following instructions listed in the **PLATE WASHING** section.

15. Add 150  $\mu$ L of Sheath Fluid (or Drive Fluid if using MAGPIX<sup>®</sup>) to all wells. Resuspend the beads on a plate shaker for 5 minutes.

16. Run plate on Luminex<sup>®</sup> 200™, HTS, FLEXMAP 3D<sup>®</sup> or MAGPIX<sup>®</sup> with xPONENT<sup>®</sup> software

17. Save and analyze the Median Fluorescent Intensity (MFI) data using a 5-parameter logistic or spline curve-fitting method for calculating cytokine/chemokines concentrations in samples. (Note: For diluted samples, multiply the calculated concentration by the dilution factor.)

Incubate 1 hour at RT.  
Do not aspirate.



Add 25  $\mu$ L Streptavidin-Phycoerythrin per well



Incubate for 30 minutes at RT



Remove well contents and wash 2X with 200  $\mu$ L Wash Buffer

Add 150  $\mu$ L Sheath Fluid or Drive Fluid per well

Read on Luminex<sup>®</sup> (100  $\mu$ L, 50 beads per bead set)

## PLATE WASHING

### 1.) Solid Plate

**If using a solid plate, use either a handheld magnet or magnetic plate washer.**

- A.) Handheld magnet (EMD Millipore Catalog # 40-285) - Rest plate on magnet for 60 seconds to allow complete settling of magnetic beads. Remove well contents by gently decanting the plate in an appropriate waste receptacle and gently tapping on absorbent pads to remove residual liquid. Wash plate with 200  $\mu$ L of Wash Buffer by removing plate from magnet, adding Wash Buffer, shaking for 30 seconds, reattaching to magnet, letting beads settle for 60 seconds and removing well contents as previously described after each wash. Repeat wash steps as recommended in Assay Procedure.
- B.) Magnetic plate washer (EMD Millipore Catalog # 40-094, # 40-095, # 40-096 and # 40-097) - Please refer to specific automatic plate washer manual for appropriate equipment settings. Please note that after the final aspiration, there will be approximately 25  $\mu$ L of residual wash buffer in each well. This is expected when using the BioTek plate washer and this volume does not need to be aspirated from the plate.

**If using an automatic plate washer other than BioTek® 405 LS or 405 TS, please refer to the manufacturer's recommendations for programming instructions.**

### 2.) Filter Plate (EMD Millipore Catalog # MX-PLATE)

If using a filter plate, use a vacuum filtration manifold to remove well contents. Wash plate with 200  $\mu$ L/well of Wash Buffer, removing Wash Buffer by vacuum filtration after each wash. Repeat wash steps as recommended in the Assay Procedure.

## EQUIPMENT SETTINGS

Luminex® 200™, HTS, FLEXMAP 3D®, and MAGPIX® with xPONENT® software:

These specifications are for the Luminex® 200™, Luminex® HTS, Luminex® FLEXMAP 3D®, and Luminex® MAGPIX® with xPONENT® software. Luminex® instruments with other software (e.g. MasterPlex®, StarStation, LiquiChip, Bio-Plex® Manager™, LABScan™ 100) would need to follow instrument instructions for gate settings and additional specifications from the vendors for reading Luminex® magnetic beads.

For magnetic bead assays, the Luminex® 200™ and HTS instruments must be calibrated with the xPONENT® 3.1 compatible Calibration Kit (EMD Millipore Catalog # 40-275) and performance verified with the Performance Verification Kit (EMD Millipore Catalog # 40-276). The Luminex® FLEXMAP 3D® instrument must be calibrated with the FLEXMAP 3D® Calibrator Kit (EMD Millipore Catalog # 40-028) and performance verified with the FLEXMAP 3D® Performance Verification Kit (EMD Millipore Catalog # 40-029). The Luminex® MAGPIX® instrument must be calibrated with the MAGPIX® Calibration Kit (EMD Millipore Catalog # 40-049) and performance verified with the MAGPIX® Performance Verification Kit (EMD Millipore Catalog # 40-050).

**NOTE: When setting up a Protocol using the xPONENT® software, you must select MagPlex® as the Bead Type in the Acquisition settings.**

**NOTE: These assays cannot be run on any instruments using Luminex® IS 2.3 or Luminex® 1.7 software.**



## EQUIPMENT SETTINGS (continued)

The Luminex® probe height must be adjusted to the plate provided in the kit. Please use Catalog #MAG-PLATE, if additional plates are required for this purpose.

Events:	50, per bead	
Sample Size:	100 µL	
Gate Settings	8,000 to 15,000	
Time Out	60 seconds	
Bead Set:	Customizable 23-Plex Beads	23-Plex Premix Beads
Eotaxin-2	12	12
MCP-2	13	13
BCA-1	15	15
MCP-4	18	18
I-309	19	19
IL-16	21	21
TARC	26	26
6CKine	28	28
Eotaxin-3	30	30
LIF	34	34
TPO	36	36
SCF	38	38
TSLP	43	43
IL-33	45	45
IL-20	51	51
IL-21	52	52
IL-23	54	54
TRAIL	56	56
CTACK	62	62
SDF-1α+β	64	64
ENA-78	66	66
MIP-1δ	76	76
IL-28A	77	77

## QUALITY CONTROLS

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert or can be located at the EMD Millipore website [www.emdmillipore.com](http://www.emdmillipore.com) using the catalog number as the keyword.

## ASSAY CHARACTERISTICS

### Cross-Reactivity

There was no or negligible cross-reactivity between the antibodies and any of the other analytes in this panel.

### Assay Sensitivities (minimum detectable concentrations, pg/mL)

Minimum Detectable Concentration (MinDC) is calculated using MILLIPLEX® Analyst 5.1. It measures the true limits of detection for an assay by mathematically determining what the empirical MinDC would be if an infinite number of standard concentrations were run for the assay under the same conditions.

Cytokines	Overnight Protocol (N=7)
	MinDC +2SD (pg/mL)
Eotaxin-2	4.4
MCP-2	2.2
BCA-1	1.3
MCP-4	3.4
I-309	1.4
IL-16	9.1
TARC	0.4
6CKine	45.4
Eotaxin-3	8.7
LIF	5.8
TPO	37.9
SCF	5.6
TSLP	3.1
IL-33	6.2
IL-20	53.3
IL-21	6.8
IL-23	31.5
TRAIL	3.5
CTACK	1.8
SDF-1 $\alpha$ + $\beta$	55.8
ENA-78	7.2
MIP-1 $\delta$	10.5
IL-28A	7.9

## Precision

Intra-assay precision is generated from the mean of the % CV's from 8 reportable results across two different concentration of cytokines in a single assay. Inter-assay precision is generated from the mean of the % CV's from 4-8 reportable results across two different concentrations of cytokine across 6 different assays.

Cytokines	Overnight Protocol (N=6 assays)	
	Intra-assay %CV	Inter-assay %CV
Eotaxin-2	4.5	8.1
MCP-2	4.5	7.3
BCA-1	4.8	6.2
MCP-4	3.6	6.1
I-309	7.0	11.0
IL-16	3.5	16.6
TARC	5.6	9.8
6CKine	8.6	13.7
Eotaxin-3	8.3	14.5
LIF	6.0	12.1
TPO	8.6	9.8
SCF	7.4	11.6
TSLP	6.9	13.8
IL-33	5.2	7.2
IL-20	4.6	7.6
IL-21	5.6	12.7
IL-23	4.6	7.5
TRAIL	5.3	10.2
CTACK	8.2	9.2
SDF-1 $\alpha$ + $\beta$	11.1	9.2
ENA-78	4.3	6.2
MIP-1 $\delta$	5.3	12.0
IL-28	4.4	10.1

## Accuracy

Defined as percent recovery, is generated from the mean of % recovery of 6 levels of cytokines spiked into matrix in 5 independent experiments.

Cytokines	Overnight Protocol
	% recovery in matrix
Eotaxin-2	100.4
MCP-2	106.2
BCA-1	101.3
MCP-4	104.6
I-309	103.4
IL-16	96.6
TARC	100.9
6CKine	93.3
Eotaxin-3	104.2
LIF	102.0
TPO	99.7
SCF	99.4
TSLP	109.5
IL-33	103.9
IL-20	98.5
IL-21	99.8
IL-23	99.7
TRAIL	102.9
CTACK	98.5
SDF-1 $\alpha$ + $\beta$	81.3
ENA-78	107.9
MIP-1 $\delta$	94.1

## TROUBLESHOOTING GUIDE

Problem	Probable Cause	Solution
Insufficient bead count	Plate washer aspirate height set too low	Adjust aspiration height according to manufacturers' instructions.
	Bead mix prepared inappropriately	Sonicate bead vials and vortex just prior to adding to bead mix bottle according to protocol. Agitate bead mix intermittently in reservoir while pipetting this into the plate.
	Samples cause interference due to particulate matter or viscosity	See above. Also sample probe may need to be cleaned with alcohol flush, back flushes and washes; or if needed probe should be removed and sonicated.
	Probe height not adjusted correctly	When reading the assay on Luminex® 200™, adjust probe height to the kit solid plate or to the recommended EMD Millipore filter plates using 3 alignment discs. When reading the assay on MAGPIX®, adjust probe height to the kit solid plate or to the recommended EMD Millipore filter plates using 2 alignment discs. When reading the assay on FLEXMAP 3D®, adjust probe height to the kit solid plate using 1 alignment disc.  For FLEXMAP 3D® when using the solid plate in the kit, the final resuspension should be with 150 µL Sheath Fluid in each well and 75 µL should be aspirated.
Background is too high	Background wells were contaminated	Avoid cross-well contamination by using sealer appropriately, and pipetting with multichannel pipettes without touching reagent in plate.
	Matrix used has endogenous analyte or interference	Check matrix ingredients for cross reacting components (e.g. interleukin modified tissue culture medium).
	Insufficient washes	Increase number of washes.
Beads not in region or gate	Luminex® instrument not calibrated correctly or recently	Calibrate Luminex® instrument based on manufacturer's instructions, at least once a week or if temperature has changed by >3°C.
	Gate settings not adjusted correctly	Some Luminex® instruments (e.g. Bio-Plex®) require different gate settings than those described in the kit protocol. Use instrument default settings.
	Wrong bead regions in protocol template	Check kit protocol for correct bead regions or analyte selection.
	Incorrect sample type used	Samples containing organic solvents or if highly viscous should be diluted or dialyzed as required.
	Instrument not washed or primed	Prime the Luminex® instrument 4 times to rid it of air bubbles, wash 4 times with Sheath Fluid or water if there is any remnant alcohol or sanitizing liquid.
	Beads were exposed to light	Keep plate and bead mix covered with dark lid or aluminum foil during all incubation steps.

<b>Problem</b>	<b>Probable Cause</b>	<b>Solution</b>
Signal for whole plate is same as background	<p>Incorrect or no Detection Antibody was added</p> <p>Streptavidin-Phycoerythrin was not added</p>	<p>Add appropriate Detection Antibody and continue.</p> <p>Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.</p>
Low signal for standard curve	<p>Detection Antibody may have been removed prior to adding Streptavidin Phycoerythrin</p> <p>Incubations done at inappropriate temperatures, timings or agitation</p>	<p>May need to repeat assay if desired sensitivity not achieved.</p> <p>Assay conditions need to be checked.</p>
Signals too high, standard curves are saturated	<p>Calibration target value set too high</p> <p>Plate incubation was too long with standard curve and samples</p>	<p>With some Luminex® instruments (e.g. Bio-Plex®) default target setting for RP1 calibrator is set at high PMT. Use low target value for calibration and reanalyze plate.</p> <p>Use shorter incubation time.</p>
Sample readings are out of range	<p>Samples contain no or below detectable levels of analyte</p> <p>Samples contain analyte concentrations higher than highest standard point</p> <p>Standard curve was saturated at higher end of curve</p>	<p>If below detectable levels, it may be possible to use higher sample volume. Check with technical support for appropriate protocol modifications.</p> <p>Samples may require dilution and reanalysis for just that particular analyte.</p> <p>See above.</p>
High variation in samples and/or standards	<p>Multichannel pipette may not be calibrated</p> <p>Plate washing was not uniform</p> <p>Samples may have high particulate matter or other interfering substances</p> <p>Plate agitation was insufficient</p> <p>Cross-well contamination</p>	<p>Calibrate pipettes.</p> <p>Confirm all reagents are removed completely in all wash steps.</p> <p>See above.</p> <p>Plate should be agitated during all incubation steps using an orbital plate shaker at a speed where beads are in constant motion without causing splashing.</p> <p>Check when reusing plate sealer that no reagent has touched sealer.</p> <p>Care should be taken when using same pipette tips that are used for reagent additions and that pipette tip does not touch reagent in plate.</p>

<b>FOR FILTER PLATES ONLY</b>		
<b>Problem</b>	<b>Probable Cause</b>	<b>Solution</b>
Filter plate will not vacuum	<p>Vacuum pressure is insufficient</p> <p>Samples have insoluble particles</p> <p>High lipid concentration</p>	<p>Increase vacuum pressure such that 0.2mL buffer can be suctioned in 3-5 seconds.</p> <p>Centrifuge samples just prior to assay set-up and use supernatant.</p> <p>After centrifugation, remove lipid layer and use supernatant.</p>
Plate leaked	<p>Vacuum pressure too high</p> <p>Plate set directly on table or absorbent towels during incubations or reagent additions</p> <p>Insufficient blotting of filter plate bottom causing wicking</p> <p>Pipette touching plate filter during additions</p> <p>Probe height not adjusted correctly</p> <p>Sample too viscous</p>	<p>Adjust vacuum pressure such that 0.2mL buffer can be suctioned in 3-5 seconds. May need to transfer contents to a new (blocked) plate and continue.</p> <p>Set plate on plate holder or raised edge so bottom of filter is not touching any surface.</p> <p>Blot the bottom of the filter plate well with absorbent towels after each wash step.</p> <p>Pipette to the side of plate.</p> <p>Adjust probe to 3 alignment discs in well H6.</p> <p>May need to dilute sample.</p>

**REPLACEMENT REAGENTS**

	<b>Catalog #</b>
Human Cytokine Panel II Standard	MXH8062
Human Cytokine Panel II Quality Controls	MXH6062
Serum Matrix	MXHSM-2
Human Cytokine Panel II Detection Antibodies	MXH1062
Streptavidin-Phycoerythrin	L-SAPE4
Assay Buffer	L-AB
Set of two 96-Well Plates with Sealers	MAG-PLATE
10X Wash Buffer	L-WB
Bead Diluent	LBD
Human Cytokine 23 Plex Premixed Magnetic Bead Panel 2– BULK PACKAGING	HCP2MAG62KPX23BK

**Antibody-Immobilized Magnetic Beads**

<u>Analytes</u>	<u>Bead #</u>	<u>Cat. #</u>
Premix 23 plex	--	HP2PMX23-MAG
Eotaxin-2	12	HETXN2-MAG
MCP-2	13	HMCP2-MAG
BCA-1	15	HBCA1-MAG
MCP-4	18	HMCP4-MAG
I-309	19	HI309-MAG
IL-16	21	HIL16-MAG
TARC	26	HTARC-MAG
6CKine	28	H6CKINE-MAG
Eotaxin-3	30	HETXN3-MAG
LIF	34	HLIF-MAG
TPO	36	HTPO-MAG
SCF	38	HCYSCF-MAG
TSLP	43	HTSLP-MAG
IL-33	45	HIL33-MAG
IL-20	51	HIL20--MAG
IL-21	52	HIL21-MAG
IL-23	54	HIL23-MAG
TRAIL	56	HCYTRAIL-MAG
CTACK	62	HCTACK-MAG
SDF-1 $\alpha$ + $\beta$	64	HSDF1AB-MAG
ENA-78	66	HENA78-MAG
MIP-1 $\delta$	76	HMIP1D-MAG
IL-28A	77	HIL28A-MAG



## **ORDERING INFORMATION**

To place an order or to obtain additional information about our immunoassay products, please contact your Customer Service or Technical Support Specialist.

Contact information for each region can be found on our website:

[emdmillipore.com/contact](http://emdmillipore.com/contact)

### **Conditions of Sale**

For Research Use Only. Not for Use in Diagnostic Procedures.

### **Safety Data Sheets (SDS)**

Safety Data Sheets for EMD Millipore products may be downloaded through our website at [emdmillipore.com/msds](http://emdmillipore.com/msds)

## WELL MAP

	1	2	3	4	5	6	7	8	9	10	11	12
A	Standard 0 (Background)	Standard 4	QC-2 Control									
B	Standard 0 (Background)	Standard 4	QC-2 Control									
C	Standard 1	Standard 5	Sample 1									
D	Standard 1	Standard 5	Sample 1									
E	Standard 2	Standard 6	Sample 2									
F	Standard 2	Standard 6	Sample 2									
G	Standard 3	QC-1 Control	Etc.									
H	Standard 3	QC-1 Control										