

User Guide

Human Cytokine/Chemokine Panel II Serum Free

96-Well Plate Assay

HCYP2SF-62K
HCP2SF-62K-PX23

Introduction.....	2	Preparation of Antibody-Immobilized Beads	11
Principle	3	Preparation of Quality Controls ...	11
Storage Conditions Upon Receipt.....	4	Preparation of Wash Buffer	11
Reagents Supplied.....	4	Preparation of Human Cytokine Panel II Standard.....	11
Human Cytokine/Chemokine Panel II Serum-Free Antibody-Immobilized Premixed Magnetic Beads	4	Immunoassay Procedure	14
Human Cytokine/Chemokine Panel II Serum-Free Antibody- Immobilized Magnetic Beads	5	Plate Washing	16
Materials Required.....	6	Solid Plate	16
Reagents	6	Filter Plate	16
Instrumentation/Materials	6	Equipment Settings.....	16
Safety Precautions.....	6	Quality Controls.....	19
Symbol Definitions.....	7	Assay Characteristics	19
Technical Guidelines	8	Cross-Reactivity	19
Sample Collection and Storage.....	10	Troubleshooting.....	20
Preparation of Tissue Culture		Product Ordering	25
Supernatant.....	10	Well Map	27
Preparation of Reagents for Immunoassay	11	Notice	28
		Technical Assistance.....	28
		Standard Warranty	28
		Contact Information	28
		Safety Data Sheets (SDS).....	28

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

Introduction

The MILLIPLEX® portfolio 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 verification process include: cross-reactivity, dilution linearity, kit stability, and sample behavior (for example detectability and stability).

Each MILLIPLEX® 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
- 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® Human Cytokine/Chemokine Panel II Serum-Free thus enables you to focus on the therapeutic potential of cytokines as well as the modulation of cytokine expression. Coupled with the Luminex® xMAP® 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.

MILLIPLEX® Human Cytokine/Chemokine Panel II Serum-Free 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 verify and build the most comprehensive library available for protein detection and quantitation.

MILLIPLEX® kits offer you:

- The ability to select a 23-plex or premixed option or
- 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.

MILLIPLEX® Human Cytokine / Chemokine Panel II Serum-Free to be used for the simultaneous quantification of any or all of the following analytes in 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 α + β /CXCL12, ENA-78/CXCL5, MIP 1 δ /MIP 5/CCL15, IL-28A/IFN λ 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® kits are based on the Luminex® xMAP® 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 beads known as MagPlex® microspheres.

- Luminex® uses proprietary techniques to internally color-code microspheres with two fluorescent dyes. Through precise concentrations of these dyes, distinctly colored bead sets of 500 5.6 µm polystyrene microspheres or 80 6.45 µm magnetic 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.
- The following Luminex® instruments can be used to acquire and analyze data using two detection methods:
- The Luminex® analyzers Luminex® 200™, FLEXMAP 3D®, and xMAP® INTELLIFLEX are flow cytometry-based instruments that integrate key xMAP® detection components, such as lasers, optics, advanced fluidics and high-speed digital signal processors.
- The Luminex® analyzer, MAGPIX®, is a CCD-based instrument that integrates key xMAP® 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. We combine the streamlined data acquisition power of Luminex® xPONENT® acquisition software with sophisticated analysis capabilities of MILLIPLEX® Analyst 5.1, integrating data acquisition and analysis seamlessly with all Luminex® instruments.
- xMAP® INTELLIFLEX runs on INTELLIFLEX software for instrument control, run setup and generating high quality data with flexible output options. Data can be exported in xPONENT® style CSV files for compatibility with many existing analytical applications, or in the new, customizable INTELLIFLEX file format. The INTELLIFLEX file format is intended for flexibility and simplicity, allowing the user to freely select which data points to include and to reduce the time to analysis.

The capability of adding multiple conjugated beads to each sample results in the ability to obtain multiple results from each sample. Open-architecture xMAP® 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

Reagents Supplied	Volume	Quantity	Component No.
Human Cytokine/Chemokine Panel II Standard	Lyophilized	1 vial	MXH8062
Human Cytokine Panel II Quality Controls 1 and 2	Lyophilized	2 vials	MXH6062
Set of one 96-Well Plate with 2 Sealers	-----	1 plate 2 sealers	-----
Assay Buffer	30 mL	1 bottle	L-AB
Bead Diluent	3.5 mL	1 bottle	LBD
10X Wash Buffer Note: Contains 0.05% Proclin™	60 mL	1 bottle	L-WB
Human Cytokine Panel II Detection Antibodies	3.2 mL	1 bottle	MXH1062
Streptavidin-Phycoerythrin	3.2 mL	1 bottle	L-SAPE4
Mixing Bottle (not provided with premixed panel)	-----	1 bottle	-----

Human Cytokine/Chemokine Panel II Serum-Free Antibody-Immobilized Premixed Magnetic Beads

Reagents Supplied	Volume	Quantity	Component No.
Premixed 23-plex Beads	3.5 mL	1 bottle	HP2PMX23-MAG

Included Human Cytokine/Chemokine Panel II Serum-Free Antibody-Immobilized Beads are dependent on customizable selection of analytes within the panel (see next page).

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

Human Cytokine/Chemokine Panel II Serum-Free Antibody-Immobilized Magnetic Beads

Bead/Analyte Name	Luminex® Magnetic Bead	Customizable 23 Analytes (50X concentration, 90 µL)		23-Plex Magnetic Premixed Beads + HCYIFNG-MAG
	Region	Available	Component No.	
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	✓

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

Materials Required (Not provided)

Reagents

MAGPIX® Drive Fluid PLUS (Catalogue No. 40-50030), xMAP® Sheath Fluid PLUS (Catalogue No. 40-50021), or xMAP® Sheath Concentrate PLUS (Catalogue No. 40-50023).

Instrumentation/Materials








- Adjustable Pipettes with Tips capable of delivering 25 µL to 1000 µL
- Multichannel Pipettes capable of delivering 5 µL to 50 µL, or 25 µL to 200 µL
- Reagent Reservoirs
- Polypropylene Microfuge Tubes
- Rubber Bands
- Aluminum Foil
- Absorbent Pads
- Laboratory Vortex Mixer
- Sonicator (Branson Ultrasonic Cleaner Model No. B200 or equivalent)
- Titer Plate Shaker (VWR® Microplate Shaker Catalogue No. 12620-926 or equivalent)
- Luminex® 200™, HTS, FLEXMAP 3D®, MAGPIX® with xPONENT® software or xMAP® INTELLIFLEX with INTELLIFLEX software by Luminex® Corporation.
- Automatic Plate Washer for magnetic beads (BioTek® 405 LS and 405 TS, Cat. Nos. 40-094, 40-095, 40-096, 40-097 or equivalent) or Handheld Magnetic Separation Block (Catalogue No. 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 (Catalogue No. MX-PLATE) to run the assay using a Vacuum Filtration Unit (Vacuum Manifold Catalogue No. MSVMHTS00 or equivalent with Vacuum Pump Catalogue No. 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.

Symbol Definitions

Ingredient	Cat. No.	Label	Label
Streptavidin-Phycoerythrin	L-SAPE4		Warning. 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		Warning. 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		Warning. 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	 	Danger. 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	 	Danger. 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.

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

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 ≤ -20 °C for 1 month and at ≤ -80 °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 Mixing 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 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 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 µL Sheath Fluid PLUS in each well and 75 µL should be aspirated.

For xMAP® INTELLIFLEX, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.

- For cell culture supernatants or tissue extraction, use the culture or extraction medium as the matrix solution in background, standard curve and control wells. If samples are diluted in Assay Buffer, use the Assay Buffer as matrix.
- 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 aggregates. Centrifuge samples before use.
- Vortex all reagents well before adding to plate.

Sample Collection and Storage

Preparation of Tissue Culture Supernatant

- Centrifuge the sample to remove debris and assay immediately or aliquot and store samples at ≤ -20 °C.
- Avoid multiple (> 2) freeze/thaw cycles.
- Tissue culture supernatant may require a dilution with an appropriate control medium prior to assay. Tissue/cell extracts should be done in neutral buffers containing reagents and conditions that do not interfere with assay performance. Excess concentrations of detergent, salt, denaturants, high or low pH, etc. will negatively affect the assay. Organic solvents should be avoided. The tissue/cell extract samples should be free of particles such as cells or tissue debris.

Note:

- A maximum of 25 μ L per well of diluted or neat 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 anti-coagulant 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

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 μL 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 $^{\circ}\text{C}$ 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 antibody-immobilized beads, add 60 μL from each of the 5 bead vials to the Mixing Bottle. Then add 2.7 mL Bead Diluent.

Example 2: When using 9 antibody-immobilized beads, add 60 μL from each of the 9 bead vials to the Mixing Bottle. Then add 2.46 mL Bead Diluent.

Preparation of Quality Controls

Before use, reconstitute Quality Control 1 and Quality Control 2 with 250 μL deionized water. Invert the vial several times to mix and vortex. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted Quality Control 1 and Quality Control 2 into two polypropylene microfuge tubes. Unused portion may be stored at $\leq -20^{\circ}\text{C}$ for up to one month.

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 the unused portion at 2-8 $^{\circ}\text{C}$ for up to one month.

Preparation of Human Cytokine Panel II Standard

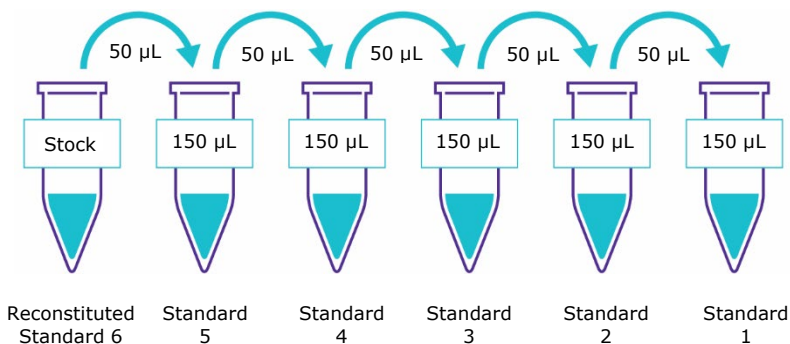
1. Prior to use, reconstitute the Human Cytokine Panel II Standard with 250 μL deionized water. Refer to table below for analyte concentrations. Invert the vial several times to mix. Vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted standard to a polypropylene microfuge tube. This will be used as "Standard 6"; the unused portion may be stored at $\leq -20^{\circ}\text{C}$ for up to one month.
2. Preparation of Working Standards
Label 5 polypropylene microfuge tubes Standard 1 through Standard 5. Add 150 μL of Assay Buffer to each of the 5 tubes. Prepare serial dilutions by adding 50 μL of the reconstituted Standard 6 to the Standard 5 tube, mix well and transfer 50 μL of Standard 5 to the Standard 4 tube, mix well and transfer 50 μL of Standard 4 to the Standard 3 tube, mix well and transfer 50 μL of Standard 3 to the Standard 2 tube, mix well and transfer 50 μL of Standard 2 to the Standard 1 tube and mix well. The Standard 0 (Background) will be Assay Buffer.

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

Standard No.	Volume of Deionized Water to Add	Volume of Standard to Add
Original (Standard 6)	250 μL	0

Standard No.	Add Assay Buffer (μL)	Add Standard (volume)
Standard 5	150 μL	50 μL of Standard 6
Standard 4	150 μL	50 μL of Standard 5
Standard 3	150 μL	50 μL of Standard 4
Standard 2	150 μL	50 μL of Standard 3
Standard 1	150 μL	50 μL of Standard 2

Preparation of Standards



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

Standard Tube No.	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)
1	1.0	2.0	4.9	9.8
2	3.9	7.8	19.5	39.1
3	15.6	31.3	78.1	156.3
4	62.5	125	312.5	625
5	250	500	1250	2500
6	1000	2000	5000	10,000

Standard Tube No.	ENA-78, 6Ckine, LIF, IL-21, IL-33 (pg/mL)	MIP-1δ, Eotaxin-3, IL-23, TPO, IL-20 (pg/mL)	SDF-1α+β, (pg/mL)
1	19.5	48.8	97.7
2	78.1	195.3	390.6
3	312.5	781.3	1562.5
4	1250	3125	6250
5	5000	12500	25,000
6	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), [Standard 1 through 6], Controls 1 and 2, and Samples on Well Map Worksheet in a vertical configuration. It is recommended to run the assay in duplicate.

Note: Most instruments will only read the 96-well plate vertically by default.

- 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.

1. Add 200 μ L of Wash Buffer into each well of the plate. Seal and mix 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 μ L of each Standard or Control into the appropriate wells. Assay Buffer should be used for 0 pg/mL standard (Background).
4. Add 25 μ L of Assay Buffer to the sample wells.
5. Add 25 μ L of appropriate control culture media solution to the background, standards, and control wells.
6. Add 25 μ L of Sample into the appropriate wells.
7. Vortex Mixing Bottle and add 25 μ 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.

Add 200 μ L Wash Buffer per well



Shake 10 min, RT
Decant

- Add 25 μ L Standard or Control to appropriate wells
- Add 25 μ L Assay Buffer to background and sample wells
- Add 25 μ L appropriate control culture media to background, standards, and control wells
- Add 25 μ L Samples to sample wells
- Add 25 μ L Beads to each well



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

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

- Add 25 μL of Detection Antibodies into each well.

Note: Allow the Detection Antibodies to warm to room temperature prior to addition.

- Seal, cover with foil and incubate with agitation on a plate shaker for 1 hour at room temperature (20-25 $^{\circ}\text{C}$). **DO NOT ASPIRATE AFTER INCUBATION.**
- Add 25 μL Streptavidin-Phycoerythrin to each well containing the 25 μL of Detection Antibodies.
- Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25 $^{\circ}\text{C}$).

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

- Add 150 μL of Sheath Fluid PLUS (or Drive Fluid PLUS if using MAGPIX[®]) to all wells. Resuspend the beads on a plate shaker for 5 minutes.

- Run plate on Luminex[®] 200™, HTS, FLEXMAP 3D[®], MAGPIX[®] with xPONENT[®] software or xMAP[®] INTELLIFLEX with INTELLIFLEX Software.

- Save and analyze the Median Fluorescent Intensity (MFI) data using a 5-parameter logistic or spline curve-fitting method for calculating analyte concentrations in samples.

Note: For diluted samples, multiply the calculated concentration by the dilution factor.



Remove well contents and wash 2X with 200 μL Wash Buffer

Add 25 μL Detection Antibodies per well



Incubate 1 hour at RT

Do Not Aspirate

Add 25 μL Streptavidin-Phycoerythrin per well



Incubate 30 minutes at RT

Remove well contents and wash 2X with 200 μL Wash Buffer

Add 150 μL Sheath Fluid PLUS or Drive Fluid PLUS per well

Read on Luminex[®] instrument (100 μL , 50 beads per bead set)

Plate Washing

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

Solid Plate

- Handheld magnet (Catalogue No. 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 μ 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.

- Magnetic plate washer (Catalogue Nos. 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 μ 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.

Filter Plate (Catalogue No. MX-PLATE)

If using a filter plate, use a vacuum filtration manifold to remove well contents. Wash plate with 200 μ 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®, MAGPIX® with xPONENT® software and xMAP® INTELLIFLEX with INTELLIFLEX software:

These specifications are for the above listed instruments and software. Luminex® instruments with other software (for example 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, each instrument must be calibrated, and performance verified with the indicated calibration and verification kits.

Instrument	Calibration Kit	Verification Kit
Luminex® 200™ and HTS	xPONENT® 3.1 compatible Calibration Kit (Catalogue No. LX2R-CAL-K25)	Performance Verification Kit (Catalogue No. LX2R-PVER-K25)
FLEXMAP 3D®	FLEXMAP 3D® Calibrator Kit (Catalogue No. F3D-CAL-K25)	FLEXMAP 3D® Performance Verification Kit (Catalogue No. F3D-PVER-K25)
xMAP® INTELLIFLEX	xMAP® INTELLIFLEX Calibration Kit (Catalogue No. IFX-CAL-K20)	xMAP® INTELLIFLEX Performance Verification Kit (Catalogue No. IFX-PVER-K20)
MAGPIX®	MAGPIX® Calibration Kit (Catalogue No. MPX-CAL-K25)	MAGPIX® Performance Verification Kit (Catalogue No. MPX-PVER-K25)

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.

The Luminex® probe height must be adjusted to the plate provided in the kit. Please use Catalogue No. MAG-PLATE, if additional plates are required for this purpose.

Events:	50, per bead	
Sample Size:	100 μ L	
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-1 α + β	64	64
ENA-78	66	66
MIP-1 δ	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 our website [SigmaAldrich.com](https://www.sigmaaldrich.com) using the catalogue number as the keyword.

Assay Characteristics

Cross-Reactivity

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

Troubleshooting

Problem	Probable Cause	Solution
	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 flushes, back flushes, and washes; or, if needed, probe should be removed and sonicated.
Insufficient bead count	Probe height not adjusted correctly	When reading the assay on Luminex® 200™, adjust probe height to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on MAGPIX®, adjust probe height to the kit solid plate or to the recommended 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 PLUS in each well and 75 µL should be aspirated.
		When reading the assay on xMAP® INTELLIFLEX, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.

Problem	Probable Cause	Solution
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 (for example, 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 (for example 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 PLUS 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.

Problem	Probable Cause	Solution
Signal for whole plate is same as background	Incorrect or no Detection Antibody was added	Add appropriate Detection Antibody and continue.
	Streptavidin-Phycoerythrin was not added	Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.
Low signal for standard curve	Detection Antibody may have been removed prior to adding Streptavidin-Phycoerythrin	May need to repeat assay if desired sensitivity not achieved.
	Incubations done at inappropriate temperatures, timings, or agitation	Assay conditions need to be checked.
Signals too high, standard curves are saturated	Calibration target value set too high	With some Luminex® instruments (for example, Bio-Plex®) default target setting for RP1 calibrator is set at high PMT. Use low target value for calibration and reanalyze plate.
	Plate incubation was too long with standard curve and samples	Use shorter incubation time.
Sample readings are out of range	Samples contain no or below detectable levels of analyte	If below detectable levels, it may be possible to use higher sample volume. Check with technical support for appropriate protocol modifications.
	Samples contain analyte concentrations higher than highest standard point	Samples may require dilution and reanalysis for just that particular analyte.
	Standard curve was saturated at higher end of curve	See above.

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

Problem	Probable Cause	Solution
High variation in samples and/or standards	Multichannel pipette may not be calibrated	Calibrate pipettes.
	Plate washing was not uniform	Confirm all reagents are removed completely in all wash steps.
	Samples may have high particulate matter or other interfering substances	See above.
	Plate agitation was insufficient	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.
	Cross-well contamination	Check when reusing plate sealer that no reagent has touched sealer. 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.

For Filter Plates Only

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

Product Ordering

Order products online at [SigmaAldrich.com](https://www.sigmaaldrich.com).

Replacement Reagents	Cat. No.
Human Cytokine Panel II Standard	MXH8062
Human Cytokine Panel II Quality Controls	MXH6062
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

Antibody-Immobilized Magnetic Beads

Analytes	Bead No.	Cat. No.
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 α + β	64	HSDF1AB-MAG
ENA-78	66	HENA78-MAG
MIP-1 δ	76	HMIP1D-MAG
IL-28A	77	HIL28A-MAG

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

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										

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

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page at SigmaAldrich.com/techservice.

Standard Warranty

The applicable warranty for the products listed in this publication may be found at SigmaAldrich.com/terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/offices.

Safety Data Sheets (SDS)

Safety Data Sheets are available on the product page at SigmaAldrich.com.

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma, Millipore, Milliplex and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2021-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

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