



Human Th17 Magnetic Bead Panel

96-Well Plate Assay

**Cat. # HTH17MAG-14K,
HT17MG-14K-PX25,
HT17MAG14PMX25BK**

The M logo is a trademark of Merck KGaA, Darmstadt, Germany.
© 2013 EMD Millipore Corporation, Billerica, MA 01821 USA.

HTH17MAG-14K
HT17MG-14K-PX25
HT17MAG14PMX25BK

MILLIPLEX[®] MAP

HUMAN Th17 MAGNETIC BEAD PANEL 96-Well Plate Assay

HTH17MAG-14K,
HT17MG-14K-PX25,
HT17MAG14PMX25BK

| <u>TABLE OF CONTENTS</u> | <u>PAGE</u> |
|---|--------------------|
| Introduction | 2 |
| Principle | 3 |
| Storage Conditions Upon Receipt | 3 |
| 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 | 13 |
| Plate Washing | 16 |
| Equipment Settings | 16 |
| Quality Controls | 18 |
| Assay Characteristics | 18 |
| Troubleshooting Guide | 21 |
| Replacement Reagents | 23 |
| Ordering Information | 25 |
| Well Map | 26 |

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 Th17 Magnetic Bead Panel

INTRODUCTION

CD4 T-helper cells are major players in adaptive immunity. They can be divided into the major subsets Th1, Th2, Th17 and Treg based on their expression profile of transcription factors and secreted cytokines. Th1 cells, induced by IL-12 and enhanced by IFN γ , secrete IFN γ for activating macrophages. Th2 cells, induced by IL-4, produce mainly IL-4, IL-5, IL-13, IL-25 and IL-17E and play a role in allergy. Treg cells, induced by TGF β , produce the anti-inflammatory cytokines IL-10 and TGF β and function to control T-cell responses to prevent autoimmune reactions. Th17 cells, a more recently discovered subset characterized by the production of IL-17, are induced by TGF β , IL-6 and IL-21 in mice or TGF β combined with IL-23 and IL-1 β instead of IL-6 and IL-21 in humans. Activated Th17 cells secrete IL-17A, IL-17F, IL-21, IL-22 and TNF α which promote tissue inflammation.

Th17 cells are involved in the clearance of extracellular bacteria and fungi. They are abundant in the intestinal lamina propria and function as a barrier against invading pathogens.

Excessive amounts of Th17 cells have been implicated in the pathogenesis of several autoimmune diseases such as multiple sclerosis, psoriasis, autoimmune uveitis, juvenile diabetes, rheumatoid arthritis, and Crohn's disease. In addition, Th17 cells play an important role in cancer development and progression, potentially playing a role in both promoting and inhibiting tumor growth.

MILLIPLEX[®] 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[®] 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[®] MAP Human Th17 Magnetic Bead Panel thus enables you to focus on the therapeutic potential of the modulation of CD4 T-helper cell cytokines. 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.

EMD Millipore's MILLIPLEX[®] MAP Human Th17 Magnetic Bead Panel is part of the most versatile system available for CD4 T-helper cell cytokine 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[®] MAP offers you:
 - The ability to select a 25-plex premixed option or
 - The ability to choose any combination of analytes from our panel of 25 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 MAP Human Th17 Magnetic Bead Panel is to be used for the simultaneous quantification of any or all of the following analytes in human serum, plasma, tissue/cell lysate and culture supernatant samples: IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-9, IL-10, IL-12p70, IL-13, IL-15, IL-17A, IL-17E/IL-25, IL-17F, IL-21, IL-22, IL-23, IL-27, IL-28A, IL-31, IL-33, GM-CSF, IFN γ , MIP-3 α , TNF α and TNF β .

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[®] MAP is 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[®]-C 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 μ m polystyrene microspheres or 80 6.45 μ m 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.
- EMD Millipore provides three Luminex[®] instruments to acquire and analyze data using two detection methods:
 - The Luminex[®] analyzers Luminex[®] 200[™] and FLEXMAP 3D[®], 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[®]), 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. EMD Millipore combines the streamlined data acquisition power of Luminex[®] xPONENT[®] acquisition software with sophisticated analysis capabilities of the new MILLIPLEX[®] Analyst 5.1, integrating data acquisition and analysis seamlessly with all Luminex[®] 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[®] 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 \leq -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 Th17 Standard | HTH17-8014 | Lyophilized | 1 vial |
| Human Th17 Quality Controls 1 and 2 | HTH17-6014 | Lyophilized | 2 vials |
| Serum Matrix Note: Contains 0.08% Sodium Azide | MXHSM | Lyophilized | 2 vials |
| Set of one 96-Well Plate with 2 sealers | ----- | ----- | 1 plate 2 sealers |
| Assay Buffer | L-AB | 30 mL | 1 bottle |
| 10X Wash Buffer Note: Contains 0.05% Proclin | L-WB | 60 mL | 1 bottle |
| Human Th17 Detection Antibodies | HTH17-1014 | 3.2 mL | 1 bottle |
| Streptavidin-Phycoerythrin | L-SAPE11 | 3.2 mL | 1 bottle |
| Mixing Bottle | ----- | ----- | 1 bottle |
| Bead Diluent | LBD | 3.5 mL | 1 bottle |

Human Th17 Antibody-Immobilized Premixed Magnetic Beads:

| | | | |
|--|--------------------------------|----------------------|------------------------|
| Premixed 24-plex Beads + IFN γ | HTH17PMX24-MAG, HCYIFNG-MAG | 3.5 mL 90 μ L | 1 bottle + 1 bead vial |
|--|--------------------------------|----------------------|------------------------|

Included Human Th17 Antibody-Immobilized Magnetic Beads are dependent on customizable selection of analytes within the panel (see nex page).

REAGENTS SUPPLIED (continued)

Human Th17 Antibody-Immobilized Magnetic Beads:

| Bead/Analyte Name | Luminex® Magnetic Bead Region | Customizable 25 Analytes (50X concentration, 90 µL) | | 24-Plex Premixed Beads + HCYIFNG-MAG |
|--------------------------------|--|--|--------------|---|
| | | Available | Cat. # | |
| Anti-Human IL-17F | 19 | ✓ | HIL17F-MAG | ✓ |
| Anti-Human GM-CSF | 20 | ✓ | HGMCSF-MAG | ✓ |
| Anti-Human IFN γ | 25 | ✓ | HCYIFNG-MAG | ✓ Single vial |
| Anti-Human IL-10 | 27 | ✓ | HCYIL10-MAG | ✓ |
| Anti-Human CCL20/MIP3 α | 28 | ✓ | HMIP3A-MAG | ✓ |
| Anti-Human IL-12P70 | 33 | ✓ | HIL12P70-MAG | ✓ |
| Anti-Human IL-13 | 35 | ✓ | HIL13-MAG | ✓ |
| Anti-Human IL-15 | 37 | ✓ | HIL15-MAG | ✓ |
| Anti-Human IL-17A | 39 | ✓ | HIL17-MAG | ✓ |
| Anti-Human IL-22 | 43 | ✓ | HIL22-MAG | ✓ |
| Anti Human IL-9 | 45 | ✓ | HIL9-MAG | ✓ |
| Anti-Human IL-1 β | 46 | ✓ | HCYIL1B-MAG | ✓ |
| Anti-Human IL-33 | 47 | ✓ | HTHIL33-MAG | ✓ |
| Anti-Human IL-2 | 48 | ✓ | HIL2-MAG | ✓ |
| Anti-Human IL-21 | 52 | ✓ | HIL21-MAG | ✓ |
| Anti-Human IL-4 | 53 | ✓ | HIL4-MAG | ✓ |
| Anti-Human IL-23 | 54 | ✓ | HIL23-MAG | ✓ |
| Anti-Human IL-5 | 55 | ✓ | HIL5-MAG | ✓ |
| Anti-Human IL-6 | 57 | ✓ | HCYIL6-MAG | ✓ |
| Anti-Human IL-17E/IL-25 | 66 | ✓ | HIL17E-MAG | ✓ |
| Anti-Human IL-27 | 72 | ✓ | HIL27-MAG | ✓ |
| Anti-Human IL-31 | 73 | ✓ | HIL31-MAG | ✓ |
| Anti-Human TNF α | 75 | ✓ | HCYTNFA-MAG | ✓ |
| Anti-Human TNF β | 76 | ✓ | HTNFB-MAG | ✓ |
| Anti-Human IL-28A | 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 | |
|--------------------------------------|------------|---|---|
| Human Th17 Detection Antibodies | HTH17-1014 |  | 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 Th17 Quality Controls 1 & 2 | HTH17-6014 |   | 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 Th17 Standard | HTH17-8014 |   | 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. |
| Streptavidin- Phycoerythrin | L-SAPE11 |  | 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 - MILLIPLEX | L-WB |  | Warning. May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water. |
| Serum Matrix | MXHSM | No Symbol Required | Harmful to aquatic life with long lasting effects. Avoid release into the environment. |

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 should be discarded.**
- 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 5-10 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 μL Sheath Fluid in each well and 75 μL should be aspirated.

TECHNICAL GUIDELINES (continued)

- 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 serum/plasma samples use the Serum Matrix provided in the kit for further dilution.
- 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

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.
- It is recommended to use neat samples in the assay.

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.
- It is recommended to use neat samples in the assay.

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. 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.

SAMPLE COLLECTION AND STORAGE (continued)

NOTE:

- A maximum of 25 μL per well of neat 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

Prepare premixed beads and individually mixed beads immediately prior to use, any unused mixed Antibody-Immobilized Beads should be discarded.

If premixed beads are used, sonicate the premixed bead bottle 30 seconds then vortex for 1 minute before use. To prepare 25-plex premixed beads, add 70 μL of HCYIFNG-MAG bead to the 24-plex premixed beads. Mix well 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.

Example 1: When using 20 antibody-immobilized beads, add 60 μL from each of the 20 bead sets to the Mixing Bottle. Then add 1.8 mL Bead Diluent.

Example 2: When using 9 antibody-immobilized beads, add 60 μL 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 μL deionized water. Invert the vial several times to mix and vortex. Allow the vial to sit for 5-10 minutes and then transfer the controls to appropriately labeled polypropylene microfuge tubes. Unused portion may be stored at $\leq 20^{\circ}\text{C}$ 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 $2-8^{\circ}\text{C}$ for up to one month.

D. Preparation of Serum Matrix

This step is required for serum or plasma samples only. Add 0.5 mL deionized water 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.

PREPARATION OF REAGENTS FOR IMMUNOASSAY (continued)

E. Preparation of Human Th17 Standard

1.) Prior to use, reconstitute the Human Th17 Standard with 250 μL 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 stock standard (Standard 7); the unused portions of this stock may be stored at $\leq -20^{\circ}\text{C}$ for up to one month.

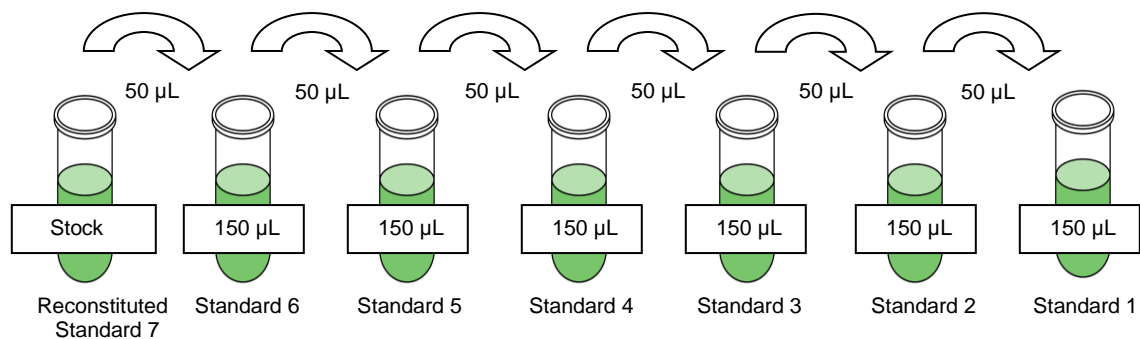
2.) Preparation of Working Standards

Label six polypropylene microfuge tubes as Standard 6, Standard 5, Standard 4, Standard 3, Standard 2 and Standard 1. Add 150 μL of Assay Buffer to each of the six tubes. Prepare 1:4 serial dilutions by adding 50 μL of the reconstituted Standard 7 to the Standard 6 tube, mix well and transfer 50 μL of the Standard 6 to the Standard 5 tube, mix well and transfer 50 μL of the Standard 5 to the Standard 4 tube, mix well and transfer 50 μL of the Standard 4 to the Standard 3 tube, mix well and transfer 50 μL of the Standard 3 to the Standard 2 tube, mix well and transfer 50 μL of the Standard 2 to the Standard 1 tube. The 0 Standard (Background) will be the Assay Buffer.

| Reconstituted Standard | Volume of Deionized Water to Add | Volume of Standard to Add |
|------------------------|----------------------------------|---------------------------|
| Standard 7 | 250 μL | 0 |

| Standard Number | Volume of Assay Buffer to Add | Volume of Standard to Add |
|-----------------|-------------------------------|--------------------------------|
| Standard 6 | 150 μL | 50 μL of Standard 7 |
| 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



PREPARATION OF REAGENTS FOR IMMUNOASSAY (continued)

After dilution, each tube has the following concentrations for each analyte:

| Standard Dilution | IL-17F & IL-4 | GMCSF & IL-27 | IFN γ | IL-10 | MIP3 α , IL12p70, IL-15, IL-1 β , IL-33 & IL-21 | IL-13 |
|-------------------|---------------|---------------|--------------|---------|--|---------|
| | (ng/mL) | (ng/mL) | (pg/mL) | (pg/mL) | (pg/mL) | (pg/mL) |
| Standard 7 | 100 | 250 | 40,000 | 5,000 | 20,000 | 30,000 |
| Standard 6 | 25 | 63 | 10,000 | 1,250 | 5,000 | 7,500 |
| Standard 5 | 6.2 | 16 | 2,500 | 313 | 1,250 | 1,875 |
| Standard 4 | 1.6 | 4 | 625 | 78 | 313 | 469 |
| Standard 3 | 0.4 | 1 | 156 | 20 | 78 | 117 |
| Standard 2 | 0.09 | 0.2 | 39 | 5 | 20 | 29 |
| Standard 1 | 0.02 | 0.06 | 10 | 1 | 5 | 7 |

| Standard Dilution | IL-17A & IL-2 | IL-22 & TNF β | IL-9 | IL-23 | IL-5 | IL-6 & TNF α |
|-------------------|---------------|---------------------|---------|---------|---------|---------------------|
| | (pg/mL) | (ng/mL) | (pg/mL) | (ng/mL) | (pg/mL) | (pg/mL) |
| Standard 7 | 50,000 | 150 | 35,000 | 1,500 | 25,000 | 10,000 |
| Standard 6 | 12,500 | 38 | 8,750 | 375 | 6,250 | 2,500 |
| Standard 5 | 3,125 | 9.4 | 2,188 | 94 | 1,563 | 625 |
| Standard 4 | 781 | 2.4 | 547 | 24 | 391 | 156 |
| Standard 3 | 195 | 0.6 | 137 | 6 | 98 | 39 |
| Standard 2 | 49 | 0.2 | 34 | 1.5 | 24 | 10 |
| Standard 1 | 12 | 0.04 | 9 | 0.3 | 6 | 2.5 |

| Standard Dilution | IL-31 & IL-28A | IL-17E/IL-25 |
|-------------------|----------------|--------------|
| | (ng/mL) | (ng/mL) |
| Standard 7 | 200 | 500 |
| Standard 6 | 50 | 125 |
| Standard 5 | 13 | 31.2 |
| Standard 4 | 3.1 | 7.81 |
| Standard 3 | 0.8 | 1.95 |
| Standard 2 | 0.2 | 0.49 |
| Standard 1 | 0.04 | 0.12 |

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 7], 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.

1. Add 200 μL of Assay 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 Assay 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 μL of each Standard or Control into the appropriate wells. Assay Buffer should be used for 0 standard (Background).
4. Add 25 μL of Assay Buffer to the sample wells.
5. Add 25 μL of appropriate matrix solution to the background, standards, and control wells. When assaying serum or plasma, use the Serum Matrix provided in the kit as the matrix solution. When assaying tissue culture or other supernatant, use proper control culture medium as the matrix solution.
6. Add 25 μL of neat Samples into the appropriate wells.
7. Vortex Mixing Bottle and add 25 μL 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 (16–18 hours) at 4°C.
9. Gently remove well contents and wash plate two times following instructions listed in the **PLATE WASHING** section.

Add 200 μL Assay Buffer per well



Shake 10 min, RT

Decant

- Add 25 μL Standard or Control to appropriate wells
- Add 25 μL Assay Buffer to background and sample wells
- Add 25 μL appropriate matrix solution to background, standards, and control wells
- Add 25 μL neat Samples to sample wells
- Add 25 μL Beads to each well



Incubate overnight (16-18 hours) at 4°C



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

10. Add 25 μ L of Detection Antibodies into each well. (Note: Allow the Detection Antibodies to warm to room temperature prior to addition.)
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 μ L Streptavidin-Phycoerythrin to each well containing the 25 μ 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 two times following instructions listed in the **PLATE WASHING** section.
15. Add 150 μ L of Sheath Fluid (or Drive Fluid if using MAGPIX[®]) to all wells. Resuspend the beads on a plate shaker for 5 minutes.
16. Run plate on Luminex[®] 200[™], HTS, FLEXMAP 3D[™] or MAGPIX[®] with xPONENT software.
17. 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.)

Add 25 μ L Detection Antibodies per well



Incubate 1 hour at RT (20-25°C)

Do Not Aspirate

Add 25 μ L Streptavidin-Phycoerythrin per well



Incubate for 30 minutes at RT



Remove well contents and wash 2X with 200 μ L Wash Buffer

Add 150 μ L Sheath Fluid or Drive Fluid per well

Read on Luminex[®] 100 μ 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 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.

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 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.

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 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®, 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 | |
| Reporter: | Default (low PMT) | |
| Time: Out: | 60 seconds | |
| Bead: Set: | Customizable 25-Plex Beads | |
| | IL-17F | 19 |
| | GM-CSF | 20 |
| | IFN γ | 25 |
| | IL-10 | 27 |
| | CCL20/MIP3 α | 28 |
| | IL-12p70 | 33 |
| | IL-13 | 35 |
| | IL-15 | 37 |
| | IL-17A | 39 |
| | IL-22 | 43 |
| | IL-9 | 45 |
| | IL-1 β | 46 |
| | IL-33 | 47 |
| | IL-2 | 48 |
| | IL-21 | 52 |
| | IL-4 | 53 |
| | IL-23 | 54 |
| | IL-5 | 55 |
| | IL-6 | 57 |
| | IL-17E/IL-25 | 66 |
| | IL-27 | 72 |
| | IL-31 | 73 |
| | TNF α | 75 |
| | TNF β | 76 |
| | IL-28A | 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 emdmillipore.com using the catalog 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.

Assay Sensitivities (minimum detectable concentrations)

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.

| Analyte | Overnight Protocol (n = 8 Assays) | |
|-----------------------------|-----------------------------------|-----------|
| | MinDC | MinDC+2SD |
| IL-17F (ng/mL) | 0.009 | 0.016 |
| GM-CSF (ng/mL) | 0.146 | 0.206 |
| IFN γ (pg/mL) | 1.8 | 2.4 |
| IL-10 (pg/mL) | 0.3 | 0.5 |
| CCL20/MIP3 α (pg/mL) | 2.2 | 3.4 |
| IL-12p70 (pg/mL) | 1.1 | 2.2 |
| IL-13 (pg/mL) | 2.4 | 3.5 |
| IL-15 (pg/mL) | 2.7 | 4.6 |
| IL-17A (pg/mL) | 2.1 | 2.8 |
| IL-22 (ng/mL) | 0.021 | 0.032 |
| IL-9 (pg/mL) | 6 | 8.7 |
| IL-1 β (pg/mL) | 2.1 | 3.6 |
| IL-33 (pg/mL) | 6.3 | 10.9 |
| IL-2 (pg/mL) | 5.1 | 9 |
| IL-21 (pg/mL) | 2 | 3.3 |
| IL-4 (ng/mL) | 0.009 | 0.014 |
| IL-23 (ng/mL) | 0.098 | 0.169 |
| IL-5 (pg/mL) | 1.2 | 1.5 |
| IL-6 (pg/mL) | 1.7 | 2.9 |
| IL-17E/IL-25 (ng/mL) | 0.099 | 0.186 |
| IL-27 (ng/mL) | 0.063 | 0.123 |
| IL-31 (ng/mL) | 0.021 | 0.041 |
| TNF α (pg/mL) | 0.9 | 1.7 |
| TNF β (ng/mL) | 0.021 | 0.035 |
| IL-28A (ng/mL) | 0.038 | 0.082 |

ASSAY CHARACTERISTICS (continued)

Precision

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

| Analyte | Overnight Protocol | |
|---------------------|--------------------|-----------------|
| | Intra-assay %CV | Inter-assay %CV |
| IL-17F | 2 | 10 |
| GM-CSF | 5 | 13 |
| IFN γ | 4 | 13 |
| IL-10 | 3 | 11 |
| CCL20/MIP3 α | 5 | 10 |
| IL-12p70 | 3 | 8 |
| IL-13 | 3 | 9 |
| IL-15 | 4 | 12 |
| IL-17A | 3 | 13 |
| IL-22 | 4 | 8 |
| IL-9 | 3 | 5 |
| IL-1 β | 4 | 11 |
| IL-33 | 5 | 11 |
| IL-2 | 3 | 11 |
| IL-21 | 3 | 11 |
| IL-4 | 4 | 11 |
| IL-23 | 3 | 9 |
| IL-5 | 4 | 7 |
| IL-6 | 5 | 7 |
| IL-17E/IL-25 | 5 | 10 |
| IL-27 | 4 | 10 |
| IL-31 | 4 | 8 |
| TNF α | 4 | 11 |
| TNF β | 3 | 7 |
| IL-28A | 3 | 13 |

ASSAY CHARACTERISTICS (continued)

Accuracy

Spike Recovery: The data represent mean percent recovery of spiked standards ranging from low, medium, and high concentration in serum matrices (n=5).

| Analyte | Overnight Protocol |
|---------------------|----------------------------|
| | % Recovery in Serum Matrix |
| IL-17F | 104 |
| GM-CSF | 105 |
| IFN γ | 107 |
| IL-10 | 96 |
| CCL20/MIP3 α | 101 |
| IL-12p70 | 96 |
| IL-13 | 96 |
| IL-15 | 94 |
| IL-17A | 101 |
| IL-22 | 101 |
| IL-9 | 100 |
| IL-1 β | 96 |
| IL-33 | 104 |
| IL-2 | 98 |
| IL-21 | 100 |
| IL-4 | 79 |
| IL-23 | 100 |
| IL-5 | 97 |
| IL-6 | 96 |
| IL-17E/IL-25 | 106 |
| IL-27 | 104 |
| IL-31 | 102 |
| TNF α | 101 |
| TNF β | 102 |
| IL-28A | 98 |

TROUBLESHOOTING GUIDE:

| Problem | Probable Cause | Solution |
|-----------------------------|--|--|
| Insufficient Bead Count | <p>Plate Washer aspirate height set too low</p> <p>Bead mix prepared inappropriately</p> <p>Samples cause interference due to particulate matter or viscosity</p> <p>Probe height not adjusted correctly</p> | <p>Adjust aspiration height according to manufacturers' instructions.</p> <p>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.</p> <p>See above. Also sample probe may need to be cleaned with Alcohol flush, Back flush and washes; or if needed probe should be removed and sonicated.</p> <p>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.</p> |
| Background is too high | <p>Background wells were contaminated</p> <p>Matrix used has endogenous analyte or interference</p> <p>Insufficient washes</p> | <p>Avoid cross-well contamination by using sealer appropriately, and pipeting with Multichannel pipets without touching reagent in plate.</p> <p>Check matrix ingredients for cross reacting components (e.g. interleukin modified tissue culture medium).</p> <p>Increase number of washes.</p> |
| Beads not in region or gate | <p>Luminex[®] not calibrated correctly or recently</p> <p>Gate Settings not adjusted correctly</p> <p>Wrong bead regions in protocol template</p> <p>Incorrect sample type used</p> <p>Instrument not washed or primed</p> <p>Beads were exposed to light</p> | <p>Calibrate Luminex[®] based on Instrument Manufacturer's instructions, at least once a week or if temperature has changed by >3°C.</p> <p>Some Luminex[®] instruments (e.g. Bioplex) require different gate settings than those described in the Kit protocol. Use Instrument default settings.</p> <p>Check kit protocol for correct bead regions or analyte selection.</p> <p>Samples containing organic solvents or if highly viscous should be diluted or dialyzed as required.</p> <p>Prime the Luminex[®] 4 times to rid of air bubbles, wash 4 times with sheath fluid or water if there is any remnant alcohol or sanitizing liquid</p> <p>Keep plate and bead mix covered with dark lid or aluminum foil during all incubation steps.</p> |

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

REPLACEMENT REAGENTS

| | |
|--|------------------|
| Human Th17 Standard | HTH17-8014 |
| Human Th17 Quality Controls 1 & 2 | HTH17-6014 |
| Human Th17 Detection Antibodies | HTH17-1014 |
| Streptavidin-Phycoerythrin | L-SAPE11 |
| Serum Matrix | MXHSM |
| Bead Diluent | LBD |
| Assay Buffer | L-AB |
| 10X Wash Buffer | L-WB |
| Set of two 96-Well plates with sealers | MAG-PLATE |
| Human Th17 25 Plex Premixed Magnetic Bead Panel – BULK PACKAGING | HT17MAG14PMX25BK |

Human Th17 Antibody Immobilized Premixed Magnetic Beads

| | |
|--------------------------------------|------------------------------|
| Premixed 24-plex beads + HCYIFNG-MAG | HTH17PMX24-MAG + HCYIFNG-MAG |
|--------------------------------------|------------------------------|

Antibody-Immobilized Magnetic Beads

| <u>Analyte</u> | <u>Bead #</u> | <u>Cat. #</u> |
|---------------------|---------------|---------------|
| IL-17F | 19 | HIL17F-MAG |
| GM-CSF | 20 | HGMCSF-MAG |
| IFN γ | 25 | HCYIFNG-MAG |
| IL-10 | 27 | HCYIL10-MAG |
| CCL20/MIP3 α | 28 | HMIP3A-MAG |
| IL-12p70 | 33 | HIL12P70-MAG |
| IL-13 | 35 | HIL13-MAG |
| IL-15 | 37 | HIL15-MAG |
| IL-17A | 39 | HIL17-MAG |
| IL-22 | 43 | HIL22-MAG |
| IL-9 | 45 | HIL9-MAG |
| IL-1 β | 46 | HCYIL1B-MAG |
| IL-33 | 47 | HTHIL33-MAG |
| IL-2 | 48 | HIL2-MAG |
| IL-21 | 52 | HIL21-MAG |
| IL-4 | 53 | HIL4-MAG |
| IL-23 | 54 | HIL23-MAG |
| IL-5 | 55 | HIL5-MAG |
| IL-6 | 57 | HCYIL6-MAG |
| IL-17E/IL-25 | 66 | HIL17E-MAG |
| IL-27 | 72 | HIL27-MAG |
| IL-31 | 73 | HIL31-MAG |
| TNF α | 75 | HCYTNFA-MAG |
| TNF β | 76 | HTNFB-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

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.

WELL MAP

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