

User Guide

MILLIPLEX® Mouse Adipokine Magnetic Bead Panel

96-Well Plate Assay

MADKMAG-71K

Introduction	2	Immunoassay Procedure	12
Principle	3	Plate Washing	14
Storage Conditions Upon Receipt	4	Equipment Settings	14
Reagents Supplied.....	4	Quality Controls	16
Materials Required.....	5	Assay Characteristics	16
Safety Precautions	6	Cross-Reactivity	16
Symbol Definitions.....	6	Assay Sensitivities	16
Technical Guidelines	7	Precision	17
Sample Collection and Storage	8	Accuracy	17
Preparation of Reagents for		Troubleshooting	18
Immunoassay	9	Product Ordering	22
Preparation of		Well Map.....	23
Antibody-Immobilized Beads	9	Notice	24
Preparation of Quality Controls	10	Technical Assistance	24
Preparation of Wash Buffer	10	Terms and Conditions of Sale	24
Preparation of Serum Matrix	10	Safety Data Sheets (SDS).....	24
Preparation of Mouse		Contact Information.....	24
Adipokine Standard	10		
Preparation of Standards	11		

Introduction

In the past, the role of white adipose tissue was thought to be limited to energy storage and internal organ protection. The discovery of leptin secretion from adipocytes in 1994 led to the recognition that white adipose tissue is involved in a variety of metabolic and physiological processes. Adipocytes secrete a number of hormones called adipokines with functions that include appetite and energy balance, insulin sensitivity and lipid metabolism. One of these adipokines, adiponectin, is involved in the regulation of lipid and glucose metabolism, influencing the body's response to insulin. Also, its anti-inflammatory effects on the cellular lining of blood vessel walls may help explain the association of high adiponectin levels with the reduced risk of heart attack.

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
- 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® Mouse Adipokine Magnetic Bead Panel thus enables you to focus on the therapeutic potential of adipocytes. 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.

The MILLIPLEX® Mouse Adipokine Magnetic Bead Panel is part of the most versatile system available for metabolic disease 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® products offer you:

- The ability to choose any combination of analytes from our panel of 7 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.

The MILLIPLEX® Mouse Adipokine Magnetic Bead Panel is a 7-plex kit to be used for the simultaneous quantification of any or all of the following analytes in mouse serum and plasma samples: IL-6, Insulin, Leptin, MCP-1, PAI-1 Total, Resistin and TNF α .

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® products 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®-C microspheres.

- Luminex® products use 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®), 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 the new 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

Store all reagents at 2-8 °C.

Reagents	Volume	Quantity	Cat. No.
Mouse Adipokine Standard	Lyophilized	1 vial	LMA-8071
Mouse Adipokine Quality Controls 1 and 2	Lyophilized	2 vials	LMA-6071
Serum Matrix	Lyophilized	1 vial	LMC-SD
Bead Diluent	3.5 mL	1 vial	LE-BD
Set of one 96-Well Plate with 2 sealers	-	1 set	-
Assay Buffer	30 mL	1 bottle	LE-ABGLP
10X Wash Buffer	60 mL	1 bottle	L-WB
Mouse Adipokine Detection Antibodies	5.5 mL	1 bottle	LMAK-1071
Streptavidin-Phycoerythrin	5.5 mL	1 bottle	L-SAPE7
Mixing Bottle	-	1 bottle	-

Included Mouse Adipokine Panel Antibody-Immobilized Beads are dependent on customizable selection of analytes within the panel.

Mouse Adipokine Panel Antibody-Immobilized Magnetic Beads

Bead/Analyte Name	Luminex® Magnetic Bead Region	Customizable 7 Analytes 20X concentration, 200 µL	
		Available	Cat. No.
Anti-IL-6	26	✓	MIL6-MAG
Anti-Insulin	37	✓	RMINS-MAG
Anti-Leptin	38	✓	RMLPTN-MAG
Anti-MCP-1	56	✓	RMMCP1-MAG
Anti-PAI-1 Total	61	✓	MTPAI1-MAG
Anti-Resistin	64	✓	MRES-MAG
Anti-TNFα	65	✓	RMTNFA-MAG

Materials Required (Not included)

Reagents

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

Instrumentation/Materials





- Adjustable pipettes with tips capable of delivering 10 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 B200 or equivalent)
- Titer plate shaker (VWR® Microplate Shaker Cat. No. 12620-926 or equivalent)
- Luminex® 200™, HTS, FLEXMAP 3D®, MAGPIX® instrument with xPONENT® software, or xMAP® INTELLIFLEX instrument with INTELLIFLEX software by Luminex® Corporation
- Automatic plate washer for magnetic beads (BioTek® 405 LS and 405 TS, Cat. No. 40-094, 40-095, 40-096, 40-097 or equivalent) or Handheld Magnetic Separation Block (Cat. 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 (Cat. No. MX-PLATE) to run the assay using a vacuum filtration unit (Vacuum Manifold Cat. No. MSVMHTS00 or equivalent with Vacuum Pump Cat. 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 has been added to some reagents as a preservative. Although the concentrations are low, Sodium azide 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	
Mouse Adipokine Quality Controls 1 & 2	LMA-6071		Warning. Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Mouse Adipokine Standard	LMA-8071		Warning. Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Serum Matrix	LMC-SD	No Symbol Required	Harmful to aquatic life with long lasting effects. Avoid release to the environment.
Streptavidin-Phycoerythrin	L-SAPE7		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.

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 should be discarded after use.
- 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 the Luminex® 200™ instrument, 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 the MAGPIX® instrument, 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 the FLEXMAP 3D® instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 1 alignment disc.
- For the FLEXMAP 3D® instrument, 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 the xMAP® INTELLIFLEX instrument, 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 serum/plasma samples that require further dilution beyond neat, use the matrix provided in the kit.
- 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 Serum Samples

- Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000 x g. Remove serum and assay immediately or aliquot and store samples at ≤ -20 °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.
- Serum samples should be tested neat. However, if dilution is required, use the Serum Matrix provided in the kit as the diluent for serum samples requiring dilution prior to assay.

Preparation of Plasma Samples

- Plasma collection using EDTA as an anti-coagulant is recommended. Centrifuge for 10 minutes at 1000 x *g* within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at ≤ -20 °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.
- Plasma samples should be tested neat. However, if dilution is required, use the Serum Matrix provided in the kit as the diluent for plasma samples requiring dilution prior to assay.

NOTE:

- A maximum of 10 μ L per well of neat serum or plasma can 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

Preparation of Antibody-Immobilized Beads

Prepare the bead mixture right before setting up the assay. Early preparation of the bead mixture should be avoided. For individual vials of beads, sonicate each antibody-bead vial for 30 seconds; vortex for 1 minute. Add 150 μ 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 of the mixed beads should be discarded after use. Left-over individual beads can 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 4 antibody-immobilized beads, add 150 μ L from each of the 4 bead vials to the Mixing Bottle. Then add 2.4 mL Bead Diluent.

Example 2: When using 7 antibody-immobilized beads, add 150 μ L from each of the 7 bead vials to the Mixing Bottle. Then add 1.95 mL Bead Diluent.

Preparation of Quality Controls

Before use, reconstitute Quality Control 1 and Quality Control 2 with 250 μ L 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 ≤ 20 $^{\circ}$ 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 unused portion at 2-8 $^{\circ}$ C for up to one month.

Preparation of Serum Matrix

This step is required for serum or plasma samples only.

Add 1 mL deionized water to the bottle containing lyophilized Serum Matrix. Mix well. Allow at least 5-10 minutes for complete reconstitution. Leftover reconstituted Serum Matrix should be stored at ≤ -20 $^{\circ}$ C for up to one month.

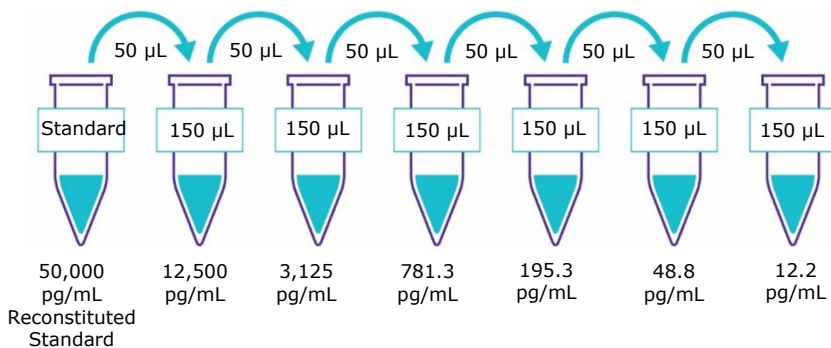
Preparation of Mouse Adipokine Standard

1. Prior to use, reconstitute the Mouse Adipokine Standard with 250 μ L deionized water to give a 50,000 pg/mL concentration of standard stock. 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. The unused portion may be stored at ≤ -20 $^{\circ}$ C for up to one month.
2. Preparation of Working Standards
Label six polypropylene microfuge tubes 12,500 pg/mL, 3,125 pg/mL, 781.3 pg/mL, 195.3 pg/mL, 48.8 pg/mL and 12.2 pg/mL. Add 150 μ L of Assay Buffer to each of the six tubes. Prepare 1:4 serial dilutions by adding 50 μ L of the 50,000 pg/mL reconstituted standard to the 12,500 pg/mL tube, mix well and transfer 50 μ L of the 12,500 pg/mL standard to the 3,125 pg/mL tube, mix well and transfer 50 μ L of the 3,125 pg/mL standard to the 781.3 pg/mL tube, mix well and transfer 50 μ L of the 781.3 pg/mL standard to the 195.3 pg/mL tube, mix well and transfer 50 μ L of the 195.3 pg/mL standard to the 48.8 pg/mL tube, mix well and transfer 50 μ L the 48.8 pg/mL standard to the 12.2 pg/mL tube and mix well. The 0 pg/mL standard (Background) will be Assay Buffer.

Standard Concentration (pg/mL)	Add Deionized Water (μL)	Add Standard (volume)
50,000	250	0

Standard Concentration (pg/mL)	Add Assay Buffer (μL)	Add Standard (volume)
12,500	150	50 μL of 50,000 pg/mL
3,125	150	50 μL of 12,500 pg/mL
781.3	150	50 μL of 3,125 pg/mL
195.3	150	50 μL of 781.3 pg/mL
48.8	150	50 μL of 195.3 pg/mL
12.2	150	50 μL of 48.8 pg/mL

Preparation of Standards



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), 12.2, 48.8, 195.3, 781.3, 3,125, 12,500, 50,000 pg/mL], 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 μ L 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 10 μ L of each Standard or Control into the appropriate wells. Assay Buffer should be used for 0 pg/mL standard (Background).
4. Add 10 μ L of Assay Buffer to the sample wells.
5. Add 10 μ L of Serum Matrix to the background, standards, and control wells.
6. Add 10 μ L of Sample (neat) into the appropriate wells.
7. Vortex Mixing Bottle and add 25 μ L of the Mixed 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 2-8 °C.

Add 200 μ L Assay Buffer per well



Shake 10 min, RT
Decant

- Add 10 μ L Standard or Control to appropriate wells
- Add 10 μ L Assay Buffer to background and sample wells
- Add 10 μ L Serum Matrix to background, standards, and control wells
- Add 10 μ L neat Samples to sample wells
- Add 25 μ L Beads to each well



Incubate overnight at 2-8 °C

9. Gently remove well contents and wash plate 3 times following instructions listed in the Plate Washing section.
10. Add 50 μL 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 30 minutes at room temperature (20-25 $^{\circ}\text{C}$). **DO NOT ASPIRATE AFTER INCUBATION.**
12. Add 50 μL Streptavidin-Phycoerythrin to each well containing the 50 μL of Detection Antibodies.
13. Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25 $^{\circ}\text{C}$).
14. Gently remove well contents and wash plate 3 times following instructions listed in the Plate Washing section.
15. Add 100 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.
16. Run plate on Luminex[®] 200™, HTS, FLEXMAP 3D[®], MAGPIX[®] instrument with xPONENT[®] software or xMAP[®] INTELLIFLEX instrument with INTELLIFLEX 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.



Remove well contents and wash 3X with 200 μL Wash Buffer

Add 50 μL Detection Antibodies per well



Incubate 30 minutes at RT
Do Not Aspirate

Add 50 μL Streptavidin-Phycoerythrin per well



Incubate 30 minutes at RT
Remove well contents and wash 3X with 200 μL Wash Buffer

Add 100 μL Sheath Fluid PLUS or Drive Fluid PLUS per well

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

Plate Washing

Solid Plate

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

- Handheld magnet (Cat. 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 (Cat. No. 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 (Cat. 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® instruments with xPONENT® software and xMAP® INTELLIFLEX instrument 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 (Cat. No. LX2RCAL-K25)	Performance Verification Kit (Cat. No. LX2RPVER-K25)
FLEXMAP 3D®	FLEXMAP 3D® Calibrator Kit (Cat. No. F3DCAL-K25)	FLEXMAP 3D® Performance Verification Kit (Cat. No. F3DPVER-K25)
xMAP® INTELLIFLEX	xMAP® INTELLIFLEX Calibration Kit (Cat. No. IFX-CAL-K20)	xMAP® INTELLIFLEX Performance Verification Kit (Cat. No. IFX-PVER-K20)
MAGPIX®	MAGPIX® Calibration Kit (Cat. No. MPXCAL-K25)	MAGPIX® Performance Verification Kit (Cat. No. MPXPVER-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 Cat. No. MAG-PLATE, if additional plates are required for this purpose.

Events	50, per bead	
Sample Size	50 µL	
Gate Settings	8,000 to 15,000	
Reporter Gain	Default (low PMT)	
Time Out	60 seconds	
Bead Set	<u>Customizable 7-Plex Beads</u>	
	IL-6	26
	Insulin	37
	Leptin	38
	MCP-1	56
	PAI-1 Total	61
	Resistin	64
	TNFα	65

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 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.

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 (n=8).

Analyte	MinDC (pg/mL)	MinDC+2SD (pg/mL)
IL-6	2.3	6.3
Insulin	13.0	27.7
Leptin	4.2	8.2
MCP-1	4.9	11.9
PAI-1 Total	4.0	10.3
Resistin	1.1	3.7
TNF α	5.3	11.0

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	Intra-assay %CV	Inter-assay %CV
IL-6	< 10%	< 20%
Insulin	< 10%	< 20%
Leptin	< 10%	< 20%
MCP-1	< 10%	< 20%
PAI-1 Total	< 10%	< 20%
Resistin	< 10%	< 20%
TNF α	< 10%	< 20%

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	% Recovery in Serum Matrix
IL-6	89%
Insulin	89%
Leptin	86%
MCP-1	83%
PAI-1 Total	92%
Resistin	77%
TNF α	85%

Troubleshooting

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 flush and washes; or if needed probe should be removed and sonicated.
	Probe height not adjusted correctly	When reading the assay on the Luminex® 200™ instrument, adjust probe height to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on the MAGPIX® instrument, adjust probe height to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on the FLEXMAP 3D® instrument, adjust probe height to the kit solid plate using 1 alignment disc. For the FLEXMAP 3D® instrument, 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 the xMAP® INTELLIFLEX instrument, 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®.
Background is too high	Background wells were contaminated	Avoid cross-well contamination by using sealer appropriately and pipetting with Multichannel pipets 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.

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

Problem	Probable Cause	Solution
Beads not in region or gate	Luminex® not calibrated correctly or recently	Calibrate Luminex® based on Instrument 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, Bioplex) 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® 4 times to rid of air bubbles, wash 4 times with sheath fluid or water if there is any remnant alcohol or sanitizing liquid.
Signal for whole plate is same as background	Beads were exposed to light	Keep plate and bead mix covered with dark lid or aluminum foil during all incubation steps.
	Incorrect or no Detection Antibody was added	Add appropriate Detection Antibody and continue.
Low signal for standard curve	Streptavidin-Phycoerythrin was not added	Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.
	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® Instrument (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.

Problem	Probable Cause	Solution
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 tech 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.
High Variation in samples and/or standards	Multichannel pipet may not be calibrated	Calibrate pipets.
	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 a vertical 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 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 setup 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

Replacement Reagents	Cat. No.
Mouse Adipokine Standard	LMA-8071
Mouse Adipokine Quality Controls	LMA-6071
Serum Matrix	LMC-SD
Mouse Adipokine Detection Antibodies	LMAK-1071
Bead Diluent	LE-BD
Streptavidin-Phycoerythrin	L-SAPE7
Assay Buffer	LE-ABGLP
Set of two 96-Well plates with sealers	MAG-PLATE
10X Wash Buffer	L-WB

Antibody-Immobilized Magnetic Beads

Analyte	Bead No.	Cat. No.
IL-6	26	MIL6-MAG
Insulin	37	RMINS-MAG
Leptin	38	RMLPTN-MAG
MCP-1	56	RMMCP1-MAG
PAI-1 Total	61	MTPAI1-MAG
Resistin	64	MRES-MAG
TNF α	65	RMTNFA-MAG

Well Map

	1	2	3	4	5	6	7	8	9	10	11	12
A	0 pg/mL Standard (Background)	781.3 pg/mL Standard	QC-1 Control	Etc.								
B	0 pg/mL Standard (Background)	781.3 pg/mL Standard	QC-1 Control	Etc.								
C	12.2 pg/mL Standard	3,125 pg/mL Standard	QC-2 Control									
D	12.2 pg/mL Standard	3,125 pg/mL Standard	QC-2 Control									
E	48.8 pg/mL Standard	12,500 pg/mL Standard	Sample 1									
F	48.8 pg/mL Standard	12,500 pg/mL Standard	Sample 1									
G	195.3 pg/mL Standard	50,000 pg/mL Standard	Sample 2									
H	195.3 pg/mL Standard	50,000 pg/mL Standard	Sample 2									

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.

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

Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://www.SigmaAldrich.com/techservice).

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://www.SigmaAldrich.com/terms).

Safety Data Sheets (SDS)

Safety Data Sheets are available on the product page at [SigmaAldrich.com](https://www.SigmaAldrich.com).

Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.SigmaAldrich.com/offices).

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

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

© 2012-2025 Merck and/or its affiliates.
All Rights Reserved.
Template 0001187FM v. 20.0

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

00003770MAN Rev 11/25

