

User Guide

MILLIPLEX® Rat Kidney Toxicity Magnetic Bead Panel 1

96 Well Plate Assay

RKTX1MAG-37K

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

Introduction

Absorption, distribution, metabolism and excretion (ADME) parameters are critical to all stages of a fully integrated drug development program and are used to augment the interpretation of toxicological findings. As toxicity has been found to be the leading cause of drug failure, this area of research is expanding in search of more sensitive, rapid methods for determining organ-specific damage as quickly as possible. Drug-induced damage to kidney cells, also known as renal toxicity, results from drug excretion. The traditional methodology for determining renal toxicity has been to measure the blood urea nitrogen (BUN) and serum creatinine. These two tests only detect kidney damage a week after it begins to occur and only show that damage has occurred somewhere in the kidney.

The MILLIPLEX® portfolio offer 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® Rat Kidney Toxicity Magnetic Bead Panel 1 thus enables you to focus on the therapeutic potential of drug induced renal toxicity. 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® Rat Kidney Toxicity Magnetic Bead Panel 1 is part of the most versatile system available for kidney toxicity 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® Rat Kidney Toxicity Magnetic Bead Panel 1 is a 7-plex kit to be used for the simultaneous quantification of any or all of the following analytes in urine samples: Clusterin, Glutathione S-Transferase alpha (GSTα), IP-10, Kidney Injury Molecule-1 (KIM-1), Osteopontin (OPN), Tissue Inhibitor of Metalloproteinase 1 (TIMP-1), and VEGF.

**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. |
|--|-------------|----------|----------------|
| Rat Kidney Toxicity Panel 1 Standard | Lyophilized | 1 vial | RKTXMG1-8037-1 |
| Rat Kidney Toxicity Panel 1 Quality Controls 1 and 2 | Lyophilized | 2 vials | RKTXMG1-6037-1 |
| Serum Matrix* | Lyophilized | 1 | HSCR-SM |
| Set of one 96-Well Plate with 2 sealers | - | 1 set | - |
| Assay Buffer | 30 mL | 1 bottle | L-AB1 |
| 10X Wash Buffer** | 60 mL | 1 bottle | L-WB |
| Bead Diluent | 4.0 mL | 1 bottle | LA-BD |
| Rat Kidney Toxicity Panel 1 Biotinylated GSTa (only supplied if GSTa is a requested analyte) | 3.2 mL | 1 bottle | RKTX1-BTGST |
| Rat Kidney Toxicity Panel 1 Detection Antibodies | 5.5 mL | 1 bottle | RKTXMG1-1037-1 |
| Streptavidin-Phycoerythrin | 5.5 mL | 1 bottle | L-SAPE7 |
| Mixing Bottle | - | 1 bottle | - |

* Contains 0.08% Sodium azide

** Contains 0.05% Proclin

Included Rat Kidney Toxicity Magnetic Bead Panel 1 Antibody-Immobilized Beads are dependent on customizable selection of analytes within the panel.

Rat Kidney Toxicity Magnetic Bead Panel 1
Antibody-Immobilized Magnetic Beads:

| Bead/Analyte Name | Luminex® Magnetic Bead Region | Customizable 7 Analytes (20X concentration, 200 µL) | |
|-------------------------|-------------------------------------|--|-------------|
| | | Available | Cat. No. |
| Anti-Rat Clusterin Bead | 46 | ✓ | RCLSTRN-MAG |
| Anti-Rat GSTα Bead | 51 | ✓ | RGSTA-MAG |
| Anti-Rat IP-10 Bead | 53 | ✓ | RKTIP10-MAG |
| Anti-Rat KIM-1 Bead | 55 | ✓ | RKIM1-MAG |
| Anti-Rat OPN Bead | 61 | ✓ | ROPN-MAG |
| Anti-Rat TIMP-1 Bead | 74 | ✓ | RTIMP1-MAG |
| Anti-Rat VEGF Bead | 76 | ✓ | RVEGF-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 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 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 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 | |
|--|----------------|---|--|
| Rat Kidney Toxicity Panel 1 Biotinylated GSTa | RKTX1-BTGST |  | 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. |
| Rat Kidney Toxicity Panel 1 Detection Antibodies | RKTXMG1-1037-1 |  | 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. |
| Rat Kidney Toxicity Panel 1 Quality Controls 1 & 2 | RKTXMG1-6037-1 |    | Danger. Harmful if swallowed. Causes serious eye damage. Toxic 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. |
| Rat Kidney Toxicity Panel 1 Standard | RKTXMG1-8037-1 |    | Danger. Harmful if swallowed. Causes serious eye damage. Toxic to aquatic life with long lasting effects. Avoid release to the environment. Wear eye protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/attention. |
| Serum Matrix | HSCR-SM | No Symbol Required | Harmful to aquatic life with long lasting effects. Avoid release to the environment. |
| Assay Buffer | L-AB1 |  | 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. |
| 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 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 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 urine samples that require further dilution beyond 1:2, use the serum matrix provided in the kit.
- Vortex all reagents well before adding to plate.

Sample Collection and Storage

Preparation of Urine Samples

- Centrifuge samples briefly to pellet debris. 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.
- Urine samples do not require dilution.

NOTE:

- A maximum of 12.5 µL per well of neat urine 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

For individual vials of beads, sonicate each antibody-bead vial for 30 seconds; vortex for 1 minute. Add 150 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 °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 3 antibody-immobilized beads, add 150 µL from each of the 3 bead vials to the Mixing Bottle. Then add 2.55 mL Bead Diluent.

Example 2: When using 7 antibody-immobilized beads, add 150 µL 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

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

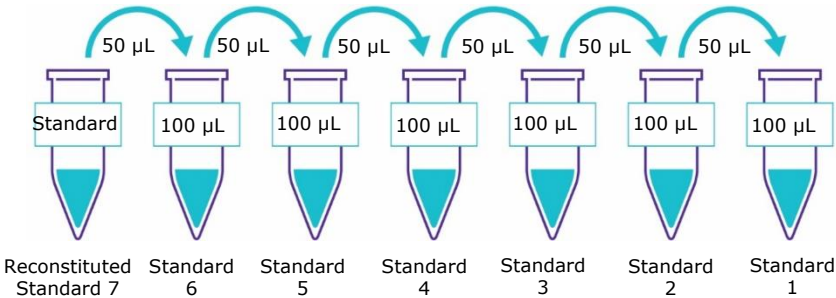
Preparation of Rat Kidney Toxicity Magnetic Bead Panel 1 Standard

1. Prior to use, reconstitute the Rat Kidney Toxicity Panel 1 Standard with 250 μ L 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 7; the unused portion may be stored at ≤ -20 $^{\circ}$ C for up to one month.
2. Preparation of Working Standards
Label 6 polypropylene microfuge tubes Standard 1 through Standard 6. Add 100 μ L of Assay Buffer to each of the 6 tubes. Prepare serial dilutions by adding 50 μ L of the reconstituted Standard 7 to the Standard 6 tube, mix well and transfer 50 μ L of Standard 6 to the Standard 5 tube, mix well and transfer 50 μ L of Standard 5 to the Standard 4 tube, mix well and transfer 50 μ L of Standard 4 to the Standard 3 tube, mix well and transfer 50 μ L of Standard 3 to the Standard 2 tube, mix well and transfer 50 μ L of Standard 2 to the Standard 1 tube and mix well. The 0 ng/mL standard (Background) will be Assay Buffer.

| Standard Tube No. | Add Deionized Water (μ L) | Add Standard (volume) |
|-------------------|--------------------------------|-----------------------|
| Standard 7 | 250 | 0 |

| Standard Tube No. | Add Assay Buffer (μ L) | Add Standard (volume) |
|-------------------|-----------------------------|--------------------------|
| Standard 6 | 100 | 50 μ L of Standard 7 |
| Standard 5 | 100 | 50 μ L of Standard 6 |
| Standard 4 | 100 | 50 μ L of Standard 5 |
| Standard 3 | 100 | 50 μ L of Standard 4 |
| Standard 2 | 100 | 50 μ L of Standard 3 |
| Standard 1 | 100 | 50 μ L of Standard 2 |

Preparation of Standards



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

| Standard | Clusterin (ng/mL) | GSTα (ng/mL) |
|------------|-------------------|--------------|
| Standard 1 | 2.06 | 4.12 |
| Standard 2 | 6.17 | 12.35 |
| Standard 3 | 18.52 | 37.04 |
| Standard 4 | 55.56 | 111.11 |
| Standard 5 | 166.67 | 333.33 |
| Standard 6 | 500.00 | 1,000.00 |
| Standard 7 | 1,500.00 | 3,000.00 |

| Standard | IP-10, KIM-1, OPN, VEGF (ng/mL) | TIMP-1 (ng/mL) |
|------------|---------------------------------|----------------|
| Standard 1 | 0.03 | 0.27 |
| Standard 2 | 0.08 | 0.82 |
| Standard 3 | 0.25 | 2.47 |
| Standard 4 | 0.74 | 7.41 |
| Standard 5 | 2.22 | 22.22 |
| Standard 6 | 6.67 | 66.67 |
| Standard 7 | 20.00 | 200.00 |

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), Standards 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 ng/mL standard (Background).
 4. Add 25 μL of Serum Matrix to the background, standards, and control wells.
 5. Add 25 μL of Assay Buffer to the sample wells.
 6. Add 12.5 μL of neat Sample to the sample wells.
 7. Add 12.5 μL of Serum Matrix into the sample wells.
 8. If measuring GST α , add 25 μL of Biotinylated GST α to each well. If GST α is not one of the selected analytes, add 25 μL of Assay Buffer instead to each well.
 9. Vortex Mixing Bottle and add 25 μL of the Mixed Beads to each well.
(**Note:** During addition of Beads, shake bead bottle intermittently to avoid settling.)

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 Serum Matrix to background, standards, and control wells
- Add 12.5 μL neat Samples to sample wells
- Add 12.5 μL Serum Matrix to sample wells
- Add 25 μL Biotinylated GST α (if GST α is a selected analyte) or 25 μL Assay Buffer (if GST α is excluded) to each well
- Add 25 μL Beads to each well

10. 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.
11. Gently remove well contents and wash plate 2 times following instructions listed in the Plate Washing section.
12. Add 50 µL of Detection Antibodies into each well.
(**Note:** Allow the Detection Antibodies to warm to room temperature prior to addition.)
13. 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.**
14. Add 50 µL Streptavidin-Phycoerythrin to each well containing the 50 µL of Detection Antibodies.
15. Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25 °C).
16. Gently remove well contents and wash plate 2 times following instructions listed in the Plate Washing section.
17. Add 125 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.
18. Run plate on Luminex® 200™, HTS, FLEXMAP 3D®, MAGPIX® instrument with xPONENT® software or xMAP® INTELLIFLEX instrument with INTELLIFLEX software.
19. 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.)



Incubate overnight
at 4 °C



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

Add 50 µL Detection
Antibodies per well



Incubate 1 hour at
RT

Do Not Aspirate

Add 50 µL
Streptavidin-Phycoerythrin
per well



Incubate for
30 minutes at RT

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

Add 125 µL Sheath Fluid
PLUS or Drive Fluid PLUS
per well

Read on Luminex® (100 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. LX2R-CAL-K25) | Performance Verification Kit (Cat. No. LX2R-PVER-K25) |
| FLEXMAP 3D® | FLEXMAP 3D® Calibrator Kit (Cat. No. F3D-CAL-K25) | FLEXMAP 3D® Performance Verification Kit (Cat. No. F3D-PVER-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. MPX-CAL-K25) | MAGPIX® Performance Verification Kit (Cat. 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 Cat. No. 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 Gain | Default (low PMT) |
| Time Out | 60 seconds |
| Bead Set | Customizable 7-Plex Beads |
| | Clusterin 46 |
| | GSTα 51 |
| | IP-10 53 |
| | KIM-1 55 |
| | OPN 61 |
| | TIMP-1 74 |
| | VEGF 76 |

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.

Assay Sensitivities (minimum detectable concentrations, ng/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.

Overnight Protocol (n = 13 Assays) 2 Hour Protocol (n = 5 Assays)

| Analyte | MinDC (ng/mL) | MinDC+2SD (ng/mL) | MinDC (ng/mL) | MinDC+2SD (ng/mL) |
|-----------|------------------|----------------------|------------------|----------------------|
| Clusterin | 0.307 | 0.845 | 0.425 | 1.559 |
| GSTα | 6.137 | 24.177 | 6.306 | 17.533 |
| IP-10 | 0.006 | 0.015 | 0.002 | 0.003 |
| KIM-1 | 0.007 | 0.015 | 0.007 | 0.012 |
| OPN | 0.008 | 0.017 | 0.007 | 0.015 |
| TIMP-1 | 0.140 | 0.216 | 0.131 | 0.214 |
| VEGF | 0.009 | 0.020 | 0.005 | 0.008 |

Precision

Intra-assay precision is generated from the mean of the %CVs from 8 reportable results across two different concentrations of analytes in a single assay. Inter-assay precision is generated from the mean of the %CVs across two different concentrations of analytes across 13 different assays.

| Analyte | Overnight Protocol | | 2 Hour Protocol |
|-----------|--------------------|-----------------|-----------------|
| | Intra-assay %CV | Inter-assay %CV | Intra-assay %CV |
| Clusterin | < 10 | < 15 | < 10 |
| GSTa | < 20 | < 30 | < 25 |
| IP-10 | < 10 | < 15 | < 10 |
| KIM-1 | < 10 | < 15 | < 10 |
| OPN | < 10 | < 15 | < 10 |
| TIMP-1 | < 10 | < 15 | < 10 |
| VEGF | < 10 | < 15 | < 10 |

Accuracy

Spike Recovery: The data represent mean percent recovery of spiked standards ranging from low, medium, and high concentration in urine samples (n=8).

| Analyte | Overnight Protocol | 2 Hour Protocol |
|-----------|-----------------------------|-----------------------------|
| | % Recovery in Urine Samples | % Recovery in Urine Samples |
| Clusterin | 123 | 126 |
| GSTa | 104 | 81 |
| IP-10 | 84 | 98 |
| KIM-1 | 102 | 110 |
| OPN | 81 | 92 |
| TIMP-1 | 121 | 128 |
| VEGF | 101 | 106 |

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 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. |

| Problem | Probable Cause | Solution |
|---|---|---|
| Beads not in region or gate | Luminex® instrument 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. |
| | Beads were exposed to light | Keep plate and bead mix covered with dark lid or aluminum foil during all incubation steps. |
| 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. |

| 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 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 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 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 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. |
|---|-----------------|
| Rat Kidney Toxicity Panel 1 Standard | RKTXMG1-8037-1 |
| Rat Kidney Toxicity Panel 1 Quality Controls 1 and 2 | RKTXMG1-6037-1 |
| Serum Matrix | HSCR-SM |
| Bead Diluent | LA-BD |
| Rat Kidney Toxicity Panel 1 Biotinylated GSTα (only supplied if GSTα is a requested analyte) | RKTX1-BTGST |
| Rat Kidney Toxicity Panel 1 Detection Antibodies | RKTXMG1-1037-1 |
| Streptavidin-Phycoerythrin | L-SAPE7 |
| Assay Buffer | L-AB1 |
| Set of two 96-Well plates with sealers | MAG-PLATE |
| 10X Wash Buffer | L-WB |

Antibody-Immobilized Magnetic Beads

| Analyte | Bead No. | Cat. No. |
|----------------|-----------------|-----------------|
| Clusterin | 46 | RCLSTRN-MAG |
| GSTα | 51 | RGSTA-MAG |
| IP-10 | 53 | RKTIP10-MAG |
| KIM-1 | 55 | RKIM1-MAG |
| OPN | 61 | R0PN-MAG |
| TIMP-1 | 74 | RTIMP1-MAG |
| VEGF | 76 | RVEGF-MAG |

Well Map

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|-------------------------------------|------------|-----------------|------|---|---|---|---|---|----|----|----|
| A | 0 ng/mL Standard (Background) | Standard 4 | QC-1 Control | Etc. | | | | | | | | |
| B | 0 ng/mL Standard (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 | | | | | | | | | |

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®.

Contact Information

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

Technical Assistance

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

Standard Warranty

The applicable warranty for the products listed in this publication 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).

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 Merck KGaA, Darmstadt, Germany and/or its affiliates.
All Rights Reserved.

RKTX1MAG37K Rev 08/21

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