

MILLIPLEX™ MAP

RAT BONE PANEL 3 KIT 96 Well Plate Assay

#RBN3-31K

<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	5
Safety Precautions	5
Technical Guidelines	5
Sample Collection And Storage	7
Preparation of Reagents for Immunoassay	8
Immunoassay Procedure	10
Equipment Settings	12
Quality Controls	12
Assay Characteristics	13
Troubleshooting Guide	14
Replacement Reagents	16
Ordering Information	17
References	18
Well Map	19

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}.

INTRODUCTION

Bone metabolism is the dynamic process of ongoing bone deposition and resorption, controlled by osteoblasts, osteocytes, and osteoclasts. While osteoblasts and osteocytes (osteoblasts surrounded by matrix) are responsible for bone deposition, osteoclasts are responsible for bone resorption. Both are required to maintain bone structure, as well as an adequate supply of calcium. To maintain this metabolic balance these cells rely on complex signaling pathways involving hormones and cytokines to achieve the appropriate rates of growth and differentiation. The disruption of bone metabolism results in such disease as osteoporosis, osteoarthritis, rheumatoid arthritis, chronic kidney disease and bone metastases.

Millipore recognizes the need to better understand the role that bone metabolism biomarkers play both in preserving normal bone structure and in the development of disease. Therefore, we are proud to announce the launch of our MILLIPLEX MAP Rat Bone Metabolism Panel 3. The panel provides biomedical researchers using rat models with quality tools for the study of bone metabolism and its related diseases.

Millipore's MILLIPLEX[™] Rat Bone Panel 3 is a versatile system for bone metabolism research.

- MILLIPLEX[™] MAP offers you the ability to:
 - Choose any combination of analytes from our panel of 3 analytes to design a custom kit that better meets your needs.
- A convenient "all-in-one" box format gives you the assurance that you will have all the necessary reagents you need to run your assay.

Millipore's MILLIPLEX[™] Rat Bone Panel 3 kit is to be used for the simultaneous quantification of the following 3 analytes: PTH, Osteopontin and Osteocalcin. This kit may be used for the analysis of all above analytes in rat tissue, serum, plasma, and cell / tissue culture supernatant samples.

This kit is for research purposes only.

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 beads known as microspheres.

- Luminex® uses proprietary techniques to internally color-code microspheres with two fluorescent dyes. Through precise concentrations of these dyes, 100 distinctly colored bead sets 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 microspheres are allowed to pass rapidly through a laser which excites the internal dyes marking the microsphere set. A second laser excites PE, the fluorescent dye on the reporter molecule.
- Finally, high-speed digital-signal processors identify each individual microsphere and quantify the result of its bioassay based on fluorescent reporter signals.

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.
- **DO NOT STORE RECONSTITUTED STANDARDS OR CONTROLS IN GLASS VIALS.** 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 Antibodies, and Streptavidin-Phycoerythrin.**

REAGENTS SUPPLIED

Note: Store all reagents at 2 – 8 °C

REAGENTS SUPPLIED	CATALOG NUMBER	VOLUME	QUANTITY
Rat Bone Panel 3 Standard	LRBN80313	lyophilized	1 vial
Rat Bone Panel 3 Quality Controls	LRBN60313	lyophilized	2 vials
Set of one 96-Well Filter Plate with 2 Sealers	MX-PLATE	-----	1 plate 2 sealers
Assay Buffer	L-AB1	30 mL	2 bottles
10X Wash Buffer Note: Contains 0.05% Proclin	L-WB	30 mL	1 bottle
Rat Bone Panel 3 Detection Antibody	LRBN10313	3.2 mL	1 bottle
Streptavidin-Phycoerythrin	L-SAPE	5.5 mL	1 bottle
Mixing Bottle	-----	-----	1 bottle

Included Rat Bone Panel 3 Antibody-Immobilized Beads are dependent on customizable selection of analytes within the panel.

Rat Bone Panel 3 Antibody-Immobilized Beads:

Bead/Analyte Name	Luminex Bead Region	(20X concentration, 200 µL)	
		Available	Cat. #
Anti - PTH Bead	14	<input type="checkbox"/>	RPTH
Anti - Osteopontin Bead	18	<input type="checkbox"/>	ROPN
Anti - Osteocalcin Bead	25	<input type="checkbox"/>	ROC

MATERIALS REQUIRED BUT NOT PROVIDED

Reagents

1. Luminex Sheath Fluid (Luminex Catalogue #40-50000)

Instrumentation / Materials

1. Adjustable Pipettes with Tips capable of delivering 25 μ L to 1000 μ L
2. Multichannel Pipettes capable of delivering 5 μ L to 50 μ L or 25 μ L to 200 μ L
3. Reagent Reservoirs
4. Polypropylene Microfuge Tubes
5. Rubber Bands
6. Absorbent Pads
7. Laboratory Vortex Mixer
8. Sonicator (Branson Ultrasonic Cleaner Model #B200 or equivalent)
9. Titer Plate Shaker (Lab-Line Instruments Model #4625 or equivalent)
10. Vacuum Filtration Unit (Millipore Vacuum Manifold Catalog #MSVMHTS00 or equivalent with Millipore Vacuum Pump Catalog #WP6111560 or equivalent)
11. Luminex 100™ IS, 200™, or HTS by Luminex Corporation
12. Plate Stand (Millipore Catalog # MX-STAND)

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. On disposal, flush with a large volume of water to prevent azide build up.

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.

- The bottom of the Microtiter Filter Plate should not be in direct contact with any surface during assay set-up or incubation times. The plate can be set on a plate stand or on the non-flat side of the plate cover or any other plate holder to raise the plate from the surface. A plate stand can be purchased separately from Millipore (Millipore Catalog #MX-STAND).
- Incomplete washing can adversely affect the assay outcome. All washing must be performed with the Wash Buffer provided.
- After the wash steps, keep the bottom of the Microtiter Filter Plate clean by blotting on paper towels or absorbent pads to prevent any leakage due to capillary action.
- Keep the vacuum suction on the plate as low as possible. It is recommended to have a vacuum setting that will remove 200 μ L of buffer in ≥ 5 seconds (equivalent to < 100 mmHg).
- The Standards prepared by serial dilution must be used within 1 hour of preparation. Discard any unused standards except the standard stock which may be stored at $\leq -20^{\circ}\text{C}$ for 1 month and at $\leq -80^{\circ}\text{C}$ for greater than one month.
- If samples fall outside the dynamic range of the assay, further dilute the samples with the appropriate diluent and repeat the assay.
- Any unused mixed Antibody-Immobilized Beads may be stored in the Mixing Bottle at $2-8^{\circ}\text{C}$ 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 $2-8^{\circ}\text{C}$ 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. Adjust probe height according to the protocols recommended by Luminex to the kit filter plate using 3 alignment discs prior to reading an assay.
- 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 the assay buffer diluted serum/plasma samples, use the Assay Buffer as matrix.
- For cell/tissue homogenate, the final cell or tissue homogenate should be prepared in a buffer that has a neutral pH, contains minimal detergents or strong denaturing detergents, and has an ionic strength close to physiological concentration. Avoid debris, lipids, and cell/tissue aggregates. Centrifuge samples before use.
- Vortex all reagents well before adding to plate.

SAMPLE COLLECTION AND STORAGE

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.
- Customers may need to determine the optimal dilution factors for their samples depending on the expected biological range. Dilution of rat serum samples in Assay Buffer at 1:20 dilution is recommended prior to the assay. The Assay Buffer provided in the kit should be used as the sample diluent.

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.
- Customers may need to determine the optimal dilution factors for their samples depending on the expected biological range. Dilution of plasma samples in Assay Buffer at 1:20 dilution is recommended prior to the assay. The Assay Buffer provided in the kit should be used as the sample diluent.

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.
- Customers may need to determine the optimal dilution factors for their samples depending on the expected biological range.

NOTE:

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

PREPARATION OF REAGENTS FOR IMMUNOASSAY

A. Preparation of Antibody-Immobilized Beads

Sonicate each antibody-bead vial for 30 seconds then 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 Assay Buffer. Vortex the mixed beads well. Unused portion may be stored at 2-8°C for up to one month.

Example 1: When using 2 rat bone antibody-immobilized beads, add 150 μ L from each of the 2 bead sets to the Mixing Bottle. Then add 2.70 mL Assay Buffer.

Example 2: When using 3 rat bone antibody-immobilized beads, add 150 μ L from each of the 3 bead sets to the Mixing Bottle. Then add 2.55 mL Assay Buffer.

B. 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. Unused portion may be stored at \leq -20°C for up to one month.

C. Preparation of Wash Buffer

Bring the 10X Wash Buffer to room temperature and mix to bring all salts into solution. Dilute 30 mL of 10X Wash Buffer with 270 mL deionized water. Store unused portion at 2-8°C for up to one month.

D. Preparation of Rat Bone Panel 3 Standard

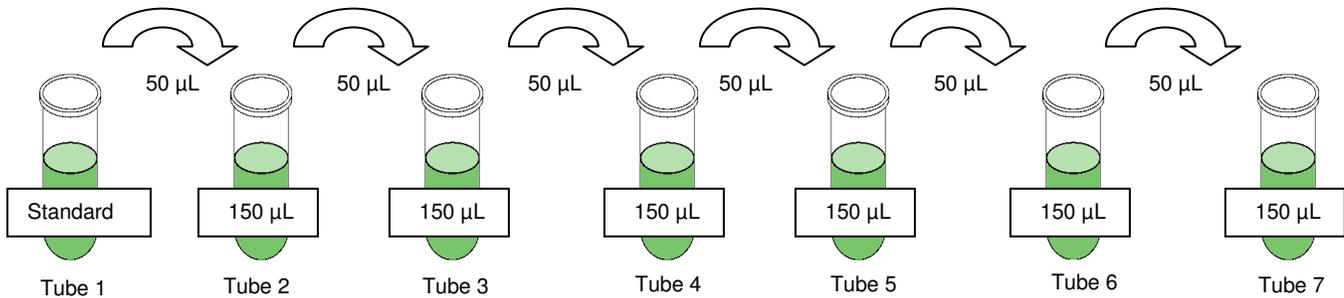
1.) Prior to use, reconstitute the Rat Bone Panel 3 Standard with 250 μ L Deionized Water to give a concentration of 5,000 pg/mL for PTH, 20,000 pg/mL for Osteopontin, and 20,000 pg/mL for Osteocalcin. Invert the vial several times to mix. Allow the vial to set for 5-10 minutes to make sure that the standards are completely reconstituted. This will be used as the Standard 1. The unused portions of this stock may be stored at \leq -20°C for up to one month.

2.) Preparation of Working Standards

Label six polypropylene microfuge tubes as Standard 2, Standard 3, Standard 4, Standard 5, Standard 6, and Standard 7. 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 reconstituted Standard 1 to the Standard 2 tube, mix well and transfer 50 μ L of the Standard 2 to the Standard 3 tube, mix well and transfer 50 μ L of the Standard 3 to the Standard 4 tube, mix well and transfer 50 μ L of the Standard 4 to Standard 5 tube, mix well and transfer 50 μ L of the Standard 5 to the Standard 6 tube, mix well and transfer 50 μ L of the Standard 6 to the Standard 7 tube and mix well. The 0 pg/mL standard (Background) will be the Assay Buffer.

Preparation of Rat Bone Panel 3 Standard

Tube Number	Standard Dilution	Volume of Deionized Water to Add	Volume of Standard to Add
1	Original	250 μ L	0
Tube Number	Standard Dilution	Volume of Assay Buffer to Add	Volume of Standard to Add
2	1:4	150 μ L	50 μ L of Standard Tube 1 (Original Standard)
3	1:16	150 μ L	50 μ L of Standard Tube 2 (1:4 Standard)
4	1:64	150 μ L	50 μ L of Standard Tube 3 (1:16 Standard)
5	1:256	150 μ L	50 μ L of Standard Tube 4 (1:64 Standard)
6	1:1024	150 μ L	50 μ L of Standard Tube 5 (1:256 Standard)
7	1:4096	150 μ L	50 μ L of Standard Tube 6 (1:1024 Standard)



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

Tube Number	Standard Dilution	PTH (pg/mL)	Osteopontin Osteocalcin (pg/mL)
	Background	0	0
7	1:4096	1.22	4.88
6	1:1024	4.88	19.53
5	1:256	19.53	78.13
4	1:64	78.13	312.5
3	1:16	312.5	1,250
2	1:4	1,250	5,000
1	Original	5,000	20,000

IMMUNOASSAY PROCEDURE

- Prior to beginning this assay, it is imperative to read this protocol completely and to thoroughly understand the Technical Guidelines.
- Allow all reagents to warm to room temperature (20-25°C) before use in the assay.
- Diagram the placement of Standards [0 (Background), Tube 7, 6, 5, 4, 3, 2, 1], 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.
- 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. Prewet the filter plate by pipetting 200 μ L of Assay Buffer into each well of the microtiter plate. Seal and mix on a plate shaker for 10 minutes at room temperature (20-25°C).
2. Remove Assay Buffer by vacuum. (**NOTE: DO NOT INVERT PLATE.**) Blot excess Assay Buffer from the bottom of the plate with an absorbent pad or paper towels.
3. Add 25 μ L of each Standard or Control into the appropriate wells. Assay Buffer should be used for the 0 pg/mL standard (Background).
4. Add 25 μ L of Assay Buffer to the sample wells.
5. Add 25 μ L of appropriate matrix solution to the background, standards, and control wells.
 - a. When tissue culture supernatant samples are tested in the analysis, add 25 μ L of control tissue culture medium to the background, standards and control wells.
 - b. When 1:20 Assay Buffer-diluted serum/plasma samples are tested in the analysis, add 25 μ L of Assay Buffer to the background, standards and control wells.
6. Add 25 μ L of Sample (tissue culture or diluted serum/plasma) 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. Add 25 μ L of Rat Bone Panel 3 Detection Antibody into each well.

Add 200 μ L Assay Buffer per well



Shake 10 min, RT

Vacuum

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

9. Seal the plate with a plate sealer then cover it with the opaque lid. Wrap a rubber band around the plate holder, plate and lid and incubate with agitation on a plate shaker overnight (16-20 hours) at 4°C
10. Gently remove fluid by vacuum. **(NOTE: DO NOT INVERT PLATE.)**
11. Wash plate 3 times with 200 µL/well of Wash Buffer, removing Wash Buffer by vacuum filtration between each wash. Blot excess Wash Buffer from the bottom the plate by with an absorbent pad or paper towels.
12. Add 50 µL of Assay Buffer to each well.
13. Add 50 µL Streptavidin-Phycoerythrin to each well containing the 50 µL Assay Buffer.
14. Seal, cover with lid and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25°C).
15. Gently remove all contents by vacuum. **(NOTE: DO NOT INVERT PLATE.)**
16. Wash plate 3 times with 200 µL/well Wash Buffer, removing Wash Buffer by vacuum filtration between each wash. Wipe any excess buffer on the bottom of the plate with a tissue.
17. Add 100 µL of Sheath Fluid to all wells. Resuspend the beads on a plate shaker for 5 minutes.
18. Run plate on Luminex 100™ IS, 200™, or HTS .
19. Save the data and analyze the Median Fluorescent Intensity (MFI) using a 5-parameter logistic or spline curve-fitting method to calculate analyte concentrations in samples.

Incubate overnight at 4°C with shaking



Vacuum and wash 3X with 200 µL Wash Buffer

- Add 50 µL Assay Buffer per well
- Add 50 µL Streptavidin-Phycoerythrin per well

Incubate for 30 minutes at RT



Vacuum and wash 3X with 200 µL Wash Buffer

Add 100 µL Sheath Fluid per well

Read on Luminex (50 µL, 50 beads per bead set)

EQUIPMENT SETTINGS

These specifications are for the Luminex^{100™} IS v.1.7 or Luminex^{100™} IS v2.1/2.2, Luminex^{200™} v2.3, xPONENT, and Luminex HTS. Luminex instruments with other software (e.g. MasterPlex, StarStation, LiquiChip, Bio-Plex, LABScan100) would need to follow instrument instructions for gate settings and additional specifications from the vendors.

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:	3-Plex Beads	
	PTH	14
	Osteopontin	18
	Osteocalcin	25

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 MILLIPORE website www.millipore.com/techlibrary/index.do using the catalog number as the keyword.

ASSAY CHARACTERISTICS

Assay Sensitivities (minimum detectable concentrations, pg/mL)

Minimum Detectable Concentration (MinDC) is calculated by the StatLIA® Immunoassay Analysis Software from Brendan Technologies. 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	MinDC (pg/mL)
PTH	0.3
Osteopontin	0.8
Osteocalcin	1.6

Precision

Intra-assay precision is generated from the mean of the %CV from 16 reportable results across two different concentrations of analyte in a single assay. Inter-assay precision is generated from the mean of the %CV from 2 reportable results across 2 different concentrations of analyte across 6 different assays.

Analyte	Intra-Assay (CV%)	Inter-Assay (CV%)
PTH	3.5	8.1
Osteopontin	3.1	5.4
Osteocalcin	2.9	4.6

Accuracy

Accuracy, defined as the percentage of measured analyte in samples spiked with known concentrations of analyte, was determined by taking the average recovery of 3 levels of analyte in eight independent 1:20 Assay Buffer diluted serum samples.

Analyte	% Recovery
PTH	77.7
Osteopontin	84.8
Osteocalcin	100.0

Cross-reactivity

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

TROUBLESHOOTING GUIDE

Problem	Probable Cause	Solution
Filter plate will not vacuum	Vacuum pressure is insufficient	Increase vacuum pressure such that 0.2mL buffer can be suctioned in 3-5 seconds.
	Samples have insoluble particles	Centrifuge samples just prior to assay set-up and use supernatant. If high lipid concentration, after centrifugation, remove lipid layer and use supernatant.
Insufficient bead count	Sample too viscous	May need to dilute sample.
	Vacuum pressure too high	Adjust vacuum pressure such that 0.2mL buffer can be suctioned in 3-5 seconds.
	Bead mix prepared incorrectly	Sonicate bead vials and vortex just prior to adding to bead mix bottle according to protocol. Agitate bead mix intermittently in reservoir while pipetting into the plate.
Plate leaked	Samples cause interference due to particulate matter or viscosity	See above. Also sample probe may need to be cleaned with alcohol flush, backflush and washes; or, if needed, probe should be removed and sonicated.
	Probe height not adjusted correctly	Adjust probe to 3 alignment discs in well H6.
	Vacuum pressure too high	Adjust vacuum pressure such that 0.2mL buffer can be suctioned in 3-5 seconds. May need to transfer contents to a new (prewetted) plate and continue.
	Plate set directly on table or absorbent towels during incubations or reagent additions	Set plate on plate stand or raised edge so bottom of filter is not touching any surface.
Background is too high	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 well.
	Probe height not adjusted correctly	Adjust probe to 3 alignment discs in well H6.
	Background wells were contaminated	Avoid cross-well contamination by using sealer appropriately and by pipeting with multichannel pipets without touching reagent in plate.
Background is too high	Matrix used has endogenous analyte or interference	Check matrix ingredients for crossreacting components (e.g. interleukin modified tissue culture medium).
	Insufficient washes	Increase number of washes.

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 (e.g. Bio-Plex) require different gate settings than those described in the kit protocol. Use instrument default settings.
	Wrong bead regions in protocol template	Check kit protocol for correct bead regions or analyte selection.
	Incorrect sample type used	Samples containing organic solvents or if highly viscous should be diluted or dialyzed as required.
	Instrument not washed or primed	Prime the Luminex 4 times to eliminate 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
Low signal for standard curve	Incubations done at incorrect 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 (e.g. Bio-Plex) default target setting for RP1 calibrator is set at High PMT. Use low target value for calibration and reanalyze plate.
	Plate incubation was too long with standard curve and samples	Use shorter incubation time.
Sample readings are out of range	Samples contain no or below detectable levels of analyte	If below detectable levels, it may be possible to use higher sample volume. Check with tech support for appropriate protocol modifications.
	Samples contain analyte concentrations higher than highest standard point	Samples may require dilution and reanalysis for 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 vacuumed out 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 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.

REPLACEMENT REAGENTS

Catalog

Rat Bone Panel 3 Standard	LRBN80313
Rat Bone Panel 3 Quality Controls	LRBN60313
Rat Bone Panel 3 Detection Antibody	LRBN10313
Streptavidin-Phycoerythrin	L-SAPE
Assay Buffer	L-AB1
Set of two 96-Well Filter Plates with Sealers	MX-PLATE
10X Wash Buffer	L-WB

Antibody-Immobilized Bead

<u>Analyte</u>	<u>Bead #</u>	<u>Catalog #</u>
PTH Bead	14	RPTH
Osteopontin Bead	18	R0PN
Osteocalcin Bead	25	ROC

ORDERING INFORMATION

To place an order:

To assure the clarity of your custom kit order, please FAX the following information to our customer service department:

- Your name, telephone and/or fax number
- Customer account number
- Shipping and billing address
- Purchase order number
- Catalog number and description of product
- Quantity of kits

FAX: (636) 441-8050

Toll-Free US: (866) 441-8400

(636) 441-8400

Mail Orders: Millipore Corp.
6 Research Park Drive
St. Charles, Missouri 63304 U.S.A.

For International Customers:

To best serve our international customers in placing an order or obtaining additional information about MILLIPLEX™ MAP products, please contact your multiplex specialist or sales representative or email our European Customer Service at customerserviceEU@Millipore.com.

Conditions of Sale

All products are for research use only. They are not intended for use in clinical diagnosis or for administration to humans or animals. All products are intended for *in vitro* use only.

Material Safety Data Sheets (MSDS)

Material Safety Data Sheets for Millipore products may be ordered by fax or phone or through our website at www.millipore.com/techlibrary/index.do

REFERENCES

Ducy P, Desbois C, Boyce B, Pinero G, Story B, Dunstan C, Smith E, Bonadio J, Goldstein S, Gundberg C, Bradley A, Karsenty G. Increased bone formation in osteocalcin-deficient mice. *Nature* 382:448-452 (1996)

Frolik CA, Bryant HU, Black EC, Magee DE, and Chandrasekhar S. Time-dependent changes in biochemical bone markers and serum cholesterol in ovariectomized rats: effects of raloxifene HCl, tamoxifen, estrogen, and alendronate. *Bone* 18: 621-627 (1996)

Khosla S and Kleetkoper M. Biochemical markers of bone turnover. In: Favus MJ Ed "Primer on the Metabolic Bone Diseases and Disorders of Mineral Metabolism. Published by the American Society for Bone and Mineral Research, Washington, DC" 166-172 (2003)

Wada C, Kataoka M, Seto H, Hayashi N, Kido J, Shinohara Y, Nagata T. High-turnover osteoporosis is induced by cyclosporine A in rats. *J Bone Miner Metab* 24: 199-205 (2006)

Noda M and Rodan GA. Transcriptional regulation of osteopontin production in rat osteoblast-like cells by parathyroid hormone. *J Cell Biol* 108:713-718 (1989)

WELL MAP

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