

User Guide

MILLIPLEX® AAV Serotyping Magnetic Bead Panel

96-Well Plate Assay

AAV SERG-230K

Introduction.....	2	Preparation of Antigen-Immobilized Beads.....	10
Principle	3	Preparation of Wash Buffer	10
Storage Conditions Upon Receipt.....	4	Immunoassay Procedure	11
Reagents Supplied.....	4	Plate Washing	13
AAV Serotyping Assay Control		Solid Plate	13
Magnetic Beads.....	4	Filter Plate	13
AAV Serotyping Antigen-Immobilized Magnetic Beads	5	Equipment Settings.....	14
Materials Required.....	6	Assay Characteristics	16
Reagents.....	6	Cross-Reactivity	16
Instrumentation/Materials	6	Precision.....	16
Safety Precautions.....	6	Troubleshooting.....	17
Symbol Definitions.....	7	Product Ordering	21
Technical Guidelines	8	Well Map	22
Sample Collection and Storage.....	9	Notice	23
Preparation of Serum Samples	9	Technical Assistance.....	23
Preparation of Plasma Samples	9	Standard Warranty	23
Preparation of Reagents for Immunoassay	10	Contact Information	23
		Safety Data Sheets (SDS).....	23

Introduction

Adeno-associated virus (AAV) is recognized as a preferred vector for gene therapy due to its infectivity, versatility, and favorable safety profile. The various AAV serotypes exhibit distinct tissue specificities, transduction efficiencies, and immunogenicity, allowing for tailored approaches to gene therapy. While AAV generally has low immunogenicity, exposure to these viruses is common which leads to a high rate of seroprevalence in the general population. Pre-existing antibodies against AAV can challenge a gene therapy's safety and efficacy making it essential to assess seroprevalence and immunogenicity in descriptive preclinical testing.

The MILLIPLEX® AAV Serotyping Magnetic Bead Panel is a research use only kit designed to measure antibodies against six common AAV serotypes to support the critical preclinical assessments needed to advance applications in gene therapy, vaccine development, and disease research.

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, each panel and kit meet stringent manufacturing criteria to ensure batch-to-batch reproducibility. The MILLIPLEX® AAV Serotyping Magnetic Bead Panel thus enables you to focus on the emerging potential of cell and gene therapies. 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® AAV Serotyping Magnetic Bead Panel is part of the most versatile system available for cell and gene therapy 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 anti-AAV antibody detection.

MILLIPLEX® kits offer you:

The ability to choose any combination of analytes from our panel of 6 analytes to design a custom kit that better meets your needs.

A convenient "all-in-one" box format that gives you the assurance that you will have all the necessary reagents you need to run your assay.

MILLIPLEX® AAV Serotyping Magnetic Bead Panel is a 6-plex kit to be used for the simultaneous detection of any or all of the following antibodies in serum and plasma samples: anti-AAV1, anti-AAV2, anti-AAV5, anti-AAV6, anti-AAV8, anti-AAV9.

For this panel, 4 assay control beads are included as part of the base format.

For Research Use Only. Not for Use in Diagnostic Procedures.

Please read entire protocol before use.

It is important to use same assay incubation conditions throughout your study.

Principle

MILLIPLEX® kits are based on the Luminex® xMAP® technology — one of the fastest growing and most respected multiplex technologies offering applications throughout the life sciences and capable of performing a variety of bioassays including immunoassays on the surface of fluorescent-coded magnetic beads known as MagPlex® microspheres.

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

The capability of adding multiple conjugated beads to each sample results in the ability to obtain multiple results from each sample. Open-architecture xMAP® technology enables multiplexing of many types of bioassays reducing time, labor and costs over traditional methods.

Storage Conditions Upon Receipt

- Recommended storage for kit components is 2-8 °C.
- DO NOT FREEZE Antigen-Immobilized Beads and PE-IgG Conjugate.

Reagents Supplied

Reagents	Volume	Quantity	Catalogue No.
Set of one 96-Well Plate with 2 sealers	1 plate 2 sealers	-	-
Assay Buffer	1 bottle	30 mL	L-AB
10X Wash Buffer	1 bottle	60 mL	L-WB
PE-IgG Conjugate	1 bottle	5.5 mL	AAV-PEIGG
Mixing Bottle	1 bottle	-	-

AAV Serotyping Assay Control Magnetic Beads

Bead/Analyte Name	Luminex® Magnetic Bead Region	Control Beads (50X concentration, 90 µL)	
		Available	Catalogue No.
Control 1 Bead	12	✓	CB1-MAG
Control 2 Bead	13	✓	CB2-MAG
Control 3 Bead	14	✓	CB3-MAG
Negative Control Bead	15	✓	NCB-MAG

The 4 Control Beads are to be combined with the 6 Antigen-Immobilized Beads to make up to a 10-plex.

Included AAV Serotyping Antigen-Immobilized Beads are dependent on customizable selection of analytes within the panel (see next page).

AAV Serotyping Antigen-Immobilized Magnetic Beads

Bead/Analyte Name	Luminex® Magnetic Bead Region	Customizable 6 Analytes (50X concentration, 90 µL)	
		Available	Catalogue No.
AAV1 Bead	38	✓	AAV1-MAG
AAV2 Bead	39	✓	AAV2-MAG
AAV5 Bead	51	✓	AAV5-MAG
AAV6 Bead	52	✓	AAV6-MAG
AAV8 Bead	55	✓	AAV8-MAG
AAV9 Bead	56	✓	AAV9-MAG

Materials Required (Not provided)

Reagents

MAGPIX® Drive Fluid PLUS (40-50030), xMAP® Sheath Fluid PLUS (40-50021), or xMAP® Sheath Concentrate PLUS (40-50023).

Instrumentation/Materials





- Adjustable pipettes with tips capable of delivering 25 µL to 1000 µL
- Multichannel pipettes capable of delivering 5 µL to 50 µL, or 25 µL to 200 µL
- Reagent reservoirs
- Polypropylene microfuge tubes
- Rubber bands
- Aluminum foil
- Absorbent pads
- Laboratory vortex mixer
- Sonicator (Branson Ultrasonic Cleaner Model No. B200 or equivalent)
- Titer plate shaker (VWR® Microplate Shaker, 12620-926 or equivalent)
- Luminex® 200™, HTS, FLEXMAP 3D®, MAGPIX® with xPONENT® software or xMAP® INTELLIFLEX with INTELLIFLEX software by Luminex® Corporation
- Automatic plate washer for magnetic beads (BioTek® 405 LS and 405, TS 40-094, 40-095, 40-096, 40-097 or equivalent) or Handheld Magnetic Separation Block (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 (MX-PLATE) to run the assay using a Vacuum Filtration Unit (Vacuum Manifold, MSVMHTS00 or equivalent with Vacuum Pump, 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	
PE-anti-Human IgG Conjugate	AAV-PEIGG	 	<p>Warning. Causes serious eye irritation. May cause damage to organs Respiratory Tract through prolonged or repeated exposure. Do not breathe mist or vapours. Wash skin thoroughly after handling. Wear eye protection/face 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 if you feel unwell. If eye irritation persists: Get medical advice/ attention. Dispose of contents/container to an approved waste disposal plant.</p>
Assay Buffer	L-AB		<p>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.</p>
10X Wash Buffer	L-WB		<p>Warning. May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.</p>

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 Antigen-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.
- Any unused mixed Antigen-Immobilized Beads may be stored in the Mixing Bottle at 2-8 °C for up to one month.
- The plate should be read immediately after the assay is finished. If, however, the plate cannot be read immediately, seal the plate, cover with aluminum foil or an opaque lid, and store the plate at 2-8 °C for up to 24 hours. Prior to reading, agitate the plate on the plate shaker at room temperature for 10 minutes. Delay in reading a plate may result in decreased sensitivity for some analytes.
- The titer plate shaker should be set at a speed to provide maximum orbital mixing without splashing of liquid outside the wells. For the recommended plate shaker, this would be a setting of 5-7 which is approximately 500-800 rpm.
- Ensure that the needle probe is clean. This may be achieved by sonication and/or alcohol flushes.
- When reading the assay on Luminex® 200™, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on MAGPIX®, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on FLEXMAP 3D®, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 1 alignment disc.
For FLEXMAP 3D®, when using the solid plate in the kit, the final resuspension should be with 150 µL Sheath Fluid PLUS in each well and 75 µL should be aspirated.
For xMAP® INTELLIFLEX, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.
- For serum/plasma samples that require further dilution beyond 1:100, use the Assay Buffer provided in the kit.
- 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 $1000 \times g$. Remove serum and assay immediately or aliquot and store samples at $\leq -20 \text{ }^{\circ}\text{C}$.
- 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.
- Serum samples should be diluted 1:100 in Assay Buffer provided in the kit. For example, in a tube, $10 \text{ }\mu\text{L}$ of serum may be combined with $90 \text{ }\mu\text{L}$ of Assay Buffer to make a 1:10 dilution and then $10 \text{ }\mu\text{L}$ of the 1:10 dilution may be combined with $90 \text{ }\mu\text{L}$ of Assay Buffer to make a 1:100 dilution. When further dilution beyond 1:100 is required, use Assay Buffer as the diluent.

Preparation of Plasma Samples

- Plasma collection using EDTA as an anti-coagulant is recommended. Centrifuge for 10 minutes at $1000 \times g$ within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at $\leq -20 \text{ }^{\circ}\text{C}$.
- 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.
- Plasma samples should be diluted 1:100 in Assay Buffer provided in the kit. For example, in a tube, $10 \text{ }\mu\text{L}$ of plasma may be combined with $90 \text{ }\mu\text{L}$ of Assay Buffer to make a 1:10 dilution and then $10 \text{ }\mu\text{L}$ of the 1:10 dilution may be combined with $90 \text{ }\mu\text{L}$ of Assay Buffer to make a 1:100 dilution. When further dilution beyond 1:100 is required, use Assay Buffer as the diluent.

Note:

- A maximum of $25 \text{ }\mu\text{L}$ per well of diluted 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 anti-coagulant since an excess of heparin will provide falsely high values. Use no more than 10 IU heparin per mL of blood collected.

Preparation of Reagents for Immunoassay

Preparation of Antigen-Immobilized Beads

For **individual vials of beads**, sonicate each Antigen-Immobilized Bead vial for 30 seconds; vortex for 1 minute. Add 60 μ L from each Antigen-Immobilized 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.

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:

- When using 3 Antigen-Immobilized Beads and the 4 control beads, add 60 μ L from each of the 7 bead vials to the Mixing Bottle. Then add 2.58 mL Assay Buffer.
- When using 6 Antigen-Immobilized Beads and the 4 control beads, add 60 μ L from each of the 10 bead vials to the Mixing Bottle. Then add 2.40 mL Assay Buffer.

Preparation of Wash Buffer

Bring the 10X Wash Buffer to room temperature and mix to bring all salts into solution. Dilute 60 mL of 10X Wash Buffer with 540 mL deionized water. Store the unused portion at 2-8 °C for up to one month.

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 Background (Assay Buffer) and Samples on Well Map Worksheet in a vertical configuration. It is recommended to run the assay in duplicate.

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

- If using a filter plate, set the filter plate on a plate holder at all times during reagent dispensing and incubation steps so that the bottom of the plate does not touch any surface.

1. Add 200 μ L of Wash Buffer into each well of the plate. Seal and mix on a plate shaker for 10 minutes at room temperature (20-25 °C).

2. Decant Wash Buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times.

3. Add 25 μ L of Assay Buffer to all wells.

4. Add 25 μ L of Sample (diluted) into the appropriate wells.

5. Add an additional 25 μ L of Assay Buffer to the background wells.

6. 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.

7. 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. Alternatively, incubate for 2 hours at room temperature (20-25 °C).

Add 200 μ L Wash Buffer per well



Shake 10 min, RT
Decant

- Add 25 μ L Assay Buffer to all wells
- Add 25 μ L diluted Samples to sample wells
- Add an additional 25 μ L Assay Buffer to background wells
- Add 25 μ L Beads to each well



Incubate overnight (16-18 hours) at 2-8 °C

8. Gently remove well contents and wash plate 3 times following instructions listed in the Plate Washing section.

9. Add 50 μL of AAV PE-IgG Conjugate into each well.

Note: Allow the PE-IgG Conjugate to warm to room temperature prior to addition.


10. Seal, cover with foil and incubate with agitation on a plate shaker for 1 hour at room temperature (20-25 $^{\circ}\text{C}$).

11. Gently remove well contents and wash plate 3 times following instructions listed in the Plate Washing section.

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


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

14. Save and analyze the Median Fluorescent Intensity (MFI) data. A working cut-off for a positive result can be established using MFI data from known negative sample population.



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

Add 50 μL AAV PE-IgG Conjugate per well



Incubate 1 hour at RT

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

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

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

Plate Washing

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

Solid Plate

- Handheld magnet (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 (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 (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™, FLEXMAP 3D®, MAGPIX® with xPONENT® software and xMAP® INTELLIFLEX with INTELLIFLEX software:

These specifications are for the above listed instruments and software. Luminex® instruments with other software (for example, MasterPlex®, StarStation, LiquiChip, Bio-Plex® Manager™, LABScan™100) would need to follow instrument instructions for gate settings and additional specifications from the vendors for reading Luminex® magnetic beads.

For magnetic bead assays, each instrument must be calibrated, and performance verified with the indicated calibration and verification kits.

Instrument	Calibration Kit	Verification Kit
Luminex® 200™	xPONENT® 3.1 compatible Calibration Kit (LX2R-CAL-K25)	Performance Verification Kit (LX2R-PVER-K25)
FLEXMAP 3D®	FLEXMAP 3D® Calibrator Kit (F3D-CAL-K25)	FLEXMAP 3D® Performance Verification Kit (F3D-PVER-K25)
xMAP® INTELLIFLEX	xMAP® INTELLIFLEX Calibration Kit (IFX-CAL-K20)	xMAP® INTELLIFLEX Performance Verification Kit (IFX-PVER-K20)
MAGPIX®	MAGPIX® Calibration Kit (MPX-CAL-K25)	MAGPIX® Performance Verification Kit (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.
- 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 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 6-plex Beads + Control Beads
	<hr/>
Control 1	12
Control 2	13
Control 3	14
Negative Control	15
AAV1	38
AAV2	39
AAV5	51
AAV6	52
AAV8	55
AAV9	56

Assay Characteristics

Cross-Reactivity

Biologically, individuals with pre-existing antibodies against one serotype may exhibit varying degrees of cross-reactivity with other serotypes due to homology and shared epitopes.

For the purposes of assay verification, where possible, control antibodies directed against specific AAV serotypes were used to establish negligible cross-reactivity on the non-specific beads. The lone exception was the AAV1 control antibody used for assay verification. This antibody has known cross-reactivity to AAV6, which was also observed in this kit.

Precision

Intra-assay precision was generated from the mean of the % CV's from 4 samples in a single assay (overnight protocol). Inter-assay precision was generated from the mean of the % CV's across 16 samples across 3 different assays (overnight protocol).

Analyte	Intra-assay %CV	Inter-assay %CV
AAV1	<10	<15
AAV2	<10	<15
AAV5	<10	<15
AAV6	<10	<15
AAV8	<10	<15
AAV9	<10	<15

Troubleshooting

Problem	Probable Cause	Solution
	Plate washer aspirate height set too low	Adjust aspiration height according to manufacturers' instructions.
	Bead mix prepared inappropriately	Sonicate bead vials and vortex just prior to adding to bead mix bottle according to protocol. Agitate bead mix intermittently in reservoir while pipetting this into the plate.
	Samples cause interference due to particulate matter or viscosity	See above. Also sample probe may need to be cleaned with alcohol flushes, back flushes, and washes; or, if needed, probe should be removed and sonicated.
Insufficient bead count	Probe height not adjusted correctly	<p>When reading the assay on Luminex[®] 200™, adjust probe height to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on MAGPIX[®], adjust probe height to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on FLEXMAP 3D[®], adjust probe height to the kit solid plate using 1 alignment disc. For FLEXMAP 3D[®] when using the solid plate in the kit, the final resuspension should be with 150 µL Sheath Fluid PLUS in each well and 75 µL should be aspirated.</p> <p>When reading the assay on xMAP[®] INTELLIFLEX, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex[®].</p>

Problem	Probable Cause	Solution
Background is too high	Background wells were contaminated	Avoid cross-well contamination by using sealer appropriately and pipetting with multichannel pipettes without touching reagent in plate.
	Insufficient washes	Increase number of washes.
Beads not in region or gate	Luminex® instrument not calibrated correctly or recently	Calibrate Luminex® instrument based on manufacturer's instructions, at least once a week or if temperature has changed by >3 °C.
	Gate settings not adjusted correctly	Some Luminex® instruments (for example Bio-Plex®) require different gate settings than those described in the kit protocol. Use instrument default settings.
	Wrong bead regions in protocol template	Check kit protocol for correct bead regions or analyte selection.
	Incorrect sample type used	Samples containing organic solvents or if highly viscous should be diluted or dialyzed as required.
	Instrument not washed or primed	Prime the Luminex® instrument 4 times to rid it of air bubbles, wash 4 times with Sheath Fluid PLUS or water if there is any remnant alcohol or sanitizing liquid.
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 PE-IgG conjugate was added	Add PE-IgG conjugate according to the protocol.

Problem	Probable Cause	Solution
High variation in samples	Multichannel pipette may not be calibrated	Calibrate pipettes.
	Plate washing was not uniform	Confirm all reagents are removed completely in all wash steps.
	Samples may have high particulate matter or other interfering substances	See above.
	Plate agitation was insufficient	Plate should be agitated during all incubation steps using an orbital plate shaker at a speed where beads are in constant motion without causing splashing.
	Cross-well contamination	Check when reusing plate sealer that no reagent has touched sealer. Care should be taken when using same pipette tips that are used for reagent additions and that pipette tip does not touch reagent in plate.
For Filter Plates Only		
Filter plate will not vacuum	Vacuum pressure is insufficient	Increase vacuum pressure such that 0.2 mL buffer can be suctioned in 3-5 seconds.
	Samples have insoluble particles	Centrifuge samples just prior to assay set-up and use supernatant.
	High lipid concentration	After centrifugation, remove lipid layer and use supernatant.

Problem	Probable Cause	Solution
	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.
Plate leaked	Insufficient blotting of filter plate bottom causing wicking	Blot the bottom of the filter plate well with absorbent towels after each wash step.
	Pipette touching plate filter during additions	Pipette to the side of plate.
	Probe height not adjusted correctly	Adjust probe to 3 alignment discs in well H6.
	Sample too viscous	May need to dilute sample.

Product Ordering

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

Replacement Reagents	Catalogue No.
Assay Buffer	L-AB
PE-IgG Conjugate	AAV-PEIGG
Control 1 Bead	CB1-MAG
Control 2 Bead	CB2-MAG
Control 3 Bead	CB3-MAG
Negative Control Bead	NCB-MAG
Set of two 96-Well plates with sealers	MAG-PLATE
10X Wash Buffer	L-WB

Antigen-Immobilized Magnetic Beads

Bead/Analyte Name	Bead No.	Catalogue No.
AAV1 Bead	38	AAV1-MAG
AAV2 Bead	39	AAV2-MAG
AAV5 Bead	51	AAV5-MAG
AAV6 Bead	52	AAV6-MAG
AAV8 Bead	55	AAV8-MAG
AAV9 Bead	56	AAV9-MAG

Well Map

	1	2	3	4	5	6	7	8	9	10	11	12
A	Background	Sample 4										
B	Background	Sample 4										
C	Sample 1	Etc.										
D	Sample 1	Etc.										
E	Sample 2											
F	Sample 2											
G	Sample 3											
H	Sample 3											

Notice

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

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

Technical Assistance

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

Standard Warranty

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

Contact Information

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

Safety Data Sheets (SDS)

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

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

Merck, 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.

© 2025 Merck KGaA, Darmstadt, Germany and/or its affiliates.
All Rights Reserved.
Template 00011187FM v. 19.0

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

23225172 Rev 06/25

The Merck logo is displayed in a bold, blue, sans-serif font. The letters are closely spaced and have a slightly irregular, blocky appearance.