

User Guide

MILLIPLEX[®] Non-Human Primate Pituitary Magnetic Bead Panel

96 Well Plate Assay

NHPPT1MG-46K

Introduction	2	Equipment Settings	14
Principle	3	Quality Controls	15
Storage Conditions Upon Receipt ...	4	Assay Characteristics	16
Reagents Supplied	4	Assay Sensitivities	16
Materials Required	5	Precision	16
Safety Precautions	6	Accuracy	17
Symbol Definitions	6	Cross-Reactivity	17
Technical Guidelines.....	7	Troubleshooting	18
Sample Collection and Storage	8	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

Pituitary hormones and other brain-derived proteins, such as hypothalamus neuropeptides, play very important roles in the regulation of various functions including metabolism, growth, and reproduction. Accurate measurement of these proteins to understand their new biological functions and molecular mechanisms of the functions are crucial. Traditional laboratory methods, such as RIA and ELISA are not able to measure multiple proteins with a small sample volume.

We recognize the integral role that these proteins play and the MILLIPLEX® Non-Human Primate Pituitary Magnetic Bead Panel 1 can be used for the simultaneous measurement of the following analytes in any combination: ACTH, AGRP, CNTF, FSH, GH, LH, and TSH. This multiplex assay can analyze these 7 proteins simultaneously and uses a small sample volume 25 µL.

The MILLIPLEX® Non-Human Primate Pituitary Magnetic Bead Panel 1 provides biomedical researchers with quality tools for the study of reproduction, growth, metabolic homeostasis, and pituitary-related diseases such as acromegaly, growth hormone deficiency, diabetes insipidus and pituitary tumors.

The MILLIPLEX® portfolio offers the broadest selection of analytes across a wide range of disease states and species. Once the analytes of interest have been identified, you can rely on the quality that we build into each kit to produce results you can trust. In addition to the assay characteristics listed in the protocol, other performance criteria evaluated during the verification process include: cross-reactivity, dilution linearity, kit stability, and sample behavior (for example, detectability and stability).

Each MILLIPLEX® panel and kit includes:

- Quality controls (QCs) provided to qualify assay performance
- Comparison of standard (calibrator) and QC lots to a reference lot to ensure lot-to-lot consistency
- Optimized serum matrix to mimic native analyte environment
- Detection antibody cocktails designed to yield consistent analyte profiles within panel

In addition, each panel and kit meets stringent manufacturing criteria to ensure batch-to-batch reproducibility. The MILLIPLEX® Non-Human Primate Pituitary Magnetic Bead Panel 1 thus enables you to focus on the therapeutic potential of pituitary hormones. 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® Non-Human Primate Pituitary Magnetic Bead Panel 1 is part of the most versatile system available for metabolism and endocrinology 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® Non-Human Primate Pituitary 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 non-human primate serum, plasma, and tissue culture samples: ACTH, AGRP, CNTF, FSH, GH, LH, and TSH.

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 Supplied	Volume	Quantity	Cat. No.
Pituitary Standard	Lyophilized	1 vial	PIT-8046
Pituitary Quality Controls 1 and 2	Lyophilized	2 vials	PIT-6046
Set of one 96-Well black Plate with 2 sealers	-	1 set	-
Assay Buffer	30 mL	1 bottle	L-AB
Serum Matrix*	1.5 mL	1 bottle	PIT-SM
Bead Diluent	3.5 mL	1 bottle	LBD
10X Wash Buffer**	60 mL	1 bottle	L-WB
Pituitary Detection Antibodies	3.2 mL	1 bottle	PIT-1046
Streptavidin-Phycoerythrin	5.5 mL	1 bottle	L-SAPE7
Mixing Bottle	-	1 bottle	-

* Contains 0.08% Sodium azide

** Contains 0.05% Proclin

Non-Human Primate Pituitary Hormone Antibody Immobilized Magnetic Beads:

Bead/Analyte Name	Luminex® Magnetic Bead Region	Customizable 7 Analytes (20X concentration, 200 µL)	
		Available	Cat. No.
ACTH	12	✓	HACTH-MAG
AGRP	15	✓	HAGRP-MAG
CNTF	20	✓	HCNTF-MAG
FSH	26	✓	RFSH-MAG
GH	29	✓	HGH-MAG
LH	33	✓	HLH-MAG
TSH	61	✓	HTSH-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 (Lab-Line Instruments Model No. 4625 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. 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	
Pituitary Detection Antibodies	PIT-1046		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.
Pituitary Quality Controls 1 & 2	PIT-6046	 	Warning. Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Pituitary Standard	PIT-8046	 	Warning. Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Streptavidin-Phycoerythrin	L-SAPE7		Warning. Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
10X Wash Buffer	L-WB		Warning. May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.

Technical Guidelines

To obtain reliable and reproducible results, the operator should carefully read this entire manual and fully understand all aspects of each assay step before running the assay. The following notes should be reviewed and understood before the assay is set-up.

- FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.
- Do not use beyond the expiration date on the label.
- Do not mix or substitute reagents with those from other lots or sources.
- The Antibody-Immobilized Beads are light sensitive and must be protected from light at all times. Cover the assay plate containing beads with opaque plate lid or aluminum foil during all incubation steps.
- It is important to allow all reagents to warm to room temperature (20-25 °C) before use in the assay.
- Incomplete washing can adversely affect the assay outcome. All washing must be performed with the Wash Buffer provided.
- The standards prepared by serial dilution must be used within 1 hour of preparation. Discard any unused standards except the standard stock which may be stored at ≤ -20 °C for 1 month and at ≤ -80 °C for greater than one month.
- If samples fall outside the dynamic range of the assay, further dilute the samples with Assay Buffer or the appropriate diluent and repeat the assay.
- Any unused mixed Antibody-Immobilized Beads may be stored in the Bead Mix 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 using 4 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. When reading the assay on the MAGPIX®

instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 2 alignment discs.

- For the xMAP® INTELLIFLEX instrument, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.
- For cell culture supernatants or tissue extraction, use the culture or extraction medium as the matrix solution in blank, standard curve and controls
- For serum/plasma samples, use the matrix provided in the kit in the kit as the matrix solution in blank, standard curve and controls.
- 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 chunks. Centrifuge samples before use.
- Vortex all reagents well before adding to plate.

Sample Collection and Storage

Preparation of Serum Samples

- Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000 x g. Remove serum and assay immediately or aliquot and store samples at ≤ -20 °C.
- Avoid multiple (> 2) freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.

Preparation of Plasma Samples

- Plasma collection using EDTA as an anticoagulant is recommended. Centrifuge for 10 minutes at 1000 x g within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at ≤ -20 °C. If measuring ACTH, store plasma samples at -70 °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.

Preparation of Tissue Culture Supernatant

- Centrifuge the sample to remove debris and assay immediately or aliquot and store samples at ≤ -20 °C.
- Avoid multiple (> 2) freeze/thaw cycles.
- Tissue Culture Supernatant may require a dilution with an appropriate control medium prior to assay.

Note:

- 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

Sonicate each antibody-bead vial for 30 seconds; vortex for 1 minute. Add 150 μ L from each antibody bead vial to the Mixing Bottle and bring final volume to 3.0 mL with Bead Diluent. Vortex the mixed beads well. Unused portions may be stored at 2-8 $^{\circ}$ 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 7 antibody-immobilized beads, add 150 μ L from each of the 7 bead sets to the Mixing Bottle. Then add 1.95 mL Bead Diluent.

Example 2: When using 3 antibody-immobilized beads, add 150 μ L from each of the 3 bead sets to the Mixing Bottle. Then add 2.55 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 portions 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 portions at 2-8 $^{\circ}$ C for up to one month.

Preparation of Serum Matrix

This step is required for serum or plasma samples only.

Thaw Serum Matrix completely. Mix well. Dilute Serum Matrix 1:2 by adding 1.5 mL of Assay Buffer and mix well. Leftover Serum Matrix should be stored at ≤ -20 $^{\circ}$ C for up to one month.

Preparation of Pituitary Standard

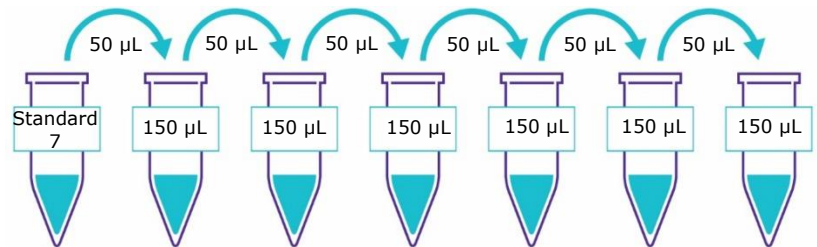
1. Prior to use, reconstitute the Pituitary Standard with 250 μL Deionized Water. Invert the vial several times to mix. Vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted standard to a polypropylene microfuge tube. This will be used as Standard 7.
2. Preparation of Working Standards
Label 6 polypropylene microfuge tubes Standard 6, Standard 5, Standard 4, Standard 3, Standard 2, and Standard 1. Add 150 μL Assay Buffer to each of the six tubes. Perform 4 times serial dilutions by adding 50 μL of the Standard 7 to the Standard 6 tube, mix well and transfer 50 μL of the Standard 6 to the Standard 5 tube, mix well and transfer 50 μL of the Standard 5 to Standard 4 tube, mix well and transfer 50 μL of the Standard 4 to the Standard 3, mix well and transfer 50 μL of the Standard 3 to the Standard 2 tube, mix well and transfer 50 μL of the Standard 2 to the Standard 1. The 0 Standard (background) will be Assay Buffer.

Reconstituted Standard	Add Deionized Water (μL)	Add Standard (volume)
-------------------------------	---	------------------------------

Standard 7	250	0
------------	-----	---

Standard	Add Assay Buffer (μL)	Add Standard (volume)
Standard 6	150	50 μL of Standard 7
Standard 5	150	50 μL of Standard 6
Standard 4	150	50 μL of Standard 5
Standard 3	150	50 μL of Standard 4
Standard 2	150	50 μL of Standard 3
Standard 1	150	50 μL of Standard 2

Preparation of Working Standards



Reconstituted Standard 7 Standard 6 Standard 5 Standard 4 Standard 3 Standard 2 Standard 1

After serial dilution, the tubes should have the following concentrations for constructing standard curves.

Standard Tube No.	AGRP (pg/mL)	FSH (mIU/mL)	GH (pg/mL)	LH (mIU/mL)	TSH (µIU/mL)	ACTH (pg/mL)	CNTF (pg/mL)
1	2.4	0.024	2.4	0.049	0.039	3	122
2	10	0.098	10	0.195	0.156	12	488
3	39	0.39	39	0.781	0.625	49	1,953
4	156	1.56	156	3.125	2.5	195	7,813
5	625	6.25	625	12.5	10	781	31,250
6	2,500	25	2,500	50	40	3,125	125,000
7	10,000	100	10,000	200	160	12,500	500,000

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), Std 1, Std 2, Std 3, Std 4, Std 5, Std 6, and Std 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 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 1:2 diluted Matrix Solution (when measuring serum or plasma samples) or appropriate culture media (when measuring culture samples) in Background, Standards, and Quality Control wells.
4. Add 25 μL Assay Buffer in Sample wells
5. Add 25 μL of Assay Buffer to the Background wells. Add 25 μL of each Standard or Control into the appropriate wells.
6. Add 25 μL of samples to the Sample wells.
7. Vortex Bead Bottle and add 25 μL of the prepared Beads to each well.
(Note: during addition of the Beads, shake beads intermittently to avoid settling.)
8. Seal the plate with a plate sealer (wrap the plate with foil if not using foil pate sealer) and incubate with agitation on a plate shaker for overnight incubation at 4 °C (16-18 hr).

Add 200 μL Assay Buffer per well



Shake 10 min, RT
Decant

- Add 25 μL Matrix Solution (or appropriate media) to Background, Standard and Quality Controls.
- Add 25 μL Assay Buffer to Sample wells.
- Add 25 μL Assay Buffer to background wells and 25 μL Standard, and 25 μL Controls to Standard and Control wells, respectively.
- Add 25 μL samples to Sample Wells.
- Add 25 μL Beads to each well



Incubate overnight at 4 °C with agitation on a plate shaker

9. Gently remove well contents and wash plate 2 times following instructions listed in the Plate Washing section.
10. Add 25 μL of Detection Antibodies into each well.
(**Note:** allow the Detection Antibodies to warm to room temperature prior to addition.)
11. Seal, cover with foil, and incubate with agitation on a plate shaker for 60 minutes at room temperature (20-25 $^{\circ}\text{C}$) **DO NOT ASPIRATE AFTER INCUBATION.**
12. Add 25 μL Streptavidin-Phycoerythrin to each well containing the 25 μL of Detection Antibodies.
13. Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25 $^{\circ}\text{C}$).
14. Gently remove well contents and wash plate 2 times following instructions listed in the Plate Washing section.
15. Add 100 μL of Sheath Fluid PLUS (or Drive Fluid PLUS if using MAGPIX[®]) to all wells. Resuspend the beads on a plate shaker for 5 minutes.
16. Run plate on Luminex[®] 200[™], HTS, FLEXMAP 3D[®], MAGPIX[®] instrument with xPONENT[®] software or xMAP[®] INTELLIFLEX instrument with INTELLIFLEX software.
17. Save and analyze the median Fluorescent Intensity (MFI) data using a weighted 5-parameter logistic or spline curve-fitting method for calculating analyte concentrations in samples.



Remove well contents.
Wash 2X with 200 μL
Wash Buffer

Add 25 μL Detection Antibody
per well



Incubate 60 min at RT
Do Not Aspirate

Add 25 μL
Streptavidin-Phycoerythrin
per well



Incubate for 30
minutes at RT

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

Add 100 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	7-Plex Beads

ACTH Beads	12
AGRP Beads	15
CNTF Bead	20
FSH Bead	26
GH Bead	29
LH Bead	33
TSH Bead	61

Quality Controls

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert or can be located at our website SigmaAldrich.com using the catalogue number as the keyword.

Assay Characteristics

Assay Sensitivities (minimum detectable concentrations)

Minimum Detectable Concentration (MinDC) defines the reportable range of the assay. It is a measure of the true limits of detection for an assay and is mathematically determined.

Overnight Protocol (N = 6 assays)		
Analyte	Mean MinDC	Mean MinDc + 2SD
ACTH (pg/mL)	0.79	1.42
AGRP (pg/mL)	0.80	1.81
CNTF (pg/mL)	37.42	100.94
FSH (mIU/mL)	0.012	0.02
GH (pg/mL)	2.03	4.03
LH (mIU/mL)	0.013	0.02
TSH (μ IU/mL)	0.005	0.01

Precision

Intra-assay precision is generated from the mean of the %CV's from 16 reportable results across two different concentration of analytes in one experiment. Inter-assay precision is generated from the mean of the %CV's from two reportable results each for two different concentrations of cytokine different experiments.

Analyte	Intra-Assay CV %	Inter-Assay CV %
ACTH	< 10	< 15
AGRP	< 10	< 15
CNTF	< 10	< 15
FSH	< 10	< 15
GH	< 10	< 15
LH	< 10	< 15
TSH	< 10	< 15

Accuracy

Spike Recovery: The data represents mean recovery of 3 levels of spiked standards using 5 independent serum samples.

Analyte	Spike and Recovery %
ACTH	112
AGRP	93
CNTF	71
FSH	98
GH	106
LH	70
TSH	113

Cross-Reactivity

The antibody pairs in the panel are specific only to the desired analyte and exhibit no or negligible (< 2%) cross-reactivity with other analytes in the panel.

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 flush, Back flush and washes; or if needed probe should be removed and sonicated.
Insufficient Bead Count	Probe height not adjusted correctly	When reading the assay on the Luminex® 200™ instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 4 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. When reading the assay on the MAGPIX® instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 2 alignment discs. 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 wells were contaminated	Avoid cross-well contamination by using sealer appropriately, and pipetting with Multichannel pipets without touching reagent in plate.
Background is too high	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® not calibrated correctly or recently	Calibrate Luminex® based on Instrument Manufacturer's instructions, at least once a week or if temperature has changed by > 3 °C.
	Gate Settings not adjusted correctly	Some Luminex® instruments (for example, Bioplex®) require different gate settings than those described in the Kit protocol. Use Instrument default settings.
	Wrong bead regions in protocol template	Check kit protocol for correct bead regions or analyte selection.
	Incorrect sample type used	Samples containing organic solvents or if highly viscous should be diluted or dialyzed as required.
	Instrument not washed or primed	Prime the Luminex® 4 times to rid of air bubbles, wash 4 times with sheath fluid or water if there is any remnant alcohol or sanitizing liquid.
Signal for whole plate is same as background	Beads were exposed to light	Keep plate and bead mix covered with dark lid or aluminum foil during all incubation steps.
	Incorrect or no Detection Antibody was added	Add appropriate Detection Antibody and continue.
Low signal for standard curve	Streptavidin-Phycoerythrin was not added	Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.
	Detection Antibody may have been removed prior to adding Streptavidin Phycoerythrin	May need to repeat assay if desired sensitivity not achieved.
Signals too high, standard curves are saturated	Incubations done at inappropriate temperatures, timings or agitation	Assay conditions need to be checked.
	Calibration target value set too high	With some Luminex® Instrument (for example, Bio-plex®) Default target setting for RP1 calibrator is set at High PMT. Use low target value for calibration and reanalyze plate.

Problem	Probable Cause	Solution
	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 just that particular analyte.
	Standard curve was saturated at higher end of curve.	See above.
High Variation in samples and/or standards	Multichannel pipet may not be calibrated	Calibrate pipets.
	Plate washing was not uniform	Confirm all reagents are removed completely in all wash steps.
	Samples may have high particulate matter or other interfering substances	See above.
	Plate agitation was insufficient	Plate should be agitated during all incubation steps using a vertical plate shaker at a speed where beads are in constant motion without causing splashing.
	Cross well contamination	Check when reusing plate sealer that no reagent has touched sealer. Care should be taken when using same pipet tips that are used for reagent additions and that pipet tip does not touch reagent in plate.

FOR FILTER PLATES ONLY

Problem	Probable Cause	Solution
Filter plate will not vacuum	Vacuum pressure is insufficient	Increase vacuum pressure such that 0.2 mL buffer can be suctioned in 3-5 seconds.
	Samples have insoluble particles	Centrifuge samples just prior to assay setup and use supernatant.
	High lipid concentration	After centrifugation, remove lipid layer and use supernatant.
Plate leaked	Vacuum Pressure too high	Adjust vacuum pressure such that 0.2 mL buffer can be suctioned in 3-5 seconds. May need to transfer contents to a new (blocked) plate and continue.
	Plate set directly on table or absorbent towels during incubations or reagent additions	Set plate on plate holder or raised edge so bottom of filter is not touching any surface.
	Insufficient blotting of filter plate bottom causing wicking	Blot the bottom of the filter plate well with absorbent towels after each wash step.
	Pipette touching plate filter during additions	Pipette to the side of plate.
	Probe height not adjusted correctly	Adjust probe to 3 alignment discs in well H6.
	Sample too viscous	May need to dilute sample.

Product Ordering

REPLACEMENT REAGENTS	Cat. No.
Pituitary Standard	PIT-8046
Pituitary Quality Controls	PIT-6046
Pituitary Detection Antibodies	PIT-1046
Serum Matrix	PIT-SM
Bead Diluent	LBD
Assay Buffer	L-AB
Streptavidin-Phycoerythrin	L-SAPE7
Set of two 96-Well Black plates with 4 sealers	MAG-PLATE
10X Wash Buffer	L-WB

Antibody-Immobilized Magnetic Beads

Analyte	Bead No.	Cat. No.
ACTH	12	HACTH-MAG
AGRP	15	HAGRP-MAG
CNTF	20	HCNTF-MAG
FSH	26	RFSH-MAG
GH	29	HGH-MAG
LH	33	HLH-MAG
TSH	61	HTSH-MAG

Well Map

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

NHPPT1MG46K Rev 08/21

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

The logo for MilliporeSigma, featuring the word "MILLIPORE" in a blue, sans-serif font above the word "SIGMA" in a larger, bold, blue, sans-serif font.