

## User Guide

# MILLIPLEX® Non-Human Primate Cytokine Magnetic Bead Panel II

## 96-Well Plate Assay

**PRCYT2MAG40K**  
**PRCYTMAG40K-PX24**  
**PRCYTMAG40K-PX25**  
**PRCY2MG40PMX24BK**  
**PRCY2MG40PMX25BK**

Introduction .....	2	Equipment Settings .....	19
Principle .....	4	Quality Controls .....	21
Storage Conditions Upon Receipt ...	5	Assay Characteristics .....	21
Reagents Supplied .....	5	Cross-Reactivity .....	21
Materials Required .....	8	Assay Sensitivities .....	22
Safety Precautions .....	9	Precision .....	23
Symbol Definitions .....	9	Accuracy .....	24
Technical Guidelines .....	10	Troubleshooting .....	25
Sample Collection and Storage ....	11	Product Ordering .....	29
Preparation of Reagents for Immunoassay .....	13	Well Map .....	31
Immunoassay Procedure .....	17	Notice .....	32
Plate Washing .....	19	Contact Information .....	32
		Technical Assistance .....	32
		Standard Warranty .....	32
		Safety Data Sheets (SDS) .....	32

---

## Introduction

Cytokines, immunomodulatory polypeptides that play key roles in both adaptive and innate immune responses, are one of the regulatory mechanisms of the immune system. Acting at the recognition, activation or effector phases of an immune response, cytokines modulate the development and functional activities of the subtypes of T cells, B cells and myeloid cells. Since chronic inflammation and aberrant immune response contribute to many disease states, research involving cytokines and chemokines plays a significant role in achieving a deeper understanding of the etiology and progression of diseases such as infectious disease, autoimmune and chronic inflammatory disease, cardiovascular disease, metabolic syndrome, neurological disorders and cancer.

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 Cytokine Magnetic Bead Panel II thus enables you to focus on the therapeutic potential of cytokines and chemokines, as well as the modulation of their expression. 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 Cytokine Magnetic Bead Panel II is part of the most versatile system available for immunology 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 select a 24-plex (for serum/plasma) or 25-plex (for cell culture) premixed kit or
- The ability to choose any combination of analytes from our panel of 25 analytes to design a custom kit that better meets your needs.  
(**Note:** RANTES is for tissue culture samples only.)
- 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 Cytokine Magnetic Bead Panel II is a 25 plex kit to be used for the simultaneous quantification of any or all of the following analytes in monkey serum, plasma, tissue/cell lysate and culture supernatant samples: sCD137, Eotaxin, sFASL, FGF-2, Fractalkine, Granzyme A, Granzyme B, IL-1 $\alpha$ , IL-2, IL-4, IL-6, IL-16, IL-17A, IL-17E/IL-25, IL-21, IL-22, IL-23, IL-28A, IL-31, IL-33, IP-10, MIP-3 $\alpha$ , Perforin, RANTES and TNF $\beta$ .

**Note: RANTES is for tissue culture samples only and cannot be combined with other analytes in this panel when measuring serum or plasma due to different dilution needs.**

**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 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  $\leq -20$  °C. Avoid multiple ( $> 2$ ) freeze/thaw cycles.
- **DO NOT FREEZE** Antibody-Immobilized Beads, Detection Antibody, and Streptavidin-Phycoerythrin.

## Reagents Supplied

Store all reagents at 2–8 °C

Reagents	Volume	Quantity	Cat. No.
Non-Human Primate Cytokine Panel II Standard	Lyophilized	1 vial	PRCYT-8040
Non-Human Primate Cytokine Panel II Quality Controls 1 and 2	Lyophilized	2 vials	PRCYT-6040
Serum Matrix*	Lyophilized	1 bottle	MXHSM-8
Set of one 96-Well Plate with 2 sealers	-	1 set	-
Assay Buffer	30 mL	2 bottles	L-AB
Bead Diluent	3.5 mL	1 bottle	LBD
10X Wash Buffer**	60 mL	1 bottle	L-WB
Non-Human Primate Cytokine Panel II Detection Antibodies	3.2 mL	1 bottle	PRCYT-1040
Streptavidin-Phycoerythrin	3.2 mL	1 bottle	L-SAPE4
Mixing Bottle (not provided with premixed panel)	-	1 bottle	-

\* Contains 0.08% Sodium azide

\*\* Contains 0.05% Proclin

## Non-Human Primate Cytokine Panel II Antibody-Immobilized Premixed Magnetic Beads

	Volume	Quantity	Cat. No.
Premixed 24-plex Beads	3.5 mL	1 bottle	PRCY2PMX24-MAG
Premixed 25-plex Beads (Premixed 24 plex + RANTES)	3.5 mL	1 bottle + 1 vial	PRCY2PMX24-MAG + HCYRNTS-MAG

Included Non-Human Primate Cytokine Panel II Antibody-Immobilized Magnetic Beads are dependent on customizable selection of analytes within the panel.

## Non-Human Primate Cytokine Panel II Antibody-Immobilized Magnetic Beads

Bead/Analyte Name	Luminex® Magnetic Bead Region	Customizable 25 Analytes (50X concentration, 90 µL)		24-Plex Magnetic Premixed Beads
		Available	Cat. No.	
Anti-FGF-2 Bead	13	✓	HCYFGF2-MAG	✓
Anti-Eotaxin Bead	14	✓	HETXN-MAG	✓
Anti-IL-16 Bead	18	✓	PRIL16-MAG	✓
Anti-Fractalkine Bead	21	✓	HFKN-MAG	✓
Anti-sCD137 Bead	22	✓	HSCD137-MAG	✓
Anti-MIP-3α Bead	28	✓	HMIP3A-MAG	✓
Anti-Granzyme A Bead	29	✓	HGRNZMA-MAG	✓
Anti-IL-2 Bead	33	✓	PRIL2-MAG	✓
Anti-Granzyme B Bead	37	✓	HCDGRNZMB-MAG	✓
Anti-IL-17A Bead	39	✓	HIL17-MAG	✓
Anti-IL-22 Bead	43	✓	HIL22-MAG	✓
Anti-IL-1α Bead	44	✓	HIL1A-MAG	✓
Anti-IL-33 Bead	45	✓	HIL33-MAG	✓
Anti-IL-21 Bead	52	✓	HIL21-MAG	✓
Anti-IL-4 Bead	53	✓	HIL4-MAG	✓
Anti-IL-23 Bead	54	✓	HIL23-MAG	✓
Anti-IL-6 Bead	57	✓	HCYIL6-MAG	✓
Anti-IP-10 Bead	65	✓	HIP10-MAG	✓
Anti-IL-17E/IL-25 Bead	66	✓	HIL17E-MAG	✓
Anti-sFasL Bead	67	✓	HCDSFASLG-MAG	✓
Anti-IL-31 Bead	73	✓	HIL31-MAG	✓
Anti-RANTES Bead	74	✓	HCYRNTS-MAG	
Anti-TNFβ Bead	76	✓	HTNFB-MAG	✓
Anti-IL-28A Bead	77	✓	HIL28A-MAG	✓
Anti-Perforin Bead	78	✓	HPRFRN-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 No. B200 or equivalent)
- Titer Plate Shaker (VWR® Microplate Shaker Cat. No. 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, Cat. No. 40-094, 40-095, 40-096, 40-097 or equivalent) or Handheld Magnetic Separation Block (Cat. No. 40-285 or equivalent).







**Note:** If a plate washer or handheld magnetic separation block for magnetic beads is not available, one can use a microtiter filter plate (Cat. No. MX-PLATE) to run the assay using a vacuum filtration unit (Vacuum Manifold, Cat. No. MSVMHTS00 or equivalent with Vacuum Pump, Cat. No. WP6111560 or equivalent).



## Safety Precautions

- All blood components and biological materials should be handled as potentially hazardous. Follow universal precautions as established by the Centers for Disease Control and Prevention and by the Occupational Safety and Health Administration when handling and disposing of infectious agents.
- Sodium azide or Proclin has been added to some reagents as a preservative. Although the concentrations are low, Sodium azide and Proclin may react with lead and copper plumbing to form highly explosive metal azides. Dispose of unused contents and waste in accordance with international, federal, state, and local regulations.

## Symbol Definitions

Ingredient	Cat. No.	Label	
Streptavidin-Phycoerythrin	L-SAPE4		<b>Warning.</b> 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		<b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.
Serum Matrix	MXHSM-8	No Symbol Required	Harmful to aquatic life with long lasting effects. Avoid release to the environment.
Non-Human Primate Cytokine Panel II Quality Control 1 & 2	PRCYT-6040	 	<b>Warning.</b> Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Non-Human Primate Cytokine Panel II Standard	PRCYT-8040	 	<b>Warning.</b> Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.

---

## 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  $\leq -20$  °C for 1 month and at  $\leq -80$  °C for greater than one month.
- If samples fall outside the dynamic range of the assay, further dilute the samples with the appropriate diluent and repeat the assay.
- Any unused mixed Antibody-Immobilized Beads may be stored in the Mixing Bottle at 2-8 °C for up to 1 month.
- During the preparation of the standard curve, make certain to mix the higher concentration well before making the next dilution. Use a new tip with each dilution.
- The plate should be read immediately after the assay is finished. If, however, the plate cannot be read immediately, seal the plate, cover with aluminum foil or an opaque lid, and store the plate at 2-8 °C for up to 24 hours. Prior to reading, agitate the plate on the plate shaker at room temperature for 10 minutes. Delay in reading a plate may result in decreased sensitivity for some analytes.
- The titer plate shaker should be set at a speed to provide maximum orbital mixing without splashing of liquid outside the wells. For the recommended plate shaker, this would be a setting of 5-7 which is approximately 500-800 rpm.
- Ensure that the needle probe is clean. This may be achieved by sonication and/or alcohol flushes.

- When reading the assay on the Luminex® 200™ instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on the MAGPIX® instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on the FLEXMAP 3D® instrument, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 1 alignment disc.
- For the FLEXMAP 3D® instrument, when using the solid plate in the kit, the final resuspension should be with 150 µL Sheath Fluid PLUS in each well and 75 µL should be aspirated.
- For 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 cell culture supernatants or tissue extraction, use the culture or extraction medium as the matrix solution in background, standard curve and control wells. If samples are diluted in assay buffer, use the assay buffer as matrix.
- For serum/plasma samples that require further dilution beyond 1:2, use the assay buffer provided in the kit.
- For cell/tissue homogenate, the final cell or tissue homogenate should be prepared in a buffer that has a neutral pH, contains minimal detergents or strong denaturing detergents, and has an ionic strength close to physiological concentration. Avoid debris, lipids, and cell/tissue aggregates. Centrifuge samples before use.
- Vortex all reagents well before adding to plate.

## Sample Collection and Storage

### Preparation of Serum Samples

- Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000 x g. Remove serum and assay immediately or aliquot and store samples at ≤ -20 °C.
- Avoid multiple > 2 freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.
- Serum samples should be diluted 1:2 in the Assay Buffer provided in the kit. For example, in a tube, 35 µL of serum may be combined with 35 µL of Assay Buffer. When further dilution beyond 1:2 is required, use Assay Buffer as the diluent.
- **(Note: RANTES is for tissue culture samples only and cannot be combined with other analytes in this panel when measuring serum or plasma.)**

## Preparation of Plasma Samples

- Plasma collection using EDTA as an anti-coagulant is recommended. Centrifuge for 10 minutes at 1000 x g within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at  $\leq -20^{\circ}\text{C}$ .
- 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.
- Plasma samples should be diluted 1:2 in the assay buffer provided in the kit. For example, in a tube, 35  $\mu\text{L}$  of serum may be combined with 35  $\mu\text{L}$  of Assay Buffer. When further dilution beyond 1:2 is required, use assay buffer as the diluent.
- **(Note: RANTES is for tissue culture samples only and cannot be combined with other analytes in this panel when measuring serum or plasma.)**

## 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.
- Tissue culture supernatant may require a dilution with an appropriate control medium prior to assay. Tissue/cell extracts should be done in neutral buffers containing reagents and conditions that do not interfere with assay performance. Excess concentrations of detergent, salt, denaturants, high or low pH, etc. will negatively affect the assay. Organic solvents should be avoided. The tissue/cell extract samples should be free of particles such as cells or tissue debris.

### NOTE:

- A maximum of 25  $\mu\text{L}$  per well of 1:2 diluted serum or plasma can be used. Tissue culture or other media may also be used.
- **Note: RANTES is for tissue culture samples only and cannot be combined with other analytes in this panel when measuring serum or plasma.**
- 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 Antibody-Immobilized Beads

- If premixed beads are used, sonicate the premixed bead bottle 30 seconds and then vortex for 1 minute before use.

To prepare 25-plex premixed beads, add 70  $\mu\text{L}$  of RANTES to the 24-plex premixed bead bottle. Mix well before use.

- For individual vials of beads, sonicate each antibody-bead vial for 30 seconds; vortex for 1 minute. Add 60  $\mu\text{L}$  from each antibody-bead vial to the Mixing Bottle and bring final volume to 3.0 mL with LBD. Vortex the mixed beads well. Unused portion may be stored at 2-8 °C for up to one month.

**(Note:** Due to the composition of magnetic beads, you may notice a slight color in the bead solution. This does not affect the performance of the beads or the kit.)

Example 1: When using 3 antibody-immobilized beads, add 60  $\mu\text{L}$  from each of the 3 bead vials to the Mixing Bottle. Then add 2.82 mL LBD.

Example 2: When using 20 antibody-immobilized beads, add 60  $\mu\text{L}$  from each of the 20 bead vials to the Mixing Bottle. Then add 1.8 mL LBD.

### Preparation of Quality Controls

Before use, reconstitute Quality Control 1 and Quality Control 2 with 250  $\mu\text{L}$  deionized water. Invert the vial several times to mix and vortex. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted Quality Control 1 and Quality Control 2 into two polypropylene microfuge tubes. Unused portion may be stored at  $\leq -20$  °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 the unused portion at 2-8 °C for up to one month.

### Preparation of Serum Matrix

**This step is required for serum or plasma samples only.**

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

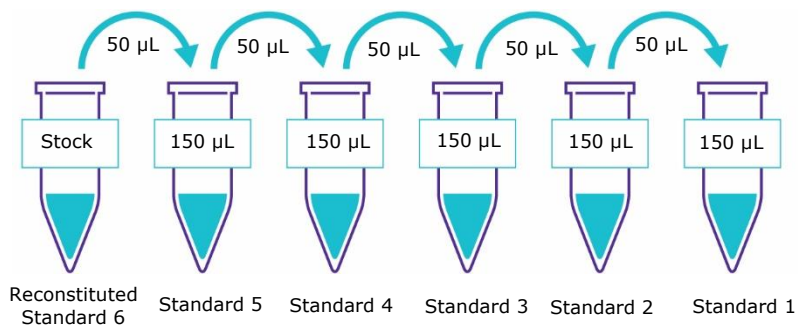
**Preparation of Non-Human Primate Cytokine Panel II Standard**

- 1. Prior to use, reconstitute the Non-Human Primate Cytokine Panel II Standard with 250 µL deionized water to give a stock concentration of standard for all analytes [see table below for analyte concentrations]. Invert the vial several times to mix. Vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted standard to a polypropylene microfuge tube. This will be used as the standard 6; the unused portion may be stored at ≤ -20 °C for up to one month.
- 2. Preparation of Working Standards  
Label five polypropylene microfuge tubes as Standard 5, Standard 4, Standard 3, Standard 2 and Standard 1. Add 150 µL of Assay Buffer to each of the five tubes. Prepare 1:4 serial dilutions by adding 50 µL of the reconstituted Standard 6 to the Standard 5 tube, mix well and transfer 50 µL of the Standard 5 to the Standard 4 tube, mix well and transfer 50 µL of the Standard 4 to the Standard 3 tube, 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 tube. The 0 Standard (Background) will be the Assay Buffer.

Standard	Add Deionized Water (µL)	Add Standard (volume)
Standard 6	250	0

Standard No.	Add Assay Buffer (µL)	Add Standard (volume)
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 Standards



Standard	FGF2, IL-33 (pg/mL)	Eotaxin (pg/mL)	IL-16 (pg/mL)	Fractalkine, IL-22, TNF $\beta$ , Perforin (pg/mL)	sCD137, IP-10, IL-2 (pg/mL)
Standard 1	20	3	73	98	10
Standard 2	78	12	293	391	39
Standard 3	313	47	1172	1563	156
Standard 4	1250	188	4688	6250	625
Standard 5	5000	750	18750	25000	2500
Standard 6	20000	3000	75000	100000	10000

<b>Standard</b>	<b>MIP3α (pg/mL)</b>	<b>Granzyme A (pg/mL)</b>	<b>IL-1α IL-28A (pg/mL)</b>	<b>Granzyme B (pg/mL)</b>	<b>IL-17A (pg/mL)</b>
Standard 1	7	78	29	4	1
Standard 2	27	313	117	16	6
Standard 3	109	1250	469	63	23
Standard 4	438	5000	1875	250	94
Standard 5	1750	20000	7500	1000	375
Standard 6	7000	80000	30000	4000	1500

<b>Standard</b>	<b>IL-21, IL-31 (pg/mL)</b>	<b>IL-4 (pg/mL)</b>	<b>IL-23 (pg/mL)</b>	<b>IL-6 (pg/mL)</b>	<b>IL-17E/IL-25 (pg/mL)</b>
Standard 1	15	146	24	3	293
Standard 2	59	586	98	12	1172
Standard 3	234	2344	391	47	4688
Standard 4	938	9375	1563	188	18750
Standard 5	3750	37500	6250	750	75000
Standard 6	15000	150000	25000	3000	300000

<b>Standard</b>	<b>sFASL (pg/mL)</b>	<b>RANTES (pg/mL)</b>
Standard 1	8	1
Standard 2	31	3
Standard 3	125	12
Standard 4	500	47
Standard 5	2000	188
Standard 6	8000	750



## 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), Standard 1 through 6], Controls 1 and 2, and Samples on Well Map Worksheet in a vertical configuration. (**Note:** Most instruments will only read the 96-well plate vertically by default.)  
It is recommended to run the assay in duplicate.
  - If using a filter plate, set the filter plate on a plate holder at all times during reagent dispensing and incubation steps so that the bottom of the plate does not touch any surface.
1. Add 200 µL of Assay Buffer into each well of the plate. Seal and mix on a plate shaker for 10 minutes at room temperature (20-25 °C).
  2. Decant Assay Buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times.
  3. Add 25 µL of each Standard or Control into the appropriate wells. Assay Buffer should be used for 0 pg/mL standard (Background).
  4. Add 25 µL of Assay Buffer to the sample wells.
  5. Add 25 µL of appropriate matrix solution to the background, standards, and control wells. When assaying serum or plasma, use the Serum Matrix provided in the kit as the matrix solution. When assaying tissue culture or other supernatant, use proper control culture medium as the matrix solution.
  6. Add 25 µL of 1:2 diluted Samples into the appropriate wells.
  7. Vortex Mixing Bottle and add 25 µL of the Mixed or Premixed Beads to each well. (**Note:** During addition of Beads, shake bead bottle intermittently to avoid settling.)
  8. Seal the plate with a plate sealer. Wrap the plate with foil and incubate with agitation on a plate shaker overnight (16–18 hours) at 4 °C or same day 2 hrs at RT.

Add 200 µL Assay Buffer per well



Shake 10 min, RT  
Decant

- Add 25 µL Standard or Control to appropriate wells
- Add 25 µL Assay Buffer to background wells and sample wells
- Add 25 µL appropriate matrix solution to background, standards, and control wells
- Add 25 µL 1:2 diluted Samples to sample wells
- Add 25 µL Beads to each well



Incubate overnight (16-18 hours) at 4 °C or same day

9. Gently remove well contents and wash plate three times following instructions listed in the Plate Washing section.
10. Add 25  $\mu$ L 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 1 hour at room temperature (20-25  $^{\circ}$ C). **DO NOT ASPIRATE AFTER INCUBATION.**
12. Add 25  $\mu$ L Streptavidin-Phycoerythrin to each well containing the 25  $\mu$ L of Detection Antibodies.
13. Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25  $^{\circ}$ C).
14. Gently remove well contents and wash plate three times following instructions listed in the Plate Washing section.
15. Add 150  $\mu$ L of Sheath Fluid PLUS (or Drive Fluid PLUS if using MAGPIX<sup>®</sup>) to all wells. Resuspend the beads on a plate shaker for 5 minutes.
16. Run plate on Luminex<sup>®</sup> 200™, HTS, FLEXMAP 3D<sup>®</sup>, MAGPIX<sup>®</sup> with xPONENT<sup>®</sup> software, or xMAP<sup>®</sup> INTELLIFLEX with INTELLIFLEX Software.
17. Save and analyze the Median Fluorescent Intensity (MFI) data using a 5-parameter logistic or spline curve-fitting method for calculating analyte concentrations in samples.  
(**Note:** For diluted samples, multiply the calculated concentration by the dilution factor.)



Remove well contents and wash 3X with 200  $\mu$ L Wash Buffer

Add 25  $\mu$ L Detection Antibodies per well



Incubate 1 hour at RT (20-25  $^{\circ}$ C) at RT

Do Not Aspirate

Add 25  $\mu$ L Streptavidin-Phycoerythrin per well



Incubate for 30 minutes at RT

Remove well contents and wash 3X with 200  $\mu$ L Wash Buffer

Add 150  $\mu$ L Sheath Fluid PLUS or Drive Fluid PLUS per well

Read on Luminex<sup>®</sup> 100  $\mu$ L, 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  $\mu$ 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  $\mu$ 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  $\mu$ 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® 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™ and HTS	xPONENT® 3.1 compatible Calibration Kit (Cat. No. LX2R-CAL-K25)	Performance Verification Kit (Cat. No. LX2R-PVER-K25)
FLEXMAP 3D®	FLEXMAP 3D® Calibrator Kit (Cat. No. F3D-CAL-K25)	FLEXMAP 3D® Performance Verification Kit (Cat. No. F3D-PVER-K25)
xMAP® INTELLIFLEX	xMAP® INTELLIFLEX Calibration Kit (Cat. No. IFX-CAL-K20)	xMAP® INTELLIFLEX Performance Verification Kit (Cat. No. IFX-PVER-K20)
MAGPIX®	MAGPIX® Calibration Kit (Cat. No. MPX-CAL-K25)	MAGPIX® Performance Verification Kit (Cat. No. MPX-PVER-K25)

**NOTE:** When setting up a Protocol using the xPONENT® software, you must select MagPlex® as the Bead Type in the Acquisition settings.

**NOTE:** These assays cannot be run on any instruments using Luminex® IS 2.3 or Luminex® 1.7 software.

The Luminex® probe height must be adjusted to the plate provided in the kit. Please use Cat. No. MAG-PLATE, if additional plates are required for this purpose.

Events	50, per bead
Sample Size	100 µL
Gate Settings	8,000 to 15,000
Reporter Gain	Default (low PMT)
Time Out	60 seconds
Bead Set	Customizable 25-Plex Beads
	FGF-2 13
	Eotaxin 14
	IL-16 18
	Fractalkine 21
	sCD137 22
	MIP-3α 28
	Granzyme A 29
	IL-2 33
	Granzyme B 37
	IL-17A 39
	IL-22 43
	IL-1α 44
	IL-33 45
	IL-21 52
	IL-4 53
	IL-23 54
	IL-6 57
	IP-10 65
	IL-17E/IL-25 66
	sFasL 67
	IL-31 73
	RANTES 74
	TNFβ 76
	IL-28A 77
	Perforin 78

## Quality Controls

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert or can be located at our website [SigmaAldrich.com](https://www.sigmaaldrich.com) using the catalogue number as the keyword.

## Assay Characteristics

### Cross-Reactivity

There was no or negligible cross-reactivity between the antibodies for an analyte and any of the other analytes in this panel.

## Assay Sensitivities (minimum detectable concentrations, pg/mL)

Minimum Detectable Concentration (MinDC) is calculated using MILLIPLEX® Analyst 5.1. It measures the true limits of detection for an assay by mathematically determining what the empirical MinDC would be if an infinite number of standard concentrations were run for the assay under the same conditions.

Analyte	Overnight Protocol (n = 9 Assays)		2 Hour Protocol (n = 4 Assays)	
	MinDC (pg/mL)	MinDC+2SD (pg/mL)	MinDC (pg/mL)	MinDC+2SD (pg/mL)
FGF-2	12.0	18.4	11.3	18.0
Eotaxin	2.5	4.5	2.5	3.1
IL-16	36.9	84.4	64.3	73.2
Fractalkine	32.6	69.2	27.1	65.1
sCD137	5.3	10.0	3.2	10.4
MIP-3α	4.3	8.5	3.6	8.5
Granzyme A	13.6	24.3	10.2	18.8
IL-2	4.4	9.2	5.5	10.6
Granzyme B	1.9	3.4	1.6	2.9
IL-17A	1.8	3.1	1.7	3.6
IL-22	92.0	100.1	76.4	142.1
IL-1α	23.7	40.2	14.3	39.7
IL-33	6.2	14.2	4.0	12.5
IL-21	5.9	11.9	6.3	14.9
IL-4	38.5	136.1	61.9	153.6
IL-23	7.5	11.4	7.1	17.9
IL-6	7.1	17.8	1.4	3.5
IP-10	5.6	10.3	5.9	9.8
IL-17E/IL-25	252.9	366.0	266.6	296.6
sFasL	5.5	11.3	5.9	12.6
IL-31	7.2	14.5	6.8	16.6
RANTES	0.7	2.0	1.7	2.2
TNFβ	43.9	106.2	39.0	92.5
IL-28A	14.0	30.8	12.8	33.8
Perforin	35.8	98.3	38.9	61.2

## Precision

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

Analyte	Overnight Protocol		2 Hour Protocol
	Intra-assay %CV	Inter-assay %CV	Intra-assay %CV
FGF-2	< 10	< 15	< 10
Eotaxin	< 10	< 15	< 10
IL-16	< 10	< 15	< 10
Fractalkine	< 10	< 15	< 10
sCD137	< 10	< 15	< 10
MIP-3 $\alpha$	< 10	< 15	< 10
Granzyme A	< 10	< 15	< 10
IL-2	< 10	< 15	< 10
Granzyme B	< 10	< 15	< 10
IL-17A	< 10	< 15	< 10
IL-22	< 10	< 15	< 10
IL-1 $\alpha$	< 10	< 15	< 10
IL-33	< 10	< 15	< 10
IL-21	< 10	< 15	< 10
IL-4	< 10	< 15	< 10
IL-23	< 10	< 15	< 10
IL-6	< 10	< 15	< 10
IP-10	< 10	< 15	< 10
IL-17E/IL-25	< 10	< 15	< 10
sFasL	< 10	< 15	< 10
IL-31	< 10	< 15	< 10
RANTES	< 10	< 15	< 10
TNF $\beta$	< 10	< 15	< 10
IL-28A	< 10	< 15	< 10
Perforin	< 10	< 15	< 10

## Accuracy

Spike Recovery: The data represent mean percent recovery of spiked standards ranging from low, medium, and high concentration in serum matrices (n=6).

Analyte	Overnight Protocol	2-Hour Protocol
	% Recovery in Serum Matrix	% Recovery in Serum Matrix
FGF-2	99	77
Eotaxin	94	87
IL-16	89	83
Fractalkine	101	87
sCD137	100	87
MIP-3 $\alpha$	104	88
Granzyme A	97	86
IL-2	93	78
Granzyme B	102	85
IL-17A	104	81
IL-22	96	78
IL-1 $\alpha$	98	82
IL-33	106	87
IL-21	92	90
IL-4	104	80
IL-23	95	95
IL-6	88	74
IP-10	99	86
IL-17E/IL-25	109	92
sFasL	86	77
IL-31	101	86
RANTES	80	80
TNF $\beta$	98	87
IL-28A	97	91
Perforin	99	86



# Troubleshooting

Problem	Probable Cause	Solution
Insufficient bead count	Plate washer aspirate height set too low	Adjust aspiration height according to manufacturers' instructions.
	Bead mix prepared inappropriately	Sonicate bead vials and vortex just prior to adding to bead mix bottle according to protocol. Agitate bead mix intermittently in reservoir while pipetting this into the plate.
	Samples cause interference due to particulate matter or viscosity	See above. Also sample probe may need to be cleaned with alcohol flushes, back flushes and washes; or if needed, probe should be removed and sonicated.
	Probe height not adjusted correctly	When reading the assay on the Luminex® 200™ instrument, adjust probe height to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on the MAGPIX® instrument, adjust probe height to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on the FLEXMAP 3D® instrument, adjust probe height to the kit solid plate using 1 alignment disc. For the FLEXMAP 3D® instrument, when using the solid plate in the kit, the final resuspension should be with 150 µL Sheath Fluid PLUS in each well and 75 µL should be aspirated. When reading the assay on the xMAP® INTELLIFLEX instrument, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.
Background is too high	Background wells were contaminated	Avoid cross-well contamination by using sealer appropriately and pipetting with multichannel pipettes without touching reagent in plate.
	Matrix used has endogenous analyte or interference	Check matrix ingredients for cross-reacting components (for example, interleukin modified tissue culture medium).
	Insufficient washes	Increase number of washes.

<b>Problem</b>	<b>Probable Cause</b>	<b>Solution</b>
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 Detection Antibody was added	Add appropriate Detection Antibody and continue.
Low signal for standard curve	Streptavidin-Phycoerythrin was not added	Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.
	Detection Antibody may have been removed prior to adding Streptavidin-Phycoerythrin	May need to repeat assay if desired sensitivity not achieved.
	Incubations done at inappropriate temperatures, timings or agitation	Assay conditions need to be checked.
Signals too high, standard curves are saturated	Calibration target value set too high	With some Luminex® instruments (for example, Bio-Plex®) default target setting for RP1 calibrator is set at high PMT. Use low target value for calibration and reanalyze plate.
	Plate incubation was too long with standard curve and samples	Use shorter incubation time.

<b>Problem</b>	<b>Probable Cause</b>	<b>Solution</b>
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 technical support for appropriate protocol modifications.
	Samples contain analyte concentrations higher than highest standard point	Samples may require dilution and reanalysis for just that particular analyte.
	Standard curve was saturated at higher end of curve	See above.
High variation in samples and/or standards	Multichannel pipette may not be calibrated	Calibrate pipettes.
	Plate washing was not uniform	Confirm all reagents are removed completely in all wash steps.
	Samples may have high particulate matter or other interfering substances	See above.
	Plate agitation was insufficient	Plate should be agitated during all incubation steps using 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

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 set-up 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.
Non-Human Primate Cytokine Panel II Standard	PRCYT-8040
Non-Human Primate Cytokine Panel II Quality Controls 1 and 2	PRCYT-6040
Serum Matrix	MXHSM-8
Non-Human Primate Cytokine Panel II Detection Antibodies	PRCYT-1040
Streptavidin-Phycoerythrin	L-SAPE4
Assay Buffer	L-AB
Set of two 96-Well plates with sealers	MAG-PLATE
10X Wash Buffer	L-WB
Bead Diluent	LBD
Non-Human Primate Cytokine Panel II 24 Plex Premixed Magnetic Bead Kit – BULK PACKAGED	PRCY2MG40PMX24BK
Non-Human Primate Cytokine Panel II 25 Plex Premixed Magnetic Bead Kit – BULK PACKAGED	PRCY2MG40PMX25BK

### Non-Human Primate Cytokine Panel II Antibody-Immobilized Premixed Magnetic Beads:

Premixed 24-plex Beads	PRCY2PMX24-MAG	3.5 mL
------------------------	----------------	--------

## Antibody-Immobilized Magnetic Beads

Analyte	Bead No.	Cat. No.
FGF-2	13	HCYFGF2-MAG
Eotaxin	14	HETXN-MAG
IL-16	18	PRIL16-MAG
Fractalkine	21	HFKN-MAG
sCD137	22	HSCD137-MAG
MIP-3 $\alpha$	28	HMIP3A-MAG
Granzyme A	29	HGRNZMA-MAG
IL-2	33	PRIL2-MAG
Granzyme B	37	HCDGRNZMB-MAG
IL-17A	39	HIL17-MAG
IL-22	43	HIL22-MAG
IL-1 $\alpha$	44	HIL1A-MAG
IL-33	45	HIL33-MAG
IL-21	52	HIL21-MAG
IL-4	53	HIL4-MAG
IL-23	54	HIL23-MAG
IL-6	57	HCYIL6-MAG
IP-10	65	HIP10-MAG
IL-17E/IL-25	66	HIL17E-MAG
sFasL	67	HCDSFASLG-MAG
IL-31	73	HIL31-MAG
RANTES	74	HCYRNTS-MAG
TNF $\beta$	76	HTNFB-MAG
IL-28A	77	HIL28A-MAG
Perforin	78	HPRFRN-MAG

## Well Map

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

---

## 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.

PRCYT2MAG-40K Rev 11/21

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