

Luminex® Performance Assay

Human High Sensitivity Cytokine Base Kit B

Catalog Number LBHS000

For the simultaneous quantitative determination of multiple human cytokine concentrations in cell culture supernates, serum, and plasma.

This package insert must be read in its entirety before using this product.
For research use only. Not for use in diagnostic procedures.

TABLE OF CONTENTS

SECTION	PAGE
INTRODUCTION	1
PRINCIPLE OF THE ASSAY	2
LIMITATIONS OF THE PROCEDURE	2
MATERIALS PROVIDED & STORAGE CONDITIONS	3
OTHER SUPPLIES REQUIRED	4
PRECAUTIONS.....	4
TECHNICAL HINTS.....	4
SAMPLE COLLECTION & STORAGE	5
SAMPLE PREPARATION.....	5
REAGENT PREPARATION	6
DILUTED MICROPARTICLE COCKTAIL PREPARATION.....	7
DILUTED BIOTIN-ANTIBODY COCKTAIL PREPARATION	7
INSTRUMENT SETTINGS	8
ASSAY PROCEDURE	9
ASSAY PROCEDURE SUMMARY	10
CALCULATION OF RESULTS	11
CALIBRATION	11
SPECIFICITY.....	12
PLATE LAYOUT	14

Manufactured and Distributed by:

USA R&D Systems, Inc.

614 McKinley Place NE, Minneapolis, MN 55413

TEL: 800 343 7475 612 379 2956

FAX: 612 656 4400

E-MAIL: info@bio-techne.com

Distributed by:

Europe | Middle East | Africa Bio-Techne Ltd.

19 Barton Lane, Abingdon Science Park

Abingdon OX14 3NB, UK

TEL: +44 (0)1235 529449

FAX: +44 (0)1235 533420

E-MAIL: info.emea@bio-techne.com

China Bio-Techne China Co., Ltd.

Unit 1901, Tower 3, Raffles City Changning Office,

1193 Changning Road, Shanghai PRC 200051

TEL: +86 (21) 52380373 (400) 821-3475

FAX: +86 (21) 52371001

E-MAIL: info.cn@bio-techne.com

INTRODUCTION

Cytokines are intercellular signaling proteins released from a wide variety of cells and tissues. They play an integral role in regulating growth and cellular proliferation as well as modulating host response to infection, injury, and inflammation. Cytokines also influence reproduction and bone remodeling. A large number of cytokines are pleiotropic and share similar functions. In addition, many cytokines influence the production of other cytokines. Analysis and quantification of cytokines in biological fluids has thus become increasingly important. Methods such as bioassay, enzyme-linked immunosorbent assay (ELISA), intracellular staining, ribonuclease protection assay (RPA) and polymerase chain reaction (PCR) have all been used for quantifying cytokines, however, each of these techniques has limitations associated with it. These techniques are not capable of measuring multiple cytokines simultaneously in a limited sample volume.

Assessing the levels of multiple cytokines may be more revealing than analyzing a single protein. Quantifying multiple cytokines on an individual level can be time consuming and expensive. When combined with separately available analyte-specific microparticle sets, this kit is an excellent tool for simultaneously assessing the levels of multiple human cytokines in a single sample.

Any combination of the following microparticle sets are suitable for use with this base kit:

Analyte	Catalog Number	Microparticle Region
GM-CSF	LHSCM215	29
IFN- γ	LHSCM285B	30
IL-1 β	LHSCM201	20
IL-2	LHSCM202	19
IL-5	LHSCM205	22
IL-6	LHSCM206	25
IL-7	LBHS207	33
IL-13	LBHS213	34
IL-15	LBHS247	35
IL-17A	LBHS317	36
IL-17F	LBHS1335	37
IL-22	LBHS5782	38
IL-23	LBHS1716	39
IL-31	LBHS2824	42
IL-33	LBHS3625	18
IL-36 β	LBHS1099	15
TNF- α	LHSCM210	12

PRINCIPLE OF THE ASSAY

Luminex® Performance Assay multiplex kits are designed for use with any Luminex analyzer including the MAGPIX®, Luminex® 100/200™, FLEXMAP 3D®, xMAP INTELLIFLEX®, or Bio-Rad® Bio-Plex®, dual laser, flow-based sorting and detection platforms.

Analyte-specific antibodies are pre-coated onto magnetic microparticles embedded with fluorophores at set ratios for each unique microparticle region. Microparticles, standards and samples are pipetted into wells and the immobilized antibodies bind the analytes of interest. After washing away any unbound substances, a biotinylated antibody cocktail specific to the analytes of interest is added to each well. Following a wash to remove any unbound biotinylated antibody, streptavidin-phycoerythrin conjugate (Streptavidin-PE), which binds to the biotinylated antibody, is added to each well. Final washes remove unbound Streptavidin-PE, the microparticles are resuspended in buffer and read using the MAGPIX®. A magnet in the analyzer captures and holds the superparamagnetic microparticles in a monolayer. Two spectrally distinct Light Emitting Diodes (LEDs) illuminate the microparticles. One LED excites the dyes inside each microparticle to identify the region and the second LED excites the PE to measure the amount of analyte bound to the microparticle. A sample from each well is imaged with a CCD camera with a set of filters to differentiate excitation levels.

Analysis with the Luminex® 100/200™, FLEXMAP 3D®, xMAP INTELLIFLEX®, or Bio-Rad® Bio-Plex® uses one laser to excite the dyes inside each microparticle to identify the microparticle region and the second laser to excite the PE to measure the amount of analyte bound to the microparticle. All excitation emitted as each microparticle passes through the flow cell is then analyzed to differentiate excitation levels using a Photomultiplier Tube (PMT) and an Avalanche Photodiode.

LIMITATIONS OF THE PROCEDURE

- FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.
- The kit should not be used beyond the expiration date on the kit label.
- Do not mix or substitute reagents with those from other lots or sources.
- If samples fall outside the dynamic range of the assay, further dilute the samples with the appropriate calibrator diluent and repeat the assay.
- Any variation in diluent, operator, pipetting technique, washing technique, incubation time or temperature, and kit age can cause variation in binding.
- Variations in sample collection, processing, and storage may cause sample value differences.
- This assay is designed to eliminate interference by other factors present in biological samples. Until these factors have been tested in the Luminex® Performance Assay, the possibility of interference cannot be excluded.
- Luminex® Performance Assays afford the user the benefit of multi-analyte analysis of biomarkers in a single complex sample. For each sample type, a single multipurpose diluent is used to optimize recovery, linearity, and reproducibility. Such a multipurpose diluent may not optimize any single analyte to the same degree that a unique diluent selected for analysis of that analyte can optimize conditions, therefore, some performance characteristics may be more variable than those for assays designed specifically for single analyte analysis.
- **Only the analytes listed on the Standard Value Card can be measured with this base kit.**

MATERIALS PROVIDED & STORAGE CONDITIONS

Store the unopened kit at 2-8 °C. Do not use past the kit expiration date.

PART	PART #	DESCRIPTION	STORAGE OF OPENED, DILUTED, OR RECONSTITUTED MATERIAL
Standard Cocktail Human HS Cytokine Panel B	894489	2 vials of recombinant human cytokines in a buffered protein base with preservatives; lyophilized.	Use a fresh standard and control for each assay. Discard after use.
Diluent RD2-5	895982	2 vials (6 mL/vial) of a buffered protein base with preservatives.	May be stored for up to 1 month at 2-8 °C.* <i>Prepare fresh 1X solutions at the time of assay. Discard after use.</i>
Calibrator Diluent RD5K	895119	21 mL of a buffered protein base with preservatives. <i>For cell culture supernate samples.</i>	May be stored for up to 1 month at 2-8 °C.*
Calibrator Diluent RD6-65	895098	21 mL of a buffered protein base with preservatives. <i>For serum/plasma samples.</i>	
Streptavidin-PE	895613	5.5 mL of a 1X streptavidin-phycoerythrin conjugate with preservatives.	
Wash Buffer Concentrate	895003	21 mL of a 25-fold concentrated solution of buffered surfactant with preservative. <i>May turn yellow over time.</i>	
Microplate	641385	1 flat-bottomed 96-well microplate used as a vessel for the assay.	
Mixing Bottles	895505	2 empty 8 mL bottles used for mixing microparticles with Diluent RD2-5.	
Plate Sealers	640445	4 adhesive foil strips.	
Standard Value Card	749810	1 card listing the standard cocktail reconstitution volume and working standard concentrations for this lot of base kit.	

*Provided this is within the expiration date of the kit.

Additional wash buffer ([#WA126](#)) and plates ([#LYX010](#)) are available for purchase.

OTHER SUPPLIES REQUIRED

- **Luminex® Performance Assay analyte-specific kit(s) (see page 1)**
- MAGPIX®, Luminex® 100/200™, FLEXMAP 3D®, xMAP INTELLIFLEX®, or Bio-Rad® Bio-Plex® analyzer with X-Y platform
- Hand-held microplate magnet or platewasher with a magnetic platform
- Pipettes and pipette tips
- Deionized or distilled water
- Multi-channel pipette, manifold dispenser, or automated dispensing unit
- 500 mL graduated cylinder
- Horizontal orbital microplate shaker (0.12" orbit) capable of maintaining a speed of 800 ± 50 rpm
- Microcentrifuge
- **Polypropylene** test tubes for dilution of standards and samples
- Luminex® Performance Assay Control (optional; [R&D Systems®](#), # QC19)

PRECAUTIONS

Some components in this kit contain a preservative which may cause an allergic skin reaction. Avoid breathing mist.

Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.

TECHNICAL HINTS

- When mixing or reconstituting protein solutions, always avoid foaming.
- To avoid cross-contamination, change pipette tips between additions of each standard level, between sample additions, and between reagent additions. Also, use separate reservoirs for each reagent.
- To ensure accurate results, proper adhesion of plate sealers during incubation steps is necessary.
- Protect microparticles and Streptavidin-PE from light at all times to prevent photobleaching.

SAMPLE COLLECTION & STORAGE

The sample collection and storage conditions listed below are intended as general guidelines. Sample stability has not been evaluated.

Cell Culture Supernates - Remove particulates by centrifugation and assay immediately or aliquot and store samples at ≤ -20 °C. Avoid repeated freeze-thaw cycles.

Serum - Use a serum separator tube (SST) and allow samples to clot for 30 minutes at room temperature before centrifuging for 15 minutes at 1000 x g. Remove serum and assay immediately or aliquot and store samples at ≤ -20 °C. Avoid repeated freeze-thaw cycles.

Plasma - Collect plasma using EDTA or heparin as an anticoagulant. Centrifuge for 15 minutes at 1000 x g within 30 minutes of collection. Assay immediately or aliquot and store samples at ≤ -20 °C. Avoid repeated freeze-thaw cycles.

Note: *Citrate plasma has not been validated for use in this assay.*

Hemolyzed, icteric, and lipemic sample are not suitable for use in this assay.

SAMPLE PREPARATION

Use polypropylene tubes.

Note: *On the day of the assay, ALL fresh and previously frozen serum and plasma samples require centrifugation at 16,000 x g for 4 minutes immediately prior to use or dilution.*

Note: *Human IL-17A cannot be multiplexed with human IL-17F (R&D Systems® Catalog #'s LBHS317 and LBHS1335, respectively).*

Note: *Serum and plasma samples are not recommended for measurement of human IL-22 (R&D Systems, Catalog # LBHS5782).*

Note: *Cell culture supernate samples are not suitable for use in the IL-23 (R&D Systems, Catalog # LBHS1716) and IL-31 (R&D Systems, Catalog # LBHS2824) assays.*

Serum and plasma samples require a 2-fold dilution. A suggested 2-fold dilution is 150 μ L of sample + 150 μ L of Calibrator Diluent RD6-65. Mix thoroughly.

REAGENT PREPARATION

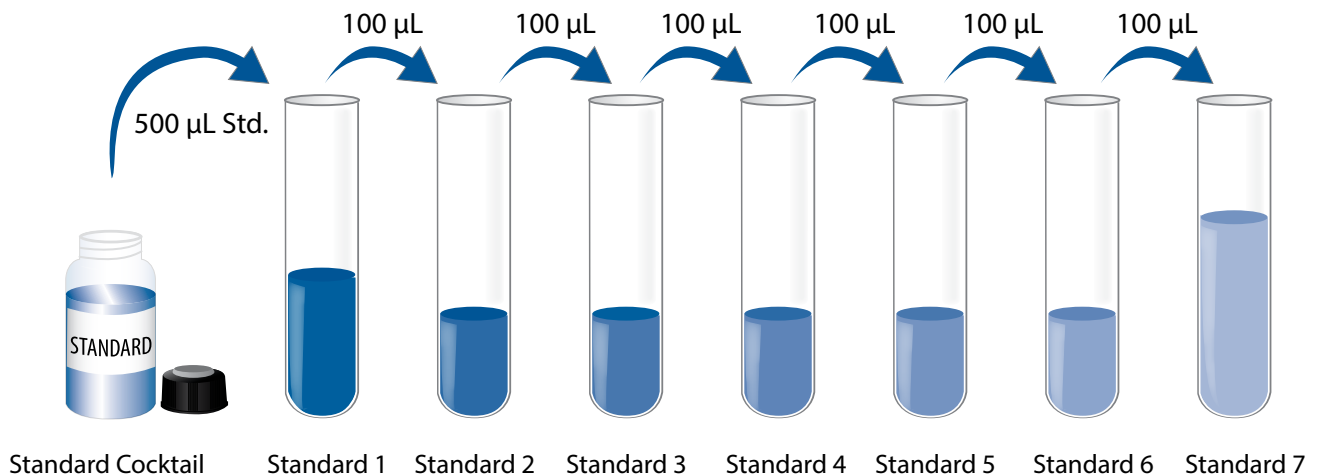
Bring all reagents to room temperature before use.

Wash Buffer - If crystals have formed in the concentrate, warm to room temperature and mix gently until the crystals have completely dissolved. Add 20 mL of Wash Buffer Concentrate into 480 mL of deionized or distilled water to prepare 500 mL of Wash Buffer.

Standard - Refer to the **Standard Value Card for the reconstitution volume and assigned values**. Reconstitute the Standard Cocktail with Calibrator Diluent RD5K (*for cell culture supernate samples*) or Calibrator Diluent RD6-65 (*for serum/plasma samples*). Allow the standard to sit for a minimum of 15 minutes with gentle agitation prior to making dilutions.

Note: For cell culture supernate samples, only calculate sample values between standard points two through seven.

Use polypropylene tubes. Pipette 500 μL of the reconstituted standard into the Standard 1 tube. Pipette 300 μL of the appropriate calibrator diluent into the remaining tubes. Use Standard 1 to produce a 4-fold dilution series (below). Refer to *analyte specific datasheet for details*. Mix each tube thoroughly before the next transfer. Standard 1 serves as the high standard. The appropriate calibrator diluent serves as the blank.



DILUTED MICROPARTICLE COCKTAIL PREPARATION

1. Centrifuge each Microparticle Concentrate vial for 30 seconds at 1000 x g prior to removing the cap.
2. Gently vortex the vials to resuspend the microparticles, taking precautions not to invert the vials.
3. Dilute the Microparticle Concentrates in the mixing bottle provided. The volume of the Microparticle Concentrate listed in the table below is for each analyte (e.g. if measuring a full plate of IL-1 β and IL-6, add 50 μ L of IL-1 β Microparticle Concentrate and 50 μ L of IL-6 Microparticle Concentrate to 2.5 mL of Diluent RD2-5).

Number of Wells Used	Microparticle Concentrate	+	Diluent RD2-5
96	50.0 μ L	+	2.50 mL
72	37.5 μ L	+	1.875 mL
48	25.0 μ L	+	1.25 mL
24	12.5 μ L	+	0.625 mL

Note: Protect microparticles from light during handling. Diluted microparticles cannot be stored. Prepare microparticles within 30 minutes of use.

DILUTED BIOTIN-ANTIBODY COCKTAIL PREPARATION

1. Centrifuge each Biotin-Antibody Concentrate vial for 30 seconds at 1000 x g prior to removing the cap.
2. Gently vortex the vials, taking precautions not to invert the vials.
3. Dilute the Biotin-Antibody Concentrates in Diluent RD2-5. The volume of the Biotin-Antibody listed in the table below is for each analyte (e.g. if measuring a full plate, add 50 μ L of each Biotin-Antibody to 5 mL of RD2-5). Mix gently.

Number of Wells Used	Biotin-Antibody Concentrate	+	RD2-5
96	50.0 μ L	+	5.00 mL
72	37.5 μ L	+	3.75 mL
48	25.0 μ L	+	2.50 mL
24	12.5 μ L	+	1.25 mL

INSTRUMENT SETTINGS

Note: Adjust the probe height setting on the analyzer to avoid puncturing the plate. Calibrate the analyzer using the proper reagents for superparamagnetic microparticles (refer to instrument manual).

MAGPIX® analyzer:

- a) Sample volume: 50 µL
- b) Assign the microparticle region for each analyte being measured (see page 1)
- c) 50 count/region
- d) Collect Median Fluorescence Intensity (MFI)

Luminex® 100/200™, FLEXMAP 3D®, xMAP INTELLIFLEX®, and Bio-Rad® Bio-Plex® analyzers:

Note: Ensure that the instrument flow rate is set to the default of 60 µL/minute (fast) for all flow based analyzers.

- a) Sample volume: 50 µL
- b) Bead Type:
 - i. Luminex® 100/200™, FLEXMAP 3D®, and xMAP INTELLIFLEX® select MagPlex
 - ii. Bio-Rad® Bio-Plex® Manager use Bio-Plex® MagPlex Beads (Magnetic)
- c) Doublet Discriminator gates:
 - i. Luminex® 100/200™ and FLEXMAP 3D®, set at 8000 and 16,500
 - ii. xMAP INTELLIFLEX® set at 7000 and 17,000
 - iii. Bio-Rad® Bio-Plex® Manager set at 8000 and 23,000
- d) Reporter Gain Setting:
 - i. Luminex® 100/200™ use Default setting
 - ii. FLEXMAP 3D® use Standard PMT setting
 - iii. xMAP INTELLIFLEX® use Luminex® 200™ Operating Mode on Low PMT setting
 - iv. Bio-Rad® Bio-Plex® Manager use the low RP1 target value for the CAL2 setting
- e) Assign the microparticle region for each analyte being measured (see page 1)
- f) 50 count/region
- g) Collect MFI

ASSAY PROCEDURE

Bring all reagents and samples to room temperature before use. It is recommended that all standards, controls, and samples be assayed in duplicate.

Note: *Protect microparticles and Streptavidin-PE from light at all times.*

1. Prepare all reagents, working standards, and samples as directed in the previous sections.
2. Add 100 μL of standard, control, or sample* per well. A plate layout is provided to record standards and samples assayed.
3. Resuspend the diluted Microparticle Cocktail by inversion or vortexing. Add 25 μL of the microparticle cocktail to each well of the microplate. Securely cover with a foil plate sealer. Incubate for 3 hours at room temperature on a horizontal orbital microplate shaker (0.12" orbit) set at 800 ± 50 rpm.
4. Using a magnetic device designed to accommodate a microplate, wash by applying the magnet to the bottom of the microplate, allow 1 minute before removing the liquid, filling each well with Wash Buffer (100 μL) and allow 1 minute before removing the liquid again. Uniform removal of liquid is essential for good performance. **Note: Do NOT blot; this may cause a loss of microparticles.** Perform the wash procedure three times.
5. Add 50 μL of diluted Biotin-Antibody Cocktail to each well. Securely cover with a foil plate sealer and incubate for 1 hour at room temperature on the shaker set at 800 ± 50 rpm.
6. Repeat the wash as in step 4.
7. Add 50 μL of Streptavidin-PE to each well. Securely cover with a foil plate sealer and incubate for 30 minutes at room temperature on the shaker set at 800 ± 50 rpm.
8. Repeat the wash as in step 4.
9. Resuspend the microparticles by adding 100 μL of Wash Buffer to each well. Incubate for 2 minutes at room temperature on the shaker set at 800 ± 50 rpm
10. Read within 90 minutes using a Luminex[®] or Bio-Rad[®] Analyzer.
Note: *Resuspend microparticles immediately prior to reading by shaking the plate for 2 minutes on the plate shaker at 800 ± 50 rpm.*

*Samples may require dilution. See Sample Preparation section.

ASSAY PROCEDURE SUMMARY

Note: Protect microparticles and Streptavidin-PE from light at all times.

- ① Prepare all reagents as instructed.
- ↓
- ② Add 100 μ L of standard, control, or sample* to each well.
- ↓
- ③ Add 25 μ L of diluted Microparticle Cocktail to each well.
Incubate for 3 hours at RT on a shaker at 800 rpm.
- ↓
- ④ Wash by removing the liquid from each well, filling with 100 μ L Wash Buffer, and removing the liquid again.
Perform the wash 3 times.
- ↓
- ⑤ Add 50 μ L of diluted Biotin-Antibody Cocktail to each well.
Cover and incubate for 1 hour at RT on the shaker at 800 rpm.
- ↓
- ⑥ Repeat the wash as in step 4.
- ↓
- ⑦ Add 50 μ L of Streptavidin-PE to each well.
Incubate for 30 minutes at RT on the shaker at 800 rpm.
- ↓
- ⑧ Repeat the wash as in step 4.
- ↓
- ⑨ Add 100 μ L of Wash Buffer to each well.
Incubate for 2 minutes at RT on the shaker at 800 rpm.
- ↓
- ⑩ Read within 90 minutes using a Luminex® or Bio-Rad® analyzer
Note: Resuspend microparticles immediately prior to reading.

*Samples require dilution. See Sample Preparation section.

CALCULATION OF RESULTS

Use the standard concentrations on the Standard Value Card and calculate 4-fold dilutions for the remaining levels. Average the duplicate readings for each standard and sample and subtract the average blank Median Fluorescence Intensity (MFI).

Create a standard curve for each analyte by reducing the data using computer software, such as [Bio-Techne® Quantist™](#), capable of generating a five parameter logistic (5-PL) curve-fit.

If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

Note: *For cell culture supernate samples only calculate sample values between standard points two through seven (see page 6).*

CALIBRATION

This assay is calibrated against highly purified recombinant human cytokines produced at R&D Systems®.

SPECIFICITY

The assay was tested for cross-reactivity and interference with the following factors. Less than 0.5% cross-reactivity and interference was observed unless otherwise noted on the analyte specific datasheet.

Note: Human IL-17A cannot be multiplexed with human IL-17F (R&D Systems® Catalog #'s LBHS317 and LBHS1335, respectively).

Recombinant human:

4-1BB Ligand	IL-2 R α	LIF R α
Amphiregulin	IL-3	LIGHT
Angiogenin	IL-3 R α	M-CSF
Angiopoietin-2	IL-4	MCP-1
APRIL	IL-4 R α	Midkine
BAFF/BLyS	IL-5 R α	MIF
BDNF	IL-6 R α	MIP-1 α
Cardiotropin-1	IL-7 R α	MIP-1 β
CD4	IL-8	β -NGF
CD40	IL-9	NT-3
CD40 Ligand	IL-10	NT-4
CNTF	IL-11	Oncostatin M (OSM)
EDA-A2	IL-12 p70	Osteoprotegerin
EDA	IL-12 p40	OX40 Ligand
β -ECGF	IL-17B	PD-ECGF
EGF	IL-17C	PDGF-AA
Epo	IL-17D	PDGF-AB
Fas	IL-17E	PDGF-BB
Fas Ligand	IL-18	Pleiotrophin/PTN
FGF acidic	IL-19	RANTES
FGF basic	IL-20	SCF
FGF-4	IL-21	SLPI
FGF-5	IL-22 R α 1	SMDF
FGF-6	IL-23 R	Lymphotoxin α 1/ β 2
G-CSF	IL-24	Lymphotoxin α 2/ β 1
GDNF	IL-26	TGF- α
GITR	IL-27	TGF- β 1
GITR Ligand	IL-28A/IFN- λ 2	TGF- β 2
GRO- α	IL-29/IFN- λ 1	TGF- β 3
GRO- β	IL-32 α	TGF- β 5
GRO- γ	IL-32 β	TGF- β RII
HB-EGF	IL-32 γ	TNF- β
HGF	IL-33 (Pro)	TNF RI
IGF-I	IL-36 α	TNF RII
IL-1 α	IL-36 β	TRAIL
IL-37/IL1F7	IL-36 γ /IL-1F9	TRANCE
IL-1ra	IL-36Ra/IL-1F5	TWEAK
IL-1 RI	Leptin	VEGF
IL-1 RII	LIF	VEGI

SPECIFICITY *CONTINUED*

Recombinant mouse:

GM-CSF
IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-7
IL-10
IL-13
IL-15
IL-17A
IL-17F
IL-22
IL-31
IL-33
IL-32 β
TNF- α

Recombinant rat:

GM-CSF
IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-10
IL-13
IL-17F
IL-22
IL-23
TNF- α

Recombinant porcine:

GM-CSF
IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-10
TNF- α

Recombinant cotton rat:

IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-10
IL-13
TNF- α

Recombinant feline:

GM-CSF
IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-10
TNF- α

Recombinant canine:

GM-CSF
IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-10
IL-13
IL-17
TNF- α

Other recombinants:

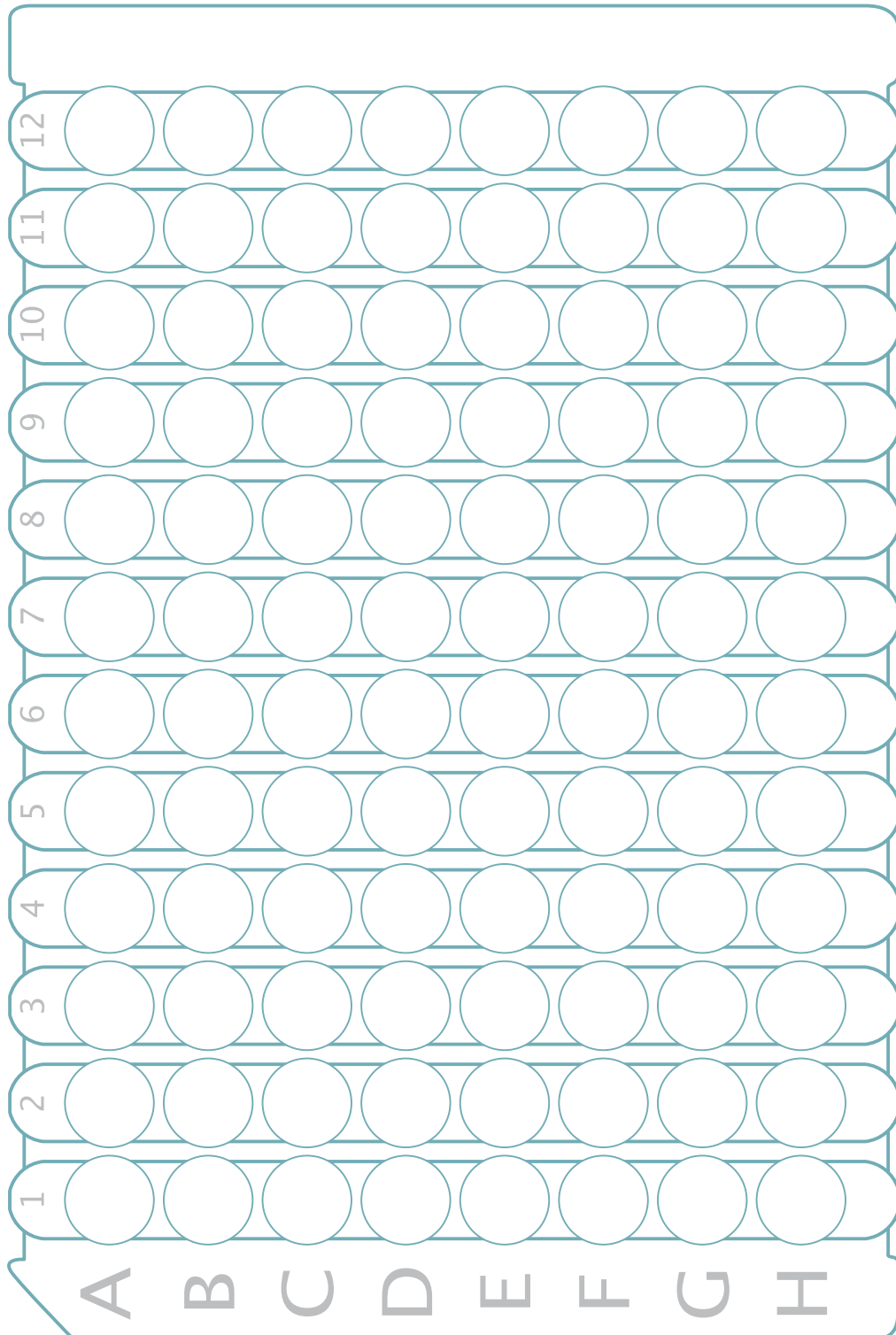
rhesus macaque IL-1 β
rhesus macaque IL-5
rhesus macaque IFN- γ
rhesus macaque TNF- α
rabbit IL-2
rabbit IL-6
rabbit TNF- α
guinea pig IL-1 β
guinea pig IL-10
bovine IFN- γ
bovine IL-2
bovine IL-5
bovine TNF- α

Recombinant human multiplex partners:

GM-CSF
IFN- γ
IL-1 β
IL-2
IL-5
IL-6
IL-7
IL-13
IL-15
IL-17A
IL-17F
IL-22
IL-23
IL-31
IL-33
IL-36 β
TNF- α

PLATE LAYOUT

Use this plate layout to record standards and samples assayed.



All trademarks and registered trademarks are the property of their respective owners.

Luminex®, MAGPIX®, Luminex® 100/200™, FLEXMAP 3D®, and xMAP INTELLIFLEX® are trademarks of Luminex Corporation.

©2025 R&D Systems®, Inc.