

Quantikine[®] QuickKit[™] ELISA

Human IL-1 β /IL-1F2 Immunoassay

Catalog Number QK201

For the quantitative determination of human Interleukin 1 beta (IL-1 β) 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
TECHNICAL HINTS.....	2
MATERIALS PROVIDED & STORAGE CONDITIONS	3
OTHER SUPPLIES REQUIRED	3
PRECAUTIONS.....	4
SAMPLE COLLECTION & STORAGE.....	4
SAMPLE PREPARATION.....	4
REAGENT PREPARATION	5
ASSAY PROCEDURE	6
CALCULATION OF RESULTS.....	7
TYPICAL DATA.....	7
PRECISION	8
RECOVERY.....	8
SENSITIVITY	8
LINEARITY.....	9
CALIBRATION	9
SAMPLE VALUES.....	9
SPECIFICITY.....	10
REFERENCES	11
PLATE LAYOUT	12

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

The Interleukin 1 (IL-1) family of proteins consists of the classic members IL-1 α , IL-1 β , and IL-1ra, plus IL-18, IL-33 and IL-1F5-F10. IL-1 α and IL-1 β bind to the same cell surface receptors and share biological functions (1). IL-1 is not produced by unstimulated cells of healthy individuals with the exception of skin keratinocytes, some epithelial cells, and certain cells of the central nervous system. However, in response to inflammatory agents, infections, or microbial endotoxins, a dramatic increase in the production of IL-1 by macrophages and various other cell types is observed. IL-1 β plays a central role in immune and inflammatory responses, bone remodeling, fever, carbohydrate metabolism, and GH/IGF-I physiology. Inappropriate or prolonged production of IL-1 has been implicated in a variety of pathological conditions including sepsis, rheumatoid arthritis, inflammatory bowel disease, acute and chronic myelogenous leukemia, insulin-dependent diabetes mellitus, atherosclerosis, neuronal injury, and aging-related diseases (2-5).

IL-1 α and IL-1 β are structurally related polypeptides that show approximately 25% homology at the amino acid level. Both are synthesized as 31 kDa precursors that are subsequently cleaved into mature proteins of approximately 17.5 kDa (6, 7). Cleavage of the IL-1 β precursor by Caspase-1/ICE is a key step in the inflammatory response (2, 8). Neither IL-1 α nor IL-1 β contains a typical hydrophobic signal peptide (9-11), but evidence suggests that these factors can be secreted by non-classical pathways (12, 13). A portion of unprocessed IL-1 α can be presented on the cell membrane and may retain biological activity (14). The precursor form of IL-1 β , unlike the IL-1 α precursor, shows little or no biological activity in comparison to the processed form (13, 15). Both unprocessed and mature forms of IL-1 β are exported from the cell.

IL-1 α and IL-1 β exert their effects through immunoglobulin superfamily receptors that additionally bind IL-1ra. The 80 kDa transmembrane type I receptor (IL-1 RI) is expressed on T cells, fibroblasts, keratinocytes, endothelial cells, synovial lining cells, chondrocytes, and hepatocytes (16, 17). The 68 kDa transmembrane type II receptor (IL-1 RII) is expressed on B cells, neutrophils, and bone marrow cells (18). The two IL-1 receptor types show approximately 28% homology in their extracellular domains but differ significantly in that the type II receptor has a cytoplasmic domain of only 29 amino acids (aa), whereas the type I receptor has a 213 aa cytoplasmic domain. IL-1 RII does not appear to signal in response to IL-1 and may function as a decoy receptor that attenuates IL-1 function (19). The IL-1 receptor accessory protein (IL-1 RAcP) associates with IL-1 RI and is required for IL-1 RI signal transduction (20). IL-1ra is a secreted molecule that functions as a competitive inhibitor of IL-1 (21, 22). Soluble forms of both IL-1 RI and IL-1 RII have been detected in human plasma, synovial fluids, and the conditioned media of several human cell lines (23, 24). In addition, IL-1 binding proteins that resemble soluble IL-1 RII are encoded by vaccinia and cowpox viruses (25).

The Quantikine® QuicKit™ Human IL-1 β /IL-1F2 Immunoassay is a one step, 80-minute solid phase ELISA designed to measure human IL-1 β in cell culture supernates, serum, and plasma. It contains *E. coli*-expressed recombinant human IL-1 β and antibodies raised against the recombinant protein. Results obtained for naturally occurring human IL-1 β showed linear curves that were parallel to the standard curves obtained using the recombinant QuicKit™ standards. These results indicate that this kit can be used to determine relative mass values for natural human IL-1 β .

PRINCIPLE OF THE ASSAY

This assay employs the quantitative sandwich enzyme immunoassay technique. An anti-tag antibody has been pre-coated onto a microplate. Standards and samples are pipetted into the wells followed by an antibody cocktail. The antibody cocktail consists of an affinity tag labeled monoclonal capture antibody and an enzyme-linked polyclonal detection antibody, specific for human IL-1 β . After washing away any unbound substances, a substrate solution is added to the wells and color develops in proportion to the amount of IL-1 β bound. The color development is stopped and the intensity of the color is measured.

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 generate values higher than the highest standard, further dilute the samples with the 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 all factors have been tested in the Quantikine[®] QuickKit[™] Immunoassay, the possibility of interference cannot be excluded.

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.
- Ensure reagent addition to plate wells is uninterrupted.
- To ensure accurate results, proper adhesion of the plate sealer during the incubation step is necessary.
- When using an automated plate washer, adding a 30 second soak period following the addition of Wash Buffer, and/or rotating the plate 180 degrees between wash steps may improve assay precision.
- Substrate Solution should remain colorless until added to the plate. Keep Substrate Solution protected from light. Substrate Solution should change from colorless to gradations of blue.
- Stop Solution should be added to the plate in the same order as the Substrate Solution. The color developed in the wells will turn from blue to yellow upon addition of the Stop Solution. Wells that are green in color indicate that the Stop Solution has not mixed thoroughly with the Substrate Solution.

MATERIALS PROVIDED & STORAGE CONDITIONS

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

PART	PART #	DESCRIPTION	STORAGE OF OPENED/ RECONSTITUTED MATERIAL
QuickKit™ Coated Microplate	899063	96 well polystyrene microplate (12 strips of 8 wells) coated with an anti-tag monoclonal antibody.	Return unused wells to the foil pouch containing the desiccant pack. Reseal along entire edge of the zip-seal. May be stored for up to 1 month at 2-8 °C.*
Human IL-1β Standard	899156	2 vials of recombinant human IL-1β in a buffered protein solution with preservatives; lyophilized. <i>Refer to the vial label for reconstitution volume.</i>	Use a new standard for each assay. Discard after use.
Human IL-1β Capture Ab Concentrate	899154	Lyophilized tagged monoclonal antibody specific for human IL-1β. <i>Refer to the vial label for reconstitution volume.</i>	May be stored for up to 1 month at 2-8 °C.*
Human IL-1β Detection Ab Concentrate	899155	400 μL of a polyclonal antibody specific for human IL-1β conjugated to horseradish peroxidase with preservatives.	
Assay Diluent RD1-132	896275	11 mL of a buffered protein base with preservatives.	
Calibrator Diluent RD5-17	895512	21 mL of a buffered protein base 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>	
Color Reagent A	895000	12 mL of stabilized hydrogen peroxide.	
Color Reagent B	895001	12 mL of stabilized chromogen (tetramethylbenzidine).	
Stop Solution	895032	6 mL of 2 N sulfuric acid.	
Plate Sealers	N/A	4 adhesive strips.	

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

OTHER SUPPLIES REQUIRED

- Microplate reader capable of measuring absorbance at 450 nm, with the correction wavelength set at 540 nm or 570 nm
- Pipettes and pipette tips
- Deionized or distilled water
- 500 mL and 100 mL graduated cylinders
- Squirt bottle, manifold dispenser, or automated microplate washer
- Horizontal orbital microplate shaker (0.12" orbit) capable of maintaining a speed of 500 ± 50 rpm
- Test tubes for dilution of standards and samples
- Human IL-1β Controls (optional; R&D Systems®, Catalog # QC266)

PRECAUTIONS

IL-1 β is detectable in saliva. Take precautionary measures to prevent contamination of kit reagents while running this assay.

The Stop Solution provided with this kit is an acid solution.

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

Color Reagent B may cause skin, eye, and respiratory irritation. Avoid breathing fumes.

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

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 centrifugation 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.*

SAMPLE PREPARATION

Cell culture supernate sample may require dilution due to high endogenous levels.

REAGENT PREPARATION

Bring all reagents to room temperature before use.

Note: *IL-1 β is detectable in saliva and sweat. Take precautionary measures to prevent contamination of the kit reagents while running the assay.*

Human IL-1 β Capture Ab Concentrate - Refer to the vial label for reconstitution volume.

Reconstitute the Human IL-1 β Capture Ab Concentrate with Assay Diluent RD1-132. This reconstitution produces a 20X Capture Antibody stock. Allow the capture antibody to sit for a minimum of 5 minutes with gentle agitation prior to diluting. Once reconstituted, the 20X Capture Antibody stock can be stored for up to 4 weeks at 2-8 °C.

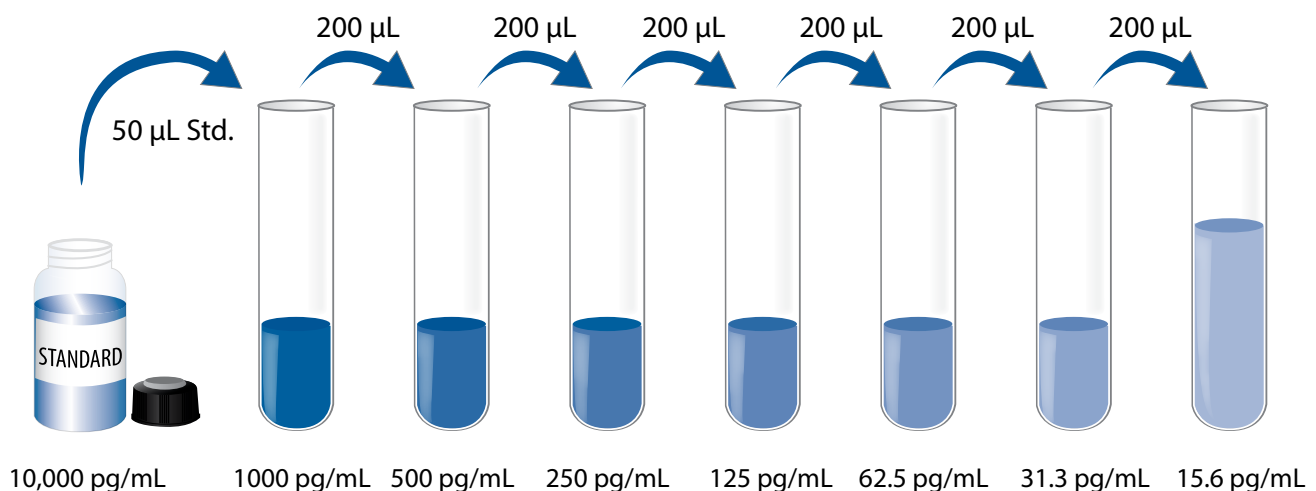
Antibody Cocktail - Dilute the reconstituted Capture Ab stock and the Detection Ab Concentrate 20-fold in Assay Diluent RD1-132. For a full plate, add 300 μ L of reconstituted Human IL-1 β Capture Ab stock and 300 μ L of Human IL-1 β Detection Ab Concentrate to 5.4 mL of Assay Diluent RD1-132 to get 6 mL of Human IL-1 β Antibody Cocktail.

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 to 480 mL of deionized or distilled water to prepare 500 mL of Wash Buffer.

Substrate Solution - Color Reagents A and B should be mixed together in equal volumes within 15 minutes of use. Protect from light. 100 μ L of the resultant mixture is required per well.

Human IL-1 β Standard - Refer to the vial label for reconstitution volume. Reconstitute the Human IL-1 β Standard with distilled or deionized water. This reconstitution produces a stock solution of 10,000 pg/mL. Allow the standard to sit for a minimum of 15 minutes with gentle agitation prior to making dilutions.

Pipette 450 μ L of Calibrator Diluent RD5-17 into the 1000 pg/mL tube. Pipette 200 μ L into each remaining tube. Use the stock solution to produce a dilution series (below). Mix each tube thoroughly before the next transfer. The 1000 pg/mL standard serves as the high standard. The Calibrator Diluent RD5-17 serves as the zero standard (0 pg/mL).



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: *IL-1 β is detectable in saliva and sweat. Take precautionary measures to prevent contamination of the kit reagents while running the assay.*

1. Prepare all reagents, samples, and working standards as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, and reseal.
3. Add 50 μ L of standard, control, or sample per well.
4. Add 50 μ L Antibody Cocktail to each well. Cover with the adhesive strip provided. Incubate for 1 hour at room temperature on a horizontal orbital microplate shaker (0.12" orbit) set at 500 \pm 50 rpm. A plate layout is provided to record standards and samples assayed.
5. Aspirate each well and wash, repeating the process twice for a total of three washes. Wash by filling each well with Wash Buffer (400 μ L) using a squirt bottle, manifold dispenser, or autowasher. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
6. Add 100 μ L of Substrate Solution to each well. Incubate for 20 minutes at room temperature **on the benchtop. Protect from light.**
7. Add 50 μ L of Stop Solution to each well. The color in the wells should change from blue to yellow. If the color in the wells is green or the color change does not appear uniform, gently tap the plate to ensure thorough mixing.
8. Determine the optical density of each well within 30 minutes, using a microplate reader set to 450 nm. If wavelength correction is available, set to 540 nm or 570 nm. If wavelength correction is not available, subtract readings at 540 nm or 570 nm from the readings at 450 nm. This subtraction will correct for optical imperfections in the plate. Readings made directly at 450 nm without correction may be higher and less accurate.

CALCULATION OF RESULTS

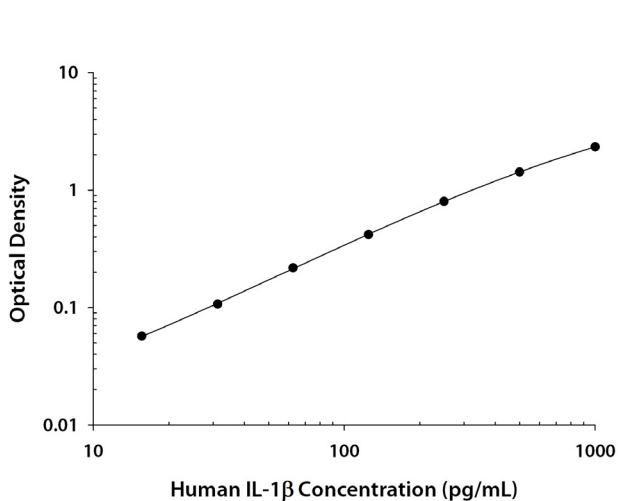
Average the duplicate readings for each standard, control, and sample and subtract the average zero standard optical density (O.D.).

Create a standard curve by reducing the data using computer software capable of generating a four parameter logistic (4-PL) curve fit. As an alternative, construct a standard curve by plotting the mean absorbance for each standard on the y-axis against the concentration on the x-axis and draw a best fit curve through the points on the graph. The data may be linearized by plotting the log of the human IL-1 β concentrations versus the log of the O.D. and the best fit line can be determined by regression analysis. This procedure will produce an adequate but less precise fit of the data.

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

TYPICAL DATA

This standard curve is provided for demonstration only. A standard curve should be generated for each set of samples assayed.



(pg/mL)	O.D.	Average	Corrected
0	0.015 0.017	0.016	—
15.6	0.072 0.073	0.073	0.057
31.3	0.122 0.123	0.123	0.107
62.5	0.233 0.234	0.234	0.218
125	0.432 0.435	0.434	0.418
250	0.812 0.823	0.818	0.802
500	1.441 1.447	1.444	1.428
1000	2.349 2.356	2.353	2.337

PRECISION

Intra-Assay Precision (Precision within an assay)

Two samples of known concentration were tested twenty times on one plate to assess intra-assay precision.

Inter-Assay Precision (Precision between assays)

Two samples of known concentration were tested in ten separate assays to assess inter-assay precision. Assays were performed by at least three technicians.

Sample	Intra-Assay Precision		Inter-Assay Precision	
	1	2	1	2
n	20	20	10	10
Mean (pg/mL)	102	526	103	528
Standard deviation	1.93	7.77	4.22	16.4
CV (%)	1.9	1.5	4.1	3.1

RECOVERY

The recovery of human IL-1 β spiked to three different levels in samples throughout the range of the assay in various matrices was evaluated.

Sample Type	Average % Recovery	Range
Cell culture media (n=4)	105	99-108%
Serum (n=2)	98	94-104%
EDTA plasma (n=2)	110	106-118%
Heparin plasma (n=2)	105	99-113%

SENSITIVITY

Fourteen assays were evaluated and the minimum detectable dose (MDD) of human IL-1 β ranged from 0.792-1.95 pg/mL. The mean MDD was 1.28 pg/mL.

The MDD was determined by adding two standard deviations to the mean O.D. value of twenty zero standard replicates and calculating the corresponding concentration.

LINEARITY

To assess the linearity of the assay, samples containing and/or spiked with high concentrations of human IL-1 β in various matrices and diluted to produce samples with values within the dynamic range of the assay. Samples were diluted prior to assay.

		Cell culture supernates* (n=6)	Serum (n=2)	EDTA plasma (n=2)	Heparin plasma (n=2)
1:2	Average % of Expected	98	98	95	94
	Range (%)	96-101	97-99	94-96	93-94
1:4	Average % of Expected	98	104	96	96
	Range (%)	94-101	102-105	95-97	95-97
1:8	Average % of Expected	97	112	97	98
	Range (%)	92-100	111-113	95-98	96-100
1:16	Average % of Expected	97	112	96	99
	Range (%)	92-94	110-115	95-97	96-103

*Samples were diluted prior to assay.

CALIBRATION

This immunoassay is calibrated against a highly purified *E. coli*-expressed recombinant human IL-1 β produced at R&D Systems®.

SAMPLE VALUES

Serum/Plasma - Ten serum and plasma samples from apparently healthy volunteers were evaluated for the presence of human IL-1 β in this assay. All samples measured less than 15.6 pg/mL. No medical histories were available for the donors used in this study.

Cell Culture Supernates - Human peripheral blood mononuclear cells (PBMCs) (1×10^6 cells/mL) were cultured in RPMI supplemented with 10% fetal bovine serum, 2 mM L-glutamine, 100 U/mL penicillin, and 100 μ g/mL streptomycin sulfate. Cells were left unstimulated or stimulated with 10 μ g/mL PHA for 24 hours. Aliquots of the culture supernates were removed, assayed for levels of human IL-1 β , and measured at 37.2 pg/mL or 2,273 pg/mL, respectively.

PBMCs (1×10^6 cells/mL) were cultured as above. Cells were left unstimulated or stimulated with 1.0 μ g/mL LPS for 24 hours. Aliquots of the cell culture supernates were removed, assayed for levels of human IL-1 β , and were not detectable or measured 25,396 pg/mL, respectively.

SPECIFICITY

This assay recognizes natural and recombinant human IL-1 β .

The factors listed below were prepared at 50 ng/mL in calibrator diluent and assayed for cross-reactivity. Preparations of the following factors at 50 ng/mL in a low level recombinant human IL-1 β control were assayed for interference. No significant cross-reactivity or interference was observed.

Recombinant human:

IL-1 Ra

IL-1 RI

IL-1 RII

ITGA-5/ITGB-1 heterodimer

Recombinant rhesus macaque IL-1 β cross-reacts approximately 177% and interferes at concentrations > 31.3 pg/mL in this assay.

Recombinant human IL-1 β precursor cross-reacts approximately 49.3% and interferes at concentrations > 62.5 pg/mL in this assay.

Recombinant rat IL-1 β cross-reacts approximately 2.0% and interferes at concentrations > 31.3 ng/mL in this assay.

Recombinant cotton rat IL-1 β cross-reacts approximately 0.56% and interferes at concentrations > 12.5 ng/mL in this assay.

Recombinant mouse IL-1 β does not interfere but does cross-react approximately 0.21% in this assay.

REFERENCES

1. Sims, J.E and D.E. Smith (2010) *Nat. Rev. Immunol.* **10**:89.
2. Martinon, F. and J. Tschopp (2007) *Cell Death Differ.* **14**:10.
3. Isoda, K. and F. Ohsuzu (2006) *J. Atheroscler. Thromb.* **13**:21.
4. Allan, S.M. *et al.* (2005) *Nat. Rev. Immunol.* **5**:629.
5. Kornman, K.S. (2006) *Am. J. Clin. Nutr.* **83**:475S.
6. Giri, J.G. *et al.* (1985) *J. Immunol.* **134**:343.
7. Hazuda, D.J. *et al.* (1988) *J. Biol. Chem.* **265**:6318.
8. Cerretti, D.P. *et al.* (1992) *Science* **256**:97.
9. Lomedico, P.T. *et al.* (1984) *Nature* **312**:458.
10. Auron, P.E. *et al.* (1987) *J. Immunol.* **138**:1447.
11. March, C.J. *et al.* (1985) *Nature* **315**:641.
12. Rubartelli, A. *et al.* (1990) *EMBO J.* **9**:1503.
13. Rubartelli, A. *et al.* (1993) *Cytokine* **5**:117.
14. Kurt-Jones, E.A. *et al.* (1985) *Proc. Natl. Acad. Sci. USA* **82**:1204.
15. Hazuda, D. *et al.* (1989) *J. Biol. Chem.* **264**:1689.
16. Urdal, D.L. *et al.* (1988) *J. Biol. Chem.* **263**:2870.
17. Sims, J.E. *et al.* (1988) *Science* **241**:585.
18. McMahan, C.J. *et al.* (1991) *EMBO J.* **10**:2821.
19. Slack, J. *et al.* (1993) *J. Biol. Chem.* **268**:2513.
20. Greenfeder, S.J. *et al.* (1995) *J. Biol. Chem.* **270**:13757.
21. Eisenberg, S.P. *et al.* (1990) *Nature* **343**:341.
22. Carter, D.B. *et al.* (1990) *Nature* **344**:633.
23. Dayer, J-M. and D. Burger (1994) *Eur. Cytokine Netw.* **5**:563.
24. Svenson, M. *et al.* (1993) *Cytokine* **5**:427.
25. Sims, J.E. and S.K. Dower (1994) *Eur. Cytokine Netw.* **5**:539.

PLATE LAYOUT

Use this plate layout to record standards and samples assayed.

The diagram shows a 12x8 microplate layout. The rows are numbered 1 through 12 on the left side, and the columns are labeled A through H at the bottom. Each well is represented by a circle. The layout is as follows:

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

NOTES

NOTES

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

©2020 R&D Systems®, Inc.