

Quantikine[®] QuickKit[™] ELISA

Human IL-2 Immunoassay

Catalog Number QK202

For the quantitative determination of human Interleukin 2 (IL-2) 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
REAGENT PREPARATION	5
ASSAY PROCEDURE	6
CALCULATION OF RESULTS	7
TYPICAL DATA	7
PRECISION	8
RECOVERY	8
SENSITIVITY	8
CALIBRATION	8
LINEARITY	9
SAMPLE VALUES	9
SPECIFICITY	10
REFERENCES	10

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

Interleukin 2 (IL-2), also known as T cell growth factor (TCGF), is a 15-18 kDa variably glycosylated α -helical polypeptide that is a member of the Common gamma Chain (γ c) cytokine family (1-4). It exists as a monomer and has a notably short half-life (< 30 minutes) (1). Human IL-2 is synthesized as a 153 amino acid (aa) precursor that contains a 20 aa signal sequence plus a 133 aa mature region (5, 6). The mature region is α -helical in nature, and contains one utilized O-linked glycosylation site at Thr3 plus three cysteines, two of which form an intrachain disulfide bond that is essential for activity (7). Mature human IL-2 shares 73%, 66%, 78%, and 97% aa identity with canine, rat, feline and rhesus monkey IL-2, respectively. Although human IL-2 shares only approximately 60% aa identity with the highly polymorphic mouse IL-2, human IL-2 is known to be active on mouse IL-2 responsive cells. Cells reported to secrete IL-2 include $\gamma\delta$ T cells (8), activated conventional CD4⁺ and CD8⁺ T cells (1, 9), neurons (10, 11), microglia (12), and hematopoietic stem cells (13).

The receptor for IL-2 (IL-2 R) is composed of three subunits, the 55 kDa CD25/IL-2 R α chain, the 70 kDa IL-2 R β chain, and the 65 kDa Common gamma Chain (1, 3). IL-2 first binds to CD25, the binary complex then recruits IL-2 R β and γ c to form the quaternary signaling complex (1, 14). In addition to IL-2, IL-2 R β is used by IL-15 in its quaternary signaling complex. γ c also serves as a signaling receptor for IL-4, -7, -9, -15, and -21 (1, 3).

In vitro studies have shown an important role for IL-2 in T cell activation and expansion. *In vivo*, IL-2 is critical for the development, maintenance and function of regulatory T cells (Treg) which provide protection against autoimmune disease. On the other hand, IL-2 can also promote autoimmune inflammation in target organs through its roles in regulating the expression of T cell trafficking genes, and production of Th2 cytokines. Within the CD8⁺ T cell subset, IL-2 is essential for optimal primary responses and differentiation into terminal effector cells. IL-2 also promotes the development of activated CD8⁺ T cells into memory cells. (1).

The Quantikine® QuickKit™ Human IL-2 Immunoassay is a one step, 80-minute solid phase ELISA designed to measure human IL-2 levels in cell culture supernates, serum, and plasma. It contains *E. coli*-expressed recombinant human IL-2 and antibodies raised against the recombinant protein. Results obtained for naturally occurring human IL-2 showed linear curves that were parallel to the standard curves obtained using the recombinant QuickKit™ standards. These results indicate that this kit can be used to determine relative mass values for natural IL-2.

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-2. After washing away any unbound substances, a substrate solution is added to the wells and color develops in proportion to the amount of IL-2 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 from other lots or sources.
- If samples generate values higher than the highest standard, dilute the samples with the 1X Kit 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® QuicKit™ 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 zip-seal. May be stored for up to 1 month at 2-8 °C.*
Human IL-2 Standard	899075	2 vials of recombinant human IL-2 in a buffered protein base with preservatives; lyophilized. <i>Refer to the vial label for the reconstitution volume.</i>	Use a new standard for each assay. Discard after use.
Human IL-2 Capture Ab Concentrate	899073	Lyophilized tagged monoclonal antibody specific for human IL-2.	May be stored for up to 1 month at 2-8 °C.*
Human IL-2 Detection Ab Concentrate	899074	400 µL of a polyclonal antibody specific for human IL-2 conjugated to horseradish peroxidase with preservatives.	
Assay Diluent RD1-76	895812	11 mL of a buffered protein base with preservatives.	
Calibrator Diluent RD5P	895151	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.
- Squirt bottle, manifold dispenser, or automated microplate washer.
- 500 mL graduated cylinder.
- Horizontal orbital microplate shaker (0.12" orbit) capable of maintaining a speed of 500 ± 50 rpm.
- Test tubes for dilution of standards.
- Human IL-2 Controls (optional; R&D Systems®, Catalog # QC261).

PRECAUTIONS

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.*

REAGENT PREPARATION

Bring all reagents to room temperature before use.

1X Kit Diluent - Prepare the 1X Kit Diluent by mixing deionized or distilled water, Assay Diluent RD1-76, and Calibrator Diluent RD5P following the volumes indicated in the table below:

	15 mL (for half plate)	30 mL (for full plate)
Deionized or distilled water	8 mL	16 mL
Assay Diluent RD1-76	5 mL	10 mL
Calibrator Diluent RD5P	2 mL	4 mL

Human IL-2 Capture Ab Concentrate - Refer to the vial label for reconstitution volume.

Reconstitute the Human IL-2 Capture Ab Concentrate with 1X Kit Diluent. 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 4 weeks at 2-8 °C.

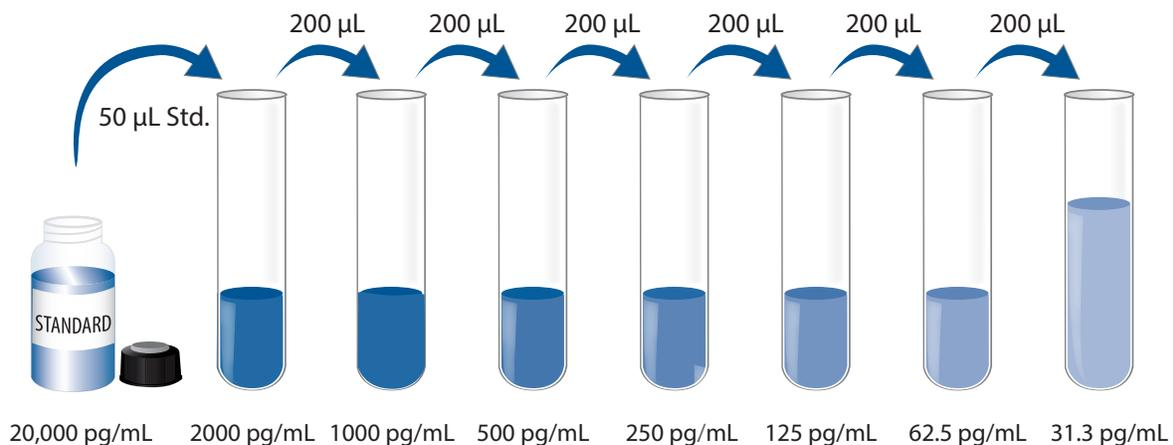
Antibody Cocktail - Dilute the reconstituted Capture Ab stock and the Detection Ab Concentrate 20-fold in 1X Kit Diluent. For a full plate, add 300 µL of reconstituted Human IL-2 Capture Ab stock and 300 µL of Human IL-2 Detection Ab Concentrate to 5.4 mL of 1X Kit Diluent.

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 µL of the resultant mixture is required per well.

Human IL-2 Standard - Refer to the vial label for reconstitution volume. Reconstitute the Human IL-2 Standard with 1X Kit Diluent. This reconstitution produces a stock solution of 20,000 pg/mL. Allow the standard to sit for a minimum of 5 minutes with gentle agitation prior to making dilutions. **Use standard stock within 45 minutes.**

Pipette 450 µL of 1X Kit Diluent into the 2000 pg/mL tube. Pipette 200 µL into the remaining tubes. Use the stock solution to produce a dilution series (below). Mix each tube thoroughly before the next transfer. The 2000 pg/mL standard serves as the high standard. The 1X Kit Diluent 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.

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 μL of standard, control, or sample per well.
4. Add 50 μL 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 ± 50 rpm.
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 μL) 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 μL of Substrate Solution to each well. Incubate for 20 minutes at room temperature **on the benchtop. Protect from light.**
7. Add 50 μL 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

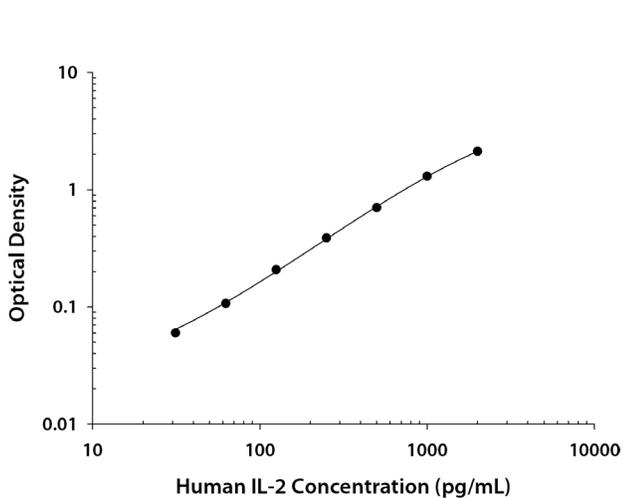
Average the duplicate readings for each standard, control, and sample and subtract the average zero standard optical density (O.D.).

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-2 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	0.007		—
	0.007	0.007	
31.3	0.058		
	0.062	0.060	0.053
62.5	0.105		
	0.108	0.107	0.100
125	0.206		
	0.209	0.208	0.201
250	0.386		
	0.389	0.388	0.381
500	0.671		
	0.734	0.703	0.696
1000	1.289		
	1.313	1.301	1.294
2000	2.083		
	2.156	2.120	2.113

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)	217	1231	213	1331
Standard deviation	5.70	33.1	17.1	67.7
CV (%)	2.6	2.7	8.0	5.1

RECOVERY

The recovery of human IL-2 spiked to three levels in samples throughout the range of the assay was evaluated.

Sample Type	Average % Recovery	Range
Cell culture media (n=4)	112	103-119%
Serum (n=2)	91	73-105%
EDTA plasma (n=2)	85	72-96%
Heparin plasma (n=2)	85	74-93%

SENSITIVITY

Fourteen assays were evaluated and the minimum detectable dose (MDD) of human IL-2 ranged from 0.684-2.01 pg/mL. The mean MDD was 1.13 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.

CALIBRATION

This immunoassay is calibrated against a highly purified *E. coli*-expressed recombinant human IL-2 produced at R&D Systems®.

LINEARITY

To assess the linearity of the assay, samples were spiked with high concentrations of human IL-2 in various matrices and diluted with 1X Kit Diluent to produce samples with values within the dynamic range of the assay.

		Cell culture media (n=4)	Serum (n=2)	EDTA plasma (n=2)	Heparin plasma (n=2)
1:2	Average % of Expected	88	93	97	94
	Range (%)	83-91	88-99	92-101	90-98
1:4	Average % of Expected	83	91	95	92
	Range (%)	77-86	88-94	92-99	89-95
1:8	Average % of Expected	80	87	92	90
	Range (%)	77-85	80-94	86-99	86-95
1:16	Average % of Expected	81	89	92	90
	Range (%)	78-85	82-97	87-98	86-94

SAMPLE VALUES

Serum/Plasma - Ten serum and plasma samples from apparently healthy volunteers were evaluated for the presence of human IL-2 in this assay. All samples measured less than the lowest standard, 31.3 pg/mL. No medical histories were available for the donors used in this study.

Cell Culture Supernates - Human peripheral blood mononuclear (PBMCs) cells were cultured in RPMI supplemented with 10% fetal bovine serum, 2 mM L-glutamine, 100 U/mL penicillin, and 100 µg/mL streptomycin sulfate. The cells were cultured unstimulated or stimulated with 10 µg/mL PHA for 24 hours. Aliquots of the cell culture supernates were removed, assayed for levels of human IL-2, and were undetectable or measured 390 pg/mL, respectively.

CD4⁺ T cells were isolated from PBMCs and cultured in RPMI supplemented with 10% fetal bovine serum, 2 mM L-glutamine, 100 U/mL penicillin, and 100 µg/mL streptomycin sulfate. Cells were stimulated with plate-bound Mouse Anti-human CD3 (Catalog # MAB100) and soluble Mouse anti-human CD28 (Catalog # MAB342) for 5 days, followed by 10 ng/mL PMA and 500 ng/mL Calcium Ionomycin for 24 hours. An aliquot of the cell culture supernate was removed, assayed for human IL-2, and measured 84,350 pg/mL.

SPECIFICITY

This assay recognizes natural and recombinant human IL-2.

The factors listed below were prepared at 50 ng/mL in 1X Kit Diluent and assayed for cross-reactivity. Preparations of the following factors at 50 ng/mL in a low level recombinant human IL-2 control were assayed for interference. No significant cross-reactivity or interference was observed.

Recombinant human:

IL-1 α
IL-1 β
IL-2R α
IL-2R β
IL-2R gamma Chain
IL-4
IL-7
IL-9
IL-15
IL-21

Other recombinants:

mouse IL-2
rat IL-2

REFERENCES

1. Malek, T.R and I. Castro (2010) *Immunity* **33**:153.
2. Waldman, T.A. (2006) *Nat. Rev. Immunol.* **6**:595.
3. Malek, T.R. (2008) *Annu. Rev. Immunol.* **26**:453.
4. Conradt, H.S. *et al.* (1989) *J. Biol. Chem.* **264**:17368.
5. Taniguchi, T. *et al.* (1983) *Nature* **302**:305.
6. Rosenberg, S.A. *et al.* (1984) *Science* **223**:1412.
7. Smith, K.A. (1984) *Annu. Rev. Immunol.* **2**:319.
8. Tsukaguchi, K. *et al.* (1995) *J. Immunol.* **154**:1786.
9. Conlon, K. *et al.* (1995) *Eur. J. Immunol.* **25**:644.
10. Hwang, I.K. *et al.* (2006) *Brain Res.* **1106**:197.
11. Giestal de Araujo, E. *et al.* (2009) *Ann. N.Y. Acad. Sci.* **1153**:57.
12. Kowalski, J. *et al.* (2004) *Pol. J. Pharmacol.* **56**:563.
13. Eguizabal, C. *et al.* (2007) *Int. J. Dev. Biol.* **51**:731.
14. Stauber, D.J. *et al.* (2006) *Proc. Natl. Acad. Sci. USA* **103**:2788.

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