

BD OptEIA™

Human IL-2 ELISA KIT II

Instruction Manual

Cat. No. 550611



BD flow cytometers are class I (1) laser products

©2005 Becton, Dickinson and Company. All rights reserved. No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from BD Biosciences.

For research use only. Not for use in diagnostic or therapeutic procedures. Purchase does not include or carry any right to resell or transfer this product either as a stand-alone product or as a component of another product. Any use of this product other than the permitted use without the express written authorization of Becton Dickinson and Company is strictly prohibited.

BD, BD Logo and all other trademarks are the property of Becton, Dickinson and Company. ©2005 BD

BD OptEIA™ ELISA Kits and Sets available from BD Biosciences – Pharmingen:

BD OptEIA ELISA Kits

| Human | Cat. No. |
|----------------------|----------|
| C3a | 550499 |
| C4a | 550947 |
| C5a | 550500 |
| C5a Kit II | 557965 |
| IFN- γ Kit II | 550612 |
| IL-1 β | 559111 |
| IL-1 β Kit II | 557966 |
| IL-2 Kit II | 550611 |
| IL-4 Kit II | 550614 |
| IL-6 Kit II | 550799 |
| IL-8 Kit II | 550999 |
| IL-10 Kit II | 550613 |
| IL-12 p40 Kit II | 551116 |
| IL-12 p70 | 559258 |
| MCP-1 | 559017 |
| TNF Kit II | 550610 |

Mouse

| | |
|---------------|--------|
| IFN- γ | 550582 |
| IL-6 | 550950 |
| TNF | 559732 |

Rat

| | |
|-----|--------|
| TNF | 550734 |
|-----|--------|

BD OptEIA ELISA Sets

(Capture antibody, Detection antibody, Streptavidin-HRP, and Standard for 5 or 20 ELISA plates).

Human

| | |
|------------------------------|---------|
| Active Caspase-3 (5 plate) | inquire |
| Cleaved PARP (5 plate) | 552592 |
| Eotaxin | 555175 |
| GM-CSF | 555126 |
| IFN- γ | 555142 |
| IL-1 β | 558848 |
| IL-1 β Set II | 557953 |
| IL-2 | 555190 |
| IL-2 sR α | 559104 |
| IL-3 | 558979 |
| IL-4 | 555194 |
| IL-5 | 555202 |
| IL-6 | 555220 |
| IL-8 | 555244 |
| IL-10 | 555157 |
| IL-12 p40 | 555171 |
| IL-12 p70 | 555183 |
| IL-15 | 559268 |
| IP-10 | 550926 |
| LT- α (TNF- β) | 550995 |
| MCP-1 | 555179 |

BD OptEIA ELISA Sets *continued*

| Human | |
|--------------|--------|
| MIG | 550998 |
| FAS | 555224 |
| sICAM-1 | 551424 |
| TNFRI | 550996 |
| TGF- β | 559119 |
| TNF | 555212 |
| TRAIL | 550948 |

Mouse

| | |
|-----------------------|--------|
| GM-CSF | 555167 |
| IFN- γ | 555138 |
| IFN- γ (AN-18) | 551866 |
| IgE | 555248 |
| IL-1 α | 550347 |
| IL-1 β | 559603 |
| IL-2 | 555148 |
| IL-3 | 555228 |
| IL-4 | 555232 |
| IL-5 | 555236 |
| IL-6 | 555240 |
| IL-10 | 555252 |
| IL-12 p40 | 555165 |
| IL-12 p70 | 555256 |
| MCP-1 | 555260 |
| TNF (Mono/Mono) | 555268 |
| TNF (Mono/Polys) | 558874 |

Rat

| | |
|---------------|--------|
| IFN- γ | 558861 |
| IL-4 | 555198 |
| IL-6 | 550319 |
| IL-10 | 555134 |
| MCP-1 | 555130 |
| TNF | 558870 |

Monkey

| | |
|---------------|--------|
| IFN- γ | 551492 |
| IL-2 | 551494 |

BD ELISA Kits

| | |
|---------------------------------|--------|
| Canine C-Reactive Protein (CRP) | 557826 |
| Rat C-Reactive Protein (CRP) | 557825 |

Table of Contents

| | |
|---|----|
| Introduction | 5 |
| Intended Use | 5 |
| Principle of the Test | 5 |
| Reagents Provided | 6 |
| Materials Required but not Provided | 6 |
| Storage Information | 7 |
| Warnings and Precautions. | 7 |
| Specimen Collection and Handling | 8 |
| Reagent Preparation | 8 |
| Assay Procedure | 9 |
| Assay Procedure Summary | 10 |
| Calculation of Results. | 10 |
| Typical Data | 11 |
| Limitations of the Procedure. | 11 |
| Performance Characteristics | 12 |
| Sensitivity | 12 |
| Recovery | 12 |
| Linearity. | 12 |
| Specificity | 13 |
| Precision | 13 |
| Intra-assay | 13 |
| Inter-assay | 13 |
| Standardization | 14 |
| Expected Values. | 14 |
| Serum/Plasma | 14 |
| Cell Culture Supernatant | 14 |
| Troubleshooting | 15 |
| References. | 15 |
| Plate Templates. | 16 |

Introduction

Interleukin-2 (IL-2), originally called T cell growth factor (TCGF), is a multi-functional cytokine that can stimulate the differentiation and proliferation of T lymphocytes and other cell types including B lymphocytes, NK cells, LAK cells, and monocytes/macrophages.¹⁻³ IL-2 exerts its biological effects by binding to specific receptors expressed by various target cells. Human IL-2 is a 15 kDa protein containing 133 amino acid residues.

The BD OptEIA™ ELISA Kit II format was developed for superior accuracy with serum and plasma specimens. The data that demonstrates this enhancement can be located in the Performance Characteristics “Recovery” and “Linearity” sections.

Intended Use

The BD OptEIA™ Human IL-2 ELISA Kit II is for the quantitative determination of human IL-2 in serum, plasma, and cell culture supernatant.

Principle of the Test

The BD OptEIA™ test is a solid phase sandwich ELISA (Enzyme-Linked Immunosorbent Assay). It utilizes a monoclonal antibody specific for IL-2 coated on a 96-well plate. Standards and samples are added to the wells, and any IL-2 present binds to the immobilized antibody. The wells are washed and streptavidin-horseradish peroxidase conjugate mixed with biotinylated anti-human IL-2 antibody is added, producing an antibody-antigen-antibody “sandwich”. The wells are again washed and TMB substrate solution is added, which produces a blue color in direct proportion to the amount of IL-2 present in the initial sample. The Stop Solution changes the color from blue to yellow, and the microwell absorbances are read at 450 nm.

Reagents Provided

| | |
|---------------------------------|--|
| Antibody Coated Wells: | 2 plates of 96 breakable wells (12 strips × 8 wells) coated with anti-human IL-2 monoclonal antibody |
| Detection Antibody: | 30 ml of biotinylated anti-human IL-2 monoclonal antibody with 0.15% ProClin™-150 as preservative |
| Standards: | 4 vials lyophilized recombinant human IL-2 |
| Enzyme Concentrate (250×): | 150 µl of 250× concentrated Streptavidin-horseradish peroxidase conjugate with BSA* and ProClin™-300 as preservative |
| Standard/Sample Diluent: | 30 ml of animal serum* with 0.09% sodium azide as preservative |
| ELISA Diluent: | 12 ml of a buffered protein base with 0.09% sodium azide as preservative |
| Wash Concentrate (20×): | 100 ml of 20× concentrated detergent solution with ProClin™-150 as preservative |
| TMB One-Step Substrate Reagent: | 30 ml of 3,3',5,5'-tetramethylbenzidine (TMB) in buffered solution |
| Stop Solution: | 13 ml of 1 M phosphoric acid |
| Plate Sealers: | 4 sheets with adhesive backing |

*Source of all serum proteins is from USDA inspected abattoirs located in the United States

Materials Required but not Provided

- Microplate reader capable of measuring absorbance at 450 nm
- Precision pipettes to deliver 50 µl and 100 µl volumes
- Adjustable 1 ml, 5 ml, 10 ml, 25 ml pipettes for reagent preparation
- Deionized or distilled water
- Wash bottle or automated microplate washer
- Log-log graph paper or computer and software for ELISA data analysis
- Tubes to prepare standard dilutions
- Laboratory timer
- Absorbent paper

Storage Information

1. Store kit at 2 – 8°C. Do not use kit after expiration date.
2. Before use, bring all reagents to room temperature (18 – 25°C). Immediately after use, return to proper storage conditions.
3. Lyophilized standards are stable until kit expiration date. After reconstitution, use freshly reconstituted standard within 12 hours (stored at 2 – 8°C).

Warnings and Precautions

1. Reagents that contain preservatives may be toxic if they are ingested, inhaled, or come in contact with skin.
2. Avoid contact of skin, eyes, or clothing with Stop Solution or Substrate Reagents.
3. Handle all serum and plasma specimens in accordance with NCCLS guidelines for preventing transmission of blood-borne infections.
4. TMB One-Step Substrate Reagent is toxic if inhaled or swallowed. Avoid contact with skin. Keep container tightly closed. In case of an accident or if you feel unwell, seek medical advice immediately.
5. Standard/Sample Diluent and ELISA Diluent contain less than 0.1% sodium azide. Sodium azide yields highly toxic hydrazoic acid under acidic conditions. Dilute azide compounds in running water before discarding to avoid accumulation of potentially explosive deposits in plumbing.
6. The Wash Concentrate contains 16% Sodium Chloride and is an irritant.
 - R36/37/38 Irritating to eyes, respiratory system, and skin.
 - S7 Keep container tightly closed.
 - S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.
 - S36/37 Wear suitable protective clothing and gloves.
 - S60 This material and its container must be disposed of as hazardous waste.
7. The Stop Solution contains 11.5% Phosphoric Acid and is a corrosive solution.
 - R34 Causes burns.
 - R36/38 Irritating to eyes and skin.
 - S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.
 - S37 Wear suitable gloves.
 - S45 In case of accident or if you feel unwell, seek medical advice immediately.
 - S60 This material and its container must be disposed of as hazardous waste.

Specimen Collection and Handling

Specimens should be clear, non-hemolyzed and non-lipemic. Samples with expected values higher than the top standard, 500 pg/ml, should be diluted with Standard/Sample Diluent prior to running the assay.

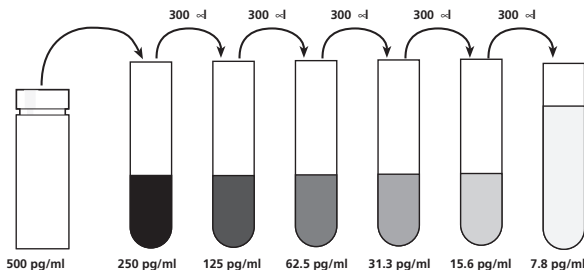
Cell culture supernatants: Remove any particulate material by centrifugation and assay immediately or store samples at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles.

Serum: Use a serum tube (eg, BD Vacutainer® Cat. No. 366430) and allow samples to clot for 30 minutes, then centrifuge for 10 minutes at $1000 \times g$. Remove serum and assay immediately or store samples at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles.

Plasma: Collect plasma using citrate, EDTA, or heparin as anticoagulant. Centrifuge for 10 minutes at $1000 \times g$ within 30 minutes of collection. Assay immediately or store samples at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles.

Reagent Preparation

1. Bring all reagents to room temperature ($18 - 25^{\circ}\text{C}$) before use.
2. Standards
 - a. Reconstitute 1 vial lyophilized Standard with required volume (noted on vial label) of Standard/Sample Diluent to prepare a 500 pg/ml stock standard. Allow the standard to equilibrate for at least 15 minutes before making dilutions. Gently vortex to mix.
 - b. Add 300 μl Standard/Sample Diluent to 6 tubes. Label as 250 pg/ml, 125 pg/ml, 62.5 pg/ml, 31.3 pg/ml, 15.6 pg/ml, and 7.8 pg/ml.
 - c. Perform serial dilutions by adding 300 μl of each standard to the next tube and vortexing between each transfer (*see figure below*). The undiluted standard serves as the high standard (500 pg/ml). The Standard/Sample Diluent serves as the zero standard (0 pg/ml).



3. Working Detector

Note: One-step incubation of Biotin/Avidin reagents.
See Assay Procedure, step 5.

4. Wash Buffer

Note: If the Wash Concentrate contains visible crystals, warm to room temperature and mix gently until dissolved. Dilute required quantity of 20× Wash Concentrate with deionized or distilled water, mix. (To prepare 2.0 L, add 100 ml Wash Concentrate to 1900 ml water. At least 500 ml solution should be prepared for a full 96-well plate).

5. TMB One-Step Substrate Reagent

No more than 15 minutes prior to use, add required volume of TMB One-Step Substrate Reagent to a clean tube or reservoir. To prevent contamination, pipette out from the tube/reservoir instead of directly from bottle. Avoid prolonged exposure to light or contact with metal, air, or extreme temperature as color may develop.

Assay Procedure

1. Bring all reagents and samples to room temperature (18 – 25°C) prior to use. It is recommended that all standards and samples be run in duplicate. A standard curve is required in each assay run.

2. Remove required quantity of test strips/wells, place in well holder.

Note: Wells are provided in breakable 8-well strips. Strips may be “broken” into individual wells, replaced in well holder, and assayed. Return any unused wells to sealed pouch for 2 – 8°C storage.

3. Pipette 50 µl of ELISA Diluent into each well.

4. Pipette 100 µl of each standard (*see Reagent Preparation, step 2*) and sample into appropriate wells. Gently shake/tap the plate for 5 seconds to mix. Cover wells with Plate Sealer and incubate for 2 hours at room temperature.

5. Prepare Working Detector. Within 15 minutes prior to use, pipette required volume of Detection Antibody into a clean tube or flask. Add in required quantity of Enzyme Concentrate (250×), vortex or mix well. For a full 96-well plate, add 48 µl of Enzyme Concentrate into 12 ml of Detection Antibody.

6. Decant or aspirate contents of wells. Wash wells by filling with at least 300 µl/well prepared Wash Buffer (*see Reagent Preparation, step 4*), followed by decanting/aspirating. Repeat wash 4 times for a total of 5 washes. After the last wash, blot plate on absorbent paper to remove any residual buffer. Complete removal of liquid is required for proper performance.

7. Add 100 µl of prepared Working Detector (*see step 5 above*) to each well. Cover wells with Plate Sealer and incubate for 1 hour at room temperature.

8. Wash wells as in Step 6, but a total of 7 times.

Note: In this final wash step, soak wells in wash buffer for 30 seconds to 1 minute for each wash. Thorough washing at this step is very important.

9. Add 100 μl of TMB One-Step Substrate Reagent to each well. Incubate plate (without Plate Sealer) for 30 minutes at room temperature in the dark.
10. Add 50 μl of Stop Solution to each well.
11. Read absorbance at 450 nm within 30 minutes of stopping reaction. If wavelength correction is available, subtract the optical density readings at 570 nm from readings at 450 nm.

Assay Procedure Summary

1. Add 50 μl ELISA Diluent to each well.
2. Add 100 μl standard or sample to each well.
Incubate 2 hours at room temperature.
3. Aspirate and wash 5 times.
4. Add 100 μl prepared Working Detector to each well.
Incubate 1 hour at room temperature.
5. Aspirate and wash/soak 7 times.
6. Add 100 μl TMB One-Step Substrate Reagent to each well.
Incubate 30 minutes at room temperature.
7. Add 50 μl Stop Solution to each well.
Read at 450 nm within 30 minutes.
 λ correction 570 nm.

Calculation of Results

Calculate the mean absorbance for each set of duplicate standards, controls and samples. Subtract the mean zero standard absorbance (ie, plate background) from each.

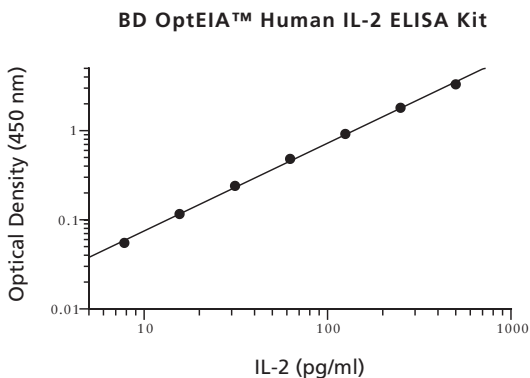
Plot the standard curve on log-log graph paper, with IL-2 concentration on the x-axis and absorbance on the y-axis. Draw the best fit straight line through the standard points.

To determine the IL-2 concentration of the unknowns, find the unknowns' mean absorbance value on the y-axis and draw a horizontal line to the standard curve. At the point of intersection, draw a vertical line to the x-axis and read the IL-2 concentration. If samples were diluted, multiply the interpolated IL-2 concentration by the dilution factor.

Computer-based curve-fitting statistical software may also be employed.

Typical Data

This standard curve is for demonstration only. A standard curve must be run with each assay.



| Concentration (pg/ml) | OD1 | OD2 | Mean | Zero Standard Subtracted |
|-----------------------|-------|-------|-------|--------------------------|
| 0 | 0.028 | 0.030 | 0.029 | 0.000 |
| 7.8 | 0.083 | 0.086 | 0.085 | 0.055 |
| 15.6 | 0.144 | 0.146 | 0.145 | 0.116 |
| 31.3 | 0.273 | 0.266 | 0.270 | 0.240 |
| 62.5 | 0.515 | 0.507 | 0.511 | 0.482 |
| 125 | 0.997 | 0.898 | 0.948 | 0.918 |
| 250 | 1.856 | 1.812 | 1.834 | 1.805 |
| 500 | 3.307 | 3.367 | 3.337 | 3.308 |

Limitations of the Procedure

1. This kit is intended for use as an integral unit. Do not mix reagents from different kit lots. Reagents from other manufacturers/other available antibody clones should not be used in this kit.
2. Interference by drug metabolites, soluble receptors, or other binding proteins in specimens has not been thoroughly investigated. The possibility of interference cannot be excluded.

Performance Characteristics

Sensitivity

The minimum detectable dose of IL-2 was determined to be 1 pg/ml. This is defined as two standard deviations above the mean optical density of 20 replicates of the zero standard.

Recovery

Three different levels of IL-2 were spiked into samples of various matrices. Results are compared with the same amounts of IL-2 spiked into Standard/Sample Diluent, as follows:

| | Spike Concentration (pg/ml) | Average % Recovery | Range |
|---------------------------------------|--|-------------------------------|--------------|
| Serum (n = 9) | 250 | 88 | 84 – 92 |
| | 125 | 86 | 70 – 107 |
| | 62.5 | 95 | 85 – 110 |
| Plasma (n = 5) | 250 | 83 | 77 – 91 |
| | 125 | 82 | 73 – 91 |
| | 62.5 | 90 | 78 – 100 |
| Cell culture media (n = 3) | 250 | 105 | 101 – 108 |
| | 125 | 116 | 110 – 120 |
| | 62.5 | 119 | 108 – 127 |

Linearity

Samples spiked with high concentrations of IL-2 were serially diluted with Standard/Sample Diluent and run in the BD OptEIA™ Human IL-2 ELISA Kit II. Results are as follows:

| Dilution | | Serum (n = 9) | Plasma (n = 5) | Cell culture media (n = 3) |
|-----------------|-----------------------|--------------------------|---------------------------|---------------------------------------|
| 1:2 | Average % of Expected | 106 | 110 | 95 |
| | Range | 101 – 111 | 105 – 116 | 90 – 100 |
| 1:4 | Average % of Expected | 107 | 108 | 96 |
| | Range | 101 – 111 | 102 – 122 | 93 – 98 |
| 1:8 | Average % of Expected | 100 | 108 | 95 |
| | Range | 93 – 106 | 101 – 119 | 92 – 101 |
| 1:16 | Average % of Expected | 100 | 109 | 91 |
| | Range | 94 – 111 | 100 – 128 | 87 – 94 |

Specificity

Cross-Reactivity: The factors listed below were spiked in Standard/Sample Diluent at 50 ng/ml to test for any cross-reactivity with the BD OptEIA™ Human IL-2 assay. No cross reactivity was identified.

Recombinant Human

sCD23, Eotaxin, Fas, GM-CSF, GRO- α , GRO- β , GRO- γ , I-309, IFN- γ , IL-3, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12 (p40), IL-12(p70), IL-15, IL-16, IP-10, MCP-1, MCP-2, MCP-3, MIG, MIP-1 α , MIP-1 β , NAP-2, PF-4, SDF-1 α , TNF- α , TNF- β

Recombinant Mouse

IFN- γ , IL-2, IL-3, IL-4, IL-5, IL-6, IL-9, IL-10, IL-12 (p40), IL-12 (p70), GM-CSF, MCP-1, MIG, MIP-1 α , MIP-1 β , TCA-3, TNF- α

Recombinant Rat

IL-2, IL-10, GM-CSF, MCP-1, RANTES

Interfering Substances: The following substances at levels > 10 mg/ml were added to Standard/Sample Diluent spiked with 100 pg/ml IL-2. No effect on assay results was observed.

- Bilirubin
- Human hemoglobin
- Human transferrin
- Triglycerides
- Heparin (600 units/ml)
- Sodium Citrate
- EDTA

Precision

Intra-assay: Twenty-four replicates each of three different levels of IL-2 were tested in one plate. The following results were observed:

| | | | |
|----------------------|-------------|-------------|------------|
| Number of replicates | 24 | 24 | 24 |
| Mean Concentration | 233.7 pg/ml | 111.0 pg/ml | 56.4 pg/ml |
| SD | 7.3 | 6.3 | 2.8 |
| %CV | 3.1 | 5.7 | 5.0 |

Inter-assay: Three different levels of IL-2 were tested in four different plates. The following results were observed:

| | | | |
|----------------------|-------------|-------------|------------|
| Number of replicates | 32 | 32 | 32 |
| Mean Concentration | 238.2 pg/ml | 119.9 pg/ml | 59.6 pg/ml |
| SD | 9.8 | 4.0 | 2.7 |
| %CV | 4.1 | 3.3 | 4.5 |

Standardization

This immunoassay is calibrated against baculovirus-expressed recombinant human IL-2 produced at BD Biosciences Pharmingen.

The NIBSC/WHO First International Standard 86/504 (natural human IL-2) was evaluated in this kit. The conversion factor for NIBSC material is as follows:

NIBSC (86/504) equivalent value (IU/ml) = 0.0096 × BD OptEIA™ IL-2 value (pg/ml)

∴ 1 µg NIBSC IL-2 = 1.37 µg BD OptEIA™ recombinant human IL-2

Expected Values

Serum: Twenty-six apparently healthy normal donors' serum samples were tested in this assay. Nineteen samples measured less than the lowest standard, 7.8 pg/ml. Seven samples measured between 8.1 – 20.8 pg/ml.

Plasma: Eight apparently healthy donors' EDTA plasma samples were tested in this assay. All eight samples measured less than the lowest standard, 7.8 pg/ml.

Cell culture supernatants: Human peripheral blood mononuclear cells or CD8⁺ cells from five apparently healthy, normal donors were stimulated with immobilized anti-human CD3, soluble anti-human CD28, recombinant human IL-2, and recombinant human IL-4 for 2 days. Cells were harvested and recultured in complete tissue culture medium containing rhIL-2 and rhIL-4 for 3 days. Finally, the *in-vitro* primed cells were harvested, washed, and restimulated with PMA and ionomycin for 4 hours, 24 hours, or 32 hours (as noted). Culture supernatants were collected and quantified for IL-2 using a BD OptEIA™ human IL-2 Kit II. The results are as follows:

| Donor No. | IL-2 (pg/ml) | |
|-----------|--------------|---|
| 1 | 7,268.6 | PBMC, 4 hr PMA/ionomycin stimulation |
| 2 | 8,658.8 | PBMC, 4 hr PMA/ionomycin stimulation |
| 3 | 10,598.1 | PBMC, 4 hr PMA/ionomycin stimulation |
| 4 | 20,757.3 | PBMC, 24 hr PMA/ionomycin stimulation |
| 5 | 23,963.3 | PBMC, 24 hr PMA/ionomycin stimulation |
| 5 | 90,070.8 | CD8 ⁺ cells, 32 hr PMA/ionomycin stimulation |

Troubleshooting

| Problem | Possible Source | Corrective Action |
|----------------------------|--|--|
| Poor Precision | <ul style="list-style-type: none"> • Inadequate washing / aspiration of wells • Inadequate mixing of reagents • Imprecise / inaccurate pipetting • Imprecise sealing of plate | <ul style="list-style-type: none"> • Check function of washing system • Ensure adequate mixing • Check / calibrate pipettes • Ensure complete sealing of plate |
| Poor Standard Curve | <ul style="list-style-type: none"> • Improper standard handling / dilution • Incomplete washing / aspiration of wells • Imprecise / inaccurate pipetting | <ul style="list-style-type: none"> • Ensure correct preparation of standards • Check function of washing system • Check / calibrate pipettes |
| Low Signal | <ul style="list-style-type: none"> • Inadequate reagent volumes added to wells • Incorrect incubation times / temperature • Overly high wash / aspiration pressure from automated plate-washer. | <ul style="list-style-type: none"> • Check / calibrate pipettes • Ensure sufficient incubation times / reagents warmed to room temperature • Utilize manual washing |

References:

1. Taniguchi, T., *et al.* 1983. Structure and expression of a cloned cDNA for human interleukin-2. *Nature*. 302:305-307.
2. Smith, K.A. 1988. Interleukin-2: inception, impact, and implications. *Science* 240:1169-1176.
3. Stern, A.S., *et al.* 1984. Purification to homogeneity and partial characterization of interleukin-2 from a human T-cell leukemia. *Proc. Natl. Acad. Sci. USA* 81:871-875.
4. Gillis, S., *et al.* 1978. T cell growth factor: Parameters of production and a quantitative microassay for activity. *J. Immunol.* 120:2027-2032.
5. Prussin, C., and D. D. Metcalfe. 1995. Detection of intracytoplasmic cytokine using flow cytometry and directly conjugated anti-cytokine antibodies. *J. Immunol. Meth.* 188:117-28.

Plate Templates

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

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

United States

877.232.8995

Canada

888.259.0187

Europe

32.53.720.211

Japan

0120.8555.90

Asia/Pacific

65.6861.0633

Latin America/Caribbean

55.11.5185.9995



BD Biosciences

10975 Torreyana Road

San Diego, CA 92121

Customer/Technical Service

Tel 877.232.8995 (US)

Fax 800.325.9637

www.bdbiosciences.com