

Human IL-1 β ELISA Kit

EH2IL1B EH2IL1B2 EH2IL1B5

1372.7

Number	Description
EH2IL1B	Human Interleukin-1 beta (IL-1β) ELISA Kit , sufficient reagents for 96 determinations
EH2IL1B2	Human Interleukin-1 beta (IL-1β) ELISA Kit , sufficient reagents for 2 \times 96 determinations
EH2IL1B5	Human Interleukin-1 beta (IL-1β) ELISA Kit , sufficient reagents for 5 \times 96 determinations

Kit Contents	EH2IL1B	EH2IL1B2	EH2IL1B5
Anti-human IL-1 β Precoated 96-well Strip Plate	1 each	2 each	5 each
Lyophilized Recombinant Human IL-1 β	2 vials	4 vials	5 vials
Standard Diluent	12mL	2 \times 12mL	75mL
Biotinylated Antibody Reagent	8mL	2 \times 8mL	35mL
30X Wash Buffer	50mL	2 \times 50mL	200mL
Streptavidin-HRP Concentrate	75 μ L	2 \times 75 μ L	250 μ L
Streptavidin-HRP Dilution Buffer	14mL	2 \times 14mL	70mL
TMB Substrate	13mL	2 \times 13mL	5 \times 13mL
Stop Solution, contains 0.16M sulfuric acid	13mL	2 \times 13mL	55mL
Adhesive plate covers	6 each	12 each	30 each

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

Storage: For maximum stability, store in a non-defrosting -20 $^{\circ}$ C freezer and refer to the expiration date for frozen storage on the label. Alternatively, store at 2-8 $^{\circ}$ C and refer to the expiration date for refrigerated storage. Once thawed, store at 4 $^{\circ}$ C until the expiration date for refrigerated storage. Kit is shipped on dry ice.












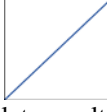
Table of Contents

Introduction	1
Procedure Summary.....	2
Additional Materials Required.....	2
Precautions.....	2
Sample Preparation.....	3
Reagent Preparation.....	3
Assay Procedure	4
Performance Characteristics	6
Cited Reference	7
Data Templates	8

Introduction

The Thermo Scientific Human Interleukin-1 beta (IL-1 β) ELISA is an *in vitro* enzyme-linked immunosorbent assay for measuring human IL-1 β in culture supernatants, plasma (heparin and sodium citrate), serum and urine.

Procedure Summary

- | | | | |
|--|---|---|--|
| 
1. Add 50µL of standards or samples to each well in duplicate. | 
2. Add 50µL of Biotinylated Antibody Reagent to each well. | 
3. Cover plate and incubate at room temperature (20-25°C) for 3 hours. | 
4. Wash plate THREE times. |
| 
5. Add 100µL of prepared Streptavidin-HRP Solution to each well. | 
6. Cover plate and incubate at room temperature for 30 minutes. | 
7. Wash plate THREE times. | 
8. Add 100µL of TMB Substrate to each well. |
| 
9. Develop plate in the dark at room temperature for 30 minutes. | 
10. Add 100µL of Stop Solution to each well. | 
11. Measure absorbance on a plate reader at 450nm minus 550nm. | 
12. Calculate results using graph paper or curve-fitting statistical software. |

Additional Materials Required

- Precision pipettors with disposable plastic tips to deliver 5-1000µL and plastic pipettes to deliver 5-15mL
- A glass or plastic two-liter container to prepare Wash Buffer
- A squirt wash bottle or an automated 96-well plate washer
- 1.5mL polypropylene or polyethylene tubes to prepare standards – do not use polystyrene, polycarbonate or glass tubes
- Disposable reagent reservoirs (Product No. 15075)
- 15mL plastic tube to prepare Streptavidin-HRP Solution
- A standard ELISA reader for measuring absorbance at 450nm and 550nm. If a 550nm filter is not available, the absorbance can be measured at 450nm only. Refer to the instruction manual supplied with the instrument being used.
- Graph paper or a computerized curve-fitting statistical software package

Precautions

- All samples and reagents must be at room temperature (20-25°C) before use in the assay.
- Review all instructions carefully and verify components against the Kit Contents list (page 1) before beginning the assay.
- Do not use a water bath to thaw samples. Thaw samples at room temperature.
- When preparing standard curve and sample dilutions in culture medium, use the same medium used to culture the cells. For example, if RPMI with 10% fetal calf serum (FCS) was used to culture cells, then use RPMI with 10% FCS to dilute the standard and samples. Do NOT use RPMI without serum supplement.
- If using a multichannel pipettor, always use a new disposable reagent reservoir.
- Use new disposable pipette tips for each transfer to avoid cross-contamination.
- Use a new adhesive plate cover for each incubation step.
- Once reagents have been added to the plate, take care NOT to let plate DRY at any time during the assay.
- Avoid microbial contamination of reagents.
- Vigorous plate washing is essential.
- Avoid exposing reagents to excessive heat or light during storage and incubation.

-
- Discard unused kit components. Do not mix reagents from different kit lots.
 - Do not use glass pipettes to measure TMB Substrate. Take care not to contaminate the solution. If the solution is blue before use, **DO NOT USE IT**.
 - Individual components may contain antibiotics and preservatives. Wear gloves while performing the assay to avoid contact with samples and reagents. Please follow proper disposal procedures.

Additional Precautions for the 5-plate kits

- Dispense and equilibrate to room temperature only reagent volumes required for the number of plates being used. Do not combine leftover reagents with those reserved for additional plates.
- Use only one bottle of the TMB Substrate per 96-well plate. Do not combine leftover substrate with that reserved for other plates.
- Use only one vial of standard per 96-well plate.

Sample Preparation

Sample Handling

- Serum, heparin and sodium citrate plasma, urine, or culture supernatants may be tested in this ELISA.
Note: EDTA interferes with human IL-1 β in this ELISA. Use only plasma collected with sodium citrate or heparin.
- 50 μ L per well of serum, plasma, urine or culture supernatant are required.
- Store samples to be assayed within 24 hours at 2-8°C. For long-term storage, aliquot and freeze samples at -70°C. Avoid repeated freeze-thaw cycles when storing samples.
- Samples and standards must be assayed in duplicate each time the ELISA is performed.
- Gradually equilibrate samples to room temperature before beginning assay. Do not use a heated water bath to thaw or warm samples.
- Mix samples by gently inverting tubes.
- If samples are clotted, grossly hemolyzed, lipemic or contaminated, or if there is any question about the integrity of a sample, make a note on the template and interpret results with caution.

Sample Dilution

- If the human IL-1 β concentration possibly exceeds the highest point of the standard curve (i.e., 400pg/mL), prepare one or more 10-fold dilutions of the sample. When testing culture supernatants, prepare serial dilutions using the culture medium. When testing serum, plasma or urine, prepare serial dilutions using the Standard Diluent provided. A 10-fold dilution is prepared by adding 50 μ L of sample to 450 μ L of appropriate diluent. Mix thoroughly between dilutions.

Reagent Preparation

For procedural differences when using partial plates, look for **(PP)** throughout these instructions.

Note: When using the 5-plate kit, only one standard per plate is supplied, therefore, partial plates cannot be used.

Wash Buffer

Note: Wash Buffer must be at room temperature before use in the assay. Do not use Wash Buffer if it becomes visibly contaminated during storage.

1. Label a clean glass or plastic 2L container "Wash Buffer." The 30X Wash Buffer may have a cloudy appearance.
2. If using a 5-plate kit, add 30mL Wash Buffer to 870mL water for each plate used, otherwise, add the entire contents of the 30X Wash Buffer bottle (50mL) to the container. Dilute buffer to a final volume of 1.5L with ultrapure water and mix thoroughly.

(PP) When using partial plates, store the reconstituted Wash Buffer at 2-8°C.

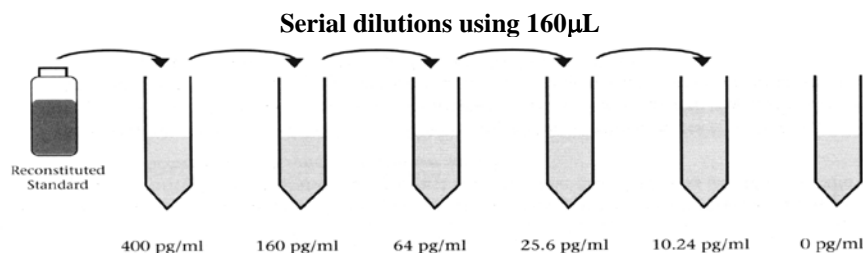
Standards

- **(PP)** Reconstitute and use one vial of the lyophilized standard per partial plate.
 - Prepare standards just before use and use within one hour of reconstitution. Do not store reconstituted standards.
1. When testing **culture supernatant samples**, reconstitute standard with ultrapure water. Reconstitution volume is stated on the standard vial label. The standard will dissolve in approximately 1 minute. Mix by gently inverting vial. Use the sample culture medium to prepare standard curve serial dilutions.

When testing **serum, plasma or urine samples**, reconstitute standard with ultrapure water. Reconstitution volume is stated on the standard vial label. The standard will dissolve in approximately 1 minute. Mix by gently inverting vial. Use the Standard Diluent provided to prepare standard curve serial dilutions.

When testing **serum, plasma or urine and cell culture supernatant samples on the same plate**, validate the medium to establish if the same standard curve can be used for the different sample types. Prepare a standard curve (including a zero/blank) using culture medium to reconstitute and dilute the standard. Use medium containing serum or other protein to maximize stability of the human IL-1 β . Perform this curve in parallel with a standard curve reconstituted in ultrapure water and diluted in the Standard Diluent provided. If OD values are within 10% of the mean for both curves, then the assay can be performed with Standard Diluent, whether testing culture supernatant, plasma, urine or serum samples.

2. Label six tubes, one for each standard curve point: 400, 160, 64, 25.6, 10.24, and 0pg/mL, then prepare 1:2.5 serial dilutions for the standard curve as follows:
3. Pipette 240 μ L of appropriate diluent into each tube.
4. Pipette 160 μ L of the reconstituted standard into the first tube (i.e., 400pg/mL) and mix.
5. Pipette 160 μ L of this dilution into the second tube (i.e., 160pg/mL) and mix.
6. Repeat the serial dilutions (using 160 μ L) three more times to complete the standard curve points.



Assay Procedure

A. Sample and Biotinylated Antibody Reagent Incubation

- **(PP)** Determine number of strips required and leave these strips in the plate frame. Tightly seal the remaining unused strips in the provided foil pouch with desiccant and store at 2-8°C. After completing assay, retain plate frame for second partial plate. When using the second partial plate, place strips securely in the plate frame.
 - Use the Data Template provided to record locations of the zero standard (blank or negative control), human IL-1 β standards and samples. Perform five standard points and one blank in duplicate with each series of unknown samples.
 - If using a multichannel pipettor, use a new reagent reservoir to add the Biotinylated Antibody Reagent. Remove from the vial only the amount required for the number of strips being used. Take care not to touch the samples in wells with the pipette tip when adding the Biotinylated Antibody Reagent.
1. Add 50 μ L of reconstituted standards or test samples in duplicate to each well.

Note: If the human IL-1 β concentration in any test sample possibly exceeds the highest point on the standard curve, 400pg/mL, see Sample Preparation – Sample Dilution section.

2. Add 50 μ L of Biotinylated Antibody Reagent to each well. Mix well by gently tapping the plate several times.
3. Add 50 μ L of Standard Diluent to all wells that do not contain standards or samples.

- Carefully cover plate with an adhesive plate cover. Ensure all edges and strips are sealed tightly by running your thumb over edges and down each strip. Incubate for three (3) hours at room temperature, 20-25°C.
- Carefully remove adhesive plate cover. Wash plate as described in the Plate Washing Section (section B).

B. Plate Washing

- Gently squeeze the long sides of the plate frame before washing to ensure all strips securely remain in the frame.
- Discard plate contents. Use a squirt wash bottle to vigorously fill each well completely with Wash Buffer, then discard plate contents. Repeat procedure four additional times for a total of FIVE washes. Blot plate onto paper towels or other absorbent material.

Note: Automated plate washers may produce sub-optimal results. For best results, perform a manual wash the first time the assay is performed. Automated plate washing may require validation to determine the number of washes necessary to achieve optimal results; typically, three washes are sufficient. For automated washing, aspirate all wells and wash by overfilling wells with Wash Buffer. Blot plate onto paper towels or other absorbent material.

C. Streptavidin-HRP Solution Preparation and Incubation

- Prepare Streptavidin-HRP Solution **immediately before use**. Do not store prepared Streptavidin-HRP Solution and do not prepare more Streptavidin-HRP Solution than required.
- Use a 15mL plastic tube to prepare Streptavidin-HRP Solution.

Note: If using a multichannel pipettor, use new reagent reservoir and pipette tips when adding the prepared Streptavidin-HRP Solution.

- Briefly centrifuge the Streptavidin-HRP Concentrate to force entire vial contents to the bottom.
- (PP)** Use only the Streptavidin-HRP Solution amount required for the number of strips being used. For each strip, mix 2.5µL of Streptavidin-HRP Concentrate with 1mL of Streptavidin-HRP Dilution Buffer. Store Streptavidin-HRP Concentrate reserved for additional strips at 2-8°C.

For one complete 96-well plate, add 30µL of Streptavidin-HRP Concentrate to 12mL of Streptavidin-HRP Dilution Buffer and mix gently.

- Add 100µL of prepared Streptavidin-HRP Solution to each well.
- Carefully attach a new adhesive plate cover, ensuring all edges and strips are tightly sealed. Incubate plate for 30 minutes at room temperature, 20-25°C.
- Carefully remove the plate cover and discard plate contents. Wash plate as described in the Plate Washing Section.

D. Substrate Incubation and Stop Step

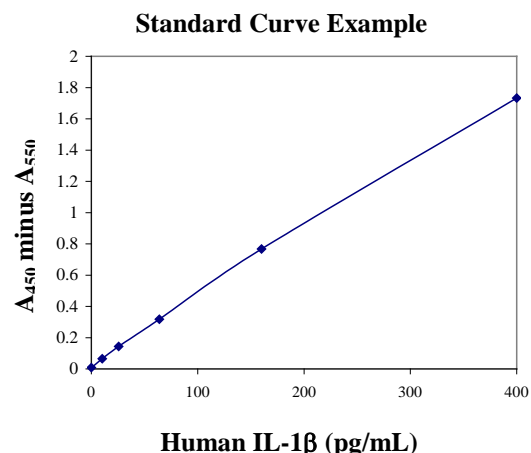
- Use new disposable reagent reservoirs when adding TMB Substrate and Stop Solution.
 - Dispense from bottle **ONLY** amount required, 100µL per well, for the number of wells being used. Use a plastic pipette (i.e., do not use glass) to measure the TMB Substrate.
 - (PP)** Do not combine leftover substrate with that reserved for the second partial plate. Take care not to contaminate remaining TMB Substrate.
- Pipette 100µL of TMB Substrate into each well.
 - Allow color reaction to develop at room temperature in the dark for 30 minutes. Do not cover plate with aluminum foil or a plate sealer. The substrate reaction yields a blue solution that turns yellow when Stop Solution is added.
 - After 30 minutes, stop the reaction by adding 100µL of Stop Solution to each well.

E. Absorbance Measurement

- Evaluate plate within 30 minutes of stopping the reaction. Measure absorbance on an ELISA plate reader set at 450nm and 550nm. Subtract 550nm values from 450nm values to correct for optical imperfections in the microplate. If 550nm is not available, measure absorbance at 450nm only.
- **Note:** Omitting the 550nm measurement will result in greater absorbance values.

F. Results Calculation

- Generate the standard curve by plotting the average absorbance obtained for each standard concentration on the vertical (Y) axis vs. the corresponding human IL-1 β concentration (pg/mL) on the horizontal (X) axis.
- Calculate results using graph paper or curve-fitting statistical software. Determine the IL-1 β amount in each sample by interpolating from the absorbance value (Y axis) to IL-1 β concentration (X axis) using the standard curve.
- Multiply the interpolated value obtained from the standard curve by the dilution factor to calculate amount of human IL-1 β in the sample.
- Absorbance values obtained for duplicates should be within 10% of the mean value. Carefully consider duplicate values that differ from the mean by greater than 10%.



Performance Characteristics

Sensitivity: <1pg/mL

The sensitivity or Lower Limit of Detection (LLD)¹ is determined by assaying replicates of zero and the standard curve. The mean signal of zero + 2 standard deviations read in dose from the standard curve is the LLD. This value is the smallest dose that is not zero with 95% confidence.

Specificity: This ELISA is specific for the measurement of natural and recombinant human IL-1 β . This ELISA does not cross-react with human IL-1 α , IL-1RA, IL-2, IL-3, IL-4, IL-6, IL-7, IL-8, TNF α , IFN α , IFN γ or mouse IL-1 β .

Assay Range: 10.24-400pg/mL

Suggested standard curve points are 400, 160, 64, 25.6, 10.24 and 0pg/mL.

Reproducibility:

Intra-assay CV: < 10%

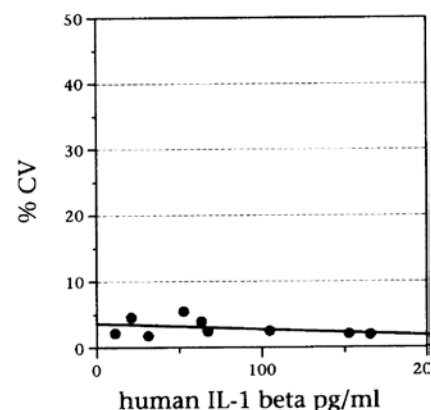
Inter-assay CV: < 10%

Calibration: The standards in this assay have been re-calibrated to the NIBSC reference standard lot 861/680. One (1) pg of Endogen standard = 2 NIBSC pg = 0.2 NIBSC units.

Precision: The intra-assay coefficient of variation is plotted against IL-1 β concentration (pg/mL). The points represent samples evaluated in replicates of four in six different kit lots.

Expected Values: The average level of human IL-1 β obtained in normal serum, plasma and urine samples is 0pg/mL.

Precision Profile of the hIL-1 β ELISA

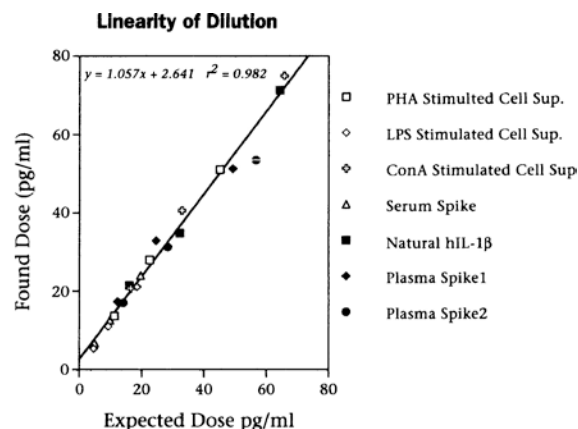


<u>Sample Type</u>	<u>Average</u>	<u>Range</u>
Urine (n=9)	1.1pg/mL	0-7.1pg/mL
Serum (n=9)	< 1pg/mL	< 1pg/mL
Plasma (n=9)	< 1pg/mL	< 1pg/mL

Recovery: Recovery was determined by spiking three different levels of recombinant human IL-1 β into human serum, sodium citrate plasma and urine samples collected from apparently healthy individuals. Mean recoveries are as follows:

<u>Spike Level</u>	<u>15pg/mL</u>	<u>40pg/mL</u>	<u>80pg/mL</u>
Serum Recovery	81%	82%	82%
Plasma Recovery	69%	74%	72%
Urine Recovery	87%	86%	85%

Linearity of Dilution: Linearity of Dilution is determined by serially diluting seven different positive samples. The dilutions are evaluated in the ELISA and the “found” doses are plotted against the “expected” doses.



Cited Reference

1. *Immunoassay: A Practical Guide*, Chan and Perlstein, Eds., 1987, Academic Press: New York, p71.

This product (“Product”) is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts (“Documentation”) and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product (“Buyer”).

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer’s exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.

Data 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												