

# Human IL-1 $\beta$ ELISA Reagent Kit

ESS0008

2473.0

**Number**

ESS0008

**Description**

**Human IL-1 $\beta$  ELISA Reagent Kit**, pre-titered coating and detection antibodies, recommended buffers and specific assay protocol optimized for the quantitative measurement of human IL-1 $\beta$  in cell culture supernatants

Kit provides sufficient reagents for approximately five 96-well plates, provided the Human IL-1 $\beta$  ELISA Reagent Kit Protocol is followed.

Kit Contents	Size	Assay Dilution
Anti-human IL-1 $\beta$ Coating Antibody	0.625mL	1:100
Lyophilized Recombinant Human IL-1 $\beta$ Standard	5 vials	See vial label
Anti-human IL-1 $\beta$ Detection Antibody	0.625mL	1:100
Streptavidin-HRP	0.25mL	1:400
Substrate Solution	55mL	Ready to use
Stop Solution, 0.16M Sulfuric Acid	55mL	Ready to use

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

**Storage:** Immediately upon receipt, aliquot and freeze the coating and detecting antibodies at  $\leq -20^{\circ}\text{C}$  in a manual defrost freezer (125 $\mu\text{L}$ /tube). Avoid repeated freeze-thaw cycles. Store all other components at 2-8 $^{\circ}\text{C}$ . Kit is shipped on dry ice.

**Table of Contents**

Introduction .....	1
ELISA Reagent Kit Buffers .....	1
Additional Materials Required .....	1
Assay Protocol .....	2
Performance Characteristics .....	3
General References .....	3

**Introduction**

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

**Materials Required**

- 8-well strip plates, clear, corner-notched (Product No. 15031)
- Plate sealers for 96-well plates (Product No. 15036)
- Reagent reservoir, sterile, 50mL capacity, 40pk (Product No. 15075)

---

## ELISA Reagent Kit Buffers

- D-PBS: 0.008M sodium phosphate, 0.002M potassium phosphate, 0.14M sodium chloride, 0.01M potassium chloride, pH 7.4, 0.2µm filtered (e.g., Thermo Scientific™ BupH™ Modified Dulbecco's Phosphate Buffered Saline Packs, Product No. 28374)
- Blocking Buffer: 4% bovine serum albumin (BSA), 5% sucrose in D-PBS, 0.2µm filtered OR ELISA Blocker Blocking Buffer, Product No. N502
- Reagent Diluent: 4% BSA in D-PBS (pH 7.4), 0.2µm filtered OR Thermo Scientific™ StartingBlock™ (PBS) Blocking Buffer, Product No. 37538
- Wash Buffer: 0.05% Tween™-20 Detergent (e.g., 0.5% Thermo Scientific™ Surfact-Amps™ 20 Detergent Solution, Product No. 28320) in D-PBS, pH 7.4 OR ELISA Wash Buffer (30X), Product No. N503

**Note:** Mix new solution daily.

## Assay Protocol

Kit components are titrated to give optimal results using the Human IL-1β ELISA Reagent Kit Protocol for cell culture supernatants. Any change, including component concentration, volumes, incubation times or temperatures, buffer content or number of wash steps may significantly affect the ELISA results and require optimization to give the best results.

**Note:** Allow all reagents and buffers to equilibrate to room temperature (22-25°C) before use. Thaw one aliquot of coating and detecting antibody for each plate. Do not use a water bath.

### A. Plate Preparation

1. Dilute the Coating Antibody 1:100 in D-PBS buffer by adding 110µL of Coating Antibody to 10.89mL of D-PBS buffer.
2. Add 100µL of diluted Coating Antibody to each well. Cover plate with plate sealer and incubate overnight at room temperature.
3. Aspirate Coating Antibody solution and add 300µL of Blocking Buffer to each well. Cover plate with plate sealer and incubate for 1 hour at room temperature.
4. Aspirate Blocking Buffer and proceed to assay or allow to dry overnight at room temperature. When sealed with dessicant, plates can be stored at 2-8°C for 6 months.

### B. Assay Procedure

1. Reconstitute standard with Reagent Diluent with volume stated on vial label. The concentration of the reconstituted standard is 1000pg/mL.
2. Dilute reconstituted standard 1:2 in Reagent Diluent to prepare top Standard (500pg/mL). Using Reagent Diluent, prepare 1:2 serial dilutions of top Standard and dilute any supernatant expected to read above the top standard. Add 100µL of sample or Standard to each well. Cover plate with plate sealer and incubate for 1 hour at room temperature.
3. Aspirate and wash three times with Wash Buffer using 300µL per well.
4. Dilute the Detection Antibody 1:100 in Reagent Diluent by adding 110µL of Detection Antibody to 10.89mL of Reagent Diluent.
5. Add 100µL of Detection Antibody to each well. Cover plate with plate sealer and incubate for 1 hour at room temperature.
6. Aspirate and wash three times with Wash Buffer, using 300µL per well.
7. Dilute Streptavidin-HRP 1:400 in Reagent Diluent by adding 30µL of Streptavidin-HRP to 12mL of Reagent Diluent.
8. Add 100µL of diluted Streptavidin-HRP reagent to each well. Cover plate with plate sealer and incubate for 30 minutes at room temperature.
9. Aspirate and wash three times with Wash Buffer, using 300µL per well.
10. Add 100µL of Substrate Solution to each well. Cover plate with plate sealer and incubate in the dark for 20 minutes at room temperature.

11. Stop the reaction by adding 100µL of Stop Solution to each well.
12. Measure the absorbance at  $A_{450}$  minus  $A_{550}$ .

### C. Absorbance Measurement

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 an absorbance at 550nm is not available, measure the absorbance at 450nm only.

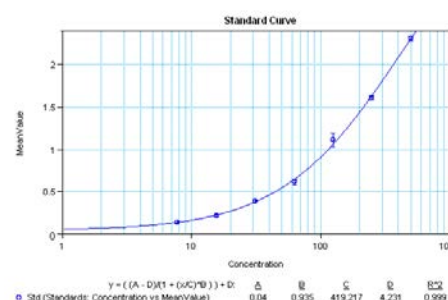
**Note:** When the 550nm measurement is omitted, absorbance values will be higher.

**Note:** Evaluate the plate within 30 minutes of stopping the reaction.

### D. Calculation of Results

- The standard curve is used to determine human IL-1β amount in an unknown sample. Generate the standard curve by plotting the average absorbance obtained for each Standard concentration on the vertical (Y) axis vs. the corresponding human IL-1β concentration (pg/mL) on the horizontal (X) axis.
- Calculate results using graph paper or curve-fitting statistical software. Determine the human IL-1β amount in each sample by interpolating from the absorbance value (Y-axis) to human IL-1β concentration (X-axis) using the standard curve.
- If the test sample was diluted, multiply the interpolated value obtained from the standard curve by the dilution factor to calculate pg/mL of human IL-1β 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%.

**Standard Curve Example**



**Standard curve based on data obtained using the Human IL-1β ELISA Reagent Kit Protocol.**

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

## Performance Characteristics

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

## General References

*Immunoassay: A Practical Guide.* Chan and Perlstein, Eds. (1987). Academic Press: New York. p.71.

Products are 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"). 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 Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

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 REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR 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.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

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

© 2013 Thermo Fisher Scientific Inc. All rights reserved. Tween is a trademark of Croda International PLC. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.