

Version 8b Last updated 3 July 2020

ab46052

Human IL-1 beta ELISA

Kit

For the qualitative and quantitative measurement of IL-1 beta in supernatants, buffered solutions or serum and plasma samples.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

| | |
|-------------------------------------|----|
| 1. Overview | 1 |
| 2. Protocol Summary | 2 |
| 3. Precautions | 3 |
| 4. Storage and Stability | 3 |
| 5. Limitations | 4 |
| 6. Materials Supplied | 4 |
| 7. Materials Required, Not Supplied | 5 |
| 8. Technical Hints | 6 |
| 9. Reagent Preparation | 8 |
| 10. Standard Preparation | 10 |
| 11. Sample Preparation | 11 |
| 12. Plate Preparation | 12 |
| 13. Assay Procedure | 13 |
| 14. Calculations | 15 |
| 15. Typical Data | 16 |
| 16. Typical Sample Values | 17 |
| 17. Assay Specificity | 18 |
| 18. Troubleshooting | 19 |
| 19. Notes | 21 |

1. Overview

Abcam's Human IL-1 beta *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-1 beta in Human serum, plasma, buffered solutions or cell culture medium.

A monoclonal antibody specific for IL-1 beta has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-1 beta concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-1 beta are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-1 beta present in the samples.

This kit will recognize both endogenous and recombinant Human IL-1 beta.

2. Protocol Summary

Remove appropriate number of antibody coated well strips



Equilibrate all reagents to room temperature



Prepare all the reagents, samples, and standards as instructed



Add standard or sample to each well used



Add prepared Biotinylated labeled detector antibody. Incubate at room temperature



Aspirate and wash each well. Add prepared Streptavidin-HRP mix to each well. Incubate at room temperature



Aspirate and wash each well. Add the TMB Solution to each well until color develops and then add the Stop Solution. Immediately begin recording the color development

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

4. Storage and Stability

Store kit at +4°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

| Item | Quantity | | Storage Condition |
|---|--------------|--------------|-------------------|
| | 1 x 96 tests | 2 x 96 tests | |
| IL-1 beta Microplate (12 x 8 well strips) | 96 wells | 2 x 96 wells | +2-8°C |
| IL-1 beta Standard (Lyophilized) | 2 vials | 4 vials | +2-8°C |
| 10X Standard Diluent Buffer | 15 mL | 25 mL | +2-8°C |
| Biotinylated anti-IL-1 beta | 400 µL | 2 x 400 µL | +2-8°C |
| Biotinylated Antibody Diluent | 7.5 mL | 13 mL | +2-8°C |
| Streptavidin-HRP | 2 x 5 µL | 4 x 5 µL | +2-8°C |
| HRP Diluent | 12 mL | 23 mL | +2-8°C |
| 200X Wash Buffer | 10 mL | 2 x 10 mL | +2-8°C |
| Chromogen TMB Substrate Solution | 11 mL | 24 mL | +2-8°C |
| Stop Reagent | 11 mL | 2 x 11 mL | +2-8°C |
| Control | 2 vials | 4 vials | +2-8°C |
| Standard Diluent (Serum) | 7 mL | 2 x 7 mL | +2-8°C |

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 2 μ L to 1 mL volumes.
- Adjustable 1-25 mL pipettes for reagent preparation.
- 100 mL and 1 liter graduated cylinders.
- Absorbent paper.
- Distilled or deionized water.
- Tubes to prepare standard or sample dilutions.
- Log-log graph paper or computer and software for ELISA data analysis.

8. Technical Hints

- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.
- Since exact conditions may vary from assay to assay, a standard curve must be established for every assay performed.
- Bacterial or fungal contamination of either samples or reagents or cross-contamination between reagents may cause erroneous results.
- Disposable pipette tips, flasks or glassware are preferred, reusable glassware must be washed and thoroughly rinsed of all detergents before use.
- Improper or insufficient washing at any stage of the procedure will result in either false positive or false negative results. Completely empty wells before dispensing fresh 1X Wash Buffer. Do not allow wells to sit uncovered or dry for extended periods.
- Kit components should be stored as indicated. All the reagents should be equilibrated to room temperature before use. Reconstituted standards should be discarded after use.
- Once the desired number of strips has been removed, immediately reseal the bag to protect the remaining strips from degradation.
- Use a clean disposable plastic pipette tip for each reagent, standard, or specimen addition in order to avoid cross-contamination; for the dispensing of the Stop Solution and substrate solution, avoid pipettes with metal parts.
- Thoroughly mix the reagents and samples before use by agitation or swirling.
- All residual washing liquid must be drained from the wells by efficient aspiration or by decantation followed by tapping the plate forcefully on absorbent paper. Never insert absorbent paper directly into the wells.
- The TMB solution is light sensitive. Avoid prolonged exposure to light. Also, avoid contact of the TMB solution with metal to prevent color development. Warning TMB is toxic avoid direct contact with hands. Dispose of properly.

- If a dark blue color develops within a few minutes after preparation, this indicates that the TMB solution has been contaminated and must be discarded. Read absorbances within 1 hour after completion of the assay.
- When pipetting reagents, maintain a consistent order of addition from well-to-well. This will ensure equal incubation times for all wells.
- Dispense the TMB solution within 15 minutes following the washing of the microtiter plate.

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells.
- Prepare only as much reagent as is needed on the day of the experiment.

9.1 1X Standard Diluent Buffer

Dilute the 10X Standard Diluent Buffer 10-fold in distilled water before use.

9.2 1X Wash Buffer

Dilute the 200X Wash Buffer Concentrate 200-fold in distilled water before use. Mix gently to avoid foaming.

The 1X Wash Buffer can be prepared as needed according to the following table:

| Number of well strips used | Volume of 200X Wash Buffer Concentrate (mL) | Volume of distilled water (mL) |
|----------------------------|---|--------------------------------|
| 1-6 | 5 | 995 |
| 1-12 | 10 | 1,990 |

9.3 1X Control Solution

Lyophilized Control vials must be reconstituted with the most appropriate diluent for your test samples:
For serum and plasma samples: Use Standard Diluent (Serum).
For cell culture supernatants: Use 1X Standard Diluent Buffer.

Control vials must be reconstituted with the volume of 1X Standard Diluent Buffer that is indicated on the vial. Reconstitution of the lyophilized material with the indicated volume will yield a solution for which the IL-1 beta concentration is stated on the vial. Allow the reconstituted 1X Control Solution to stand for 5 minutes with gentle swirling prior to use in the assay procedure. **Do not store the 1X Control Solution after reconstitution.**

9.4 1X Biotinylated anti-IL-1 beta

Prepare the 1X Biotinylated anti-IL-1 beta immediately prior to use. According to the table below, dilute the Biotinylated anti-IL-1 beta with the Biotinylated Antibody Diluent based on the number of wells being used in the assay procedure:

| Number of well strips used | Volume of Biotinylated anti-IL-1 beta (μL) | Volume of Biotinylated Antibody Diluent (μL) |
|----------------------------|---|---|
| 2 | 40 | 1,060 |
| 3 | 60 | 1,590 |
| 4 | 80 | 2,120 |
| 6 | 120 | 3,180 |
| 12 | 240 | 6,360 |

9.5 1X Streptavidin-HRP Solution

Add 500 μL of HRP-Diluent to the Streptavidin-HRP vial prior to use to create a Streptavidin-HRP Concentrate. Do not keep this solution for further experiments.

Subsequently, prior to use in the assay procedure, prepare the 1X Streptavidin-HRP Solution by further diluting the Streptavidin-HRP Concentrate with HRP-Diluent. Use the table below to determine the volumes of each solution required to prepare the final 1X Streptavidin-HRP Solution:

| Number of well strips used | Volume of Streptavidin HRP (μL) | Volume of HRP Diluent (mL) |
|----------------------------|--|----------------------------|
| 2 | 30 | 2 |
| 3 | 45 | 3 |
| 4 | 60 | 4 |
| 6 | 75 | 5 |
| 12 | 150 | 10 |

10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

10.1 Standard vials must be reconstituted with the appropriate diluent for your samples.

For serum and plasma samples: Use Standard Diluent (Serum).

For cell culture supernatants: Use 1X Standard Diluent Buffer.

Prepare a 500 pg/mL **Standard #1** by reconstituting with the volume indicated on the vial using the appropriate standard diluent from above.

10.2 Label tubes #2-6 and add 100 µL of the appropriate standard diluent into each tube.

10.3 Prepare **Standard #2** by adding 100 µL of Standard #1 to tube #2 and mix thoroughly.

10.4 Prepare **Standard #3** by adding 100 µL of Standard #2 to tube #3 and mix thoroughly.

10.5 Using the table below as a guide, prepare further serial dilutions.

10.6 The appropriate standard diluent serves as the zero standard (0 pg/mL).

| Standard # | Volume to dilute (µL) | Diluent (µL) | Total Volume (µL) | Starting conc. (pg/mL) | Final conc. (pg/mL) |
|------------|-----------------------|--------------|-------------------|------------------------|---------------------|
| 1 | - | - | - | 500 | 500 |
| 2 | 100 | 100 | 200 | 500 | 250 |
| 3 | 100 | 100 | 200 | 250 | 125 |
| 4 | 100 | 100 | 200 | 125 | 62.5 |
| 5 | 100 | 100 | 200 | 62.5 | 31.25 |
| 6 | 100 | 100 | 200 | 31.25 | 15.6 |

11. Sample Preparation

11.1 Preparation of Cell culture Supernatants:

Centrifuge cell culture media at 1,000 x g for 10 minutes to remove debris. Collect supernatants and assay. Store samples at -20°C or below. Avoid repeated freeze-thaw cycles.

11.2 Preparation of Plasma Samples:

Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 1,000 x g for 30 minutes. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.3 Preparation of Serum Samples:

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 1,000 x g for 10 minutes and collect serum. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

12. Plate Preparation

- The 96 well plate strips included with this kit is supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused well strips should be returned to the plate packet and stored at 4°C.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

13. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
- We recommend that you assay all standards, controls and samples in duplicate.

13.1 Prior to use, mix all reagents thoroughly taking care not to create any foam within the vials.

13.2 Determine the number of microplate strips required to test the desired number of samples, plus appropriate number of wells needed for controls and standards. Remove sufficient microplate strips from the pouch.

13.3 Add 100 μL of each standard (see Section 10), including blank controls to the appropriate wells.

13.4 Add 100 μL of sample and 1X Control Solution to the appropriate wells.

13.5 Add 50 μL of 1X Biotinylated anti-IL-1 beta to all wells (see Section 9).

13.6 Cover and incubate for 3 hours at room temperature (18-25°C).

13.7 Remove the cover and wash the plate as follows:

13.7.1 Aspirate the liquid from each well.

13.7.2 Add 300 μL of 1X Wash Buffer into each well.

13.7.3 Aspirate the liquid from each well.

13.7.4 Repeat for a total of 3 washes.

13.8 Add 100 μL of 1X Streptavidin-HRP solution into all wells, including the blank wells. Re-cover and incubate at room temperature for 30 minutes.

13.9 Wash as described in Step 13.7.

13.10 Add 100 μL of Chromogen TMB substrate solution into each well and incubate in the dark for 10-20 minutes at room temperature. Avoid direct exposure to light by wrapping the plate in aluminum foil.

Δ Note: Incubation time of the substrate solution is usually determined by the microplate reader performances: many microplate readers record absorbance only up to 2.0 O.D. The O.D. values of the plate should be monitored and the substrate reaction stopped before positive wells are no longer accurately readable (maximum ~20 minutes).

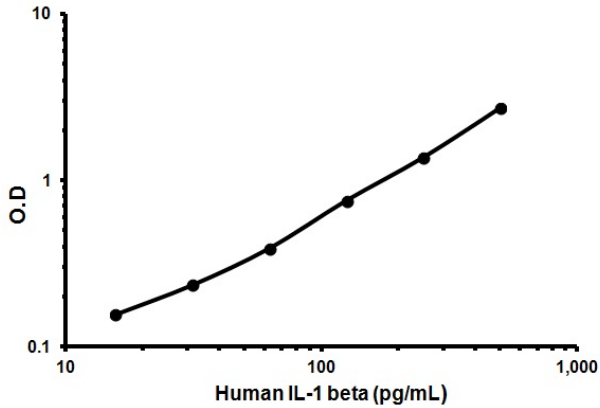
- 13.11** Add 100 μ L of Stop Reagent into each well. Results must be taken immediately after the addition of Stop Reagent, or within one hour, if the microplate is stored at 2-8°C in the dark.
- 13.12** Read absorbance of each well on a spectrophotometer using 450 nm as the primary wavelength and optionally 620 nm (610 nm to 650 nm is acceptable) as the reference wavelength.

14. Calculations

Calculate the mean absorbance for each set of duplicate standards, controls and samples, and subtract the average zero standard optical density. Plot the standard curve on log-log graph paper, with standard concentration on the x-axis and absorbance on the y-axis. Draw the best-fit straight line through the standard points.

15. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



| Conc. (pg/mL) | O.D. |
|---------------|-------|
| 500 | 2.708 |
| 250 | 1.368 |
| 125 | 0.754 |
| 62.5 | 0.390 |
| 31.25 | 0.235 |
| 15.6 | 0.156 |
| 0 | 0.069 |

Figure 1. Example of Human IL-1 beta standard curve.

16. Typical Sample Values

EXPECTED SERUM VALUES –

A panel of 20 Human sera and 20 Plasma samples were tested for IL-1 beta. All were evaluated below the detection level (6.5 pg/mL).

SENSITIVITY –

The sensitivity or minimum detectable dose of IL-1 beta using this IL-1 beta ELISA kit was found to be 6.5 pg/mL. This was determined by adding 2 standard deviations to the mean OD obtained when the zero standard was assayed in 6 independent experiments.

PRECISION –

| | Intra-assay Precision | Inter-Assay Precision |
|--------|-----------------------|-----------------------|
| n= | 6 | 6 |
| CV (%) | 4.5 | 8.7 |

DILUTION PARALLELISM –

Three spiked Human serum samples with different levels of IL-1 beta were analysed at different serial two fold dilutions (1:2 to 1:16) with two replicates each. Recoveries ranged from 83 to 111% with an overall **mean recovery of 96%**.

SPIKE RECOVERY -

The spike recovery was evaluated by spiking in Human serum and culture medium samples different concentrations of IL-1 beta in 3 separate experiments. Recoveries ranged from 72 to 105% with an overall **mean recovery of 89%**.

17. Assay Specificity

The assay recognizes both natural and recombinant Human IL-1 beta. To define the specificity of this ELISA several proteins were tested for cross reactivity. There was no cross reactivity observed for any protein tested (IL-1 α , IL-2, IL-10, IL-12, IL-17A, IL-23, IFN γ , Gp130, IL-33, TNF α).

Please contact our Technical Support team for more information.

18. Troubleshooting

| Problem | Reason | Solution |
|--------------------------|---|---|
| Poor standard curve | Inaccurate pipetting | Check pipettes |
| | Improper standards dilution | Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing |
| Low Signal | Incubation times too brief | Ensure sufficient incubation times; change to overnight standard/sample incubation |
| | Inadequate reagent volumes or improper dilution | Check pipettes and ensure correct preparation |
| Low Precision / Large CV | Use of expired components | Check the expiration date listed before use. Do not interchange components from different lots |
| | Improper wash step | Check that the correct wash buffer is being used. Check that all wells are empty after aspiration. Check that the microplate washer is dispensing properly. If washing by pipette, check for proper pipetting technique |
| | Contaminated wash buffer | Prepare fresh wash buffer |
| | Splashing of reagents while loading wells | Pipette properly in a controlled and careful manner |

| | | |
|--|---|--|
| <p>Low Precision / Large CV</p> | <p>Inconsistent volumes loaded into wells</p> | <p>Pipette properly in a controlled and careful manner. Check pipette calibration. Check pipette for proper performance</p> |
| | <p>Insufficient mixing of reagent dilutions</p> | <p>Thoroughly agitate the lyophilized components after reconstitution. Thoroughly mix dilutions</p> |
| | <p>Improperly sealed microplate</p> | <p>Check the microplate pouch for proper sealing. Check that the microplate pouch has no punctures. Check that three desiccants are inside the microplate pouch prior to sealing</p> |
| <p>Low sensitivity</p> | <p>Improper storage of the ELISA kit</p> | <p>Store the reconstituted protein at -80°C, all other assay components 4°C. Keep substrate solution protected from light.</p> |

19. Notes

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)