

Human IL-10 ELISA Kit High Sensitivity ab46059

[5 References](#) [1 Image](#)

Overview

Product name Human IL-10 ELISA Kit High Sensitivity

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|------|
| Overall | 3 | | | 7.8% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | 3 | | | 10.2% |

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 1.3 pg/ml

Range 1.56 pg/ml - 50 pg/ml

Recovery 87 %

Sample specific recovery

| Sample type | Average % | Range |
|--------------------------|-----------|-----------|
| Cell culture supernatant | 94 | % - % |
| Serum | 80 | 59% - 93% |

Assay time 3h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's IL-10 Human High Sensitivity *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 in Human sera, plasmas, buffered solutions and cell culture media.

A monoclonal antibody specific for IL-10 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-10 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-10 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-10 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-10.

Get results in 90 minutes with Human IL-10 ELISA Kit ([ab185986](#)) from our SimpleStep ELISA® range.

Platform Microplate

Properties

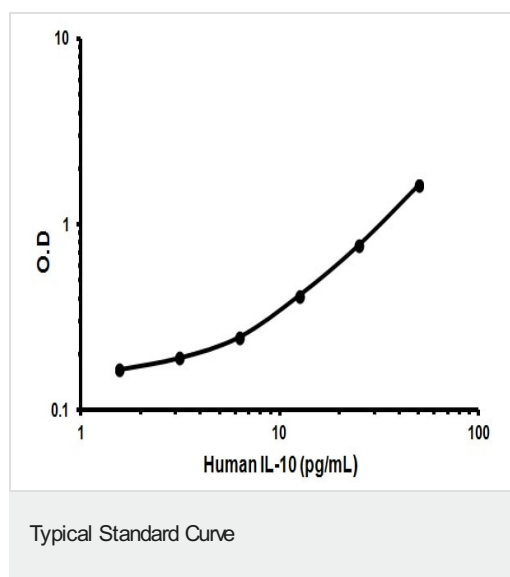
Storage instructions Store at +4°C. Please refer to protocols.

| Components | Identifier | 1 x 96 tests |
|---------------------------------------|-------------------|--------------|
| 10X Standard Diluent Buffer | Black | 1 x 15ml |
| 200X Wash Buffer | White | 1 x 10ml |
| Amplification Diluent | Brown & blue spot | 1 x 15ml |
| Amplifier | Yellow | 1 x 200µl |
| Biotinylated anti IL-10 | Red | 1 x 400µl |
| Biotinylated Antibody Diluent | Red | 1 x 7.5ml |
| Chromogen TMB Substrate Solution | | 1 x 11ml |
| HRP Diluent | Red | 1 x 23ml |
| IL-10 Microplate (12 x 8 well strips) | | 1 unit |
| IL-10 Standard (Lyophilized) | Yellow | 2 vials |
| Stop Reagent | Black | 1 x 11ml |

| Components | Identifier | 1 x 96 tests |
|------------------|------------|--------------|
| Streptavidin-HRP | | 2 x 5µl |

| | |
|------------------------------|---|
| Function | Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells. |
| Tissue specificity | Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types. |
| Sequence similarities | Belongs to the IL-10 family. |
| Cellular localization | Secreted. |

Images



Representative Standard Curve using ab46059

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors