

Human IL-10 ELISA Kit ab108871

1 Image

Overview

Product name Human IL-10 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				4.6%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				9.6%

Sample type Cell culture supernatant, Serum, Plasma, Tissue

Assay type Sandwich (quantitative)

Sensitivity > 60 pg/ml

Range 0.5 ng/ml - 4 ng/ml

Recovery 97 %

Assay time 5h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's IL-10 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 levels in cell culture supernatants, serum, plasma and tissue extracts.

An IL-10 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an IL-10 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of IL-10 captured in plate. This assay recognizes both natural and recombinant Human IL-10.

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

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Platform Microplate

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human IL-10 Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
IL-10 Microplate (12 x 8 well strips)	1 unit
IL-10 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

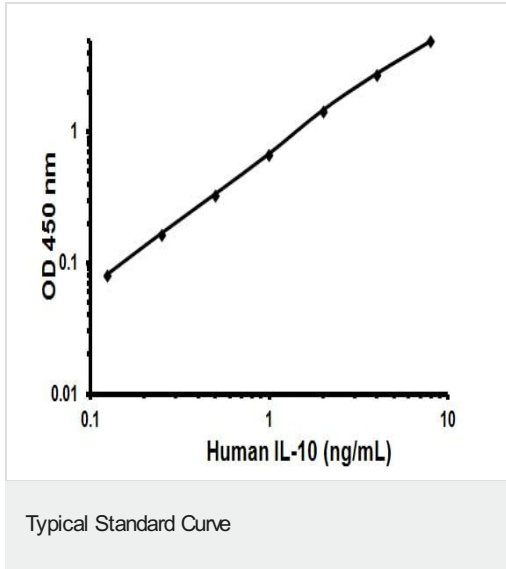
Function Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

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 ab108871.

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