

Mouse IL-10 ELISA Kit ab100697

★★★★★ [1 Abreviews](#) [19 References](#) [2 Images](#)

Overview

Product name	Mouse IL-10 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 45 pg/ml
Range	78.13 pg/ml - 5000 pg/ml
Recovery	93 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	103.7	79% - 120%
Serum	85.97	77% - 93%
Plasma	92.2	71% - 134%

Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Mouse

**Product overview**  
Abcam’s Mouse IL-10 ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of IL-10 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse IL-10 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IL-10 antibody is added. After washing away unbound biotinylated antibody, HRP conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IL-10 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform	Microplate
----------	------------

## Properties

### Storage instructions

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-Mouse IL-10	2 vials
IL-10 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-10 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### 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.

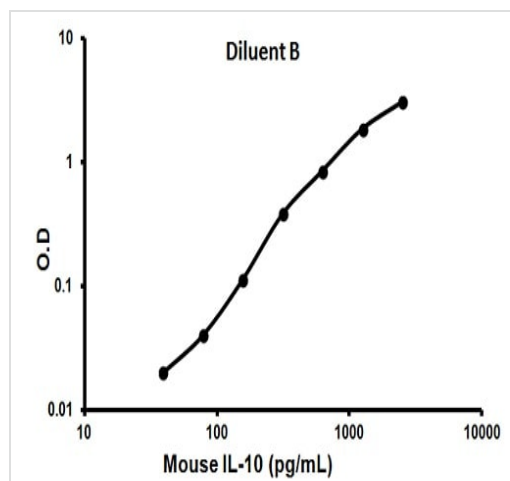
### Sequence similarities

Belongs to the IL-10 family.

### Cellular localization

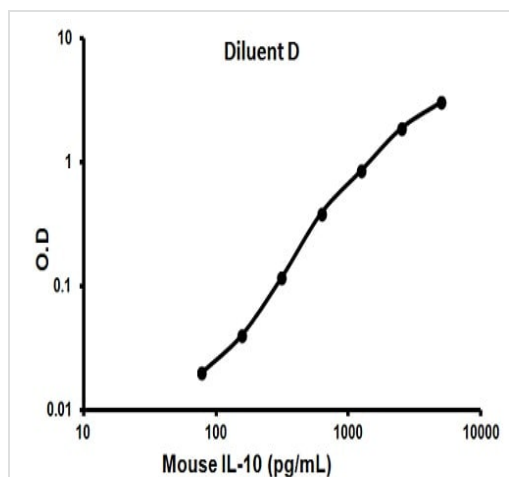
Secreted.

## Images



Representative standard curve using ab100697

Typical standard curve



Representative standard curve using ab100697

Typical standard curve

**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