

Human IL-10RB ELISA Kit ab100550

[2 Images](#)

Overview

Product name	Human IL-10RB ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 5 pg/ml
Range	4.12 pg/ml - 3000 pg/ml
Recovery	> 100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	129.1	120% - 140%
Serum	121.9	110% - 129%
Plasma	104.5	92% - 112%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

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

This assay employs an antibody specific for Human IL-10RB coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-10RB present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-10RB 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-10RB bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples

Platform Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-10RB	2 vials
IL-10RB Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-10RB Standard (lyophilized)	2 x 1ml
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

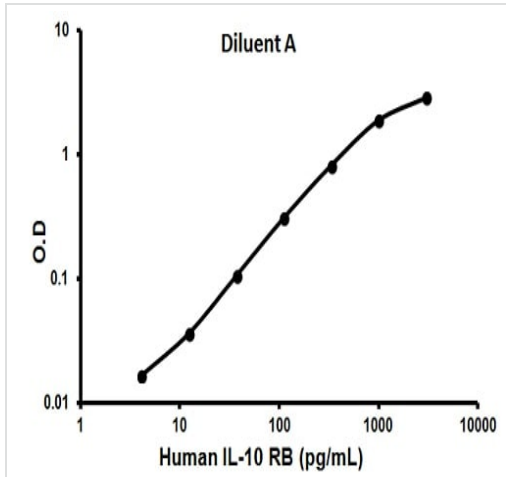
Function Shared cell surface receptor required for the activation of five class 2 cytokines: IL10, IL22, IL26, IL28, and IL29.

Involvement in disease Defects in IL10RB are the cause of inflammatory bowel disease type 25 (IBD25) [MIM:612567]. It is a chronic, relapsing inflammation of the gastrointestinal tract with a complex etiology. It is subdivided into Crohn disease and ulcerative colitis phenotypes. Crohn disease may affect any part of the gastrointestinal tract from the mouth to the anus, but most frequently it involves the terminal ileum and colon. Bowel inflammation is transmural and discontinuous; it may contain granulomas or be associated with intestinal or perianal fistulas. In contrast, in ulcerative colitis, the inflammation is continuous and limited to rectal and colonic mucosal layers; fistulas and granulomas are not observed. Both diseases include extraintestinal inflammation of the skin, eyes, or joints.

Sequence similarities Belongs to the type II cytokine receptor family.
Contains 2 fibronectin type-III domains.

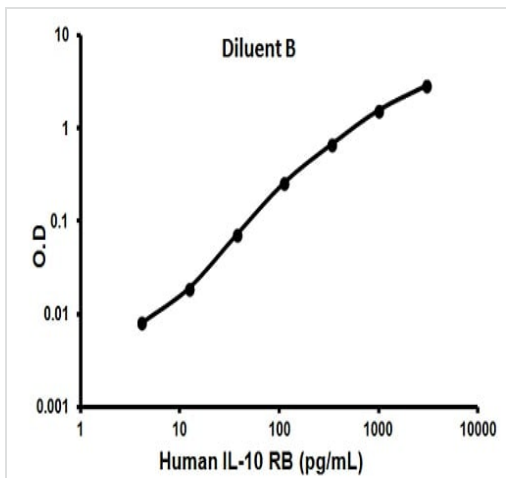
Cellular localization Membrane.

Images



Representative standard curve using ab100550

Typical standard curve



Representative standard curve using ab100550

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