

Human IL-17A Receptor ELISA Kit α b100558

2 Images

Overview

Product name	Human IL-17A Receptor ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	12.35 pg/ml - 3000 pg/ml
Recovery	81 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	81.54	70% - 92%
Serum	87.64	83% - 104%
Plasma	73.27	67% - 87%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	Abcam's IL-17A Receptor Human ELISA (Enzyme Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-17A in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-17A coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-17A present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-17A 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-17A 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

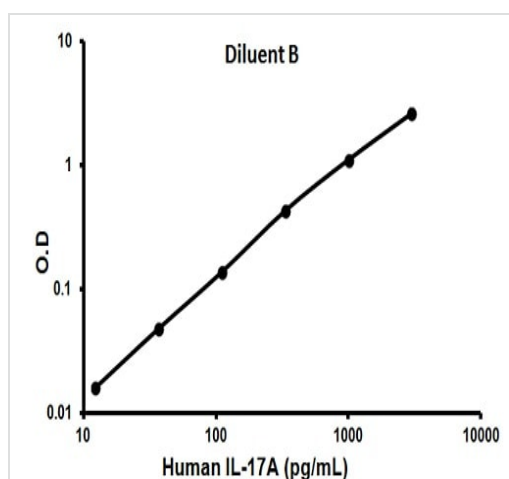
Properties

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Human IL-17A	2 vials
IL-17A Microplate (12 x 8 wells)	1 x 200µl
Recombinant Human IL-17A Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

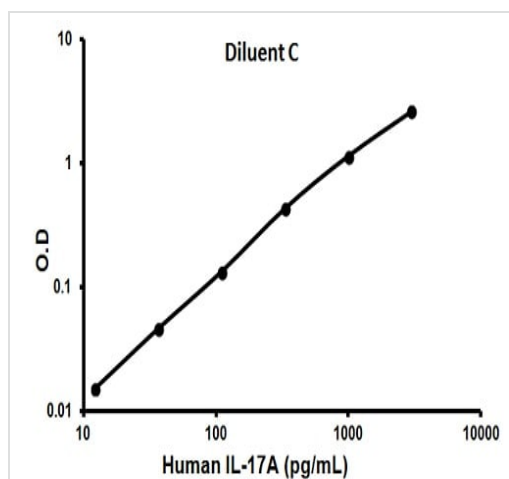
Function	Receptor for IL17A and IL17F. Binds its IL17A ligand with low affinity, suggesting that additional components are involved in IL17A-induced signaling.
Tissue specificity	Widely expressed.
Involvement in disease	Defects in IL17RA are the cause of familial candidiasis type 5 (CANDF5) [MIM:613953]. CANDF5 is a rare disorder with altered immune responses and impaired clearance of fungal infections, selective against Candida. It is characterized by persistent and/or recurrent infections of the skin, nails and mucous membranes caused by organisms of the genus Candida, mainly Candida albicans.
Sequence similarities	Contains 1 SEFIR domain.
Post-translational modifications	Glycosylated.
Cellular localization	Membrane.

Images



Representative standard curve using ab100558

Typical standard curve



Representative standard curve using ab100558

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