

Product datasheet

Human IL-17F ELISA Kit (Interleukin-17F) ab100557

[1 Image](#)

Overview

Product name	Human IL-17F ELISA Kit (Interleukin-17F)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	13.72 pg/ml - 10000 pg/ml
Recovery	93.5 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	90.63	81% - 99%
Serum	95.7	87% - 103%
Plasma	94.25	81% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's IL-17F (Interleukin 17F) Human ELISA (Enzyme- Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-17F in serum, plasma, and cell culture supernatants.

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

Tested applications **Suitable for:** Sandwich ELISA

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 Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Human IL-17F (lyophilized)	2 vials
IL-17F Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-17F Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function	Stimulates the production of other cytokines such as IL-6, IL-8 and granulocyte colony-stimulating factor, and can regulate cartilage matrix turnover. Stimulates PBMC and T-cell proliferation. Inhibits angiogenesis.
Tissue specificity	Expressed in activated, but not resting, CD4+ T-cells and activated monocytes.
Sequence similarities	Belongs to the IL-17 family.
Cellular localization	Secreted.

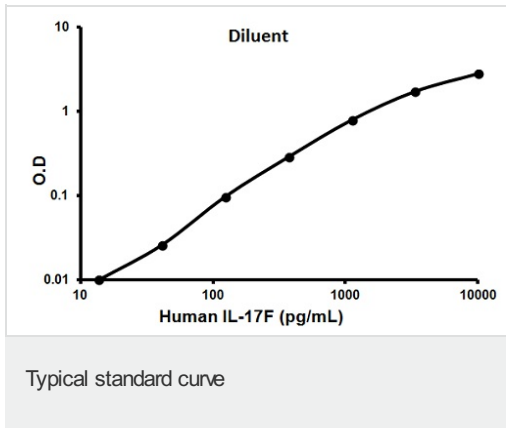
Applications

Our [Abpromise guarantee](#) covers the use of **ab100557** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative standard curve using
ab100557

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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