

### Human IL-17F ELISA Set (without plates) ab105233

[1 Image](#)

#### Overview

<b>Product name</b>	Human IL-17F ELISA Set (without plates)
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 6 pg/ml
<b>Range</b>	15.6 pg/ml - 500 pg/ml
<b>Assay time</b>	3h 45m
<b>Assay duration</b>	Multiple steps standard assay
<b>Product overview</b>	<p>Abcam's IL-17F Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-17F in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>ab105233 includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>ab105233 is intended to measure human IL-17F in cell culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>The assay recognizes natural and recombinant human IL-17A.</p> <p><b>RANGE:</b> 15.6pg/ml - 500pg/ml  <b>SENSITIVITY:</b> &lt; 6 pg/ml</p>
<b>Tested applications</b>	<b>Suitable for:</b> Sandwich ELISA
<b>Platform</b>	Reagents

#### Properties

<b>Storage instructions</b>	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	5 x 96 tests	10 x 96 tests
Detection biotinylated anti-IL-17F antibody	1 vial	2 vials
IL-17F Capture Antibody	1 x 0.5ml	2 x 0.5ml
IL-17F Standard	5 vials	10 vials
Ready-to-use TMB	2 vials	4 vials
Streptavidin-HRP	1 vial	2 vials

## Function

Ligand for IL17RA and IL17RC (PubMed:17911633). The heterodimer formed by IL17A and IL17F is a ligand for the heterodimeric complex formed by IL17RA and IL17RC (PubMed:18684971). Involved in stimulating the production of other cytokines such as IL6, IL8 and CSF2, and in regulation of cartilage matrix turnover (PubMed:11591732, PubMed:11591768, PubMed:11574464). Also involved in stimulating the proliferation of peripheral blood mononuclear cells and T-cells and in inhibition of angiogenesis (PubMed:11591732). Plays a role in the induction of neutrophilia in the lungs and in the exacerbation of antigen-induced pulmonary allergic inflammation.

## Tissue specificity

Expressed in activated, but not resting, CD4+ T-cells and activated monocytes.

## Involvement in disease

Candidiasis, familial, 6

## Sequence similarities

Belongs to the IL-17 family.

## Cellular localization

Secreted.

## Applications

### The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab105233 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
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

## Images

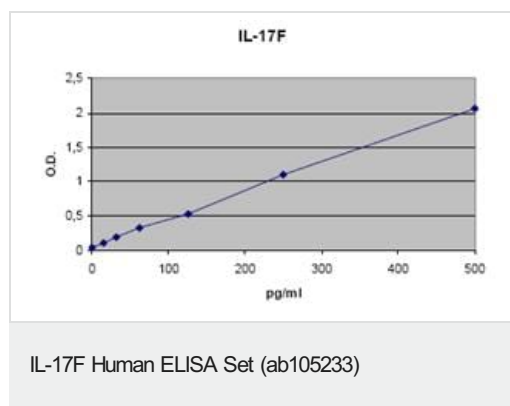


Figure 2: IL-17F Standard curve ranging from 15.6 to 500 pg/mL obtained using ab105233

**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