

## Product datasheet

# Human CXCL11 ELISA Kit (I-TAC) ab100580

★★★★★ 1 Abreviews 1 References 2 Images

### Overview

<b>Product name</b>	Human CXCL11 ELISA Kit (I-TAC)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 2 pg/ml
<b>Range</b>	2.5 pg/ml - 600 pg/ml
<b>Recovery</b>	93 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	92.82	80% - 105%
Serum	95.28	84% - 106%
Plasma	93.47	83% - 104%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

### Product overview

Abcam's Human CXCL11 ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CXCL11 in serum, plasma (heparin or EDTA), and cell culture supernatants.

This assay employs an antibody specific for Human CXCL11 coated on a 96-well plate. Standards and samples are pipetted into the wells and CXCL11 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CXCL11 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 CXCL11 C bound. The stop solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get higher sensitivity in only 90 minutes with Human CXCL11 ELISA Kit (I-TAC) ([ab187392](#)) from our SimpleStep ELISA® range.

**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 A	1 x 30ml
Biotinylated anti-Human CXCL11	2 vials
CXCL11 Microplate (12 strips x 8 wells)	1 unit
Recombinant Human CXCL11 Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Chemotactic for interleukin-activated T-cells but not unstimulated T-cells, neutrophils or monocytes. Induces calcium release in activated T-cells. Binds to CXCR3. May play an important role in CNS diseases which involve T-cell recruitment. May play a role in skin immune responses.

**Tissue specificity** High levels in peripheral blood leukocytes, pancreas and liver astrocytes. Moderate levels in thymus, spleen and lung. Low levels in placenta, prostate and small intestine. Also found in epidermal basal layer keratinocytes in skin disorders.

**Sequence similarities** Belongs to the intercrine alpha (chemokine CxC) family.

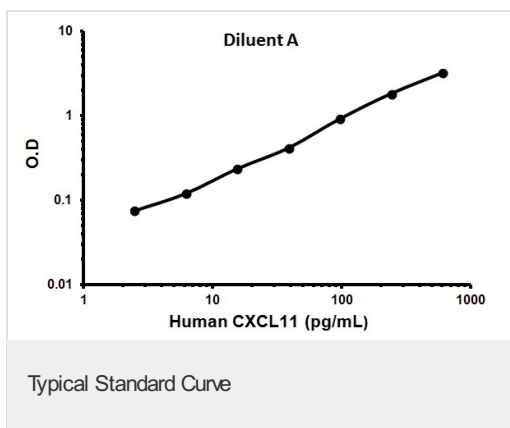
**Cellular localization** Secreted.

## Applications

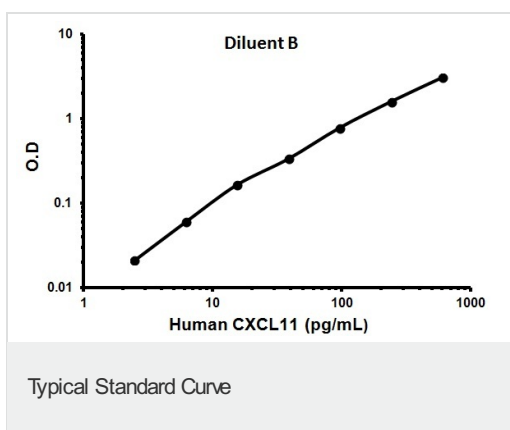
Our [Abpromise guarantee](#) covers the use of **ab100580** in the following tested applications.

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.



Representative Standard Curve using ab100580.



Representative Standard Curve using ab100580.

**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 <http://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