

## Product datasheet

# Human LEC ELISA Kit (CCL16) ab100532

★★★★★ 1 Abreviews 2 Images

### Overview

<b>Product name</b>	Human LEC ELISA Kit (CCL16)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 8 pg/ml
<b>Range</b>	8.23 pg/ml - 6000 pg/ml
<b>Recovery</b>	95 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	96.47	85% - 104%
Serum	95.29	84% - 103%
Plasma	94.68	83% - 103%

**Assay duration** Multiple steps standard assay

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

**Product overview** Abcam's LEC (CCL16) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human LEC in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human LEC coated on a 96-well plate. Standards and samples are pipetted into the wells and LEC present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human LEC 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 LEC 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
20X Wash Buffer Concentrate	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 LEC	2 vials
LEC Microplate (12 x 8 wells)	1 unit
Recombinant Human LEC Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Shows chemotactic activity for lymphocytes and monocytes but not neutrophils. Also shows potent myelosuppressive activity, suppresses proliferation of myeloid progenitor cells. Recombinant SCYA16 shows chemotactic activity for monocytes and THP-1 monocytes, but not for resting lymphocytes and neutrophils. Induces a calcium flux in THP-1 cells that were desensitized by prior expression to RANTES.

**Tissue specificity** Mainly expressed in liver, also found in spleen and thymus. Highly expressed in LPS- and IFN-gamma-activated monocytes, weakly in some lymphocytes, including natural killer cells, gamma-delta T-cells, and some T-cell clones.

**Sequence similarities** Belongs to the intercrine beta (chemokine CC) family.

**Cellular localization** Secreted.

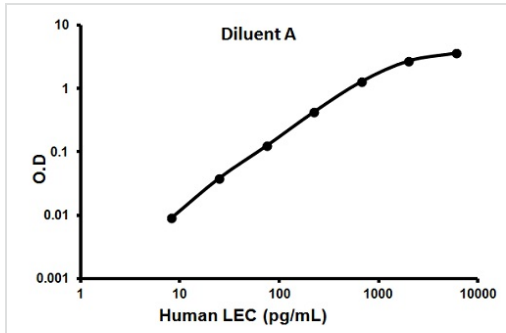
## Applications

Our [Abpromise guarantee](#) covers the use of **ab100532** in the following tested applications.

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

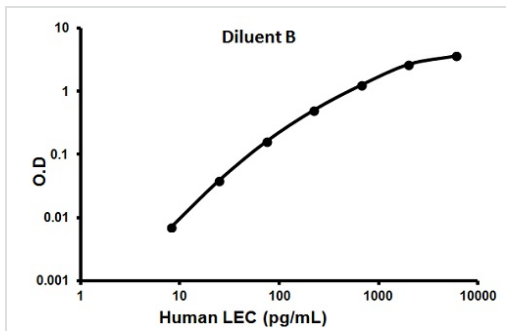
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



Representative standard curve using ab100532

Typical standard curve



Representative standard curve using ab100532

Typical standard curve

**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