

Human CCL23 ELISA Kit **ab100611**

[2 References](#) [2 Images](#)

Overview

Product name	Human CCL23 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 7 pg/ml
Range	8.23 pg/ml - 6000 pg/ml
Recovery	93 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	93.48	81% - 105%
Serum	92.78	82% - 105%
Plasma	94.16	83% - 106%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's CCL23 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CCL23 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human CCL23 coated on a 96-well plate. Standards and samples are pipetted into the wells and CCL23 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CCL23 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 CCL23 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 CCL23 ELISA Kit (**ab216169**) from our SimpleStep ELISA® range.

Notes Optimisation 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 A	1 x 30ml
Biotinylated anti-Human CCL23	2 vials
CCL23 Microplate (12 x 8 wells)	1 unit
Recombinant Human CCL23 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

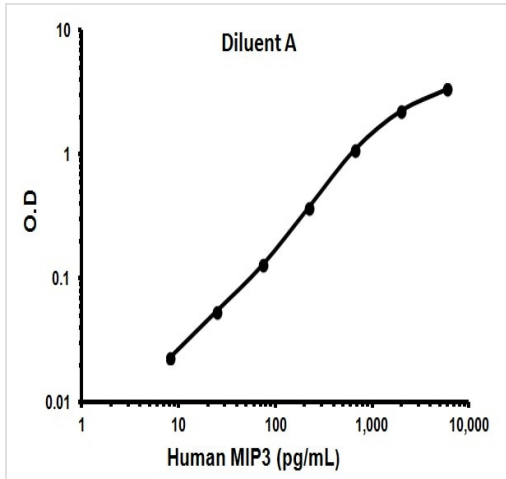
Function Shows chemotactic activity for monocytes, resting T-lymphocytes, and neutrophils, but not for activated lymphocytes. Inhibits proliferation of myeloid progenitor cells in colony formation assays. This protein can bind heparin. Binds CCR1. CCL23(19-99), CCL23(22-99), CCL23(27-99), CCL23(30-99) are more potent chemoattractants than the small-inducible cytokine A23.

Tissue specificity High levels in adult lung, liver, skeletal muscle and pancreas. Moderate levels in fetal liver, adult bone marrow and placenta. The short form is the major species and the longer form was detected only in very low abundance. CCL23(19-99), CCL23(22-99), CCL23(27-99), CCL23(30-99) are found in high levels in synovial fluids from rheumatoid patients.

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

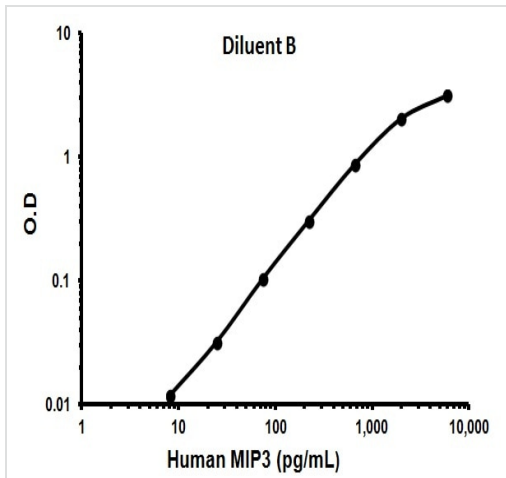
Cellular localization Secreted.

Images



Representative Standard Curve using ab100611.

Typical Standard Curve



Representative Standard Curve using ab100611.

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