

Rat CCL3 ELISA Kit ab242242

Recombinant SimpleStep ELISA

[5 Images](#)

Overview

Product name Rat CCL3 ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
supernatant	3			3.5%

Inter-assay

Sample	n	Mean	SD	CV%
supernatant	8			3.5%

Sample type Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 1.96 pg/ml

Range 7.81 pg/ml - 500 pg/ml

Recovery Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	91	90% - 93%
Serum	84	82% - 85%
Hep Plasma	85	81% - 89%
EDTA Plasma	75	72% - 78%
Cit plasma	78	77% - 78%

Assay time 1h 30m

Assay duration One step assay

Species reactivity**Reacts with:** Rat**Product overview**

Rat CCL3 ELISA Kit (ab242242) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CCL3 protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Rat CCL3 with 1.96 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties**Storage instructions**

Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Rat CCL3 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Plate Seals	1 unit
Rat CCL3 Capture Antibody (Lyophilized)	1 vial
Rat CCL3 Lyophilized Recombinant Protein	2 vials
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function

Monokine with inflammatory and chemokinetic properties. Binds to CCR1, CCR4 and CCR5. One of the major HIV-suppressive factors produced by CD8+ T-cells. Recombinant MIP-1-alpha

induces a dose-dependent inhibition of different strains of HIV-1, HIV-2, and simian immunodeficiency virus (SIV).

Sequence similarities

Belongs to the intercrine beta (chemokine CC) family.

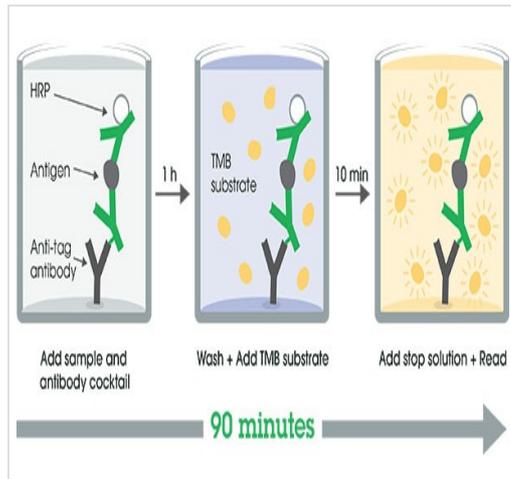
Post-translational modifications

N-terminal processed form LD78-alpha(4-69) is produced by proteolytic cleavage after secretion from HTLV1-transformed T-cells.

Cellular localization

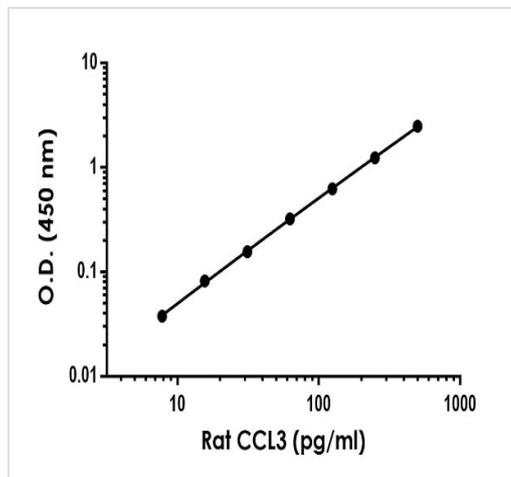
Secreted.

Images



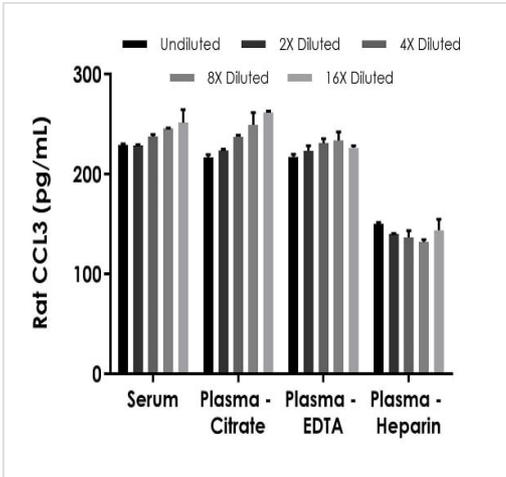
SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Other - Rat CCL3 ELISA Kit (ab242242)



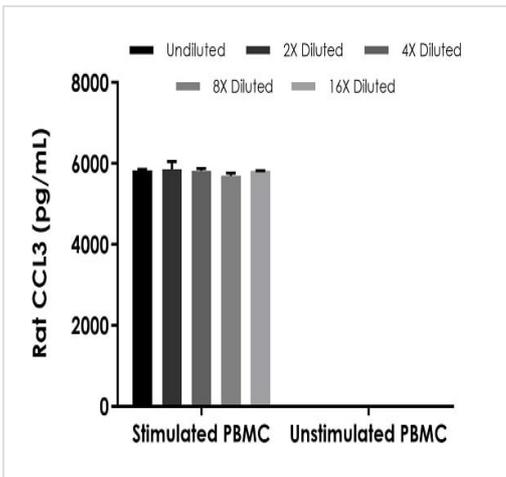
The CCL3 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of rat CCL3 standard curve in Sample Diluent NS.



Interpolated concentrations of spiked CCL3 in rat serum and plasma samples.

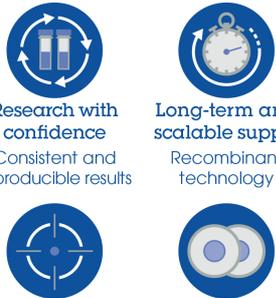
The concentrations of CCL3 were measured in duplicates, interpolated from the CCL3 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 25%, plasma (EDTA) 50% and plasma (heparin) 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).



Interpolated concentrations of native CCL3 in stimulated and unstimulated rat PBMC supernatant samples.

The concentrations of CCL3 were measured in duplicates, interpolated from the CCL3 standard curves and corrected for sample dilution. Undiluted samples are as follows: stimulated rat PBMC supernatant 5% and unstimulated rat PBMC supernatant 5%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CCL3 concentration was determined to be 5,805 pg/mL in stimulated rat PBMC supernatant and was undetectable in unstimulated rat PBMC supernatant.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results

Long-term and scalable supply
Recombinant technology

Success from the first experiment
Confirmed specificity

Ethical standards compliant
Animal-free production

Sandwich ELISA - Rat CCL3 ELISA Kit (ab242242)

To learn more about the advantages of recombinant antibodies see [here](#).

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