

Mouse CCL6 ELISA Kit ab277715

Recombinant SimpleStep ELISA

[7 Images](#)

Overview

Product name Mouse CCL6 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum	8			3.2%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			5.8%

Sample type

Cell culture supernatant, Serum, Cell culture extracts, Tissue Extracts, Hep Plasma, EDTA Plasma, Cit plasma

Assay type

Sandwich (quantitative)

Sensitivity

0.97 pg/ml

Range

6.25 pg/ml - 400 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	110	104% - 116%
Serum	115	112% - 116%
Tissue Extracts	94	92% - 97%
Hep Plasma	114	110% - 117%
EDTA Plasma	112	106% - 118%
Cit plasma	110	106% - 115%

Assay time 1h 30m

Assay duration One step assay

Species reactivity **Reacts with:** Mouse

Product overview Mouse CCL6 ELISA kit (ab277715) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Mouse CCL6 protein in mouse serum, plasma, cell culture supernatant, and cell and tissue extract samples. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse CCL6 with .97 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.

Notes CCL6, identified only in rodents, is a small cytokine that belongs to the CC chemokine family in which plays important roles in the immune system such as wound healing, cellular differentiation, macrophage chemoattractant, angiogenesis, and inflammation. Studies have shown CCL6 to promote innate immunity through cell activation and recruitment as well potentially having a role in tumorigenic and metastatic behavior.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

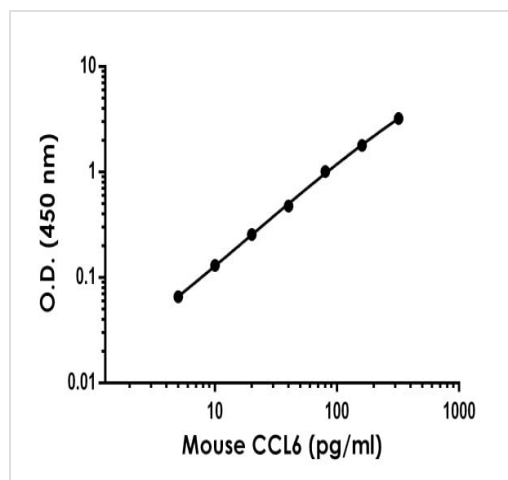
Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Mouse CCL6 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml

Components	1 x 96 tests
Antibody Diluent 4BR	1 x 6ml
Mouse CCL6 Capture Antibody (Lyophilized)	1 x 600µl
Mouse CCL6 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
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

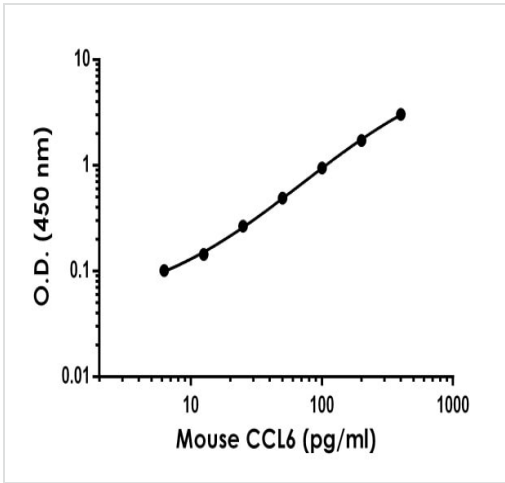
Function	CCL6(22-95) and CCL6(23-95) are potent chemoattractants.
Tissue specificity	Expressed in myelopoietic bone marrow cultures stimulated by GM-CSF.
Sequence similarities	Belongs to the intercrine beta (chemokine CC) family.
Cellular localization	Secreted.

Images



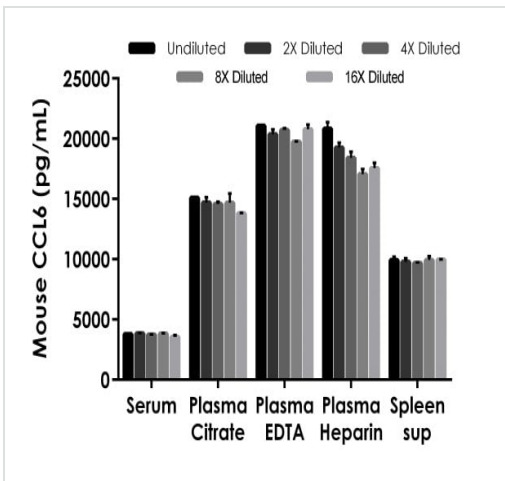
Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse CCL6 standard curve in Sample Diluent NS. The CCL6 standard curve was prepared as described in Section 10.



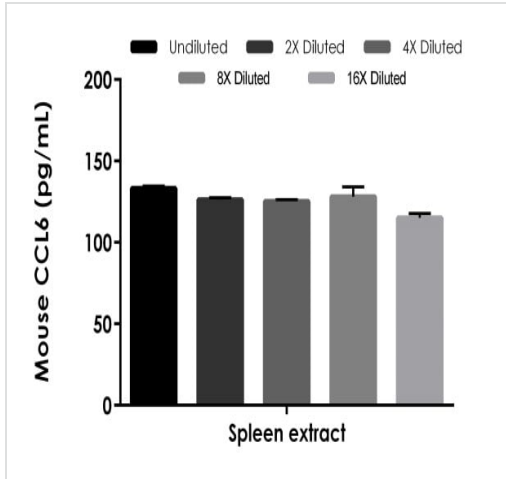
Example of mouse CCL6 standard curve in 1X Cell Extraction Buffer PTR.

The CCL6 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.



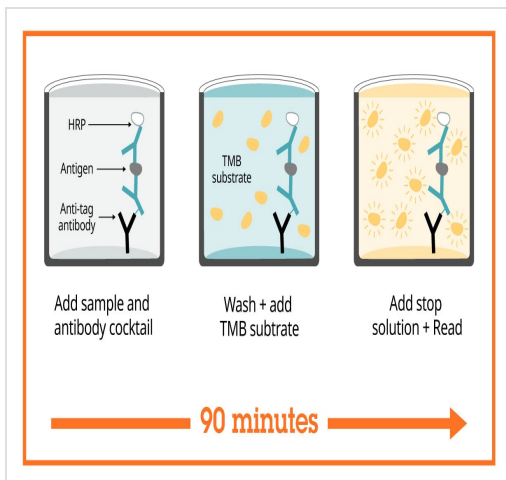
Interpolated concentrations of native or spike CCL6 in mouse serum, plasma and cell culture supernatant samples.

The concentrations of CCL6 were measured in duplicates, interpolated from the CCL6 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1:20, plasma (citrate) 1:80, plasma (EDTA) 1:80, plasma (heparin) 1:80, and spleen supernatant 1:40. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CCL6 concentration was determined to be 3,747 pg/mL in serum, 14,568 pg/mL in plasma (citrate), 20,533 pg/mL in plasma (EDTA), 18,616 pg/mL in plasma (heparin), and 9,852 pg/mL in spleen supernatant.



Interpolated concentrations of native CCL6 in mouse spleen extract sample based on a 400 µg/mL extract load.

The concentrations of CCL6 were measured in duplicate and interpolated from the CCL6 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CCL6 concentration was determined to be 125.6 pg/mL in spleen extract.



Sandwich ELISA - Mouse CCL6 ELISA Kit (ab277715)

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.

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

Recombinant Antibody Benefits

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

Get more done with
SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Mouse CCL6 ELISA Kit
(ab277715)

To learn more about the advantages of SimpleStep ELISA[®] kits 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