

## Product datasheet

# Mouse Fractalkine ELISA Kit (CX3CL1) ab100683

[2 Images](#)

### Overview

<b>Product name</b>	Mouse Fractalkine ELISA Kit (CX3CL1)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 0.3 ng/ml
<b>Range</b>	0.274 ng/ml - 200 ng/ml
<b>Recovery</b>	78 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	74.84	67% - 81%
Serum	74.56	67% - 82%
Plasma	84.97	68% - 95%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

### Product overview

Abcam's CX3CL1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse CX3CL1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse CX3CL1 coated on a 96- well plate. Standards and samples are pipetted into the wells and CX3CL1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse CX3CL1 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 CX3CL1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Tested applications** **Suitable for:** Sandwich ELISA

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
160X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse CX3CL1 (lyophilized)	2 vials
CX3CL1 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse CX3CL1 Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## Relevance

CX3CL1 is a member of the delta chemokine subfamily that contains a unique CX3C cysteine motif near the N-terminal. Unlike other known chemokines, it is a type 1 membrane protein containing a chemokine domain tethered on a long mucin-like stalk. CX3CL1, a leukocyte chemoattractant, is expressed in various tissues including the brain, heart, lung, kidney, skeletal muscle, and testis. The expression is reported to be up-regulated in endothelial cells and microglia by inflammatory signals. CX3CR1, a specific receptor for fractalkine, mediates both leukocyte migration and adhesion.

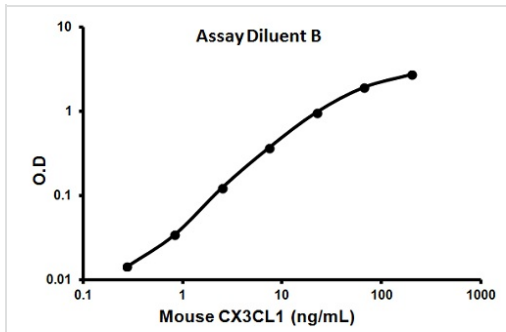
## Applications

Our [Abpromise guarantee](#) covers the use of **ab100683** in the following tested applications.

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

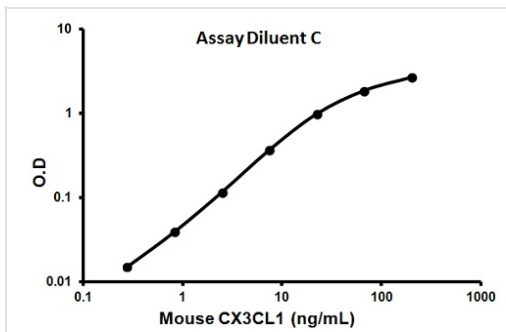
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



Representative standard curve using  
ab100683

Typical Standard Curve



Representative standard curve using  
ab100683

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 <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