

Mouse Fractalkine ELISA Kit (CX3CL1) ab100683

7 References 2 Images

Overview

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

Notes The standard is affinity purified.

Platform Microplate

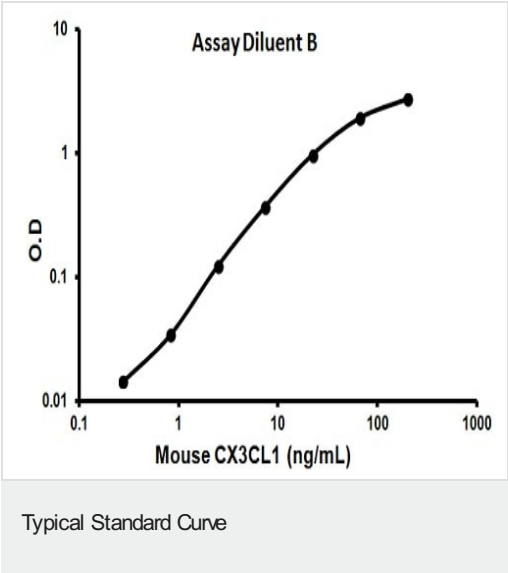
Properties

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

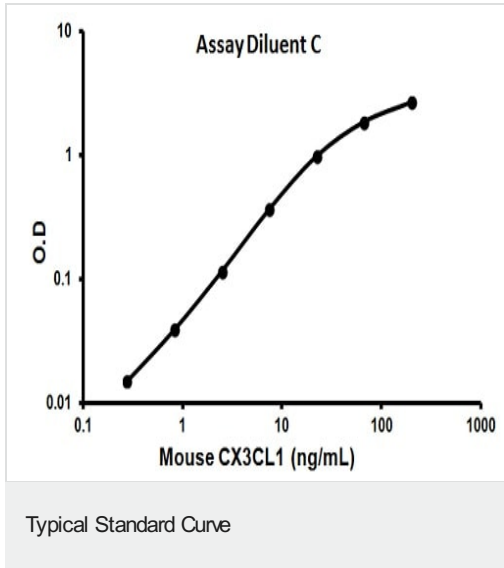
Components	1 x 96 tests
160X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse CX3CL1	2 vials
CX3CL1 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse CX3CL1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml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

Images



Representative standard curve using ab100683



Representative standard curve using ab100683

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