

Mouse CXCL5 ELISA Kit (GCP-2) ab100719

4 References 1 Image

Overview

Product name	Mouse CXCL5 ELISA Kit (GCP-2)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 35 pg/ml
Range	31.25 pg/ml - 2000 pg/ml
Recovery	91 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	83.94	74% - 94%
Serum	98.05	86% - 106%
Plasma	93.11	83% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's CXCL5 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse CXCL5 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for CXCL5 coated on a 96-well plate. Standards and samples are pipetted into the wells and CXCL5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse CXCL5 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 CXCL5 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

Storage instructions

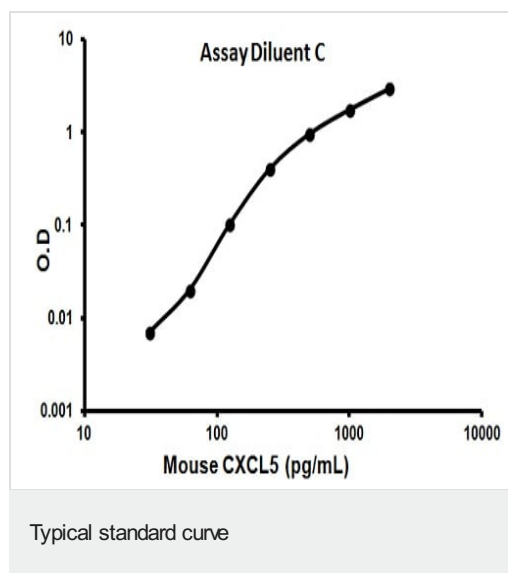
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse CXCL5	2 vials
CXCL5 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse CXCL5 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

CXCL5 is a 9.9kD inflammatory chemokine that belongs to the CXC chemokine family. CXCL5 is expressed by fibroblasts, is induced by bacterial lipopolysaccharides, and is chemotactic for neutrophils. This chemokine is produced concomitantly with Interleukin 8 (IL8) in response to stimulation with either Interleukin 1 (IL1) or Tumor Necrosis Factor alpha (TNFA). This chemokine is a potent chemotaxin involved in neutrophil activation.

Images



Representative standard curve using ab100719

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