

Mouse CXCL1 ELISA Kit (GRO alpha) ab100717

1 Image

Overview

Product name	Mouse CXCL1 ELISA Kit (GRO alpha)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1 pg/ml
Range	0.614 pg/ml - 150 pg/ml
Recovery	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	123.5	115% - 139%
Serum	89.3	78% - 97%
Plasma	101.4	88% - 113%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse
Product overview	Abcam's Mouse CXCL1 ELISA (Enzyme-Linked Immunosorbent Assay) kit (GRO alpha) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse CXCL1 (GRO alpha) in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse CXCL1 coated on a 96-well plate. Standards and samples are pipetted into the wells and CXCL1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse CXCL1 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 CXCL1 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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated anti-mouse GRO alpha	2 vials
GRO alpha Microplate (12 x 8 wells)	1 unit
Recombinant mouse GRO alpha Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function

Has chemotactic activity for neutrophils. May play a role in inflammation and exerts its effects on endothelial cells in an autocrine fashion. In vitro, the processed forms GRO-alpha(4-73), GRO-alpha(5-73) and GRO-alpha(6-73) show a 30-fold higher chemotactic activity.

Sequence similarities

Belongs to the intercrine alpha (chemokine CxC) family.

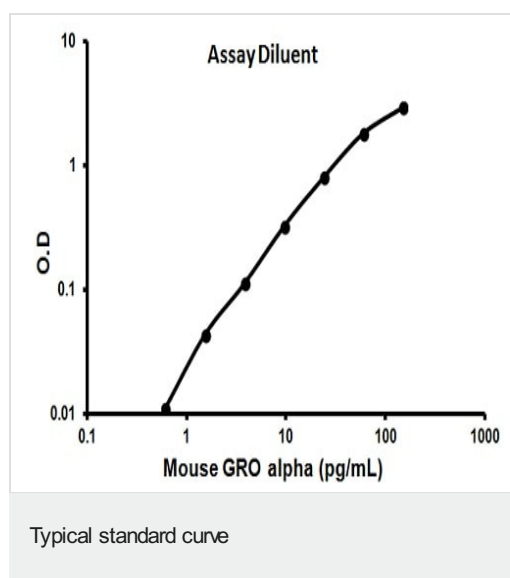
Post-translational modifications

N-terminal processed forms GRO-alpha(4-73), GRO-alpha(5-73) and GRO-alpha(6-73) are produced by proteolytic cleavage after secretion from peripheral blood monocytes.

Cellular localization

Secreted.

Images



Representative standard curve using ab100717

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