

Mouse Eotaxin-1 ELISA Kit ab100680

3 References 2 Images

Overview

Product name	Mouse Eotaxin-1 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 1 pg/ml				
Range	1.02 pg/ml - 100 pg/ml				
Recovery	94 %				
	Sample specific recovery				
	Sample type		Average %		Range
	Cell culture supernatant		95.25		83% - 104%
	Serum		94.95		83% - 104%
	Plasma		92.54		84% - 103%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's Eotaxin-1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse Eotaxin in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for mouse Eotaxin-1 coated on a 96-well plate. Standards and samples are pipetted into the wells and Eotaxin-1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse Eotaxin-1 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 Eotaxin-1 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
5X Assay Diluent B	1 x 15ml
80x HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse Eotaxin-1	2 vials
Eotaxin-1 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse Eotaxin-1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

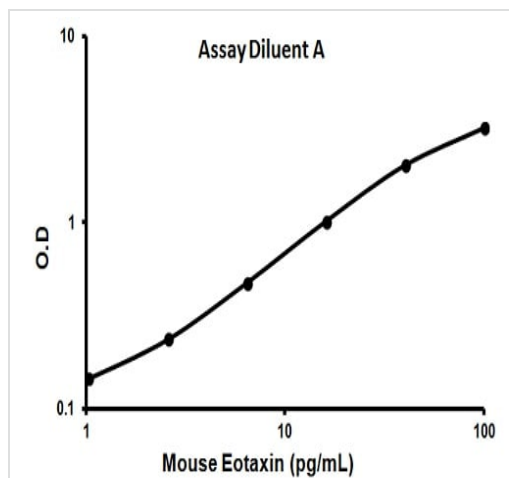
Function In response to the presence of allergens, this protein directly promotes the accumulation of eosinophils, a prominent feature of allergic inflammatory reactions. Binds to CCR3.

Sequence similarities Belongs to the intercrine beta (chemokine CC) family.

Post-translational modifications O-linked glycan consists of a Gal-GalNAc disaccharide which is modified with up to 2 sialic acid residues.

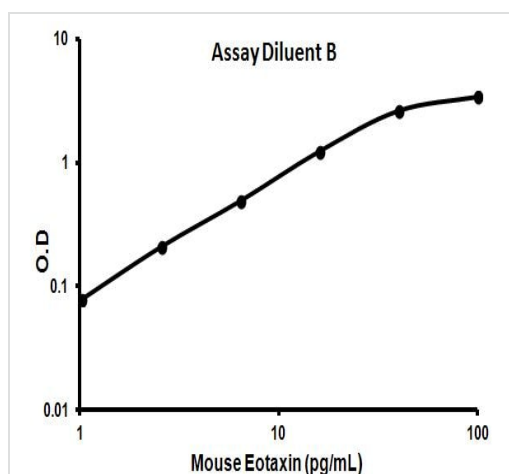
Cellular localization Secreted.

Images



Representative standard curve using ab100680

Typical Standard Curve



Representative standard curve using ab100680

Typical Standard Curve

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