

## Product datasheet

# Mouse Eotaxin-2 ELISA Kit (CCL24) ab100681

[1 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Mouse Eotaxin-2 ELISA Kit (CCL24)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 4 pg/ml
<b>Range</b>	3.28 pg/ml - 800 pg/ml
<b>Recovery</b>	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	96.65	89% - 102%
Serum	103.1	93% - 112%
Plasma	113.1	79% - 137%

**Assay duration** Multiple steps standard assay

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

**Product overview** Abcam's Eotaxin-2 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse Eotaxin-2 in serum, plasma and cell culture supernatants.

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

### Relevance

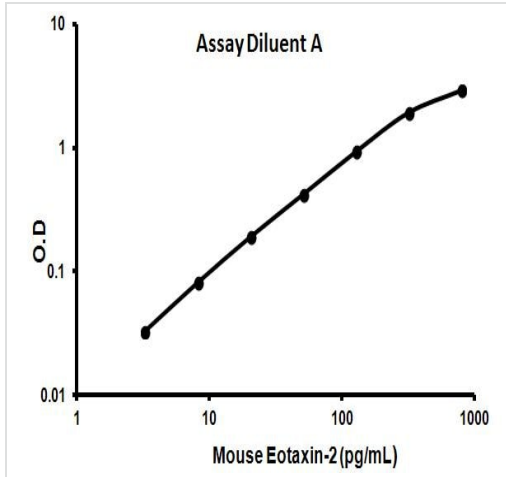
Eotaxin 2 is a member of the CC chemokine family, based on the presence of the CC motif and homology with other known CC chemokines. Eotaxin 2 cDNA encodes a 119 amino acid residue precursor protein with a 26 amino acid residue signal peptide that is cleaved to generate a mature protein predicted to contain 93 amino acid residues with an N glycosylation site. Mature human eotaxin 2 has a predicted molecular mass of approximately 10.6 kDa. Compared to other CC chemokines, eotaxin 2 exhibits 40 %, 42 %, and 39 % amino acid identity to MCP3, MIP1 alpha, and eotaxin, respectively. Human CC chemokine eotaxin 2 maps to chromosome 7q11.23. Both eotaxin and eotaxin 2 activate and attract eosinophils and basophils. A receptor for human eotaxin has been identified and found to be the third numbered receptor in the C-C chemokine subfamily of receptors (CCR3. On eosinophils, the effects of eotaxin 2 is inhibited by an CCR3 antibody and cross-desensitized by eotaxin and MCP4, suggesting that all three CC chemokines act through CCR3. Eotaxin 2 mRNA is weakly expressed in activated monocytes and T lymphocytes. Recombinant eotaxin 2 induces chemotaxis of eosinophils, basophils, and resting T lymphocytes but not monocytes and activated T lymphocytes. Eotaxin 2 inhibits colony formation in myeloid progenitor cells.

### Cellular localization

Secreted

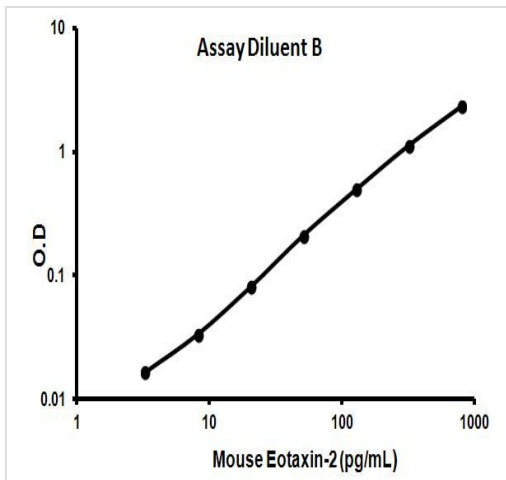
### Images

---



Representative Standard Curve using ab100681

Typical Standard Curve



Representative Standard Curve using ab100681

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