

Human Eotaxin 2 ELISA Kit (CCL24) ab100509

1 References 画像数 2

製品の概要

製品名	Human Eotaxin 2 ELISA Kit (CCL24)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 2 pg/ml
検出範囲	1.37 pg/ml - 1000 pg/ml
添加回収試験	95 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	95.77	82% - 103%
Serum	94.86	83% - 103%
Plasma	96.74	85% - 104%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	Abcam's Eotaxin 2 (CCL24) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human Eotaxin 2 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human 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-Human 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.

Get results in 90 minutes with Human Eotaxin-2 ELISA Kit (CCL24) ([ab179882](#)) from our

	SimpleStep ELISA [®] range.
特記事項	Optimization may be required with urine samples.
試験プラットフォーム	Microplate

製品の特性

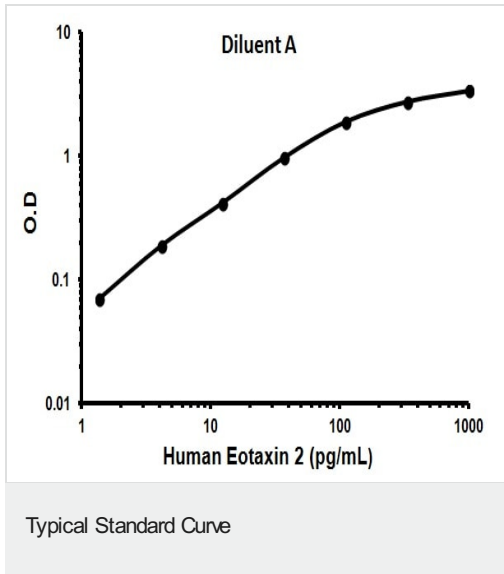
保存方法 Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Eotaxin 2	2 vials
Eotaxin 2 Microplate (12 x 8 well strips)	1 unit
Recombinant Human Eotaxin 2 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

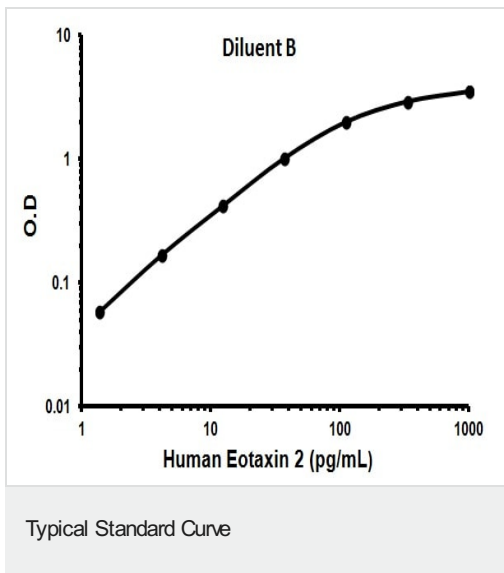
関連性 Eotaxin 2 is a member of the CC chemokine family, based on the presence of the CC motif and homology with other known CC chemokines. Eotaxin 2 cDNA encodes a 119 amino acid residue precursor protein with a 26 amino acid residue signal peptide that is cleaved to generate a mature protein predicted 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.

細胞内局在 Secreted

画像



Representative Standard Curve using ab100509.



Representative Standard Curve using ab100509.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors