

Rat Fractalkine ELISA Kit (CX3CL1) ab100761

1 References 画像数 1

製品の概要

| | |
|---------|--|
| 製品名 | Rat Fractalkine ELISA Kit (CX3CL1) |
| 検出方法 | Colorimetric |
| サンプルの種類 | Cell culture extracts, Tissue Extracts |
| アッセイタイプ | Sandwich (quantitative) |
| 検出感度 | < 5 pg/ml |
| 検出範囲 | 2.74 pg/ml - 2000 pg/ml |
| 添加回収試験 | 93 % |

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 extracts | 93.54 | 84% - 104% |
| Tissue Extracts | 92.87 | 83% - 103% |

| | |
|------|-------------------------------|
| ステップ | Multiple steps standard assay |
| 種交差性 | 交差種: Rat |

製品の概要
Abcam's Fractalkine Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat Fractalkine in cell lysates and tissue lysates.

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

特記事項
The standard is affinity purified.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.
It is the responsibility of our customers to check the necessity of application of REACH

Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム

Microplate

製品の特性

保存方法

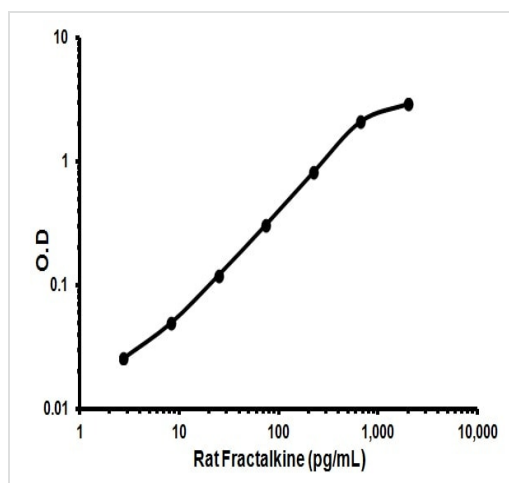
Store at -20°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 240X HRP-Streptavidin Concentrate | 1 x 200µl |
| 2X Cell Lysis Buffer | 1 x 5ml |
| 5X Assay Diluent | 1 x 15ml |
| 5X Sample Diluent Buffer | 1 x 10ml |
| Biotinylated anti-Rat Fractalkine | 2 vials |
| Fractalkine Microplate (12 x 8 wells) | 1 unit |
| Recombinant Rat Fractalkine Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

関連性

CX3CL1 is a member of the delta chemokine subfamily that contains a unique CX3C cysteine motif near the N-terminal. Unlike other known chemokines, it is a type 1 membrane protein containing a chemokine domain tethered on a long mucin-like stalk. CX3CL1, a leukocyte chemoattractant, is expressed in various tissues including the brain, heart, lung, kidney, skeletal muscle, and testis. The expression is reported to be up-regulated in endothelial cells and microglia by inflammatory signals. CX3CR1, a specific receptor for fractalkine, mediates both leukocyte migration and adhesion.

画像



Representative Standard Curve using ab100761

Representative 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.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