

Product datasheet

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.

アプリケーション 適用あり: Sandwich ELISA

試験プラットフォーム Microplate

法規制情報

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 Concentrate	1 x 25ml
240X HRP-Streptavidin Concentrate	1 x 200µl
2x Cell lysate 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	2 vials
Stop Solution	1 x 8ml
TMB 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.

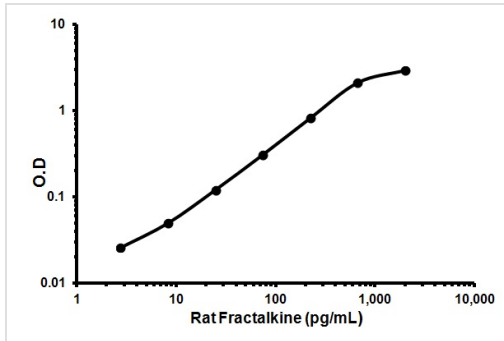
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab100761** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



Representative Standard Curve using
ab100761

Representative Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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