

Rat Fractalkine ELISA Kit (CX3CL1) ab100761

1 References 1 图像

概述

产品名称	大鼠Fractalkine ELISA试剂盒(CX3CL1)
检测方法	Colorimetric
样品类型	Cell culture extracts, Tissue Extracts
检测类型	Sandwich (quantitative)
灵敏度	< 5 pg/ml
范围	2.74 pg/ml - 2000 pg/ml
回收率	93 %

特定样本回收率

样品类型	平均%	范围
Cell culture extracts	93.54	84% - 104%
Tissue Extracts	92.87	83% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Rat
产品概述	Abcam's Fractalkine Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> 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.

说明 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.

平台

Microplate

性能

存放说明

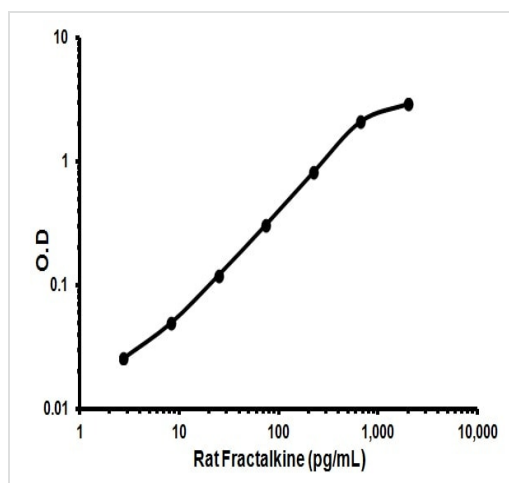
Store at -20°C. Please refer to protocols.

组件	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

相关性

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.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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