

Rat CXCL5 ELISA Kit ab100774

2 图像

概述

产品名称	大鼠CXCL5 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 20 pg/ml
范围	24.69 pg/ml - 6000 pg/ml
回收率	90 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	93.53	85% - 103%
Serum	88.63	83% - 102%
Plasma	89.45	84% - 103%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 Abcam's CXCL5 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of rat CXCL5 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for rat CXCL5 coated on a 96-well plate. Standards and samples are pipetted into the wells and CXCL5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-rat CXCL5 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 CXCL5 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

平台 Microplate

性能

存放说明

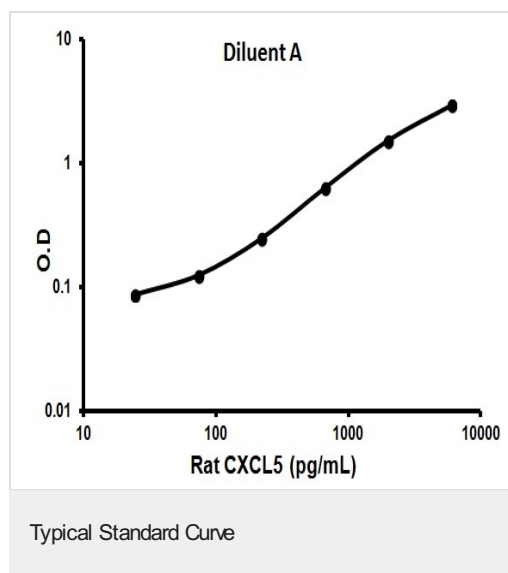
Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Rat CXCL5	2 vials
CXCL5 Microplate (12 x 8 wells)	1 unit
Recombinant Rat CXCL5 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

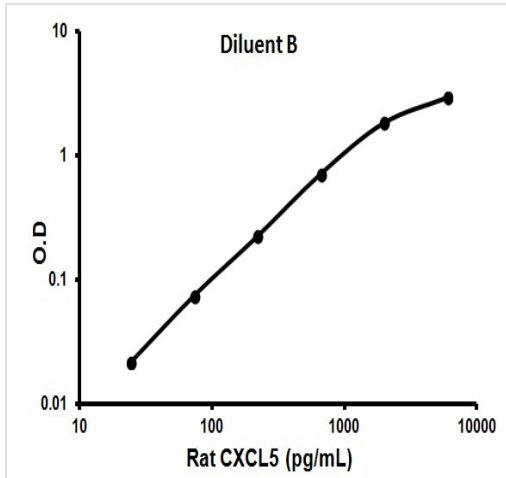
相关性

CXCL5 is a 9.9kD inflammatory chemokine that belongs to the CXC chemokine family. CXCL5 is expressed by fibroblasts, is induced by bacterial lipopolysaccharides, and is chemotactic for neutrophils. This chemokine is produced concomitantly with Interleukin 8 (IL8) in response to stimulation with either Interleukin 1 (IL1) or Tumor Necrosis Factor alpha (TNFA). This chemokine is a potent chemotaxin involved in neutrophil activation.

图片



Representative Standard Curve using ab100774



Representative Standard Curve using ab100774

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.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