

Mouse CXCL16 ELISA Kit ab100677

2 References 2 图像

概述

产品名称 小鼠CXCL16 ELISA试剂盒

检测方法 Colorimetric

精确度	批次内				
	样品	n	Mean	SD	CV%
	Overall				< 10%

	批次间				
	样品	n	Mean	SD	CV%
	Overall				< 12%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 1.3 pg/ml

范围 1.25 pg/ml - 80 pg/ml

回收率 94 %

特定样本回收率		
样品类型	平均%	范围
Cell culture supernatant	96.71	84% - 103%
Serum	94.34	83% - 103%
Plasma	93.49	82% - 102%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's CXCL16 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse CXCL16 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse CXCL16 coated on a 96-well plate.

Standards and samples are pipetted into the wells and CXCL16 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse CXCL16 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 CXCL16 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

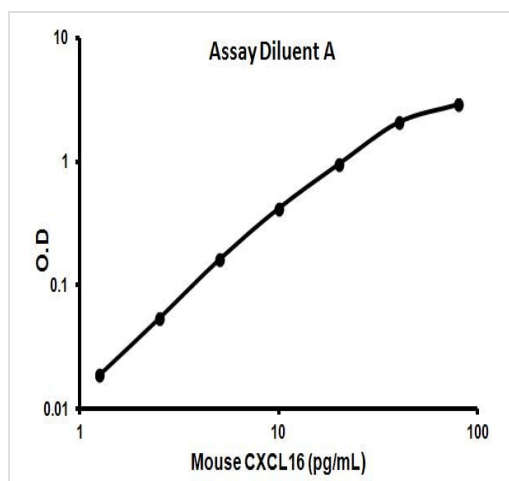
存放说明 Store at -20°C. Please refer to protocols.

组 件	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse CXCL16	2 vials
CXCL16 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse CXCL16 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

相关性 CXCL16 acts as a scavenger receptor on macrophages, which specifically binds to OxLDL (oxidized low density lipoprotein), suggesting that it may be involved in pathophysiology such as atherogenesis. It induces a strong chemotactic response and calcium mobilization. It binds to CXCR6/Bonzo.

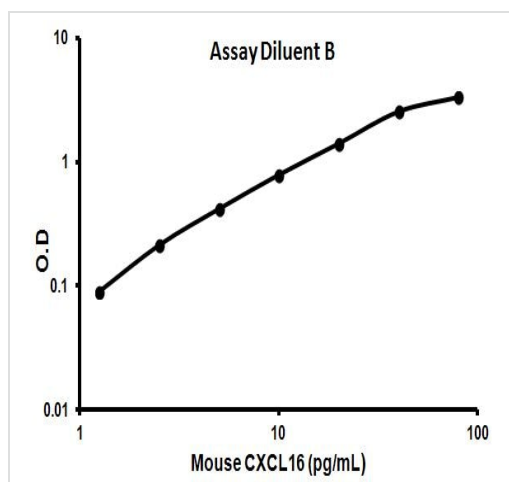
细胞定位 Cell Membrane and Secreted

图片



Typical Standard Curve

Representative standard curve using ab100677



Typical Standard Curve

Representative standard curve using ab100677

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