

Mouse CXCL16 ELISA Kit ab100677

2 References 画像数 2

製品の概要

製品名	Mouse CXCL16 ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				< 10%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				< 12%
サンプルの種類	Cell culture supernatant, Serum, Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 1.3 pg/ml			
検出範囲	1.25 pg/ml - 80 pg/ml			
添加回収試験	94 %			

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 supernatant	96.71	84% - 103%
Serum	94.34	83% - 103%
Plasma	93.49	82% - 102%

ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	Abcam's CXCL16 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> 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.

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
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

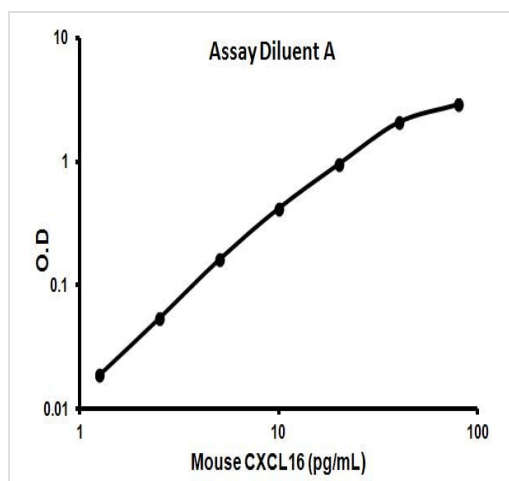
関連性

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.

細胞内局在

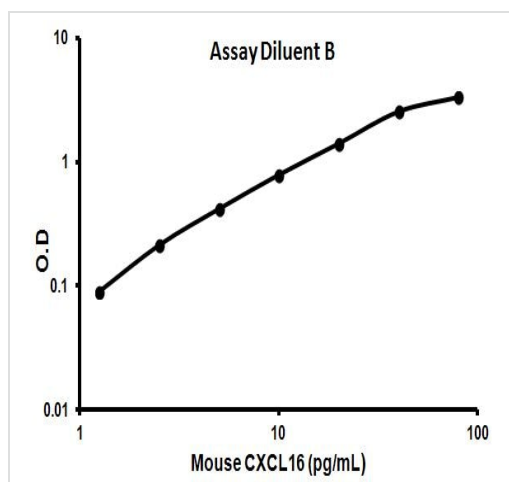
Cell Membrane and Secreted

画像



Representative standard curve using ab100677

Typical Standard Curve



Representative standard curve using ab100677

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.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