

Mouse MCP5 ELISA Kit ab100723

画像数 2

製品の概要

製品名	Mouse MCP5 ELISA Kit
検出方法	Colorimetric
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	< 0.5 pg/ml
検出範囲	0.41 pg/ml - 100 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 supernatant	95.34	84% - 105%
Serum	92.42	81% - 103%
Plasma	94.14	82% - 104%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要

Abcam's MCP5 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse MCP5 in serum, plasma and cell culture supernates.

This assay employs an antibody specific for mouse MCP5 coated on a 96-well plate. Standards and samples are pipetted into the wells and MCP5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse MCP5 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 MCP5 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

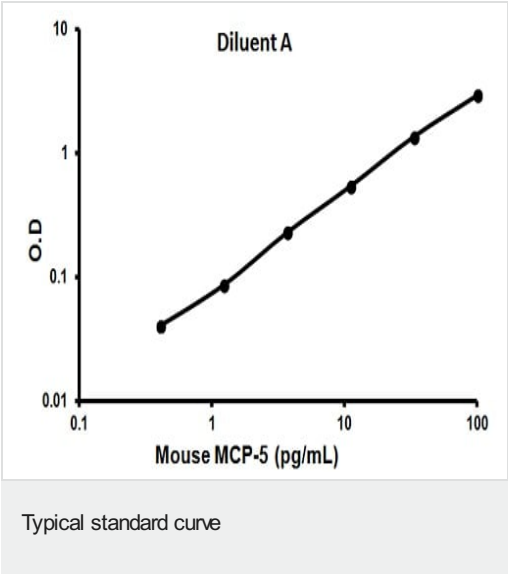
保存方法

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

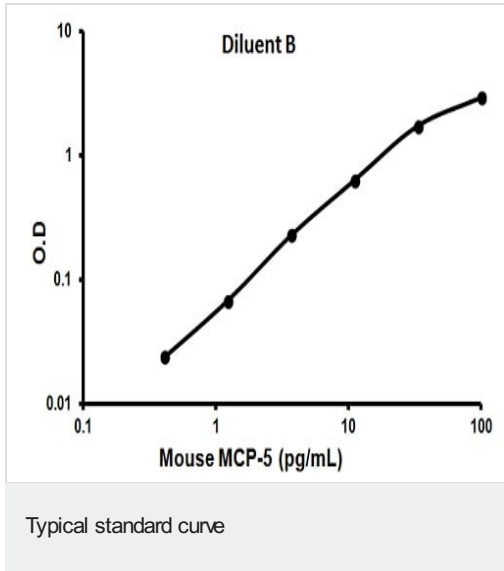
内容	1 x 96 tests
160X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-mouse MCP5	2 vials
MCP5 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse MCP5 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

関連性 MCP5 is a member of the b-chemokine family (C-C) of cytokines. This chemokine is strongly chemotactic for human monocytes and murine peritoneal macrophages, but is not active on mouse neutrophils and shows only a weak response on mouse eosinophils at high concentrations

画像



Representative standard curve using ab100723



Representative standard curve using ab100723

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