

Mouse GM-CSF ELISA Kit ab46078

2 References [画像数 1](#)

製品の概要

製品名 Mouse GM-CSF 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	CV%
A	8	279.2pg/ml	14.7	5.2%
B	8	62.5pg/ml	2.52	4%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
A	7	271.9pg/ml	10.8	3.9%
B	7	65.4pg/ml	4.3	6.5%

サンプルの種類 Cell culture supernatant, Serum

アッセイタイプ Sandwich (quantitative)

検出感度 < 9 pg/ml

検出範囲 15.6 pg/ml - 500 pg/ml

全工程の試験時間 3h 45m

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Abcam's GM-CSF mouse *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of GM-CSF in mouse serum, buffered solutions and cell culture medium.

A monoclonal antibody specific for GM-CSF has been coated onto the wells of the microtiter strips provided. Samples, including standards of known GM-CSF concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for GM-CSF are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is

added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of GM-CSF present in the samples.

This kit will recognize both endogenous and recombinant mouse GM-CSF.

Get higher sensitivity in only 90 minutes with Mouse GM-CSF ELISA Kit (CSF2) ([ab201276](#)) from our SimpleStep ELISA[®] range.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
Biotinylated anti-GM-CSF	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
GM-CSF Microplate (12 x 8 well strips)		1 unit	2 units
GM-CSF Standard (Lyophilized)	Yellow	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

機能

Cytokine that stimulates the growth and differentiation of hematopoietic precursor cells from various lineages, including granulocytes, macrophages, eosinophils and erythrocytes.

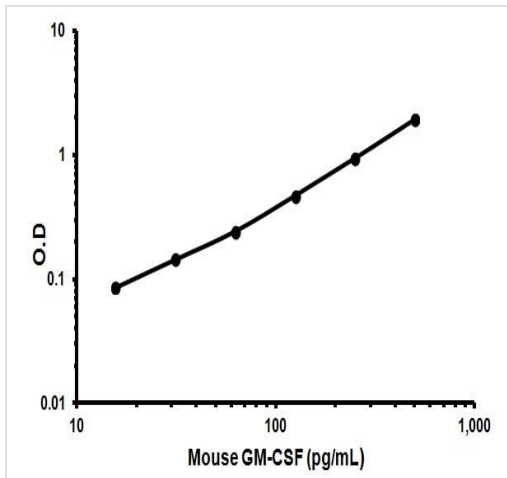
配列類似性

Belongs to the GM-CSF family.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab46078

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