

Rat GM-CSF ELISA Kit ab100762

画像数 2

製品の概要

製品名	Rat GM-CSF 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	< 10 pg/ml
検出範囲	12.35 pg/ml - 3000 pg/ml
添加回収試験	100 %

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	90.53	81% - 101%
Serum	136.4	128% - 143%
Plasma	118.8	88% - 136%

ステップ Multiple steps standard assay

種交差性 交差種: Rat

製品の概要

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

This assay employs an antibody specific for rat GM-CSF coated on a 96-well plate. Standards and samples are pipetted into the wells and GM-CSF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti- rat GM-CSF 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 GM-CSF 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
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Rat GM-CSF	2 vials
GM-CSF Microplate (12 x 8 wells)	1 unit
Recombinant Rat GM-CSF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能

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

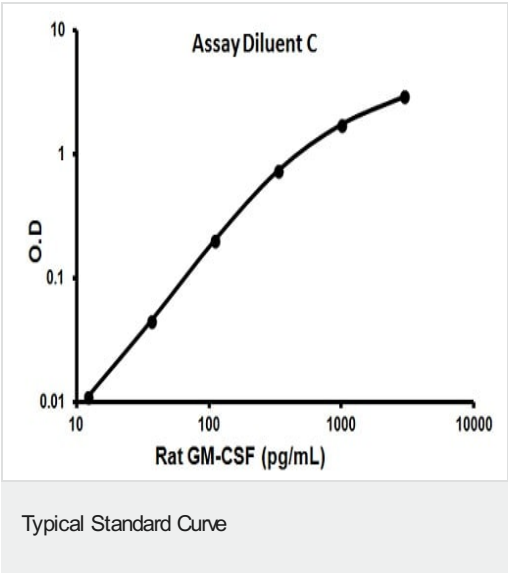
配列類似性

Belongs to the GM-CSF family.

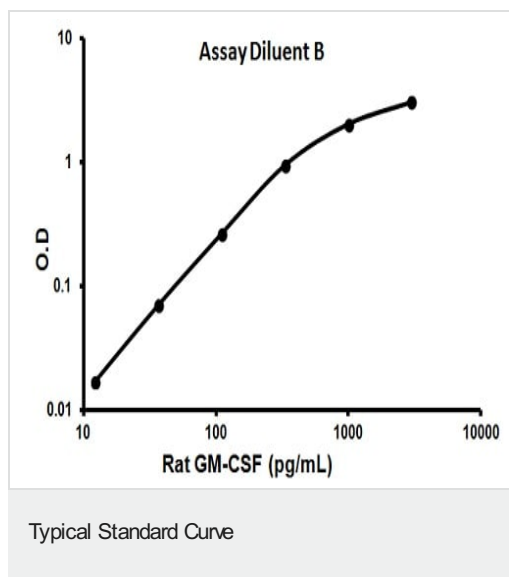
細胞内局在

Secreted.

画像



Representative standard curve using ab100762



Representative standard curve using ab100762

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