

Product datasheet

Mouse RANTES ELISA Kit (CCL5) ab100739

1 References [画像数 2](#)

製品の概要

製品名	Mouse RANTES ELISA Kit (CCL5)
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	< 1 pg/ml
検出範囲	0.82 pg/ml - 200 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	114.7	86% - 122%
Serum	109.9	98% - 119%
Plasma	75.41	68% - 87%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要

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

This assay employs an antibody specific for mouse RANTES coated on a 96-well plate. Standards and samples are pipetted into the wells and RANTES present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse RANTES 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 RANTES bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Mouse RANTES ELISA Kit ([ab215537](#)) from our SimpleStep

ELISA[™] range.

アプリケーション 適用あり: Sandwich ELISA
試験プラットフォーム Microplate

法規制情報

医薬用外毒物

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 Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Assay Diluent C	1 x 30ml
Biotinylated anti-mouse RANTES (lyophilized)	2 vials
RANTES Microplate (12 x 8 wells)	1 unit
Recombinant Mouse RANTES Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能 Chemoattractant for blood monocytes, memory T-helper cells and eosinophils. Causes the release of histamine from basophils and activates eosinophils. Binds to CCR1, CCR3, CCR4 and CCR5. One of the major HIV-suppressive factors produced by CD8+ T-cells. Recombinant RANTES protein induces a dose-dependent inhibition of different strains of HIV-1, HIV-2, and simian immunodeficiency virus (SIV). The processed form RANTES(3-68) acts as a natural chemotaxis inhibitor and is a more potent inhibitor of HIV-1-infection. The second processed form RANTES(4-68) exhibits reduced chemotactic and HIV-suppressive activity compared with RANTES(1-68) and RANTES(3-68) and is generated by an unidentified enzyme associated with monocytes and neutrophils.

組織特異性 T-cell and macrophage specific.

配列類似性 Belongs to the intercrine beta (chemokine CC) family.

翻訳後修飾 N-terminal processed form RANTES(3-68) is produced by proteolytic cleavage, probably by DPP4, after secretion from peripheral blood leukocytes and cultured sarcoma cells.

The identity of the O-linked saccharides at Ser-27 and Ser-28 are not reported in PubMed:1380064. They are assigned by similarity.

細胞内局在

Secreted.

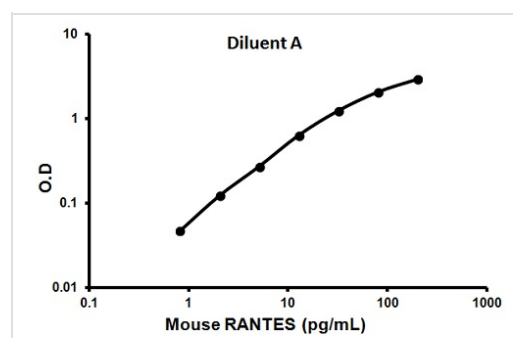
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab100739** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

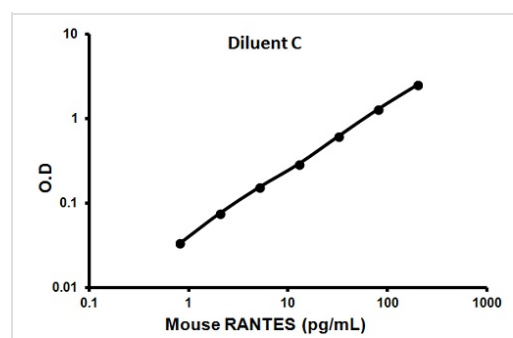
アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



Representative standard curve using
ab100739

Typical standard curve



Representative standard curve using
ab100739

Typical standard curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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