

## Product datasheet

# Mouse Exodus 2 ELISA Kit ab100667

### 画像数 2

#### 製品の概要

製品名	Mouse Exodus 2 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	< 20 pg/ml
検出範囲	31.25 pg/ml - 1000 pg/ml
添加回収試験	95 %

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	95.88	86% - 107%
Serum	95.51	84% - 105%
Plasma	94.29	83% - 102%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

#### 製品の概要

Abcam's Exodus 2 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse Exodus 2 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse Exodus 2 coated on a 96-well plate. Standards and samples are pipetted into the wells and Exodus 2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse Exodus 2 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 Exodus 2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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

## 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
Biotinylated anti-Mouse Exodus 2	2 vials
Exodus 2 Microplate (12 strips x 8 wells)	1 unit
Recombinant Mouse Exodus 2 Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### 関連性

Exodus 2 is one of several CC cytokine genes clustered on the p arm of chromosome 9. Cytokines are a family of secreted proteins involved in immunoregulatory and inflammatory processes. The CC cytokines are proteins characterised by two adjacent cysteines. Similar to other chemokines the protein encoded by this gene inhibits hemopoiesis and stimulates chemotaxis. This protein is chemotactic in vitro for thymocytes and activated T cells, but not for B cells, macrophages, or neutrophils. The cytokine encoded by this gene may also play a role in mediating homing of lymphocytes to secondary lymphoid organs. It is a high affinity functional ligand for chemokine receptor 7 (CCR7) that is expressed on T and B lymphocytes and a known receptor for another member of the cytokine family (small inducible cytokine A19).

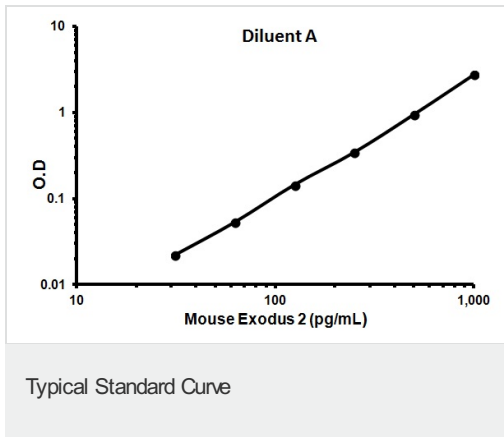
### アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab100667** in the following tested applications.

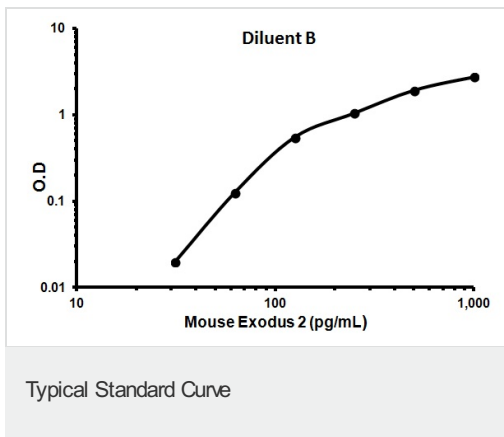
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

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 ab100667.



Representative Standard Curve using ab100667.

**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