

## Product datasheet

# Mouse GRO alpha ELISA Kit (CXCL1) ab100717

### 画像数 1

#### 製品の概要

製品名	Mouse GRO alpha ELISA Kit (CXCL1)
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	0.614 pg/ml - 150 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	123.5	115% - 139%
Serum	89.3	78% - 97%
Plasma	101.4	88% - 113%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

#### 製品の概要

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

This assay employs an antibody specific for mouse GRO alpha coated on a 96-well plate. Standards and samples are pipetted into the wells and GRO alpha present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse GRO alpha 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 GRO alpha 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	1 x 15ml
Biotinylated anti-mouse GRO alpha	2 vials
GRO alpha Microplate (12 x 8 wells)	1 unit
Recombinant mouse GRO alpha Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### 関連性

The GRO gene was originally identified by subtractive hybridization studies between normal and tumorigenic Chinese hamster embryo fibroblasts. The hamster cDNA was cloned and used as a probe for cloning of the human GRO cDNA. The GROalpha gene initially cloned from T24 cells and the gene in melanoma cells encoding melanoma growth stimulating protein (MGSA) are identical. Human cells contain three closely related, but distinct GRO genes: GRO alpha, GRO beta, and GRO gamma. GRO beta and GRO gamma share 93% and 82% identity, respectively, with GRO alpha at the nucleotide level. GROs are members of the chemokine alpha family that is characterized by the separation with one amino acid of the first two cysteine residues, C-X-C, in the amino acid sequence. The GRO gene has been mapped to chromosome 4q21. In normal cells, human mRNA GRO expression is found in foreskin fibroblasts, synovial fibroblasts, chondrocytes and osteocytes. Additionally, GRO mRNA has been detected in mammary fibroblasts, mammary epithelial cells, endothelial cells, activated monocytes, macrophages, and neutrophils. Characterization of the GROalpha receptor indicates the presence of low and high affinity receptors on human neutrophils.

### 細胞内局在

Secreted

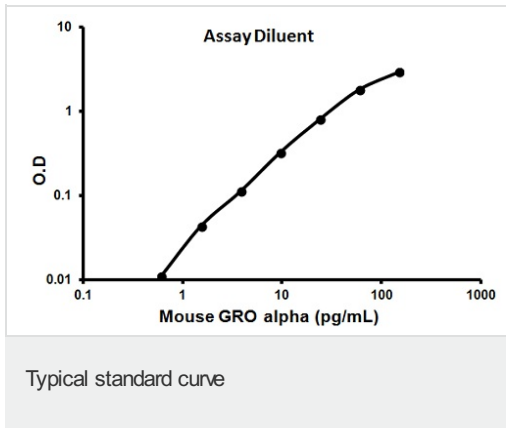
## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab100717** 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  
ab100717

**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