

Rat CXCL5 ELISA Kit ab100775

画像数 1

製品の概要

| | |
|---------|--|
| 製品名 | Rat CXCL5 ELISA Kit |
| 検出方法 | Colorimetric |
| サンプルの種類 | Cell culture extracts, Tissue Extracts |
| アッセイタイプ | Sandwich (quantitative) |
| 検出感度 | < 15 pg/ml |
| 検出範囲 | 8.23 pg/ml - 6000 pg/ml |
| 添加回収試験 | 92 % |

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 extracts | 91.25 | 82% - 102% |
| Tissue Extracts | 93.47 | 84% - 103% |

| | |
|------|-------------------------------|
| ステップ | Multiple steps standard assay |
| 種交差性 | 交差種: Rat |

製品の概要

Abcam's CXCL5 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat CXCL5 cell lysate and tissue lysate.

This assay employs an antibody specific for Rat CXCL5 coated on a 96- well plate. Standards and samples are pipetted into the wells and CXCL5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat CXCL5 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 CXCL5 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

特記事項

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

| | |
|------------|------------|
| 試験プラットフォーム | Microplate |
|------------|------------|

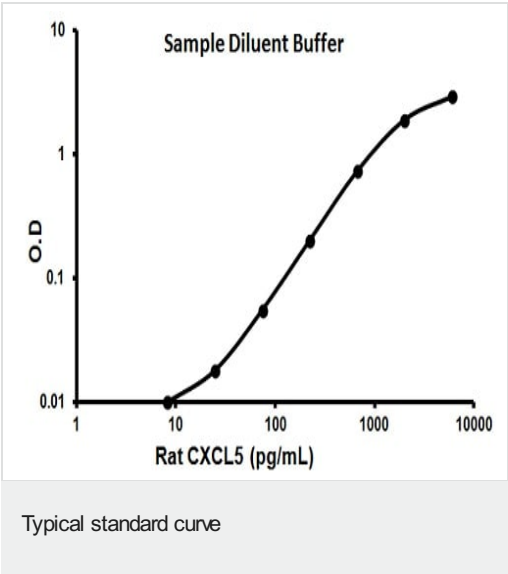
製品の特性

保存方法Store at -20°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 2X Cell Lysis Buffer | 1 x 5ml |
| 400X HRP-Streptavidin Concentrate | 1 x 200µl |
| 5X Assay Diluent | 1 x 15ml |
| 5X Sample Diluent Buffer | 1 x 10ml |
| Biotinylated anti-Rat CXCL5 | 2 vials |
| CXCL5 Microplate (12 x 8 wells) | 1 unit |
| Recombinant Rat CXCL5 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

関連性CXCL5 is a 9.9kD inflammatory chemokine that belongs to the CXC chemokine family. CXCL5 is expressed by fibroblasts, is induced by bacterial lipopolysaccharides, and is chemotactic for neutrophils. This chemokine is produced concomitantly with Interleukin 8 (IL8) in response to stimulation with either Interleukin 1 (IL1) or Tumor Necrosis Factor alpha (TNFA). This chemokine is a potent chemotaxin involved in neutrophil activation.

画像



Representative standard curve using ab100775

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