

Rabbit IL-4 ELISA Kit ab273244

画像数 1

製品の概要

| | | | | |
|------|-----------------------|--|--|--|
| 製品名 | Rabbit IL-4 ELISA Kit | | | |
| 検出方法 | Colorimetric | | | |
| 再現性 | Intra-Assay (同時再現性) | | | |

| サンプル | N | 平均値 | SD | CV% |
|---------|---|-----|----|-------|
| Overall | | | | < 10% |

| Inter-Assay (日差再現性) | | | | |
|---------------------|---|-----|----|-------|
| サンプル | N | 平均値 | SD | CV% |
| Overall | | | | < 12% |

| | |
|---------|---|
| 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 | 13.7 pg/ml |
| 検出範囲 | 13.72 pg/ml - 10000 pg/ml |

| 添加回収試験 | 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 |
| Serum | 132.2 | 121% - 142% |
| Plasma | 115.6 | 104% - 126% |
| Cell culture media | 120.4 | 102% - 133% |

| | |
|------|-------------------------------|
| ステップ | Multiple steps standard assay |
|------|-------------------------------|

| | |
|------|-------------|
| 種交差性 | 交差種: Rabbit |
|------|-------------|

| | |
|-------|---|
| 製品の概要 | Rabbit IL-4 ELISA Kit (ab273244) is an <i>in-vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Rabbit IL-4 in serum (Rabbit IL-4 concentration is low in normal serum/plasma, and may not be detectable in this assay), plasma and cell culture supernatants. |
|-------|---|

This assay employs an antibody specific for Rabbit IL-4 coated on a 96-well plate. Standards and

samples are pipetted into the wells and IL-4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed, and biotinylated anti-Rabbit IL-4 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 IL-4 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

Pre-coated microplate (12 x 8 well strips)

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 |
| 5X Assay Diluent B | 1 x 15ml |
| 800X HRP-Streptavidin Concentrate | 1 x 200µl |
| Anti-Rabbit IL-4 coated Microplate (12 x 8 wells) | 1 unit |
| Assay Diluent C | 1 x 30ml |
| Biotinylated Anti-Rabbit IL-4 Antibody | 2 vials |
| Rabbit IL-4 Standard (Lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB Substrate Solution | 1 x 12ml |

機能

Participates in at least several B-cell activation processes as well as of other cell types. It is a costimulator of DNA-synthesis. It induces the expression of class II MHC molecules on resting B-cells. It enhances both secretion and cell surface expression of IgE and IgG1. It also regulates the expression of the low affinity Fc receptor for IgE (CD23) on both lymphocytes and monocytes.

関連疾患

Genetic variations in IL4 may be a cause of susceptibility to ischemic stroke (ISCHSTR) [MIM:601367]; also known as cerebrovascular accident or cerebral infarction. A stroke is an acute neurologic event leading to death of neural tissue of the brain and resulting in loss of motor, sensory and/or cognitive function. Ischemic strokes, resulting from vascular occlusion, is considered to be a highly complex disease consisting of a group of heterogeneous disorders with multiple genetic and environmental risk factors.

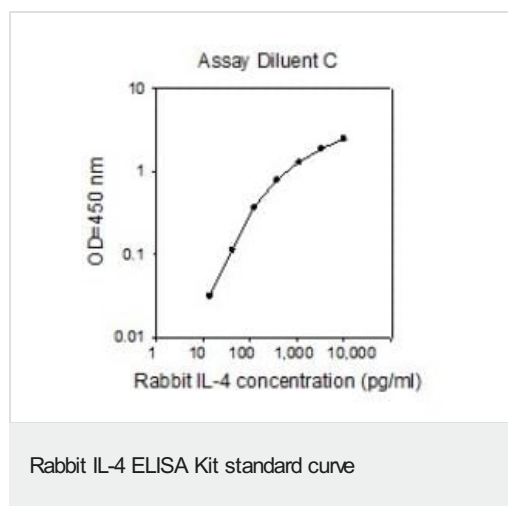
配列類似性

Belongs to the IL-4/IL-13 family.

細胞内局在

Secreted.

画像



This standard curve is for demonstration only. A standard curve must be run with each assay.

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