

Rat IL-4 ELISA Kit ab100771

4 References 1 Image

Overview

|                  |  |
|------------------|--|
| Product name     | Rat IL-4 ELISA Kit                     |
| Detection method | Colorimetric                           |
| Sample type      | Cell culture extracts, Tissue Extracts |
| Assay type       | Sandwich (quantitative)                |
| Sensitivity      | < 1.5 pg/ml                            |
| Range            | 0.66 pg/ml - 160 pg/ml                 |
| Recovery         | 108 %                                  |

Sample specific recovery

| Sample type           | Average % | Range       |
|-----------------------|-----------|-------------|
| Cell culture extracts | 93.49     | 85% - 105%  |
| Tissue Extracts       | 123.9     | 114% - 133% |

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Rat

Product overview Abcam's IL-4 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat IL-4 in cell lysates and tissue lysates.

This assay employs an antibody specific for Rat 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-Rat 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.

Notes 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.

Platform Microplate

## Properties

### Storage instructions

Store at -20°C. Please refer to protocols.

| Components                                  | 1 x 96 tests |
|---|--------------|
| 20X Wash Buffer                             | 1 x 25ml     |
| 2X Cell Lysis Buffer                        | 1 x 5ml      |
| 5X Assay Diluent                            | 1 x 15ml     |
| 5X Sample Diluent Buffer                    | 1 x 10ml     |
| 600X HRP-Streptavidin Concentrate           | 1 x 200µl    |
| Biotinylated anti-Rat IL-4                  | 2 vials      |
| IL-4 Microplate (12 x 8 wells)              | 1 unit       |
| Recombinant rat IL-4 Standard (lyophilized) | 2 vials      |
| Stop Solution                               | 1 x 8ml      |
| TMB One-Step Substrate Reagent              | 1 x 12ml     |

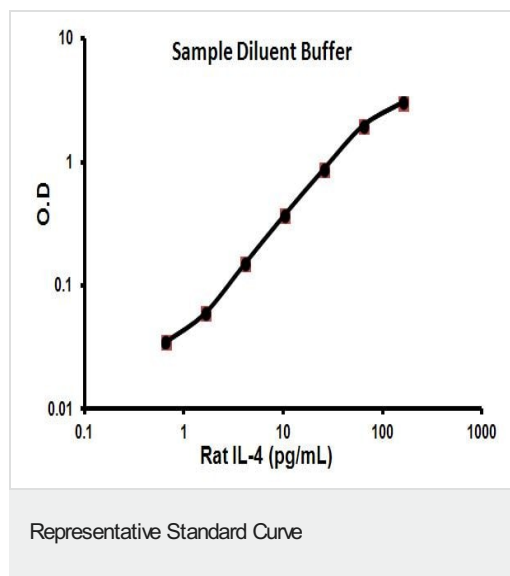
### Relevance

IL-4 is a pleiotropic cytokine produced by activated T cells. It is a ligand for interleukin 4 receptor. The interleukin 4 receptor also binds to IL13, which may contribute to many overlapping functions of this cytokine and IL13. STAT6, a signal transducer and activator of transcription, has been shown to play a central role in mediating the immune regulatory signal of IL-4. IL-4, IL3, IL5, IL13, and CSF2 form a cytokine gene cluster on chromosome 5q, with this gene particularly close to IL13. IL-4, IL13 and IL5 are found to be regulated coordinately by several long-range regulatory elements in an over 120 kilobase range on the chromosome. Two alternatively spliced transcript variants of IL-4 encoding distinct isoforms have been reported.

### Cellular localization

Secreted

## Images



Representative Standard Curve using ab100771

**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.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors