

Product datasheet

Rat IL-4 ELISA Kit (Interleukin-4) ab100771

[1 Image](#)

Overview

Product name	Rat IL-4 ELISA Kit (Interleukin-4)
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.

Tested applications **Suitable for:** Sandwich ELISA

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
2x Cell lysate 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 (lyophilized)	2 vials
IL-4 Microplate (12 x 8 wells)	1 unit
Recombinant rat IL-4 Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

IL4 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 IL4. IL4, IL3, IL5, IL13, and CSF2 form a cytokine gene cluster on chromosome 5q, with this gene particularly close to IL13. IL4, 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 IL4 encoding distinct isoforms have been reported.

Cellular localization

Secreted

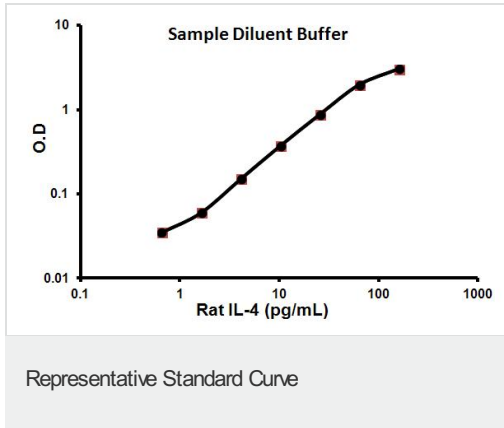
Applications

Our [Abpromise guarantee](#) covers the use of **ab100771** in the following tested applications.

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

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative Standard Curve using
ab100771

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.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