

Rat IL-13 ELISA Kit ab100766

★★★★★ 1 Abreviews 5 References 1 Image

Overview

Product name	Rat IL-13 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			

Sample	n	Mean	SD	CV%
Overall				< 10%

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type	Cell culture extracts, Tissue Extracts
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	27.43 pg/ml - 20000 pg/ml
Recovery	114 %

Sample type	Average %	Range
Cell culture extracts	120.4	80% - 139%
Tissue Extracts	107.8	81% - 135%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Rat
Product overview	<p>Abcam's IL-13 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Rat IL-13 cell lysate and tissue lysate.</p> <p>This assay employs an antibody specific for Rat IL-13 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-13 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat IL-13 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the</p>

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-13 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
700X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated anti-rat IL-13	2 vials
IL-13 Microplate (12 x 8 wells)	1 unit
Recombinant Rat IL-13 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

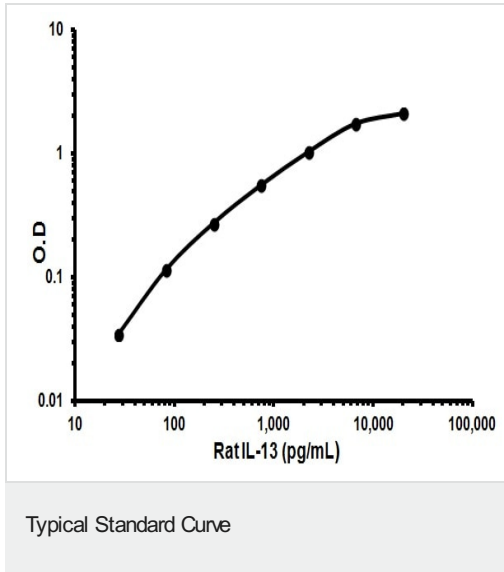
Function Cytokine (PubMed:8096327, PubMed:8097324). Inhibits inflammatory cytokine production (PubMed:8096327). Synergizes with IL2 in regulating interferon-gamma synthesis (PubMed:8096327). May be critical in regulating inflammatory and immune responses (PubMed:8096327, PubMed:8097324). Positively regulates IL31RA expression in macrophages.

Involvement in disease Allergic rhinitis

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

Cellular localization Secreted.

Images



Representative standard curve using ab100766

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