

Human IL-4 Receptor ELISA Kit ab46022

1 References [1 图像](#)

概述

产品名称 人IL-4 Receptor ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Sample A	8	563.7pg/ml	17.8	3.1%
Sample B	8	140.6pg/ml	9.6	6.8%

批次间

样品	n	Mean	SD	CV%
Sample A	8	530.5pg/ml	35.1	6.6%
Sample B	8	106.5pg/ml	7.58	7.1%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 11.5 pg/ml

范围 31.25 pg/ml - 1000 pg/ml

检测时间 1h 45m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Human IL-4 Receptor (IL-4R) in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of soluble IL-4 receptor molecule (CD124) in Human serum, plasma, buffered solutions or cell culture medium.

A monoclonal antibody specific for IL-4R has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-4R concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-4R are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and

washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-4R present in the samples.

This kit will recognize both endogenous and recombinant Human IL-4R.

说明 The standard is affinity purified.

平台 Microplate

性能

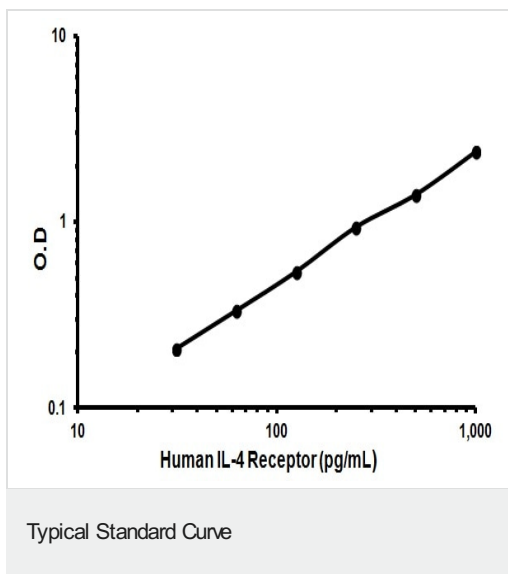
存放说明 Store at +4°C. Please refer to protocols.

组件	标识符	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
Biotinylated anti-IL-4R	Red	1 x 0.4ml	2 x 0.4ml
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-4R Microplate (12 x 8 well strips)		1 unit	2 units
IL-4R Standard (Lyophilized)	Orange	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

相关性 IL-4R binds interleukin 4 and interleukin 13 to regulate IgE production. It also can bind interleukin 4 to promote differentiation of Th2 cells. A soluble form of the IL-4R can be produced by an alternate splice variant or by proteolysis of the membrane-bound protein, and this soluble form can inhibit IL4-mediated cell proliferation and IL5 upregulation by T-cells. Allelic variations in this gene have been associated with atopy, a condition that can manifest itself as allergic rhinitis, sinusitis, asthma, or eczema. Two transcript variants encoding different isoforms, a membrane-bound and a soluble form, have been found for this gene.

细胞定位 Cell membrane; single-pass type I membrane protein. Isoform 2: Secreted protein

图片



Representative Standard Curve using ab46022

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.cn/abpromise> or contact our technical team.

Terms and conditions

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