

Rat IL-4 ELISA Set (without plates) ab48423

1 图像

概述

产品名称	大鼠IL-4 ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 15 pg/ml
范围	31.25 pg/ml - 1000 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Rat
产品概述	<p>Abcam's IL-4 Rat <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Rat IL-4 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The rat IL-4 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>This ELISA Set is intended to measure rat IL-4 in cell -culture supernatants, as well as in serum samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>Recognizes both natural and recombinant Rat IL-4 Range: 31.25 pg/ml - 1000 pg/ml Sensitivity < 15 pg/ml</p>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

性能

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

组件	标识符	15 x 96 tests	10 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution		6 vials	4 vials	2 vials

组件	标识符	15 x 96 tests	10 x 96 tests	5 x 96 tests
Detection biotinylated anti-Rat IL-4 antibody		3 x 0.55ml	2 x 0.55ml	1 x 0.55ml
Rat IL-4 Capture antibody		3 x 0.5ml	2 x 0.5ml	1 x 0.5ml
Rat IL-4 Standard (Lyophilized)	Yellow	15 vials	10 vials	5 vials
Streptavidin-HRP		3 vials	2 vials	1 vial

相关性 IL-4 is a pleiotropic cytokine produced by activated T cells, mast cells and basophils. 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. IL-4 elicits many different biological responses, but has two dominant functions. The first is regulating differentiation of naïve CD4+ T cell to the Th2 type. Th2 cells produce IL-4, IL5, IL10 and IL13, which tend to favor a humoral immune response while suppressing a cell mediated immune response controlled by Th1 cells. STAT6, a signal transducer and activator of transcription, has been shown to play a central role in mediating the immune regulatory signal of this cytokine. The second is regulating IgE and IgG1 production by B cells. Two alternatively spliced transcript variants of this gene encoding distinct isoforms have been reported.

细胞定位 Secreted

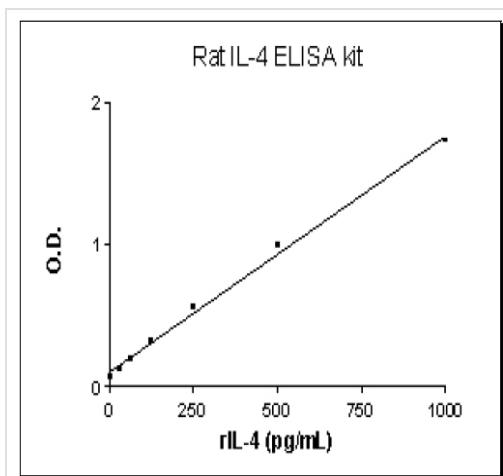
应用

The Abpromise guarantee Abpromise™承诺保证使用ab48423于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



Rat IL-4 Standard curve ranging from 31.25 to 1000 pg/mL

Rat IL-4 ELISA Set (ab48423)

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