

画像数 1

製品名	Rat IL-4 ELISA Set (without plates)
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 15 pg/ml
検出範囲	31.25 pg/ml - 1000 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Rat
製品の概要	<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>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	15 x 96 tests	10 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution		6 vials	4 vials	2 vials

内容	ラベル	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

関連性 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.

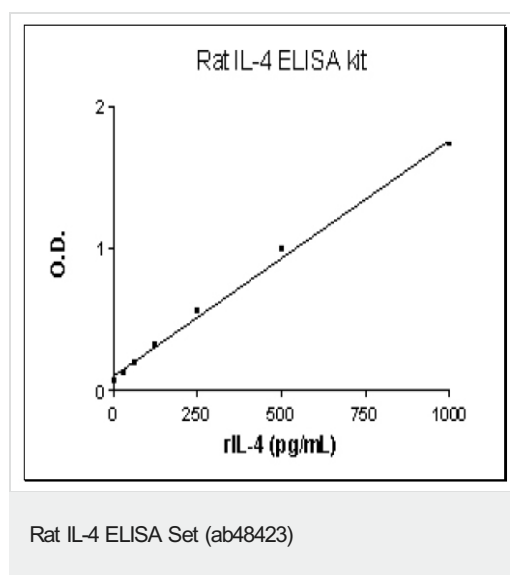
細胞内局在 Secreted

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab48423の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



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

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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