

Product datasheet

Rat IL-1 beta ELISA Kit (Interleukin-1 beta) ab100767

3 References 2 图像

概述

产品名称	大鼠IL-1 beta ELISA试剂盒(Interleukin-1 beta)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 80 pg/ml
范围	68.59 pg/ml - 50000 pg/ml
复检	99 %

特定样本复检结果

样品类型	平均%	范围
Cell culture supernatant	92.89	81% - 109%
Serum	107.2	95% - 115%
Plasma	97.55	89% - 108%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 Abcam's IL-1 beta (Interleukin-1 beta) Rat ELISA (Enzyme-Linked Immunosorbent Assay) Kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of rat IL-1 beta in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for IL-1 beta coated on a 96-well plate. Standards and samples are pipetted into the wells and IL- 1 beta present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-rat IL-1 beta 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 colour develops in proportion to the amount of IL-1 beta bound. The Stop Solution changes the colour from blue to yellow, and the intensity of the colour is measured at 450 nm.

经测试应用 适用于: Sandwich ELISA

平台 Microplate

性能

存放说明

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

组件	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-rat IL-1 beta	2 vials
IL-1 beta Microplate (12 x 8 wells)	1 unit
Recombinant rat IL-1 beta Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

功能

Produced by activated macrophages, IL-1 stimulates thymocyte proliferation by inducing IL-2 release, B-cell maturation and proliferation, and fibroblast growth factor activity. IL-1 proteins are involved in the inflammatory response, being identified as endogenous pyrogens, and are reported to stimulate the release of prostaglandin and collagenase from synovial cells.

序列相似性

Belongs to the IL-1 family.

结构域

The similarity among the IL-1 precursors suggests that the amino ends of these proteins serve some as yet undefined function.

细胞定位

Secreted. The lack of a specific hydrophobic segment in the precursor sequence suggests that IL-1 is released by damaged cells or is secreted by a mechanism differing from that used for other secretory proteins.

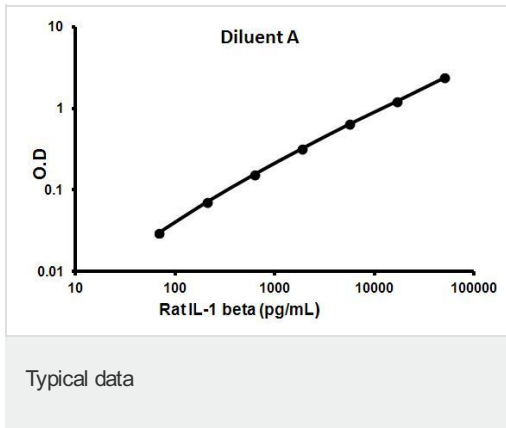
应用

Our [Abpromise guarantee](#) covers the use of **ab100767** in the following tested applications.

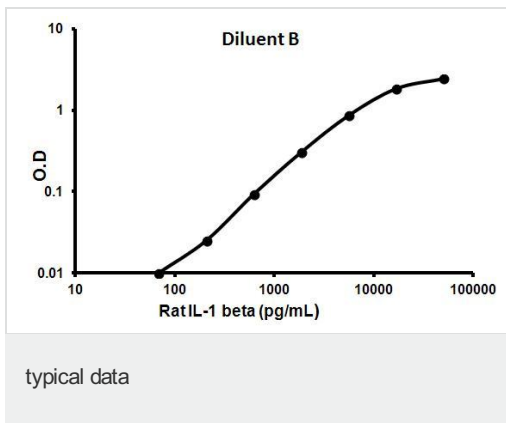
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

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

图片



Representative standard curve using ab100767



Representative standard curve using ab100767

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.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