

## 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
回收率	99 %

特定样本回收率

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.

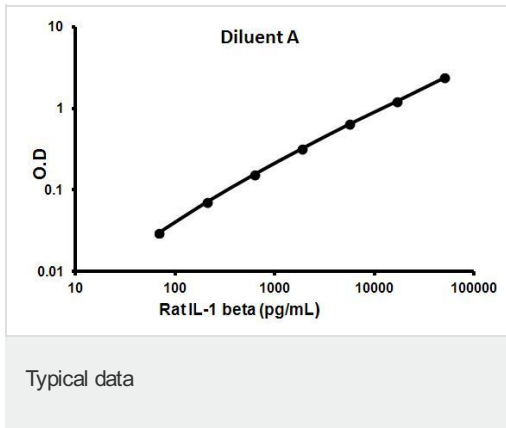
## 应用

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

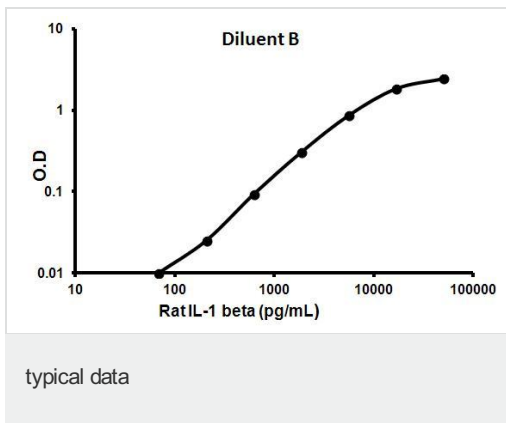
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

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