

Rat IL-1 alpha ELISA Kit ab113350

[3 References](#) [1 Image](#)

Overview

| | |
|-------------------------|---|
| Product name | Rat IL-1 alpha ELISA Kit |
| Detection method | Colorimetric |
| Sample type | Cell culture supernatant, Serum, Plasma |
| Assay type | Sandwich (quantitative) |
| Sensitivity | < 15 pg/ml |
| Range | 11 pg/ml - 8000 pg/ml |
| Recovery | 81 % |

Sample specific recovery

| Sample type | Average % | Range |
|--------------------------|-----------|-----------|
| Cell culture supernatant | 87.48 | 76% - 95% |
| Serum | 79.95 | 69% - 88% |
| Plasma | 75.73 | 68% - 84% |

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat

Product overview Abcam's IL-1 alpha Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat IL-1 alpha in serum, plasma and cell culture supernatants.

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

Platform Microplate

Properties

Storage instructions

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

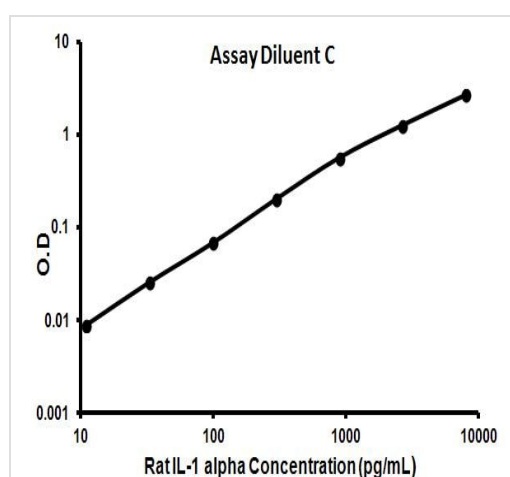
| Components | 1 x 96 tests |
|---|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 5X Assay Diluent B | 1 x 15ml |
| 500X HRP-Streptavidin Concentrate | 1 x 200µl |
| Assay Diluent C | 1 x 30ml |
| Biotinylated anti-Rat IL-1 alpha | 2 vials |
| IL-1 alpha Microplate (12 x 8 wells) | 1 unit |
| Recombinant rat IL-1 alpha Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

Relevance

IL-1 alpha is a member of the interleukin 1 cytokine family. It is a pleiotropic cytokine involved in various immune responses, inflammatory processes, and hematopoiesis. IL-1 alpha is produced by monocytes and macrophages as a proprotein, which is proteolytically processed and released in response to cell injury, and thus induces apoptosis. This gene and eight other interleukin 1 family genes form a cytokine gene cluster on chromosome 2. It has been suggested that the polymorphism of these genes is associated with rheumatoid arthritis and Alzheimer's disease.

Cellular localization

Secreted

Images

Representative standard curve using ab113350

Typical Data

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.com/abpromise> or contact our technical team.

Terms and conditions

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