

### Human IL-1 alpha ELISA Kit ab46028

#### 1 References [画像数 1](#)

#### 製品の概要

**製品名** Human IL-1 alpha ELISA Kit

**検出方法** Colorimetric

**再現性** Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Sample A	22	459.8pg/ml	13	2.8%
Sample B	21	72.6pg/ml	3.9	4.3%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Sample A	8	259.9pg/ml	15	5.8%
Sample B	8	58.6pg/ml	4.3	7.3%

**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** < 10 pg/ml

**検出範囲** 31.25 pg/ml - 1000 pg/ml

**添加回収試験** 95.3 %

**全工程の試験時間** 3h 45m

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Human

**製品の概要** Abcam's IL-1 alpha Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-1 alpha in cell culture supernatants, buffered solutions, serum, plasma and other body fluids. This assay will recognize both natural and recombinant Human IL-1a.

A monoclonal antibody specific for IL-1 alpha has been coated to the wells of the microtiter strip plate provided. Samples, including standards of known IL-1 alpha concentrations and unknown are pipetted into these wells. During incubation, binding of IL-1 alpha to the capture antibodies is

completed and then any excess unbound analyte is removed by washing. A biotinylated monoclonal antibody specific for IL-1 alpha is then incubated. After washing, the enzyme (streptavidin-peroxydase) is added. After incubation and washing to remove all the unbound enzyme, a substrate solution which is acting on the bound enzyme is added to induce a coloured reaction product. The intensity of this coloured product is directly proportional to the concentration of IL-1 alpha present in the sample.

This kit will recognize both endogenous and recombinant Human IL-1 alpha.

Get higher sensitivity in only 90 minutes with Human IL-1 alpha ELISA Kit ([ab178008](#)) from our SimpleStep ELISA® range.

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Biotinylated anti-IL-1 alpha	Red	2 x 0.4ml	1 x 0.4ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
IL-1 alpha Control	Silver	4 vials	2 vials
IL-1 alpha Microplate (12 x 8 well strips)		2 units	1 unit
IL-1 alpha Standard (Lyophilized)	Yellow	4 vials	2 vials
Standard Diluent (Serum)		2 x 7ml	1 x 7ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

#### 関連性

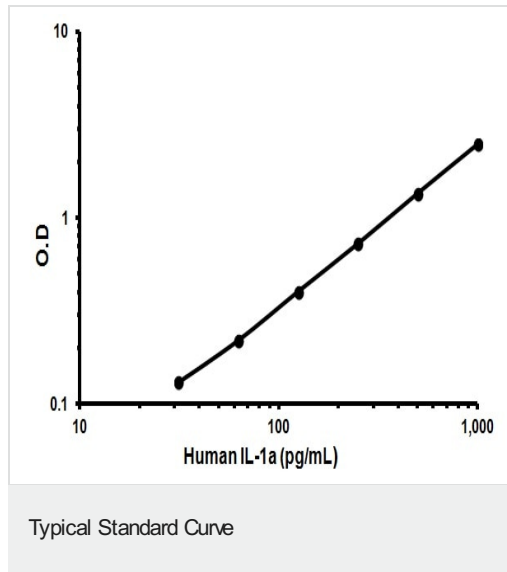
IL-1 alpha is a member of the interleukin 1 cytokine family. It is a pleiotropic cytokine involved in various immune responses, inflammatory processes, and hematopoiesis. IL-1 alpha is produced by monocytes and macrophages as a proprotein, which is proteolytically processed and released in response to 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.

細胞内局在

Secreted

画像



Representative Standard Curve using ab46028

**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