

Human IL-1 alpha ELISA Kit ab46028

1 References 1 Image

Overview

Product name	Human IL-1 alpha ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	22	459.8pg/ml	13	2.8%
	Sample B	21	72.6pg/ml	3.9	4.3%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	259.9pg/ml	15	5.8%
	Sample B	8	58.6pg/ml	4.3	7.3%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 10 pg/ml				
Range	31.25 pg/ml - 1000 pg/ml				
Recovery	95.3 %				
Assay time	3h 45m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's IL-1 alpha Human <i>in vitro</i> 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.

**Platform** Microplate

## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	Identifier	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

## Relevance

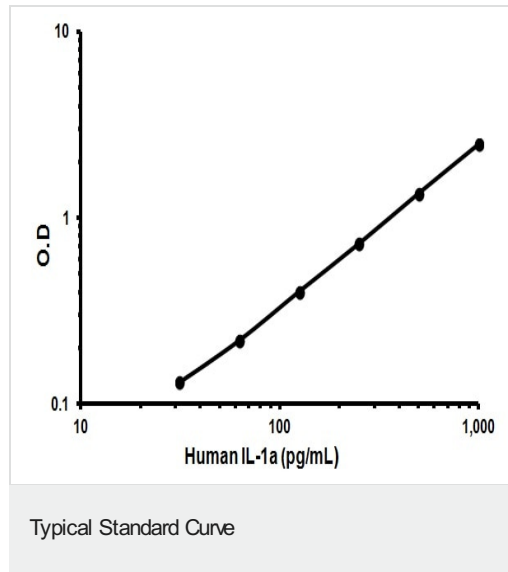
IL-1 alpha is a member of the interleukin 1 cytokine family. It is a pleiotropic cytokine involved in various immune responses, inflammatory processes, and hematopoiesis. IL-1 alpha is produced by monocytes and macrophages as a proprotein, which is proteolytically processed and released in response to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