

Human IL-1 alpha ELISA Set (without plates) ab47350

1 Image

Overview

Product name	Human IL-1 alpha ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	31.25 pg/ml - 1000 pg/ml
Assay time	3h 45m
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's IL-1α Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-1α in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA set is intended to measure human IL-1 alpha in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>The IL-1 alpha ELISA set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>Recognizes both natural and recombinant human IL-1a.</p>
Tested applications	Suitable for: Sandwich ELISA
Platform	Reagents

Properties

Storage instructionsStore at +4°C. Please refer to protocols.

Components	Identifier	5 x 96 tests	10 x 96 tests	15 x 96 tests
Chromogen TMB Substrate Solution		2 vials	4 vials	6 vials

Components	Identifier	5 x 96 tests	10 x 96 tests	15 x 96 tests
Detection biotinylated anti-IL-1 alpha antibody		1 x 0.55ml	2 x 0.55ml	3 x 0.55ml
IL-1 alpha Capture antibody		1 x 0.1ml	2 x 0.1ml	3 x 0.1ml
IL-1 alpha Standard (Lyophilized)	Yellow	5 vials	10 vials	15 vials
Streptavidin-HRP		1 vial	2 vials	3 vials

Relevance

IL-1 is a member of the interleukin 1 cytokine family. This cytokine is a pleiotropic cytokine involved in various immune responses, inflammatory processes, and hematopoiesis. This cytokine 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.

Applications

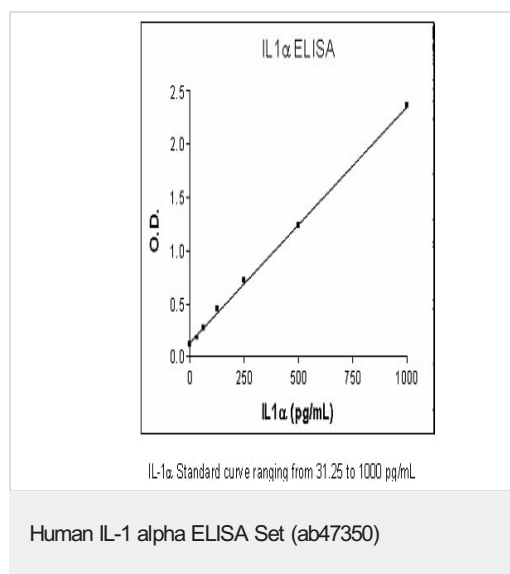
The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab47350 in the following tested applications.

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

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



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