

Product datasheet

Human IL-5 ELISA Set (without plates) ab47356

1 Image

Overview

Product name	Human IL-5 ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 5 pg/ml
Range	7.8 pg/ml - 250 pg/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	Abcam's IL-5 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-5 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

This ELISA Set is intended to measure human IL-5 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.

These reagents are intended for research use only, not for diagnostic purposes.

The IL-5 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.

Recognizes both natural and recombinant human IL-5.

Tested applications	Suitable for: Sandwich ELISA
Platform	Reagents

Properties

Storage instructions

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

Components	Identifier	5 x 96 tests	10 x 96 tests	15 x 96 tests
Detection biotinylated anti-IL-5 antibody		1 x 0.55ml	2 x 0.55ml	3 x 0.55ml
IL-5 Capture antibody		1 x 0.5ml	2 x 0.5ml	3 x 0.5ml
IL-5 standard	Yellow	5 vials	10 vials	15 vials
Chromogen TMB Substrate Solution		2 vials	4 vials	6 vials
Streptavidin-HRP		1 vial	2 vials	3 vials

Relevance

Interleukin 5 (IL-5) is a cytokine produced primarily by activated T lymphocytes. It exists as an antiparallel disulfide linked homodimeric glycoprotein with 115 amino acid residues in each chain. Known also as EDF (eosinophil differentiating factor), functions predominantly as an eosinophilopoietic factor. Analysis of its crystal structure reveals a novel two domain structure, with each domain showing significant homology to the cytokine fold in GM-CSF, MCSF, IL2, IL4 and growth hormone. Human and mouse IL5 have 70% amino acid sequence homology.

Cellular localization

Secreted

Applications

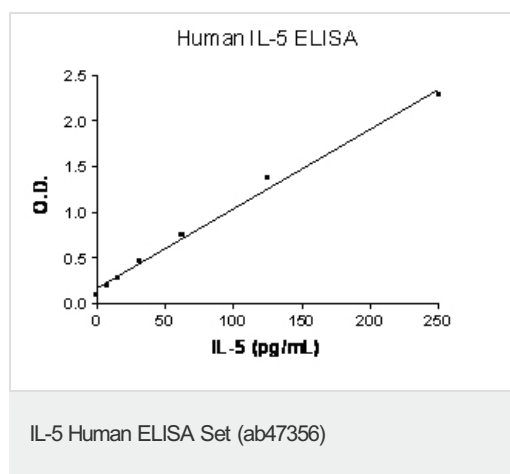
The Abpromise guarantee

Our [Abpromise guarantee](#) covers the use of ab47356 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



IL-5 Standard curve ranging from 7.8 to 250 pg/mL

IL-5 Human ELISA Set (ab47356)

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