

Human IL-7 ELISA Kit ab46031

1 References 1 Image

Overview

Product name Human IL-7 ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	198.7pg/ml	1.667	0.83%
	Sample B	8	28.69pg/ml	1.118	3.86%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Sample A	29	198.9pg/ml	3.772	1.89%
	Sample B	29	25.68pg/ml	1.505	5.84%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 3 pg/ml

Range 6.25 pg/ml - 200 pg/ml

Recovery 97.6 %

Assay time 2h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

**Product overview**  
Abcam's IL-7 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-7 in Human serum, plasma, buffered solutions and cell culture medium.

A monoclonal antibody specific for IL-7 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-7 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-7 are simultaneously incubated. After washing,

the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-7 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-7.

Get higher sensitivity in only 90 minutes with Human IL-7 ELISA Kit ([ab185984](#)) from our SimpleStep ELISA® range.

## Platform

Microplate

## Properties

### Storage instructions

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

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
Biotinylated anti-IL-7	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 23ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-7 Control	Silver	2 vials	4 vials
IL-7 Microplate (12 x 8 well strip)		1 unit	2 units
IL-7 standard	Yellow	2 vials	4 vials
Standard Diluent (Serum)		1 x 7ml	2 x 7ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

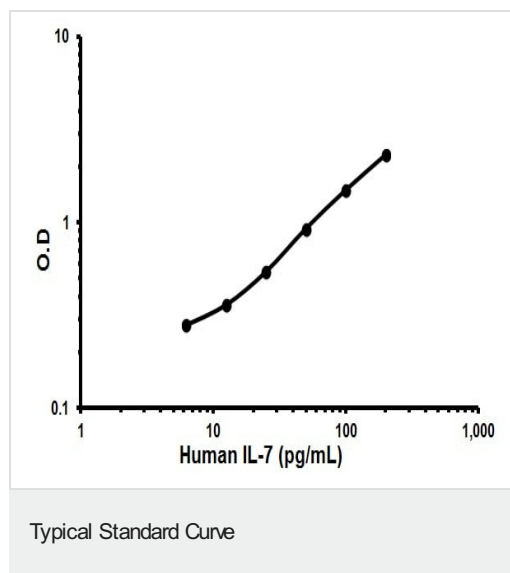
## Relevance

IL-7 is a cytokine important for B and T cell development. It and the hepatocyte growth factor form a heterodimer that functions as a pre-pro-B cell growth-stimulating factor. IL-7 is found to be a cofactor for V(D)J rearrangement of the T cell receptor beta (TCRB) during early T cell development. It can be produced locally by intestinal epithelial and epithelial goblet cells, and may serve as a regulatory factor for intestinal mucosal lymphocytes. Knockout studies in mice suggested that IL-7 plays an essential role in lymphoid cell survival.

## Cellular localization

Secreted

## Images



Representative Standard Curve using ab46031

**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