

Product datasheet

Mouse IL-2 ELISA Kit (Interleukin-2) ab46096

[1 References](#) [1 Image](#)

Overview

Product name Mouse IL-2 ELISA Kit (Interleukin-2)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
A	10	170pg/ml	13	7.8%

Inter-assay

Sample	n	Mean	SD	CV%
B	8	137pg/ml	8.9	6.4%

Sample type Cell culture supernatant, Serum

Assay type Sandwich (quantitative)

Sensitivity < 5 pg/ml

Range 15.6 pg/ml - 500 pg/ml

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview

Abcam's IL-2 (Interleukin-2) mouse *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-2 in mouse serum, buffered solutions and cell culture medium.

A monoclonal antibody specific for IL-2 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-2 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-2 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-2 present in the samples.

This kit will recognize both endogenous and recombinant mouse IL-2.

Tested applications

Suitable for: Sandwich ELISA

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 25ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
IL2 Biotinylated Detection antibody		1 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 23ml	1 x 23ml
IL-2 Microplate (12 x 8 well strips)		1 unit	2 units
Murine IL2 standard (Lyophilized)	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

Function Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

Involvement in disease Note=A chromosomal aberration involving IL2 is found in a form of T-cell acute lymphoblastic leukemia (T-ALL). Translocation t(4;16)(q26;p13) with involves TNFRSF17.

Sequence similarities Belongs to the IL-2 family.

Cellular localization Secreted.

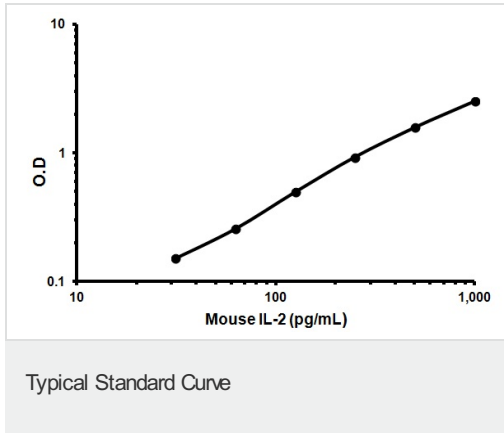
Applications

Our [Abpromise guarantee](#) covers the use of **ab46096** 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



Representative Standard Curve using ab46096.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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