

Product datasheet

Mouse IL-2 ELISA Set (without plates) ab47591

1 Image

Overview

Product name	Mouse IL-2 ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 5 pg/ml
Range	15.6 pg/ml - 500 pg/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse
Product overview	Abcam's IL-2 Murine <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Murine IL-2 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

This ELISA Set is intended to measure murine IL-2 in cell -culture supernatants, as well as in serum samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.

The murine IL-2 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 Murine IL-2

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-murine IL-2 antibody		1 x 0.55ml	2 x 0.55ml	3 x 0.55ml
Murine IL-2 Capture antibody		1 x 0.1ml	2 x 0.1ml	3 x 0.1ml
Murine IL-2 standard (Lyophilized)	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

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.

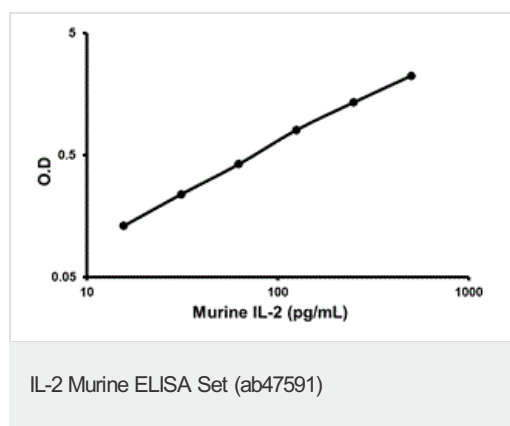
Applications

Our [Abpromise guarantee](#) covers the use of **ab47591** 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