

Mouse IL-2 ELISA Kit ab46096

3 References 1 图像

概述

产品名称

检测方法

精确度

小鼠IL-2 ELISA试剂盒

Colorimetric

批次内

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

批次间

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

样品类型

检测类型

灵敏度

范围

检测时间

实验步骤

种属反应性

产品概述

Cell culture supernatant, Serum

Sandwich (quantitative)

< 3.4 pg/ml

15.6 pg/ml - 500 pg/ml

3h 45m

Multiple steps standard assay

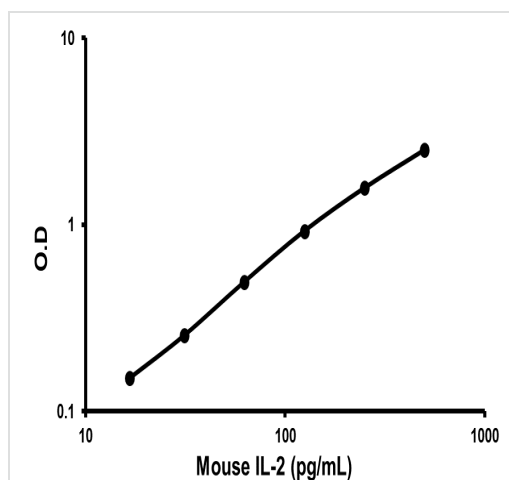
与反应: Mouse

Abcam's IL-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.

平台	Microplate		
性能			
存放说明	Store at +4°C. Please refer to protocols.		
组件	标识符	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
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-2 Biotinylated Detection antibody		1 x 400µl	2 x 400µl
IL-2 Microplate (12 x 8 well strips)		1 unit	2 units
Murine IL-2 standard (Lyophilized)	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl
功能	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.		
疾病相关	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.		
序列相似性	Belongs to the IL-2 family.		
细胞定位	Secreted.		
图片			



Representative Standard Curve using ab46096.

Typical Standard Curve

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors