

Human IL-2 ELISA Kit ab46033

1 图像

概述

产品名称	人IL-2 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Sample	6			3.4%
	批次间				
	样品	n	Mean	SD	CV%
	Sample	6			9.4%
样品类型	Cell culture supernatant, Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 10 pg/ml				
范围	31.2 pg/ml - 1000 pg/ml				
回收率	94 %				
检测时间	1h 50m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's IL-2 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for				

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 Human IL-2.

Get results in 90 minutes with Human IL-2 ELISA Kit ([ab174444](#)) from our SimpleStep ELISA® range.

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 anti IL-2	Red	1 x 400µl	2 x 400µl
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
Control	Silver	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-2 Microplate (12 x 8 well strips)		1 unit	2 units
IL-2 Standard (Lyophilized)	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

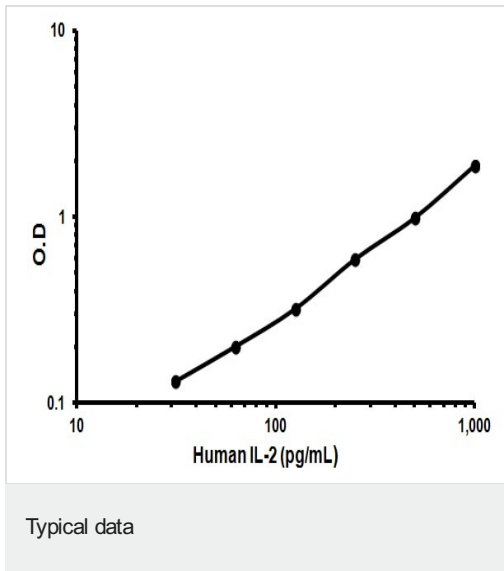
功能 Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

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 ab46033

**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