

Human IL-2 ELISA Kit ab100566

2 图像

概述

产品名称	人IL-2 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 4 pg/ml
范围	2.07 pg/ml - 1500 pg/ml
回收率	95 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	96.27	84% - 103%
Plasma	95.52	83% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Human IL-2 ELISA kit is designed for the quantitative measurement of Human IL-2 in plasma, and cell culture supernatants. (Human IL-2 concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for Human IL-2 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-2 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IL-2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Human IL-2 in normal serum with ab100566, only in serum spiked with Human IL-2.

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

range.

**说明** Optimization may be required with urine samples.

**平台** Microplate

## 性能

**存放说明** Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-2	2 vials
IL-2 Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-2 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

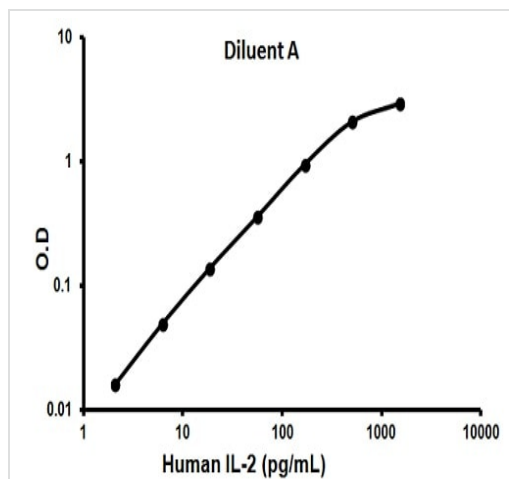
**功能** Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

**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.

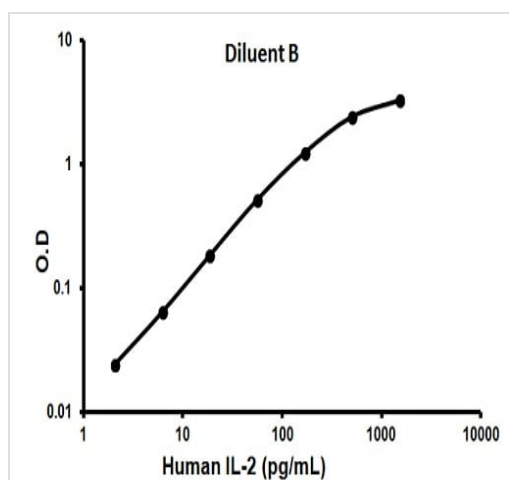
**细胞定位** Secreted.

## 图片



Representative standard curve using ab100566

Typical standard curve



Representative standard curve using ab100566

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