

# Human IL-2 ELISPOT Kit ab62926

## 概述

产品名称	人IL-2 ELISPOT试剂盒
检测方法	Colorimetric
样品类型	Suspension cells
检测类型	Sandwich (qualitative)
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>The ELISPOT assay is designed to enumerate cytokine producing cells in a single cell suspension. This method has the advantage of requiring a minimum of in-vitro manipulations allowing cytokine production analysis as close as possible to in-vivo conditions in a highly specific way. This technique is designed to determine the frequency of cytokine producing cells under a given stimulation, and the follow-up of such frequency during a treatment and/or a pathological state. Elispot assay constitutes an ideal tool in the TH1 / TH2 response, vaccine development, viral infection monitoring and treatment, cancerology, infectious diseases, autoimmune diseases and transplantation.</p> <p>Abcam Elispot assay is based on sandwich immuno-enzyme technology. Cell secreted cytokines or soluble molecules are captured by coated antibodies avoiding diffusion in supernatant, protease degradation or binding on soluble membrane receptors. After cell removal, the captured cytokines are revealed by tracer antibodies and appropriate conjugates.</p> <p><b>Principal:</b></p> <p>After cell stimulation, locally produced cytokines are captured by a specific monoclonal antibody. After cell lysis, trapped cytokine molecules are revealed by a secondary biotinylated detection antibody, which is in turn recognised by streptavidin conjugated to alkaline phosphatase. PVDF-bottomed-well plates are then incubated with BCIP/NBT substrate. Colored "purple" spots indicate cytokine production by individual cells.</p> <p>Recognizes natural human IL-2</p>
经测试应用	适用于: ELISpot
平台	Microplate

## 性能

## 存放说明

Store at +4°C. Please refer to protocols.

组件	1 x 96 tests	5 x 96 tests
Bovine Serum Albumin	1 x 200mg	1 x 1g
Human IL-2 precoated 96 well plates	1 unit	5 units
IL-2 Biotinylated detection antibody	1 vial	1 vial
Ready-to-use BCIP/NBT substrate buffer	1 x 10ml	1 x 50ml
Streptavidin - Alkaline Phosphatase conjugated	1 x 10µl	1 x 50µl

## 相关性

Human interleukin 2 (IL-2) is a secreted cytokine that is important for the proliferation of T and B lymphocytes. The receptor of this cytokine is a heterotrimeric protein complex whose gamma chain is also shared by interleukin 4 (IL-4) and interleukin 7 (IL-7). The expression of this gene in mature thymocytes is monoallelic, which represents an unusual regulatory mode for controlling the precise expression of a single gene. The targeted disruption of a similar gene in mice leads to ulcerative colitis like disease, which suggests an essential role of this gene in the immune response to antigenic stimuli. IL-2 has been shown to have antitumor effects in some studies. This is probably mediated by cytotoxic effector cells.

## 细胞定位

Secreted

## 应用

### The Abpromise guarantee

**Abpromise™**承诺保证使用ab62926于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
ELISpot		Use at an assay dependent dilution.

**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