

# Human IL2 ELISPOT Kit ab62926

## 製品の概要

---

<b>製品名</b>	Human IL2 ELISPOT Kit
<b>検出方法</b>	Colorimetric
<b>サンプルの種類</b>	Cell culture supernatant, Cell culture extracts
<b>アッセイタイプ</b>	Sandwich (qualitative)
<b>ステップ</b>	Multiple steps standard assay
<b>種交差性</b>	<b>交差種:</b> Human
<b>製品の概要</b>	<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 IL2</p>
<b>アプリケーション</b>	<b>適用あり:</b> ELISpot
<b>試験プラットフォーム</b>	Microplate

## 製品の特性

---

<b>保存方法</b>	Store at +4°C. Please refer to protocols.
-------------	---

内容	1 x 96 tests	5 x 96 tests
Bovine Serum Albumin	1 x 200mg	1 x 1g
Human IL2 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

**関連性** Human interleukin 2 (IL2) 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 (IL4) and interleukin 7 (IL7). 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. IL2 has been shown to have antitumor effects in some studies. This is probably mediated by cytotoxic effector cells.

**細胞内局在** Secreted

## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab62926** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
ELISpot		Use at an assay dependent dilution.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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 <http://www.abcam.co.jp/abpromise> or contact our technical team.

## Terms and conditions

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