

### Human IL-2 ELISPOT Kit (with un-coated plates) ab46554

#### 製品の概要

製品名	Human IL-2 ELISPOT Kit (with un-coated plates)
検出方法	Colorimetric
サンプルの種類	Suspension cells
アッセイタイプ	Sandwich (qualitative)
ステップ	Multiple steps standard assay
種交差性	交差種: Human, Monkey
製品の概要	<p>The Human IL-2 ELISPOT Kit (with un-coated plates) 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>This 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>
アプリケーション	適用あり: ELISpot
試験プラットフォーム	Microplate

## 製品の特徴

### 保存方法

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

内容	5 x 96 tests	10 x 96 tests
96 PVDF-bottomed-well plates.	5 units	10 units
Bovine Serum Albumin	1 x 2g	2 x 2g
Human IL-2 Capture antibody	1 x 500µl	2 x 500µl
IL-2 Biotinylated detection antibody	1 vial	2 vials
Ready-to-use BCIP/NBT substrate buffer	1 x 50ml	2 x 50ml
Streptavidin - Alkaline Phosphatase conjugated	1 x 50µl	2 x 50µl

### 関連性

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.

### 細胞内局在

Secreted

## アプリケーション

### The Abpromise guarantee

**Abpromise保証は、次のテスト済みアプリケーションにおけるab46554の使用に適用されます**

アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

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. 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.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