

# Human IL-17 ELISPOT Kit ab83723

## 製品の概要

製品名	Human IL-17 ELISPOT Kit
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

## 製品の概要

### Intended Use

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. 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.

### Principle of Method

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.

アプリケーション	適用あり: ELISpot
試験プラットフォーム	Microplate

## 製品の特性

保存方法	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 IL-17 Biotinylated detection antibody	1 vial	1 vial
Human IL-17 Pre-coated 96-well PDVF-bottomed plates	1 unit	5 units
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** IL-17, also known as CTLA8, is a T cell derived hematopoietic cytokine. It was originally cloned from a T cell hybridoma produced by fusion of a mouse cytotoxic T cell clone and a rat T lymphoma. IL-17 exhibits multiple biological activities on a variety of cells including: the induction of IL-6, IL-8 and GCSF production in fibroblasts; the enhancement of surface expression of ICAM 1 in fibroblasts; activation of NF kappaB and costimulation of T cell proliferation. IL-17 is an approximately 16 kDa polypeptide of 136 amino acids. The precursor form of IL-17 consists of 155 amino acids. To generate the mature IL-17 (136 amino acids), the precursor cleaves a 19 amino acid signal peptide. Human IL-17 shows approximately 62.5% amino acid homology to mouse IL-17 and 58% amino acid homology to rat IL-17.

**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ab83723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 concentration.

**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