

Human IL17 ELISPOT Kit ab83723

概述

产品名称	人IL17 ELISPOT试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Cell culture extracts
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

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.

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 IL17 Biotinylated detection antibody	1 vial	1 vial
Human IL17 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

相关性 IL17, 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. IL17 exhibits multiple biological activities on a variety of cells including: the induction of IL6, IL8 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. IL17 is an approximately 16 kDa polypeptide of 136 amino acids. The precursor form of IL17 consists of 155 amino acids. To generate the mature IL17 (136 amino acids), the precursor cleaves a 19 amino acid signal peptide. Human IL17 shows approximately 62.5% amino acid homology to mouse IL17 and 58% amino acid homology to rat IL17.

细胞定位 Secreted

应用

Our [Abpromise guarantee](#) covers the use of **ab83723** in the following tested applications.

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

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

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.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