

Human IL-17F ELISPOT Kit (with un-coated plates) ab105242

1 图像

概述

产品名称	人IL-17F ELISPOT试剂盒(with un-coated plates)
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
产品概述	<p>ab105242 is a highly specific immunoassay for the analysis of cytokine and other soluble molecule production and secretion from T-cells at a single cell level in conditions closely comparable to the in-vivo environment with minimal cell manipulation. This technique is designed to determine the frequency of cytokine producing cells under a given stimulation and the comparison of such frequency against a specific treatment or pathological state. The ELISPOT assay constitutes an ideal tool in the investigation of Th1 / Th2 responses, vaccine development, viral infection monitoring and treatment, oncology, infectious disease, autoimmune diseases and transplantation.</p> <p>Utilising sandwich immune-enzyme technology, ELISPOT assays can detect both secreted cytokines and single cells that simultaneously produce multiple cytokines. 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>A capture antibody highly specific for the molecule of interest is coated to the wells of a PVDF bottomed 96 well microtiter plate either during kit manufacture or in the laboratory. The plate is then blocked to minimise any non-antibody dependent unspecific binding and washed. Cell suspension and stimulant are added and the plate incubated allowing the specific antibodies to bind any molecules produced. Cells are then removed by washing prior to the addition of Biotinylated detection antibodies which bind to the previously captured molecule. Enzyme conjugated streptavidin is then added binding to the detection antibodies. Following incubation and washing substrate is then applied to the wells resulting in coloured spots which can be quantified using appropriate analysis software or manually using a microscope. See Figure 1 for visual scheme.</p> <p>The assay recognizes natural human IL-17F. After testing, no cross reactivity was observed for IL17A, IL17B, IL17D, IL17E, IL5, IL23 and Perforin. The antibody pair shows cross reactivity with the human IL17A + IL-17F heterodimer.</p>

说明

Store kit reagents between 4°C.

Uncoated plates should be stored at RT.

Immediately after use remaining reagents should be returned to cold storage (4°C).

All reagents should be warmed to room temperature before use.

BCIP/NBT buffer is potentially carcinogenic and should be disposed of appropriately, caution should be taken when handling this reagent, always wear gloves.

Capture Antibody is supplied sterile: once opened keep the vial sterile or aliquot and store at -20°C. For optimal performance prepare the Capture Antibody dilution immediately before use.

If not used within a short period of time, reconstituted Detection Antibody should be aliquoted and stored at -20°C. In these conditions the reagent is stable for at least one year.

For optimal performance prepare the Streptavidin-AP dilution immediately prior to use. Do not keep this solution for further experiments.

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	5 x 96 tests	10 x 96 tests
Bovine Serum Albumin	1 x 1g	2 x 1g
Human IL-17F Biotinylated detection antibody	1 vial	2 vials
Human IL-17F Capture antibody	1 x 500µl	2 x 500µl
Ready-to-use BCIP/NBT substrate buffer	1 x 50ml	2 x 50ml
Sterile 96 PVDF-bottomed-well plates	5 units	10 units
Streptavidin - Alkaline Phosphatase conjugated	1 x 50µl	2 x 50µl

功能

Ligand for IL17RA and IL17RC (PubMed:17911633). The heterodimer formed by IL17A and IL17F is a ligand for the heterodimeric complex formed by IL17RA and IL17RC (PubMed:18684971). Involved in stimulating the production of other cytokines such as IL6, IL8 and CSF2, and in regulation of cartilage matrix turnover (PubMed:11591732, PubMed:11591768, PubMed:11574464). Also involved in stimulating the proliferation of peripheral blood mononuclear cells and T-cells and in inhibition of angiogenesis (PubMed:11591732). Plays a role in the induction of neutrophilia in the lungs and in the exacerbation of antigen-induced pulmonary allergic inflammation.

组织特异性	Expressed in activated, but not resting, CD4+ T-cells and activated monocytes.
疾病相关	Candidiasis, familial, 6
序列相似性	Belongs to the IL-17 family.
细胞定位	Secreted.

应用

The Abpromise guarantee

Abpromise™承诺保证使用ab105242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.

图片

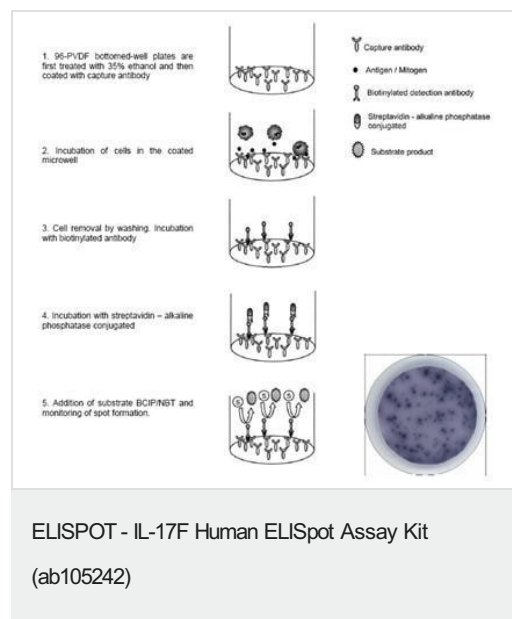


Figure 1: principle of assay

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