

Human IL12 p70 ELISPOT Kit (without plates) ab46552

概述

产品名称	人IL12 p70 ELISPOT试剂盒(without plates)
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	<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>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>Principal:</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 IL12 Specific to p70 heterodimer</p>
经测试应用	适用于: ELISpot
平台	Reagents

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	5 x 96 tests	10 x 96 tests	15 x 96 tests	20 x 96 tests
Bovine Serum Albumin	1 x 1g	2 x 1g	3 x 1g	4 x 1g
Dry Skimmed milk	1 vial	2 vials	3 vials	4 vials
IL12 p70 Biotinylated Detection Antibody	1 x 0.55ml	2 x 0.55ml	3 x 0.55ml	4 x 0.55ml
IL12 p70 Capture antibody (clone B-T21)	1 x 0.5ml	2 x 0.5ml	3 x 0.5ml	4 x 0.5ml
Ready-to-use BCIP/NBT substrate buffer	1 x 50ml	2 x 50ml	3 x 50ml	4 x 50ml
Streptavidin - Alkaline Phosphatase conjugated	1 x 50µl	2 x 50µl	3 x 50µl	4 x 50µl

相关性 Interleukin 12 is a cytokine that acts on T and natural killer cells, and has a broad array of biological activities. Interleukin 12 is a disulfide-linked heterodimer composed of the 35 kD cytokine receptor like subunit encoded by IL12A gene, and a 40 kD subunit encoded by IL12B. This cytokine is expressed by activated macrophages that serve as an essential inducer of Th1 cells development. It is required for the T-cell-independent induction of interferon (IFN)-gamma, and is important for the differentiation of both Th1 and Th2 cells. This cytokine has been found to be important for sustaining a sufficient number of memory/effector Th1 cells to mediate long-term protection to an intracellular pathogen. The responses of lymphocytes to this cytokine are mediated by the activator of transcription protein STAT4. Nitric oxide synthase 2A (NOS2A/NOS2) is found to be required for the signaling process of this cytokine in innate immunity. Overexpression of this gene was observed in the central nervous system of patients with multiple sclerosis (MS), suggesting a role of this cytokine in the pathogenesis of the disease. The promoter polymorphism of this gene has been reported to be associated with the severity of atopic and non-atopic asthma in children.

细胞定位 Secreted

应用

Our [Abpromise guarantee](#) covers the use of **ab46552** 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 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 <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