

Human IL-12 p70 ELISPOT Kit ab62939

概述

| | |
|-------|---|
| 产品名称 | 人IL-12 p70 ELISPOT试剂盒 |
| 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 |
| 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>Abcam 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 IL-12, specific to p70 heterodimer.</p> |
| 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 0.2g | 1 x 1g |
| IL-12 p70 Biotinylated Detection Antibody | 1 x 0.1ml | 1 x 0.55ml |
| Human IL-12 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

IL12A is a subunit of a cytokine that acts on T and natural killer cells, and has a broad array of biological activities. The cytokine is a disulfide-linked heterodimer composed of the 35-kD subunit encoded by this gene, and a 40-kD subunit that is a member of the cytokine receptor family. This cytokine is required for the T-cell-independent induction of interferon (IFN)-gamma, and is important for the differentiation of both Th1 and Th2 cells. 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. IL12B is a cytokine that can act as a growth factor for activated T and NK cells, enhance the lytic activity of NK/lymphokine-activated killer cells, and stimulate the production of IFN-gamma by resting PBMC.

细胞定位

Secreted

应用

The Abpromise guarantee

Abpromise™ 承诺保证使用ab62939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. |

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