

Human Perforin ELISPOT Kit ab62943

概述

产品名称	人Perforin 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>This ELISPOT assay is designed to enumerate Perforin producing cells in a single cell suspension. This method has the advantage of requiring a minimum of in-vitro manipulations allowing perforin production analysis as close as possible to in-vivo conditions in a highly specific way. This technique is designed to determine the frequency of perforin 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 or molecule are captured by a specific monoclonal antibody. After cell lysis, trapped specific soluble 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 Perforin</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 1g	1 x 1g
Human Perforin Pre-coated 96-well PDVF-bottomed plates	1 unit	1 x 5 units
Perforin Biotinylated Detection Antibody	1 vial	1 vial
Ready-to-use BCIP/NBT substrate buffer	1 x 27ml	1 x 50ml
Streptavidin - Alkaline Phosphatase conjugated	1 x 10µl	1 x 50µl

相关性

Perforin has structural and functional similarities to complement component 9 (C9). Like C9, this protein creates transmembrane tubules and is capable of lysing non-specifically a variety of target cells. Perforin is one of the main cytolytic proteins of cytolytic granules, and it is known to be a key effector molecule for T-cell- and natural killer-cell-mediated cytotoxicity. Defects in this gene cause familial hemophagocytic lymphohistiocytosis type 2 (HPLH2), a rare and lethal autosomal recessive disorder of early childhood.

细胞定位

Cell Membrane and Cytoplasmic

应用

The Abpromise guarantee

Abpromise™承诺保证使用ab62943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