

Human Interferon Gamma + IL2 FLUOROSPOT Kit ab48452

1 References

概述

产品名称	人Interferon Gamma + IL2 FLUOROSPOT试剂盒
检测方法	Fluorescent
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, oncology, 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>
经测试应用	适用于: FluroSpot
平台	Microplate

性能

存放说明 Please refer to protocols.

组件	1 x 96 tests	5 x 96 tests	10 x 96 tests
96 well PVDF bottomed Microplate (non-sterile)	1 unit	5 units	10 units
Anti-FITC antibody green fluorescence conjugate.	1 vial	1 vial	2 vials
Biotinylated detection antibody for IL-2 (resuspend in 0.55mL).	1 vial	1 vial	2 vials

组件	1 x 96 tests	5 x 96 tests	10 x 96 tests
Bovine Serum Albumin	1 x 2g	1 x 2g	2 x 2g
FITC conjugated detection antibody for IFN γ (resuspend in 0.55mL)	1 vial	1 vial	2 vials
Human IFN Gamma Capture antibody	1 x 100 μ l	1 x 500 μ l	2 x 500 μ l
Human IL-2 Capture antibody	1 x 100 μ l	1 x 500 μ l	2 x 500 μ l
Streptavidin-phycoerythrin conjugate.	1 vial	1 vial	2 vials

相关性 IFN-gamma is produced by lymphocytes activated by specific antigens or mitogens. IFN-gamma, in addition to having antiviral activity, has important immunoregulatory functions. It is a potent activator of macrophages, it has antiproliferative effects on transformed cells and it can potentiate the antiviral and antitumor effects of the type I interferons.

细胞定位 Secreted

应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab48452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
FluroSpot		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