

Product datasheet

Human Interferon Gamma + Granzyme B ELISPOT Kit (without plates) ab48463

製品の概要

製品名 Human Interferon Gamma + Granzyme B ELISPOT Kit (without plates)

サンプルの種類 Cell culture supernatant, Cell culture extracts

アッセイタイプ Sandwich (qualitative)

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

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.

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.

The dual colour Elispot allows you to monitor the production of two cytokines simultaneously in the same well.

Principle of the Method

After cell stimulation, locally produced cytokines are captured by IFN gamma and Granzyme B specific monoclonal antibodies. After cell lysis, trapped cytokine molecules are revealed by a secondary anti-IFN gamma FITC conjugated antibody and a biotinylated anti-Granzyme B antibody. Those are in turn recognised by anti-FITC HRP and streptavidin-AP conjugates. PVDF-bottomed-well plates are then incubated first with AEC substrate buffer, washed and subsequently incubated with BCIP/NBT. Coloured red/brownish spots indicate IFN gamma production while Granzyme B is revealed by blue/purple spots.

アプリケーション 適用あり: ELISpot

試験プラットフォーム Reagents

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	5 x 96 tests	10 x 96 tests	15 x 96 tests	20 x 96 tests
10 x concentrate buffer for the preparation of AEC buffer	1 x 5ml	2 x 5ml	3 x 5ml	4 x 5ml
50 x concentrate AEC substrate buffer	1 x 1ml	2 x 1ml	3 x 1ml	4 x 1ml
Anti-FITC antibody HRP conjugate	1 x 100µl	2 x 100µl	3 x 100µl	4 x 100µl
Bovine Serum Albumin	1 x 1g	2 x 1g	3 x 1g	4 x 1g
Dry Skimmed milk	1 x 1g	2 x 1g	3 x 1g	4 x 1g
Granzyme B Biotinylated detection antibody	1 vial	2 vials	3 vials	4 vials
Granzyme B Capture Antibody	1 x 500µl	2 x 500µl	3 x 500µl	4 x 500µl
Human IFN γ Capture antibody	1 x 500µl	2 x 500µl	3 x 500µl	4 x 500µl
IFN γ FITC conjugated detection antibody	1 vial	2 vials	3 vials	4 vials
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

関連性

Interferon 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. Granzyme B is necessary for target cell lysis in cell-mediated immune responses. It cleaves after Asp. Seems to be linked to an activation cascade of caspases (aspartate-specific cysteine proteases) responsible for apoptosis execution. Cleaves caspase-3, -7, -9 and 10 to give rise to active enzymes mediating apoptosis.

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab48463** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
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 <http://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors