

Human IL10 + IL2 ELISPOT Kit (without plates) ab48455

製品の概要

製品名 Human IL10 + IL2 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, cancerology, 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

After cell stimulation, locally produced cytokines are captured by IL10 and IL2 specific monoclonal antibodies. After cell lysis, trapped cytokine molecules are revealed by a secondary anti-IL10 FITC conjugated antibody and a biotinylated anti-IL2 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 IL10 production while IL2 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 |
| Human IL-10 Capture antibody | 1 x 500µl | 2 x 500µl | 3 x 500µl | 4 x 500µl |
| Human IL-2 Capture antibody | 1 x 500µl | 2 x 500µl | 3 x 500µl | 4 x 500µl |
| IL10 FITC conjugated detection antibody | 1 vial | 2 vials | 3 vials | 4 vials |
| IL-2 Biotinylated 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 IL2 is produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells. IL10 inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

細胞内局在 Secreted

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab48455** 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