

Product datasheet

Human Interferon Gamma + IL2 ELISPOT Kit ab48448

Overview

Product name	Human Interferon Gamma + IL2 ELISPOT Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Cell culture extracts
Assay type	Sandwich (qualitative)
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human

Product overview

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.

Tested applications	Suitable for: ELISpot
Platform	Microplate

Properties

Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	5 x 96 tests	10 x 96 tests
10 x concentrate buffer for the preparation of AEC buffer	1 x 5ml	2 x 5ml
50 x concentrate AEC substrate buffer	1 x 1ml	2 x 1ml

Components	5 x 96 tests	10 x 96 tests
96 PVDF-bottomed-well plates.	5 units	10 units
Anti-FITC antibody HRP conjugate	1 x 120µl	2 x 100µl
Bovine Serum Albumin	1 x 1g	2 x 1g
Human IFN γ Capture antibody	1 x 500µl	2 x 500µl
Human IL-2 Capture antibody	1 x 500µl	2 x 500µl
IFN γ FITC conjugated detection antibody	1 vial	2 vials
IL-2 Biotinylated detection antibody	1 vial	2 vials
Ready-to-use BCIP/NBT substrate buffer	2 x 27ml	2 x 50ml
Streptavidin - Alkaline Phosphatase conjugated	1 x 50µl	2 x 50µl

Relevance 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.

Cellular localization Secreted

Applications

Our [Abpromise guarantee](#) covers the use of **ab48448** in the following tested applications.

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

Application	Abreviews	Notes
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.com/abpromise> or contact our technical team.

Terms and conditions

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