

Human IL-6 ELISPOT Kit ab62936

Overview

Product name Human IL-6 ELISPOT Kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|----|------|----|-------|
| Overall | 12 | | | 5.9% |
| Overall | 12 | | | 3.9% |
| Overall | 12 | | | 4.7% |
| Overall | 12 | | | 6.3% |
| Overall | 12 | | | 10.4% |

Sample type Suspension cells

Assay type Sandwich (qualitative)

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human IL-6 ELISpot (ab62936) is a highly specific immunoassay for the analysis of IL-6 production and secretion from T-cells at a single cell level in conditions closely comparable to the in-vivo environment with minimal cell manipulation.

A capture antibody highly specific for the analyte of interest is coated to the wells of a PVDF bottomed 96 well microtiter plate either during kit manufacture or in the laboratory. The plate is then blocked to minimise any non-antibody dependent unspecific binding and washed. Cell suspension and stimulant are added and the plate incubated allowing the specific antibodies to bind any analytes produced. Cells are then removed by washing prior to the addition of Biotinylated detection antibodies which bind to the previously captured analyte. Enzyme conjugated streptavidin is then added binding to the detection antibodies. Following incubation and washing substrate is then applied to the wells resulting in coloured spots which can be quantified using appropriate analysis software or manually using a microscope.

For the quantitative determination of the frequency of Human IL-6 -producing cells.

For research use only - not intended for diagnostic use.

Tested applications **Suitable for:** ELISpot
Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

| Components | 1 x 96 tests |
|--|--------------|
| Biotinylated detection antibody, clone B-E4: | 1 x 0.55ml |
| Bovine Serum Albumin | 1 x 2g |
| Human IL-6 Pre-coated 96-well PDVF-bottomed plates | 1 unit |
| Ready-to-use BCIP/NBT substrate buffer | 1 x 50ml |
| Streptavidin - Alkaline Phosphatase conjugated | 1 x 50µl |

Function Cytokine with a wide variety of biological functions. It is a potent inducer of the acute phase response. Plays an essential role in the final differentiation of B-cells into Ig-secreting cells Involved in lymphocyte and monocyte differentiation. It induces myeloma and plasmacytoma growth and induces nerve cells differentiation Acts on B-cells, T-cells, hepatocytes, hematopoietic progenitor cells and cells of the CNS. Also acts as a myokine. It is discharged into the bloodstream after muscle contraction and acts to increase the breakdown of fats and to improve insulin resistance.

Involvement in disease Genetic variations in IL6 are associated with susceptibility to rheumatoid arthritis systemic juvenile (RASJ) [MIM:604302]. An inflammatory articular disorder with systemic-onset beginning before the age of 16. It represents a subgroup of juvenile arthritis associated with severe extraarticular features and occasionally fatal complications. During active phases of the disorder, patients display a typical daily spiking fever, an evanescent macular rash, lymphadenopathy, hepatosplenomegaly, serositis, myalgia and arthritis.
Note=A IL6 promoter polymorphism is associated with a lifetime risk of development of Kaposi sarcoma in HIV-infected men.

Sequence similarities Belongs to the IL-6 superfamily.

Post-translational modifications N- and O-glycosylated.

Cellular localization Secreted.

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab62936 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. 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.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