

Human IL-6 ELISPOT Kit ab62936

概述

产品名称 人IL-6 ELISPOT试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall	12			5.9%
Overall	12			3.9%
Overall	12			4.7%
Overall	12			6.3%
Overall	12			10.4%

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 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.

经测试应用

适用于: ELISpot

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

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

组件	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

功能

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.

疾病相关

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.

序列相似性

Belongs to the IL-6 superfamily.

翻译后修饰

N- and O-glycosylated.

细胞定位

Secreted.

应用

The Abpromise guarantee

Abpromise™ 承诺保证使用ab62936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

应用	Ab评论	说明
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.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