

Human IL-21 ELISA Kit ab119542

5 References 1 图像

概述

产品名称

检测方法

精确度

样品类型

检测类型

灵敏度

范围

回收率

实验步骤

种属反应性

产品概述

人IL-21 ELISA试剂盒

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall	8			5.8%

批次间

样品	n	Mean	SD	CV%
Overall	8			7.7%

Cell culture supernatant, Serum, Plasma

Sandwich (quantitative)

20 pg/ml

78 pg/ml - 5000 pg/ml

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	95	% - %
Serum	71	% - %

Multiple steps standard assay

与反应: Human

Abcam's IL-21 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human IL-21 concentrations in Cell culture supernatant, serum and plasma (EDTA, citrate).

IL-21 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated IL-21 detection antibody then incubated at room temperature. Following washing, a Streptavidin-HRP conjugate is added to each well,

incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of IL-21 captured on the plate.

**平台** Microplate

## 性能

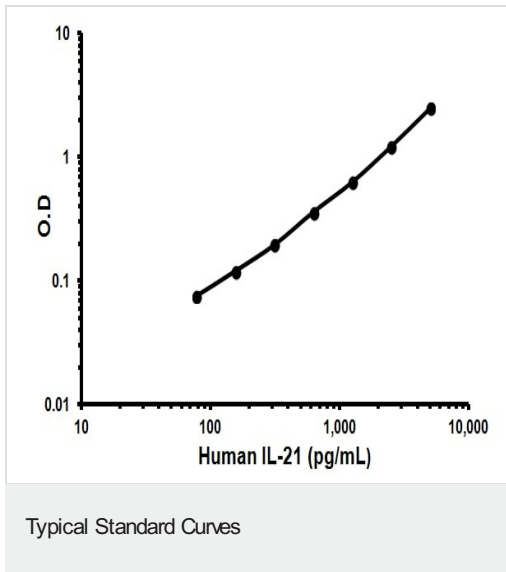
**存放说明** Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-human IL-21 monoclonal antibody	1 x 70µl
Human IL-21 Standard lyophilized (10 ng/ml upon reconstitution)	2 vials
Microplate coated with monoclonal antibody to Human IL-21	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

**相关性** A novel cytokine related to IL-2 and IL-15 was recently identified and designated IL-21. IL-21 has been found to be a powerful growth factor for naive B cells. The receptor for IL-21 (IL-21R, also termed NILR for novel Interleukin receptor) is a new member of the class I cytokine receptor family. IL-21R forms a complex with the common cytokine receptor  $\gamma$  chain,  $\gamma c$ , and mediates IL-21 signaling. Both IL-21R and the  $\gamma c$  are necessary for the IL-21 function. IL-21 and its receptor activate JAK/STAT signaling pathway. IL-21R is expressed in spleen, thymus, natural killer (NK), T and B cell lines. IL-21 plays a role in the proliferation and maturation of NK, B and T cell populations.

**细胞定位** Secreted

## 图片



Representative Standard Curve using ab119542.

**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