

Product datasheet

Human IL-21 ELISA Kit (Interleukin-21) ab119542

画像数 1

製品の概要

製品名 Human IL-21 ELISA Kit (Interleukin-21)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall	8			5.8%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	8			7.7%

サンプルの種類 Cell culture supernatant, Serum, Plasma

アッセイタイプ Sandwich (quantitative)

検出感度 20 pg/ml

検出範囲 78 pg/ml - 5000 pg/ml

添加回収試験

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	95	% - %
Serum	71	% - %

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Abcam's IL-21 (Interleukin-21) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human IL-21 (Interleukin-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.

アプリケーション

適用あり: Sandwich ELISA

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	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 IL21 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

関連性

A novel cytokine related to IL2 and IL15 was recently identified and designated IL21. IL21 has been found to be a powerful growth factor for naive B cells. The receptor for IL21 (IL21R, also termed NILR for novel Interleukin receptor) is a new member of the class I cytokine receptor family. IL21R forms a complex with the common cytokine receptor γ chain, γc , and mediates IL21 signaling. Both IL21R and the γc are necessary for the IL21 function. IL21 and its receptor activate JAK/STAT signaling pathway. IL21R is expressed in spleen, thymus, natural killer (NK), T and B cell lines. IL21 plays a role in the proliferation and maturation of NK, B and T cell populations.

細胞内局在

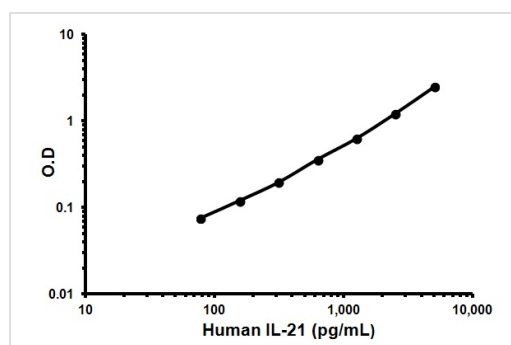
Secreted

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab119542** 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
Sandwich ELISA		Use at an assay dependent concentration.



Representative Standard Curve using ab119542.

Typical Standard Curves

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