

Human IL-2 ELISA Kit High Sensitivity ab46054

3 References 画像数 1

製品の概要

| | | | | | |
|----------|---|---|------|-------|------|
| 製品名 | Human IL-2 ELISA Kit High Sensitivity | | | | |
| 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 | 5 | | | 4.2% |
| | Inter-Assay (日差再現性) | | | | |
| | サンプル | N | 平均値 | SD | CV% |
| | Overall | 5 | | | 9% |
| 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 | < 0.97 pg/ml | | | | |
| 検出範囲 | 1.87 pg/ml - 60 pg/ml | | | | |
| 添加回収試験 | 100 % | | | | |
| | 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 | | 100 | % - % | |
| 全工程の試験時間 | 2h 00m | | | | |
| ステップ | Multiple steps standard assay | | | | |
| 種交差性 | 交差種: Human | | | | |
| 製品の概要 | Abcam's Human High Sensitivity IL-2 ELISA is to be used for the <i>in-vitro</i> quantitative determination interleukin-2 (IL-2) in Human sera, plasmas, buffered solutions or cell culture media. | | | | |

A monoclonal antibody specific for IL-2 has been coated onto the wells of the microtiter strips

provided. Samples, including standards of known IL-2 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-2 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-2 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-2.

Get results in 90 minutes with Human IL-2 ELISA Kit ([ab174444](#)) from our SimpleStep ELISA® range.

試験プラットフォーム

Microplate

製品の特長

保存方法

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

| 内容 | ラベル | 1 x 96 tests | 1 x 96 tests |
|--------------------------------------|-------------------|--------------|--------------|
| 10X Standard Diluent Buffer | Black | 1 x 15ml | 1 x 15ml |
| 200X Wash Buffer | White | 1 x 10ml | 1 x 10ml |
| Amplification Diluent | Brown & blue spot | 1 x 15ml | 1 x 15ml |
| Amplifier | Yellow | 1 x 200µl | 1 x 200µl |
| Biotinylated anti IL-2 | Red | 1 x 400µl | 1 x 400µl |
| Biotinylated Antibody Diluent | Red | 1 x 7.5ml | 1 x 7.5ml |
| Chromogen TMB Substrate Solution | | 1 x 11ml | 1 x 11ml |
| HRP Diluent | Red | 1 x 23ml | 1 x 23ml |
| IL-2 Microplate (12 x 8 well strips) | | 1 unit | 1 unit |
| IL-2 Standard (Lyophilized) | Yellow | 2 vials | 2 vials |
| Standard Diluent (Serum) | | 1 x 7ml | 1 x 7ml |
| Stop Reagent | Black | 1 x 11ml | 1 x 11ml |
| Streptavidin-HRP | | 2 x 5µl | 2 x 5µl |

機能

Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

関連疾患

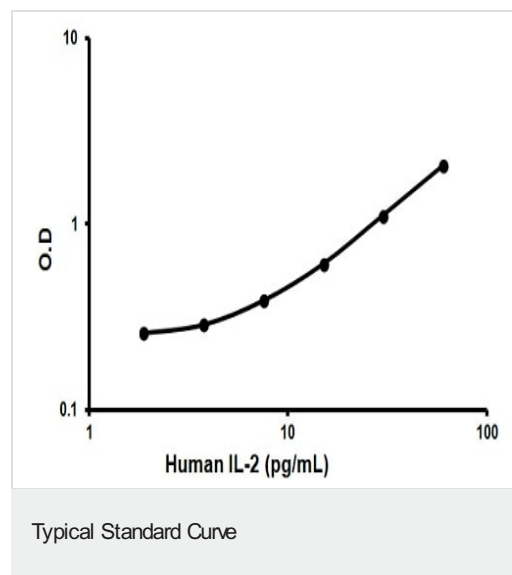
Note=A chromosomal aberration involving IL2 is found in a form of T-cell acute lymphoblastic leukemia (T-ALL). Translocation t(4;16)(q26;p13) with involves TNFRSF17.

配列類似性

Belongs to the IL-2 family.

細胞内局在

Secreted.

画像

Representative standard curve using ab46054

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.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