

Human IL-2 ELISA Kit ab46033

画像数 1

製品の概要

| 製品名 | Human IL-2 ELISA Kit | | | | | | | | | | |
|----------|---|------|----|------|----|-----|--------|---|--|--|------|
| 検出方法 | Colorimetric | | | | | | | | | | |
| 再現性 | Intra-Assay (同時再現性) | | | | | | | | | | |
| | <table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Sample</td> <td>6</td> <td></td> <td></td> <td>3.4%</td> </tr> </tbody> </table> | サンプル | N | 平均値 | SD | CV% | Sample | 6 | | | 3.4% |
| サンプル | N | 平均値 | SD | CV% | | | | | | | |
| Sample | 6 | | | 3.4% | | | | | | | |
| | Inter-Assay (日差再現性) | | | | | | | | | | |
| | <table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Sample</td> <td>6</td> <td></td> <td></td> <td>9.4%</td> </tr> </tbody> </table> | サンプル | N | 平均値 | SD | CV% | Sample | 6 | | | 9.4% |
| サンプル | N | 平均値 | SD | CV% | | | | | | | |
| Sample | 6 | | | 9.4% | | | | | | | |
| 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 | < 10 pg/ml | | | | | | | | | | |
| 検出範囲 | 31.2 pg/ml - 1000 pg/ml | | | | | | | | | | |
| 添加回収試験 | 94 % | | | | | | | | | | |
| 全工程の試験時間 | 1h 50m | | | | | | | | | | |
| ステップ | Multiple steps standard assay | | | | | | | | | | |
| 種交差性 | 交差種: Human | | | | | | | | | | |
| 製品の概要 | <p>Abcam's IL-2 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-2 in supernatants, buffered solutions, serum and plasma samples.</p> <p>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.</p> | | | | | | | | | | |

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

製品の特性

保存方法

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

| 内容 | ラベル | 1 x 96 tests | 2 x 96 tests |
|--------------------------------------|--------|--------------|--------------|
| 10X Standard Diluent Buffer | Black | 1 x 15ml | 1 x 25ml |
| 200X Wash Buffer | White | 1 x 10ml | 2 x 10ml |
| Biotinylated anti IL-2 | Red | 1 x 400µl | 2 x 400µl |
| Biotinylated Antibody Diluent | Red | 1 x 7ml | 1 x 13ml |
| Chromogen TMB Substrate Solution | | 1 x 11ml | 1 x 24ml |
| Control | Silver | 2 vials | 4 vials |
| HRP Diluent | Red | 1 x 12ml | 1 x 23ml |
| IL-2 Microplate (12 x 8 well strips) | | 1 unit | 2 units |
| IL-2 Standard (Lyophilized) | Yellow | 2 vials | 4 vials |
| Standard Diluent (Serum) | | 1 x 7ml | 2 x 7ml |
| Stop Reagent | Black | 1 x 11ml | 2 x 11ml |
| Streptavidin-HRP | | 2 x 5µl | 4 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.

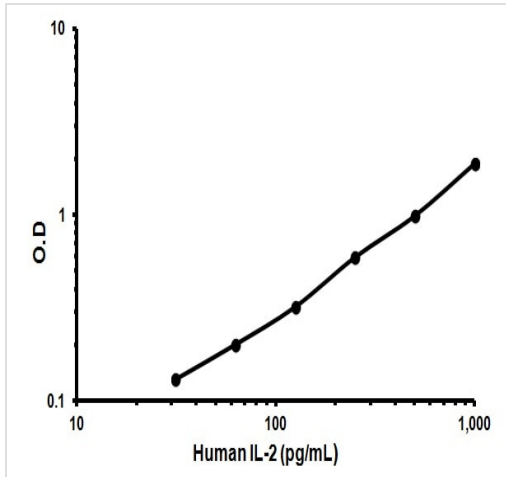
配列類似性

Belongs to the IL-2 family.

細胞内局在

Secreted.

画像



Representative standard curve using ab46033

Typical data

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