

Human IL-2 ELISA Set (without plates) ab48471

画像数 1

製品の概要

製品名	Human IL-2 ELISA Set (without plates)
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	31.25 pg/ml - 1000 pg/ml
全工程の試験時間	1h 45m
ステップ	Multiple steps standard assay
種交差性	交差種: Human

製品の概要

Abcam's IL-2 Human *in vitro* ELISA Set is designed for the quantitative measurement of IL-2 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

The IL-2 ELISA set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.

This ELISA set is intended to measure human IL-2 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.

アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	5 x 96 tests	10 x 96 tests	15 x 96 tests
Detection biotinylated anti-IL-2 antibody		1 x 0.55ml	2 x 0.55ml	3 x 0.55ml
IL-2 Capture antibody		1 x 0.5ml	2 x 0.5ml	3 x 0.5ml

内容	ラベル	5 x 96 tests	10 x 96 tests	15 x 96 tests
IL-2 Standard (Lyophilized)	Yellow	5 vials	10 vials	15 vials
Ready-to-Use TMB		2 vials	4 vials	6 vials
Streptavidin-HRP		1 vial	2 vials	3 vials

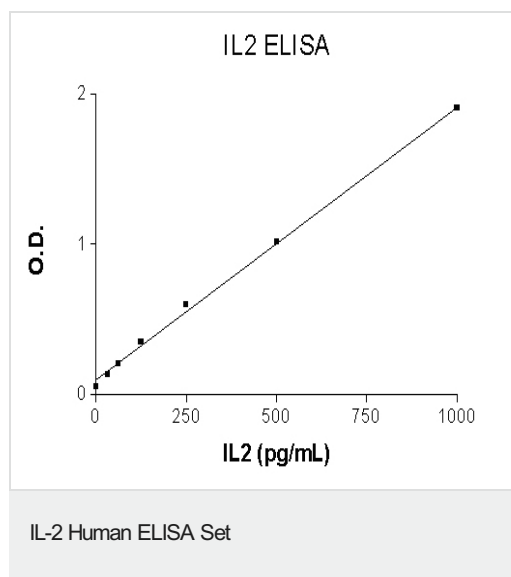
機能	Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.
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

アプリケーション

The Abpromise guarantee **Abpromise保証は、次のテスト済みアプリケーションにおけるab48471の使用に適用されます**
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



IL-2 Standard curve ranging from 31.25 to 1000 pg/mL

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