

Mouse IL-2 ELISA Set (without plates) ab47591

2 References 画像数 1

製品の概要

製品名	Mouse 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	< 5 pg/ml
検出範囲	15.6 pg/ml - 500 pg/ml
全工程の試験時間	3h 45m
ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	<p>Abcam's IL-2 Murine <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Murine IL-2 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure murine IL-2 in cell -culture supernatants, as well as in serum samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>The murine 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.</p> <p>Recognizes both natural and recombinant Murine IL-2</p>
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	10 x 96 tests	5 x 96 tests	15 x 96 tests
Chromogen TMB Substrate Solution		4 vials	2 vials	6 vials

内容	ラベル	10 x 96 tests	5 x 96 tests	15 x 96 tests
Detection biotinylated anti-murine IL-2 antibody		2 x 0.55ml	1 x 0.55ml	3 x 0.55ml
Murine IL-2 Capture antibody		2 x 0.1ml	1 x 0.1ml	3 x 0.1ml
Murine IL-2 standard (Lyophilized)	Yellow	10 vials	5 vials	15 vials
Streptavidin-HRP		2 vials	1 vial	3 vials

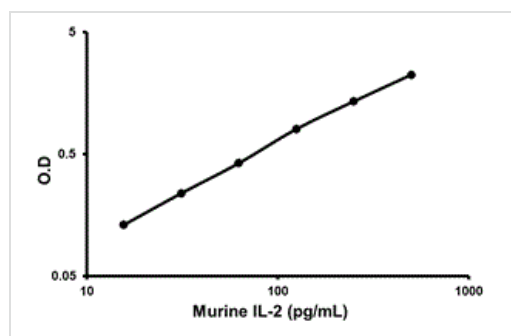
機能	Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.
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ab47591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 Murine ELISA Set (ab47591)

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