

Product datasheet

Human TNF alpha ELISA Set (without plates) ab47347

画像数 1

製品の概要

製品名	Human TNF alpha 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	25 pg/ml - 800 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human

製品の概要
 Abcam's TNF α Human *in vitro* ELISA Set is designed for the quantitative measurement of TNF α in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

The TNF alpha Elipair 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 Elipair set is intended to measure human TNF alpha 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.

Recognizes both natural and recombinant human TNF alpha.

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	5 x 96 tests	10 x 96 tests
Detection Biotinylated anti-TNF alpha Antibody	1 x 0.55ml	2 x 0.55ml
Ready-to-use TMB	2 vials	4 vials

Streptavidin-HRP	1 vial	2 vials
TNF alpha Capture Antibody	1 x 0.5ml	2 x 0.5ml
TNF alpha Standard	5 vials	10 vials

関連性 TNF-alpha is a cytokine that binds to TNFRSF1A/TNFR1 and TNFRSF1B/TNFR. It is mainly secreted by macrophages and can induce cell death of certain tumor cell lines. TNF-alpha is a potent pyrogen causing fever by direct action or by stimulation of interleukin 1 secretion and is implicated in the induction of cachexia, under certain conditions it can stimulate cell proliferation and induce cell differentiation.

細胞内局在 Cell membrane; single-pass type II membrane protein. Processed form: Secreted protein.

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab47347** 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
----------	-----------	------

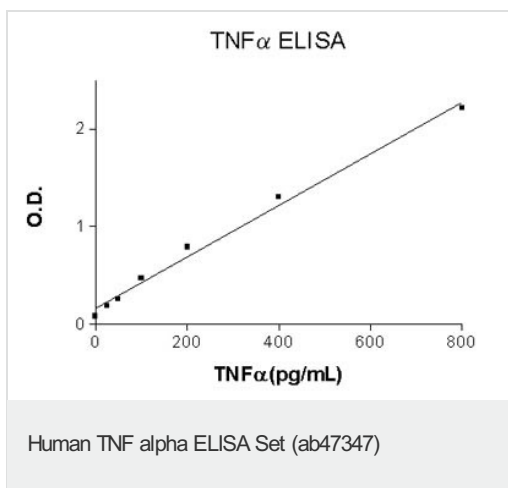
Sandwich ELISA

追加情報 ELISA: Use at an assay dependent dilution.

Not yet tested in other applications.

Optimal dilutions/concentrations should be determined by the end user.

Human TNF alpha ELISA Set (without plates) 画像



TNF alpha Standard curve ranging from 25 to 800 pg/mL

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