

Product datasheet

Human TNF alpha ELISA Set (without plates) ab47347

1 图像

概述

产品名称	人TNF alpha ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 10 pg/ml
范围	25 pg/ml - 800 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's TNFα Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of TNFα in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>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.</p> <p>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.</p> <p>Recognizes both natural and recombinant human TNF alpha.</p>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	5 x 96 tests	10 x 96 tests
Detection Biotinylated anti-TNF alpha Antibody	1 x 0.55ml	2 x 0.55ml

组件	5 x 96 tests	10 x 96 tests
Chromogen TMB Substrate Solution	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

相关性 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.

细胞定位 Cell membrane; single-pass type II membrane protein. Processed form: Secreted protein.

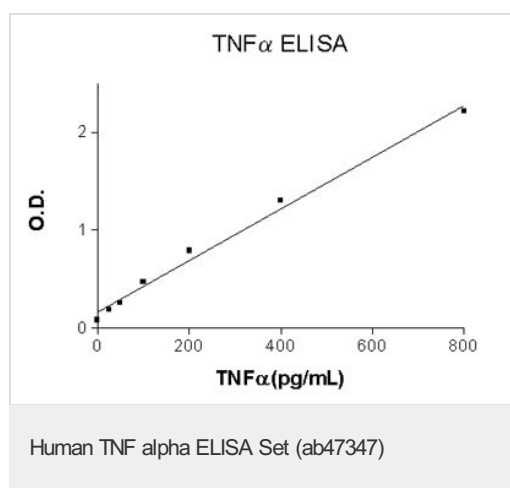
应用

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.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors