

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
检测时间	2h 45m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's TNF<math>\alpha</math> Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of TNF<math>\alpha</math> 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	10 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution	4 vials	2 vials

组件	10 x 96 tests	5 x 96 tests
Detection Biotinylated anti-TNF alpha Antibody	2 x 0.55ml	1 x 0.55ml
Streptavidin-HRP	2 vials	1 vial
TNF alpha Capture Antibody	2 x 0.5ml	1 x 0.5ml
TNF alpha Standard	10 vials	5 vials

功能	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. It is 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.
疾病相关	Genetic variations in TNF are a cause of susceptibility psoriatic arthritis (PSORAS) [MIM:607507]. PSORAS is an inflammatory, seronegative arthritis associated with psoriasis. It is a heterogeneous disorder ranging from a mild, non-destructive disease to a severe, progressive, erosive arthropathy. Five types of psoriatic arthritis have been defined: asymmetrical oligoarthritis characterized by primary involvement of the small joints of the fingers or toes; asymmetrical arthritis which involves the joints of the extremities; symmetrical polyarthritis characterized by a rheumatoidlike pattern that can involve hands, wrists, ankles, and feet; arthritis mutilans, which is a rare but deforming and destructive condition; arthritis of the sacroiliac joints and spine (psoriatic spondylitis).
序列相似性	Belongs to the tumor necrosis factor family.
翻译后修饰	The soluble form derives from the membrane form by proteolytic processing. The membrane form, but not the soluble form, is phosphorylated on serine residues. Dephosphorylation of the membrane form occurs by binding to soluble TNFRSF1A/TNFR1. O-glycosylated; glycans contain galactose, N-acetylgalactosamine and N-acetylneuraminic acid.
细胞定位	Secreted and Cell membrane.

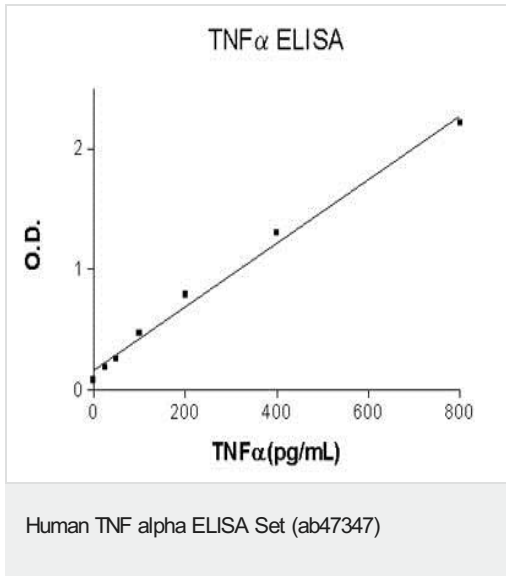
应用

**The Abpromise guarantee**      **Abpromise™**承诺保证使用ab47347于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

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. 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.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