

Human TGF alpha ELISA Kit ab100646

★★★★★ [1 Abreviews](#) [3 References](#) [2 图像](#)

概述

产品名称	人TGF alpha ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 3 pg/ml
范围	1.37 pg/ml - 1000 pg/ml
回收率	93 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	94.49	84% - 105%
Serum	92.34	82% - 104%
Plasma	93.46	83% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's TGF alpha Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human TGF alpha in serum, plasma, cell culture supernatants</p> <p>This assay employs an antibody specific for Human TGF alpha coated on a 96-well plate. Standards and samples are pipetted into the wells and TGF alpha present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TGF alpha antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of TGF alpha bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
说明	Optimization may be required with urine samples.
平台	Microplate

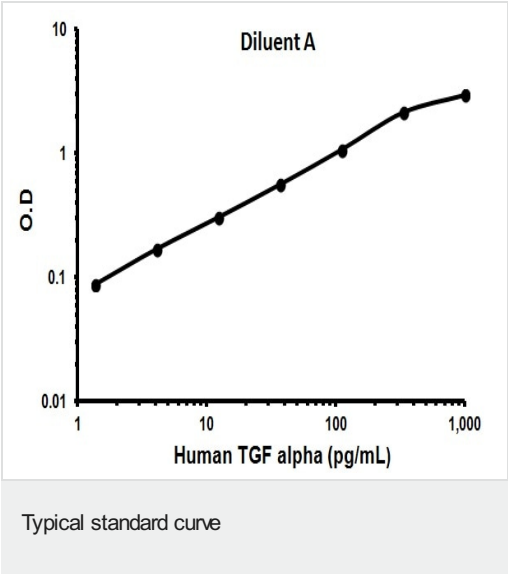
性能

存放说明Store at -20°C. Please refer to protocols.

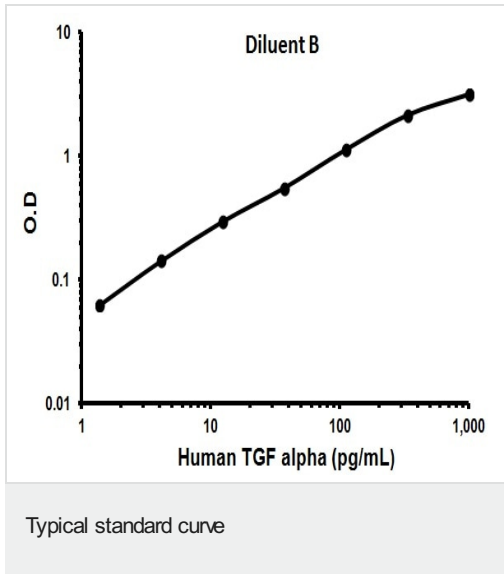
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human TGF alpha	2 vials
Recombinant Human TGF alpha Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TGF alpha Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

相关性Transforming growth factor alpha (TGF alpha) is an acid- and heat-stable originally discovered in rodents and humans. TGF alpha protien is 33% homologous to epidermal growth factor (EGF). TGF alpha binds to the EGF receptor, mediating tyrosine phosphorylation of the receptor, and promotes anchorage-independent growth of normal rat fibroblasts in soft agar in the presence of transforming growth factor-beta. TGF alpha is secreted by a variety of transformed cells and tumors, embryonic cells and some normal adult cells . TGF alpha bioactivity has also been found in the urine of cancer patients. It has been suggested that it may act as an autocrine growth factor for the induction or maintenance of malignancy.

图片



Representative standard curve using ab100646



Representative standard curve using ab100646

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