

Human Soluble TNF Receptor II ELISA Kit ab100643

[1 References](#) [2 图像](#)

概述

产品名称	人Soluble TNF Receptor II 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	< 5 pg/ml
范围	2.74 pg/ml - 2000 pg/ml
回收率	100 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	102.65	91% - 112%
Serum	104.92	95% - 115%
Plasma	105.49	94% - 118%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Soluble TNF Receptor II ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Soluble TNF Receptor II in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human Soluble TNF Receptor II coated on a 96-well plate. Standards and samples are pipetted into the wells and Soluble TNF Receptor II present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Soluble TNF Receptor II 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 Soluble TNF Receptor II bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明 Optimization may be required with urine samples.

平台 Microplate

性能

存放说明

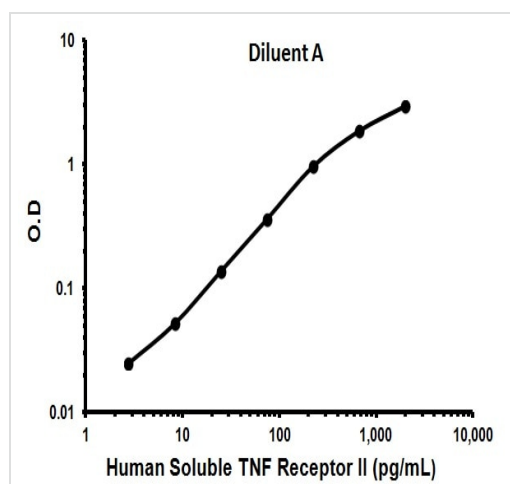
Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Soluble TNF Receptor II	2 vials
Recombinant Human Soluble TNF Receptor II Standard (lyophilized)	2 vials
Soluble TNF Receptor II Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

相关性

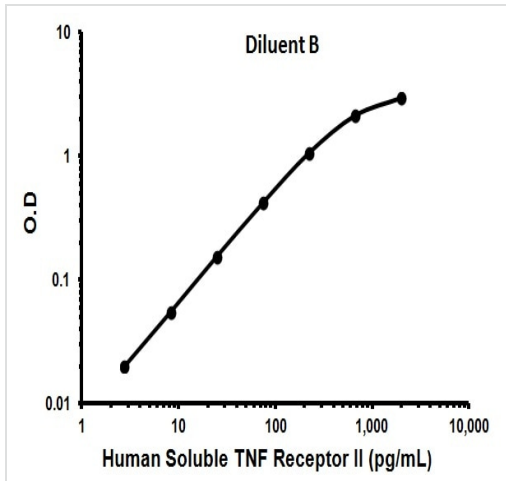
TNFR II (TNF Receptor II) is a member of the TNFR family of transmembrane proteins, and is expressed in immune cells and certain endothelial cells. It is a high affinity receptor for TNF alpha but manifests a lower affinity to TNF beta. Signaling through this receptor regulates various biological processes including cell proliferation, differentiation, apoptosis, lipid metabolism, coagulation, and neurotransmission. TNFR II consists of a 240 amino acid residue extracellular ligand binding portion, a 27 amino acid residue transmembrane segment and a 173 amino acid residue cytoplasmic domain. Soluble TNFR II (Soluble TNF Receptor II) is a 268 amino acid isoform containing the extracellular ligand-binding portion of TNFR II. It binds to TNF alpha and is capable of inhibiting TNF alpha induced activities by acting as a decoy receptor.

图片



Representative standard curve using ab100643

Typical standard curve



Representative standard curve using ab100643

Typical standard curve

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