

Product datasheet

Human Soluble TNF Receptor II ELISA Kit ab100643

画像数 2

製品の概要

製品名	Human Soluble TNF Receptor II ELISA Kit
検出方法	Colorimetric
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	< 5 pg/ml
検出範囲	2.74 pg/ml - 2000 pg/ml
添加回収試験	100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	102.65	91% - 112%
Serum	104.92	95% - 115%
Plasma	105.49	94% - 118%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

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.

特記事項 Optimization may be required with urine samples.

アプリケーション 適用あり: Sandwich ELISA

製品の特性

保存方法

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

内容	1 x 96 tests
20X Wash Buffer Concentrate	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 (lyophilized)	2 vials
Recombinant Human Soluble TNF Receptor II Standard	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

関連性

TNFRII (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. TNFRII 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 TNFRII (Soluble TNF Receptor II) is a 268 amino acid isoform containing the extracellular ligand-binding portion of TNFRII. It binds to TNF alpha and is capable of inhibiting TNF alpha induced activities by acting as a decoy receptor.

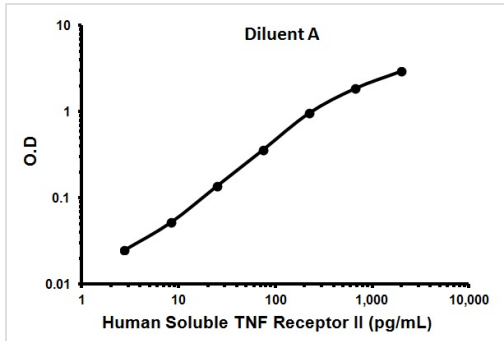
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab100643** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

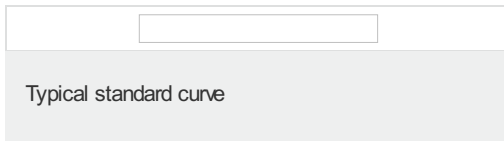
アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



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 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