

Human Soluble TNF Receptor I ELISA Kit ab100642

3 References 1 Image

Overview

Product name	Human Soluble TNF Receptor I ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1 pg/ml
Range	0.55 pg/ml - 400 pg/ml
Recovery	80 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	80.02	71% - 90%
Serum	80.43	72% - 91%
Plasma	81.36	73% - 91%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's Human STNF-R1 (soluble tumor necrosis factor receptor I) ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human STNF-R1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human STNF-R1 coated on a 96-well plate. Standards and samples are pipetted into the wells and STNF-R1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human STNF-R1 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 STNF-R1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Platform Microplate

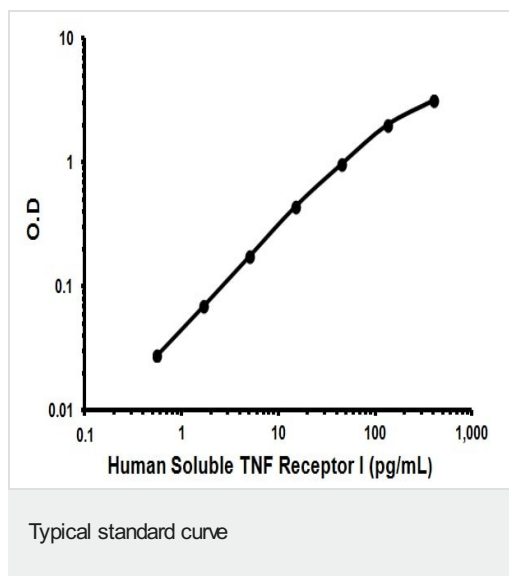
Properties

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

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

Function	Receptor for TNFSF2/TNF-alpha and homotrimeric TNFSF1/lymphotoxin-alpha. The adapter molecule FADD recruits caspase-8 to the activated receptor. The resulting death-inducing signaling complex (DISC) performs caspase-8 proteolytic activation which initiates the subsequent cascade of caspases (aspartate-specific cysteine proteases) mediating apoptosis. Contributes to the induction of non-cytocidal TNF effects including anti-viral state and activation of the acid sphingomyelinase.
Involvement in disease	Familial hibernian fever Multiple sclerosis 5
Sequence similarities	Contains 1 death domain. Contains 4 TNFR-Cys repeats.
Domain	The domain that induces A-SMASE is probably identical to the death domain. The N-SMASE activation domain (NSD) is both necessary and sufficient for activation of N-SMASE. Both the cytoplasmic membrane-proximal region and the C-terminal region containing the death domain are involved in the interaction with TRPC4AP.
Post-translational modifications	The soluble form is produced from the membrane form by proteolytic processing.
Cellular localization	Cell membrane. Golgi apparatus membrane. Secreted. A secreted form is produced through proteolytic processing and Secreted. Lacks a Golgi-retention motif, is not membrane bound and therefore is secreted.

Images



Representative standard curve using ab100642

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.com/abpromise> or contact our technical team.

Terms and conditions

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