

Product datasheet

Human TNFSF4 ELISA Kit (OX40L) ab213842

[2 References](#) [1 Image](#)

Overview

Product name Human TNFSF4 ELISA Kit (OX40L)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	209pg/ml	8.56	= 4.1%
2	16	1225pg/ml	51.45	= 4.2%
3	16	4641pg/ml	305.7	= 6.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	228pg/ml	12.54	= 5.5%
2	24	1277pg/ml	72.78	= 5.7%
3	24	4885pg/ml	365.77	= 7.5%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 156 pg/ml - 10000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview The Human TNFSF4 Enzyme-Linked Immunosorbent Assay (ELISA) kit (OX40L) (ab213842) is designed for the quantitative measurement of Human TNFSF4 in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for TNFSF4/OX40L has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: Q51-L183) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for TNFSF4/OX40L is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human TNFSF4/OX40L amount of sample captured in plate.

Notes OX40L is the ligand for CD134 and is expressed on such cells as DC2s enabling amplification of Th2 cell differentiation. OX40L has also been designated CD252. By analysis of an interspecific backcross, the mouse *Tnfsf4* gene is mapped to chromosome 1. Using fluorescence *in situ* hybridization, they localized the human TNFSF4 gene to 1q25, a region sharing homology of synteny with the portion of mouse chromosome 1 containing the *Tnfsf4* gene. OX40L strongly inhibited the generation of IL-10-producing Tr1 cells induced by two physiologic stimuli, the inducible costimulatory ligand and immature dendritic cells. In addition, OX40L strongly inhibited IL-10 production and suppressive function of differentiated IL-10-producing Tr1 cells. *Tnfsf4* underlies Ath1 in mice and polymorphisms in its human homolog TNFSF4 increase the risk of myocardial infarction in humans.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human TNFSF4 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human TNFSF4 antibody		1 x 100µl
Lyophilized recombinant Human TNFSF4 standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml

Function Cytokine that binds to TNFRSF4. Co-stimulates T-cell proliferation and cytokine production.

Involvement in disease Genetic variations in TNFSF4 influence susceptibility to systemic lupus erythematosus (SLE) [MIM:152700]. SLE is a chronic, inflammatory and often febrile multisystemic disorder of connective tissue. It affects principally the skin, joints, kidneys and serosal membranes. It is

thought to represent a failure of the regulatory mechanisms of the autoimmune system. Note=The upstream region of TNFSF4 contains a single risk haplotype for SLE, which is correlated with increased expression of both cell-surface TNFSF4 and TNFSF4 transcripts. Increased levels of TNFSF4 are thought to augment T cell-APC interaction and the functional consequences of T cell activation, thereby destabilizing peripheral tolerance.

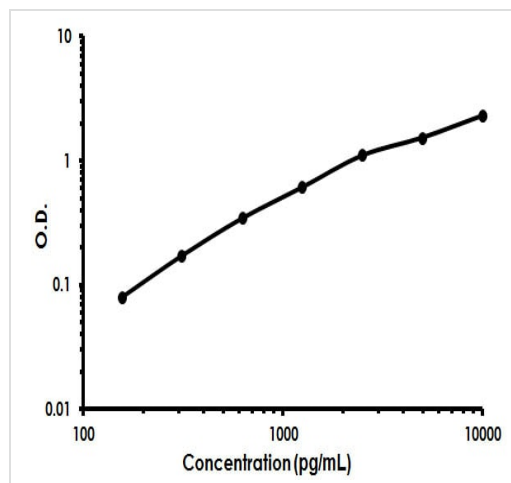
Sequence similarities

Belongs to the tumor necrosis factor family.

Cellular localization

Membrane.

Images



Human TNFSF4 ELISA Kit (OX40L) (ab213842) Standard Curve.

Human TNFSF4 ELISA Kit (OX40L) (ab213842)
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.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