

Mouse TNFRSF7 ELISA Kit (CD27) ab213895

1 Image

Overview

Product name Mouse TNFRSF7 ELISA Kit (CD27)

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	1	16	533pg/ml	22.4	4.2%
	2	16	1626pg/ml	56.91	3.5%
	3	16	2563pg/ml	138.4	5.4%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	1	24	658pg/ml	39.5	6%
	2	24	1856pg/ml	89.1	4.8%
	3	24	2731pg/ml	180.3	6.6%

Sample type Cell culture supernatant, Serum

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 62.5 pg/ml - 4000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview The Mouse TNFRSF7 Enzyme-Linked Immunosorbent Assay (ELISA) kit (CD27) (ab213895) is designed for the quantitative detection of Mouse TNFRSF7 in cell culture supernatants and serum.

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from rat specific for TNFRSF7/CD27 has been precoated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: T21-R182) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for TNFRSF7/CD27 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was 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 Mouse TNFRSF7/CD27 amount of sample captured in plate.

## Notes

CD27 antigen, also called CD27 or TNFRSF7 is a tumor necrosis factor receptor. This gene is mapped to 12p13.31. The protein encoded by this gene is a member of the TNF-receptor superfamily. This receptor is required for generation and long-term maintenance of T cell immunity. It binds to ligand CD70, and plays a key role in regulating B-cell activation and immunoglobulin synthesis. This receptor transduces signals that lead to the activation of NF-kappaB and MAPK8/JNK. Adaptor proteins TRAF2 and TRAF5 have been shown to mediate the signaling process of this receptor. CD27-binding protein (SIVA), a proapoptotic protein, can bind to this receptor and is thought to play an important role in the apoptosis induced by this receptor.

## 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-Mouse TNFRSF7 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Mouse TNFRSF7 antibody		1 x 100µl
Lyophilized recombinant Mouse TNFRSF7 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
Wash Buffer (25X)		1 x 20ml

## Function

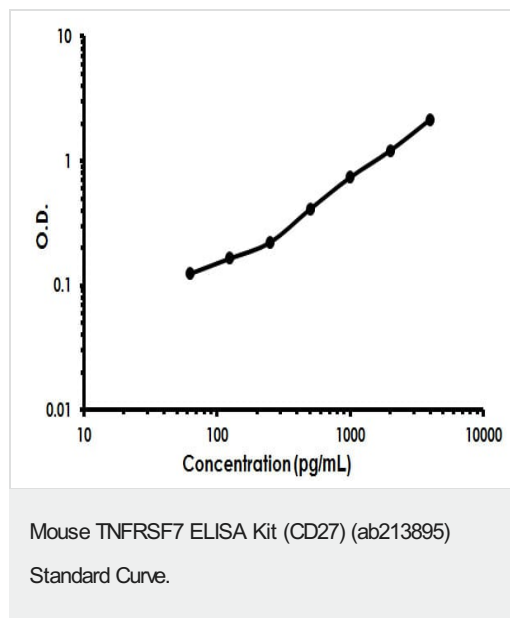
Receptor for CD70/CD27L. May play a role in survival of activated T-cells. May play a role in apoptosis through association with SIVA1.

## Tissue specificity

Found in most T-lymphocytes.

<b>Sequence similarities</b>	Contains 3 TNFR-Cys repeats.
<b>Post-translational modifications</b>	Phosphorylated and O-glycosylated.
<b>Cellular localization</b>	Membrane.

## Images



Mouse TNFRSF7 ELISA Kit (CD27) (ab213895) 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