

Product datasheet

Human TNFSF18 / GITRL ELISA Kit ab113323

[1 Image](#)

Overview

Product name	Human TNFSF18 / GITRL ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 25 pg/ml
Range	41 pg/ml - 10000 pg/ml
Recovery	97 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	87.83	67% - 100%
Serum	116.9	108% - 123%
Plasma	87.69	67% - 100%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human TNFSF18 / GITRL ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of human TNFSF18 / GITRL in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for human TNFSF18 / GITRL coated on a 96-well plate. Standards and samples are pipetted into the wells and TNFSF18 / GITRL present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TNFSF18 / GITRL 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 TNFSF18 / GITRL 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	Identifier	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 GITR Ligand	Item F	2 vials
TNFSF18 Ligand Microplate (12 x 8 wells)		1 unit
Recombinant Human GITR Ligand Standard (lyophilized)		2 vials
Stop Solution		1 x 8ml
TMB One-Step Substrate Reagent		1 x 12ml

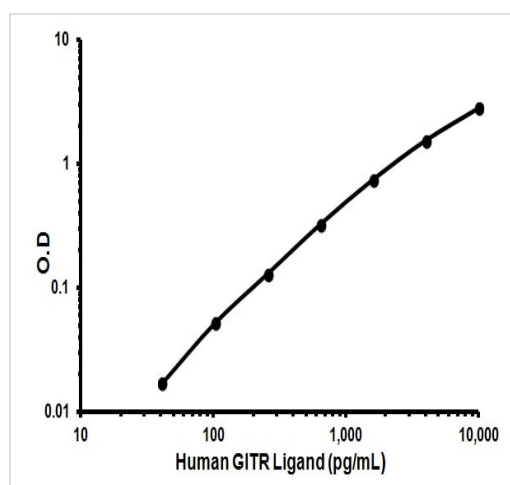
Function Cytokine that binds to TNFRSF18/AITR/GITR. Important for interactions between activated T-lymphocytes and endothelial cells and may modulate T-lymphocyte survival in peripheral tissues.

Tissue specificity Expressed at high levels in the small intestine, ovary, testis, kidney and endothelial cells.

Sequence similarities Belongs to the tumor necrosis factor family.

Cellular localization Membrane.

Images



Representative standard curve using ab113323.

Typical data

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