

Human TNFSF18 / GITRL ELISA Kit ab113323

1 图像

概述

| | |
|------|---|
| 产品名称 | 人TNFSF18 / GITRL ELISA试剂盒 |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Serum, Plasma |
| 检测类型 | Sandwich (quantitative) |
| 灵敏度 | < 25 pg/ml |
| 范围 | 41 pg/ml - 10000 pg/ml |
| 回收率 | 97 % |

特定样本回收率

| 样品类型 | 平均 % | 范围 |
|--------------------------|-------|-------------|
| Cell culture supernatant | 87.83 | 67% - 100% |
| Serum | 116.9 | 108% - 123% |
| Plasma | 87.69 | 67% - 100% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 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.

说明 Optimization may be required with urine samples.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

| 组件 | 标识符 | 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 |

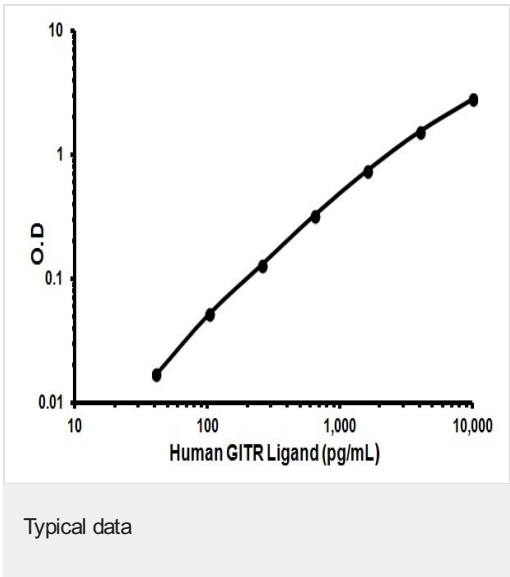
功能 Cytokine that binds to TNFRSF18/AITR/GITR. Important for interactions between activated T-lymphocytes and endothelial cells and may modulate T-lymphocyte survival in peripheral tissues.

组织特异性 Expressed at high levels in the small intestine, ovary, testis, kidney and endothelial cells.

序列相似性 Belongs to the tumor necrosis factor family.

细胞定位 Membrane.

图片



Representative standard curve using ab113323.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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