

Product datasheet

Human TNFRSF18 ELISA Kit (GITR) ab100528

1 图像

概述

产品名称	人TNFRSF18 ELISA试剂盒(GITR)
检测方法	Colorimetric
样品类型	Cell culture extracts, Tissue Extracts
检测类型	Sandwich (quantitative)
灵敏度	< 15 pg/ml
范围	13.7 pg/ml - 10000 pg/ml
复检	94 %

特定样本复检结果

样品类型	平均%	范围
Cell culture extracts	93.49	83% - 103%
Tissue Extracts	94.92	82% - 102%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's TNFRSF18 (GITR) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human TNFRSF18 in cell lysates and tissue lysates.

This assay employs an antibody specific for Human TNFRSF18 coated on a 96-well plate. Standards and samples are pipetted into the wells and TNFRSF18 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TNFRSF18 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 TNFRSF18 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

经测试应用 适用于: Sandwich ELISA

平台 Microplate

性能

存放说明

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

组件	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
2x Cell lysate buffer	1 x 5ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
5X Sample Diluent Buffer	1 x 10ml
Biotinylated anti-Human TNFRSF18	2 vials
Recombinant Human TNFRSF18 Standard	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml
TNFRSF18 Microplate (12 x 8 wells)	1 unit

相关性

TNFRSF18 or GITR (glucocorticoid-induced TNF receptor family-regulated gene) is a 25 kD TNF receptor superfamily member (also known as AITR and TNFRSF18) originally identified in the mouse by comparing untreated and dexamethasone-treated murine T cell hybridoma cells. Human GITR was subsequently identified by searching an expressed sequence tag database. GITR is expressed on activated lymphocytes and is upregulated by T cell receptor engagement. The cytoplasmic domain of GITR is homologous to CD40, 4-1BB and CD27 and has been shown to interact with TRAF 1-3, but not TRAF 5 or 6. GITR signaling has been shown to regulate T cell proliferation and TCR-mediated apoptosis, and to break immunological self-tolerance. GITR appears to be highly expressed on CD4+CD25+ T regulatory cells and has been shown to induce NF-κB activation through TRAF2/NIK signaling. GITR has been proposed to be involved in the development of regulatory T cells and to regulate the activity of Th1 subsets. GITR binds GITRL, a TNF superfamily ligand expressed on endothelial cells, dendritic cells, macrophages, and B cells.

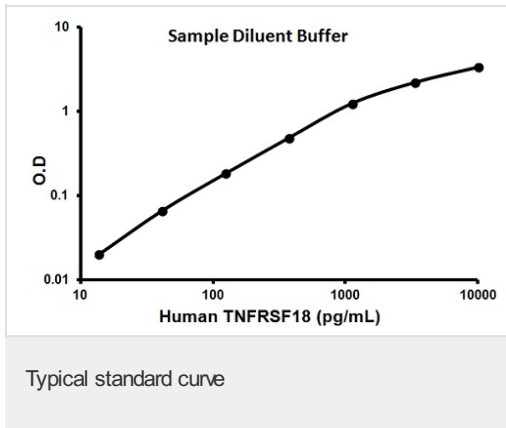
应用

Our [Abpromise guarantee](#) covers the use of **ab100528** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



Representative standard curve using
ab100528

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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