

Human TIE1 ELISA Kit ab100649

2 图像

概述

产品名称	人TIE1 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant
检测类型	Sandwich (quantitative)
灵敏度	< 0.6 ng/ml
范围	0.68 ng/ml - 500 ng/ml
回收率	100 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	105.9	93% - 115%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's TIE1 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human TIE1 in cell culture supernatants.

This assay employs an antibody specific for Human TIE1 coated on a 96-well plate. Standards and samples are pipetted into the wells and TIE1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TIE1 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 TIE1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Human Tie-1 in normal serum with ab100649, only in serum spiked with Human Tie-1. Plasma is not recommended.

说明 Serum has been validated as a sample matrix with this kit. However, TIE1 concentrations are low in normal serum and therefore may not be detected by this assay.

Plasma is not recommended.

Optimization may be required with urine samples.

平台

Microplate

性能

存放说明

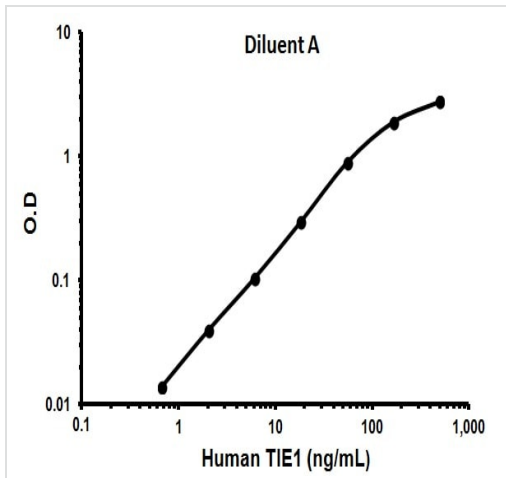
Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human TIE1	2 vials
Recombinant Human TIE1 Standard (lyophilized)	2 vials
Stop Solution	1 x 25ml
TIE1 Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

相关性

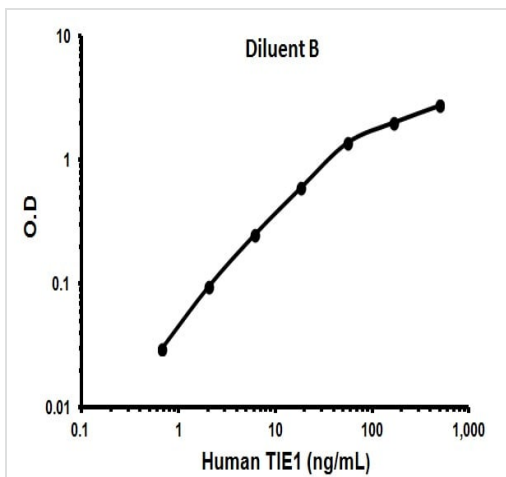
TIE1/TIE (tyrosine kinase with Ig and EGF homology domains 1) and TIE2/Tek define a new class of the receptor tyrosine kinase (RTK) subfamily with unique structural characteristics: two immunoglobulin like domains flanking three epidermal growth factor (EGF) like domains followed by three fibronectin type III like repeats in the extracellular region and a split tyrosine kinase domain in the cytoplasmic region. Human TIE1 cDNA encodes a 1138 amino acid residue precursor protein with a putative signal peptide, an extracellular domain, and a cytoplasmic domain. Human TIE1/Fc, a disulfide linked homodimeric protein, has a calculated molecular mass of approximately 107 kDa. Due to glycosylation, the protein migrates to approximately 160 kDa in SDS PAGE under reducing conditions. TIE1 and TIE2, expressed primarily on endothelial and hematopoietic progenitor cells, play important roles in angiogenesis, vasculogenesis, and hematopoiesis. In developing vascular endothelial cells, TIE1 and TIE2 are specifically expressed. Two ligands that bind TIE have been identified, angiopoietin 1 and angiopoietin 2. Based on gene targeting studies, the in vivo functions of TIE1 are related to endothelial cell differentiation. The receptor tyrosine kinase TIE also plays a role in the survival and integrity of the endothelium.

图片



Representative standard curve using ab100649

Typical standard curve



Representative standard curve using ab100649

Typical 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.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