

Human TIE1 ELISA Kit ab100649

画像数 2

製品の概要

製品名	Human TIE1 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant
アッセイタイプ	Sandwich (quantitative)
検出感度	< 0.6 ng/ml
検出範囲	0.68 ng/ml - 500 ng/ml
添加回収試験	100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	105.9	93% - 115%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

特記事項 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.

試験プラットフォーム

Microplate

製品の特性

保存方法

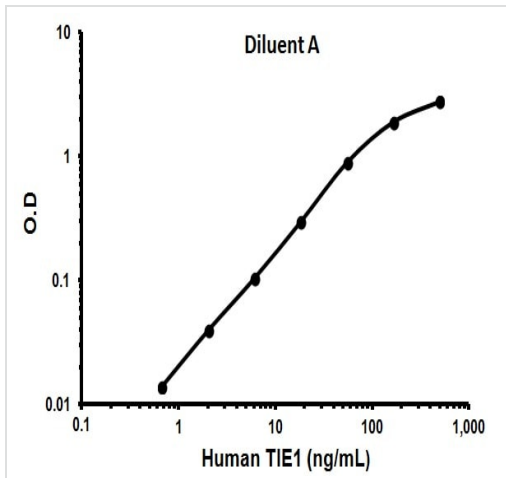
Store at -20°C. Please refer to protocols.

内容	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

関連性

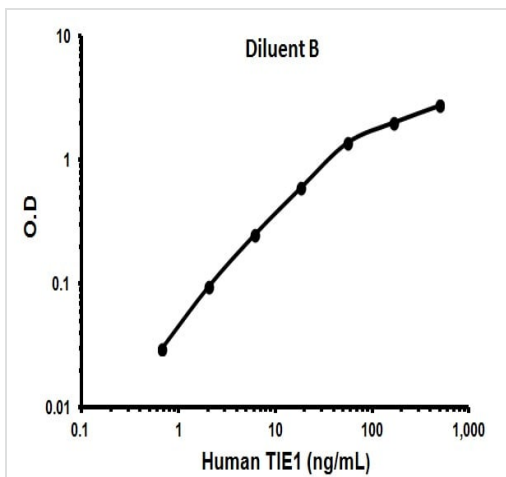
TIE1/TIE (tyrosine kinase with Ig and EGF homology domains 1) and TIE2/Tek define a new class of the receptor tyrosine kinase (RTK) subfamily with unique structural characteristics: two immunoglobulin like domains flanking three epidermal growth factor (EGF) like domains followed by three fibronectin 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.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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