

Human TIMP4 ELISA Kit ab113328

1 References 画像数 2

製品の概要

製品名	Human TIMP4 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 25 pg/ml
検出範囲	28.67 pg/ml - 7000 pg/ml
添加回収試験	80 %

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	96.11	89% - 104%
Serum	73.49	67% - 82%
Plasma	72.83	67% - 81%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's TIMP4 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human TIMP4 in serum, plasma (collect plasma using EDTA and citrate as an anticoagulant. Heparin is not recommended), cell culture supernatants.

This assay employs an antibody specific for Human TIMP4 coated on a 96- well plate. Standards and samples are pipetted into the wells and TIMP4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TIMP4 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 TIMP4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

特記事項 Optimization may be required with urine samples.

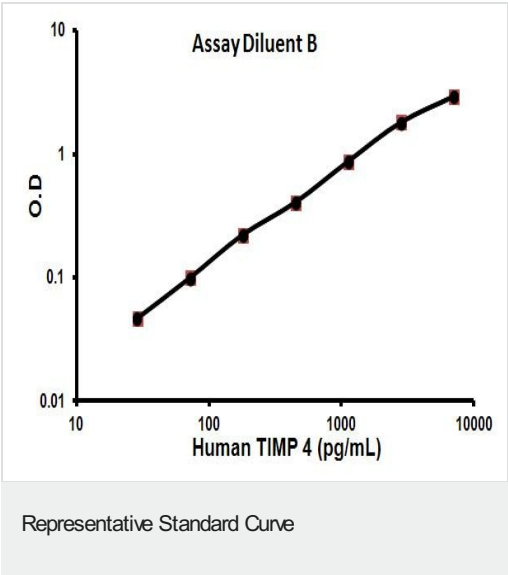
The standard is affinity purified.

製品の特Ⓕ性

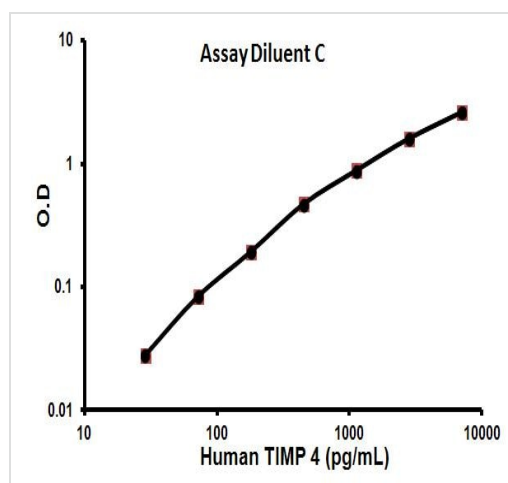
保存方法Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200μl
Assay Diluent C	1 x 30ml
Biotinylated anti-Human TIMP4	2 vials
Recombinant Human TIMP4 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TIMP4 Microplate (12 strips x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

画像



Representative Standard Curve using ab113328



Representative Standard Curve using ab113328

Representative 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