

3 References 画像数 2

製品名	Human Endostatin 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	< 10 pg/ml
検出範囲	13.7 pg/ml - 10000 pg/ml
添加回収試験	> 90 %

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	95.32	84% - 108%
Serum	88.47	78% - 102%
Plasma	92.31	80% - 104%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Abcam's Human Endostatin ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Endostatin in serum, plasma, and cell culture supernatants.

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

Get results in 90 minutes with Human Endostatin ELISA Kit (**ab184864**) from our SimpleStep

ELISA® range.

特記事項

Optimisation 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

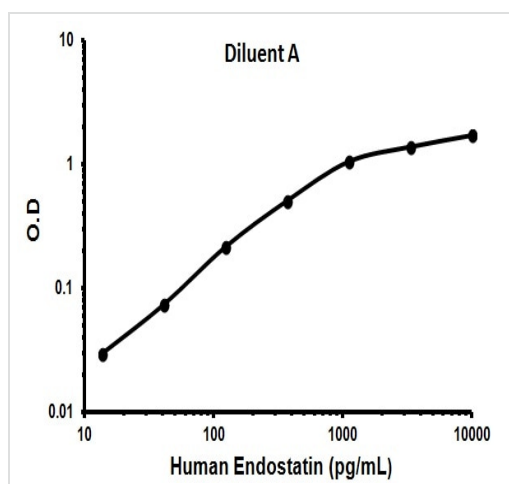
Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Detection Antibody Endostatin	2 units
Endostatin Microplate (12 strips x 8 wells)	1 unit
Endostatin Standard Human(recombinant) (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

関連性

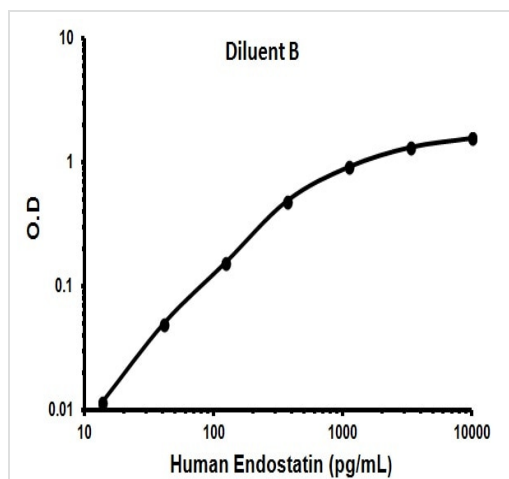
Endostatin (20-22 kDa), a cleaved product of the C-terminal domain of collagen XVIII, is an endogenous angiogenesis inhibitor. It was originally found in conditioned media from a murine endothelial tumor cell line, hemangioendothelioma. Endostatin inhibits endothelial cell migration (in vivo and in vitro) and induces endothelial cell apoptosis. It inhibits tumor growth and impairs blood vessel maturation in wound healing. Endostatin has an important role in endothelial cell adhesion and cytoskeletal organization. Endostatin can be found in vessel walls (elastic fibers) and basement membranes.

画像



Representative Standard Curve using ab100508

Typical Standard Curve



Representative Standard Curve using ab100508

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