

# Human Antithrombin III ELISA Kit ab108801

画像数 7

## 製品の概要

**製品名** Human Antithrombin III ELISA Kit

**検出方法** Colorimetric

**再現性** Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall				4.8%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				7.1%

**サンプルの種類** Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Cell culture media, Cerebral Spinal Fluid

**アッセイタイプ** Sandwich (quantitative)

**検出感度** = 3 ng/ml

**検出範囲** 3.13 ng/ml - 200 ng/ml

**添加回収試験** 98.5 %

**全工程の試験時間** 4h 0m

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Human

**製品の概要** Abcam's Antithrombin III Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Antithrombin III in urine, plasma, milk, saliva, cerebrospinal fluid, serum, and cell culture supernatants.

An antithrombin III specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an antithrombin III specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the

amount of antithrombin III captured in plate.

#### アプリケーション

**適用あり:** Sandwich ELISA

#### 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
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Antithrombin III Antibody	1 x 120µl
Antithrombin III Microplate (12 x 8 well strips)	1 unit
Antithrombin III Standard	1 vial
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml

#### 機能

Most important serine protease inhibitor in plasma that regulates the blood coagulation cascade. AT-III inhibits thrombin, matriptase-3/TMPRSS7, as well as factors IXa, Xa and XIa. Its inhibitory activity is greatly enhanced in the presence of heparin.

#### 組織特異性

Found in plasma.

#### 関連疾患

Antithrombin III deficiency

#### 配列類似性

Belongs to the serpin family.

#### 翻訳後修飾

Phosphorylation sites are present in the extracellular medium.

#### 細胞内局在

Secreted > extracellular space.

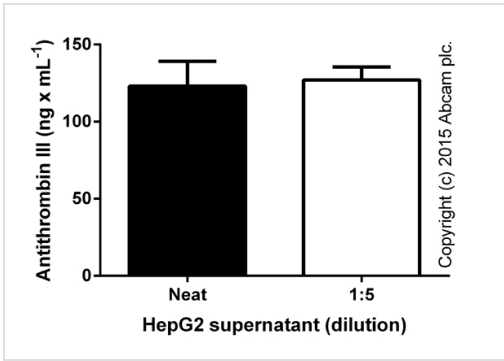
### アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab108801** in the following tested applications.

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

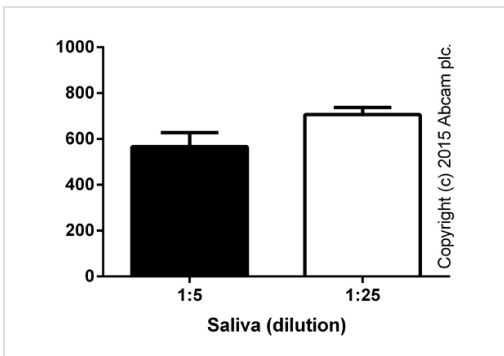
アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

### 画像



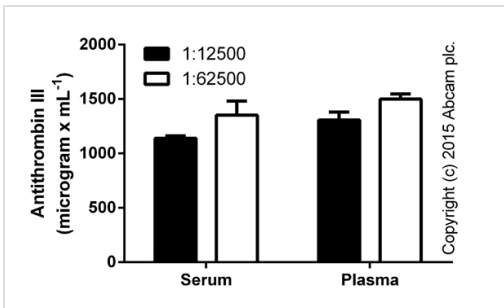
Linearity: Antithrombin III in samples at different dilutions, corrected for the factor of dilution (duplicates +/- SD)

ELISA: Antithrombin III Human ELISA Kit



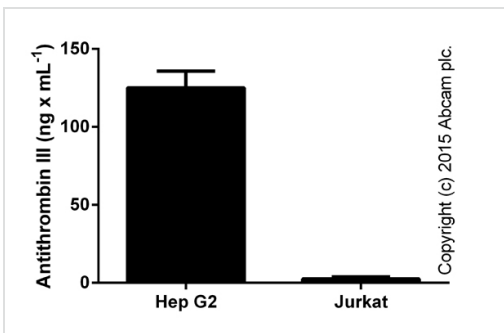
Linearity: Antithrombin III in samples at different dilutions, corrected for the factor of dilution (duplicates +/- SD), showing ng x mL<sup>-1</sup> on y-axis.

ELISA: Antithrombin III Human ELISA Kit



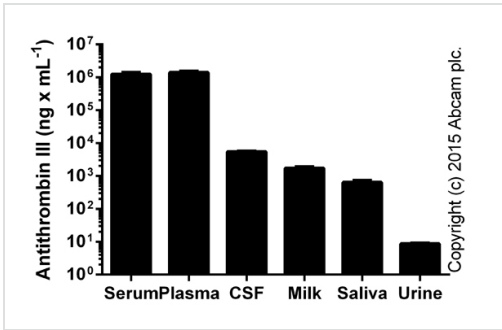
Linearity: Antithrombin III in samples at different dilutions, corrected for the factor of dilution (duplicates +/- SD)

ELISA: Antithrombin III Human ELISA Kit



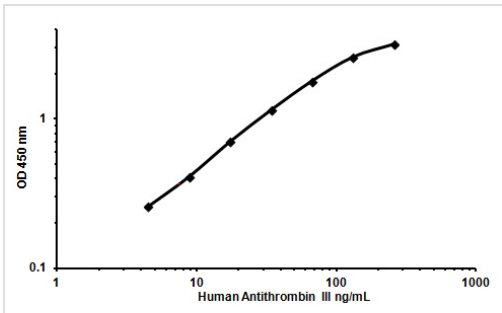
Antithrombin III in cell culture supernatants (duplicates +/- SD), tested in the dilution range of 1:1-1:100.

ELISA: Antithrombin III Human ELISA Kit

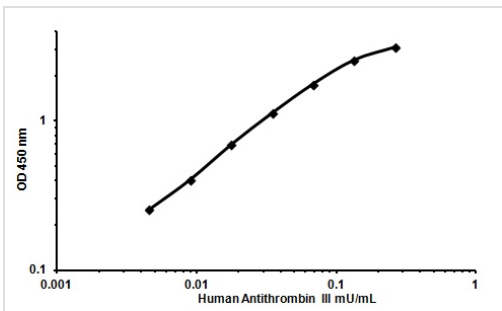


Antithrombin III in human biological fluids (duplicates +/- SD). The samples were differently diluted (serum and Li-heparine plasma: 1e4-1e5; CSF: 1e2-1e3; milk: 10-100; saliva: 5-100; urine: neat)

ELISA: Antithrombin III Human ELISA Kit



Typical Standard Curve



Representative standard curve using ab108801.

Typical Standard Curve

**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 <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