

### Mouse Prothrombin / Thrombin Total ELISA Kit ab157527

3 References [画像数 1](#)

#### 製品の概要

**製品名** Mouse Prothrombin / Thrombin Total 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%
1	20	2.81ng/ml	0.267	9.51%
2	20	14.1ng/ml	0.56	3.98%
3	20	39.5ng/ml	1.4	3.53%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
1	10	2.42ng/ml	0.241	9.96%
2	10	13.8ng/ml	1.2	8.72%
3	10	145ng/ml	7.18	4.95%

**サンプルの種類** Serum, Plasma, Cell culture media

**アッセイタイプ** Sandwich (quantitative)

**検出感度** 0.52 ng/ml

**検出範囲** 1 ng/ml - 500 ng/ml

**添加回収試験** 95.18 %

**全工程の試験時間** 1h 50m

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Mouse

**製品の概要** Abcam's Total Prothrombin/Thrombin ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for is for the quantitative determination of total Prothrombin and Thrombin in mouse plasma, serum and cell culture media.

Mouse Prothrombin, Thrombin and thrombin-antithrombin complex will bind to the capture antibody coated on the microtiter plate. After appropriate washing steps, biotinylated primary antibody binds to the captured protein. Excess primary antibody is washed away and bound antibody is reacted with horseradish peroxidase conjugated streptavidin. TMB substrate is used for color development at 450nm. A standard calibration curve is prepared along with the samples to be measured using dilutions of Prothrombin. The amount of color development is directly proportional to the concentration of Prothrombin in the sample.

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Wash Buffer	1 x 50ml
5X Diluent	1 x 50ml
Anti-mouse Prothrombin primary antibody	1 vial
Anti-prothrombin coated Microtiter Plate (8 x 12 wells)	1 unit
HRP-conjugated Streptavidin	1 vial
Prothrombin Standard	1 vial
TMB Substrate Solution	1 x 10ml

#### 機能

Thrombin, which cleaves bonds after Arg and Lys, converts fibrinogen to fibrin and activates factors V, VII, VIII, XIII, and, in complex with thrombomodulin, protein C. Functions in blood homeostasis, inflammation and wound healing.

#### 組織特異性

Expressed by the liver and secreted in plasma.

#### 関連疾患

Defects in F2 are the cause of factor II deficiency (FA2D) [MIM:613679]. It is a very rare blood coagulation disorder characterized by mucocutaneous bleeding symptoms. The severity of the bleeding manifestations correlates with blood factor II levels.

Genetic variations in F2 may be a cause of susceptibility to ischemic stroke (ISCHSTR) [MIM:601367]; also known as cerebrovascular accident or cerebral infarction. A stroke is an acute neurologic event leading to death of neural tissue of the brain and resulting in loss of motor, sensory and/or cognitive function. Ischemic strokes, resulting from vascular occlusion, is considered to be a highly complex disease consisting of a group of heterogeneous disorders with multiple genetic and environmental risk factors.

Defects in F2 are a cause of susceptibility to thrombosis (THR) [MIM:188050]. It is a multifactorial disorder of hemostasis characterized by abnormal platelet aggregation in response to various agents and recurrent thrombi formation. Note=A common genetic variation in the 3-prime untranslated region of the prothrombin gene is associated with elevated plasma prothrombin levels and an increased risk of venous thrombosis.

#### 配列類似性

Belongs to the peptidase S1 family.

Contains 1 Gla (gamma-carboxy-glutamate) domain.

Contains 2 kringle domains.

Contains 1 peptidase S1 domain.

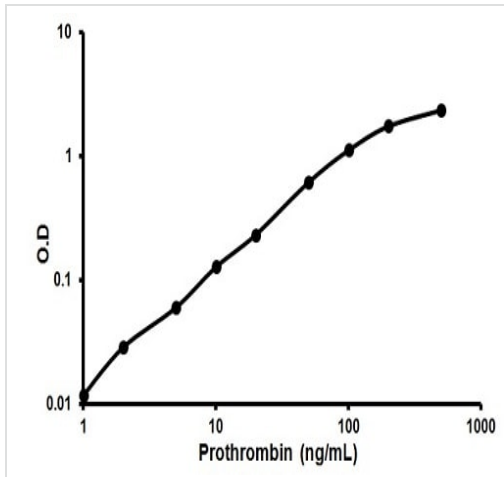
#### 翻訳後修飾

The gamma-carboxyglutamyl residues, which bind calcium ions, result from the carboxylation of glutamyl residues by a microsomal enzyme, the vitamin K-dependent carboxylase. The modified residues are necessary for the calcium-dependent interaction with a negatively charged phospholipid surface, which is essential for the conversion of prothrombin to thrombin.

#### 細胞内局在

Secreted > extracellular space.

#### 画像



Example Standard Curve for ab157527

**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