

### Human Factor XI ELISA Kit ab108834

1 References [画像数 1](#)

#### 製品の概要

| 製品名      | Human Factor XI ELISA Kit  |      |    |      |    |     |         |  |  |  |      |
|----------|--|------|----|------|----|-----|---------|--|--|--|------|
| 検出方法     | Colorimetric   |      |    |      |    |     |         |  |  |  |      |
| 再現性      | Intra-Assay (同時再現性)  |      |    |      |    |     |         |  |  |  |      |
|          | <table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>4.9%</td> </tr> </tbody> </table>  | サンプル | N  | 平均値  | SD | CV% | Overall |  |  |  | 4.9% |
| サンプル     | N  | 平均値  | SD | CV%  |    |     |         |  |  |  |      |
| Overall  |  |      |    | 4.9% |    |     |         |  |  |  |      |
|          | Inter-Assay (日差再現性)  |      |    |      |    |     |         |  |  |  |      |
|          | <table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>10%</td> </tr> </tbody> </table>   | サンプル | N  | 平均値  | SD | CV% | Overall |  |  |  | 10%  |
| サンプル     | N  | 平均値  | SD | CV%  |    |     |         |  |  |  |      |
| Overall  |  |      |    | 10%  |    |     |         |  |  |  |      |
| サンプルの種類  | Serum, Plasma  |      |    |      |    |     |         |  |  |  |      |
| アッセイタイプ  | Sandwich (quantitative)  |      |    |      |    |     |         |  |  |  |      |
| 検出感度     | 0.69 ng/ml   |      |    |      |    |     |         |  |  |  |      |
| 検出範囲     | 1.563 ng/ml - 50 ng/ml   |      |    |      |    |     |         |  |  |  |      |
| 添加回収試験   | 96 %   |      |    |      |    |     |         |  |  |  |      |
| 全工程の試験時間 | 4h 00m   |      |    |      |    |     |         |  |  |  |      |
| ステップ     | Multiple steps standard assay  |      |    |      |    |     |         |  |  |  |      |
| 種交差性     | 交差種: Human   |      |    |      |    |     |         |  |  |  |      |
| 製品の概要    | <p>Abcam's Factor XI Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor XI levels in human plasma and serum samples.</p> <p>A Factor XI specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor XI 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 Factor XI captured in plate.</p> |      |    |      |    |     |         |  |  |  |      |

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

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 M Concentrate                 | 1 x 30ml     |
| 20X Wash Buffer Concentrate               | 2 x 30ml     |
| 50X Biotinylated Human Factor XI Antibody | 1 x 120µl    |
| Chromogen Substrate                       | 1 x 7ml      |
| Factor XI Microplate (12 x 8 well strips) | 1 unit       |
| Factor XI Standard                        | 1 vial       |
| Sealing Tapes                             | 3 units      |
| Stop Solution                             | 1 x 11ml     |

### 機能

Factor XI triggers the middle phase of the intrinsic pathway of blood coagulation by activating factor IX.

### 組織特異性

Isoform 2 is produced by platelets and megakaryocytes but absent from other blood cells.

### 関連疾患

Defects in F11 are the cause of factor XI deficiency (FA11D) [MIM:612416]; also known as plasma thromboplastin antecedent deficiency or Rosenthal syndrome. It is a hemorrhagic disease characterized by reduced levels and activity of factor XI resulting in moderate bleeding symptoms, usually occurring after trauma or surgery. Patients usually do not present spontaneous bleeding but women can present with menorrhagia. Hemorrhages are usually moderate.

### 配列類似性

Belongs to the peptidase S1 family. Plasma kallikrein subfamily.

Contains 4 apple domains.

Contains 1 peptidase S1 domain.

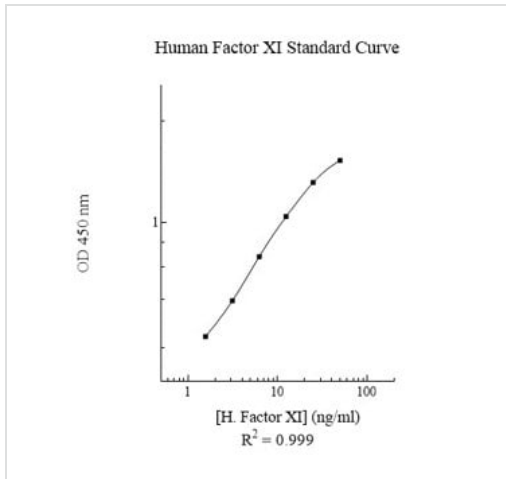
### 翻訳後修飾

Activated by factor XIIa (or XII), which cleaves each polypeptide after Arg-387 into the light chain, which contains the active site, and the heavy chain, which associates with high molecular weight (HMW) kininogen.

### 細胞内局在

Secreted.

## 画像



Representative Standard Curve using ab108834.

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