

## Product datasheet

# Human Factor XI ELISA Kit (with plasma controls) ab168547

★★★★★ [1 Abreviews](#) [1 References](#) [1 Image](#)

### Overview

**Product name** Human Factor XI ELISA Kit (with plasma controls)

**Detection method** Colorimetric

**Precision**

Intra-assay

| Sample  | n | Mean | SD | CV%  |
|---------|---|------|----|------|
| Overall |   |      |    | 4.9% |

Inter-assay

| Sample  | n | Mean | SD | CV% |
|---------|---|------|----|-----|
| Overall |   |      |    | 10% |

**Sample type** Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** > 0.69 ng/ml

**Range** 3.13 ng/ml - 25 ng/ml

**Recovery** = 96 %

Sample specific recovery

| Sample type | Average % | Range      |
|-------------|-----------|------------|
| Serum       |           | 95% - 106% |
| Plasma      |           | 94% - 105% |

**Assay time** 4h 00m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's Factor XI Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor XI concentrations in serum and plasma.

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 Conjugate 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.

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

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

| Components                                  | 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       |
| Negative control (Depleted Human Plasma)    | 1 vial       |
| Positive control (Reference Plasma Control) | 1 vial       |
| Sealing Tapes                               | 1 x 3 units  |
| Stop Solution                               | 1 x 11ml     |

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

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

**Involvement in disease** 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.

**Sequence similarities** Belongs to the peptidase S1 family. Plasma kallikrein subfamily. Contains 4 apple domains.

Contains 1 peptidase S1 domain.

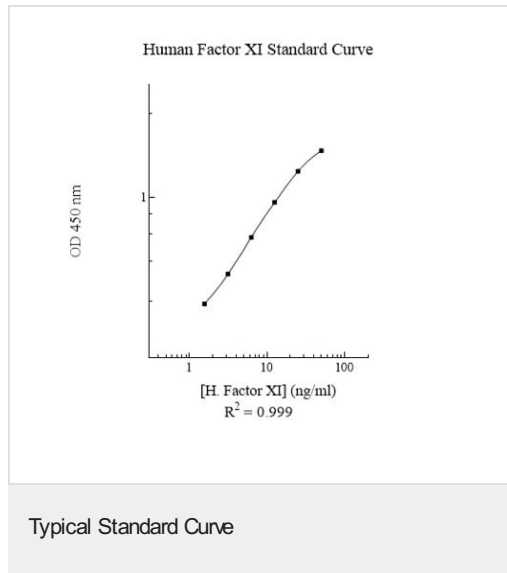
**Post-translational modifications**

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.

**Cellular localization**

Secreted.

**Images**



Representative Standard Curve using ab168547

**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.com/abpromise> or contact our technical team.

**Terms and conditions**

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors