

### Human FX ELISA kit (total FX antigen) ab272773

1 Image

#### Overview

**Product name** Human FX ELISA kit (total FX antigen)

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Sample 1	20	0.71ng/ml	0.039	5.44%
Sample 2	20	2.97ng/ml	0.155	5.23%
Sample 3	20	12.63ng/ml	0.095	2.45%

**Sample type** EDTA Plasma, Cit plasma

**Assay type** Quantitative

**Range** 0.1 ng/ml - 50 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Spike	98	% - %

**Assay duration** Multiple steps standard assay

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

**Product overview**

Human FX ELISA kit (total FX antigen) (ab272773) is intended for the quantitative determination of total Factor X antigen in human plasma.

Human Factor X will bind to the capture antibody coated on the microtiter plate. Factor X and Xa will react with the antibody on the plate. After appropriate washing steps, polyclonal anti-human Factor X primary antibody binds to the captured protein. Excess primary antibody is washed away and bound antibody, which is proportional to the total Factor X present in the samples, is reacted with the secondary antibody. Following an additional washing step, 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 human Factor X. Color development is proportional to the concentration of Factor X in the samples.

**Factor X standard provided is calibrated against the WHO 4th International Standard.**

**Platform** Microplate

## Properties

---

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Wash Buffer	1 x 50ml
Anti-Human FX Primary Antibody Lyophilized Vial	1 vial
Anti-Rabbit HRP Secondary Reagent	1 vial
FX ELISA Plate	1 unit
Human FX Standard Lyophilized Vial	1 vial
TMB Substrate	1 x 10ml

**Function** Factor Xa is a vitamin K-dependent glycoprotein that converts prothrombin to thrombin in the presence of factor Va, calcium and phospholipid during blood clotting.

**Tissue specificity** Plasma; synthesized in the liver.

**Involvement in disease** Defects in F10 are the cause of factor X deficiency (FA10D) [MIM:227600]. A hemorrhagic disease with variable presentation. Affected individuals can manifest prolonged nasal and mucosal hemorrhage, menorrhagia, hematuria, and occasionally hemarthrosis. Some patients do not have clinical bleeding diathesis.

**Sequence similarities** Belongs to the peptidase S1 family.  
Contains 2 EGF-like domains.  
Contains 1 Gla (gamma-carboxy-glutamate) domain.  
Contains 1 peptidase S1 domain.

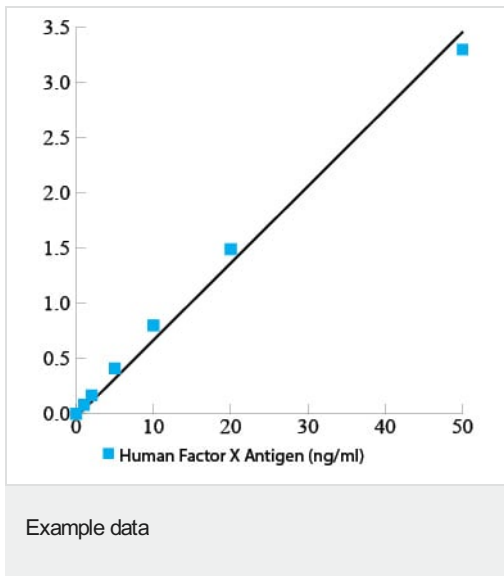
**Post-translational modifications** The vitamin K-dependent, enzymatic carboxylation of some glutamate residues allows the modified protein to bind calcium.  
N- and O-glycosylated.  
The activation peptide is cleaved by factor IXa (in the intrinsic pathway), or by factor VIIa (in the extrinsic pathway).  
The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.

**Cellular localization** Secreted.

---

## Images

---



A typical standard curve. Example only.

**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