

# Factor X Human Chromogenic Activity Assay Kit ab108833

[1 References](#) [5 Images](#)

### Overview

---

**Product name** Factor X Human Chromogenic Activity Assay Kit

**Detection method** Colorimetric

**Sample type** Cell culture supernatant, Plasma

**Assay type** Enzyme activity (quantitative)

**Sensitivity** = 0.009 µg/ml

**Range** 0.016 µg/ml - 4 µg/ml

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

**Product overview** Factor X Human Chromogenic Activity Assay Kit (ab108833) is developed to determine human Factor X activity in plasma and cell culture. The assay measures the activation of zymogen Factor X to Factor Xa by RVV. The amidolytic activity of the Factor Xa is quantitated using a highly specific Factor Xa substrate releasing a yellow para-nitroaniline (pNA) chromophore. The change in absorbance of the pNA at 405 nm is directly proportional to the Factor X enzymatic activity.

High level of FX activity samples: range of 0.125-8 µg/ml  
Low level of FX activity samples: range of 0.0156-1 µg/ml

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

**Notes** Factor X is a plasma serine protease zymogen involved in the blood coagulation cascade. Factor X is purified from plasma as a two-chain protein consisting of a 45-kDa heavy chain and a 17-kDa light chain. The Factor X heavy chain is cleaved during coagulation by several different proteases including the intrinsic Xase complex, the Factor X-activating enzyme from Russell's viper venom (RVV) and trypsin, and also by extrinsic (tissue Factor/Factor VIIa) pathway to give an active enzyme Factor Xa. Factor Xa as the activator of prothrombin occupies a central position linking the two blood coagulation pathways.

**Platform** Microplate reader

### Properties

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

Components	100 tests
Assay Diluent	1 x 5ml
Factor X Standard	1 unit
Factor Xa Substrate	2 vials
Microplate	1 unit
RVV	1 vial
Sample Diluent	1 x 5ml
Sealing Tapes	3 units

**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.

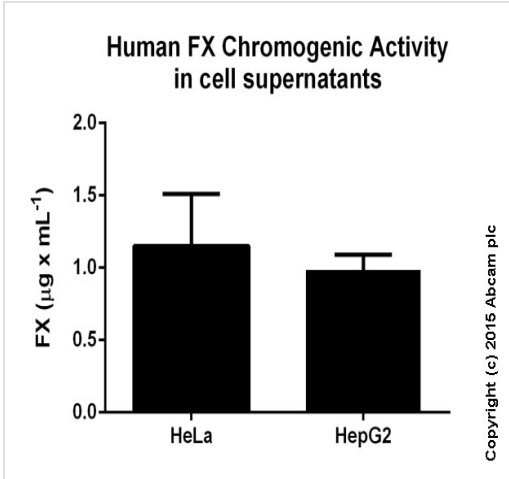
**Involvement in disease** Defects in F10 are the cause of factor X deficiency (FA10D) [MIM:227600]. A hemorrhagic disease with variable presentation. Affected individuals can manifest prolonged nasal and mucosal hemorrhage, menorrhagia, hematuria, and occasionally hemarthrosis. Some patients do not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

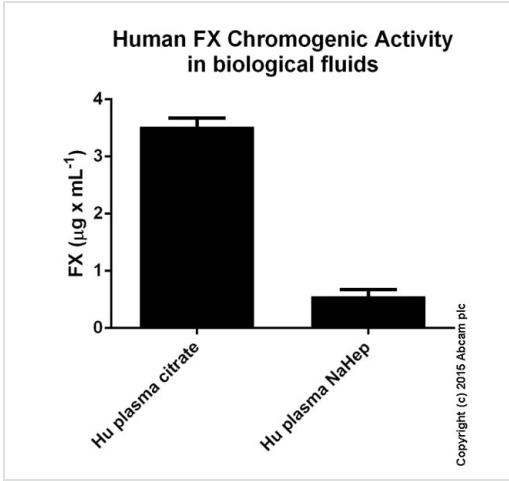
**Cellular localization** Secreted.

**Images**



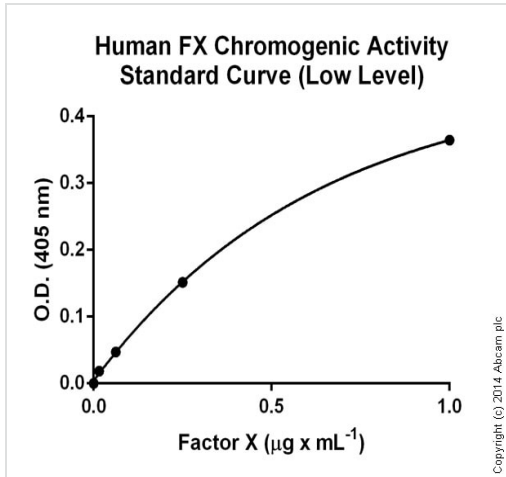
Factor X Chromogenic Activity measured in cell culture supernatants showing quantity (ug) per mL of tested sample after 40 min of incubation. Values calculated based on Low Level Standard Curve. Samples were diluted 2-10 fold.

Functional Studies - Factor X Human Chromogenic Activity Assay Kit (ab108833)



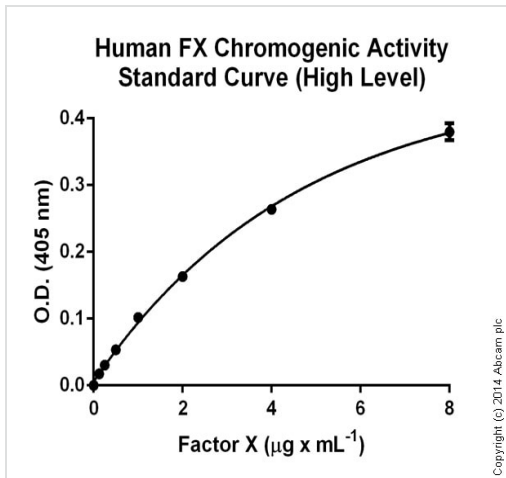
Factor X Chromogenic Activity measured in biological fluids showing quantity (ug) per mL of tested sample after 10 min of incubation. Values calculated based on High Level Standard Curve. Samples were diluted 2 fold.

Functional Studies - Factor X Human Chromogenic Activity Assay Kit (ab108833)



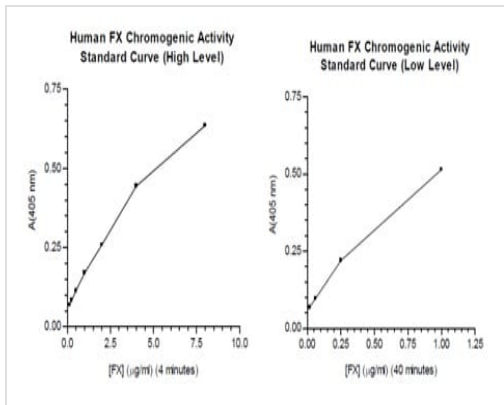
Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Functional Studies - Factor X Human Chromogenic Activity Assay Kit (ab108833)



Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Functional Studies - Factor X Human Chromogenic Activity Assay Kit (ab108833)



Functional Studies - Factor X Human Chromogenic Activity Assay Kit - 1 x 96 wells plate (ab108833)

**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