

## Product datasheet

### Mouse FX ELISA kit (total FX antigen) ab272779

[1 Image](#)

#### Overview

<b>Product name</b>	Mouse FX ELISA kit (total FX antigen)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Other biological fluids, Cit plasma
<b>Assay type</b>	Quantitative
<b>Range</b>	2.5 ng/ml - 500 ng/ml
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse
<b>Product overview</b>	Mouse FX ELISA kit (total FX antigen) (ab272779) is for the quantitative determination of total Factor X and Xa in biological fluids.

Mouse Factor X will bind to the affinity purified capture antibody coated on the microtiter plate. Factor X, Xa, and Xa in complex with inhibitors will react with the antibody on the plate. After appropriate washing steps, biotin labeled polyclonal anti-mouse Factor X primary antibody binds to the Factor X. Excess antibody is washed away and bound polyclonal antibody is then reacted with Streptavidin conjugated to horseradish peroxidase. Following an additional washing step, TMB substrate is used for color development at 450nm. The amount of color development is directly proportional to the concentration of total Factor X in the sample.

**Platform** Microplate

#### Properties

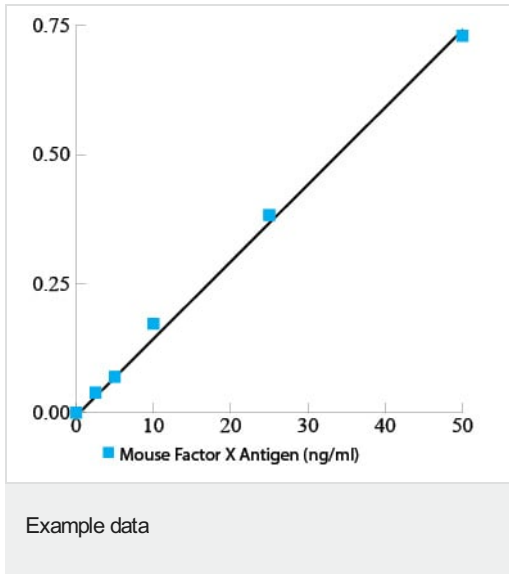
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Wash Buffer	1 x 50ml
Anti-Mouse FX Primary Antibody Lyophilized Vial	1 vial
Mouse FX ELISA Plate	1 unit
Mouse FX Standard Lyophilized Vial	1 vial

Components	1 x 96 tests
Streptavidin-HRP Secondary Reagent	1 vial
TMB Substrate	1 x 10ml

<b>Function</b>	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.
<b>Tissue specificity</b>	Plasma; synthesized in the liver.
<b>Involvement in disease</b>	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.
<b>Sequence similarities</b>	Belongs to the peptidase S1 family. Contains 2 EGF-like domains. Contains 1 Gla (gamma-carboxy-glutamate) domain. Contains 1 peptidase S1 domain.
<b>Post-translational modifications</b>	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.
<b>Cellular localization</b>	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