

Mouse FX ELISA kit (total FX antigen) ab272779

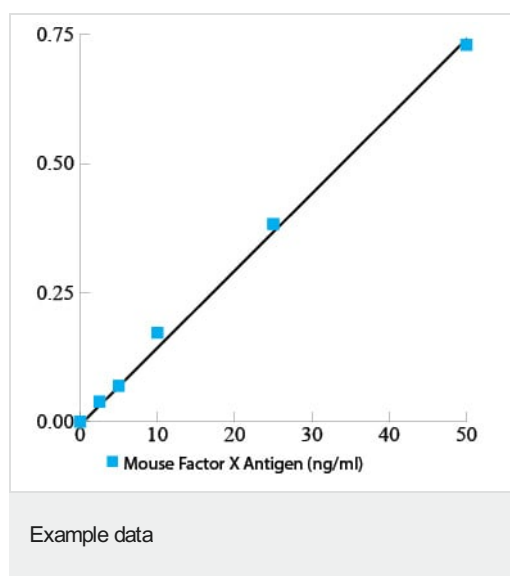
1 图像

概述											
产品名称	小鼠FX ELISA试剂盒(total FX antigen)										
检测方法	Colorimetric										
样品类型	Other biological fluids, Cit plasma										
检测类型	Quantitative										
范围	2.5 ng/ml - 500 ng/ml										
实验步骤	Multiple steps standard assay										
种属反应性	与反应: Mouse										
产品概述	<p>Mouse FX ELISA kit (total FX antigen) (ab272779) is for the quantitative determination of total Factor X and Xa in biological fluids.</p> <p>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.</p>										
平台	Microplate										
性能											
存放说明	Store at +4°C. Please refer to protocols.										
<table><tr><th>组件</th><th>1 x 96 tests</th></tr><tr><td>10X Wash Buffer</td><td>1 x 50ml</td></tr><tr><td>Anti-Mouse FX Primary Antibody Lyophilized Vial</td><td>1 vial</td></tr><tr><td>Mouse FX ELISA Plate</td><td>1 unit</td></tr><tr><td>Mouse FX Standard Lyophilized Vial</td><td>1 vial</td></tr></table>		组件	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
组件	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										

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

功能	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.
组织特异性	Plasma; synthesized in the liver.
疾病相关	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.
序列相似性	Belongs to the peptidase S1 family. Contains 2 EGF-like domains. Contains 1 Gla (gamma-carboxy-glutamate) domain. Contains 1 peptidase S1 domain.
翻译后修饰	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.
细胞定位	Secreted.

## 图片



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.cn/abpromise> or contact our technical team.

### **Terms and conditions**

---

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