

Human Factor IX ELISA Kit ab108831

4 References 3 Images

Overview

Product name	Human Factor IX ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			5.1%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			10.2%
Sample type	Cell culture supernatant, Serum, Plasma, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	0.4 ng/ml			
Range	1.56 ng/ml - 100 ng/ml			
Recovery	99 %			
Assay time	4h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Factor IX Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor IX concentrations in cell culture supernatants, serum and plasma.			

A Factor IX specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor IX 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 IX captured in plate.

Get higher sensitivity in only 90 minutes with Human Factor IX ELISA Kit ([\*\*ab188393\*\*](#)) from our SimpleStep ELISA® range.

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

**Platform** Microplate

## 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 IX Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Factor IX Microplate (12 x 8 well strips)	1 unit
Factor IX Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

**Function** Factor IX is a vitamin K-dependent plasma protein that participates in the intrinsic pathway of blood coagulation by converting factor X to its active form in the presence of Ca(2+) ions, phospholipids, and factor VIIIa.

**Tissue specificity** Synthesized primarily in the liver and secreted in plasma.

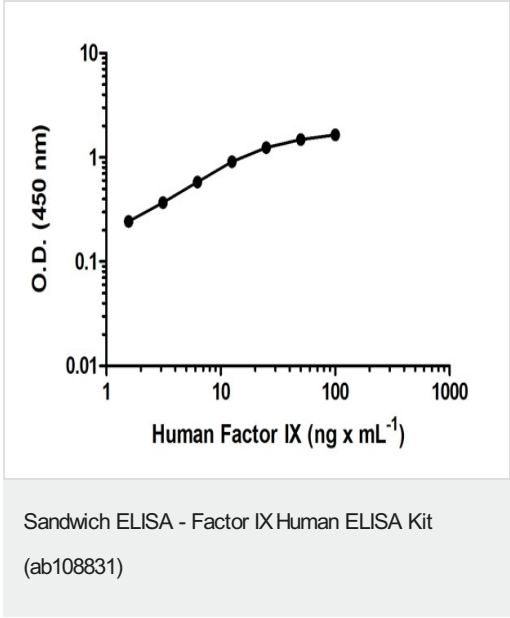
**Involvement in disease** Defects in F9 are the cause of recessive X-linked hemophilia B (HEMB) [MIM:306900]; also known as Christmas disease.  
Note=Mutations in position 43 (Oxford-3, San Dimas) and 46 (Cambridge) prevents cleavage of the propeptide, mutation in position 93 (Alabama) probably fails to bind to cell membranes, mutation in position 191 (Chapel-Hill) or in position 226 (Nagoya OR Hilo) prevent cleavage of the activation peptide.  
Defects in F9 are the cause of thrombophilia due to factor IX defect (THR-FIX) [MIM:300807]. A hemostatic disorder characterized by a tendency to thrombosis.

**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.

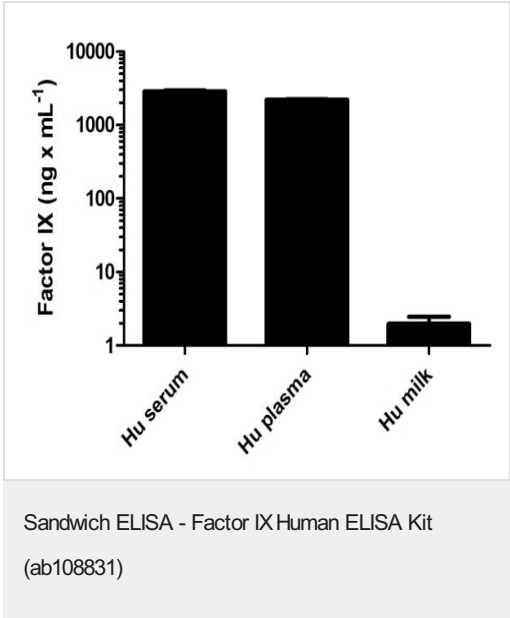
**Domain** Calcium binds to the gamma-carboxyglutamic acid (Gla) residues and, with stronger affinity, to

	another site, beyond the Gla domain.
<b>Post-translational modifications</b>	Activated by factor XIa, which excises the activation peptide. The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.
<b>Cellular localization</b>	Secreted.

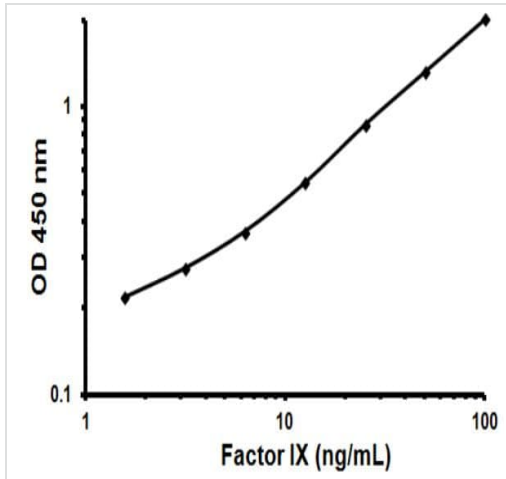
Images



Human Factor IX standard curve: mean of duplicates (+/- SD)



Human Factor IX measured in biological fluids showing quantity (ng) per mL of tested sample. Human serum and plasma were diluted 400-3200 fold. Human milk was diluted 1-8 fold.



Representative Standard Curve using ab108831

Typical Standard Curve

**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