

Human Factor IX/PTC ELISA Kit ab300307

重组 SimpleStep ELISA

★★★★★ 1 Abreviews 10 图像

概述

产品名称 人Factor IX/PTC ELISA试剂盒

检测方法 Colorimetric

精确度

批次内

样品	n	Mean	SD	CV%
Serum	8			7.4%

批次间

样品	n	Mean	SD	CV%
Serum	3			10.1%

样品类型 Serum, Hep Plasma, EDTA Plasma, Cit plasma

检测类型 Sandwich (quantitative)

灵敏度 15.884 pg/ml

范围 97.656 pg/ml - 6250 pg/ml

回收率

特定样本回收率

样品类型	平均%	范围
Serum	105	92% - 113%
Hep Plasma	112	103% - 118%
EDTA Plasma	106	95% - 115%
Cit plasma	116	111% - 119%

检测时间 1h 30m

实验步骤 One step assay

种属反应性 与反应: Human

## 产品概述

Human Factor IX/PTC SimpleStep ELISA® kit is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Factor IX/PTC protein in human serum and plasma. Quantitate Human Factor IX/PTC with 15.884 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

## 平台

Pre-coated microplate (12 x 8 well strips)

## 性能

### 存放说明

Store at +4°C. Please refer to protocols.

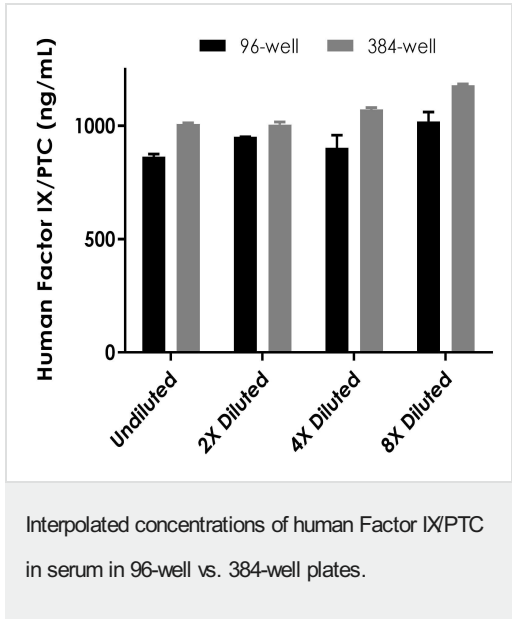
组件	10 x 96 tests	96 tests	1 x 384 tests
10X Human Factor IX/PTC Capture Antibody	1 x 6000µl	1 x 600µl	1 x 600µl
10X Human Factor IX/PTC Detector Antibody	1 x 6000µl	1 x 600µl	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 200ml	1 x 20ml	1 x 20ml
384 well CaptSure™ microplates	0 x 0 unit	0 x 0 unit	1 unit
Antibody Diluent 4BI	10 x 6ml	1 x 6ml	1 x 6ml
Human Factor IX/PTC Lyophilized Recombinant Protein	2 x 10 vials	2 vials	2 vials
Plate Seals	1 x 10 units	1 unit	1 unit
Sample Diluent NS (ab193972)	2 x 250ml	1 x 50ml	2 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 x 10 units	1 unit	0 x 0 unit
Stop Solution	1 x 120ml	1 x 12ml	2 x 12ml
TMB Development Solution	1 x 120ml	1 x 12ml	2 x 12ml

## 功能

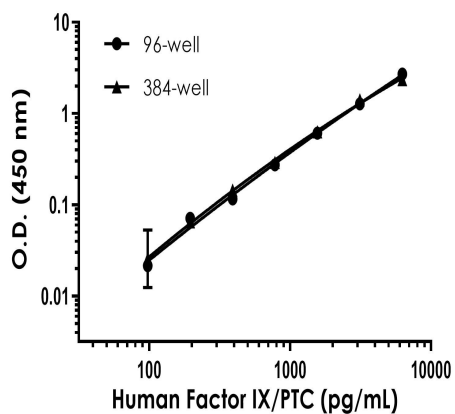
Factor IX is a vitamin K-dependent plasma protein that participates in the intrinsic pathway of blood coagulation by converting factor X to its active form in the presence of Ca(2+) ions,

	phospholipids, and factor VIIIa.
组织特异性	Synthesized primarily in the liver and secreted in plasma.
疾病相关	<p>Defects in F9 are the cause of recessive X-linked hemophilia B (HEMB) [MIM:306900]; also known as Christmas disease.</p> <p>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.</p> <p>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.</p>
序列相似性	<p>Belongs to the peptidase S1 family.</p> <p>Contains 2 EGF-like domains.</p> <p>Contains 1 Gla (gamma-carboxy-glutamate) domain.</p> <p>Contains 1 peptidase S1 domain.</p>
结构域	Calcium binds to the gamma-carboxyglutamic acid (Gla) residues and, with stronger affinity, to another site, beyond the Gla domain.
翻译后修饰	<p>Activated by factor XIa, which excises the activation peptide.</p> <p>The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.</p>
细胞定位	Secreted.

图片

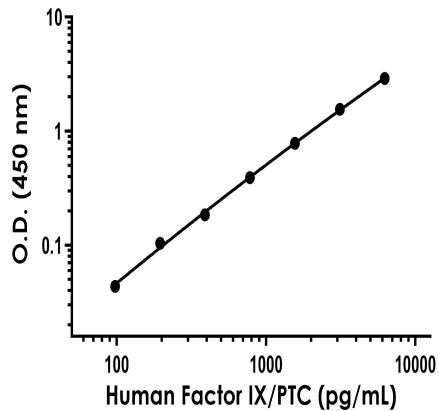


Interpolated concentration of native Factor IX/PTC was measured in duplicate at different sample concentrations in 96-well vs. 384-well plates. Undiluted samples are 1:400 serum. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). Sample dilutions are made in Sample Diluent NS.



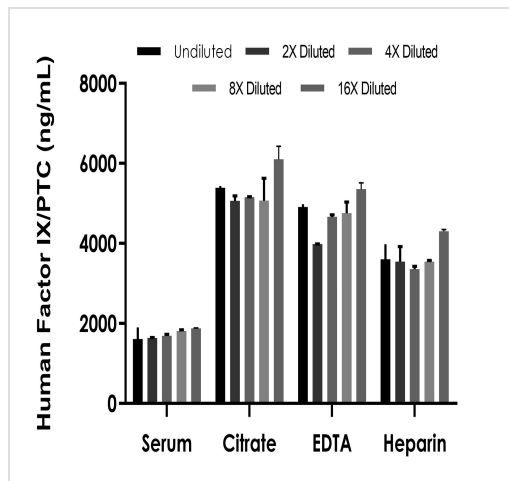
Example of human Factor IX/PTC standard curve in 96-well vs. 384-well plate. Background-subtracted data values (mean  $\pm$  SD) are graphed.

Example of human Factor IX/PTC standard curve in Sample Diluent NS in 96-well vs. 384-well plate.



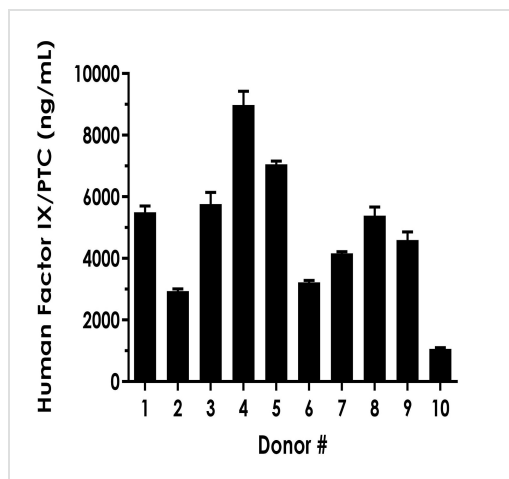
Example of human Factor IX/PTC standard curve. Background-subtracted data values (mean  $\pm$  SD) are graphed.

Example of human Factor IX/PTC standard curve in Sample Diluent NS.



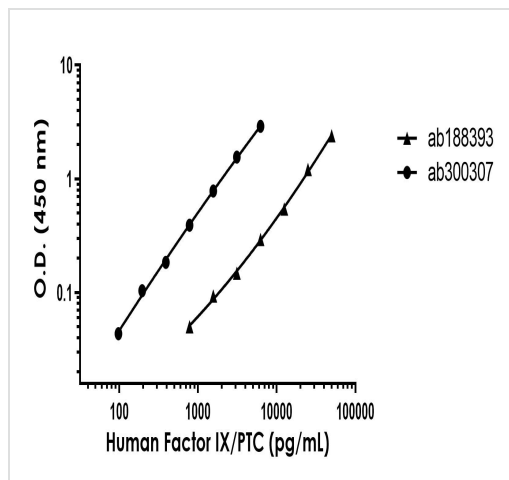
Interpolated concentrations of human Factor IX/PTC in serum, plasma (citrate), plasma (EDTA), and plasma (heparin).

Interpolated concentration of native Factor IX/PTC was measured in duplicate at different sample concentrations. Undiluted samples are as follows: serum 0.15%, plasma (citrate) 0.075%, plasma (EDTA) 0.075%, and plasma (heparin) 0.15%. The interpolated dilution factor corrected values are plotted (mean  $\pm$  SD,  $n=2$ ). Sample dilutions are made in Sample Diluent NS.



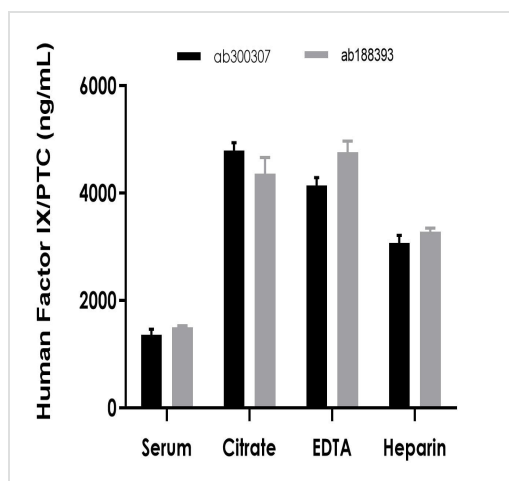
Interpolated concentrations of Factor IX/PTC in normal human serum donors.

Serum of ten individual healthy human female donors was measured in duplicate. Interpolated dilution factor corrected values are plotted (mean  $\pm$  SD,  $n=2$ ). The mean Factor IX/PTC concentration was determined to be 4,867.18 ng/mL with a range of 1,062.16 – 8,981.25 ng/mL.



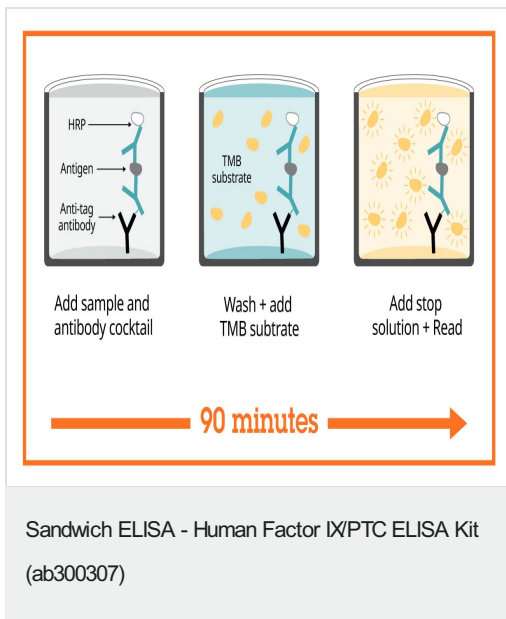
Standard curve comparison between the original Human Factor IX/PTC SimpleStep ELISA (**ab188393**) and current Human Factor IX/PTC SimpleStep ELISA (ab300307).

Human Factor IX/PTC standard curve comparison.

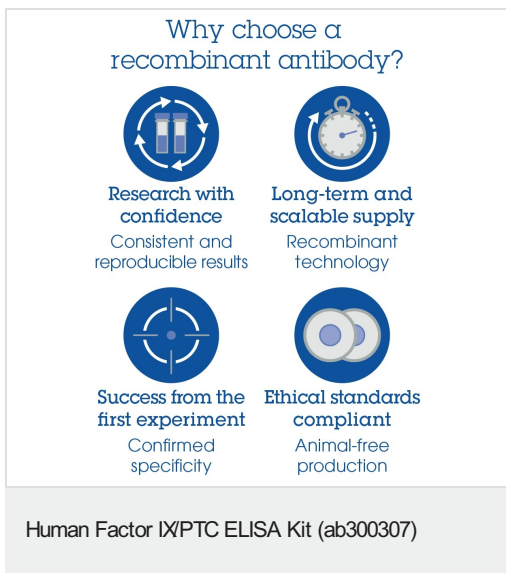


Serum and plasma comparison between the original Human Factor IX/PTC SimpleStep ELISA (**ab188393**) and current Human Factor IX/PTC SimpleStep ELISA (ab300307).

Human Factor IX/PTC serum and plasma comparison.



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



**Get more done with  
SimpleStep ELISA**



**Easy to use**  
Single-wash 90-minute  
protocol



**Flexible**  
Matched antibody pairs  
available



**Precision antibodies**  
High sensitivity, specificity  
and reproducibility



**Scalable**  
Now in 10-pack and  
384-well formats

Sandwich ELISA - Human Factor IX/PTC ELISA Kit  
(ab300307)

To learn more about the advantages of SimpleStep ELISA® kits  
see [here](#).

**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