

Human Tissue Factor Pathway Inhibitor ELISA Kit (TFPI) ab108904

6 References 2 图像

概述

产品名称	人组织Factor Pathway抑制剂ELISA试剂盒(TFPI)																				
检测方法	Colorimetric																				
精确度	<div>批次内</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>5%</td></tr></table> <div>批次间</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>9.9%</td></tr></table>	样品	n	Mean	SD	CV%	Overall				5%	样品	n	Mean	SD	CV%	Overall				9.9%
样品	n	Mean	SD	CV%																	
Overall				5%																	
样品	n	Mean	SD	CV%																	
Overall				9.9%																	
样品类型	Cell culture supernatant, Milk, Serum, Plasma, Cerebral Spinal Fluid																				
检测类型	Sandwich (quantitative)																				
灵敏度	> 0.081 ng/ml																				
范围	0.625 ng/ml - 5 ng/ml																				
回收率	98 %																				
检测时间	5h 00m																				
实验步骤	Multiple steps standard assay																				
种属反应性	与反应: Human																				
产品概述	Abcam's Tissue Factor Pathway Inhibitor (TFPI) Human <i>in vitro</i> ELISA (Enzyme-Linked																				

changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Tissue Factor Pathway Inhibitor captured in plate.

Get results in 90 minutes with Human Tissue Factor Pathway Inhibitor ELISA Kit ([ab188392](#)) 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.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Tissue Factor Pathway Inhibitor Antibody	1 x 120µl
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml
Tissue Factor Pathway Inhibitor Microplate (12 x 8 well strips)	1 x 96 units
Tissue Factor Pathway Inhibitor Standard	2 vials

功能 Inhibits factor X (X(a)) directly and, in a Xa-dependent way, inhibits VIIa/tissue factor activity, presumably by forming a quaternary Xa/LACI/VIIa/TF complex. It possesses an antithrombotic action and also the ability to associate with lipoproteins in plasma.

组织特异性 Mostly in endothelial cells.

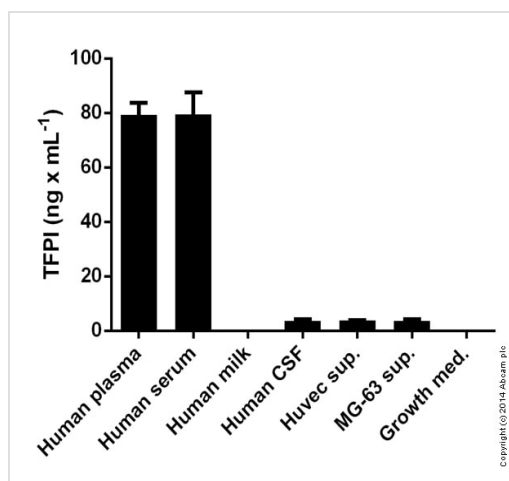
序列相似性 Contains 3 BPTI/Kunitz inhibitor domains.

结构域 This inhibitor contains three inhibitory domains. The first domain interacts with VIIa and TF, the second one with Xa.

翻译后修饰 O-glycosylated.

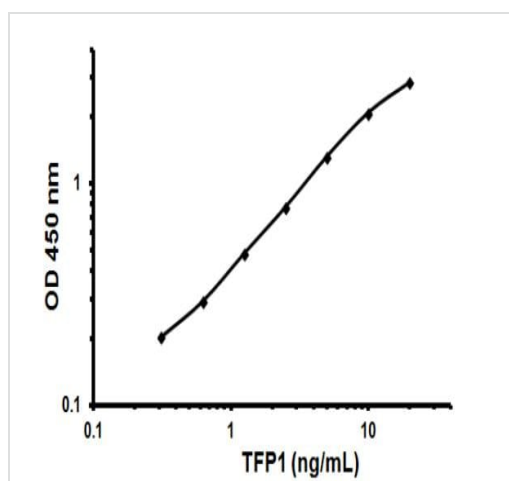
细胞定位 Secreted.

图片



TFPI measured in various samples showing quantity (ng) per mL of tested sample

Sandwich ELISA - ab108904 Tissue Factor Pathway Inhibitor (TFPI) Human ELISA Kit



Representative Standard Curve using ab108904.

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.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