

Human PAI1 + tPA ELISA Kit ab108892

2 References 1 图像

概述

产品名称	人PAI1 + tPA ELISA试剂盒																				
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>4.1%</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>10.4%</td></tr></table>	样品	n	Mean	SD	CV%	Overall				4.1%	样品	n	Mean	SD	CV%	Overall				10.4%
样品	n	Mean	SD	CV%																	
Overall				4.1%																	
样品	n	Mean	SD	CV%																	
Overall				10.4%																	
样品类型	Cell culture supernatant, Serum, Plasma																				
检测类型	Sandwich (quantitative)																				
灵敏度	= 0.039 ng/ml																				
范围	0.25 ng/ml - 2 ng/ml																				
回收率	98 %																				
检测时间	4h 00m																				
实验步骤	Multiple steps standard assay																				
种属反应性	与反应: Human																				
产品概述	Abcam's PAI1 + tPA Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is																				

A PAI1 + tPA specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a PAI1 + tPA 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 PAI1 + tPA captured in plate.

Get results in 90 minutes with Human PAI1+tPA ELISA Kit ([ab192151](#)) 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

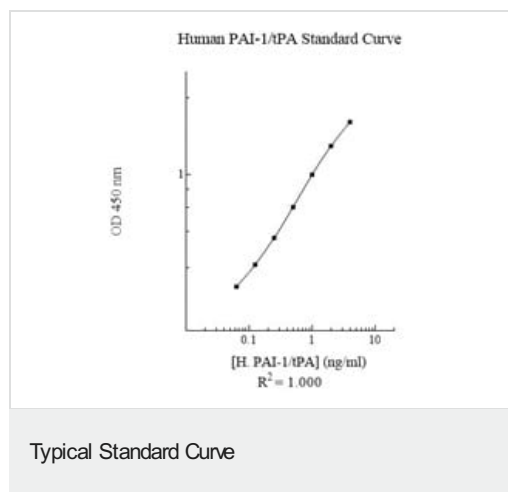
## 性能

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 M Concentrate	1 x 20ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human PAI1 + tPA Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
PAI1 + tPA Microplate (12 x 8 well strips)	1 unit
PAI1 + tPA Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

细胞定位 Secreted

## 图片



Representative Standard Curve using ab108892.

**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