

Product datasheet

Human Annexin V ELISA Kit ab119503

1 图像

概述

产品名称 人Annexin V ELISA试剂盒

检测方法 Colorimetric

精确度

批次内				
样品	n	Mean	SD	CV%
Overall				< 10%

批次间				
样品	n	Mean	SD	CV%
Overall				11%

样品类型 Cell culture supernatant

检测类型 Sandwich (quantitative)

灵敏度 0.33 ng/ml

范围 0.78 ng/ml - 50 ng/ml

复检 85 %

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Annexin V Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human Annexin V concentrations in cell culture supernatant and serum.

Annexin V specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotinylated Annexin V detection antibody then incubated at room temperature and washed. A Streptavidin-HRP conjugate is then added to each well, incubated at room temperature and washed. TMB is added and catalyzed by HRP to produce a blue color product that changes into yellow after addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of Annexin V sample captured in plate.

经测试应用 适用于: Sandwich ELISA

平台 Microplate

性能

存放说明

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

组件	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-human Annexin V monoclonal antibody	1 x 70µl
Human Annexin V Standard	2 vials
Microplate coated with monoclonal antibody to Human Annexin V (12 x 8 wells)	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

功能

This protein is an anticoagulant protein that acts as an indirect inhibitor of the thromboplastin-specific complex, which is involved in the blood coagulation cascade.

疾病相关

Pregnancy loss, recurrent, 3

序列相似性

Belongs to the annexin family.
Contains 4 annexin repeats.

结构域

The [IL]-x-C-x-x-[DE] motif is a proposed target motif for cysteine S-nitrosylation mediated by the iNOS-S100A8/A9 transnitrosylase complex.
A pair of annexin repeats may form one binding site for calcium and phospholipid.

翻译后修饰

S-nitrosylation is induced by interferon-gamma and oxidatively-modified low-density lipoprotein (LDL(ox)) possibly implicating the iNOS-S100A8/9 transnitrosylase complex.

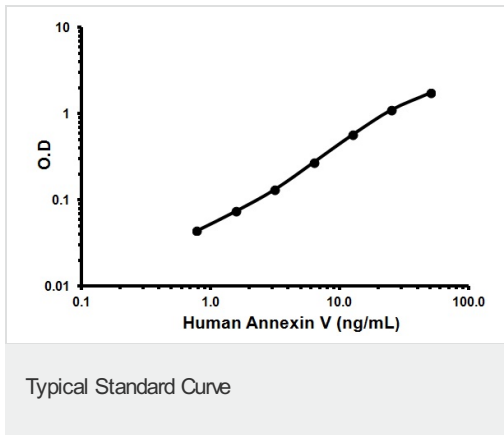
应用

Our [Abpromise guarantee](#) covers the use of **ab119503** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



Representative Standard Curve using
ab119503.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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