

Human Kininogen ELISA Kit (KNG) ab108876

[3 References](#) [1 图像](#)

概述

产品名称 人Kininogen ELISA试剂盒(KNG)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				3.5%

批次间

样品	n	Mean	SD	CV%
Overall				9.1%

样品类型 Cell culture supernatant, Saliva, Milk, Urine

检测类型 Sandwich (quantitative)

灵敏度 = 0.7 ng/ml

范围 1.563 ng/ml - 100 ng/ml

回收率 96 %

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 Kininogen (KNG) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Kininogen in urine, saliva, milk, cerebrospinal fluid and cell culture supernatants.

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

Get results in 90 minutes with Human HMW Kininogen ELISA Kit ([ab189574](#)) 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 Kininogen Antibody	1 x 120µl
Chromogen Substrate	1 x 8ml
Kininogen Microplate (12 x 8 well strips)	1 unit
Kininogen Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 12ml

功能

(1) Kininogens are inhibitors of thiol proteases; (2) HMW-kininogen plays an important role in blood coagulation by helping to position optimally prekallikrein and factor XI next to factor XII; (3) HMW-kininogen inhibits the thrombin- and plasmin-induced aggregation of thrombocytes; (4) the active peptide bradykinin that is released from HMW-kininogen shows a variety of physiological effects: (4A) influence in smooth muscle contraction, (4B) induction of hypotension, (4C) natriuresis and diuresis, (4D) decrease in blood glucose level, (4E) it is a mediator of inflammation and causes (4E1) increase in vascular permeability, (4E2) stimulation of nociceptors (4E3) release of other mediators of inflammation (e.g. prostaglandins), (4F) it has a cardioprotective effect (directly via bradykinin action, indirectly via endothelium-derived relaxing factor action); (5) LMW-kininogen inhibits the aggregation of thrombocytes; (6) LMW-kininogen is in contrast to HMW-kininogen not involved in blood clotting.

组织特异性

Secreted in plasma. T-kinin is detected in malignant ovarian, colon and breast carcinomas, but not in benign tumors.

疾病相关

Defects in KNG1 are the cause of high molecular weight kininogen deficiency (HMWK deficiency) [MIM:228960]. HMWK deficiency is an autosomal recessive coagulation defect. Patients with HMWK deficiency do not have a hemorrhagic tendency, but they exhibit abnormal surface-mediated activation of fibrinolysis.

序列相似性

Contains 3 cystatin domains.

翻译后修饰

Bradykinin is released from kininogen by plasma kallikrein.

Hydroxylation of Pro-383 occurs prior to the release of bradykinin.

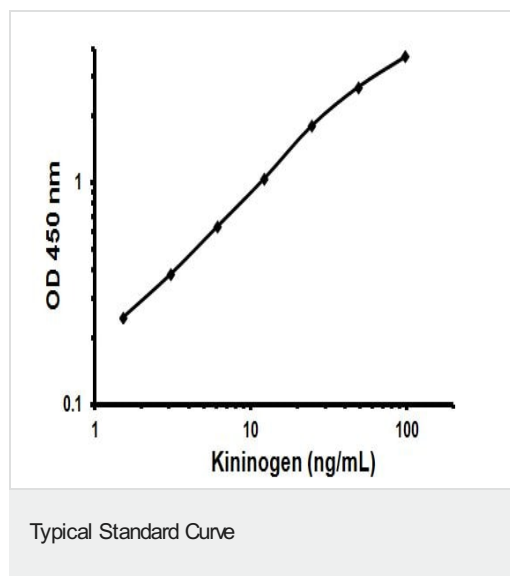
Phosphorylation sites are present in the extracellular medium.

N- and O-glycosylated. O-glycosylated with core 1 or possibly core 8 glycans.

细胞定位

Secreted > extracellular space.

图片



Representative Standard Curve Using ab108876.

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