

Human Kallikrein 12 ELISA Kit ab213802

1 图像

概述

产品名称 人Kallikrein 12 ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
1	16	1409pg/ml	73.26	5.2%
2	16	3882pg/ml	213.51	5.5%
3	16	6756pg/ml	405.36	6%

批次间

样品	n	Mean	SD	CV%
1	24	1523pg/ml	89.85	5.9%
2	24	4129pg/ml	251.86	6.1%
3	24	6526pg/ml	443.76	6.8%

样品类型 Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate

检测类型 Sandwich (quantitative)

灵敏度 < 10 pg/ml

范围 156 pg/ml - 10000 pg/ml

检测时间 3h 30m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 The Human Kallikrein 12 (KLK12) Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213802) is designed for the quantitative measurement of Human KLK12 in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for KLK12 has been pre-coated onto 96-well plates. Standards and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for KLK12 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human KLK12 amount of sample captured in plate.

说明

Kallikrein-12 is a protein that in humans is encoded by the KLK12 gene. It belongs to the kallikrein subgroup of serine proteases, which have diverse physiologic functions in many tissues. KLK12 is expressed primarily in salivary gland, stomach, uterus, trachea, prostate, thymus, lung, colon, brain, breast, and thyroid gland and at lower levels in testis, pancreas, small intestine, and spinal cord. The expression of KLK12 is downregulated at the mRNA level in breast cancer tissues and is upregulated by steroid hormones in breast and prostate cancer cell lines.

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

Store at -20°C. Please refer to protocols.

组件	标识符	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human KLK12 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated KLK12 Antibody		1 x 100µl
Human KLK12 Standard (Lyophilized)		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

相关性

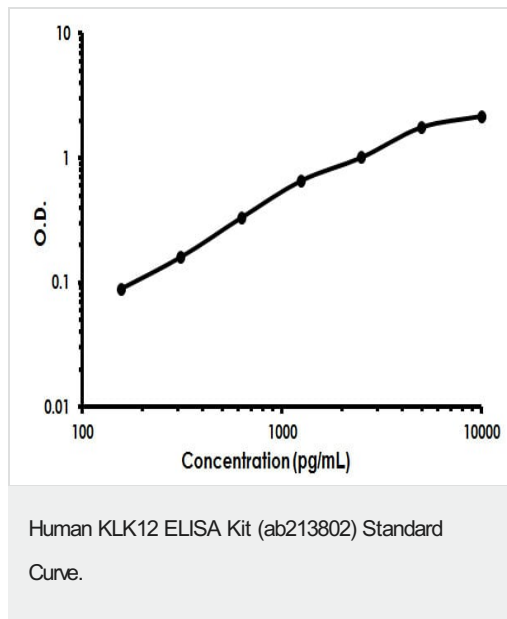
The human tissue Kallikrein gene family encodes 15 serine proteases. All Kallikreins share structural similarities including cysteine residues, a catalytic triad of His, Asp, and Ser residues, typically five coding exons, and varied intron phases. Kallikreins are predominantly secreted as inactive zymogens prior to activation by cleavage of an N-terminal peptide, and all function extracellularly. Kallikreins can be activated autocatalytically, via other Kallikreins, or additional proteases. While structurally similar, Kallikrein family members have distinct functions and have key roles in many physiological and pathological processes. Many human tissue Kallikreins also

show promise as cancer biomarkers, which may facilitate earlier detection and characterization of many forms of cancer. Kallikrein 12, also known as Kallikrein-like5 (KLK-L5), is a trypsin-like serine proteinase. It was discovered by cloning the kallikrein locus on chromosome 19, and was found to be similar to the other kallikrein family genes. It is produced by a number of tissues, including pancreas, salivary gland, breast, prostate, uterus, lung, thymus, thyroid, brain, stomach and colon. Kallikrein 12 message has been reported to be decreased in breast cancer. Different splice variants of kallikrein-12 have been identified; a 254 amino acid form, 248 amino acids form, and a truncated 111 amino acid form.

细胞定位

Secreted

图片



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