

Human Kallikrein 1 ELISA Kit (KLK1) ab119598

1 Image

Overview

Product name	Human Kallikrein 1 ELISA Kit (KLK1)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma, Cell Lysate, Tissue Homogenate
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	156 pg/ml - 10000 pg/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's Human KLK1 <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human KLK1 in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).</p> <p>A KLK1 specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat specific for KLK1 is then 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. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP 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 Human KLK1 amount of sample captured in plate.</p>
Platform	Microplate

Properties

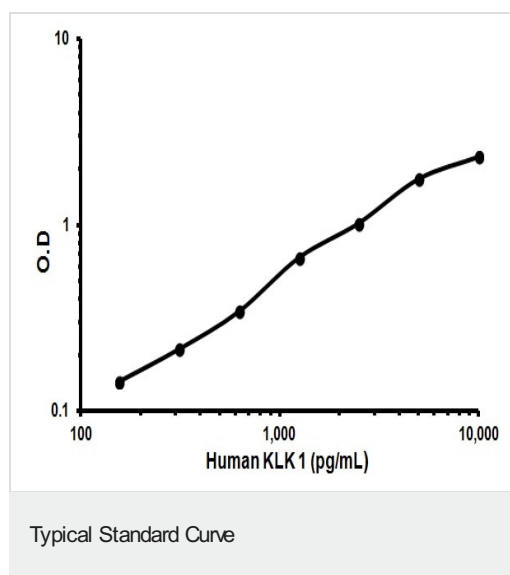
Storage instructionsStore at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Kallikrein 1 antibody Microplate (12 x 8 wells)		1 unit	1 unit

Components	Identifier	1 x 96 tests	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human Kallikrein 1 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Kallikrein 1 standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function	Glandular kallikreins cleave Met-Lys and Arg-Ser bonds in kininogen to release Lys-bradykinin.
Tissue specificity	Isoform 2 is expressed in pancreas, salivary glands, kidney, colon, prostate gland, testis, spleen and the colon adenocarcinoma cell line T84.
Sequence similarities	Belongs to the peptidase S1 family. Kallikrein subfamily. Contains 1 peptidase S1 domain.
Post-translational modifications	The O-linked polysaccharides on Ser-93, Ser-104 and Ser-167 are probably the mucin type linked to GalNAc. In PubMed:3163150, GalNAc was detected with the corresponding peptides but not located.

Images



Representative Standard Curve using ab119598

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors