

Product datasheet

Human Prostatin ELISA Kit ab213817

[1 Image](#)

Overview

Product name Human Prostatin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	3.254ng/ml	0.247	= 7.6%
2	16	12.396ng/ml	0.595	= 4.8%
3	16	41.368ng/ml	1.944	= 4.7%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	3.115ng/ml	0.267	= 8.6%
2	24	13.251ng/ml	0.689	= 5.2%
3	24	40.618ng/ml	2.599	= 6.4%

Sample type Cell culture supernatant, Urine, Serum, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 50 pg/ml

Range 1.56 ng/ml - 100 ng/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview The Human Prostatin Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213817) is designed for the quantitative measurement of Human Prostatin in cell culture supernatants, serum, plasma (heparin, EDTA) and urine.

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

polyclonal antibody from sheep specific for Prostasin has been pre-coated onto 96-well plates. Standards (Expression system for standard: CHO; Immunogen sequence: A33-G319) and test samples are added to the wells, a biotinylated detection polyclonal antibody from sheep specific for Prostasin 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 Prostasin amount of sample captured in plate.

Notes Prostasin is a protein that in humans is encoded by the PRSS8 gene. It is a single-copy gene and mapped to 16p11.2. Prostasin is likely to be the more physiologically relevant protease. This gene encodes a trypsinogen, which is a member of the trypsin family of serine proteases. Co-expression of Prostasin with either Xenopus or rat ENaC in Xenopus oocytes resulted in a 60% to 80% increase in amiloride-sensitive sodium currents, and that the addition of aprotinin, a serine protease inhibitor, completely prevented this activation. This enzyme is highly expressed in prostate epithelia and is one of several proteolytic enzymes found in seminal fluid. The proprotein is cleaved to produce a light chain and a heavy chain which are associated by a disulfide bond. It is active on peptide linkages involving the carboxyl group of lysine or arginine.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	Identifier	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 Prostasin coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human Prostasin antibody		1 x 130µl
Lyophilized recombinant Human Prostasin standard		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

Function Possesses a trypsin-like cleavage specificity.

Tissue specificity Found in prostate, liver, salivary gland, kidney, lung, pancreas, colon, bronchus and renal proximal tubular cells. In the prostate gland it may be synthesized in epithelial cells, secreted into the ducts, and excreted into the seminal fluid.

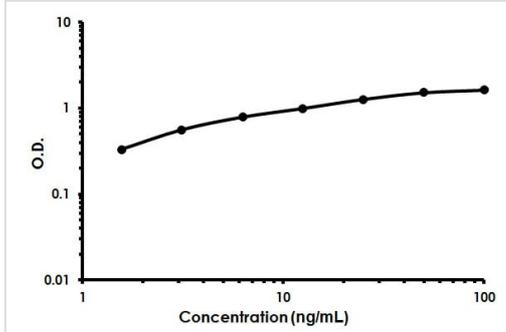
Sequence similarities Belongs to the peptidase S1 family.

Contains 1 peptidase S1 domain.

Cellular localization

Cell membrane and Secreted > extracellular space. Found in the seminal fluid. Secreted after cleavage of its C-terminus.

Images



Human Prostasin ELISA Kit (ab213817) Standard Curve.

Human Prostasin ELISA Kit (ab213817) Standard Curve.

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