

Product datasheet

Human CRISP3 ELISA Kit ab213768

[1 References](#) [1 Image](#)

Overview

Product name Human CRISP3 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	110pg/ml	4.51	= 4.1%
2	16	383pg/ml	23.75	= 6.2%
3	16	1477pg/ml	66.47	= 4.5%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	119pg/ml	5.59	= 4.7%
2	24	408pg/ml	26.11	= 6.4%
3	24	1764pg/ml	95.26	= 5.4%

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

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 62.5 pg/ml - 4000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview The Human CRISP3 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213768) is designed for the quantitative measurement of Human CRISP3 in cell culture supernatants, serum, plasma (heparin, EDTA), saliva and urine.

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

monoclonal antibody from mouse specific for CRISP3 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: N21-Y245) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for CRISP3 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 CRISP3 amount of sample captured in plate.

Notes Cysteine-rich secretory protein 3 is a cysteine-rich secretory protein that in humans is encoded by the CRISP3 gene. CRISP3 expression is found in the secretory epithelium throughout the male genital tract, with particularly high expression in the cauda epididymis and ampulla vas deferens. It is a potential biomarker for prostate cancer. CRISP3 gene can be defined as an early-transcribed gene with a putative role in the pathophysiology of sjogren's syndrome. CRISP3 is present in exocrine secretions and in secretory granules of neutrophilic granulocytes and is believed to play a role in innate immunity.

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

Properties

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

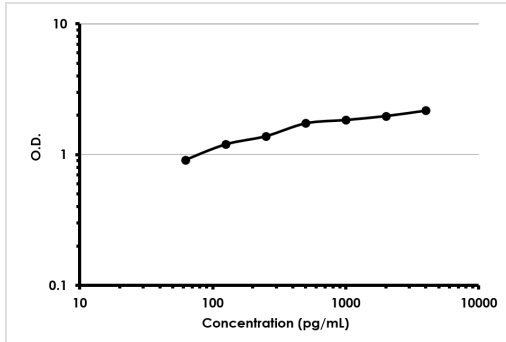
Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human CRISP3 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti- Human CRISP3 antibody		1 x 130µl
Lyophilized recombinant Human CRISP3 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

Tissue specificity Salivary gland, pancreas and prostate > epididymis, ovary, thymus and colon.

Sequence similarities Belongs to the CRISP family.

Cellular localization Secreted. In neutrophils, localized in specific granules.

Images



Human CRISP3 ELISA Kit (ab213768) Standard Curve.

Human CRISP3 ELISA Kit (ab213768) 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