

Human Cystatin C ELISA Kit ab119589

1 References 1 图像

概述

产品名称	人Cystatin C ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	1	16	0.687ng/ml	0.044	= 6.4%
	2	16	2.837ng/ml	0.207	= 7.3%
	3	16	10.374ng/ml	0.684	= 6.6%
	批次间				
	样品	n	Mean	SD	CV%
	1	24	0.647ng/ml	0.051	= 7.9%
	2	24	2.715ng/ml	0.217	= 8%
	3	24	10.112ng/ml	0.859	= 8.5%
样品类型	Cell culture supernatant, Milk, Urine, Serum, Hep Plasma, EDTA Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 10 pg/ml				
范围	0.312 ng/ml - 20 ng/ml				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's Human Cystatin C <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human Cystatin C in cell culture supernatants, serum, plasma (heparin, EDTA) saliva, urine and Human milk.				

A Cystatin C 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 Cystatin C is then added 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 Cystatin C amount of sample captured in plate.

Get results in 90 minutes with Human Cystatin C ELISA Kit ([ab179883](#)) from our SimpleStep ELISA® range.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	标识符	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 Cystatin C antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Human Cystatin C antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Cystatin C standard		2 x 20ng	2 x 20ng
Plate Seal		4 units	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

功能 As an inhibitor of cysteine proteinases, this protein is thought to serve an important physiological role as a local regulator of this enzyme activity.

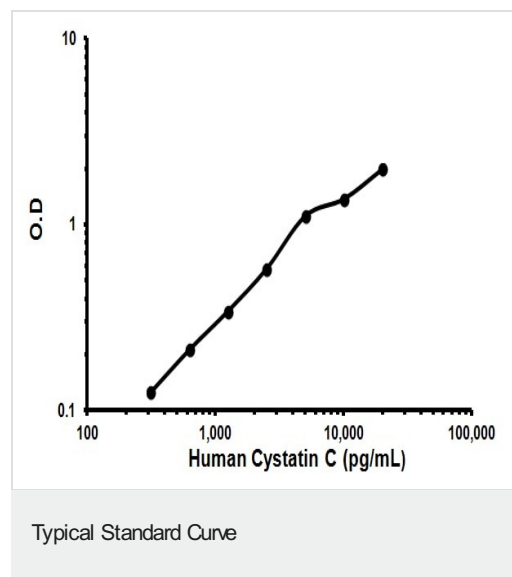
组织特异性 Expressed in submandibular and sublingual saliva but not in parotid saliva (at protein level). Expressed in various body fluids, such as the cerebrospinal fluid and plasma. Expressed in highest levels in the epididymis, vas deferens, brain, thymus, and ovary and the lowest in the submandibular gland.

疾病相关 Amyloidosis 6
Macular degeneration, age-related, 11

序列相似性 Belongs to the cystatin family.

翻译后修饰 The Thr-25 variant is O-glycosylated with a core 1 or possibly core 8 glycan. The signal peptide of the O-glycosylated Thr-25 variant is cleaved between Ala-20 and Val-21.

图片



Representative Standard Curve using 119589.

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