

Human Cystatin C ELISA Kit ab119589

1 References 画像数 1

製品の概要

製品名	Human Cystatin C ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	1	16	0.687ng/ml	0.044
	2	16	2.837ng/ml	0.207
	3	16	10.374ng/ml	0.684
				CV%
				= 6.4%
				= 7.3%
				= 6.6%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	1	24	0.647ng/ml	0.051
	2	24	2.715ng/ml	0.217
	3	24	10.112ng/ml	0.859
				CV%
				= 7.9%
				= 8%
				= 8.5%
サンプルの種類	Cell culture supernatant, Milk, Urine, Serum, Hep Plasma, EDTA Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 10 pg/ml			
検出範囲	0.312 ng/ml - 20 ng/ml			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	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.

#### 試験プラットフォーム

Microplate

#### 製品の特長

#### 保存方法

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

内容	ラベル	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

#### 機能

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

#### 組織特異性

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.

#### 関連疾患

Amyloidosis 6  
Macular degeneration, age-related, 11

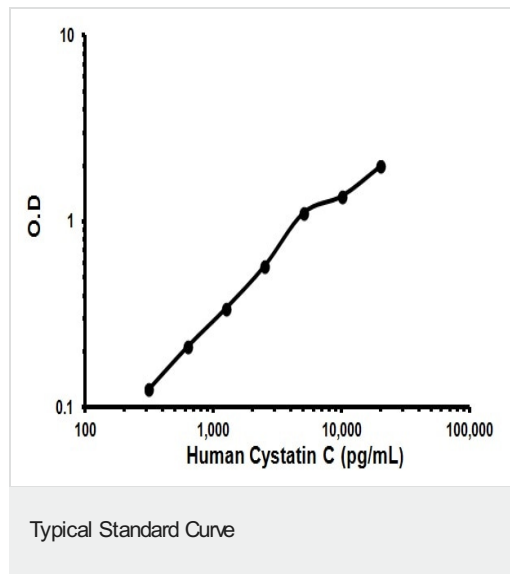
#### 配列類似性

Belongs to the cystatin family.

#### 翻訳後修飾

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.

## 画像



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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

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