

Product datasheet

Monkey Cystatin C ELISA Kit ab190542

[1 Image](#)

Overview

Product name Monkey Cystatin C ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Serum				< 10%

Sample type Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.33 ng/ml

Range 0.125 ng/ml - 4 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	% - %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Monkey

Product overview

Abcam's Cystatin C Monkey ELISA Kit (ab190542) is for the measurement of Cystatin C in monkey serum and plasma.

In this assay the Cystatin C present in samples reacts with the anti-Cystatin C antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, the Detection Antibody, biotin conjugated anti-Cystatin C, is added and complexes are formed. Following a wash step, the horseradish peroxidase (HRP) conjugated Streptavidin is added and complexes are formed. After another washing step, the complexes are assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The

quantity of bound enzyme varies directly with the concentration of Cystatin C in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Cystatin C in the test sample. The quantity of Cystatin C in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Notes Cystatin C is a small cysteine proteinase inhibitor present in body fluids. Studies have shown Cystatin C levels to be directly correlated with the glomerular filtration rate.

The calibrator used in this kit is of human origin.

Platform Microplate

Properties

Components	1 x 96 tests
1X Diluent Concentrate	1 x 60ml
20X Wash Buffer Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Cystatin C Monkey Antibody coated microwells	1 x 96 tests
Cystatin C Monkey Calibrator (lyophilized)	1 vial
Cystatin C Monkey Detection Antibody	1 vial
Cystatin C Monkey HRP Streptavidin Conjugate	1 vial
Stop Solution	1 x 12ml

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

Tissue specificity 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.

Involvement in disease Defects in CST3 are the cause of amyloidosis type 6 (AMYL6) [MIM:105150]; also known as hereditary cerebral hemorrhage with amyloidosis (HCHWA), cerebral amyloid angiopathy (CAA) or cerebroarterial amyloidosis Icelandic type. AMYL6 is a hereditary generalized amyloidosis due to cystatin C amyloid deposition. Cystatin C amyloid accumulates in the walls of arteries, arterioles, and sometimes capillaries and veins of the brain, and in various organs including lymphoid tissue, spleen, salivary glands, and seminal vesicles. Amyloid deposition in the cerebral vessels results in cerebral amyloid angiopathy, cerebral hemorrhage and premature stroke. Cystatin C levels in the cerebrospinal fluid are abnormally low. Genetic variations in CST3 are associated with age-related macular degeneration type 11 (ARMD11) [MIM:611953]. ARMD is a multifactorial eye disease and the most common cause of irreversible vision loss in the developed world. In most patients, the disease is manifest as ophthalmoscopically visible yellowish accumulations of protein and lipid that lie beneath the retinal pigment epithelium and within an elastin-containing structure known as Bruch membrane.

Sequence similarities Belongs to the cystatin family.

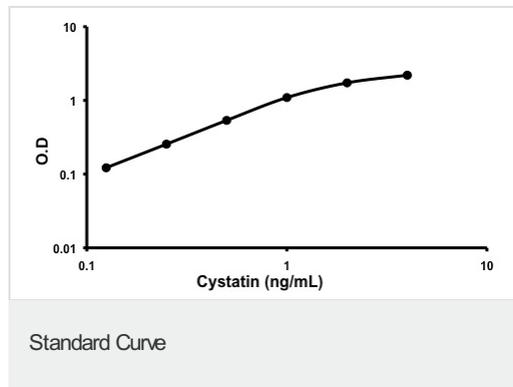
Post-translational modifications

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.

Cellular localization

Secreted.

Images



Representative standard curve using Cystatin Monkey ELISA kit (ab190542)

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