

Product datasheet

Human alpha 2 Macroglobulin ELISA Kit ab108883

★★★★★ 1 Abreviews 2 References 1 Image

Overview

Product name Human alpha 2 Macroglobulin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				3.3%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				8.9%

Sample type Cell culture supernatant, Saliva, Milk, Cerebral Spinal Fluid

Assay type Sandwich (quantitative)

Sensitivity = 0.95 ng/ml

Range 1 ng/ml - 25 ng/ml

Recovery 97 %

Assay time 4h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's alpha 2 Macroglobulin Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of alpha 2 Macroglobulin in saliva, milk, cerebrospinal fluid and cell culture supernatants.

An alpha 2 Macroglobulin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an alpha 2 Macroglobulin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase 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

amount of alpha 2 Macroglobulin captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Platform Microplate

Properties

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

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated alpha 2 Macroglobulin Antibody	1 x 120µl
alpha 2 Macroglobulin Microplate (12 x 8 well strips)	1 unit
alpha 2 Macroglobulin Standard	1 vial
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml

Function Is able to inhibit all four classes of proteinases by a unique 'trapping' mechanism. This protein has a peptide stretch, called the 'bait region' which contains specific cleavage sites for different proteinases. When a proteinase cleaves the bait region, a conformational change is induced in the protein which traps the proteinase. The entrapped enzyme remains active against low molecular weight substrates (activity against high molecular weight substrates is greatly reduced). Following cleavage in the bait region a thioester bond is hydrolyzed and mediates the covalent binding of the protein to the proteinase.

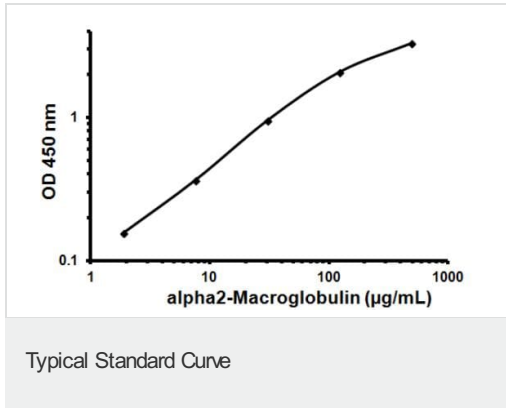
Tissue specificity Secreted in plasma.

Sequence similarities Belongs to the protease inhibitor I39 (alpha-2-macroglobulin) family.

Developmental stage Contrary to the rat protein, which is an acute phase protein, this protein is always present at high levels in circulation.

Cellular localization Secreted.

Images



Representative Standard Curve using ab108883.

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