

Human Fetuin A ELISA Kit (AHSG) ab108855

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

Overview

Product name	Human Fetuin A ELISA Kit (AHSG)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			3.6%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			8.7%
Sample type	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	1.4 ng/ml			
Range	3.125 ng/ml - 200 ng/ml			
Recovery	98 %			
Assay time	4h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Fetuin A (AHSG) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Fetuin A concentrations in Human plasma, serum, urine, milk and cell culture supernatants.			

A Fetuin A specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Fetuin A specific biotinylated detection antibody is added and then followed by washing with wash buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize 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 amount of Fetuin A captured in plate.

Get higher sensitivity in only 90 minutes with Human Fetuin A ELISA Kit (AHSG) ([ab185989](#)) from our SimpleStep ELISA[®] range.

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 N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Fetuin A Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Fetuin A Microplate (12 x 8 well strips)	1 unit
Fetuin A Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function Promotes endocytosis, possesses opsonic properties and influences the mineral phase of bone. Shows affinity for calcium and barium ions.

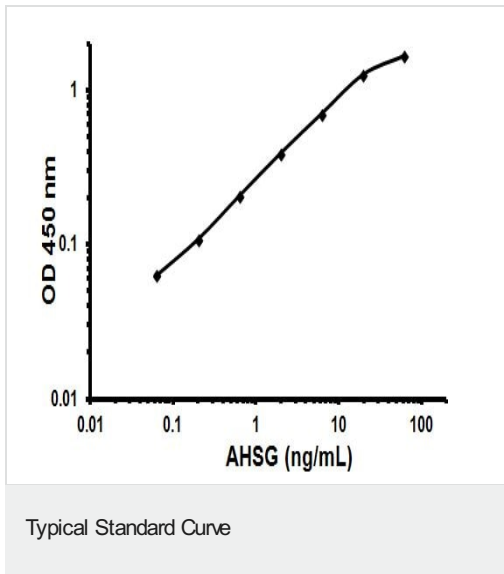
Tissue specificity Synthesized in liver and selectively concentrated in bone matrix. Secreted in plasma. It is also found in dentin in much higher quantities than other plasma proteins.

Sequence similarities Belongs to the fetuin family.
Contains 2 cystatin fetuin-A-type domains.

Post-translational modifications Phosphorylation sites are present in the extracellular medium.

Cellular localization Secreted.

Images



Representative Standard Curve using ab108855.

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