

Mouse Fetuin A ELISA Kit ab205074

1 Image

Overview

Product name	Mouse Fetuin A ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
Sample type	Urine, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.846 ng/ml				
Range	3.125 ng/ml - 100 ng/ml				
Recovery	> 85 %				
	Sample specific recovery				
	Sample type	Average %		Range	
	Serum	> 85		3.125ng/ml - 100ng/ml	
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Fetuin A in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Fetuin A protein in mouse plasma, serum and urine.				
Notes	In this assay the Fetuin present in samples reacts with the anti-Fetuin antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-Fetuin antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound Fetuin. Following another washing step, the enzyme bound to the immunosorbent is 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 Fetuin in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Fetuin in the test sample. The quantity of Fetuin in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X HRP Conjugated Enzyme Antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentration	1 x 50ml
Anti-Mouse Fetuin ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Mouse Fetuin Calibrator (lyophilized)	1 vial
Stop Solution	1 x 12ml

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

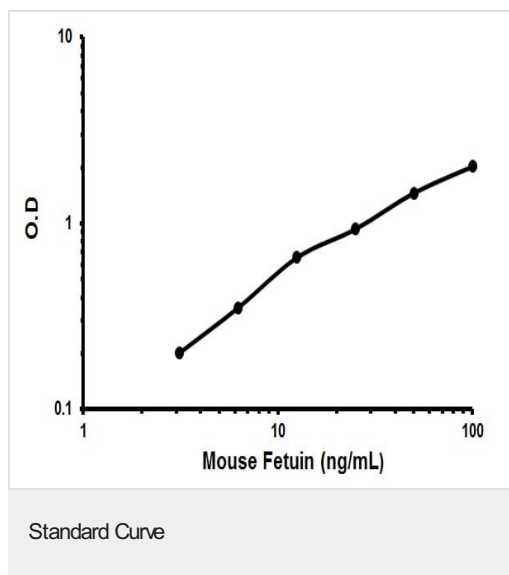
Tissue specificity Synthesized in liver and selectively concentrated in bone matrix. Secreted in plasma. It is also found in dentin in much higher quantities than other plasma proteins.

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 ab205074 Mouse Fetuin ELISA Kit.

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