

Mouse Periostin ELISA Kit ab193727

2 Images

Overview

Product name	Mouse Periostin ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma				
Assay type	Sandwich (qualitative)				
Sensitivity	0.05 ng/ml				
Range	0.05 ng/ml - 16 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		116.6		88% - 145%
	Plasma		109		84% - 134%
	Cell culture media		107		91% - 121%
Assay duration	One step assay				
Species reactivity	Reacts with: Mouse				
Product overview	<p>Abcam's Periostin Mouse ELISA Kit (ab193727) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse Periostin in serum, plasma (Plasma sample should be collected using EDTA or heparin as an anticoagulant. Citrate is not recommended.) and cell culture supernatants.</p> <p>This assay employs an antibody specific for mouse Periostin coated on a 96-well plate.</p>				

Standards and samples are pipetted into the wells and the immobilized antibody captures Periostin present in the samples. The wells are washed and biotinylated anti-mouse Periostin antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of Periostin bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated Mouse Periostin detection antibody	2 vials
mouse Periostin Standards (lyophilized)	2 vials
Pre-coated mouse Periostin Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function Binds to heparin. Induces cell attachment and spreading and plays a role in cell adhesion. May play a role in extracellular matrix mineralization.

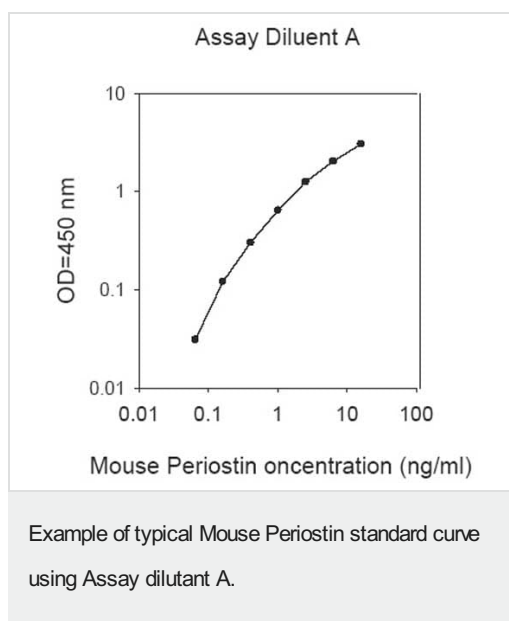
Tissue specificity Widely expressed with highest levels in aorta, stomach, lower gastrointestinal tract, placenta, uterus and breast. Up-regulated in epithelial ovarian tumors. Not expressed in normal ovaries. Also highly expressed at the tumor periphery of lung carcinoma tissue but not within the tumor. Overexpressed in breast cancers.

Sequence similarities Contains 1 EMI domain.
Contains 4 FAS1 domains.

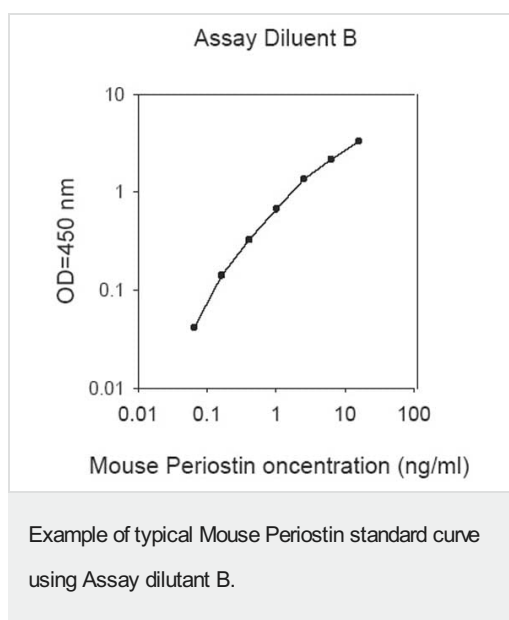
Post-translational modifications Gamma-carboxyglutamate residues are formed by vitamin K dependent carboxylation. These residues are essential for the binding of calcium.

Cellular localization Secreted > extracellular space > extracellular matrix.

Images



Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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