

Product datasheet

Mouse Osteopontin ELISA Kit ab100734

★★★★★ 2 Abreviews 3 References 1 Image

Overview

Product name	Mouse Osteopontin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 4 pg/ml
Range	2.74 pg/ml - 2000 pg/ml
Recovery	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	85.46	78% - 105%
Serum	115.6	103% - 122%
Plasma	110.4	87% - 130%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's Osteopontin Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse Osteopontin in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Osteopontin coated on a 96-well plate. Standards and samples are pipetted into the wells and Osteopontin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse Osteopontin antibody is added. After washing away unbound biotinylated antibody, HRP conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Osteopontin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse Osteopontin	2 vials
Osteopontin Microplate (12 x 8 wells)	1 unit
Recombinant Mouse Osteopontin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function

Binds tightly to hydroxyapatite. Appears to form an integral part of the mineralized matrix. Probably important to cell-matrix interaction. Acts as a cytokine involved in enhancing production of interferon-gamma and interleukin-12 and reducing production of interleukin-10 and is essential in the pathway that leads to type I immunity.

Tissue specificity

Bone. Found in plasma.

Sequence similarities

Belongs to the osteopontin family.

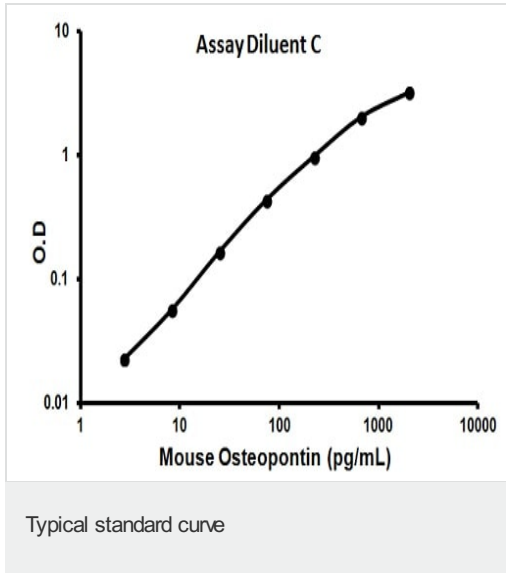
Post-translational modifications

Extensively phosphorylated on clustered serine residues. N- and O-glycosylated. Phosphorylation sites are present in the extracellular medium.

Cellular localization

Secreted.

Images



Representative standard curve using ab100734

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