

Mouse Osteopontin ELISA Kit ab100734

★★★★★ [2 Abreviews](#) [5 References](#) [1 图像](#)

概述

产品名称	小鼠Osteopontin ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 4 pg/ml
范围	2.74 pg/ml - 2000 pg/ml
回收率	100 %

特定样本回收率

样品类型	平均 %	范围
Cell culture supernatant	85.46	78% - 105%
Serum	115.6	103% - 122%
Plasma	110.4	87% - 130%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 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.

平台 Microplate

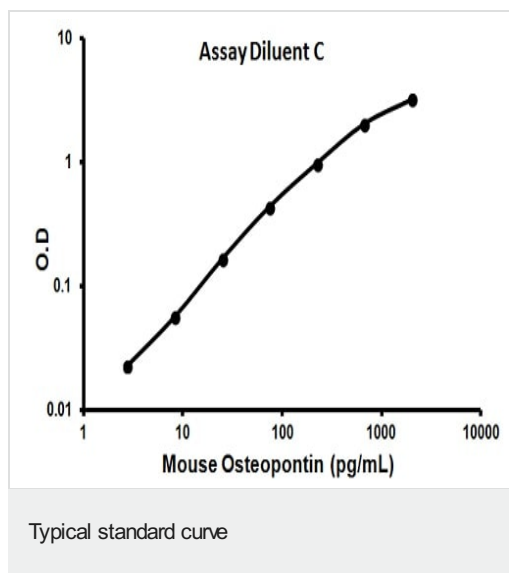
性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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

功能	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.
组织特异性	Bone. Found in plasma.
序列相似性	Belongs to the osteopontin family.
翻译后修饰	Extensively phosphorylated on clustered serine residues. N- and O-glycosylated. Phosphorylation sites are present in the extracellular medium.
细胞定位	Secreted.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors