

### Mouse Osteoactivin ELISA Kit ab267574

1 Image

#### Overview

**Product name** Mouse Osteoactivin 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** Serum, Plasma, Cell culture media

**Assay type** Sandwich (quantitative)

**Sensitivity** 80 pg/ml

**Range** 81.92 pg/ml - 20000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	96	88% - 104%
Plasma	91	71% - 116%
Cell culture media	108	97% - 122%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** Mouse Osteoactivin ELISA Kit is designed for the quantitative determination of Osteoactivin in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for mouse Osteoactivin coated on a 96-well plate. Standards and samples are pipetted into the wells and Osteoactivin present in a sample is bound

to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse Osteoactivin 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 Osteoactivin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit shows no cross-reactivity with any of the cytokines tested: Mouse CD30, L CD30, T CD40, CRG-2, CTACK, Eotaxin, Eotaxin-2, Fas Ligand, Fractalkine, GCSF, GM-CSF, IFN-gamma, IGFBP-3, IGFBP-5, IGFBP-6, IL-1 alpha, IL-1 beta, IL-2, IL-3, IL-3 Rb, IL-5, IL-6, IL-9, IL-10, IL-12 p40/p70, IL-12 p70, IL-13, IL-17, KC, Leptin R, LEPTIN(OB), LIX, L-Selectin, Lymphotactin, MCP-1, MCP-5, M-CSF, MIG, MIP-1, MIP-2, MIP-3, PF-4, P-Selectin, RANTES, SCF, SDF-1, TARC, TCA-3, TECK, TIMP-1, TNFalpha, TNF RI, TNF RII, TPO, VCAM-1, VEGF.

**Platform** Pre-coated 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	1 x 15ml
Anti-Mouse Osteoactivin coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Mouse Osteoactivin Detection Antibody	2 vials
Mouse Osteoactivin Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

**Function** Could be a melanogenic enzyme.

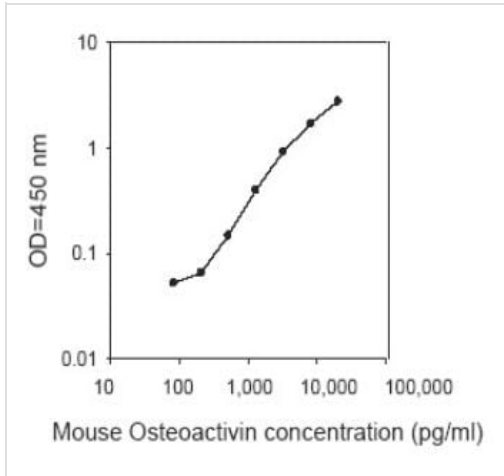
**Tissue specificity** Not restricted to the melanocytic lineage.

**Sequence similarities** Belongs to the PMEL/NMB family.  
Contains 1 PKD domain.

**Developmental stage** Expression in poorly metastatic melanoma cell lines; no expression in highly metastatic melanoma cell lines.

**Cellular localization** Membrane. Melanosome. Identified by mass spectrometry in melanosome fractions from stage I to stage IV.

**Images**



Mouse Osteoactivin ELISA kit (ab267574) Standard curve.

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

Standard curve

**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