

Human Osteoactivin ELISA Kit ab193711

[2 References](#) [2 图像](#)

概述

产品名称 人Osteoactivin ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

| 样品 | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 10% |

批次间

| 样品 | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 12% |

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 45 pg/ml

范围 45 pg/ml - 12000 pg/ml

回收率 特定样本回收率

| 样品类型 | 平均% | 范围 |
|--------------------|-------|-------------|
| Serum | 117.8 | 111% - 126% |
| Plasma | 116.5 | 107% - 126% |
| Cell culture media | 112.9 | 98% - 149% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Osteoactivin Human ELISA Kit (ab193711) is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Osteoactivin in sera, plasma and cell culture supernatants.

This assay employs an antibody specific for Human Osteoactivin coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures

Osteoactivin present in the samples. The wells are washed and biotinylated anti-Human Osteoactivin 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 Osteoactivin 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.

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

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 |
| 5X Assay Diluent B | 1 x 15ml |
| 500X HRP-Streptavidin Concentrate | 1 x 200µl |
| Assay Diluent A | 1 x 30ml |
| Biotinylated Human Osteoactivin detection antibody | 2 vials |
| Human Osteoactivin Standard (lyophilized) | 2 vials |
| Pre-coated Human Osteoactivin Microplate (12 strips x 8 wells) | 1 unit |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

功能 Could be a melanogenic enzyme.

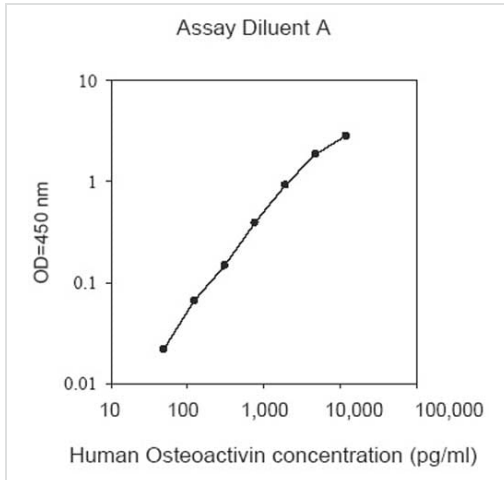
组织特异性 Not restricted to the melanocytic lineage.

序列相似性 Belongs to the PMEL/NMB family.
Contains 1 PKD domain.

发展阶段 Expression in poorly metastatic melanoma cell lines; no expression in highly metastatic melanoma cell lines.

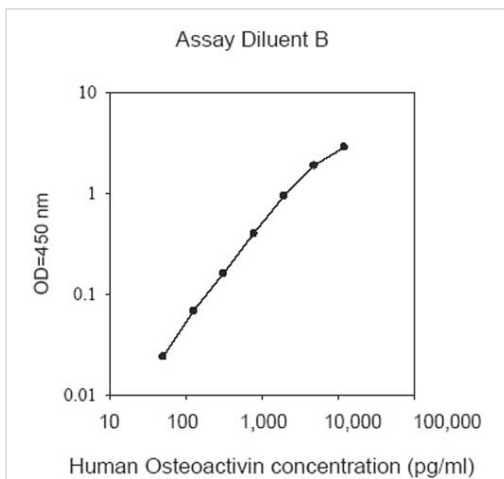
细胞定位 Membrane. Melanosome. Identified by mass spectrometry in melanosome fractions from stage I to stage IV.

图片



Typical standard curve in Assay Diluent A

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



Typical standard curve in Assay Diluent B

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.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