

Rat Osteopontin ELISA kit ab205076

1 References 1 图像

概述

产品名称 大鼠Osteopontin 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				< 10%

样品类型 Urine, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 0.02 ng/ml

范围 0.06 ng/ml - 4 ng/ml

回收率 > 85 %

			特定样本回收率
	样品类型	平均%	范围
	Serum	> 85	0.06ng/ml - 4ng/ml

检测时间 3h 20m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 Osteopontin in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Osteopontin protein in rat plasma, serum and urine.

说明 In this assay the Osteopontin present in samples reacts with the anti-Osteopontin antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, the Detection Antibody, biotin conjugated anti-Osteopontin, is added and complexes are formed. Following a wash step, horseradish peroxidase (HRP)

conjugated Streptavidin is added and complexes are formed. After another washing step, the complexes are assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of Osteopontin in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Osteopontin in the test sample. The quantity of Osteopontin in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

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

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
100X Detection Antibody Conjugate	1 x 150µl
100X HRP-Streptavidin	1 x 150µl
1X Diluent Concentrate	1 x 60ml
20X Wash Buffer Concentrate	1 x 50ml
Anti-Rat Osteopontin ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Rat Osteopontin Calibrator (Lyophilized)	1 vial
Stop Solution	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.

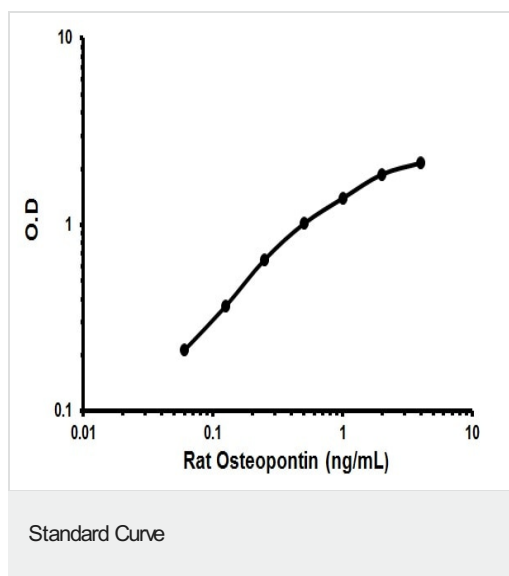
组织特异性 Bone. Found in plasma.

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 ab205076 Rat Osteopontin ELISA Kit.

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