

Product datasheet

Rat Osteopontin ELISA kit ab205076

1 Image

Overview

Product name Rat Osteopontin ELISA kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Sample type Urine, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.02 ng/ml

Range 0.06 ng/ml - 4 ng/ml

Recovery > 85 %

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	0.06ng/ml - 4ng/ml

Assay time 3h 20m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat

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

Notes 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.

Tested applications **Suitable for:** Sandwich ELISA
Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	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

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.

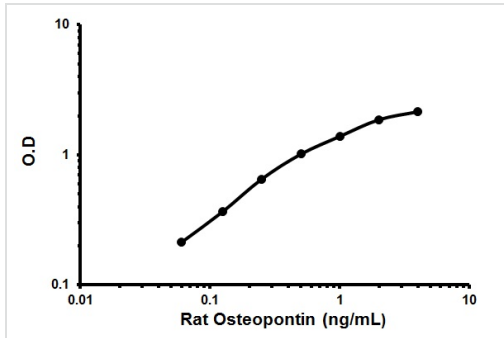
Applications

Our [Abpromise guarantee](#) covers the use of **ab205076** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative standard curve using ab205076 Rat Osteopontin ELISA Kit.

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