

# Osteopontin (human), ELISA kit

The Osteopontin (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The Osteopontin (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. Ready-to-use liquid color-coded reagents reduce error. Excellent publications in peer-reviewed publications demonstrate the utility and consistency of this kit developed with highly cited double monoclonal antibodies.

Osteopontin (OPN) is an acidic extracellular matrix cell adhesion protein that is relatively abundant not only in bone matrix, plasma, urine, and milk, but is also found in malignant and atherogenic tissues. Phosphorylation, glycosylation and calcium modifications allow intact and fragmented Osteopontin to direct a variety of diverse responses including tissue remodeling, inflammation and cell survival. Plasma Osteopontin has been shown to be a positive indicator of colon and lung cancers as well as metastatic carcinomas. The notable presence of Osteopontin in a variety of tumors is strongly correlated to pathological stage, suggesting its critical role in tumor invasiveness, progression and metastasis. In addition, Osteopontin inhibits inducible nitric oxide synthase activity, thereby protecting tumor cells from NO-mediated macrophage cytotoxic attack. Osteopontin is found in atherosclerotic plaques and may drive a number of diabetic vascular pathologies.

Citations: 14

[View Online »](#)

## Ordering Information

[Order Online »](#)

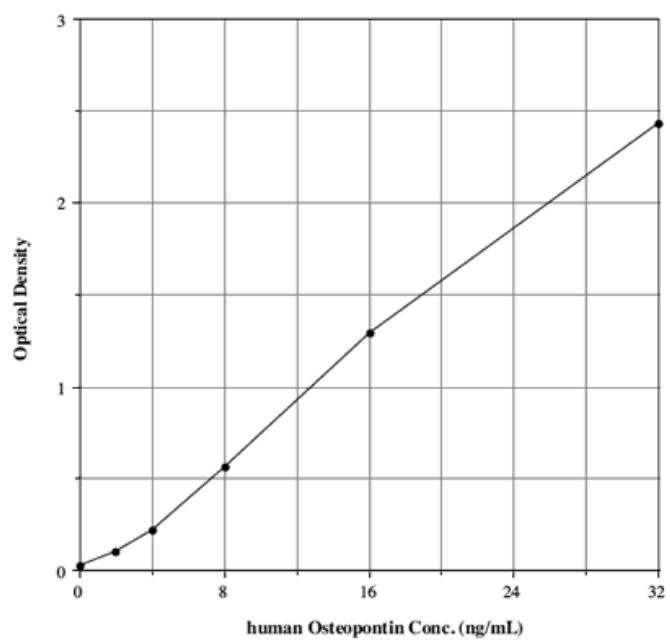
ADI-900-142

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Tested - thoroughly validated in plasma, urine, and milk
- Fast - results in just 3 hours from up to 41 samples in duplicate
- Trusted - data correlating assay results to cancer patient survival published in peer-reviewed journals
- Reliable - reproducible results day-after-day and lot-after-lot
- Convenient - ready-to-use pre-coated plates and liquid color-coded reagents reduce errors



## Handling & Storage

Use/Stability	Store all components at +4°C, except standard and conjugate at -20°C.
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	OPN, SPP1, Secreted phosphoprotein 1, Bone sialoprotein 1
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of human Osteopontin in culture supernatants, milk, plasma, and urine. Cited sample type includes serum.
Assay Time	3 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Microtiter plate, Conjugate, Antibody, Assay buffer 10 concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution
Crossreactivity	Human Osteopontin (100%), native human Osteopontin (111%) and <0.1%: native bovine Osteopontin, recombinant mouse Osteopontin
Sensitivity	0.110 ng/ml (range 2 – 32 ng/ml)
Species Reactivity	Human
UniProt ID	P10451
Wavelength	405 nm



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)