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 »

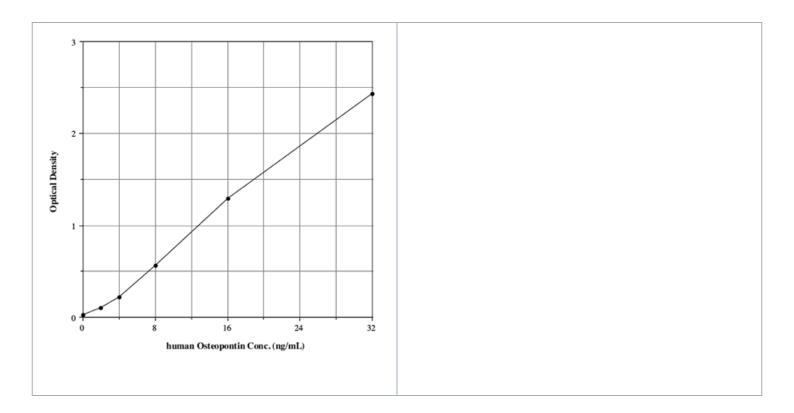
ADI-900-142

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 precoated plates and liquid colorcoded 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 NotesFor 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 Absorbance 96 Plate

Reader.

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com