

Osteopontin Antibody [#C0284]

Catalog Number: C0284-1, C0284-2, C0284

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No. : P10451

All Names: Bone sialoprotein 1, Nephropontin, OPN, OSTP, SPP-1, SPP1, Secreted phosphoprotein 1, Urinary stone protein, Uropontin

Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Storage/Stability: Store at -20°C/1 year

Immunogen: The antiserum was produced against synthesized peptide derived from human osteopontin

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity Osteopontin antibody detects endogenous levels of total osteopontin protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 ELISA: 1:20000

References: George R. Beck, J. Biol. Chem., Oct 2003; 278: 41921 - 41929.

KELLY L. HUDKINS, J. Am. Soc. Nephrol., Mar 1999; 10: 444.

Lígia R. Rodrigues, Cancer Epidemiol. Biomarkers Prev., Jun 2007; 16: 1087 - 1097.

Yu-Hua Hsieh, Cancer Res., Jul 2006; 66: 7119 - 7127.