

Osteoprotegerin (OPG) antibody (mAb)

Catalog No: 40938

RRID: AB_2793452

Clone: 98A1071

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 1.0 µg/µl

Molecular Weight: 55 kDa

Background: Osteoprotegerin (OPG) / osteoclastogenesis inhibitory factor (OCIF) is a cytokine that can inhibit the production of osteoclasts. It is a member of the tumor necrosis factor (TNF) receptor superfamily. It is a basic glycoprotein comprising 401 amino acid residues arranged into 7 structural domains. It is found as either a 60 kDa monomer or 120 kDa dimer linked by disulfide bonds. Osteoprotegerin inhibits the differentiation of osteoclast precursors into osteoclasts and also regulates the resorption of osteoclasts *in vitro* and *in vivo*. Osteoprotegerin is a RANK homolog, and works by binding to RANKL ligand on osteoblast/stromal cells, thus blocking the RANKL-RANK ligand interaction between osteoblast/stromal cells and osteoclast precursors. This has the effect of inhibiting the differentiation of the osteoclast precursor into a mature osteoclast.

Immunogen: This Osteoprotegerin (OPG) antibody was raised against a peptide corresponding to amino acid residues 20-37 of human Osteoprotegerin (OPG).

Buffer: PBS containing 0.02% gelatin and 0.05% sodium azide. Sodium azide is highly toxic.

Application Notes:

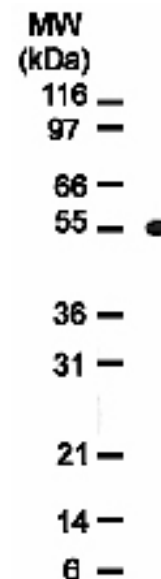
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Osteoprotegerin (OPG) mAb tested by Western blot.

Osteoprotegerin (OPG) mAb detection by Western blot. The analysis was performed using Daudi nuclear extract and Osteoprotegerin (OPG) mAb at a 2 µg/ml dilution. The immunoreactivity was detected by enhanced chemiluminescence.