Osteoprotegerin monoclonal antibody (98A1071)

Osteoprotegerin (OPG) is a secreted glycoprotein belonging to the TNF receptor (TNF-R) superfamily that functions as a decoy receptor for RANK-Ligand (RANKL/OPGL). Unlike other TNF-Rs, OPG lacks a membrane-anchoring hydrophobic domain and lacks any apparent cell-associated signaling capability. OPG secreted by osteoblasts competes with RANK (expressed on osteoblast precursors) for binding to RANKL, blocking osteoclastogenesis and increasing bone density.

This antibody is covered by our Worry-Free Guarantee.

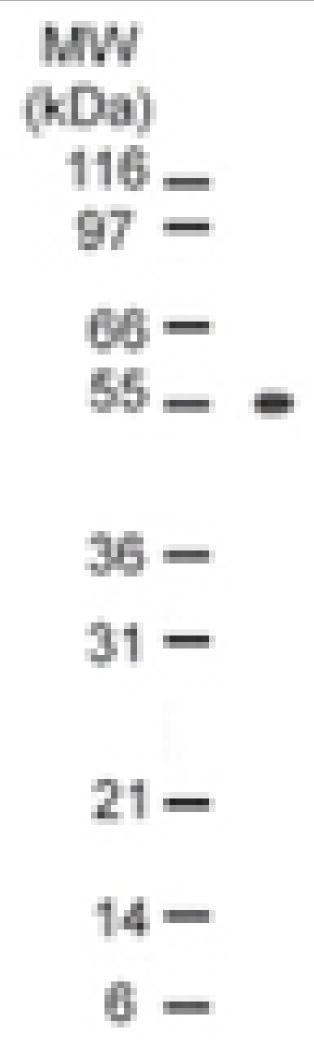
Ordering Information

Order Online »

ADI-AAM-020-E 100μg

Manuals, SDS & CofA

View Online »



Western blot analysis of Daudi cell lysate with OBG mAb at $1\mu g/ml$.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name OPG, TNFRSF 11B, Osteoclastogenesis inhibitory factor

Application WB

Application NotesDetects a band of ~55kDa & ~110kDa by Western blot.

Clone 98A1071

Formulation Liquid. In PBS containing 0.05% BSA and 0.05% sodium

azide.

Host Mouse

Immunogen Synthetic peptide corresponding to the sequence near the

N-terminus of human OPG (osteoprotegerin).

lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (1µg/ml AP, 1µg/ml ECL)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID 000300

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com