

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](#).

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Daudi cell lysate with OBG mAb at 1 µg/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 Notes Detects 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).

Isotype IgG1

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 O00300

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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)