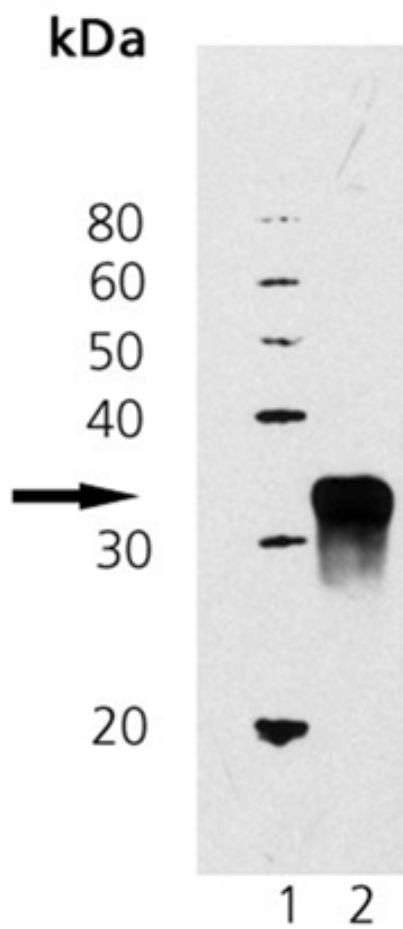


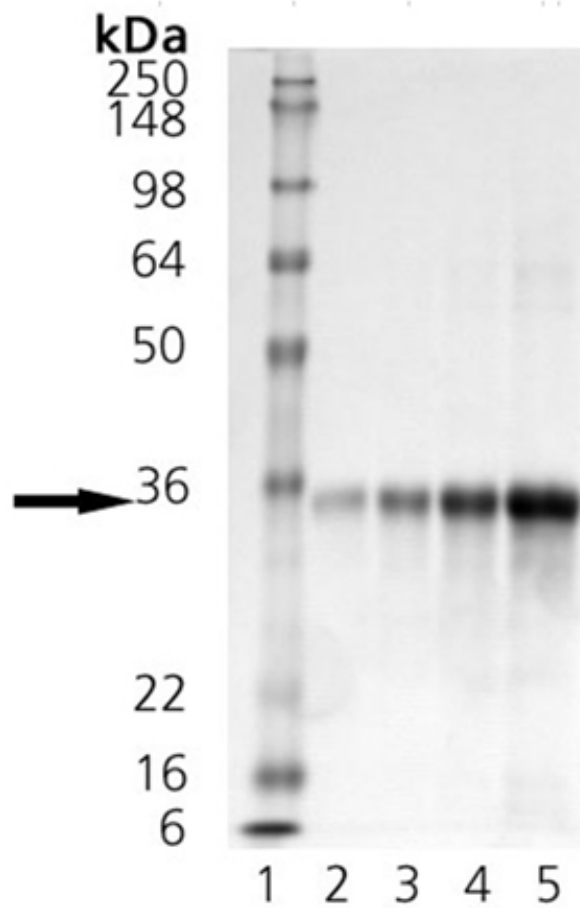
Dkk-1 (mouse), (recombinant) (His- tag)

Recombinant mouse Dkk-1 (His-tag) that inhibits Wnt/ β -catenin signaling and regulates bone remodeling.

Dkk-1 (Dickkopf or Dickkopf-1) is a secreted inhibitor of the Wnt/canonical signaling pathway which mediates a variety of cellular processes. Wnt proteins are extracellular signaling molecules that initiate activation of the canonical pathway by binding low-density lipoprotein receptor-related protein 5&6 (LRP5/6) coreceptors, triggering a cascade of events that stabilize β -catenin. Wnts mediate osteoblast development and cell-to-cell interactions during embryogenesis, and are also responsible for regulating bone mass in adults during normal bone remodeling processes. Wnts are also found to be expressed during the stationary phase of the cell cycle and are thought to control expansion of bone marrow stromal cells preventing overpopulation in marrow. There are at least two families of secreted inhibitors of Wnt signaling: the secreted frizzled-related protein family and the Dickkopf family. The mouse Dkk-1 protein consists of 272 amino acids including a 31 amino acid N-terminal signal peptide. Mature native Dkk-1 has a predicted molecular weight of 26kD and an apparent molecular weight of approximately 38-42kD by SDS-PAGE and western blot analysis. Dkk-1 disrupts the canonical pathway by binding to the LRP5/6 coreceptors, preventing interaction with Wnt and facilitating degradation of β -catenin, which ultimately results in osteolytic breakdown of bone tissue. Wnt signaling dysregulation has been implicated in cancer of the colon and stomach. Gonzalez-Sancho reported that the Wnt/ β -catenin pathway is downregulated by the induction of Dkk-1 expression, a mechanism that is lost in colon cancer, suggesting that Dkk-1 acts as a tumor suppressor gene in this cancer type. Dkk-1 proteins are expressed during log phase of the cell cycle, suggesting that dysregulated Dkk-1 could be responsible for the growth of some malignancies. The Dkk-1 gene has also been determined to be a target of p53, as Cappuccio suggests it is a necessary requirement for the execution of cell death. Tian reports that elevated Dkk-1 levels in bone marrow, plasma and peripheral blood from patients with multiple myeloma correlated with the gene-expression patterns of Dkk-1, and were associated with the presence of focal bone lesions. A study reported serum levels of Dkk-1 in patients with multiple myeloma before autologous stem cell transplantation. The range of Dkk-1 serum levels in this group of patients was 6.4-234.8 ng/mL with a mean of 63.6 ng/mL and a standard deviation of 77.3. Hall et al. suggest that Dkk-1 may act like a switch on prostate cancer cell activity, causing cells to change from osteoblastic to a highly osteolytic cell line.



Western Blot Analysis: Lane 1: MWM, Lane 2: 10 ng Dkk-1, probed with goat anti-mouse Dkk-1 pAb



SDS-PAGE Analysis: Lane 1: MWM, Lane 2: 0.5 mg, Lane 3: 1 mg, Lane 4: 2 mg, Lane 5: 5 mg Dkk-1

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Dickkopf-related protein-1
Application Notes	Western blot control.
Formulation	Liquid. In 50mM TRIS, pH 8.5, containing 250mM sodium chloride, 2mM DTT, and 1mM EDTA.
MW	~35kDa
Purity	≥95% (SDS-PAGE; Western blot)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Mouse Dkk-1 is fused at the N-terminus to a His-tag.
UniProt ID	O54908

Last modified: October 9, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com