

Dkk-1 (mouse), (recombinant) (active)

Modulator of embryogenesis and known effector of tissue development and morphogenesis

Dickkopf-related protein 1 (Dkk-1), which belongs to the dickkopf family, is a secreted protein with two cysteine rich domains and is involved in embryonic development and many human diseases. The regulation of Wnt signaling and Dkk1 inhibition has implications for bone formation and homeostasis. Dkk-1 can inhibit Wnt signaling by binding to its receptors LRP5/6 and kremen1/2 to form a complex, which can prevent Wnt from binding to LRP5/6. The predicted mouse Dkk1 protein is 272 amino acid residues with an N-terminal signal peptide. It shares 86% amino acid identity to human Dkk-1 protein.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-60002-C005	5µg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as supplied at -80°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Dickkopf-related protein-1
Biological Activity	Determined measuring the ability to inhibit the induction of the LEADING LIGHT [®] Wnt Reporter Cell Line (Prod. No. ENZ-61002) induced for at least 6 hours with 200 ng/ml Wnt3a (Prod. No. ENZ-60001). The EC50 of this inhibition is 45-55 ng/ml.
Endotoxin Content	<1.0EU/µg (LAL method)
Formulation	Lyophilized from a solution containing 1X PBS, 1% CHAPS, and 0.1% BSA.
MW	~40kDa (SDS-PAGE)
Purity	≥80% (SDS-PAGE)
Quantity	5µg
Reconstitution	Reconstitute by adding 1.5ml water to make a 3.3µg/ml solution.
Sequence	Mouse Dkk-1 amino acids 33-272, as well as a TEV recognition site and 6 histidines at the C terminus.
Source	Produced in insect cells. Produced in a baculovirus expression system.

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)