

Wnt-3a (mouse), (recombinant)

Modulator of embryogenesis, cell proliferation, stem cell differentiation, and bone, glucose and lipid metabolism

Wnt expression and signal transduction have been identified as critical components in regulating embryogenesis, cell proliferation, and bone, glucose and lipid metabolism. Wnt-3a signalling also plays a role in regulating the differentiation of stem cells. Wnt-3a is a member of the highly conserved, structurally related cysteine-rich Wnt family glycoproteins important for normal developmental processes and oncogenesis. Wnt-3a proteins are secreted and bind to receptors of the Frizzled and LRP families on the cell surface, which leads to the stabilization of β -catenin and increase of β -catenin translocation into the nucleus where it binds to TCF/LEF transcription factors to promote gene expression. The predicted mouse Wnt-3a protein of 352 amino acid residues and contains an N-terminal signal peptide for secretion. It shares 96% amino acid identity to human Wnt-3a protein. Due to glycosylation, it migrates at an apparent molecular weight of ~40 kDa by SDS-PAGE analysis under reducing conditions.

Ordering Information

[Order Online »](#)

ENZ-60001-C001	1 μ g
----------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored, as supplied, at -80°C. Following reconstitution, stable for three months when stored at or below -20°C.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Biological Activity	Wnt-3a can induce β -catenin mediated response in a TCF/LEF reporter gene luciferase assay.
Endotoxin Content	Less than 0.2 ng per μ g (2EU/ μ g)
Formulation	Lyophilized solid.
MW	~40kDa (SDS-PAGE)
Purity	>75% by SDS-PAGE and >75% by HPLC analyses
Reconstitution	Prepare 5 μ g/mL stock solution of Wnt3a by adding 200 μ L deionized water into the vial of lyophilized Wnt3a.
Source	Cell culture
UniProt ID	P27467

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)