

DLL4 (soluble) (human), (recombinant)

Notch ligand

Human sDLL4 comprises the extracellular signaling domain of DLL, a member of a structurally-related family of single-pass type I trans-membrane proteins that serve as ligands for Notch receptors. DLL4 functions to specifically activate the Notch-1 and Notch-4 receptors. The Notch signaling pathway regulates endothelial-cell differentiation, proliferation and apoptosis, and is essential for the development, maintenance and remodeling of the vascular system. Targeted deletion of the DLL4 gene in mice resulted in severe vascular defects and death before birth. Up-regulation of DLL4 expression has been implicated in the vascular development of certain tumors.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-766-0025	25µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	The lyophilized protein is stable for at least 2 years from date of receipt at -20°C. Reconstituted sDLL-4 is stable for at least 3 months when stored in working aliquots with a carrier protein at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Delta-like protein 4, Delta4
Biological Activity	The sDLL-4, when immobilized at concentrations >1.5 µg/mL will inhibit myogenesis in C2C12 cells.
Endotoxin Content	< 0.1 ng per µg of sDLL-4.
Formulation	Lyophilized.
GenBank ID	54567
MW	~54kDa
Purity	≥95% (SDS-PAGE and HPLC)
Source	Produced in HEK 293 cells. Recombinant human sDLL-4 contains 498 amino-acid residues.
UniProt ID	Q9NR61



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)