

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 »](#)

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 |

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)