

DLK1 (human), (recombinant) (His- tag)

Notch ligand

Protein Delta Homolog 1 (DLK-1) is a transmembrane protein which contains a signal peptide, an extracellular domain with six tandem epidermal growth factor (EGF)-like domains, a single pass transmembrane domain, and a short cytoplasmic tail. It is found within the stromal cells in close contact to the vascular structure of placental villi, yolk sac, fetal liver, adrenal cortex and pancreas and in the beta cells of the islets of Langerhans in the adult pancreas. In addition, it is detected in some forms of neuroendocrine lung tumor tissue. DLK-1 may have a important role in neuroendocrine differentiation.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-764-0010

10µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Lyophilized protein should be stored at < -20°C, though stable at room temperature for 3 weeks. Reconstituted protein solution can be stored at 4-7°C for 2-7 days. Aliquots of reconstituted samples are stable at < -20°C for 3 months.
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	Fetal antigen 1, Protein delta homolog 1, Preadipocyte factor 1, Pref-1, FA1, pG2
Formulation	Lyophilized. In 20mM PBS, pH 7.2, containing 150mM sodium chloride.
Purity	≥95% (SEC-HPLC and reducing SDS-PAGE)
Reconstitution	Dissolve the lyophilized protein in 1X PBS. It is not recommended to reconstitute to a concentration less than 100µg/ml.
Source	Produced by HEK 293 cells. DLK1 (Ala24-Pro297) is fused at the C-terminus to a His-tag.
UniProt ID	P80370

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)