DLK1 (human), (recombinant) (Histag)

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 improtant role in neuroendocrine differentiation.

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-201-764-0010	10µg
ALX-201-764-0050	50µg

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

Endotoxin Content <0.1ng/µg protein (1 IEU/µg).

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

