

DLK Protein

Recombinant human protein expressed in Sf9 cells

Catalog # M20-31G

Lot # S347-2

Product Description

Recombinant human DLK (1-520) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_006301](#).

Gene Aliases

MAP3K12; MUK; ZPK; ZPKP1; MEKK12; MAP3K12

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

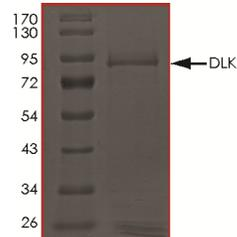
Scientific Background

DLK or mitogen-activated protein kinase kinase kinase 12 is a transmembrane protein containing six epidermal growth factor repeats and is a serine/threonine protein kinase. DLK is involved in the differentiation of several cell types, including adipocytes. DLK plays a role of a tumor suppressor protein and is predominately expressed in neuronal cells. DLK as an activator of the Jnk pathway. Overexpression of DLK results in activation of Jnk1 and the accumulation of a hyperphosphorylated form of c-Jun (1). Postnatal loss of DLK in stem cells and niche astrocytes has been shown to regulate neurogenesis (2).

References

- Hirai, S. et al: Activation of the JNK pathway by distantly related protein kinases, MEKK and MUK. *Oncogene* 12: 641-650, 1996.
- Ferron, S. R. et al: Postnatal loss of Dlk1 imprinting in stem cells and niche astrocytes regulates neurogenesis. *Nature* 475: 381-385, 2011.

Purity



The purity of DLK was determined to be **>80%** by densitometry. Approx. MW **94 kDa**.

DLK Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	M20-31G
Lot #	S347-2
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.