

LOXL3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # L263-31G

Lot # J583-2

Product Description

Recombinant human LOXL3 (526-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_032603](#).

Gene Aliases

LOXL

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.

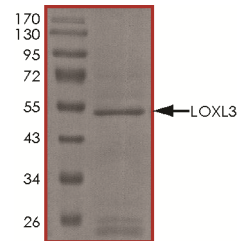
Scientific Background

LOXL3 is a member of the lysyl oxidase gene family. This family is essential to the biogenesis of connective tissue. The lysyl oxidase gene family encoding an extracellular copper-dependent amine oxidase that catalyses the first step in the formation of crosslinks in collagens and elastin (1). LOXL3 is highly expressed in central nervous system, uterus, heart, and leukocyte (1). LOXL3 plays an important role in developmental regulation, senescence, tumor suppression, cell growth control, and chemotaxis.

References

- Zhong, Z. et al: Stat3: a STAT family member activated by tyrosine phosphorylation in response to epidermal growth factor and interleukin-6. *Science*. 1994 Apr 1;264(5155):95-8.
- Tian, S. S. et al: Rapid activation of the STAT3 transcription factor by granulocyte colony-stimulating factor. *Blood*. 1994 Sep 15;84(6):1760- P53.

Purity



The purity of LOXL3 was determined to be **>90%** by densitometry. Approx. MW **54kDa**.

LOXL3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	L263-31G
Specific Lot Number	J583-2
Purity	>90%
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.