

LOXL1 Protein

Recombinant protein expressed in Sf9 cells

Catalog # L261-31G

Lot # O892-2

Product Description

Recombinant human LOXL1 (292-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC015090](#).

Gene Aliases

LOL, LOXL

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM 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.

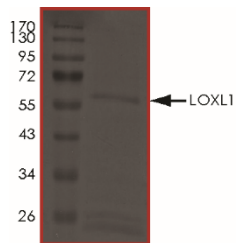
Scientific Background

LOXL1 is a member of the lysyl oxidase gene family which is essential to the biogenesis of connective tissue. LOXL1 encodes an extracellular copper-dependent amine oxidase that catalyses the first step in the formation of crosslinks in collagens and elastin. LOXL1 is responsible for catalyzing the oxidative deamination of lysine residues of tropoelastin and this deamination causes spontaneous cross-linking and formation of elastin polymer fibers (1). LOXL1 serves both as a crosslinking enzyme and an element of the scaffold to ensure spatially defined deposition of elastin (2).

References

- Hewitt, A. W.et.al: Ancestral LOXL1 variants are associated with pseudoexfoliation in Caucasian Australians but with markedly lower penetrance than in Nordic people. Hum. Molec. Genet. 17: 710-716, 2008.
- Liu, X. et.al: Elastic fiber homeostasis requires lysyl oxidase-like 1 protein. Nature Genet. 36: 178-182, 2004.

Purity



The purity of LOXL1 was determined to be **>70%** by densitometry. Approx. MW **61 kDa**.

LOXL1 Protein

Recombinant protein expressed in Sf9 cells

Catalog #	L261-31G
Lot #	O892-2
Purity	>70%
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.