

LOXL2 Protein

Recombinant protein expressed in Sf9 cells

Catalog # L262-31G

Lot # T1100-2

Product Description

Recombinant human LOXL2 (519-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC000594](#).

Gene Aliases

LOR2, WS9-14

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.

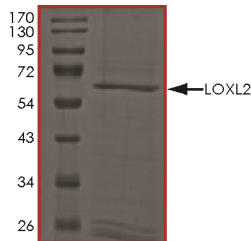
Scientific Background

LOXL2 is a member of the lysyl oxidase gene family which is essential for the biogenesis of connective tissue. LOXL2 encodes an extracellular copper-dependent amine oxidase that catalyses the first step in the formation of crosslinks in collagens and elastin. LOXL2 may play an important role in developmental regulation, senescence, tumor suppression, cell growth control, and chemotaxis to each member of the family (1). LOXL2 may mediate fibroblast activation through enzymatic crosslinking of fibrillar collagen and local matrix tension, resulting in activation of TGFβ1 signaling (2).

References

1. Jourdan-Le Saux, C. et al: The LOXL2 gene encodes a new lysyl oxidase-like protein and is expressed at high levels in reproductive tissues. *J. Biol. Chem.* 274: 12939-12944, 1999.
2. Barry-Hamilton, V. et al: Allosteric inhibition of lysyl oxidase-like-2 impedes the development of pathologic microenvironment. *Nature Med.* 16: 1009-1017, 2010.

Purity



The purity of LOXL2 was determined to be **>80%** by densitometry. Approx. MW **65 kDa**.

LOXL2 Protein

Recombinant protein expressed in Sf9 cells

Catalog #	L262-31G
Lot #	T1100-2
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.