

## LOXL4 Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # L264-31G**

Lot # Q2319-5

### Product Description

Recombinant human LOXL4 (492-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC013153](#).

### Gene Aliases

LOXC, FLJ21889

### 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.

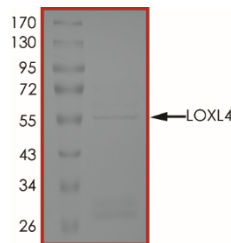
### Scientific Background

LOXL4 or lysyl oxidase-like 4 is a member of the lysyl oxidase family of amine oxidases. LOXL4 catalyzes the final enzymatic step required for the lysine-derived crosslinks essential for the formation of collagen fibrils and insoluble elastic fibers within the extracellular matrix (1). LOXL4 can also contribute to developmental regulation, senescence, tumor suppression, cell growth control, and chemotaxis. LOXL4 is highly expressed in skeletal muscle, testis, and pancreas (2).

### References

1. Asuncion, L.et.al: A novel human lysyl oxidase-like gene (LOXL4) on chromosome 10q24 has an altered scavenger receptor cysteine rich domain. *Matrix Biol.* 20: 487-491, 2001.
2. Maki, J. M. et.al: Cloning and characterization of a fifth human lysyl oxidase isoenzyme: the third member of the lysyl oxidase-related subfamily with four scavenger receptor cysteine-rich domains. *Matrix Biol.* 20: 493-496, 2001.

### Purity



The purity of LOXL4 was determined to be **>70%** by densitometry. Approx. MW **56 kDa**.

## LOXL4 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	L264-31G
Lot #	Q2319-5
Purity	>70%
Concentration	0.05µg/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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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