

LOX Protein

Recombinant human protein expressed in Sf9 cells

Catalog # L260-31G

Lot # U1751-8

Product Description

Recombinant human LOX (169-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC074872](#).

Gene Aliases

Lysyl oxidase

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.

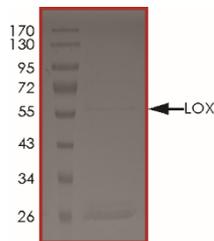
Scientific Background

LOX or lysyl oxidase is an extracellular copper enzyme that initiates the cross-linking of collagens and elastin. LOX catalyzes oxidative deamination of the epsilon-amino group in certain lysine and hydroxylysine residues of collagens and lysine residues of elastin (1). LOX plays an important role in tumor suppression (2). LOX may be required to create a niche permissive for metastatic growth and it is essential for hypoxia-induced metastasis.

References

1. Hamalainen, E.-R. et.al: Molecular cloning of human lysyl oxidase and assignment of the gene to chromosome 5q23.3-31.2. *Genomics* 11: 508-516, 1991.
2. Kaneda, A. et.al: Lysyl oxidase is a tumor suppressor gene inactivated by methylation and loss of heterozygosity in human gastric cancers. *Cancer Res.* 64: 6410-6415, 2004

Purity



The purity of LOX was determined to be **>70%** by densitometry. Approx. MW **57 kDa**.

LOX Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	L260-31G
Lot #	U1751-8
Purity	>70%
Concentration	0.05 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.