

# LOXL2 (human), (recombinant) (His- tag)

LOXL2, also known as Lysyl oxidase homolog 2, is a secreted protein that belongs to the lysyl oxidase family. LOXL2 contains four SRCR domains. The lysyl oxidase family is made up of five members that share conserved C-terminal catalytic domains that provide for lysyl oxidase or lysyl oxidase-like enzyme activity; and more divergent propeptide regions. LOXL2 is expressed by pre-hypertrophic and hypertrophic chondrocytes in vivo, and that LOXL2 expression is regulated in vitro as a function of chondrocyte differentiation. LOXL2 expression could be explored as a molecular target in the prevention of breast cancer progression.

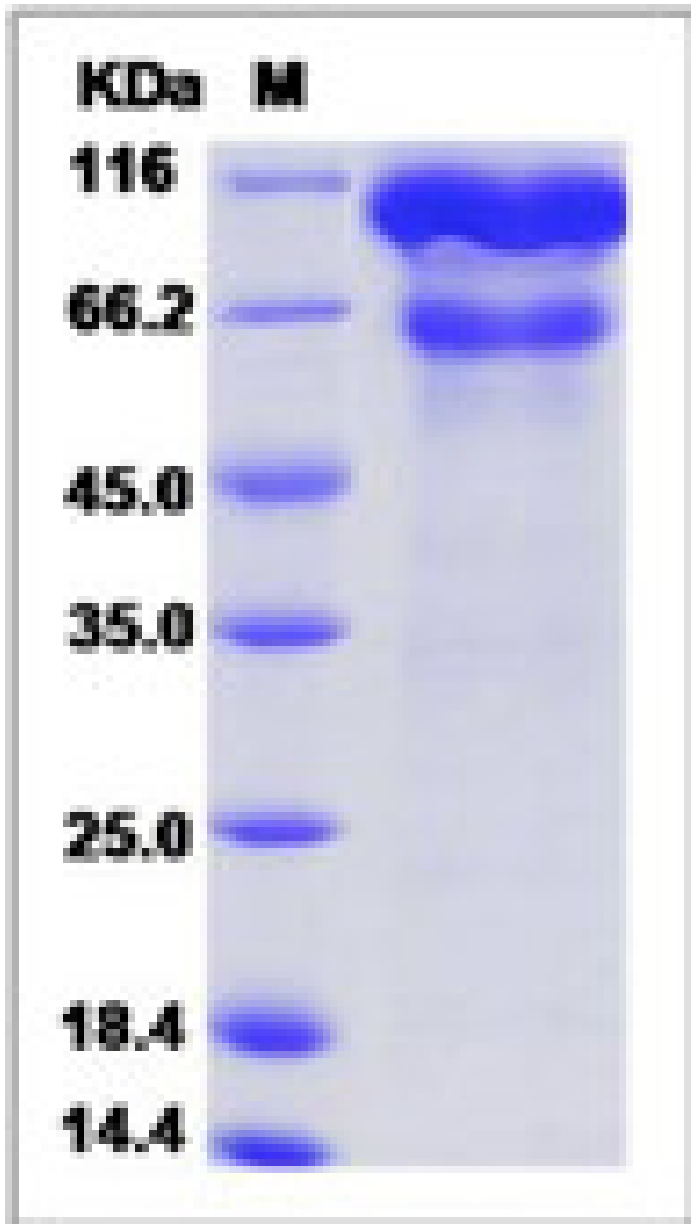
## Ordering Information

[Order Online »](#)

ENZ-PRT324-0050	50µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



As a result of glycosylation, the recombinant protein migrates as an approximately 168.3 kDa protein in SDS-PAGE under reducing conditions.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Lysyl oxidase homolog 2, Lysyl oxidase-like protein 2

**Endotoxin Content** <1 EU/µg (LAL method)

**Formulation** Lyophilized from sterile 20 mM MES, 50 mM NaCl

**MW** ~85.5 kDa

**Purity** ≥95% (SDS-PAGE)

**Reconstitution** Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

**Source** Produced in CHO Stable cells. A DNA sequence encoding the human LOXL2 (Met1-Gln774) was expressed, fused with a polyhistidine tag at the C-terminus.

**UniProt ID** Q9Y4K0

**Last modified: November 7, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)