

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.

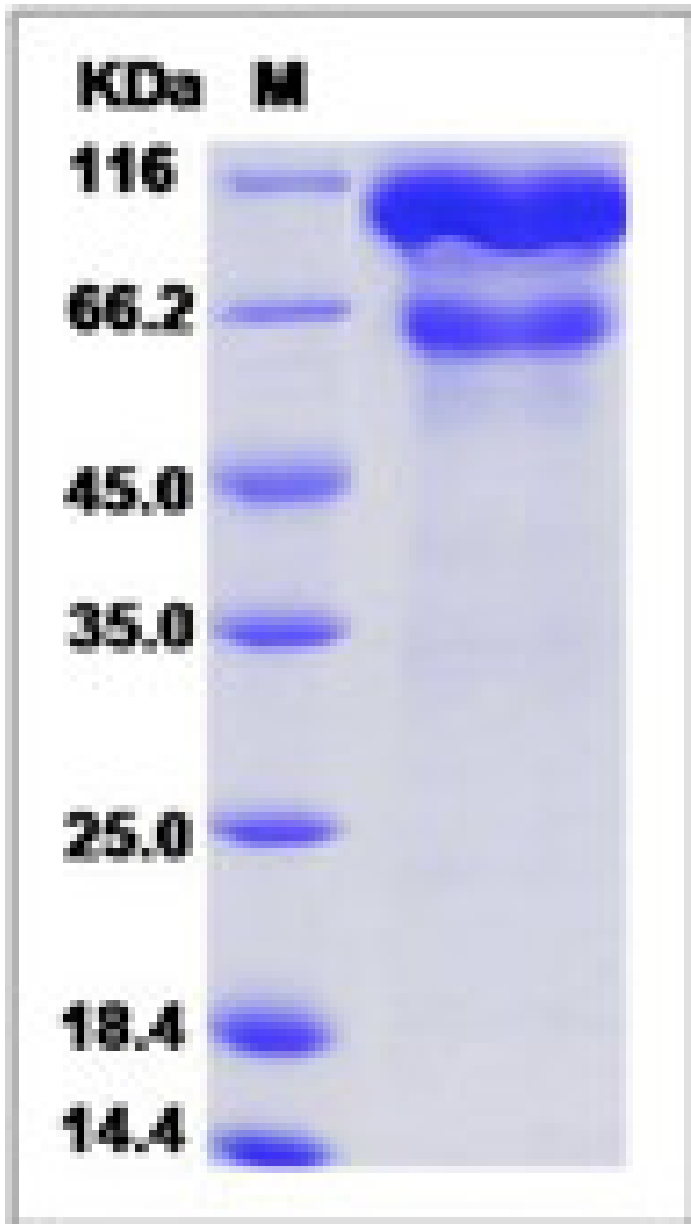
Ordering Information

[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



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)