

## Product Information

<b>Description:</b>	Recombinant human LOXL4 (Lysyl oxidase-like 4), encompassing amino acids 570-756(end). This construct contains a C-terminal His-tag. The recombinant protein was affinity purified.
<b>Background:</b>	LOXL4, also known as LOXC, belongs to the lysyl oxidase (LOX) gene family, and functions to stabilize collagen and elastin fibers via oxidative deamination. This is a gene of interest in connective tissue studies and is associated with diseases such as squamous cell carcinoma and phacoantigenic glaucoma. Further studies will reveal the mechanism of LOXL4 and other LOX genes in human cancer formation.
<b>Species:</b>	Human
<b>Construct:</b>	LOXL4 (570-756e-His)
<b>Concentration:</b>	0.33 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	9 mM Na <sub>2</sub> HPO <sub>4</sub> /KH <sub>2</sub> PO <sub>4</sub> pH 7.4, 124 mM NaCl, 2.4 mM KCl and 10% glycerol
<b>MW:</b>	23 kDa
<b>Genbank Accession:</b>	NM_032211
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Applications:</b>	Useful for SDS-PAGE.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

