

Catalog #	Aliquot Size
<b>L261-911-05</b>	<b>3 x 5 nmol</b>
<b>L261-911-20</b>	<b>3 x 20 nmol</b>
<b>L261-911-50</b>	<b>3 x 50 nmol</b>

## LOXL1 siRNA Set I

siRNA duplexes targeted against three exon regions

**Catalog # L261-911**

Lot # Z2049-71

### Specificity

LOXL1 siRNAs are designed to specifically knock-down human LOXL1 expression.

### Product Description

LOXL1 siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human LOXL1 mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [BC015090](#).

### Gene Aliases

LOL, LOXL

### Storage and Stability

The lyophilized powder is stable for at least 4 weeks at room temperature. It is recommended that the lyophilized and resuspended siRNAs are stored at or below -20°C. After resuspension, siRNA stock solutions  $\geq 2$   $\mu$ M can undergo up to 50 freeze-thaw cycles without significant degradation. For long-term storage, it is recommended that the siRNA is stored at -70°C. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

### Scientific Background

LOXL1 is a member of the lysyl oxidase gene family which is essential to the biogenesis of connective tissue. LOXL1 encodes an extracellular copper-dependent amine oxidase that catalyses the first step in the formation of crosslinks in collagens and elastin. LOXL1 is responsible for catalyzing the oxidative deamination of lysine residues of tropoelastin and this deamination causes spontaneous cross-linking and formation of elastin polymer fibers (1). LOXL1 serves both as a crosslinking enzyme and an element of the scaffold to ensure spatially defined deposition of elastin (2).

### References

- Hewitt, A. W.et.al: Ancestral LOXL1 variants are associated with pseudoexfoliation in Caucasian Australians but with markedly lower penetrance than in Nordic people. *Hum. Molec. Genet.* 17: 710-716, 2008.
- Liu, X. et.al: Elastic fiber homeostasis requires lysyl oxidase-like 1 protein. *Nature Genet.* 36: 178-182, 2004.

### Formulation

The siRNAs are supplied as a lyophilized powder and shipped at room temperature.

### Reconstitution Protocol

Briefly centrifuge the tubes (maximum RCF 4,000g) to collect lyophilized siRNA at the bottom of the tube. Resuspend the siRNA in 50  $\mu$ l of DEPC-treated water (supplied by researcher), which results in a 1x stock solution (10  $\mu$ M). Gently pipet the solution 3-5 times to mix and avoid the introduction of bubbles. Optional: aliquot 1x stock solutions for storage.

### Related Products

Product Name	Catalog Number
LOXL1 Protein	L261-31G
LOXL2 Protein	L262-31G
LOXL3 Protein	L263-31G

## LOXL1 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	L261-911
Specific Lot Number	Z2049-71
Packaging Specifications	2.5 nmol/tube for 3 x 5 nmol
Format	Lyophilized powder
Stability	1yr at -70°C from date of shipment
Storage & Shipping	The lyophilized powder is stable for at least 4 weeks at room temperature. It is recommended that the lyophilized and resuspended siRNAs are stored at or below -20°C. After resuspension, siRNA stock solutions $\geq 2$ $\mu$ M can undergo up to 50 freeze-thaw cycles without significant degradation. For long-term storage, it is recommended that the siRNA is stored at -70°C. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

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.**