

Catalog #	Aliquot Size
L263-911-05	3 x 5 nmol
L263-911-20	3 x 20 nmol
L263-911-50	3 x 50 nmol

## LOXL3 siRNA Set I

siRNA duplexes targeted against three exon regions

### Catalog # L263-911

Lot # Z2049-73

### Specificity

LOXL3 siRNAs are designed to specifically knock-down human LOXL3 expression.

### Product Description

LOXL3 siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human LOXL3 mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM\\_032603](#).

### Gene Aliases

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

LOXL3 is a member of the lysyl oxidase gene family. This family is essential to the biogenesis of connective tissue. The lysyl oxidase gene family encoding an extracellular copper-dependent amine oxidase that catalyses the first step in the formation of crosslinks in collagens and elastin (1). LOXL3 is highly expressed in central nervous system, uterus, heart, and leukocyte (1). LOXL3 plays an important role in developmental regulation, senescence, tumor suppression, cell growth control, and chemotaxis.

### References

- Zhong, Z. et al: Stat3: a STAT family member activated by tyrosine phosphorylation in response to epidermal growth factor and interleukin-6. *Science*. 1994 Apr 1;264(5155):95-8.
- Tian, S. S. et al: Rapid activation of the STAT3 transcription factor by granulocyte colony-stimulating factor. *Blood*. 1994 Sep 15;84(6):1760- P53.

### 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

## LOXL3 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	L263-911
Specific Lot Number	Z2049-73
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.**