

Catalog #	Aliquot Size
<b>A42-911-05</b>	<b>3 x 5 nmol</b>
<b>A42-911-20</b>	<b>3 x 20 nmol</b>
<b>A42-911-50</b>	<b>3 x 50 nmol</b>

## AFX siRNA Set I

siRNA duplexes targeted against three exon regions

### Catalog # A42-911

Lot # Z2013-14

### Specificity

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

### Product Description

AFX siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human AFX mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM\\_001170931](#).

### Gene Aliases

FOXO4; AFX1; MLLT7

### 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

AFX is a member of the O class of winged helix/forkhead transcription factor family. AFX proteins are involved in cell cycle regulation and inactivation of these proteins help in oncogenic transformation. AFX regulates apoptosis in part by suppressing the level of antiapoptotic BCLX through the transcriptional repressor BCL6 (1). AFX is down-regulated in patients with inflammatory bowel disease. AFX is expressed in human luteinized mural granulosa cells and may be involved in human folliculogenesis and luteinization. AFX acts as a novel anticancer agent for HER2-overexpressing cells.

### References

1. Tang, T. et al: The forkhead transcription factor AFX activates apoptosis by induction of the BCL-6 transcriptional repressor. *J. Biol. Chem.* 277:14255-14265, 2002.

### 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
Anti-AFX	A42-63R

## AFX siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	A42-911
Specific Lot Number	Z2013-14
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.**