

Catalog #	Aliquot Size
<b>A06-911B-05</b>	<b>3 x 5 nmol</b>
<b>A06-911B-20</b>	<b>3 x 20 nmol</b>
<b>A06-911B-50</b>	<b>3 x 50 nmol</b>

## Amyloid beta 42 siRNA Set I

siRNA duplexes targeted against three exon regions

### Catalog # A06-911B

Lot # Z2013-44

### Specificity

Amyloid beta 42 siRNAs are designed to specifically knock-down human Amyloid beta 42 expression.

### Product Description

Amyloid beta 42 siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human Amyloid beta 42 mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM\\_000484](#).

### Gene Aliases

### 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

Amyloid peptides, derived from amyloid precursor protein (APP), are thought to play a role in the development of the senile plaques associated with Alzheimer's Disease. The amyloid hypothesis presupposes that flaws in the processing of APP result in abnormally high levels of the longer, "stickier" forms of beta amyloid, known as Abeta42 and Abeta43, leading to aggregation of amyloid in the neuronal cell death and ultimately neuronal death. Mutations in the structure of Abeta40 and related peptides as well as in some of the enzymes involved in the processing of APP have been shown to alter the processing of APP. The sporadic (i.e., non-genetic) form of the disease, however, is far more common, caused by aging in concert with a number of both genetic and environmental risk factors.

### References

1. Esler WP, et al: Stereochemical specificity of Alzheimer's disease beta-peptide assembly. *Biopolymers*. 1999 May;49(6):505-14.
2. Roher AE, et al: beta-Amyloid-(1-42) is a major component of cerebrovascular amyloid deposits: implications for the pathology of Alzheimer disease. *Proc Natl Acad Sci U S A*. 1993 Nov 15;90 (22):10836-40.

### 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
Amyloid beta 42 Peptide	A06-58
Amyloid beta 42 (A21G), Flemish Mutant	A06-58C
Amyloid beta 42 (D23N), Iowa Mutant	A06-58F
Amyloid beta 42 (E22G), Arctic Mutant	A06-58B
Amyloid beta 42 (E22K), Italian Mutant	A06-58D
Amyloid beta 42 (E22Q), Dutch Mutant	A06-58E

## Amyloid beta 42 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	A06-911B
Specific Lot Number	Z2013-44
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.**