

Catalog #	Aliquot Size
A61-911-05	3 x 5 nmol
A61-911-20	3 x 20 nmol
A61-911-50	3 x 50 nmol

Annexin 5 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog # A61-911

Lot # Z2013-45

Specificity

Annexin 5 siRNAs are designed to specifically knock-down human Annexin 5 expression.

Product Description

Annexin 5 siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human Annexin 5 mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM_001154](#).

Gene Aliases

ANXA5, PP4, ANX5, ENX2

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 ≥ 2 μ 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

Annexin 5 is a member of the annexin family of calcium-dependent phospholipid binding proteins. Annexin 5 is a phospholipase A2 and protein kinase C inhibitory protein with calcium channel activity and a potential role in cellular signal transduction, inflammation, growth and differentiation. In vitro studies have shown that Annexin 5 is a cellular protein which plays a role in the inhibition of blood coagulation by competing for phosphatidylserine binding sites with prothrombin (1). Through affinity assays Annexin 5 has been localized on cells that expressed phosphatidylserine on the cell surface, a feature found in apoptosis as well as other forms of cell death (2).

References

1. Koopman, G. et al: Annexin V for flow cytometric detection of phosphatidylserine expression on B cells undergoing apoptosis. *Blood*, 1994; 84 (5): 1415–20.
2. Vermes, I. et al: A novel assay for apoptosis—flow cytometric detection of phosphatidylserine expression on early apoptotic cells using fluorescein labelled Annexin V. *J Immunol Methods*, 1995; 184 (1): 39.

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 μ l of DEPC-treated water (supplied by researcher), which results in a 1x stock solution (10 μ 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
Annexin 5 Protein	A61-30H

Annexin 5 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	A61-911
Specific Lot Number	Z2013-45
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 ≥ 2 μ 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
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.