

Catalog #	Aliquot Size
Y75-911-05	3 x 5 nmol
Y75-911-20	3 x 20 nmol
Y75-911-50	3 x 50 nmol

14-3-3 epsilon siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog # Y75-911

Lot # Z2013-2

Specificity

14-3-3 epsilon siRNAs are designed to specifically knock-down human 14-3-3 epsilon expression.

Product Description

14-3-3 epsilon siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human 14-3-3 epsilon mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM_006761](#).

Gene Aliases

14-3-3 epsilon, YWHAE, MDS, MDCR, KCIP-1, 14-3-3E, FLJ45465

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

14-3-3 \square also known as tyrosine 3-monooxygenase/tryptophan 5-monooxygenase activation protein is a member of the 14-3-3 family. 14-3-3 \square has been reported to be involved in heat shock transcription factor 1 (HSF1) and extracellular signal regulated protein kinase (ERK) pathways (1). 14-3-3 \square interaction with the deubiquitinating enzyme UBPY, also known as USP8, regulates cargo sorting and membrane traffic at early endosomes. Association of 14-3-3 \square , HSF1 with ERK during heat shock may modulate the amplitude of the response and control the termination of HSP expression on resumption of growth conditions (2).

References

1. Wang, X. et al: Interactions between extracellular signal-regulated protein kinase 1, 14-3-3epsilon, and heat shock factor 1 during stress. *J. Biol. Chem.* 2004; 279(47):49460-9.
2. Mizuno, E. et al: 14-3-3-dependent inhibition of the deubiquitinating activity of UBPY and its cancellation in the M phase. *Exp Cell Res.* 2007; 313(16):3624-34.

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
14-3-3 epsilon Protein	Y75-30G
14-3-3 epsilon Protein	Y75-30N

14-3-3 epsilon siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	Y75-911
Specific Lot Number	Z2013-2
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.