

Catalog #	Aliquot Size
T06-911B-05	3 x 5 nmol
T06-911B-20	3 x 20 nmol
T06-911B-50	3 x 50 nmol

TESK2 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog # T06-911B

Lot # Z2097-64

Specificity

TESK2 siRNAs are designed to specifically knock-down human TESK2 expression.

Product Description

TESK2 siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human TESK2 mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM_007170](#).

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 ≥ 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

TESK2 or testis-specific kinase 2 is a serine/threonine protein kinase that contains an N-terminal protein kinase domain and belongs to the TESK subgroup of the LIMK/TESK family of protein kinases. The kinase domain of TESK2 is structurally similar to the kinase domain of testis-specific protein kinase-1 and the LIM motif-containing protein kinases (LIMKs). TESK2 is predominantly expressed in testis and prostate and plays an important role in spermatogenesis (1). TESK2 phosphorylates cofilin specifically at Ser-3 and induces formation of actin stress fibers and focal adhesions (2).

References

- Rosok, O. et al: Identification and characterization of TESK2, a novel member of the LIMK/TESK family of protein kinases, predominantly expressed in testis. *Genomics* 61: 44-54, 1999.
- Toshima, J. et al: Cofilin phosphorylation and actin reorganization activities of testicular protein kinase 2 and its predominant expression in testicular Sertoli cells. *J Biol Chem.* 2001 Aug 17;276(33):31449-58.

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
TESK1 Protein	T05-34G
TESK2, Active	T06-10G

TESK2 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog #	T06-911B
Lot #	Z2097-64
Packaging Specifications	2.5 nmol/tube for 3 x 5nmol
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.