

Catalog #	Aliquot Size
<b>T833-911-05</b>	<b>3 x 5 nmol</b>
<b>T833-911-20</b>	<b>3 x 20 nmol</b>
<b>T833-911-50</b>	<b>3 x 50 nmol</b>

## TGFB3 siRNA Set I

siRNA duplexes targeted against three exon regions

**Catalog # T833-911**

Lot # Z2097-71

### Specificity

TGFB3 siRNAs are designed to specifically knock-down human TGFB3 expression.

### Product Description

TGFB3 siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human TGFB3 mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NP\\_003230.1](#).

### Gene Aliases

ARVD; FLJ16571; TGFB3; TGF-b3

### 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

TGFB3 or transforming growth factor, beta 3 is a member of the TGF $\beta$  family of proteins that is a ser/thr protein kinase. TGFB3 is involved in embryogenesis and cell differentiation. TGFB3 signaling activates transcription of the LEF1 gene to induce epithelial mesenchymal transformation during mouse palate development (1). TGFB3 function in normal palate and lung morphogenesis and implicates this cytokine in epithelial-mesenchymal interaction (2).

### References

1. Nawshad, A. et.al: TGF-beta-3 signaling activates transcription of the LEF1 gene to induce epithelial mesenchymal transformation during mouse palate development. *J. Cell Biol.* 163: 1291-1301, 2003.
2. Kaartinen, V. et.al: Abnormal lung development and cleft palate in mice lacking TGF-beta-3 indicates defects of epithelial-mesenchymal interaction. *Nature Genet.* 11: 415-421, 1995.

### 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
TGFB1	T831-40N
TGFB3	T833-40CN

## TGFB3 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog #	T833-911
Lot #	Z2097-71
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 $\geq 2$ $\mu$ 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.**