

Catalog #	Aliquot Size
B05-911-05	3 x 5 nmol
B05-911-20	3 x 20 nmol
B05-911-50	3 x 50 nmol

ALK6 (BMPR1B) siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog # B05-911

Lot # Z2013-29

Specificity

ALK6 (BMPR1B) siRNAs are designed to specifically knock-down human ALK6 (BMPR1B) expression.

Product Description

ALK6 (BMPR1B) siRNA is a pool of three individual synthetic siRNA duplexes designed to knock-down human ALK6 (BMPR1B) mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM_001203](#).

Gene Aliases

ALK-6; ALK6; CDw293

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

ALK6 (also known as BMPR1B) is a member of the transmembrane serine/threonine kinase that is the member of the bone morphogenetic protein (BMP) receptor, which is closely related to the activin receptors, ACVR1 and ACVR2. ALK6 is mainly involved in the endochondral bone formation and embryogenesis. ALK6 expressed in normal and cancerous prostate tissues and used in the endocrine therapy that given to the prostate cancer patients (1). ALK6 receptor trafficking also play a significant role in FOP pathogenesis and used in human T-cell differentiation (2).

References

1. Ide, H. et.al: Cloning of human bone morphogenetic protein type 1B receptor (BMPR-1B) and its expression in prostate cancer in comparison with other BMPRs. *Oncogene* 14: 1377-1382, 1997.
2. Cejalvo, T. et.al: Bone morphogenetic protein-2/4 signalling pathway components are expressed in the human thymus and inhibit early T-cell development. *Immunology* 121: 94-104, 2007.

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
ALK6 (BMPR1B), Active	B05-11G
ALK3 (BMPR1B), Active	B04-11G
BMPR2, Active	B06-11H
ALK, Active	A19-11G
ALK1, Active	A09-11G
ALK2, Active	A06-11G
ALK4, Active	A07-11G

ALK6 (BMPR1B) siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	B05-911
Specific Lot Number	Z2013-29
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.