

Catalog #	Aliquot Size
P50-911-05	3 x 5 nmol
P50-911-20	3 x 20 nmol
P50-911-50	3 x 50 nmol

AMPK B2 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog # P50-911

Lot # Z2194-4

Specificity

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

Product Description

AMPK is a pool of three individual synthetic siRNA duplexes designed to knock-down human AMPK mRNA expression. Each siRNA is 19-25 bases in length. The gene accession number is [NM_005399](#).

Gene Aliases

PRKAB2, MGC61468

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.

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
AMPK (A1/B1/G1), Active	P47-10H
AMPK (A1/B1/G2), Active	P55-10H
AMPK (A1/B1/G3), Active	P56-10H
AMPK (A1/B2/G1), Active	P50-10H
AMPK (A1/B2/G2), Active	P45-10G
AMPK (A1/B2/G3), Active	P83-10G

AMPK B2 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog Number	P50-911
Specific Lot Number	Z2194-4
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.