

| Catalog # | Aliquot Size |
|--------------------|--------------------|
| P51-911B-05 | 3 x 5 nmol |
| P51-911B-20 | 3 x 20 nmol |
| P51-911B-50 | 3 x 50 nmol |

AMPK G1 siRNA Set I

siRNA duplexes targeted against three exon regions

Catalog # P51-911B

Lot # Z2194-5

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_002733](#).

Gene Aliases

PRKAG1, AMPKG, MGC8666

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 G1 siRNA Set I

siRNA duplexes targeted against three exon regions

| | |
|--------------------------|---|
| Catalog Number | P51-911B |
| Specific Lot Number | Z2194-5 |
| 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.