

FAM Labeled Negative Control siRNA

FAM-labeled siRNA duplexes for monitoring siRNA delivery

Catalog # S03-910N

Lot # Z2207-4

Product Description

FAM Labeled Negative Control siRNA duplex is a 3'-dye-conjugated siRNA transfection efficiency control with no homology with mammalian genes. The synthetic siRNAs (5'-UUC UCC GAA CGU GUC ACG UTT-FAM-3' and 5'-ACG UGA CAC GUU CGG AGA ATT-3') enable qualitative assessment of cellular uptake, distribution and intracellular localization of siRNAs 1-4 hours post transfection. FAM Negative Control siRNA is commonly used to optimize transfection conditions. Transfection can be monitored by fluorescence microscopy or FACS.

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.

Spectral Information

Filter: FITC

Excitation max (λ_{max}): 494 nm

Emission max (λ_{max}): 520 nm

Formulation

The siRNAs are supplied as a lyophilized powder and shipped at room temperature. Package Specifications: 2.5 nmol/tube.

Reconstitution Protocol

Briefly centrifuge the tubes (maximum RCF 4,000g) to collect lyophilized siRNA at the bottom of the tube. Resuspend the siRNA in 25 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.

Purity

$\geq 90\%$ by HPLC

FAM Labeled Negative Control siRNA

FAM-labeled siRNA duplexes for monitoring siRNA delivery

Catalog #	S03-910N
Lot #	Z2207-4
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.