

siSTABLE™ control siRNA reagents

Product description

- Double-stranded, chemically synthesized oligonucleotides
- Duplexes modified with siSTABLE for extended stability and enhanced performance in nuclease-rich environments
- Mass of each strand confirmed by MALDI-TOF mass spectrometry

Shipping and storage

- Dharmacon™ siSTABLE™ Control siRNA reagents are shipped as dried pellets at room temperature (23 °C). Under these conditions, the reagents are stable for at least four weeks.
- Upon receipt, all siRNAs should be stored at –20 °C to –80 °C. Under these conditions, the siRNAs are stable for at least one year.
- siRNA should be resuspended in RNase-free solutions. We recommend 1x siRNA buffer (diluted from 5x siRNA buffer – Dharmacon Cat. #B-002000-UB-100). For *in vivo* use, RNase-free water (for short-term storage) or an RNase-free buffer (pH 7.3–7.6) such as PBS may be used for resuspension of concentrated stocks (20–100 μM).
- Upon resuspension, aliquot the siRNA into small volumes and store at –20 °C to –80 °C. For best results, limit freeze-thawing of each tube to no more than five events. Under these conditions, the siRNA is stable for at least one year.

Handling precautions

Oligonucleotides are susceptible to enzymatic degradation by nucleases and to chemical degradation by extreme pH and temperature. We recommend wearing gloves and maintaining nuclease-free conditions when handling the oligonucleotides.

Table 1. siSTABLE control siRNA reagents.

Product	Description	Cat. #
siSTABLE Non-targeting siRNA # 1	Negative control siRNA (chemically modified for increased stability) with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested. [^] MW 13,000 g/mol	D-001700-01-XX
siSTABLE Cyclophilin B siRNA (Human)	Positive silencing control for guaranteed silencing of Cyclophilin B mRNA in human cells (chemically modified for increased stability). Target NM_000942 MW 13,000 g/mol	D-001710-02-XX
siSTABLE Cyclophilin B siRNA (Mouse)	Positive silencing control for guaranteed silencing of Cyclophilin B mRNA in mouse cells (chemically modified for increased stability). Target NM_011149 MW 13,000 g/mol	D-001710-03-XX
siSTABLEG-APD siRNA (Mouse)	Positive silencing control for guaranteed silencing of GAPD mRNA in mouse cells (chemically modified for increased stability). Target NM_008084 MW 13,000 g/mol	D-001720-03-XX

[^]Non-targeting siRNA #1 may reduce EGFR mRNA by ~ 50% (NM_005228) in an assay- and cell-specific manner.

XX=05,20, or 50 for 5, 20, and 50 nmol amounts

Accompanying document

[Basic siRNA resuspension protocol.](#)

Supplemental documents

Go to dharmacon.horizondiscovery.com/protocols, to find:

- Basic and cell-line specific transfection protocols

Related products

DharmaFECT siRNA Transfection Reagents are available in four formulations that are optimized for transfecting siRNA into a wide variety of cell lines. For more information, click [here](#).

Publication reference guide

When referencing the use of siSTABLE Control siRNA Reagents, please include the following information: siSTABLE Control siRNA (type, Cat. # Dharmacon Inc, Lafayette, CO).

If you have any questions, contact

t +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

f + 44 (0)1223 655 581

w horizondiscovery.com/contact-us **or** dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2018 Horizon Discovery Group Company—All rights reserved. First published July 2014. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.