

# Lincode<sup>TM</sup> control reagents

## Product Description

- Double-stranded, chemically synthesized oligonucleotides
- Both sense and antisense strands contain UU for 3'-overhangs
- Antisense strands are modified with 5'-phosphate
- Mass of each strand confirmed by MALDI-TOF mass spectrometry
- Duplex integrity confirmed by non-denaturing polyacrylamide gel electrophoresis

## Shipping and storage

- RNAi Control reagents are shipped as dried pellets at room temperature (23 °C). Under these conditions, they are stable for at least four weeks.
- Upon receipt, RNAi Controls reagents should be stored at –20 °C to –80 °C. Under these conditions, the reagents 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). RNase-free water (for short-term storage) is also appropriate for resuspension of concentrated stocks (20-100 µM). Alternatively, an RNase-free buffer (pH 7.3-7.6) may be used such as PBS (Dharmacon Cat. #NC9826748).
- 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 6 months.

## 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.

## Accompanying documents

- [Basic siRNA Transfection Protocol](#).

## Related products

DharmaFECT<sup>TM</sup> siRNA Transfection Reagents are optimized for transfecting siRNA into a wide variety of cell lines.

For more information, go to [dharmacon.horizondiscovery.com/transfection/dharmafect-1-transfection-reagent](https://dharmacon.horizondiscovery.com/transfection/dharmafect-1-transfection-reagent).

## Publication Reference Guide

When referencing the use of DharmaFECT Duo Co-Transfection Reagents, please include the following information: DharmaFECT Duo Transfection Reagent, Dharmacon, Inc., Lafayette, CO.

**Table 1. Accession numbers for target genes of positive control siRNA.** GAS5 is a long noncoding RNA (lncRNA) expressed in a wide variety of human cell types, but its expression should be verified in each cell type prior to use as a positive control siRNA. Validated ON-TARGETplus™ positive control siRNAs targeting protein coding genes (such as Cyclophilin B) may also be used in experiments with Lincode siRNA reagents to verify transfection efficiency and successful RNAi.

Target Gene	Human
GAS5	NR_002578

**Table 2. Lincode RNAi controls.**

Product	Description	Size (nmol)	Cat. #
Lincode GAS5 Control siRNA <i>human</i>	Positive control for silencing of GAS5 lncRNA in human cells. MW ~ 13,400 g/mol	5	D-001310-01-05
		20	D-001310-01-20
		50	D-001310-01-50
Lincode GAS5 Control Pool <i>human</i>	Pool of 4 positive control siRNA for silencing of GAS5 lncRNA in human cells. Useful control for experiments with SMARTpool™ Reagents. MW ~ 13,400 g/mol	5	D-001310-10-05
		20	D-001310-10-20
		50	D-001310-10-50
Lincode Non-targeting siRNA #1 <i>human, mouse, and rat</i>	Negative control siRNA with 3 or more mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13,400 g/mol	5	D-001320-01-05
		20	D-001320-01-20
		50	D-001320-01-50
Lincode Non-targeting siRNA #2 <i>human, mouse, and rat</i>	Negative control siRNA with 3 or more mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13,400 g/mol	5	D-001320-02-05
		20	D-001320-02-20
		50	D-001320-02-50
Lincode Non-targeting siRNA #3 <i>human, mouse, and rat</i>	Negative control siRNA with 3 or more mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13,400 g/mol	5	D-001320-03-05
		20	D-001320-03-20
		50	D-001320-03-50
Lincode Non-targeting siRNA #4 <i>human, mouse, and rat</i>	Negative control siRNA with 3 or more mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13,400 g/mol	5	D-001320-04-05
		20	D-001320-04-20
		50	D-001320-04-50
Lincode Non-targeting Pool <i>human, mouse, and rat</i>	Negative control pool of four Non-targeting siRNAs. 5MW ~ 13,400 g/mol	5	D-001320-10-05
		20	D-001320-10-20
		50	D-001320-10-50

**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](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://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.