

LincodTM siRNA reagents

Product description

- Pooled and individual siRNAs designed using the SMARTselectionTM algorithm to silence long noncoding RNAs (lncRNAs)
- Lincod siRNA reagents are modified with a proprietary dual-strand modification to enhance specificity
- Annealed double-stranded RNA oligonucleotides
- 3'-UU overhangs on both strands
- 5'-Phosphate on antisense strand
- Mass of each strand confirmed by MALDI-TOF mass spectrometry

Product	Description	Cat. #
Lincod SMART-pool TM Reagent	<ul style="list-style-type: none"> • One tube containing a mixture of four SMARTselection-designed siRNAs targeting one gene • Sequence information provided 	R-XXXXXX-XX
Lincod set of 4 Upgrade Lincod individual siRNA	<ul style="list-style-type: none"> • Four individual SMARTselection—designed siRNAs from corresponding SMARTpool reagent • Sequence information provided 	RU-XXXXXX-XX N-XXXXXX-XX

Shipping and storage

- siRNA reagents are shipped as dry pellets at room temperature (23 °C). Under these conditions, they are stable for at least four weeks.
- Upon receipt, siRNA reagents should be stored at –20 °C to –80 °C. Under these conditions, they 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.
- 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.

Related products

- It is recommended to include a positive and negative control, such as Dharmacon RNAi Control Reagents, in every RNAi experiment. For more information, click [here](#).
- DharmaFECT™ siRNA Transfection Reagents are optimized for transfecting siRNA into a wide variety of cell lines. For more information, click [here](#).

Accompanying documents

- Basic siRNA resuspension protocol.

Supplemental documents

- Go to dharmacon.horizondiscovery.com to find:
- siRNA Recommended Reading List SMARTpool Journal Citations

References

References detailing the development of the SMARTselection algorithm:

1. Khvorova, A., A. Reynolds, *et al.* *Cell*, 2003. **115**(1): p. 209-216.
2. Reynolds, A., D. Leake, *et al.* *Nature Biotechnology*, 2004. **22**(3): p. 326-330.

For additional RNAi references please refer to the [siRNA Recommended Reading List](#).

Publication reference guide

When referencing the use of Dharmacon siRNA reagents, please include the following information: product name (either Lincode SMARTpool Reagent or siRNA) catalog number, 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.