

Dharmacon™ Accell™ siRNA Reagents

Please Read:

Dharmacon Accell siRNA is specially modified for use without a transfection reagent and works at a higher concentration than conventional siRNA with minimal disruption of the expression profile. When used with Accell siRNA Delivery Media, little to no delivery optimization is required.

Product Description:

- Chemically modified for stability, target specificity, and uptake by cells without a transfection reagent
- Pooled and individual siRNAs designed to silence specific target genes
- Annealed double-stranded RNA oligonucleotides
- 3'-UU overhangs
- 5'-Phosphate on antisense strand
- Mass of each strand confirmed by MALDI-TOF mass spectrometry

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:

- Accell Delivery Protocol

Supplemental Documents:

Click [here](#) to find:

- siRNA Recommended Reading List
- SMARTpool Journal Citations

Product	Description	Cat. #
Accell SMARTpool™ siRNA Reagent	<ul style="list-style-type: none"> • One tube containing a mixture of four siRNAs targeting one gene • Sequence information provided 	E-XXXXXX-XX
Accell Set of 4 siRNA Accell Upgrade Set of 4 siRNA Accell Individual siRNA	Individual siRNA(s) from corresponding SMARTpool reagent. <ul style="list-style-type: none"> • Sequence information provided 	EQ-XXXXXX-XX EU-XXXXXX-XX A-XXXXXX-XX
Accell siRNA Delivery Media	Media optimized and required for use with delivery of Accell siRNA.	B-005000-100 B-005000-500
5x siRNA buffer	For use in resuspending Accell siRNA. 300 mM KCl, 30 mM HEPES-pH 7.5, 1.0 mM MgCl ₂	B-002000-UB-100
RNase-free water	Molecular Grade RNase-free water for use in preparing 1x siRNA buffer.	B-003000-WB-100

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 – 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 nine months.

Related Products:

A positive and negative control is recommended in every RNAi experiment. For more information on Accell control products, visit gelifesciences.com/dharmacon.

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.

Publication Reference Guide:

When referencing the use of siRNA reagents, please include the following information: product name Accell SMARTpool™ or siRNA (Cat #), GE Healthcare Dharmacon, Inc., Lafayette, CO.

Each Accell siRNA product is covered by one or more of the following patents: US8252755, US8501706, US8188060 and US8415466

GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon, Inc., a General Electric company doing business as GE Healthcare. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2017 General Electric Company—All rights reserved. First version published April 2014. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK



Orders can be placed at:
gelifesciences.com/dharmacon

Customer Support: cs.dharmacon@ge.com
Technical Support: ts.dharmacon@ge.com or
1.800.235.9880; 303.604.9499 if you have any questions.