

DharmaconTM AccellTM RNAi control 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 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 positive control siRNAs designed to silence specific target genes (Table 1)
- Pooled and individual negative control siRNAs designed to not target any genes in human, mouse, or rat (Table 3)
- Accell Green and Red Control siRNAs are labeled with a fluorophore (Table 2)
- Annealed double-stranded RNA oligonucleotides
- 3'-UU overhangs
- 5'-Phosphate on antisense strand
- Mass of each strand confirmed by MALDI-TOF mass spectrometry
- The approximate molecular weight of all controls is 13,400 g/mol

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.

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

Accell siRNA Delivery media is required for use with every Accell siRNA product, and has been validated with over 40 cell types to date. For more information on Accell siRNA Delivery medium, go to dharmacon.horizondiscovery.com and search for Cat. #B-005000-100.

Accompanying documents

[Accell Delivery Protocol](#)

Publication reference guide

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

Table 1. Accession Numbers for Target Genes of Positive Control siRNAs.

Target gene	Human	Mouse	Rat
Cyclophilin B	NM_000942	NM_011149	NM_022536
GAPD	NM_002046	NM_008085	XM_575242

Table 2. Spectral Properties of Accell Green and Red Control siRNAs.

Accell Red & Green control siRNAs	Fluorophore absorption/emission Max	Extinction coefficient	Filter
Accell Green	6-FAM (494 nm /520 nm)	75,000 M ⁻¹ cm ⁻¹	FITC
Accell Red	DY-547 (557 nm /570 nm)	150,000 M ⁻¹ cm ⁻¹	Cy3, Rhodamine and PE

Table 3. Accell RNAi Negative Controls.

Product	Description	Cat. #
Accell Non-targeting siRNA #1-4	Negative control siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested.	D-001910-01-XX D-001910-02-XX D-001910-03-XX D-001910-04-XX
Accell Non-targeting Pool	Negative control pool of 4 siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested.	D-001910-10-XX
Accell Green Non-targeting siRNA Accell Red Non-targeting siRNA	Fluorescently-labeled negative control siRNA #1 with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested.	D-001950-01-XX D-001960-01-XX

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

Table 4. Accell RNAi Positive Controls.

Product	Description	Species	Cat. #
Accell Cyclophilin B Control siRNA	Positive silencing control for guaranteed silencing of Cyclophilin B mRNA.	H M R	D-001920-01-XX D-001920-02-XX D-001920-03-XX
Accell Cyclophilin B Control Pool	Pool of 4 positive silencing control siRNA for guaranteed silencing of Cyclophilin B mRNA.	H M R	D-001920-10-XX D-001920-20-XX D-001920-30-XX
Accell GAPD Control siRNA	Positive silencing control for guaranteed silencing of GAPD (GAPDH) mRNA.	H M R	D-001930-01-XX D-001930-02-XX D-001930-03-XX
Accell GAPD Control Pool	Pool of 4 positive silencing control siRNA for guaranteed silencing of GAPD (GAPDH) mRNA.	H M R	D-001930-10-XX D-001930-20-XX D-001930-30-XX
Accell eGFP Control siRNA	Positive control siRNA silencing control for guaranteed silencing of eGFP mRNA.	N/A	D-001940-01-XX
Accell eGFP Control Pool	Pool of 4 positive control siRNA silencing control for guaranteed silencing of eGFP mRNA.	N/A	D-001940-10-XX
Accell Green Cyclophilin B Control siRNA	Fluorescently-labeled positive silencing control for guaranteed silencing of Cyclophilin B mRNA.	H M R	D-001970-01-XX D-001970-02-XX D-001970-03-XX
Accell Red Cyclophilin B Control siRNA	Fluorescently-labeled positive silencing control for guaranteed silencing of Cyclophilin B mRNA.	H M R	D-001975-01-XX D-001975-02-XX D-001975-03-XX

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

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

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.

