

Dharmacon™ Accell™ siRNA Reagents

Please read

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

Product	Description	Cat.#
Accell SMARTpool™ siRNA Reagent	• 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. 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-002000UB-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.

Accompanying documents

· Accell Delivery Protocol

Knockdown guarantee

We guarantee at least 75% silencing (measured by qPCR) by the Accell SMARTpool and three of the Set of 4 individual siRNAs targeting your gene of interest. To learn about this guarantee click here.

Supplemental documents

Click here to find additional Accell siRNA resources:

- · Application notes
- · Protocols
- Recommended Reading lists
- Accell siRNA published citations

Related products

A positive and negative control is recommended in every RNAi experiment. For more information on Accell control products, visit https://dharmacon.horizondiscovery.com/rnai/controls/sirna/accell/#overview

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 Dharmacon siRNA reagents, please include the following information: Product brand name (siGENOME, ON-TARGETplus, Accell, Lincode, or custom) product format (SMARTpool or siRNA), Dharmacon catalog number, Horizon Discovery (Dharmacon), Lafayette, CO.

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

