

siRNA libraries

Product	Description	Cat. #
siGENOME™ SMARTpool™ siRNA Library ON-TARGET <i>plus</i> ™ SMARTpool siRNA Library Accell™ SMARTpool siRNA Library Lincode™ SMARTpool siRNA Library	<ul style="list-style-type: none"> Each well contains a mixture of four siRNAs targeting one gene Sequence information provided 	G-0XXXXX-XX G-1XXXXX-XX G-2XXXXX-XX G-3XXXXX-XX
siGENOME Set of 4 siRNA Library ON-TARGET <i>plus</i> Set of 4 siRNA Library Accell Set of 4 siRNA Library Lincode Set of 4 siRNA Library	<ul style="list-style-type: none"> Four individual siRNAs from corresponding SMARTpool reagent Sequence information provided 	GU-0XXXXX-XX GU-1XXXXX-XX GU-2XXXXX-XX GU-3XXXXX-XX
5x siRNA Buffer	<ul style="list-style-type: none"> Recommended for the resuspension of siRNA 300 mM KCl, 30 mM HEPES-pH 7.5, 1.0 mM MgCl₂ 	B-002000-UB-100

USB drive with detailed plate map information provided with all products

Product description

- Pooled and individual siRNAs designed to silence specific target genes
- ON-TARGET*plus* and Lincode siRNA reagents are modified with a patented chemical modification to enhance target specificity and reduce off-target effects
- Accell reagents are chemically modified for stability, target specificity, and uptake by cells without a transfection reagent
- Sequence information provided with all products
- 5'-Phosphate on antisense strand
- Mass of each strand confirmed by ESI-MS (electrospray ionization 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.

Functional guarantee

Each Accell, siGENOME and ON-TARGET*plus* SMARTpool reagent or 3 of 4 individuals is guaranteed to silence target gene expression at the mRNA level by at least 75% when used under optimized transfection and detection conditions.

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, the siRNA 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). For short term storage, RNase-free water (Cat # B-003000-WB-100) 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 (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 one year.

Supplemental documents

- Delivery protocol for Accell siRNA reagents <http://dharmacon.horizondiscovery.com/uploadedFiles/Resources/accell-delivery-protocol.pdf>
- Detailed recommendations for resuspension and handling of siRNA in multi-well plates <http://dharmacon.horizondiscovery.com/uploadedfiles/resources/dharmacon-sirna-libraries-guidelines.pdf>

Recommended experimental controls

We recommend including a positive and negative control, such as siGENOME, ON-TARGET*plus*, and Accell RNAi Control reagents, in every RNAi experiment. For more information, go to <http://dharmacon.horizondiscovery.com/dharmacon/rnaicontrols>.

Publication reference guide

When referencing the use of siRNA Libraries, please include the following information: siGENOME siRNA Library, ON-TARGET*plus* siRNA Library, or Accell siRNA Library (specific gene collection, 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.

ON-TARGET*plus* is covered by patent: US7595387.

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