

# siGLO<sup>™</sup>, siGENOME<sup>™</sup>, and ON-TARGET*plus*<sup>™</sup> control reagents

# **Product description**

- Double-stranded, chemically synthesized oligonucleotides
- Antisense strands are modified with 5'-phosphate
- siGLO reagents are labeled with a fluorophore (see Table 2) on the sense strand
- Mass of each strand confirmed by MALDI-TOF mass spectrometry
- Duplex integrity confirmed by non-denaturing polyacrylamide gel electrophoresis

# **Shipping and storage**

- RNAi Control reagents are shipped as dried pellets at room temperature (23 °C). Under these conditions, they are stable for at least four weeks.
- Upon receipt, our RNAi Controls and siGLO reagents should be stored at -20 °C to -80 °C. In addition, store siGLO siRNAs in the dark to avoid exposure to strong light. Under these conditions, the reagents 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.
- 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.

# **Accompanying document**

Basic siRNA resuspension protocol.

### **Supplemental documents**

Go to dharmacon.horizondiscovery.com to find basic transfection and resuspension protocols.

### **Related products**

DharmaFECT siRNA Transfection Reagents are available in four formulations that are optimized for transfecting siRNA into a wide variety of cell lines. For more information, click here.

# **Publication reference guide**

When referencing the use of the siRNA reagents listed in this document, please include the following information: Control type, catalog number, Horizon Discovery, Lafayette, CO.

# dharmacon.horizondiscovery.com

#### Table 1. Accession numbers for target genes of positive control siRNA

Target gene	Human	Mouse	Rat
Cyclophilin B	NM_000942	NM_011149	NM_022536
Lamin A/C	NM_005572, NM_170707, NM_170708 (Targets all three isoforms)	NM_019390, NM_001002011 (Targets both isoforms)	NM_001002016
GAPD	NM_002046	_	_

### Table 2. Spectral properties of siGLO reagents

Absorption max.	Emission max.	Extinction coefficient	Excitation filters
557 nM	570 nM	150,000 M-1 cm-1	Cy3™, Rhodamine, or PE

### Table 3. siGLO reagents

Product	Description*	Cat. #
siGLO Cyclophilin B Control siRNA human/mouse/rat	Dual purpose silencing control tar- geting, Cyclophilin B with fluorescent label for human, mouse and rat cells. MW ~ 14,000 g/mol	D-001610- 01-XX
siGLO Lamin A/C Control siRNA <i>human/mouse/rat</i>	Dual purpose silencing control targeting, Lamin A/C with fluorescent label for human, mouse and rat cells. MW ~ 14,100 g/mol	D-001620- 01-XX
	Dual purpose silencing control targeting Lamin A/C with fluorescent label for human cells. MW ~ 14,100 g/mol	D-001620- 02-XX
siGLO Lamin A/C Control siRNA human/mouse/rat	Dual purpose silencing control targeting Lamin, A/C with fluorescent label for mouse cells. MW ~ 14,100 g/mol	D-001620- 03-XX
	Dual purpose silencing control targeting Lamin, A/C with fluorescent label for rat cells. MW ~ 14,100 g/mol	D-001620- 04-XX
siGLO RISC-Free™ Control siRNA	Non-targeting siRNA with fluorescent label and impaired processing and uptake by RISC; a useful co-transfec- tion marker.	D-001600- 01-XX

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

### Table 4. siGENOME control siRNA

		<u> </u>
Product	Description*	Cat. #
siGENOME Lamin A/C Control siRNA human/mouse/rat	Positive silencing control for guaran- teed silencing of Lamin A/C mRNA in human, mouse and rat cells. MW ~ 13,400 g/mol	D-001050- 01-XX
siGENOME Cyclophilin B Control siRNA human/mouse/rat	Positive silencing control for guaranteed silencing of Cyclophilin B mRNA in human, mouse and rat cells. MW ~ 13,300 g/mol	D-001136- 01-XX
siGENOME GAPD Control siRNA human	Positive silencing control for guaranteed silencing of GAPD (GAPDH) mRNA in human cells. MW ~ 13,300 g/mol	D-001140- 01-XX
siGENOME Non- targeting siRNA #1	Negative control siRNA with at least 4 mismatches to any human, mouse or rat gene. Microarray tested.^ MW ~ 13,300 g/mol	D-001210- 01-XX
siGENOME Non- targeting siRNA #2	Negative control siRNA with at least 4 mismatches to any human, mouse or rat gene. Microarray tested.^ MW ~ 13,300 g/mol	D-001210- 02-XX
siGENOME Non- targeting siRNA #3	Negative control siRNA with at least 4 mismatches to any human, mouse or rat gene. Microarray tested.^ MW ~ 13,300 g/mol	D-001210- 03-XX
siGENOME Non- targeting siRNA #4	Negative control siRNA with at least 4 mismatches to any human, mouse or rat gene. Microarray tested.^ MW ~ 13,400 g/mol	D-001210- 04-XX
siGENOME Non- targeting siRNA #5	Negative control siRNA with at least 4 mismatches to any human, mouse or rat gene. Microarray tested.^ MW ~ 13,400 g/mol	D-001210- 05-XX
siGENOME Non-targeting siRNA Pool #1	Pool of 4 Non-Targeting siRNAs. Useful control for experiments with SMARTpool reagents.^ MW ~ 13,400 g/mol	D-001206- 13-XX
siGENOME Non-targeting siRNA Pool #2	Pool of 4 Non-Targeting siRNAs. Useful control for experiments with SMARTpool reagents.^ MW ~ 13,400 g/mol	D-001206- 14-XX
siGENOME RISC-Free Control siRNA	Non-Targeting siRNA with impaired ability for RISC interaction.	D-001220- 01-XX

^Non-Targeting siRNA #1 reduces EGFR mRNA by ~ 50% (Accession NM\_005228), Non-Targeting siRNA #2 reduces firefly luciferase mRNA by ~ 75% (Cat. #U47296), and the Non-Targeting Pool includes both Non-Targeting siRNAs. #1 and #2 and targets both genes as noted above. \*XX=05,20, or 50 for 5, 20, and 50 nmol amounts.

# dharmacon.horizondiscovery.com

Table 5. ON-TARGET			Product	Description*	Cat. #
Product ON-TARGET <i>plus</i> Cyclophilin B Control siRNA <i>human/mouse/rat</i>	Description* Positive silencing control for guaranteed silencing of Cyclophilin B mRNA in human cells. MW ~ 13,400 g/mol	Cat. # D-001820-01-XX	ON-TARGET <i>plus</i> GAPD Control Pool human/mouse/rat	Pool of 4 positive silencing control siRNA for guaranteed silencing of GAPD (GAPDH) mRNA in human cells. MW ~ 13,300 g/mol	D-001830-10-XX
	Positive silencing control for guaranteed silencing of Cyclophilin B mRNA in mouse cells. MW ~ 13,400 g/mol	D-001820-02-XX		Pool of 4 positive silencing control siRNA for guaranteed silencing of GAPD (GAPDH) mRNA in mouse cells. MW ~ 13,300 g/mol	D-001830-20-X>
	Positive silencing control for guaranteed silencing of Cyclophilin B mRNA in rat cells. MW ~ 13,400 g/mol	D-001820-03-XX		Pool of 4 positive silencing control siRNA for guaranteed silencing of GAPD (GAPDH) mRNA in rat cells. MW ~ 13,300 g/mol	D-001830-30-XX
ON-TARGET <i>plus</i> - Cyclophilin B Control Pool <i>human/mouse/rat</i>	Pool of 4 positive silencing control siRNA for guaranteed silencing of Cyclophilin B mRNA in human cells. Useful control for experiments with SMARTpool reagents.	D-001820-10-XX	ON-TARGET <i>plus</i> Non-targeting siRNA #1 <i>human/mouse/rat</i>	Negative control siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13.400 g/mol	D-001810-01-XX
	MW ~ 13,400 g/mol Pool of 4 positive silencing control siRNA for guaranteed silencing of Cyclophilin B mRNA in mouse cells. Useful control for experiments with SMARTpool reagents. MW ~ 13,400 g/mol		ON-TARGET <i>plus</i> Non-targeting siRNA #2 <i>human/mouse/rat</i>	Negative control siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13.400 g/mol	D-001810-02-XX
		D-001820-20-XX	ON-TARGET <i>plus</i> Non-targeting siRNA #3	Negative control siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested.	D-001810-03-XX
	Pool of 4 positive silencing control siRNA for guaranteed silencing of Cyclophilin B mRNA in rat cells. Useful control for experiments with SMARTpool reagents. MW ~ 13,400 g/mol	D-001820-30-XX	human/mouse/rat ON-TARGET <i>plus-</i> Non-targeting siRNA #4 human/mouse/rat	MW ~ 13.400 g/mol Negative control siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested. MW ~ 13.400 g/mol	D-001810-04-XX
ON-TARGET <i>plus</i> GAPD Control siRNA <i>human/mouse/rat</i>	Positive silencing control for guar- anteed silencing of GAPD (GAPDH) mRNA in human cells. MW ~ 13,300 g/mol	D-001830-01-XX	ON-TARGET <i>plus</i> Non-targeting Pool <i>human/mouse/rat</i>	Negative control siRNA with at least 4 mismatches to any human, mouse, or rat gene. Microarray tested. 5MW ~ 13.400 g/mol	D-001810-10-XX
	Positive silencing control for guar- anteed silencing of GAPD (GAPDH) mRNA in mouse cells. MW ~ 13,300 g/mol	D-001830-02-XX	*XX=05,20, or 50 for 5, 20, and 50 nmol amounts.		
	Positive silencing control for guar- anteed silencing of GAPD (GAPDH) mRNA in rat cells.	D-001830-03-XX			

### If you have any questions, contact

+44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA) t

MW ~ 13,300 g/mol

+ 44 (0)1223 655 581 f

**w** horizondiscovery.com/contact-us **or** dharmacon.horizondiscovery.com/service-and-support Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

horizon INSPIRED CELL SOLUTIONS

©2018 Horizon Discovery Group Company—All rights reserved. All trademarks are the property of Horizon Discovery Company unless otherwise specified. First published June 2015. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

ΤM