

Dharmacon™

RNAi, Gene Expression & Gene Editing

Decode™ Pooled Lentiviral shRNA Screening: Decode Indexing PCR and Sequencing Primer Kit

Cat. #RHS5339

Product Description:

Dharmacon Decode Pooled Lentiviral shRNA Screening Libraries are an efficient and cost-effective method for screening large numbers of shRNAs without costly automation. Decode library construction, pooling techniques, and NGS-compatible screening workflows have been experimentally validated to ensure reproducibility and accurate hit identification. Decode libraries are offered for a variety of targeted gene families as well as druggable genome and whole genome to fit many experimental designs.

The Decode Pooled shRNA screening system includes Illumina™-adapted PCR primers for identification of shRNA hairpin sequences from genomic DNA (gDNA) by high-throughput sequencing on Illumina instrumentation. Decode primer pairs are optimized for efficient amplification of shRNA while minimizing thermodynamic bias and variation in representation. In addition to efficiently amplifying shRNA hairpins from gDNA, Decode PCR primers have built-in adaptor and index sequences that allow the researcher to easily move from PCR amplification to Illumina™ highthroughput sequencing. Direct identification of shRNA passenger strand facilitates data analysis and ensures accurate target gene identification.

The Decode Indexing PCR and Sequencing Primer Kit contains the following 15 DNA oligonucleotides:

- 1 × 600 µL tube of Decode Forward PCR Primer
- 1 × 600 µL tube of each Decode Reverse Index 1-12 PCR Primers
- 1 × 100 µL tube of Decode Read 1 Sequencing Primer
- 1 × 100 µL tube of Decode Index Read Sequencing Primer

Concentration:

Cat. #	Description	Item #	Concentration of oligo solution (µM)	Volume per aliquot (µL)
PRM6178	Decode Indexing PCR and Sequencing Primer Kit			
	Decode Forward PCR Primer (600 µL)	PRM6163	50	600
	Decode Reverse Index PCR Primer 1 (600 µL)	PRM6164	50	600
	Decode Reverse Index PCR Primer 2 (600 µL)	PRM6165	50	600
	Decode Reverse Index PCR Primer 3 (600 µL)	PRM6166	50	600
	Decode Reverse Index PCR Primer 4 (600 µL)	PRM6167	50	600
	Decode Reverse Index PCR Primer 5 (600 µL)	PRM6168	50	600
	Decode Reverse Index PCR Primer 6 (600 µL)	PRM6169	50	600
	Decode Reverse Index PCR Primer 7 (600 µL)	PRM6170	50	600
	Decode Reverse Index PCR Primer 8 (600 µL)	PRM6171	50	600
	Decode Reverse Index PCR Primer 9 (600 µL)	PRM6172	50	600
	Decode Reverse Index PCR Primer 10 (600 µL)	PRM6173	50	600
	Decode Reverse Index PCR Primer 11 (600 µL)	PRM6174	50	600
	Decode Reverse Index PCR Primer 12 (600 µL)	PRM6175	50	600
	Decode Read 1 Illumina Sequencing Primer (100 µL)	PRM6176	100	100
	Decode Index Read Illumina Sequencing Primer (100 µL)	PRM6177	100	100

Shipping and Storage:

Store at -20° C. The Decode Indexing PCR and Sequencing Primer Kit is shipped at ambient temperature and should be stored in a -20° C freezer. Components provided with this kit are for single use only.

Illumina is a trademark of Illumina, Inc. GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon is a trademark of GE Healthcare companies. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2014 General Electric Company—All rights reserved. Version published June 2014. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK



GE Healthcare

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.