

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™ high-throughput 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

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

Illumina is a trademark of Illumina, Inc. All trademarks are the property of Horizon Discovery Company unless otherwise specified.
©2018 Horizon Discovery Group Company—All rights reserved. First published June 2014. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

Concentration

Table 1. Decode Indexing PCR and Sequencing Primer Kit (Cat. # PRM6178)

Description	Item #	Concentration of oligo solution (µM)	Volume per aliquot (µL)
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.