

# Dharmacon™ Edit-R™ Lentiviral sgRNA Pooled Screening Libraries:

## Edit-R Pooled sgRNA Indexing PCR and Sequencing Primer Kits

### Cat. #PRM10184 & PRM10185

#### Product Description:

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

The Edit-R Lentiviral sgRNA Pooled Library screening system includes Illumina™-adapted PCR primers for identification of sgRNA targeting sequences from genomic DNA (gDNA) by high-throughput sequencing on Illumina instrumentation. Edit-R primer pairs are optimized for efficient amplification of sgRNA while minimizing thermodynamic bias and variation in representation. In addition to efficiently amplifying sgRNA sequences from gDNA, Edit-R 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 sgRNA sequences facilitates data analysis and ensures accurate target gene identification.

The Edit-R Pooled sgRNA Indexing PCR and Sequencing Primer Kits A & B contain the following 15 DNA oligonucleotides (See tables for kit-specific part numbers):

- 1 × 600 µL tube of Edit-R Pooled sgRNA Forward PCR Primer
- 1 × 600 µL tube of each of 12 Edit-R Pooled sgRNA Reverse Index PCR Primers
- 1 × 100 µL tube of Edit-R Pooled sgRNA Read 1 Sequencing Primer
- 1 × 100 µL tube of Edit-R Pooled sgRNA Index Read Sequencing Primer

#### Concentration:

Cat. #	Description	Item #	Concentration of oligo solution (µM)	Volume per aliquot (µL)
<b>PRM10184</b>	<b>Edit-R Pooled sgRNA Indexing PCR and Sequencing Primer Kit A</b>			
	Edit-R Pooled sgRNA Forward PCR Primer	PRM10180	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 2	PRM10157	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 4	PRM10159	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 5	PRM10160	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 6	PRM10161	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 7	PRM10162	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 12	PRM10167	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 13	PRM10168	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 14	PRM10169	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 15	PRM10170	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 16	PRM10171	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 18	PRM10172	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 19	PRM10173	50	600
	Edit-R Pooled sgRNA Index Read Sequencing Primer	PRM10183	100	100
	Edit-R Pooled sgRNA Read 1 Sequencing Primer	PRM10181	100	100



Cat. #	Description	Item #	Concentration of oligo solution (μM)	Volume per aliquot (μL)
PRM10185	<b>Edit-R Pooled sgRNA Indexing PCR and Sequencing Primer Kit B</b>			
	Edit-R Pooled sgRNA Forward PCR Primer	PRM10180	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 1	PRM10156	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 3	PRM10158	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 8	PRM10163	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 9	PRM10164	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 10	PRM10165	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 11	PRM10166	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 20	PRM10174	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 21	PRM10175	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 22	PRM10176	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 23	PRM10177	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 25	PRM10178	50	600
	Edit-R Pooled sgRNA Reverse Index PCR Primer 27	PRM10179	50	600
	Edit-R Pooled sgRNA Index Read Sequencing Primer	PRM10183	100	100
	Edit-R Pooled sgRNA Read 1 Sequencing Primer	PRM10181	100	100

## Shipping and Storage:

Store at -20° C. The Edit-R Pooled sgRNA 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, Inc., a General Electric company doing business as GE Healthcare. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2016 General Electric Company—All rights reserved. Version published November 2016. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK



Orders can be placed at:  
[gelifsciences.com/dharmacon](http://gelifsciences.com/dharmacon)

Customer Support: [cs.dharmacon@ge.com](mailto:cs.dharmacon@ge.com)  
 Technical Support: [ts.dharmacon@ge.com](mailto:ts.dharmacon@ge.com)  
 1.800.235.9880; 303.604.9499 if you have any questions.