

DharmaconTM Edit-RTM Lentiviral sgRNA pooled screening libraries: Edit-R pooled sgRNA indexing PCR and sequencing primer kits

Cat. #PRM10184 & PRM10185

Product description

DharmaconTM Edit-RTM 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 IlluminaTM-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 IlluminaTM 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)
PRM10184	Edit-R Pooled sgRNA Indexing PCR and Sequencing Primer Kit A			
	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
PRM10185	Edit-R Pooled sgRNA Indexing PCR and Sequencing Primer Kit 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.

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. ©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.

horizon[™]
INSPIRED CELL SOLUTIONS

V20918