

DharmaconTM SMARTvectorTM lentiviral shRNA pooled libraries: SMARTvector indexing PCR and sequencing primer kits A & B

Cat. #PRM7668 & PRM10186

Product description

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

The SMARTvector Lentiviral shRNA Pooled Library screening system includes IlluminaTM-adapted PCR primers for identification of shRNA hairpin sequences from genomic DNA (gDNA) by high-throughput sequencing on Illumina instrumentation. SMARTvector 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, SMARTvector 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 shRNA passenger strand facilitates data analysis and ensures accurate target gene identification.

The SMARTvector Indexing PCR and Sequencing Primer Kits A & B contain the following 15 DNA oligonucleotides (See tables for kit-specific part

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

Concentration

Cat. #	Description	Item #	Concentration of oligo solution (μM)	Volume per aliquot (μL)
PRM7668	SMARTvector Indexing PCR and Sequencing Primer Kit A			
	SMARTvector Forward PCR Primer	PRM10211	50	600
	SMARTvector Reverse Index PCR Primer 2	PRM10188	50	600
	SMARTvector Reverse Index PCR Primer 4	PRM10190	50	600
	SMARTvector Reverse Index PCR Primer 5	PRM10191	50	600
	SMARTvector Reverse Index PCR Primer 6	PRM10192	50	600
	SMARTvector Reverse Index PCR Primer 7	PRM10193	50	600
	SMARTvector Reverse Index PCR Primer 12	PRM10198	50	600
	SMARTvector Reverse Index PCR Primer 13	PRM10199	50	600
	SMARTvector Reverse Index PCR Primer 14	PRM10200	50	600
	SMARTvector Reverse Index PCR Primer 15	PRM10201	50	600
	SMARTvector Reverse Index PCR Primer 16	PRM10202	50	600
	SMARTvector Reverse Index PCR Primer 18	PRM10203	50	600
	SMARTvector Reverse Index PCR Primer 19	PRM10204	50	600
	SMARTvector Index Read Sequencing Primer	PRM10214	100	100
	SMARTvector Read 1 Sequencing Primer	PRM10212	100	100
PRM10186	SMARTvector Indexing PCR and Sequencing Primer Kit B			
	SMARTvector Forward PCR Primer	PRM10211	50	600
	SMARTvector Reverse Index PCR Primer 1	PRM10187	50	600
	SMARTvector Reverse Index PCR Primer 3	PRM10189	50	600
	SMARTvector Reverse Index PCR Primer 8	PRM10194	50	600
	SMARTvector Reverse Index PCR Primer 9	PRM10195	50	600
	SMARTvector Reverse Index PCR Primer 10	PRM10196	50	600
	SMARTvector Reverse Index PCR Primer 11	PRM10197	50	600
	SMARTvector Reverse Index PCR Primer 20	PRM10205	50	600
	SMARTvector Reverse Index PCR Primer 21	PRM10206	50	600
	SMARTvector Reverse Index PCR Primer 22	PRM10207	50	600
	SMARTvector Reverse Index PCR Primer 23	PRM10208	50	600
	SMARTvector Reverse Index PCR Primer 25	PRM10209	50	600
	SMARTvector Reverse Index PCR Primer 27	PRM10210	50	600
	SMARTvector Index Read Sequencing Primer	PRM10214	100	100
	SMARTvector Read 1 Sequencing Primer	PRM10212	100	100

Shipping and storage

Store at -20 °C. The SMARTvector 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. All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2018 Horizon Discovery Group Company—All rights reserved. First published July 2014. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.