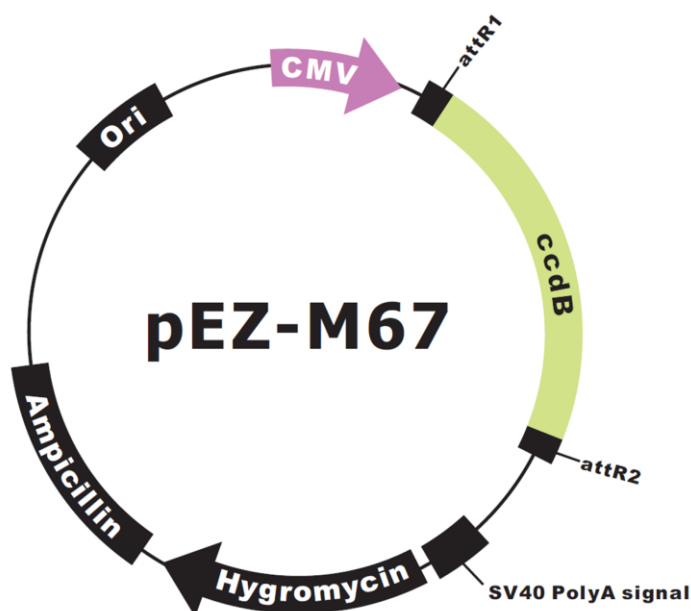


EZShuttle™ Recombination Cloning Expression Vector pEZ-M67

7662 bp, circular



Vector Features

Feature	Position
T7 promoter	41-59
SV40 polyA	358-625
f1 intergenic region	751-1005
SV40 early promoter and origin	1366-1675
Hygromycin resistance ORF	1733-2758
Synthetic polyA	2811-2859
Ampicillin resistance ORF	3269-4129
pUC Ori	4275-4948
CMV promoter	5349-5942
attR1	5966-6090
CmR	6199-6858
ccdB	7200-7505
attR2	7546-7670

Sequencing Primers

Primer	Sequence
Forward	5'-CAGCCTCCGGACTCTAGC-3'
Reverse	5'-TAATACGACTCACTATAGGG-3'

GeneCopoeia, Inc.
9620 Medical Center Drive, #101
Rockville, MD 20850
Tel: 301-762-0888 Fax: 301-762-3888
Email: inquiry@genecopoeia.com Web: www.genecopoeia.com

GeneCopoeia products are for research use only

© 2015, GeneCopoeia, Inc