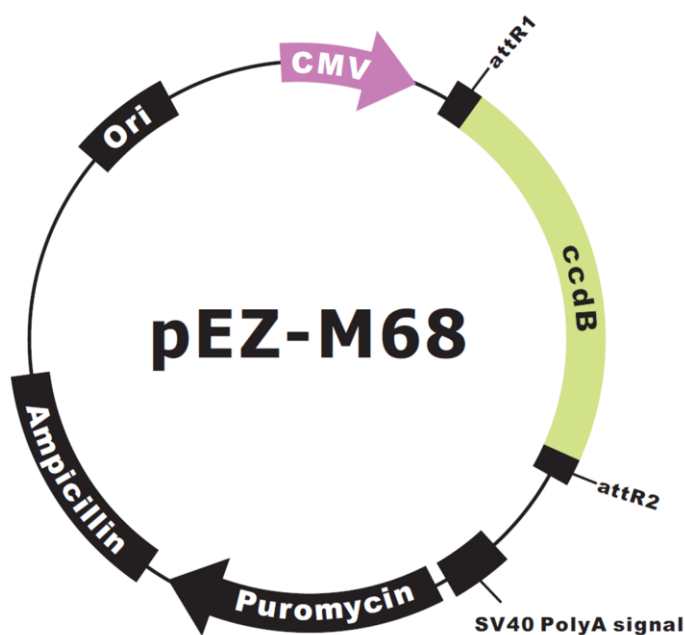


EZShuttle™ Recombination Cloning Expression Vector pEZ-M68

7238 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	1733-2332
Synthetic polyA	2387-2435
Ampicillin resistance ORF	2845-3705
pUC Ori	3851-4524
CMV promoter	4925-5518
attR1	5542-5666
CmR	5775-6434
ccdB	6776-7018
attR2	7122-7246

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