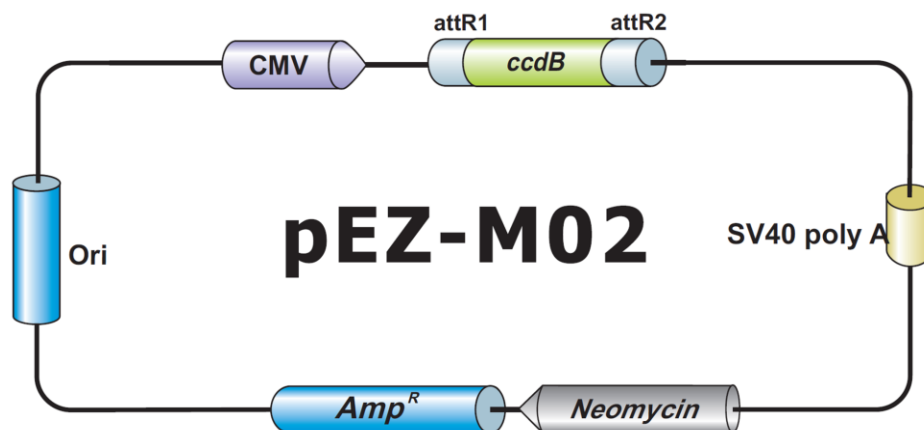


EZShuttle™ Recombination Cloning Expression Vector pEZ-M02

7445 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
Neomycin resistance gene	1733 2527
Synthetic polyA	2594 2642
Ampicillin resistance ORF	3052 3912
pUC Ori	4058 4731
CMV promoter	5132 5725
attR1	5749 5873
CmR	5982 6641
ccdB	6983 7288
attR2	7329-7453

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