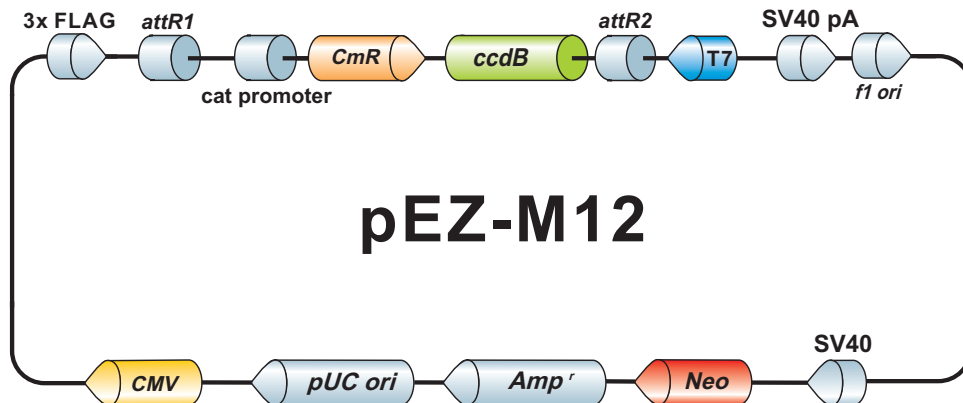


EZShuttle™ Recombination Cloning Expression Vector pEZ-M12

7660bp, circular



Vector Features

Name	Position
T7 promoter(complementary strand)	16 34
SV40 poly A	333 600
f1 ori	726 980
SV40 pomoter	1341 1650
Neomycin resistance gene	1708 2502
Ampicillin resistance gene	3027 3887
pUC ori	4033 4706
CMV promoter	5107 5700
3xFLAG	5726 5791
attR1	5798 5922
cat promoter	6069 6171
CmR	6172 6831
ccdB	7173 7478
attR2	7519 7643

Sequencing Primers

Primer	Sequence
Forward	5'-cagcctccggactctagc-3'
Reverse	5'-taatcgactcactataggg-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

© 2016, GeneCopoeia, Inc