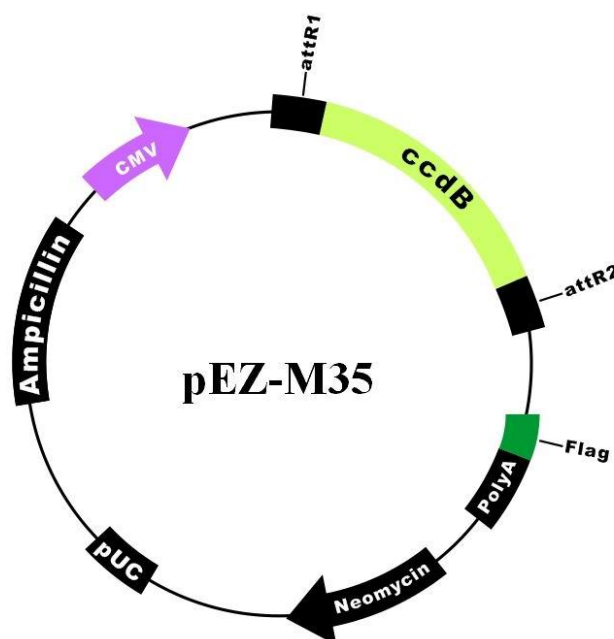


EZShuttle™ Recombination Cloning Expression Vector

pEZ-M35

7376 bp, circular



Vector Features

Feature	Position
Flag®	6 29
bGH polyA	43 267
f1 ori	313 741
SV40 promoter	755 1084
Neomycin resistance ORF	1151 1945
SV40 polyA	2121 2242
pUC Ori	2693 3281
Ampicillin resistance ORF (complementary strand)	3452 4312
CMV promoter	5199 5402
attR1	5683 5807
CmR	5916 6575
ccdB	6905 7210
attR2	7251 7374

Sequencing Primers

Primer	Sequence
Forward	5'-AGGCACTGGGCAGGTAAG-3'
Reverse	5'-GTGGCACCTTCCAGGGTC-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

© 2020, GeneCopoeia, Inc