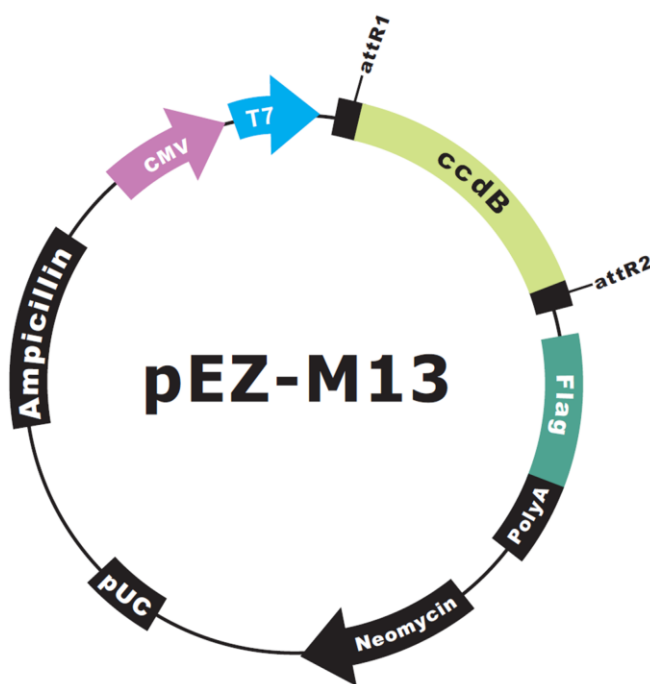


EZShuttle™ Recombination Cloning Expression Vector pEZ-M13

7080 bp, circular



Vector Features

Feature	Position
Flag®	19 54
bGH polyA	62 289
f1 ori	335 763
SV40 promoter	790 1099
Neomycin resistance ORF	1173 1967
SV40 polyA	2144 2273
pUC Ori	2656 3329
Ampicillin resistance ORF	3474 4334
CMV promoter	4702 5329
T7 promoter	5333 5352
attR1	5384 5508
CmR	5617 6276
ccdB	6618 6923
attR2	6964-7453

Sequencing Primers

Primer	Sequence
Forward	5'-GCGGTAGGCGTGTACGG-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

