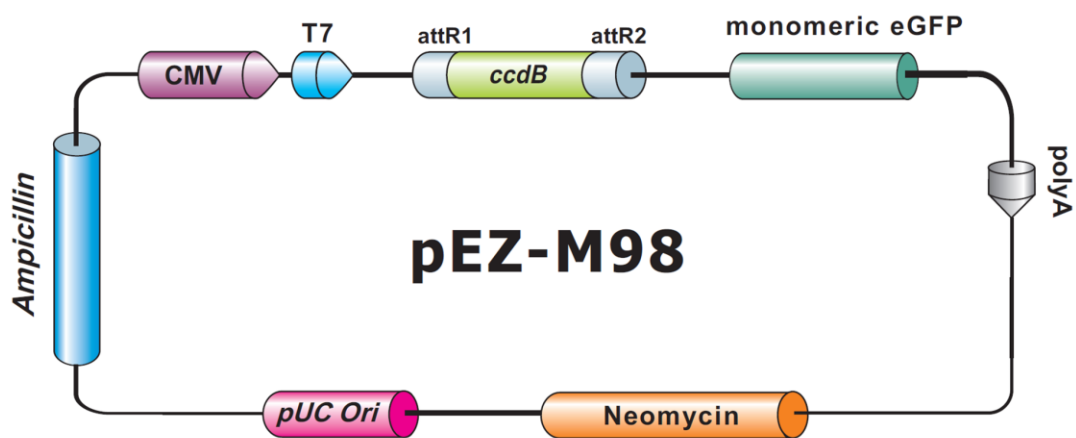


EZShuttle™ Recombination Cloning Expression Vector pEZ-M98

7445 bp, circular



Vector Features

| Feature | Position |
|--------------------------------|-----------|
| monomeric eGFP | 28-744 |
| bGH polyA | 964-1191 |
| f1 ori | 1237-1665 |
| SV40 early promoter and origin | 1692-2001 |
| Neomycin resistance gene | 2075-2869 |
| SV40 polyA | 3046-5175 |
| pUC Ori | 3558-4267 |
| Ampicillin resistance ORF | 4376-5236 |
| CMV promoter | 5604-6191 |
| T7 promoter | 6235-6254 |
| attR1 | 6286-6410 |
| CmR | 6519-7178 |
| ccdB | 7520-7825 |
| attR2 | 7866-7990 |

Sequencing Primers

| Primer | Sequence |
|---------|---------------------------|
| Forward | 5'-GCGGTAGGCGTGTACGGT-3' |
| Reverse | 5'-CCGGACACGCTGAACTTGT-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