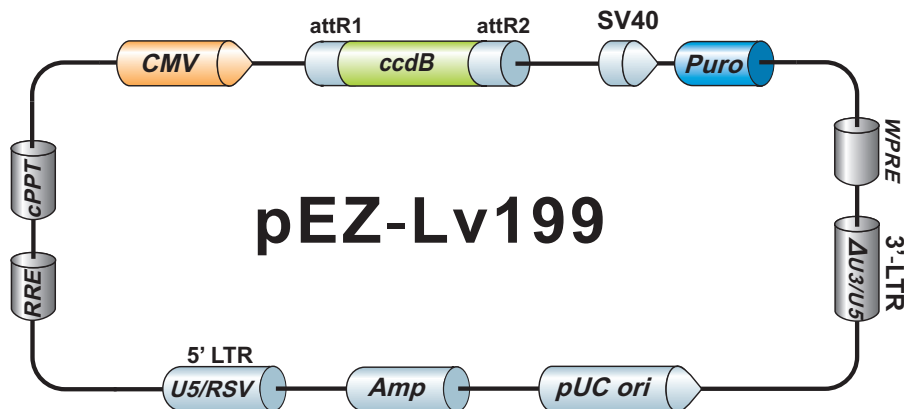


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv199

8921 bp, circular



Vector Features

| Name | Position |
|---|-----------|
| SV40 promoter | 1 368 |
| Puromycin resistance gene | 403 1002 |
| WPRE | 1016 1607 |
| 3'-LTR(delta U3/U5) | 1810 2046 |
| pUC ori (complementary strand) | 2505 3175 |
| Ampicillin resistance gene (complementary strand) | 3320 4180 |
| 5'-LTR(RSV/U5) | 4595 5006 |
| RRE | 5677 5910 |
| cPPT | 6402 6525 |
| CMV promoter | 6622 7216 |
| attR1 | 7223 7347 |
| CmR | 7456 8115 |
| ccdB | 8445 8750 |
| attR2 | 8791 8914 |

Sequencing Primers

| Primer | Sequence |
|---------|---------------------------------|
| Forward | 5'- cggtgggaggtctatataagcag -3' |
| Reverse | 5'- ctggaatagctcagaggc -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