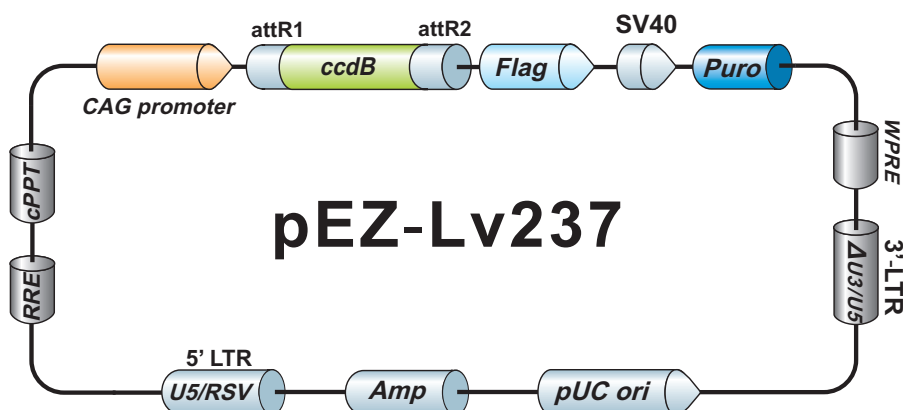


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv237

10087 bp, circular



Vector Features

Name	Position
Flag	1 24
SV40 promoter	33 400
Puromycin resistance gene	435 1034
WPRE	1048 1639
3'-LTR(delta U3/U5)	1842 2078
pUC ori (complementary strand)	2537 3207
Ampicillin resistance gene (complementary strand)	3352 4212
5'-LTR(RSV/U5)	4627 5038
RRE	5709 5942
cPPT	6434 6557
CAG promoter	6660 8376
attR1	8389 8513
CmR	8622 9281
ccdB	9611 9916
attR2	9957 10080

Sequencing Primers

Primer	Sequence
Forward	5'- ggcttctggcgtgtgacc -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