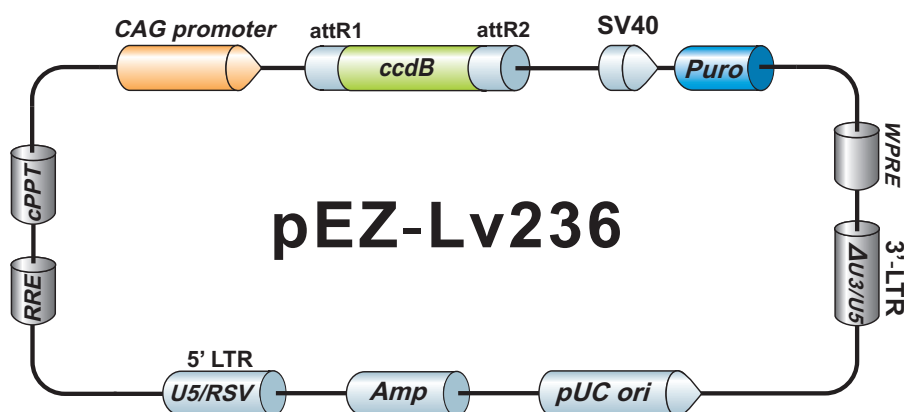


## EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv236

10055 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
CAG promoter	6628 8344
attR1	8357 8481
CmR	8590 9249
ccdB	9579 9884
attR2	9925 10048

### 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](mailto:inquiry@genecopoeia.com) Web: [www.genecopoeia.com](http://www.genecopoeia.com)

GeneCopoeia products are for research use only

© 2015, GeneCopoeia, Inc