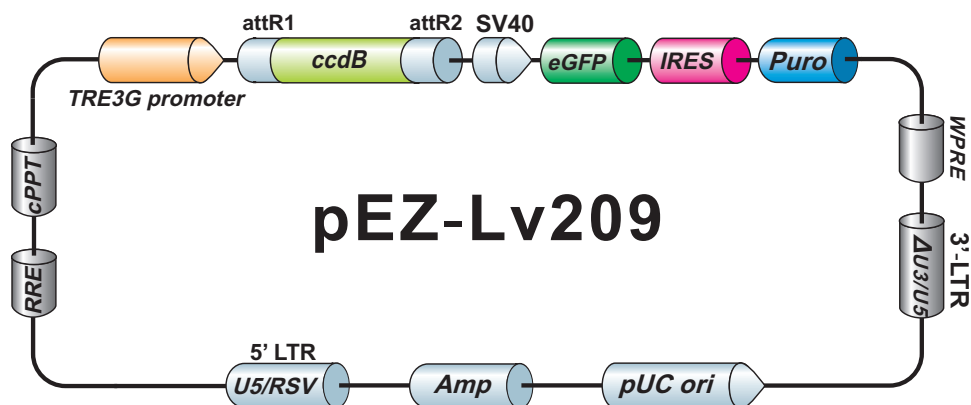


## EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv209

10059 bp, circular



### Vector Features

Name	Position
SV40 promoter	7 374
eGFP	404 1123
IRES	1137 1713
Puromycin resistance gene	1746 2345
WPRE	2359 2950
3'-LTR(delta U3/U5)	3156 3392
pUC Ori (complementary strand)	3851 4521
Ampicillin resistance gene (complementary strand)	4666 5526
5'-LTR(RSV/U5)	5941 6352
RRE	7023 7256
cPPT	7748 7871
TRE3G promoter	7968 8359
attR1	8367 8491
CmR	8600 9259
ccdB	9589 9894
attR2	9935 10058

### Sequencing Primers

Primer	Sequence
Forward	5'- GCTCGTTTAGTGAACCGTC -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