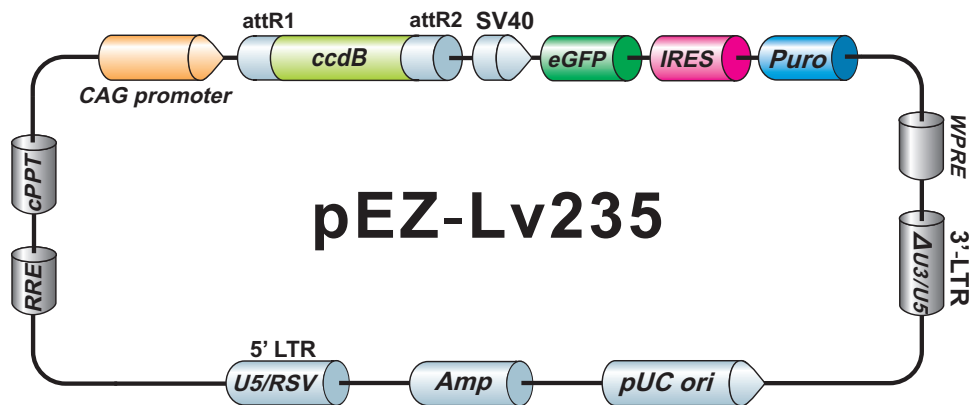


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv235

11397 bp, circular



Vector Features

Name	Position
CAG promoter	1 1717
attR1	1730 1854
CmR	1963 2622
ccdB	2952 3257
attR2	3298 3421
SV40 promoter	3429 3796
eGFP	3831 4550
IRES	4564 5140
Puromycin resistance gene	5173 5772
WPRE	5786 6377
3'-LTR(delta U3/U5)	6580 6816
pUC Ori (complementary strand)	7275 7945
Ampicillin resistance gene (complementary strand)	8090 8950
5'-LTR(RSV/U5)	9365 9776
RRE	10447 10680
cPPT	11172 11295

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