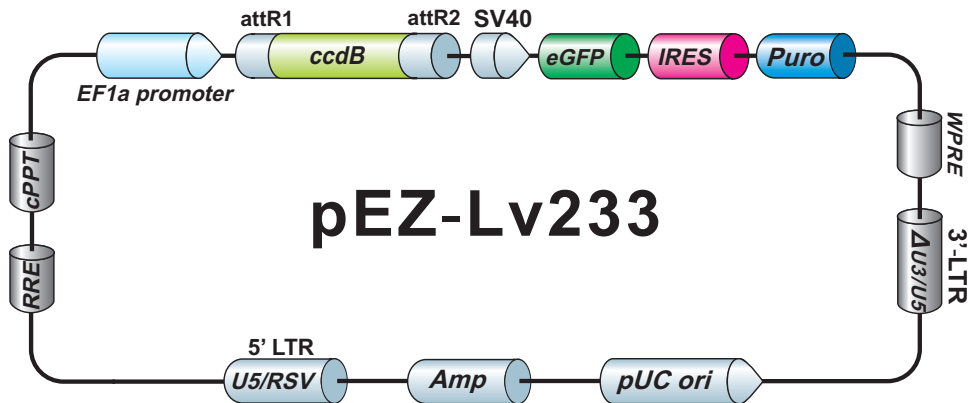


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv233

10215 bp, circular



Vector Features

Name	Position
SV40 promoter	7 374
eGFP	404 1128
IRES	1142 1718
Puromycin resistance gene	1751 2350
WPRE	2364 2955
3'-LTR(delta U3/U5)	3158 3394
pUC Ori (complementary strand)	3853 4523
Ampicillin resistance gene (complementary strand)	4668 5528
5'-LTR(RSV/U5)	5943 6354
HIV-1 RRE	7025 7258
cPPT	7750 7873
EF1a promoter	7970 8515
attR1	8523 8647
CmR	8756 9415
ccdB	9745 10050
attR2	10091 10214

Sequencing Primers

Primer	Sequence
Forward	5'- gtttcgtttctgttctgc -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