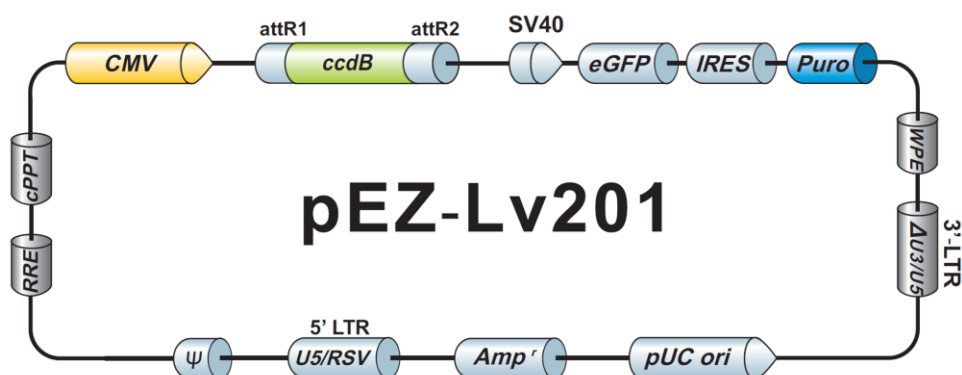


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv201

10835 bp, circular



Vector Features

Feature	Position
Sv40 promoter	7 374
eGFP	409 1128
IRES	1142 1718
Puromycin resistance gene	1751 2350
WPE	2364 2955
HIV 3' LTR	3158 3394
pUC Ori	3853 4523
Ampicillin resistance ORF	4668 5528
HIV 5' LTR	5943 6354
RNA packaging signal (Psi) sequence	6463 6507
RRE	7025 7258
cPPT	7750 7873
CMV promoter	7970 8564
attR1	8570 8694
ccdB	9123 9428
CmR	9770 10429
attR2	10710-10834

Sequencing Primers

Primer	Sequence
Forward	5'-GCGGTAGGCGTGTACGGT-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