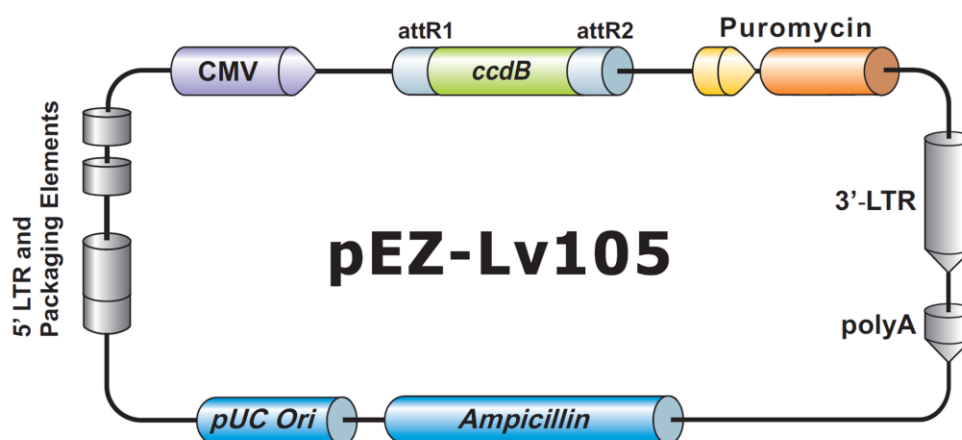


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv105

9661 bp, circular



Vector Features

Feature	Position
cPPT	78-197
hpGK promoter	244-750
Puromycin resistance gene	772-1371
3' self inactivating long terminal repeat (SIN/ 3;)	1498-1742
SV40 PolyA	1804-1934
Ampicillin resistance gene	2893-3753
pUC Ori	3954-4494
RSV promoter	5057-5285
HIV 5' LTR	5286-5466
RNA packaging signal (Psi) sequence	5577-5621
5' gag (delta 3rdG)	5624-5974
RRE	6141-6345
CMV promoter	7314-7940
attR1	7956-8080
CmR	8189-8848
ccdB	9190-9495
attR2	9536-9660

Sequencing Primers

Primer	Sequence
Forward	5'-GCGGTAGGCGTGTACGGT-3'
Reverse	5'-ATTGTGGATGAATACTGCC-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