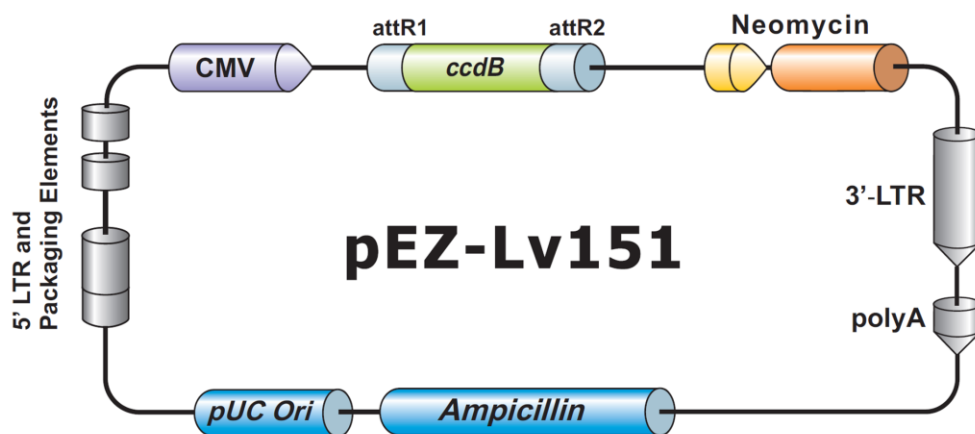


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv151

10437 bp, circular



Vector Features

Feature	Position
attR1	1-125
ccdB	554-859
CmR	1201-1860
attR2	2141-2265
cPPT	2344-2359
hpGK promoter	2510-3016
Neomycin resistance gene	3038-3841
3' self inactivating long terminal repeat (SIN/ 3')	3979-4223
SV40 PolyA	4285-4415
Ampicillin resistance ORF	5374-6234
pUC Ori	6435-6975
RSV promoter	7538-7766
HIV 5' LTR	7767-7947
RNA packaging signal (Psi) sequence	8058-8102
5' gag (delta 3rdG)	8105-8455
RRE	8622-8826
CMV promoter	9776-10422

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