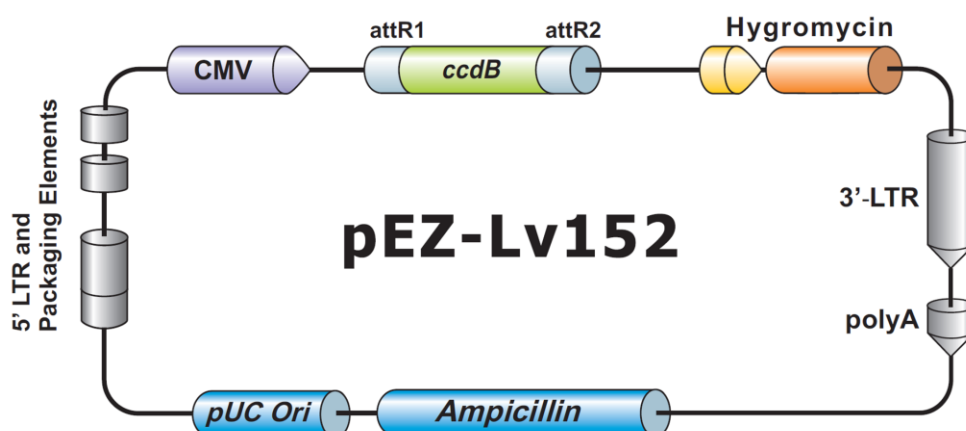


EZShuttle™ Recombination Cloning Expression Vector pEZ-Lv152

10596 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
Hygromycin resistance gene	3034-4059
3' self inactivating long terminal repeat (SIN/ 3')	4138-4382
SV40 PolyA	4444-4574
Ampicillin resistance ORF	5533-6393
pUC Ori	6594-7134
RSV promoter	7697-7925
HIV 5' LTR	7926-8106
RNA packaging signal (Psi) sequence	8217-8261
5' gag (delta 3rdG)	8264-8614
RRE	8781-8985
CMV promoter	9935-10581

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