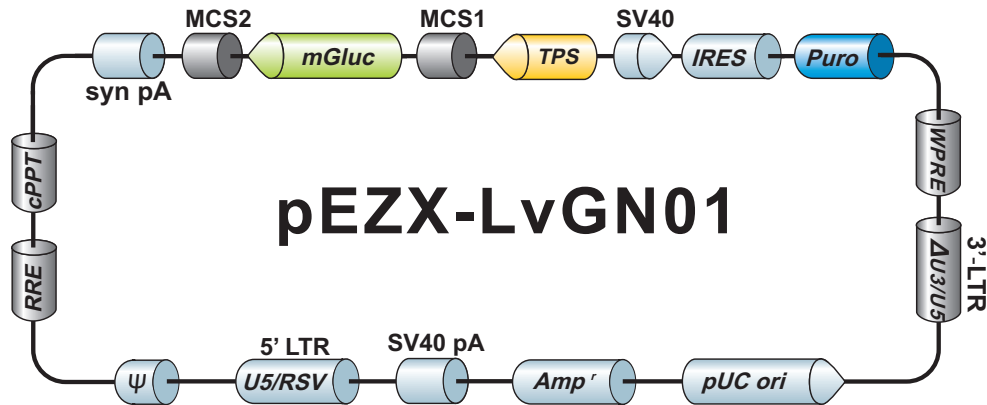


## pEZX-LvGN01

8137bp, circular



**MCS1:** Afel, AgeI, PmeI, EcoRI, MfeI, BstBI  
**MCS2:** XhoI, PspXI

### Vector Features

Name	Position
SV40 Promoter	1 368
IRES	399 972
Puromycin resistance gene	1005 1604
WPRE	1618 2206
3'LTR	2413 2646
pUC ori(complementary strand)	3166 3751
Ampicillin resistance gene(complementary strand)	3922 4782
SV40 polyA	4935 5069
5'LTR	5197 5608
HIV-1 psi pack sequence	5653 5778
RRE	6279 6512
cPPT	7004 7127
Synthesis polyA	7251 7304
MCS2	7305 7310
mGluc(complementary strand)	7311 7868
MCS1	7869 7906
TPS(complementary strand)	7907 8131

### Sequencing Primers

Primer	Sequence
Forward	5'-ttgttctcgggtgggcttggc-3'
Reverse	5'-caccctaactgacacacattccac-3'

GeneCopoeia, Inc.  
9620 Medical Center Drive, #101  
Rockville, MD 20850  
Tel: 301-762-0888 Fax: 301-762-3888  
Email: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com) Web: [www.genecopoeia.com](http://www.genecopoeia.com)

GeneCopoeia products are for research use only

© 2016, GeneCopoeia, Inc