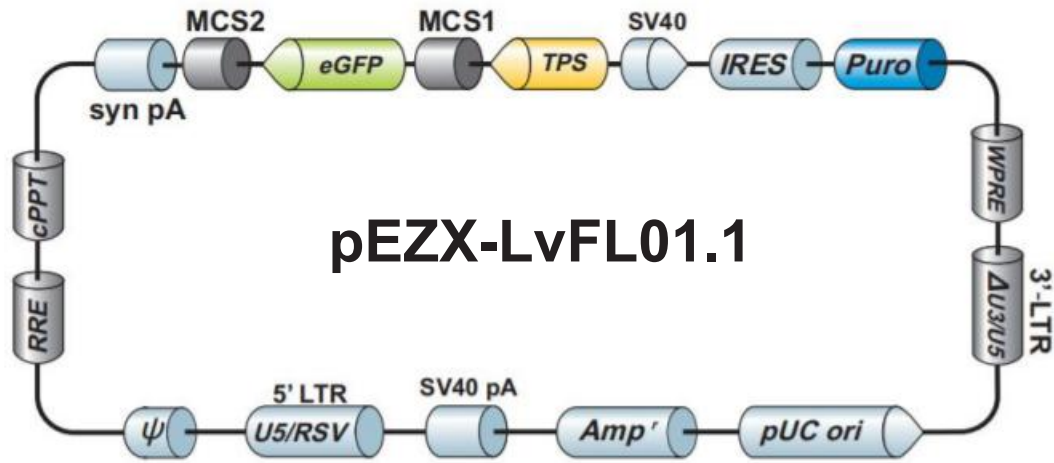


pEZX-LvFL01.1

8308 bp, circular



MCS1: XhoI

MCS2: AfeI, AgeI, PmeI, EcoRI, MfeI, BstBI

Vector Features

Name	Position
SV40 Promoter	10-339
IRES	399-972
Puromycin resistance gene	1005-1604
WPRE	1618-2206
3'LTR	2416-2649
pUC ori(complementary strand)	3169-3754
Ampicillin resistance gene(complementary strand)	3925-4785
SV 40 polyA	4938-5072
5'LTR	5429-5609
HIV-1 psi pack sequence	5656-5781
RRE	6282-6515
cPPT	7007-7122
Synthesis polyA	7254-7307
MCS2	7308-7313
eGFP(complementary strand)	7314-8033
MCS1	8039-8077
TPS(complementary strand)	8078-8302

Sequencing Primers

Primer	Sequence
Forward	5'- taatacgactcactataggg -3'
Reverse	5'- gtcattctattctggggg -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

© 2022, GeneCopoeia, Inc