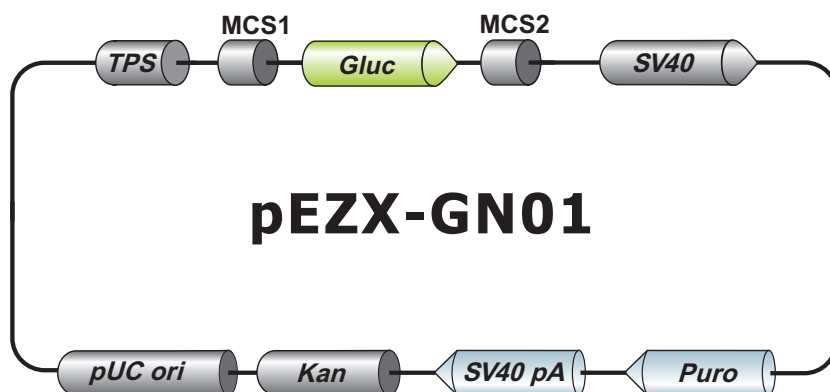


pEZX-GN01

4549bp, circular



MCS1: BstBI, BglII, EcoRI, EcoRV, HindIII
MCS2: NotI, XbaI

Vector Features

Name	Position
GLuc	7 564
MCS2	565 589
SV40 promoter	816 1145
Puromycin resistance gene	1207 1806
SV40 polyA	1906 2027
Kanamycin resistance gene	2239 3048
pUC ori	3138 3726
TPS	4249 4473
MCS1	4474 4549

Sequencing Primers

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

© 2016, GeneCopoeia, Inc