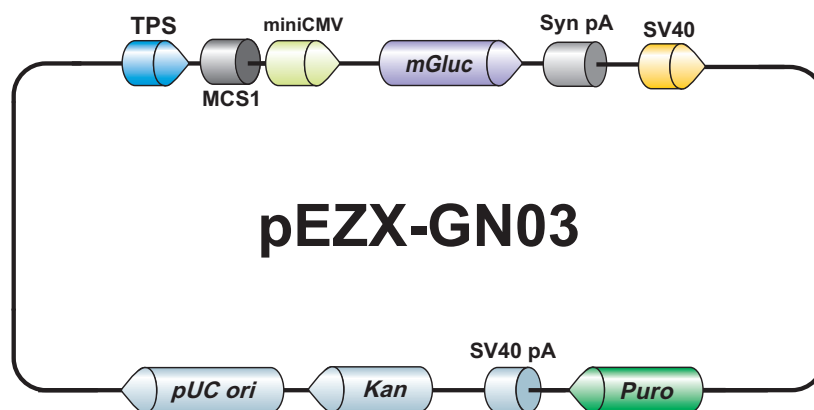


## pEZX-GN03

4367bp, circular



**MCS1:** SpeI, MluI, SmaI, HindIII

### Vector Features

Name	Position
TPS	489 649
MCS1	650 675
miniCMV promoter	676 707
mGluc	753 1310
Synthesis polyA	1324 1377
SV40 promoter	1395 1724
Puromycin resistance gene	1786 2385
SV40 polyA	2476 2697
Kanamycin resistance gene	2818 3627
pUC ori	3691 4364

### Sequencing Primers

Primer	Sequence
Forward	5'-agttacttaagctcgggcc-3'
Reverse	5'-ttgttctcgggtgggcttggc-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)