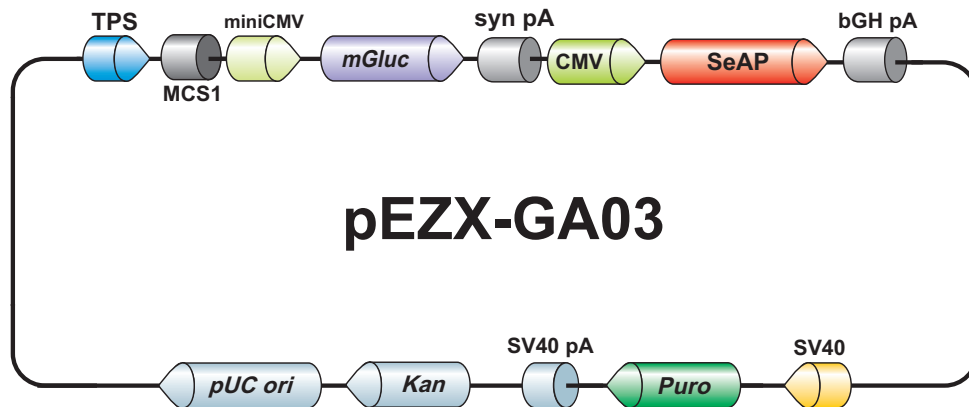


pEZX-GA03

6889bp, 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
CMV promoter	1386 1979
SeAP	2004 3545
bGH polyA	3670 3894
SV40 promoter	3917 4246
Puromycin resistance gene	4308 4907
SV40 polyA	4998 5219
Kanamycin resistance gene	5340 6149
pUC ori	6213 6886

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 Web: www.genecopoeia.com

GeneCopoeia products are for research use only

© 2016, GeneCopoeia, Inc