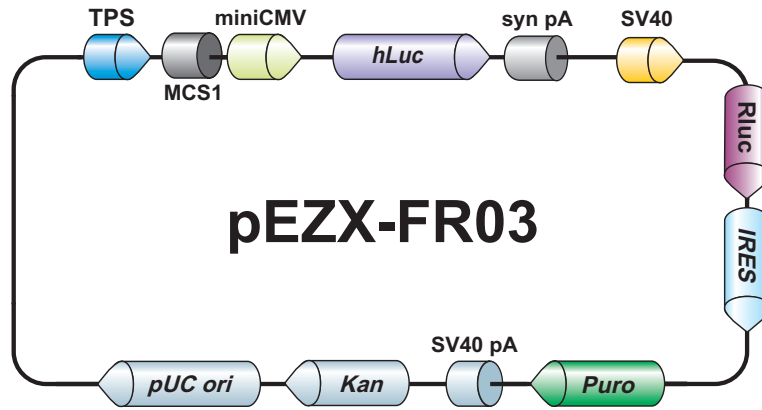


## pEZX-FR03

7026bp, circular



**MCS1**: SpeI, MluI, SmaI, Sall, PmeI, XbaI

### Vector Features

Name	Position
TPS	489 649
MCS1	650 689
miniCMV promoter	690 721
hLuc	766 2418
Synthesis polyA	2425 2478
SV40 promoter	2496 2825
Rluc	2888 3823
IRES	3839 4412
Puromycin resistance gene	4445 5044
SV40 polyA	5135 5356
Kanamycin resistance gene	5477 6286
pUC ori	6350 7023

### Sequencing Primers

Primer	Sequence
Forward	5'-agttacttaagctcgggcc-3'
Reverse	5'-atgtgggcatcgggaag-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