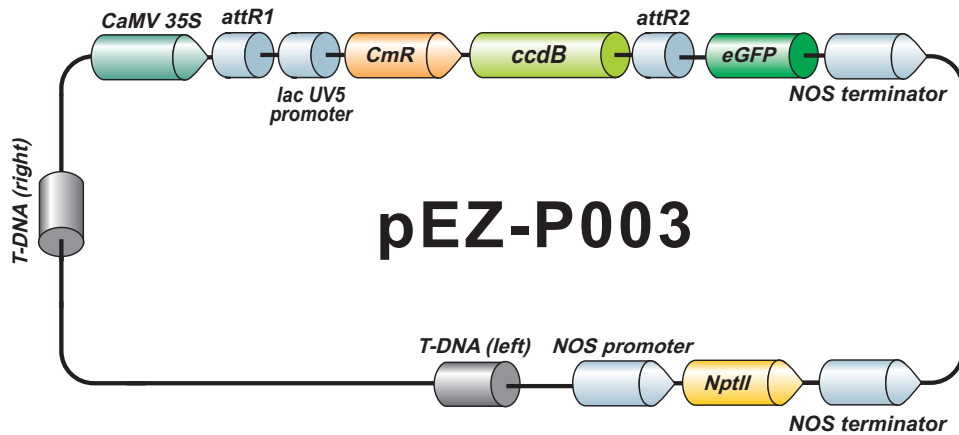


## EZShuttle™ Recombination Cloning Expression Vector pEZ-P003

12414 bp, circular



### Vector Features

Name	Position
T-DNA right border	54 78
CaMV 35S promoter	325 1159
attR1	1179 1303
lac UV5 promoter	1340 1370
CmR	1424 2083
ccdB	2425 2730
attR2	2771 2895
eGFP	2912 3631
NOS terminator	3640 4049
NOS terminator (complementary strand)	4083 4342
NptII (complementary strand)	4732 5526
NOS promoter (complementary strand)	5539 5845
T-DNA left border	5936 5960

### Sequencing Primers

Primer	Sequence
Forward	5'-CGTAAGGGATGACGCACAATC -3'
Reverse	5'-GAACAGCTCCTCGCCCTTGCTCAC -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

© 2015, GeneCopoeia, Inc