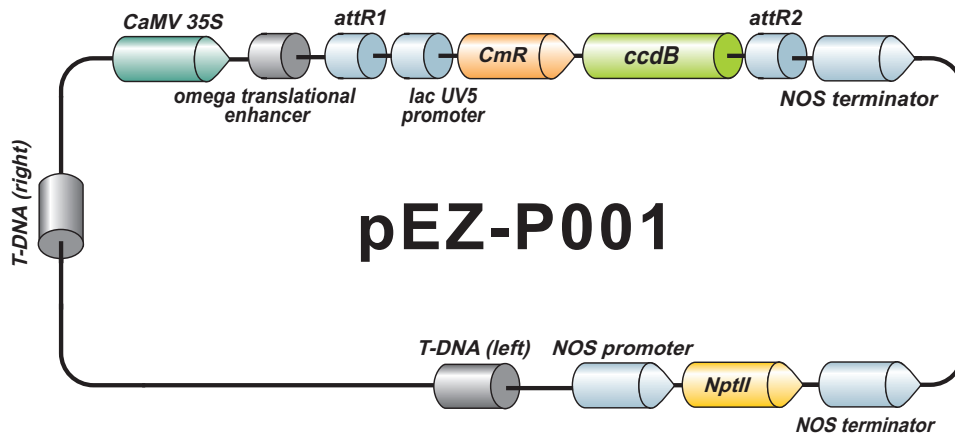


EZShuttle™ Recombination Cloning Expression Vector pEZ-P001

11678 bp, circular



Vector Features

Name	Position
T-DNA right border	54 78
CaMV 35S promoter	330 1079
omega translational enhancer	1091 1157
attR1	1170 1294
lac UV5 promoter	1333 1363
CmR	1417 2076
ccdB	2418 2723
attR2	2764 2888
NOS terminator	2904 3313
NOS terminator (complementary strand)	3342 3606
NptII (complementary strand)	3996 4790
NOS promoter (complementary strand)	4803 5109
T-DNA left border	5200 5224

Sequencing Primers

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

© 2015, GeneCopoeia, Inc