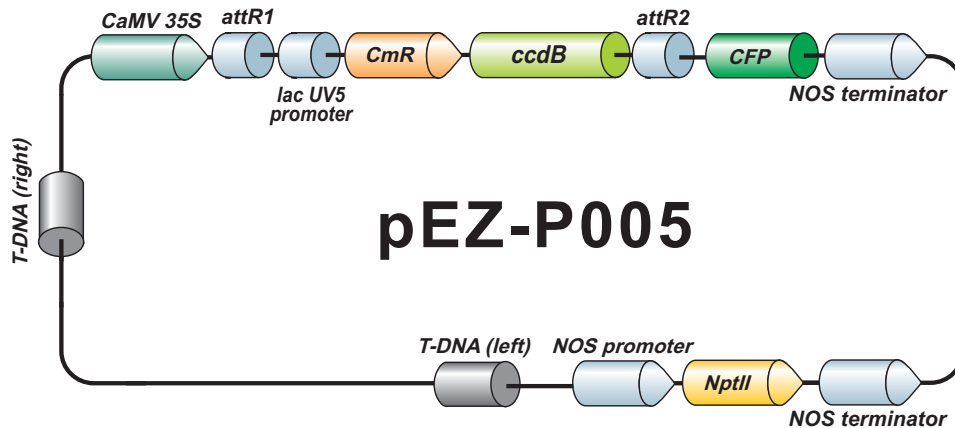


EZShuttle™ Recombination Cloning Expression Vector pEZ-P005

12402 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
CFP	2906 3622
NOS terminator	3628 4037
NOS terminator (complementary strand)	4071 4330
NptII (complementary strand)	4720 5514
NOS promoter (complementary strand)	5527 5833
T-DNA left border	5924 5948

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 Web: www.genecopoeia.com

GeneCopoeia products are for research use only

© 2015, GeneCopoeia, Inc