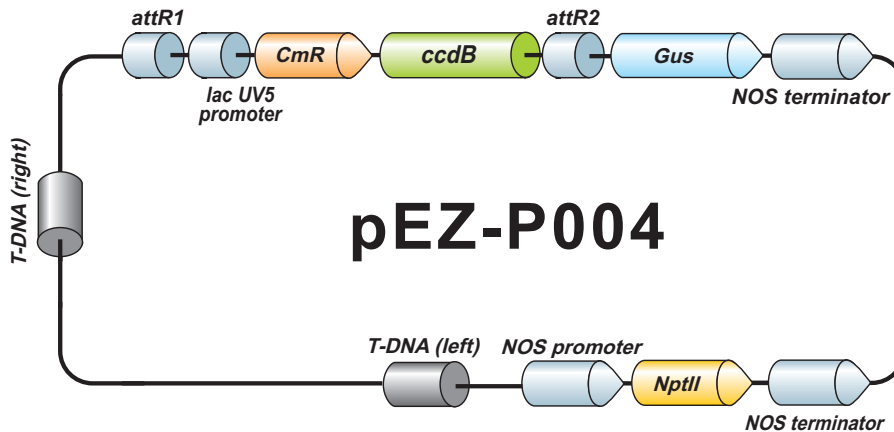


## EZShuttle™ Recombination Cloning Expression Vector pEZ-P004

12656 bp, circular



### Vector Features

Name	Position
T-DNA right border	54 78
attR1	338 462
lac UV5 promoter	499 529
CmR	583 1242
ccdB	1584 1889
attR2	1930 2054
Gus	2065 3876
NOS terminator	3882 4291
NOS terminator (complementary strand)	4328 4584
NptII (complementary strand)	4974 5768
NOS promoter (complementary strand)	5781 6087
T-DNA left border	6178 6202

### Sequencing Primers

Primer	Sequence
Forward	5'-CAGGGTTTTCCCAGTCACGAC -3'
Reverse	5'-GATCCAGACTGAATGCCAC -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