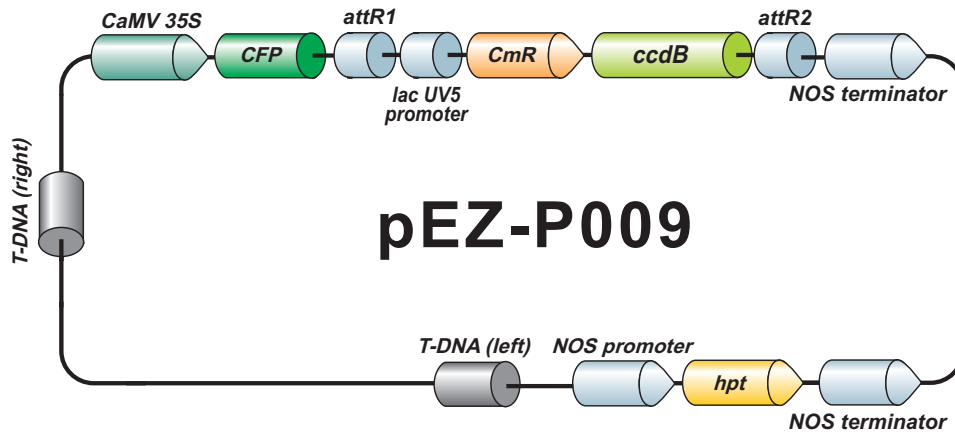


## EZShuttle™ Recombination Cloning Expression Vector pEZ-P009

12741 bp, circular



### Vector Features

Name	Position
T-DNA right border	54 78
CaMV 35S promoter	325 1159
CFP	1178 1894
attR1	1904 2028
lac UV5 promoter	2081 2111
CmR	2165 2824
ccdB	3166 3471
attR2	3521 3636
NOS terminator	3646 4055
NOS terminator (complementary strand)	4100 4365
hpt (complementary strand)	4501 5526
NOS promoter (complementary strand)	5554 5840
T-DNA left border	6263 6287

### Sequencing Primers

Primer	Sequence
Forward	5'-CCGACAACCACTACCTGA -3'
Reverse	5'-GATCGGGGAAATTCGAGC -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