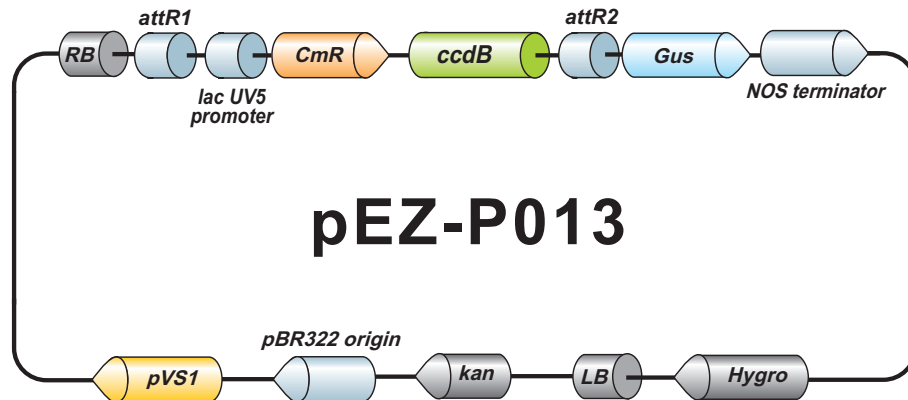


## EZShuttle™ Recombination Cloning Expression Vector pEZ-P013

12834bp, circular



### Vector Features

Name	Position
hygromycin resistance	1061 2086
LB	2377 2402
kanamycin resistance	2827 3621
pBR322 origin	3912 4192
pVS1	5002 6002
RB	8637 8662
attR1	8930 9054
lac UV5 promoter	9079 9109
CmR	9163 9822
ccdB	10164 10469
attR2	10510 10634
Gus	10693 12504
Nos terminator	12565 12822

### Sequencing Primers

Primer	Sequence
Forward	5'-TTTGATTTACGGGTTGGG -3'
Reverse	5'-CAGGGTTTTCCAGTCACGAC -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