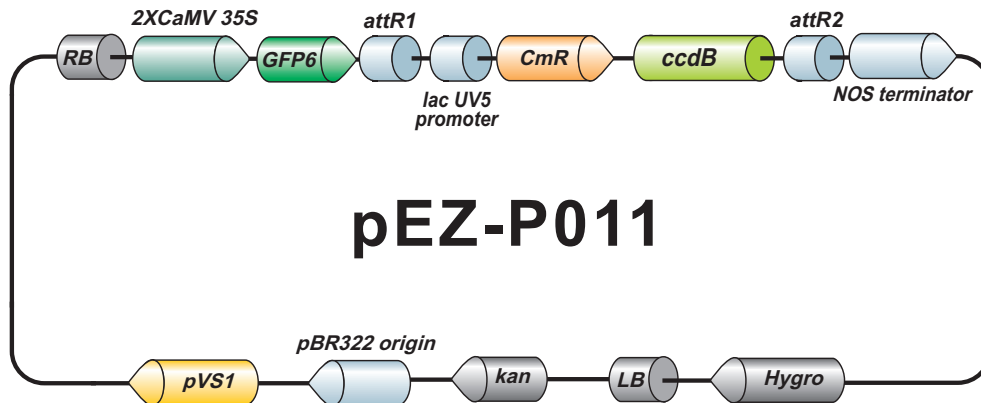


EZShuttle™ Recombination Cloning Expression Vector pEZ-P011

12460bp, circular



Vector Features

Name	Position
hygromycin resistance	1061 2086
LB	2377 2402
kanamycin resistance	2827 3621
pBR322 origin	3921 4192
pVS1	5002 6002
RB	8636 8661
2XCaMV 35S promoter	8921 9689
GFP6	9708 10429
attR1	10435 10564
lac UV5 promoter	10589 10619
CmR	10673 11332
ccdB	11674 11979
attR2	12020 12144
Nos terminator	12198 12455

Sequencing Primers

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