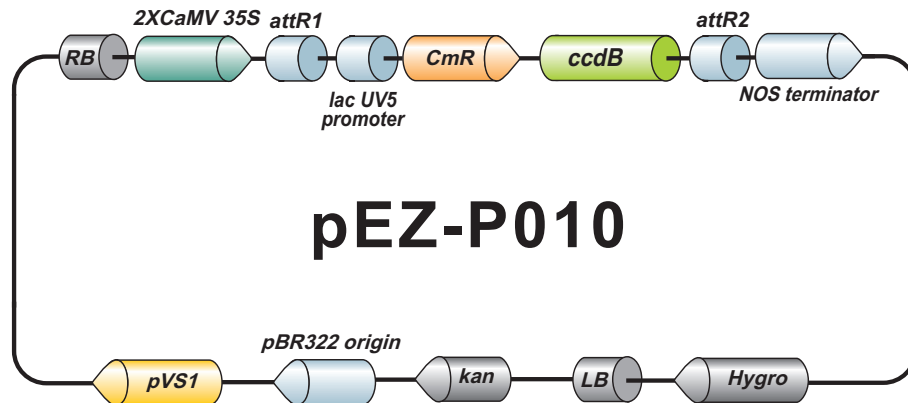


EZShuttle™ Recombination Cloning Expression Vector pEZ-P010

11752bp, circular



Vector Features

Name	Position
hygromycin resistance	1061 2086
LB	2377 2402
kanamycin resistance	2827 3621
pBR322 origin	3912 4192
pVS1	5002 6002
RB	8636 8661
2XCaMV 35S promoter	8921 9689
attR1	9725 9854
lac UV5 promoter	9879 9909
CmR	9963 10622
ccdB	10964 11269
attR2	11310 11434
Nos terminator	11490 11747

Sequencing Primers

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