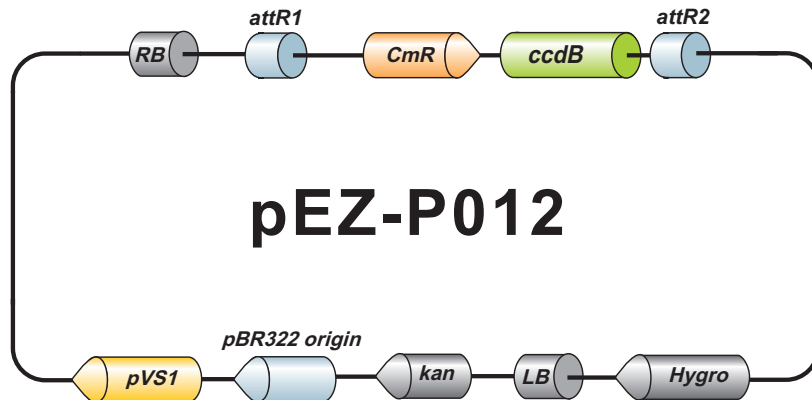


EZShuttle™ Recombination Cloning Expression Vector pEZ-P012

10742bp, circular



Vector Features

Name	Position
hygromycin resistance	1060 2085
LB	2376 2401
kanamycin resistance	2826 3620
pBR322 origin	3911 4191
pVS1	5001 6001
RB	8636 8661
attR1	8975 9104
CmR	9213 9872
ccdB	10214 10519
attR2	10560 10684

Sequencing Primers

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