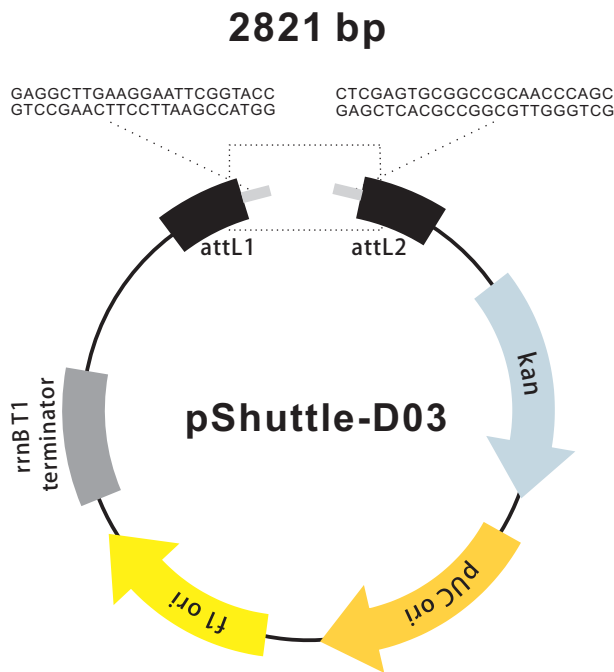


EZShuttle™ Recombination Cloning Shuttle cloning Vector pShuttle-D03



Note:Design gene specific primers

Forward Primer: 5'-^{Homologous Sequence}CAGGCTTGAAGGAATTCGGTACC^{specific sequence}→3'

Reverse Primer: 5'-^{Homologous Sequence}GCTTGGGTTGCGGCCGCACTCGAG^{specific sequence}→3'

Vector Features

Name	Position
attL2	17 116
Kanamycin resistance gene	286 1095
pUC ori	1242 1830
f1 ori	1897 2352
rrnB T1 terminator	2523 2609
attL1	2705 2804

Sequencing Primers

Primer	Sequence
Forward	5'- tgtaaacgacggccagt -3'
Reverse	5'- caggaaacagctatgacc -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