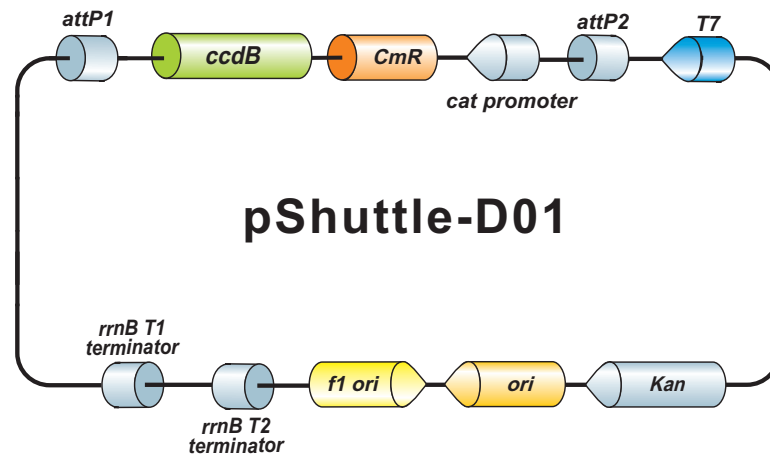


## EZShuttle™ Recombination Cloning Shuttle cloning Vector pShuttle-D01

5003 bp, circular



### Vector Features

Name	Position
f1 ori (complementary strand)	2 457
rrnB T2 terminator	513 540
rrnB T1 terminator	632 718
M13 Forward priming site	782 798
attP1	815 1046
ccdB (complementary strand)	1442 1747
CmR (complementary strand)	2089 2748
cat promoter (complementary strand)	2749 2851
attP2	2996 3227
T7 promoter (complementary strand)	3246 3264
M13 Reverse priming site	3269 3285
Kanamycin resistance gene	3398 4207
ori	4354 4942

### 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](mailto:inquiry@genecopoeia.com) Web: [www.genecopoeia.com](http://www.genecopoeia.com)

GeneCopoeia products are for research use only

© 2015, GeneCopoeia, Inc