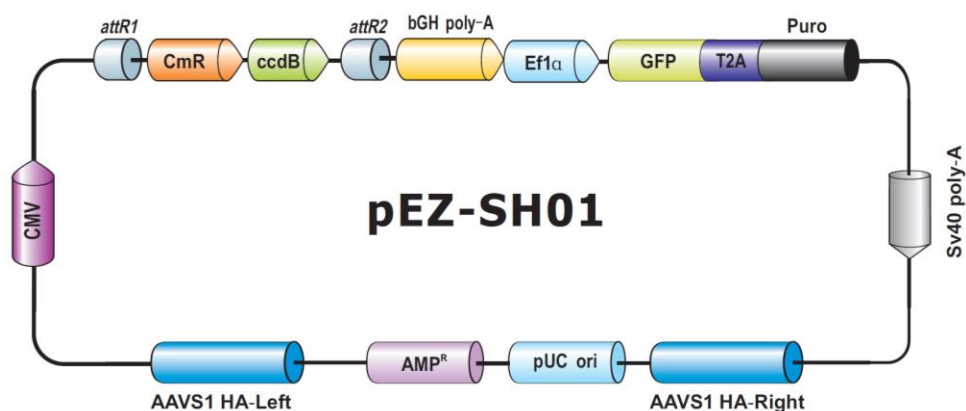


## EZShuttle™ Recombination Cloning Human AAVS1 Safe Harbor Knockin Vector pEZ-SH01

10442 bp, circular



### Vector Features

Feature	Position
AAVS1 Left Homology Arm	1402-2205
CMV promotor	2346-2549
attR1	2628-2752
CmR	2861-3520
ccdB	3862-4167
attR2	4208-4332
bGH PolyA	4360-4574
EF1alpha promoter	4589-5134
copGFP	5148-5903
T2A	5904-5957
Puromycin resistance gene	5624-6557
SV40 PolyA	6665-6786
AAVS1 Right Homology Arm	7048-7886
pUC Ori	8823-9411
Ampicillin resistance gene	10374-10442

## Sequencing Primers

Primer	Sequence
Forward	5'-CGGTGGGAGGTCTATATAAGCAG-3'
Reverse	5'-GACAGTGGGAGTGGCACCTTC-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