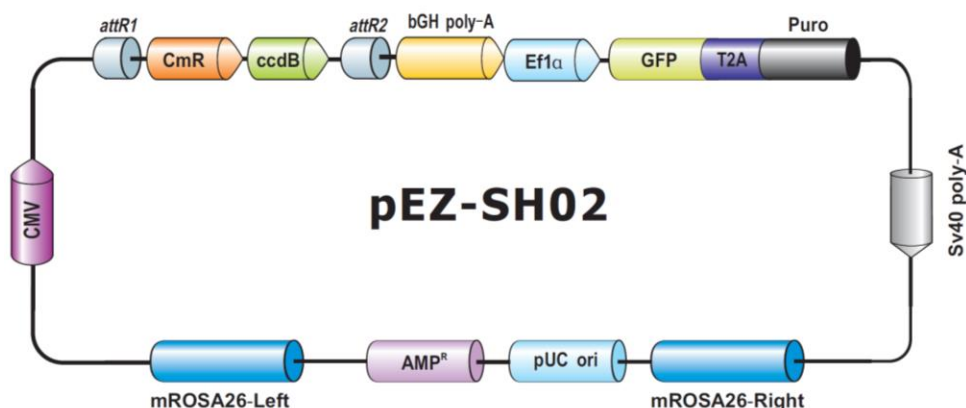


EZShuttle™ Recombination Cloning Mouse ROSA26 Safe Harbor Knockin Vector pEZ-SH02

10351 bp, circular



Vector Features

Feature	Position
ROSA26 Left Homology Arm	1365-2159
CMV promotor	2299-2502
attR1	2582-2706
CmR	2815-3474
ccdB	3804-4109
attR2	4150-4273
bGH PolyA	4287-4511
EF1alpha promoter	4526-5071
copGFP	5085-5840
T2A	5841-5894
Puromycin resistance gene	5895-6494
SV40 PolyA	6602-6723
ROSA26 Right Homology Arm	6985-7795
pUC Ori	8732-9320
Ampicillin resistance gene	9491-10351

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 Web: www.genecopoeia.com

GeneCopoeia products are for research use only

© 2015, GeneCopoeia, Inc