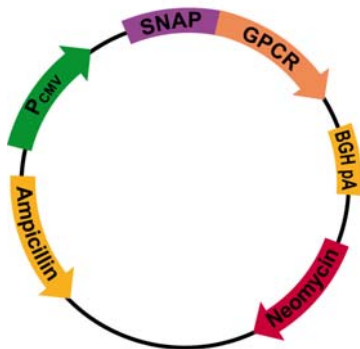


www.htrf.com

Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-C5A

Plasmid Tag-Lite™ Receptor

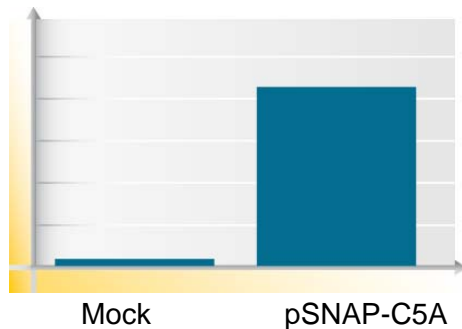
pSNAP-C5A
 Complement component 5
 receptor 1

Plasmid code
Storage temperature
Quantity

PSNAPC5A
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-C5A
Plasmid size	7094 bp
C5a ORF size	1065 bp
ORF sequence (based on NCBI nucleic sequence accession number)	NM_001736.3
Promoter	CMV
Host	Mammalian
Antibiotic resistance for E. Coli	Ampicillin
Mammalian selection marker	Neomycin

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-C5A	23

*The plasmid is transiently expressed into HEK 293. Cells are plated and labelled 24 hours after transfection with the Tag-lite SNAP-Lumi4™-Tb (SSNPTBE) according to the protocol described on the web site. Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

Sequence

Information available on: www.htrf.com

File under a txt format can be downloaded from Tag-lite section on the website.

MSDS

Information available on: www.htrf.com