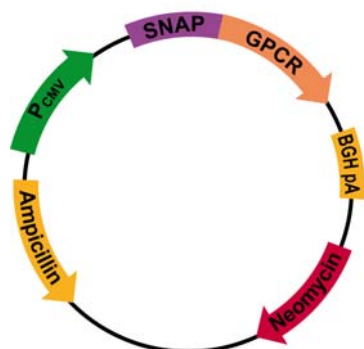


www.htrf.com

Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-V2

Plasmid Tag-Lite™ Receptor

pSNAP-V2
Arginine vasopressin receptor 2

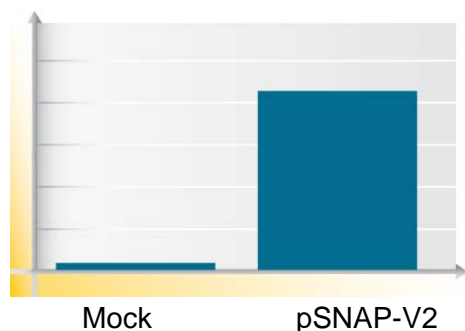
Plasmid code
Storage temperature
Quantity

PSNAPV2
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-V2
Plasmid size	7154 bp
V2 ORF size	1125 bp
ORF sequence (based on NCBI nucleic sequence accession number)	NM_000054.4

Promoter	CMV
Host	Mammalian
Antibiotic resistance for E. Coli	Ampicillin
Mammalian selection marker	Neomycin

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-V2	11

*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