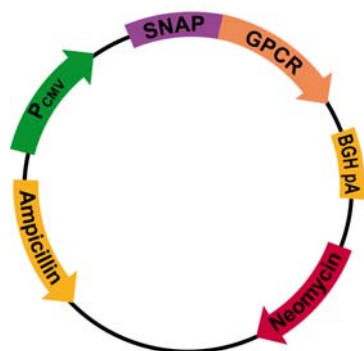


www.htrf.com

Plasmid Information

For in vitro research use only

Version June.-10-C



Tag-lite pSNAP-Alpha1A

Plasmid Tag-Lite™ Receptor

pSNAP-Alpha1A Adrenergic receptor alpha1A

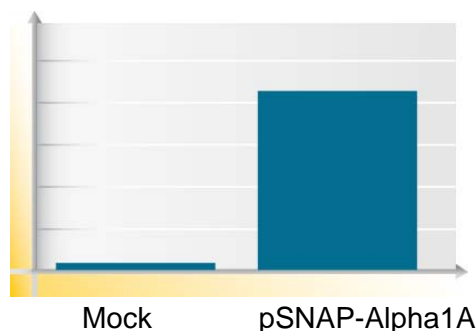
Plasmid code
Storage temperature
Quantity

PSNAPADR1
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-Alpha1A
Plasmid size	7457 bp
Alpha1A ORF size	1411 bp
ORF sequence (based on NCBI nucleic sequence accession number)	NM_000680.2

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

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-Alpha1A	37

*The plasmid is transiently expressed into HEK293 T. Cells are plated and labelled 24 hours after transfection with the Tag-lite SNAP-Lumi4™-Tb (SSNPTBC) 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