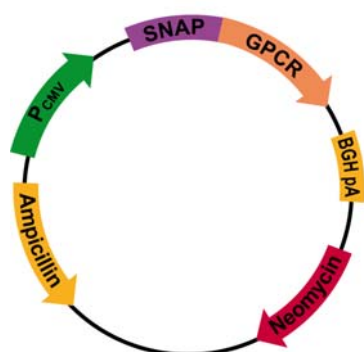


Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-NK2
Tachykinin

Plasmid Tag-Lite™ Receptor

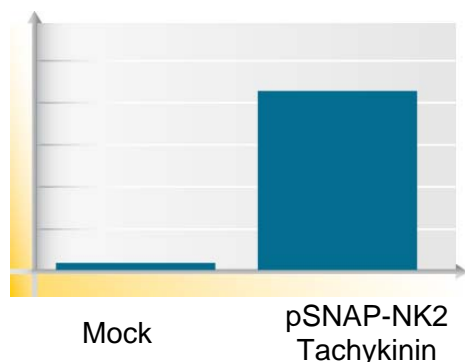
pSNAP-NK2 Tachykinin
Tachykinin/Substance K receptor

Plasmid code
Storage temperature
Quantity

PSNAPNK2
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-NK2 Tachykinin
Plasmid size	7229 bp
NK2 ORF size	1206 bp
ORF sequence (based on NCBI nucleic sequence accession number)	AY322545.1
Promoter	CMV
Host	Mammalian
Antibiotic resistance for E. Coli	Ampicillin
Mammalian selection marker	Neomycin

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-NK2 Tachykinin	92

*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