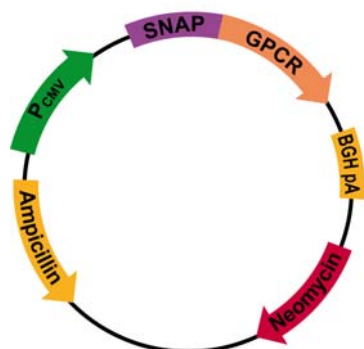


www.htrf.com

Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-CCK1

Plasmid Tag-Lite™ Receptor

pSNAP-CCK1
 Cholecystokinin receptor type A

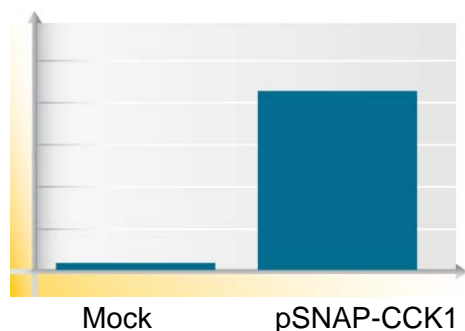
Plasmid code
Storage temperature
Quantity

PSNAPCCK1
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-CCK1
Plasmid size	7328 bp
CCK1 ORF size	1299 bp
ORF sequence (based on NCBI nucleic sequence accession number)	NM_000730.2

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

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-CCK1	10

*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