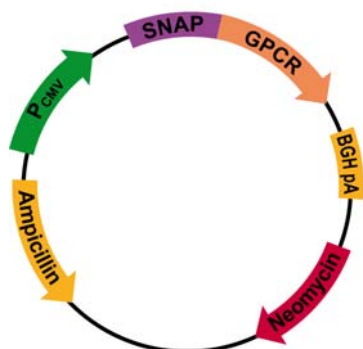


## Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-Kappa opioid

## Plasmid Tag-Lite™ Receptor

## pSNAP-Kappa opioid Kappa opioid receptor

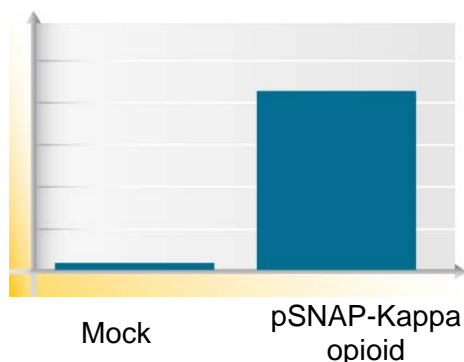
**Plasmid code**  
**Storage temperature**  
**Quantity**

PSNAPOPRK1  
 -20°C  
 10 µg/10 µL H<sub>2</sub>O

<b>Plasmid name</b>	pSNAP-Kappa opioid
<b>Plasmid size</b>	7181 bp
<b>Kappa opioid ORF size</b>	1152 bp
<b>ORF sequence</b> (based on NCBI nucleic sequence accession number)	NM_000912.3

<b>Promoter</b>	CMV
<b>Host</b>	Mammalian
<b>Antibiotic resistance for E. Coli</b>	Ampicillin
<b>Mammalian selection marker</b>	Neomycin

## Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-Kappa opioid	9

\*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](http://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](http://www.htrf.com)