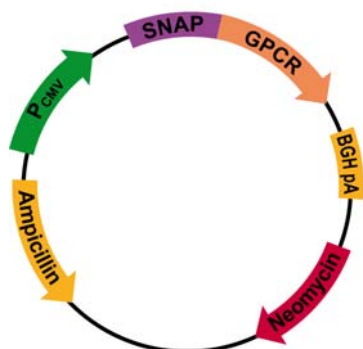


www.htrf.com

## Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-GIPR

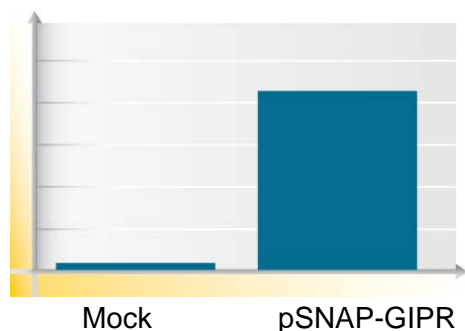
## Plasmid Tag-Lite™ Receptor

pSNAP-GIPR  
 Gastric inhibitory polypeptide  
 receptor

**Plasmid code** PSNAPGIPR  
**Storage temperature** -20°C  
**Quantity** 10 µg/10 µL H<sub>2</sub>O

<b>Plasmid name</b>	pSNAP-GIPR
<b>Plasmid size</b>	7364 bp
<b>GIPR ORF size</b>	1335 bp
<b>ORF sequence</b> (based on NCBI nucleic sequence accession number)	NM_000164.2
<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-GIPR	4

\*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)