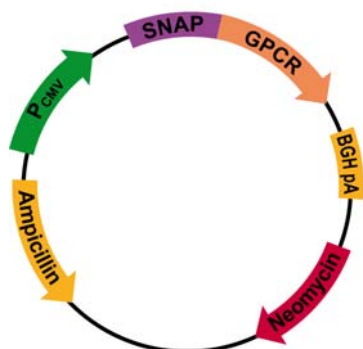


[www.htrf.com](http://www.htrf.com)

## Plasmid Information

For in vitro research use only

Version Feb.-10-A



Tag-lite pSNAP-Beta1

## Plasmid Tag-Lite™ Receptor

pSNAP-Beta1  
Adrenergic receptor beta1

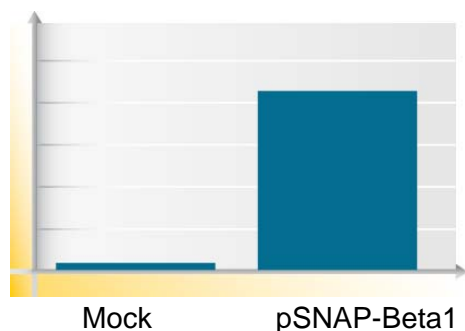
**Plasmid code**  
**Storage temperature**  
**Quantity**

PSNAPBETA1  
 -20°C  
 10 µg/10 µL H<sub>2</sub>O

<b>Plasmid name</b>	pSNAP-Beta1
<b>Plasmid size</b>	7412 bp
<b>Beta1 ORF size</b>	1389 bp
<b>ORF sequence</b> (based on NCBI nucleic sequence accession number)	NM_000684.2

<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-Beta1	8

\*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)