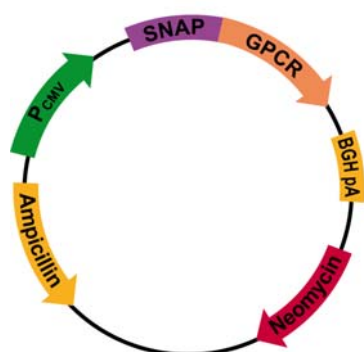


www.htrf.com

Plasmid Information

For in vitro research use only

Version June.-10-C



Tag-lite pSNAP-Beta2

Plasmid Tag-Lite™ Receptor

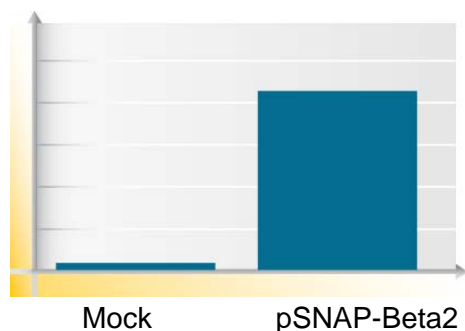
pSNAP-Beta2 Adrenergic receptor beta2

Plasmid code
Storage temperature
Quantity

PSNAPBD2
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-Beta2
Plasmid size	7292 bp
Beta2 ORF size	1253 bp
ORF sequence (based on NCBI nucleic sequence accession number)	NM_000024.4
Promoter	CMV
Host	Mammalian
Antibiotic resistance for E. Coli	Ampicillin
Mammalian selection marker	Neomycin

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-Beta2	5

*The plasmid is transiently expressed into HEK293 T. Cells are plated and labelled 24 hours after transfection with the Tag-lite SNAP-Lumi4™-Tb (SSNPTBC) 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