

R&D, Administration and Europe Office
 Cisbio Bioassays
 Phone: +33 (0)4 66 79 67 05
 Fax: +33 (0)4 66 79 19 20
 E-mail: bioassays@cisbio.com

Japan Office
 Sctei Medical Labo K.K.
 Phone: +81 (0)3 5510 2932
 Fax: +81 (0)3 5510 0130
 E-mail: reagent@sctimedilabo.co.jp

USA Office
 Cisbio US, Inc.
 Phone : 888 963 4567
 Fax : 781 687 1500
 E-mail : htrfinfo@cisbio.us

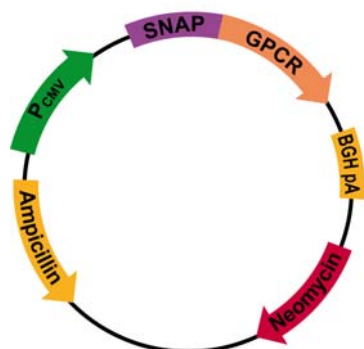


www.htrf.com

Plasmid Information

For in vitro research use only

Version Feb.-10-B



Tag-lite pSNAP-B2

Plasmid Tag-Lite™ Receptor

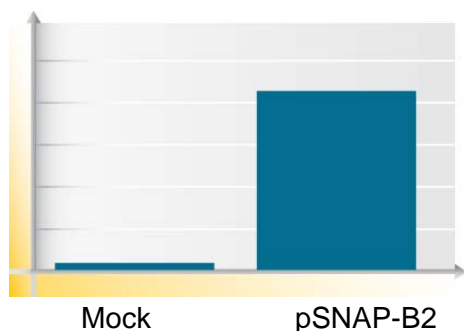
pSNAP-B2
 Bradykinin receptor

Plasmid code
Storage temperature
Quantity

PSNAPB2
 -20°C
 10 µg/10 µL H₂O

Plasmid name	pSNAP-B2
Plasmid size	7229 bp
B2 ORF size	1190 bp
ORF sequence (based on NCBI nucleic sequence accession number)	NM_000623.3
Promoter	CMV
Host	Mammalian
Antibiotic resistance for E. Coli	Ampicillin
Mammalian selection marker	Neomycin

Quality Control



Plasmid	Normalized 620 nm signal*
pSNAP-B2	39

*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