



# AdrenoceptorPLASMID TAG-LITE® pSNAP-Beta1

## Receptor: Adrenergic receptor beta1

Product Information Revision: #02 of September 2023

Part#: PSNAPBETA1

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H2O

Plasmid name: pSNAP-Beta1

Plasmid size: 7412 bp

beta1-ORF size: 2049 bp

ORF sequence\*: NM\_000684.2 (-PS)

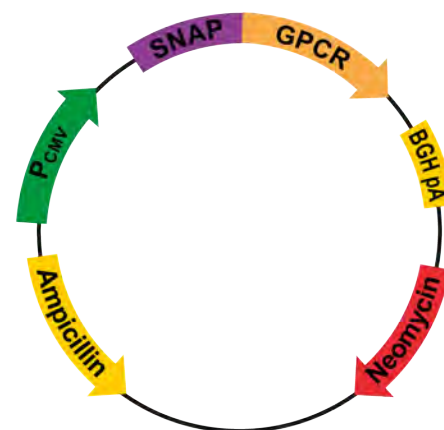
\*based on NCBI nucleic sequence accession number

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

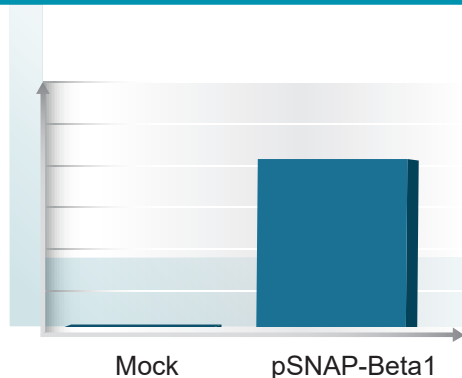
Mammalian selection marker: Neomycin



Tag-lite pSNAP-Beta1

For research use only. Not for use in diagnostic procedures.

### 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 (SSNPTBC) according to the manual described on the web site.

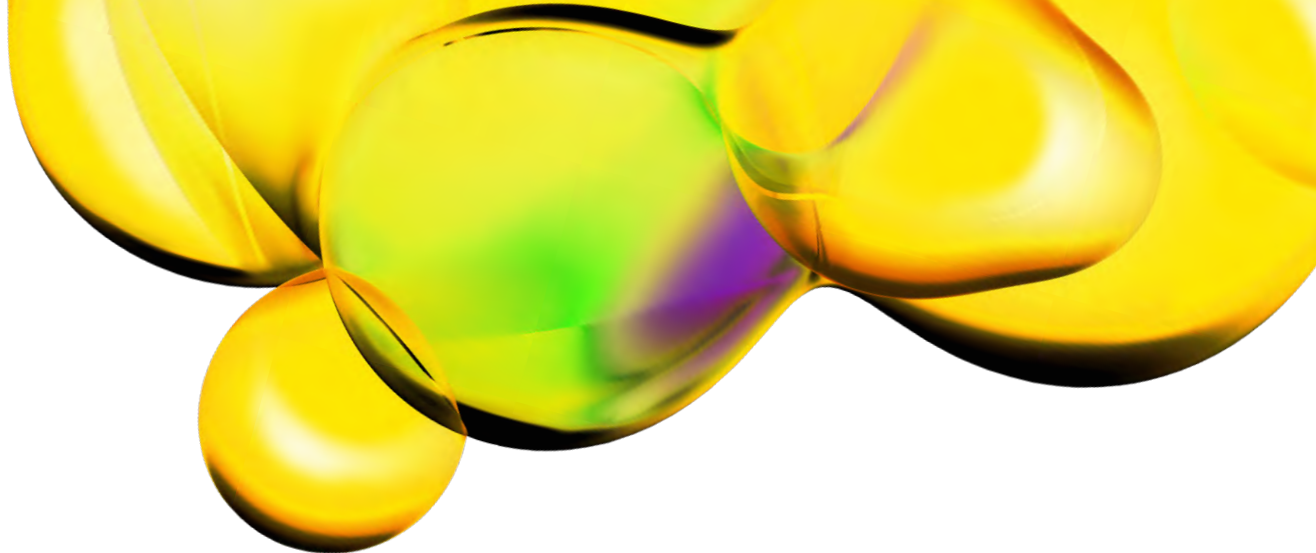
Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

## SEQUENCE & MSDS

MSDS can be downloaded from Tag-lite section on the website [www.revvy.com](http://www.revvy.com).

**Sequence:** F-Snap-N-beta1ADR GC (923bp - 2971bp, direct) 2049bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCCACCTGG
ATAGCCCTCTGGGCAAGCTGGAAGTGTCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCCGAGCCGCGTGTGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGCT
CAACGCCTACTTTACCCAGCCTGAGGCCATCGAGGAGTTCCCTGTGCCAGCCCTGCACCACCCAGTGTTCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGGAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTCATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCGCCGTGAAAACCGCCCTGAGCGGAAATCCCGTGCCCATTCCTGATCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGCTGCTGGCCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCGCACCCGCTCCCCGACGGCGCGGCCACCGCGGGCGGGCTGCTGGT
CCCGCGTCGCCGCCCGCCTCGTTGCTGCCTCCCGCCAGCGAAAGCCCCGAGCCGCTGTCTCAGCAGTGGACAGCGGGCAT
GGGTCTGCTGATGGCGCTCATCGTGTGCTCATCGTGGCGGGCAATGTGCTGGTGATCGTGGCCATCGCCAAGACGCCGC
GGCTGCAGACGCTACCAACCTCTTCATCATGTCCCTGGCCAGCGCCGACCTGGTCATGGGGCTGCTGGTGGTGCCGTTT
GGGGCCACCATCGTGGTGTGGGGCCGCTGGGAGTACGGCTCCTTCTTCTGCGAGCTGTGGACCTCAGTGGACGTGCTGTG
CGTGACGGCCAGCATCGAGACCCTGTGTGTATTGCCCTGGACCGCTACCTCGCCATCACCTCGCCCTTCCGCTACCAGA
GCCTGCTACGCGCGCGCGGGCGGGGCTCGTGTGCACCGTGTGGCCATCTCGGCCCTGGTGTCTTCCCTGCCATC
TCATGCACTGGTGGCGGGCGGAGAGCGACGAGGCGCGCCGCTGCTACAACGACCCCAAGTGCTGCGACTTCGTCACCAA
CCGGGCCTACGCCATCGCCTCGTCCGTAGTCTCCTTCTACGTGCCCTGTGCATCATGGCCTTCGTGTACCTGCGGGTGT
TCCGCGAGGCCAGAAAGCAGGTGAAGAAGATCGACAGCTGCGAGCGCCGTTTCCCTCGGCGGCCAGCGCGGCCGCCCTCG
CCCTCGCCCTCGCCCGTCCCGCGCCCGCGCCGCCCGCCGACCCCGCGCCCGCCGCCCGCCGCCACCGCCCCGCT
GGCCAACGGGCGTGCGGTAAGCGGGCGGCCCTCGCGCCTCGTGGCCCTGCGCGAGCAGAAGGCCTCAAGACGCTGGGCA
TCATCATGGGCGTCTTACGCTCTGCTGGCTGCCCTTCTTCTGGCCAACGTGGTGAAGGCCTTCCACCGCGAGCTGGTG
CCCGACCGCCTCTTCGTCTTCTTCAACTGGCTGGGCTACGCCAACTCGGCCTTCAACCCCATCATCTACTGCCGAGCCC
CGACTTCCGCAAGGCCTTCCAGGGACTGCTCTGCTGCGCGCGCAGGGCTGCCCGCCGGCGCCACGCGACCCACGGAGACC
GGCCGCGCGCCTCGGGCTGTCTGGCCCGGCCCGGACCCCGCCATCGCCGGGGCCGCTCGGACGACGACGACGACGAT
GTCGTCGGGGCCACGCCGCCCGCGCCTGCTGGAGCCCTGGGCGGGTGC AACGCGGGGCGGGCGGGACAGCGACTC
GAGCCTGGACGAGCCGTGCCGCCCGGCTTCGCTCGGAATCCAAGGTG
```



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

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

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.