

Adenosine PLASMID TAG-LITE® pSNAP-A1 Receptor: Adenosine receptor A1

Product Information Revision: #03 of September 2023

Part#: PSNAPA1

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H₂O

Plasmid name: pSNAP-A1

Plasmid size: 7019 bp

A1-ORF size: 1650 bp

ORF sequence*: NM_000674.2

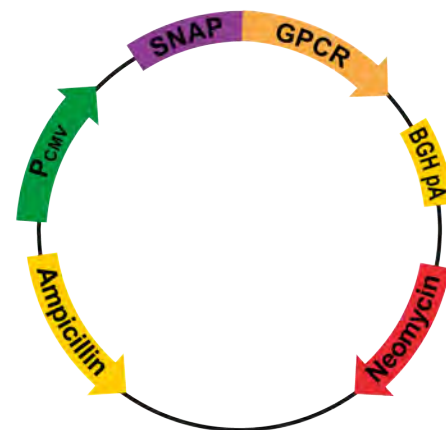
***based on NCBI nucleic sequence accession number**

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

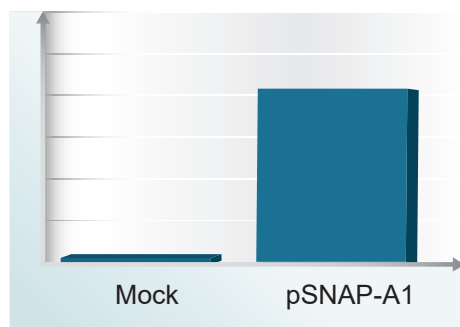
Mammalian selection marker: Neomycin



Tag-lite pSNAP-A1

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

QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pSNAP-A1	7

*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.

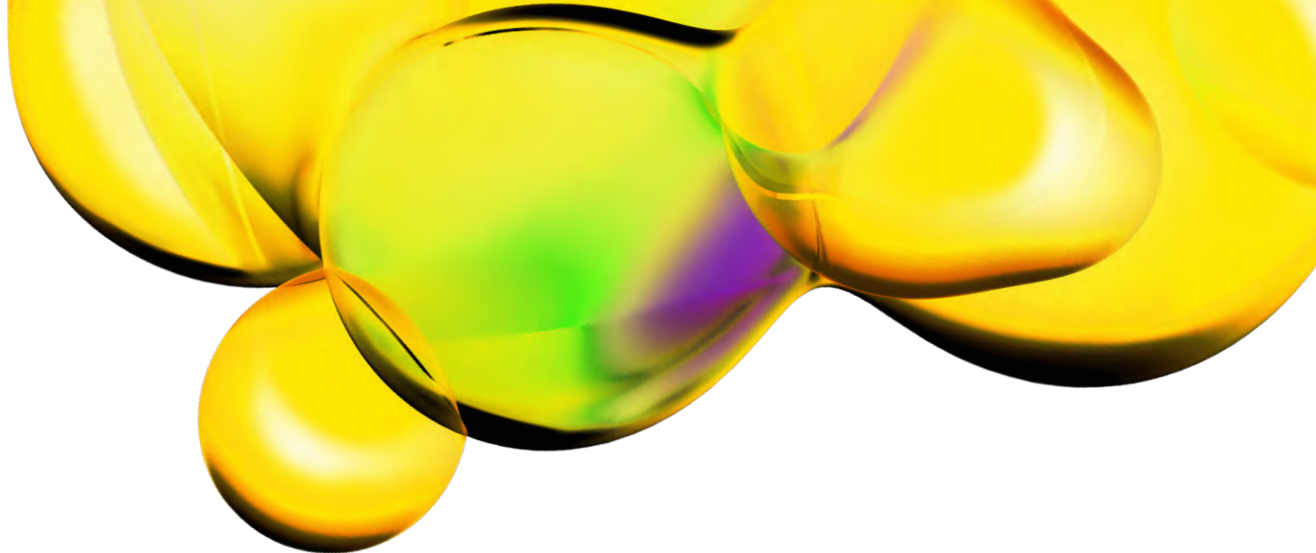
Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

SEQUENCE & MSDS

MSDS can be downloaded from Tag-lite section on the website www.revvy.com.

Sequence: N° 18 F-Snap-N-A1 GC (923bp - 2572bp, direct) 1650bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCCACCTGG
ATAGCCCTCTGGGCAAGCTGGAAGTGTCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCCCAGCCGCCGTGCTGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGT
CAACGCCTACTTTACCAGCCTGAGGCCATCGAGGAGTTCCCTGTGCCAGCCCTGCACCACCCAGTGTTCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGAAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTCATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCGCCGTGAAAACCGCCCTGAGCGGAAATCCCGTGCCCATCTGATCCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGCTGCTGGCCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCCCGCCCTCCATCTCAGCTTTCCAGGCCGCCTACATCGGCATCGAG
GTGCTCATCGCCCTGGTCTCTGTGCCCGGAACGTGCTGGTGTCTGGGCGGTGAAGGTGAACCAGGCCGCTGCGGGATGC
CACCTTCTGCTTTCATCGTGTGCTGGCGGTGGCTGATGTGGCCGTGGGTGCCCTGGTTCATCCCCCTCGCCATCCTCATCA
ACATTGGGCCACAGACCTACTTCCACACCTGCCTCATGGTTGCCCTGTCCGGTCCCTCATCCTCACCAGAGCTCCATCCTG
GCCCTGCTGGCAATTGCGGTGGACCGCTACCTCCGGGTCAAGATCCCTCTCCGGTACAAGATGGTGGTGACCCCCGGAG
GGCGGCGGTGGCCATAGCCGGCTGCTGGATCCTCTCCTTCTGGTGGGACTGACCCCTATGTTTGGCTGGAACAATCTGA
GTGCGGTGGAGCGGGCCTGGGAGCCAACGGCAGCATGGGGGAGCCCGTATCAAGTGCAGTTCGAGAAGGTCATCAGC
ATGGAGTACATGGTCTACTTCAACTTCTTTGTGTGGGTGCTGCCCCGCTTCTCCTCATGGTCCCTCATCTACCTGGAGGT
CTTCTACCTAATCCGCAAGCAGCTCAACAAGAAGGTGTGGCCTCCTCCGGCGACCCGCAGAAGTACTATGGGAAGGAGC
TGAAGATCGCCAAGTCGCTGGCCCTCATCCTCTTCTCTTTGCCCTCAGCTGGCTGCCTTTGCACATCCTCAACTGCATC
ACCCTTCTGCCCCGCTGCCACAAGCCCAGCATCCTTACCTACATTGCCATCTTCTCACGCACGGCAACTCGGCCAT
GAACCCATTGTCTATGCCTTCCGCATCCAGAAGTCCGCGTCACCTTCTTAAGATTGGAATGACCATTCCGCTGCC
AGCCTGCACCTCCATTGACGAGGATCTCCAGAAGAGAGGCCTGATGAC
```



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

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.