



# Adenosine PLASMID TAG-LITE® pSNAP-A3 Receptor: Adenosine receptor A3

PRODUCT INFORMATION - rev. 2 (Sept. 2018)

**Part#:** PSNAPA3

**Storage temperature:** -20°C

**Quantity:** 10 µg/ 10 µL H2O

**Plasmid name:** pSNAP-A3

**Plasmid size:** 6998 bp

**A3-ORF size:** 1629 bp

**ORF sequence\*:** NM\_000677.3

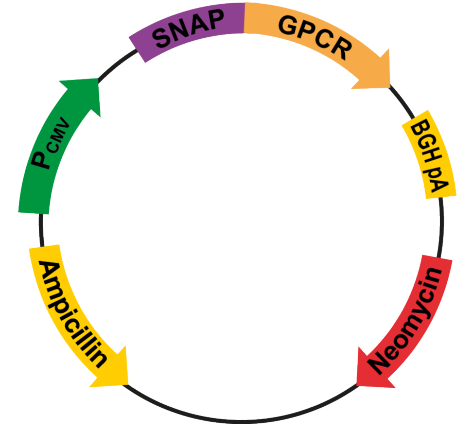
**\*based on NCBI nucleic sequence accession number**

**Promoter:** CMV

**Host:** Mammalian

**Antibiotic resistance for E. Coli:** Ampicillin

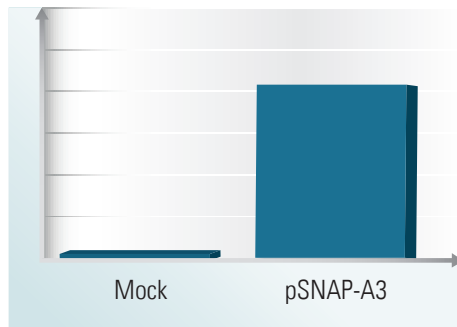
**Mammalian selection marker:** Neomycin



Tag-lite pSNAP-A3

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

## QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pSNAP-A3	10

\*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 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 & MSDS

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

**Sequence:** N° 32 F-Snap-N-A3 (923bp - 2551bp, direct) 1629bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCCCTGG
ATAGCCCTCTGGGCAAGCTGGAAGTGTCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCCCAGCCGCCGTGCTGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGCT
CAACGCCACTTTACCAGCCTGAGGCCATCGAGGAGTTCCTGTGCCAGCCCTGCACCACCCAGTGTCCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGAAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTCATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCGCCGTGAAAACCGCCCTGAGCGGAAATCCCGTGCCATTCTGATCCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGCTGCTGGCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCATGCCAACACAGCACTGCTCTGTCATTGGCCAATGTTACCTAC
ATCACCATGGAAATTTTCATTGGACTCTGCCCATAGTGGGCAACGTGCTGGTCATCTGCGTGGTCAAGCTGAACCCAG
CCTGCAGACCACCTTCTATTTTCATTGTCTCTAGCCCTGGCTGACATTGCTGTTGGGGTGGTGGTCAAGCTGACCTTTGG
CCATTGTTGTCAGCCTGGGCATCACAATCCACTTCTACAGCTGCCTTTTATGACTTGCCTACTGCTTATCTTTACCCAC
GCCTCCATCATGTCCTTGTGGCCATCGCTGTGGACCGATACTGCGGGTCAAGCTTACCGTCAGATAACAAGAGGGTCA
CACTCACAGAAGAATATGGCTGGCCCTGGGCCTTTGCTGGCTGGTGTATTCCTGGTGGGATTGACCCCATGTTTGGCT
GGAACATGAAACTGACCTCAGAGTACCACAGAAATGTCACCTTCTTTTCATGCCAATTTGTTTCCGTCATGAGAATGGAC
TACATGGTATACTTACGCTTCTCACCTGGATTTTCATCCCCTGGTGTGTCATGTGCGCCATCTATCTTGACATCTTTTA
CATCATTGGAAACAACTCAGTCTGAAGTATCTAACTCCAAAGAGACAGGTGCATTTTATGGACGGGAGTTCAAGACGG
CTAAGTCCTTGTCTGGTCTTTTCTTGTGCTGTGTCATGGCTGCCTTTATCTATCATCAACTGCATCATCTACTTT
AATGGTGAGGTACCACAGCTTGTGCTGTACATGGGCATCCTGCTGTCCCATGCCAACTCCATGATGAACCCATCGTCTA
TGCCTATAAAATAAAGAAGTTCAGGAAACCTACCTTTTGATCCTCAAAGCCTGTGTGGTCTGCCATCCCTCTGATTCTT
TGGACACAAGCATTGAGAAGAATTCTGAG
```

Copyright 2018 Cisbio Bioassays. All rights reserved. HTRF, Tag-lite, EPLigeneous and the HTRF logo are trademarks or registered trademarks of Cisbio Bioassays. CLIP-tag is a registered trademark of New England Biolabs, Inc. HaloTag is a registered trademark of Promega Corporation. SNAP-tag is a registered trademark of New England Biolabs, Inc.

### FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86-21-5018-9880 Japan +81-(0)43-306-8712

Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors

