



SerotoninPLASMID TAG-LITE® pSNAP-5HT1A Receptor: 5-Hydroxytryptamine receptor type 1A

Product Information Revision: #04 of September 2023

Part#: PSNAP5HT1A

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H₂O

Plasmid name: pSNAP-5HT1A

Plasmid size: 7307 bp

5HT1A-ORF size: 1938 bp

ORF sequence*: NM_000524.2

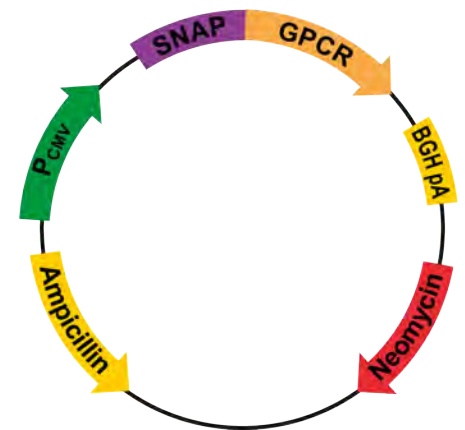
***based on NCBI nucleic sequence accession number**

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

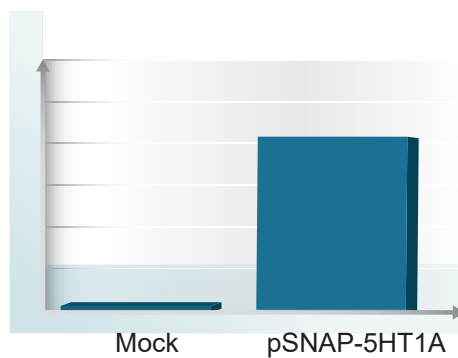
Mammalian selection marker: Neomycin



Tag-lite pSNAP-5HT1A

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

QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pSNAP-5HT1A	11

*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 manual described on the web site.

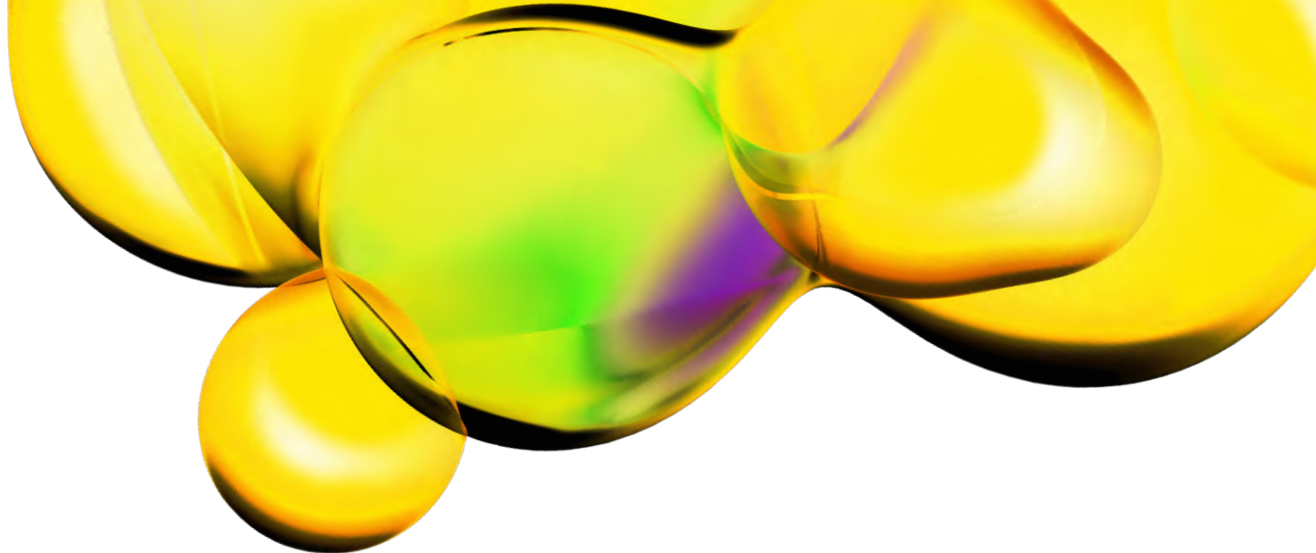
Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

SEQUENCE & MSDS

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

Sequence: F-Snap-N-5HT1A (923bp - 2860bp, direct) 1938bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCACCCTGG
ATAGCCCTCTGGGCAAGCTGGAAGTGTCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCACGCCCGCTGCTGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGCT
CAACGCCTACTTTACCAGCCTGAGGCCATCGAGGAGTTCCCTGTGCCAGCCCTGCACCACCCAGTGTTCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGAAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTCATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCCGCTGAAAACCGCCCTGAGCGGAAATCCCGTGCCCATTTGATCCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGCTGCTGGCCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCGATGTGCTCAGCCCTGGTCAGGGCAACAACACCACATCACCACCG
GCTCCCTTTGAGACCGGCCGGAACACTACTGGTATCTCCGACGTGACCGTCAGCTACCAAGTGATCACCTCTCTGCTGCT
GGGCACGCTCATCTTCTGCGCGGTGCTGGGCAATGCGTGCGTGGTGGCTGCCATCGCCTTGGAGCGCTCCCTGCAGAACG
TGGCCAATTATCTTATTGGCTCTTTGGCGGTCACCGACCTCATGGTGTGCGGTGTTGGTGTGCCATGGCCGCGCTGTAT
CAGGTGCTCAACAAGTGGACACTGGGCCAGGTAACCTGCGACCTGTTTATCGCCCTCGACGTGCTGTGCTGCACCTCATC
CATCTTGCACCTGTGCGCCATCGCGCTGGACAGGTAAGTGGGCCATCACGGACCCCATCGACTACGTGAACAAGAGGACGC
CCCGCGCGCCGCTGCGCTCATCTGCTCACTTGGCTTATTGGCTTCTCATCTATCCCGCCCATGCTGGGCTGGCGC
ACCCCGGAAGACCGCTCGGACCCCGACGCATGCACCATAGCAAGGATCATGGCTACACTATCTATTCCACCTTTGGAGC
TTTCTACATCCCGCTGCTGCTCATGCTGGTTCTCTATGGGCGCATATTCGAGCTGCGCGCTTCCGCATCCGCAAGACGG
TCAAAAAGGTGGAGAAGACCGGAGCGGACACCCGCCATGGAGCATCTCCCGCCCGCAGCCCAAGAAGAGTGTGAATGGA
GAGTCGGGGAGCAGGAAGTGGAGGCTGGGCGTGGAGAGCAAGGCTGGGGGTGCTCTGTGCGCCAATGGCGCGGTGAGGCA
AGGTGACGATGGCGCCGCCCTGGAGGTGATCGAGGTGCACCGAGTGGGCAACTCCAAAGAGCACTTGCCCTGCCCCAGCG
AGGCTGGTCTTACCCCTTGTGCCCCCGCCTCTTTGAGAGGAAAAATGAGCGCAACGCCGAGGCGAAGCGCAAGATGGCC
CTGGCCCCGAGAGAGGAAGACAGTGAAGACGCTGGGCATCATCATGGGCACCTTCATCCTCTGCTGGCTGCCCTTCTCAT
CGTGGCTCTTGTCTGCCCTTCTGCGAGAGCAGCTGCCACATGCCACCCTGTTGGGCGCCATAATCAATTGGCTGGGCT
ACTCCAACCTCTGCTTAAACCCGTCATTTACGCATACTTCAACAAGGACTTTCAAACGCGTTTAAAGAAGATCATTAAG
TGTAAGTTCTGCCGCCAG
```



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.