



# Serotonin PLASMID TAG-LITE® pSNAP-HTR4 Receptor: 5-Hydroxytryptamine receptor type 4

Product Information Revision: #03 of September 2023

**Part#:** PSNAPHTR4

**Storage temperature:** ≤-16°C

**Quantity:** 10 µg/ 10 µL H2O

**Plasmid name:** pSNAP-HTR4

**Plasmid size:** 7205 bp

**5HT4-ORF size:** 1836 bp

**ORF sequence\*:** NM\_001040172.1

\*based on NCBI nucleic sequence accession number

**Promoter:** CMV

**Host:** Mammalian

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

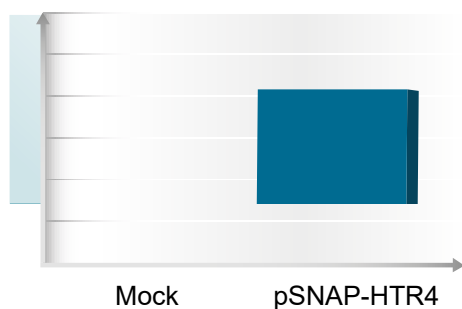
**Mammalian selection marker:** Neomycin



Tag-lite pSNAP-HTR4

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

## QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pSNAP-HTR4	11

\*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.

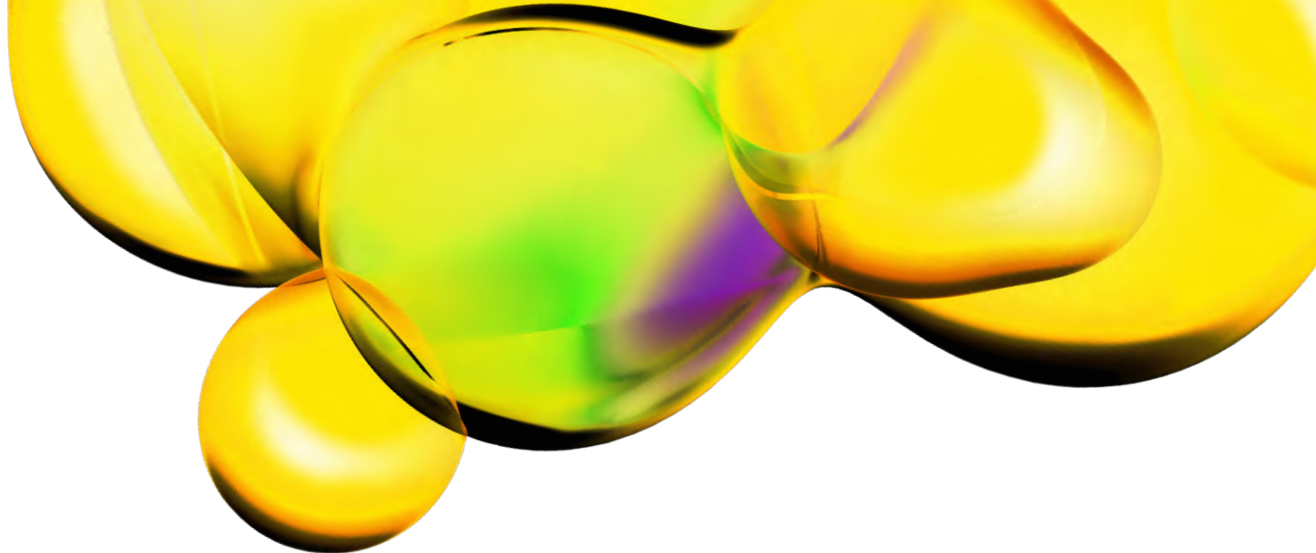
Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

## SEQUENCE & MSDS

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

**Sequence:** N°125) pT8-flag-ST26m-5HT4 (923bp - 2758bp, direct) 1836bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCACCCTGG
ATAGCCCTCTGGGCAAGCTGGAAGTGTCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCCCAGCCGCGTGTGGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGCT
CAACGCCTACTTTACCAGCCTGAGGCCATCGAGGAGTTCCCTGTGCCAGCCCTGCACCACCCAGTGTTCCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGGAAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCGCGTGAAAACCGCCCTGAGCGGAAATCCCCTGCCATTCTGATCCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGCTGCTGGCCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCGACAACTTGATGCTAATGTGAGTTCTGAGGAGGGTTTCGGGTCA
GTGGAGAAGGTGGTGTCTGCTCACGTTTCTCTCGACGGTTATCCTGATGGCCATCTTGGGGAACCTTGCTGGTGTGGTGGC
TGTGTGCTGGGACAGGCAGCTCAGGAAAATAAAAAACAATTATTTTCATTGTATCTCTTGCTTTTGCGGATCTGCTGGTTT
CGGTGCTGGTGTATGCCCTTTGGTGCCATTGAGCTGGTTCAAGACATCTGGATTTATGGGGAGGTGTTTTGTCTTGTTCGG
ACATCTCTGGACGTCCTGCTACAACGGCATCGATTTTTCACCTGTGCTGCATTTCTCTGGATAGGTATTACGCCATCTG
CTGCCAGCCTTTGGTCTATAGGAACAAGATGACCCCTCTGCGCATCGCATTAATGCTGGGAGGCTGCTGGGTGATCCCA
CGTTTATTTCTTTTCTCCCTATAATGCAAGGCTGGAATAACATTGGCATAATTGATTTGATAGAAAAGAGGAAGTTCAAC
CAGAACTCTAACTCTACGTACTGTGTCTTCATGGTCAACAAGCCCTACGCCATCACCTGCTCTGTGGTGGCCTTCTACAT
CCCATTTCTCCTCATGGTGTGGCCTATTACCGCATCTATGTCACAGCTAAGGAGCATGCCATCAGATCCAGATGTTAC
AACGGGCAGGAGCCTCCTCCGAGAGCAGGCCTCAGTCGGCAGACCAGCATAGCACTCATCGCATGAGGACAGAGACAAA
GCAGCCAAGACCCTGTGCATCATATGGGTTGCTTCTGCCTCTGCTGGGCACCATTCTTTGTACCAATATTGTGGATCC
TTTCATAGACTACACTGTCCCTGGGCAGGTGTGGACTGCTTTCCCTGCTGGCTCGGCTATATCAATTCCGGGTGAACCCTT
TTCTCTACGCCTTCTTGAATAAGTCTTTTAGACGTGCCTTCTCATCATCCTCTGCTGTGATGATGAGCGCTACCGAAGA
CCTTCCATTCTGGGCCAGACTGTCCCTTGTTCACCAACAACCATTAATGGATCCACACATGTAATAAGGGATGCAGTGGA
GTGTGGTGGCCAGTGGGAGAGTCAAGTGTCAACCCGCCAGCAACTTCTCCTTTGGTGGCTGCTCAGCCCAGTGACACT
```



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.