



OrexinPLASMID TAG-LITE® pSNAP-OX2

Receptor: Orexin receptor 2

PRODUCT INFORMATION - rev. 2 (Sept. 2018)

Part#: PSNAPOX2

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H2O

Plasmid name: pSNAP-OX2

Plasmid size: 7385 bp

OX2-ORF size: 2022 bp

ORF sequence*: NM_001526.3

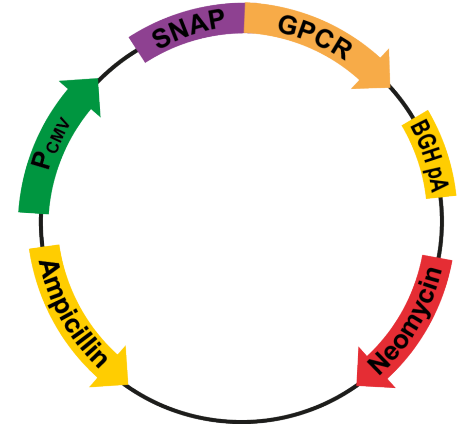
***based on NCBI nucleic sequence accession number**

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

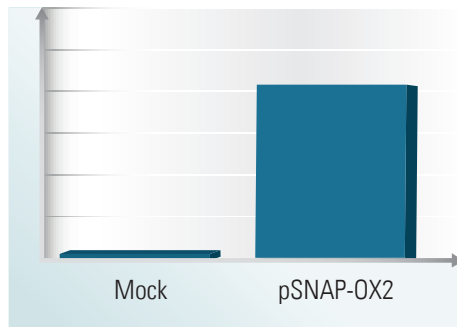
Mammalian selection marker: Neomycin



Tag-lite pSNAP-OX2

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

QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pSNAP-OX2	5

*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.

Sequence: F-Snap-N-OX2R GC (923bp - 2944bp, direct) 2022bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCACCTGG
ATAGCCCTCTGGGCAAGCTGGAACCTGCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCCCAGCCGCCGTGCTGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGCT
CAACGCCACTTTACCAGCCTGAGGCCATCGAGGAGTTCCTGTGCCAGCCCTGCACCACCCAGTGTCCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGAAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTGATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCGCCGTGAAAACCGCCCTGAGCGGAAATCCCGTGCCATTCTGATCCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGTGTGGCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCCAGCACAGTGGCGGCCGCTCCGGCACCAATGGAGGACTCCCCC
CCTTGTGCGCAACTGGTCACTGCTTCGGAGCTGAATGAAACTCAAGAGCCCTTTTAAACCCACCGACTATGACGACGA
GGAATTCCTGCGGTACCTGTGGAGGGAATACCTGCACCCGAAAGAATATGAGTGGGTCTGATCGCCGGGTACATCATCG
TGTTGCTGCTGGCTCTCATTGGGAACGCTCCTGGTTTGTGTGGCAGTGTGGAAGAACCACCACATGAGGACGGTAACCAAC
TACTTCATAGTCAATCTTCTCTGGCTGATGTGCTCGTGACCATCACCTGCCTTCCAGCCACACTGGTCTGGATATCAC
TGAGACCTGGTTTTTGGACAGTCCCTTTGCAAAGTGATTCCCTTATCTACAGACCGTGTGCGGTGCTGTGTCTGTCTCA
CACTGAGCTGTATCGCCTGGATCGGTGGTATGCAATCTGTACCCTTTGATGTTAAGAGCACAGCAAAGCGGGCCCGT
AACAGCATTGTATCATCTGGATTGCTCCTGCATTATAATGATTCCTCAGGCCATCGTATGGAGTGCAGCACCGTGT
CCCAGGCTTAGCCAATAAAACCACCTCTTACGGTGTGTGATGAGCGCTGGGGTGGTGAATTTATCCCAAGATGTACC
ACATCTGTTTCTTCTGGTGACATACATGGCACCCTGTGTCTCATGGTGTGGCTTATCTGCAAATATTCGCAAACCTC
TGGTGTGACAGATCCCTGGAACATCATCTGTAGTTCAGAGAAAATGGAAGCCCTGCAGCCTGTTTACAGCCTCGAGG
GCCAGGACAGCCAACGAAGTCCCGGATGAGCGCTGTGGCGGCTGAAATAAAGCAGATCCGAGCCAGAAGGAAAACAGCCC
GGATGTTGATGATTGTGCTTTTGGTATTGCAATTTGCTATACCAATTAGCATCCTCAATGTGCTAAAGAGAGTATTT
GGGATGTTTGGCCATACTGAAGACAGAGAGACTGTGTATGCCTGGTTTACCTTTTACACTGGCTTGTATATGCCAATAG
TGCTGCGAATCCAATATTTATAATTTTCTCAGTGGAAAATTTCCGAGAGGAATTTAAAGCTGCGTTTTCTTGCTGTTGCC
TTGGAGTTCACCATCGCCAGGAGGATCGGCTCACCAGGGGACGAACTAGCACAGAGAGCCGGAAGTCTTGACCACTCAA
ATCAGCAACTTTGATAACATATCAAACCTTTCTGAGCAAGTTGTGCTCACTAGCATAAGCACACTCCCAGCAGCCAATGG
AGCAGGACCACTTCAAACTGG
```

Copyright 2018 Cisbio Bioassays. All rights reserved. HTRF, Tag-lite, EPLgeneous 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 to find a list of our regional distributors

