



Angiotensin PLASMID TAG-LITE® pSNAP-AT2 Receptor: Angiotensin receptor type 2

Product Information Revision: #04 of September 2023

Part#: PSNAPAT2

Storage temperature: ≤-16°C

Quantity: 10 µg/ 10 µL H2O

Plasmid name: pSNAP-AT2

Plasmid size: 7133 bp

AT2-ORF size: 1764 bp

ORF sequence*: NM_000686.4

*based on NCBI nucleic sequence accession number

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

Mammalian selection marker: Neomycin



Tag-lite pSNAP-AT2

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

QUALITY CONTROL

Mock pSNAP-AT2

Plasmid	Normalized 620 nm signal*
pSNAP-AT2	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 manual described on the web site.

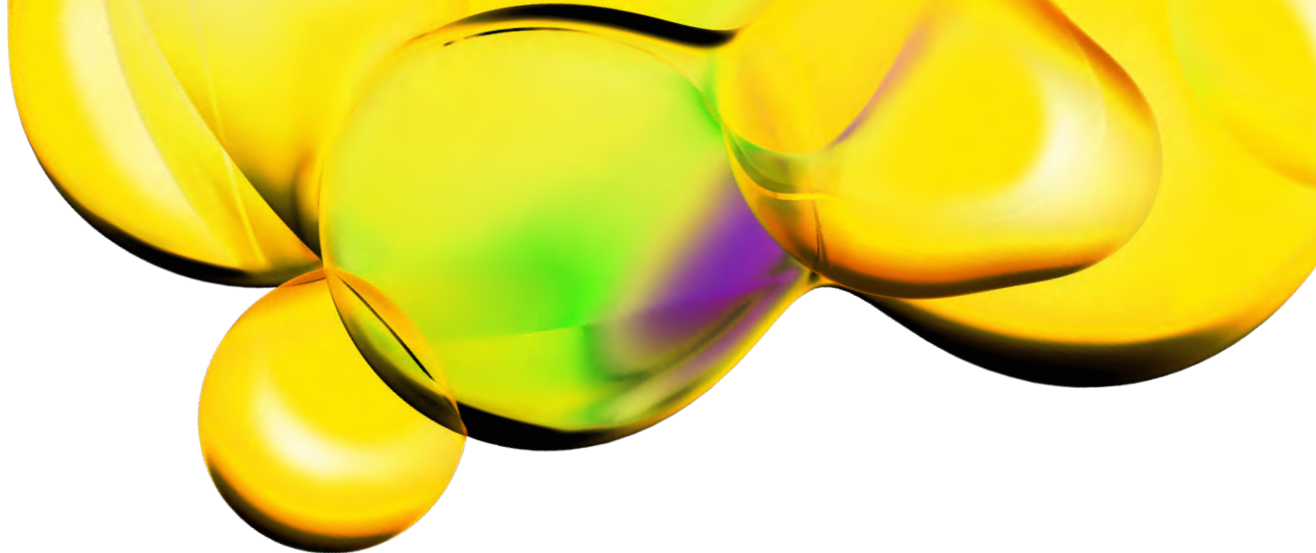
Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

SEQUENCE & MSDS

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

Sequence: N° 67 F-Snap-N-AGTR2 (923bp - 2686bp, direct) 1764bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGACTACAAGGACGACGATGACAAGGCCGGCCATCGATGCCATCATGGACAAAGACTGCGAAATGAAGCGCACCACCCTGG
ATAGCCCTCTGGGCAAGCTGGAAGTGTCTGGGTGCGAACAGGGCCTGCACGAGATCAAGCTGCTGGGCAAAGGAACATCT
GCCGCCGACGCCGTGGAAGTGCCTGCCCCAGCCGCCGTGCTGGGCGGACCAGAGCCACTGATGCAGGCCACCGCCTGGCT
CAACGCCTACTTTACCAGCCTGAGGCCATCGAGGAGTTCCCTGTGCCAGCCCTGCACCACCCAGTGTTCAGCAGGAGA
GCTTTACCCGCCAGGTGCTGTGGAAACTGCTGAAAGTGGTGAAGTTCGGAGAGGTCATCAGCTACCAGCAGCTGGCCGCC
CTGGCCGGCAATCCCGCCGCCACCGCCGCCGTGAAAACCGCCCTGAGCGGAAATCCCGTGCCCATTTCTGATCCCCTGCCA
CCGGGTGGTGTCTAGCTCTGGCGCCGTGGGGGGCTACGAGGGCGGGCTCGCCGTGAAAGAGTGGCTGCTGGCCACGAGG
GCCACAGACTGGGCAAGCCTGGGCTGGGTGATATCATGAAGGGCAACTCCACCCTTGCCACTACTAGCAAAAAACATTACC
AGCGGTCTTCACTTCGGGCTTGTGAACATCTCTGGCAACAATGAGTCTACCTTGAAGTGTTCACAGAAACCATCAGATAA
GCATTTAGATGCAATTCCTATCTTTACTACATTATATTTGTAATTGGATTTCTGGTCAATATGTCGTGGTTACTGT
TTTGTGTCAAAAAGGGTCCAAAAAGGTTTCTAGCATATACATCTTCAACCTCGCTGTGGCTGATTTACTCCTTTTGGCT
ACTCTTCTCTATGGGCAACCTATTATCTTATAGATATGACTGGCTCTTTGGACCTGTGATGTGCAAAGTTTTTGGTTC
TTTTCTTACCCTGAACATGTTTGCAAGCATTTTTTTATCACCTGCATGAGTGTGATAGGTACCAATCTGTCATCTACC
CCTTCTGTCTCAAAGAAGAAATCCCTGGCAAGCATCTTATATAGTTCCCCTTGTTGGTGTATGGCCTGTTTGCCTCA
TTGCCAACATTTATTTTCGAGACGTCAGAACCATTGAATACTTAGGAGTGAATGCTTGCATTATGGCTTTCCACCTGA
GAAATATGCCCAATGGTCAGCTGGGATTGCCTTAATGAAAAATATCCTTGGTTTTATATCCCTTTAATATTCATAGCAA
CATGCTATTTTGAATTAGAAAACCTTACTGAAGACGAATAGCTATGGGAAGAACAGGATAACCCGTGACCAAGTCCTG
AAGATGGCAGCTGCTGTTGTTCTGGCCTTCATCATTTGCTGGCTTCCCTTCCATGTTCTGACCTTCTGGATGCTCTGGC
CTGGATGGGTGTCATTAATAGCTGCGAAGTTATAGCAGTCATTGACCTGGCACTTCCTTTTGCCATCCTCTTGGGATTCA
CCAACAGCTGCGTTAATCCGTTTCTGTATTGTTTTGTTGGAAACCGGTTCCAACAGAAGCTCCGCAGTGTGTTAGGGTT
CCAATTACTTGGCTCCAAGGGAAAAGAGAGAGATATGCTTGGCGAAAAGCAGTTCTTAGAGAAATGGAGACCTTTGTGTCT
```



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.