



# PLASMID TAG-LITE® SNAP T8

Product Information Revision: #03 of September 2023

Part#: PT8SNAPHYG

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H2O

Plasmid name: Tag-lite pT8-SNAP

Plasmid size: 6231bp

T8-SNAP-ORF size from ATG to MCS: 704bp

Tag name: SNAP

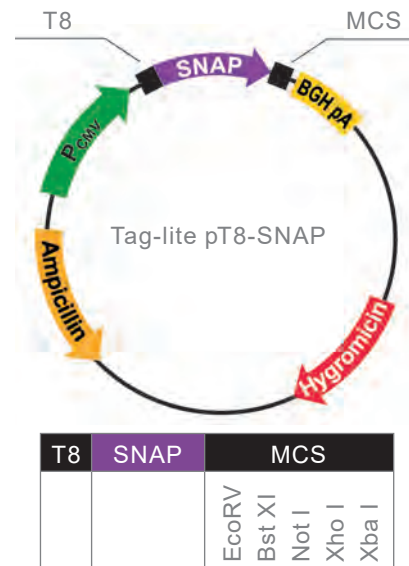
Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

Mammalian selection marker: Hygromycin

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



## T8-SNAP-ORF SEQUENCE FROM ATG TO MCS (end)

1	ATGGCCTTAC	CAGTGACCGC	CTTGCTCCTG	CCGCTGGCCT	TGCTGCTCCA	CGCCGCCAG	CGGCGCCCG	CTAGCGGCAT	CGACTACAAG	GACGACGATG
	TACCGGAATG	GTCACCTGGC	GAACGAGGAC	GGCGACCGGA	ACGACGAGGT	GGGGCGGTCC	GGCCGGCGGC	GATCGCGGTA	GCTGATGTTT	CTGCTGCTAC
		Clal								
101	ACAAGGCCGG	CATCGATGCC	ATCATGGACA	AAGACTGCGA	AATGAAGCGC	ACCACCCTGG	ATAGCCCTCT	ATAGCCCTCT	GAAGTGTCTG	GGTGCGAACA
	TGTTCCGGCC	GTAGCTACGG	TAGTACCTGT	TTCTGACGCT	TTACTTCGCG	TGGTGGGACC	TATCGGGAGA	CCCGTTGCAC	CTTGACAGAC	CCACGCTTGT
		DpnI								
201	GGGCCTGCAC	GAGATCAAGC	TGCTGGGCAA	AGGAACATCT	GCCGCCGACG	CCGTGGAAGT	GCCTGCCCA	GCCGCCGTGC	TGGGCGGACC	AGAGCCACTG
	CCCGGACGTG	CTCTAGTTCC	ACGACCCGTT	TCCTTGTAGA	CGGCGGCTGC	GGCACCTTCA	CGGACGGGGT	CGGCGGCACG	ACCCGCCTGG	ACCCGCCTGG
301	ATGCAGGCCA	CCGCCTGGCT	CAACGCCTAC	TTTCACCAGC	CTGAGGCCAT	CGAGGAGTTC	CCTGTGCCAG	CCCTGCACCA	CCCAGTGTTT	CAGCAGGAGA
	TACGTCCGGT	GGCGGACCGA	GTTGCGGATG	AAAGTGGTCG	GACTCCGGTA	GCTCCTCAAG	GGACACGGTC	GGGACGTGGT	GGGTACACAAG	GTCGTCTCTT
							PvuII			
401	GCTTTACCCG	CCAGGTGCTG	TGGAACTGCG	TGAAAGTGGT	TGAAAGTGGT	GAGGTCATCA	GCTACCAGCA	GCTGGCCGCC	CTGGCCGGCA	ATCCCGCCGC
	CGAAATGGGC	CGAAATGGGC	ACCTTTGACG	ACTTTCACCA	CTTCAAGCCT	CTCCAGTAGT	CGATGGTCGT	CGACCGGCCG	GACCGGCCGT	TAGGGCGGCC
					DpnI					
501	CACCGCCGCC	GTGAAAACCG	CCCTGAGCGG	AAATCCCGTG	CCCATTCTGA	TCCCTGCCA	CCGGTGGTG	TCTAGCTCTG	GCGCCGTGGG	GGGCTACGAG
	GTGGCGCGCG	CACCTTTGGC	GGGACTCGCC	TTTAGGGCAC	GGGTAAGACT	AGGGGACGGT	GGCCACCAC	AGATCGAGAC	CGCGGCACCC	CCCGATGCTC
							EcoRV		NotI	
										XhoI
										AvaI
601	GGCGGGCTCG	CCGTGAAAGA	GTGGCTGCTG	GCCCACGAGG	GCCACAGACT	GGGCAAGCCT	GGGCTGGGTG	ATATCCAGCA	CAGTGGCGGC	CGCTCGAGTC
	CCGCCCGAGC	GGCACTTTCT	CACCGACGAC	CGGGTGTCTC	CGGTGTCTGA	CCCGTTCGGA	CCCGACCCAC	TATAGTCTGT	GTCACCGCCG	GCGAGCTCAG
701	TAGA									
	AICT									

## MSDS

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



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.