



PLASMID TAG-LITE® SNAP T8

Product Information Revision: #03 of September 2023

Part#: PT8SNAPNEO

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H2O

Plasmid name: Tag-lite pT8-SNAP

Plasmid size: 6058bp

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

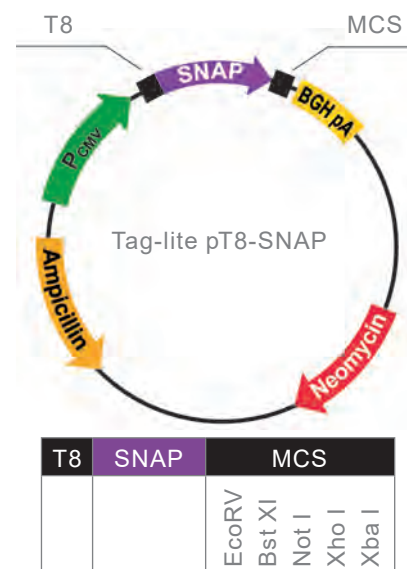
Tag name: SNAP

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

Mammalian selection marker: Neomycin



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	CCGGCCCGCG	CTAGCCGCAT	CGACTACAAG	GACGACGATG
	TACCGGAATG	GTCACCTGGC	GAACGAGGAC	GGCGACCGGA	ACGACGAGGT	GCGGCGGTCC	GGCCGGCGGC	GATCGCCGTA	GCTGATGTTT	CTGCTGCTAC
		Clal								
101	ACAAGGCCGG	CATCGATGCC	ATCATGGACA	AAGACTGCGA	AATGAAGCGC	ACCACCTGG	ATAGCCCTCT	ATAGCCCTCT	GAAGTGTCTG	GGTGCGAACA
	TGTTCCGGCC	GTAGCTACGG	TAGTACCTGT	TTCTGACGCT	TTACTTCGGG	TGGTGGGACC	TATCGGGAGA	CCCGTTCGAC	CTTGACAGAC	CCACGCTTGT
		DpnI								
201	GGGCCTGCAC	GAGATCAAGC	TGCTGGGCAA	AGGAACATCT	GCCGCCGACG	CCGTGGAAGT	GCCTGCCCA	GCCGCCGTGC	TGGCGGACC	AGAGCCACTG
	CCCGGACGTG	CTTAGTTCG	ACGACCCGTT	TCCTTGTAGA	CGGCGGCTGC	GGCACCTTCA	CGGACGGGGT	CGGCGGCACG	ACCCGCTGG	ACCCGCTGG
301	ATGCAGGCCA	CCGCCTGGCT	CAACGCCTAC	TTTCACCAGC	CTGAGGCCAT	CGAGGAGTTC	CCTGTGCCAG	CCCTGCACCA	CCAGTGTTC	CAGCAGGAGA
	TACGTCCGGT	GGCGGACCGA	GTTGCGGATG	AAAGTGGTCG	GACTCCGGTA	GCTCCTCAAG	GGACACGGTC	GGGACGTGGT	GGTCCACAAG	GTCGTCCTCT
							PvuII			
401	GCTTTACCCG	CCAGGTGCTG	TGGAAACTGC	TGAAAGTGGT	TGAAAGTGGT	GAGGTCATCA	GCTACCAGCA	GCTGGCCGCC	CTGGCCGGCA	ATCCCGCCGC
	CGAAATGGGC	CGAAATGGGC	ACCTTTGACG	ACTTTCACCA	CTTCAAGCCT	CTCCAGTAGT	CGATGGTCGT	CGACCGGCGG	GACCGGCCGT	TAGGGCGGCG
					DpnI					
501	CACCGCCGCC	GTGAAAACCG	CCCTGAGCGG	AAATCCCGTG	CCCATTCTGA	TCCCTGCCA	CCGGTGGTG	TCTAGCTCTG	GCGCCGTGGG	GGGCTACGAG
	GTGGCGGCGG	CACTTTTGGC	GGGACTCGCC	TTTAGGGCAC	GGGTAAGACT	AGGGGACGGT	GGCCACCAC	AGATCGAGAC	CGCGGCACCC	CCCAGTGCTC
							EcoRV		NotI	
									XhoI	
									AvaI	
601	GGCGGGCTCG	CCGTGAAAGA	GTGGCTGCTG	GCCCACGAGG	GCCACAGACT	GGGCAAGCCT	GGGCTGGGTG	ATATCCAGCA	CAGTGGCGGC	CGCTCGAGTC
	CCGCCCGAGC	GGCACITTTCT	CACCGACGAC	CGGGTGTCTC	CGGTGTCTGA	CCGTTTCGGA	CCCGACCCAC	TATAGTTCGT	GTCACCGCCG	GCGAGCTCAG
		XbaI						BstXI		
701	TAGA									
	AICT									

MSDS

Files can be downloaded from Tag-lite section on the website 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

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.