

MANUAL

Technology: HTRF™ Tag-lite

Plasmid Tag-lite™ SNAP T8

Part number PT8SNAPZEO

Storage temperature: -20°C

Quantity: 10µg/10µL H₂O

Version: 04

Date: October 2024

Plasmid name: Tag-lite pT8-SNAP

Plasmid size: 5645bp

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

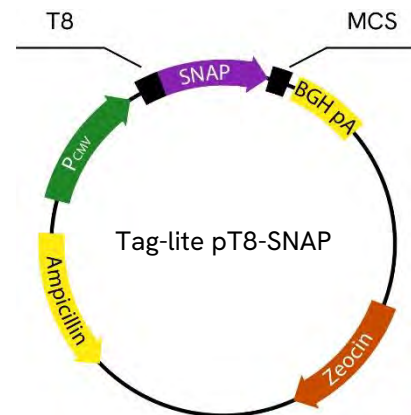
Tag name: SNAP

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

Mammalian selection marker: Zeocin



T8	SNAP	MCS				
		EcoRV	BstXI	NotI	XhoI	XbaI

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

1	AAGCTTTGAG	ACATGGCCTT	ACCAGTGACC	GCCTTGCTCC	TGCCGCTGGC	CTTGCTGCTC	CACGCCGCCA	GGCCGGCCGC	CGCTAGCGGC	ATCGACTACA
	TTCGAAACTC	TGTACCGGAA	TGTCACTGG	CGGAACGAGG	ACGGCGACCG	GAACGACGAG	GTGCGGCGGT	CCGGCCGGCG	GCGATCGCCG	TAGCTGATGT
101	AGGACGACGA	TGACAAGGCC	GGCATCGATG	CCATCATGGA	CAAAAGACTGC	GAAATGAAGC	GCACCACCCT	GGATAGCCCT	CTGGGCAAGC	TGGAACGTTT
	TCCTGCTGCT	ACTGTTCCGG	CCGTAGCTAC	GGTAGTACCT	GTTTCTGACG	CTTTACTTCG	CGTGGTGGA	CCTATCGGGA	GACCCGTTTG	ACCTTGCAAG
201	TGGGTGCGAA	CAGGGCCTGC	ACGAGATCAA	GCTGCTGGGC	AAAGGAACAT	CTGCCGCCGA	CGCCGTGGAA	GTGCTGCC	CAGCCGCCGT	GCTGGGCGGA
	ACCCACGCTT	GTCCCGGACG	TGCTCTAGTT	CGACGACCCG	TTTCCGTTGA	GACGGCGGCT	GCGGCACCTT	CACGGACGGG	GTGCGCGGCA	CGACCCGCTT
301	CCAGAGCCAC	TGATGAGGC	CACCGCCTGG	CTCAACGCCT	ACTTTCACCA	GCCTGAGGCC	ATCGAGGAGT	TCCCTGTGCC	AGCCCTGCAC	CACCCAGTGT
	GGTCTCGGTG	ACTACGTCCG	GTGGCGGACC	GAGTTGCGGA	TGAAAGTGGT	CGGACTCCGG	TAGCTCCTCA	AGGGACACGG	TGCGGACGTG	GTGGTGCACA
401	TCCAGCAGGA	GAGCTTTACC	CGCCAGGTGC	TGTGAAACT	GCTGAAAGTG	GTGAAGTTTG	GAGAGGTCAT	CAGCTACCAG	CAGCTGGCCG	CCCTGGCCGG
	AGGTGTCCT	CTCGAAATGG	GCGGTCCACG	ACACCTTTGA	CGACTTTCAC	CACTTCAAGC	CTCTCCAGTA	GTGATGGTGC	GTGACCGGCG	GGACCGGCC
501	CAATCCCGCC	GCCACCGCCG	CGTGAAATC	CGCCCTGAGC	GGAAATCCCG	TGCCATTCT	GATCCCTGTC	CACCGGGTGG	TGTCTAGCTC	TGCGCCGTG
	GTTAGGGCGG	CGGTGGCGGC	GGCACTTTTG	GCGGGACTCG	CTTTAGGGCG	ACGGGTAAGA	CTAGGGACGG	GTGGCCACCC	ACAGATCGAG	ACGCGGCAC
601	GGGGGCTACG	AGGGCGGGCT	CGCCGTGAAA	GAGTGGCTGC	TGGCCACGA	GGGCCACAGA	CTGGGCAAGC	CTGGGCTGGG	TGATATCAGG	CACAGTGGCG
	CCCCGAAGC	TCCCGCCCGG	AGCGGCACTT	TCTCACCAGC	GACCGGGTGC	TCCCGGTGTC	TGACCCGTTT	GGACCCGACC	ACTATAGGTC	GTGTCACCGC
701	GCCGCTCGAG	TCTAGAGGGC	CCGTTTAA							
	CGGCAGGCTC	AGATCTCCCG	GGCAAATT							

MSDS

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