



PLASMID TAG-LITE® CLIP T8

Product Information Revision: #03 of September 2023

Part#: PT8CLIPNEO

Storage temperature: -20°C

Quantity: 10 µg/ 10 µL H2O

Plasmid name: Tag-lite pT8-CLIP

Plasmid size: 6055bp

T8-CLIP-ORF size from ATG to MCS: 701bp

Tag name: CLIP

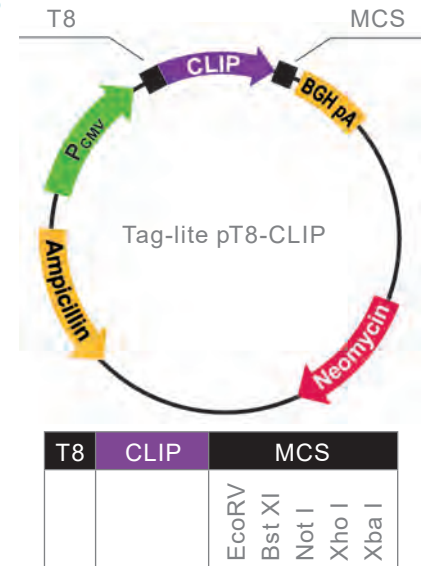
Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

Mammalian selection marker: Neomycin

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



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

								NheI		
1	ATGGCCTTAC	CAGTGACCGC	CTTGCTCCTG	CGGCTGGCCT	TGCTGTCCA	CGCCGCCAGG	CGGGCCGCGC	CTAGCGGCAT	CGACTACAAG	GACGACGATG
	TACCGGAATG	GTCACCTGGCG	GAACGAGGAC	GGCGACCGGA	ACGACGAGGT	GCGGCGGTCC	GGCCGGCGGC	GATCGCCGTA	GCTGATGTTT	CTGCTGTACT
		Clal						Alul		
101	ACAAGGCCGG	CATCGATAAC	ATGGACAAAG	ACTGCGAAAT	GAAGCGCACC	ACCCTGGATA	GCCCTCTGGG	CAAGCTGGAA	CTGTCTGGGT	GCGAACAGGG
	TGTTCCGGCC	GTAGCTATGG	TACCTGTTTC	TGACGCTTTA	CTTCGCGTGG	TGGGACCTAT	CGGGAGACCC	GTTGACCTT	GACAGACCCA	CGCTTGTCCC
201	CCTGCACGAG	ATCATCTTCC	TGGGCAAAGG	AAACATCTGC	GCCGACGCCG	TGGAAGTGCC	TGCCCCAGCC	GCCGTGCTGG	GCGGACCAGA	GCCACTGATC
	GGACGTGCTC	TAGTAGAAGG	ACCCGTTTCC	TTGTAGACGG	CGGCTGCGGC	ACCTTCACGG	ACGGGGTCGG	CGGCACGACC	CGCCTGGTCT	CGGTGACTAG
										Alul
301	CAGGCCACCG	CCTGGCTCAA	CGCCTACTTT	CACCAGCCTG	AGGCCATCGA	GGAGTTCCTT	GTGCCAGCCC	TGCACCACCC	AGTGTCCAG	CAGGAGAGCT
	GTCCGGTGGC	GGACCGAGTT	GCGGATGAAA	GTGGTCGGAC	TCCGGTAGCT	CCTCAAGGGA	CACGGTCCGG	ACGTGGTGGG	TCACAAGGTC	GTCTCTCGA
401	TTACCCGCCA	GGTGCTGTGG	AAACTGCTGA	AAAGTGGTAA	GTTTCGAGAG	GTCATCAGCG	AGAGCCACCT	GGCCGCCCTG	GTGGGCAATC	CCGCCGCCAC
	AATGGCGCGT	CCACGACACC	TTTGACGACT	TTCACCACTT	CAAGCCTCTC	CAGTAGTCGC	TCTCGGTGGA	CGGGCGGGAC	CACCCGTTAG	GGCGGCGGTG
										PspOMI
										ApaI
501	CGCCGCCGTG	AACACCGCCC	TGGACGAAA	TCCCGTCCCC	ATTCTGATCC	CCTGCCACCG	GGTGGTGCAG	GGCGACAGCG	ACGTGGGGCC	CTACCTGGGC
	GCGGGCGCAC	TTGTGGCGGG	ACCTGCCTTT	AGGGCACGGG	TAAGACTAGG	GGACGGTGGC	CCACCACGTC	CCGCTGTCGC	TGCACCCCGG	GATGGACCCG
							EcoRV		NotI	XbaI
										XhoI
										AvaI
601	GGGCTCGCCG	TGAAAGAGTG	GCTGCTGGCC	CACGAGGGCC	ACAGACTGGG	CAAGCCTGGG	CTGGGTGATA	TCCAGCACAG	TGGCGGCCCG	TCGAGTCTAG
	CCCGAGCGGC	ACTTTCTCAC	CGACGACCGG	GTGCTCCCGG	TGTTCTGACC	GTTCTGGACC	GACCCACTAT	AGGTCGTGTC	ACCGCCGGCG	AGCTCAGATC
701		A								
		T								

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.