

MANUAL

Technology: HTRF™

Tag-lite

Plasmid Tag-lite™ HaloTag T8

Part number

PT8HALONEO

Storage temperature: -20°C

Quantity: 10µg/10µL H₂O

Version: 04

Date: October 2024

Plasmid name: Tag-lite pT8-HALO

Plasmid size: 6406bp

T8- CLIP-ORF size from ATG to MCS: 1052bp

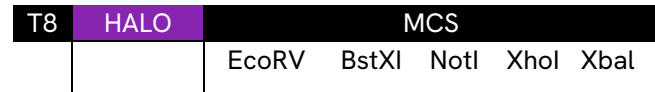
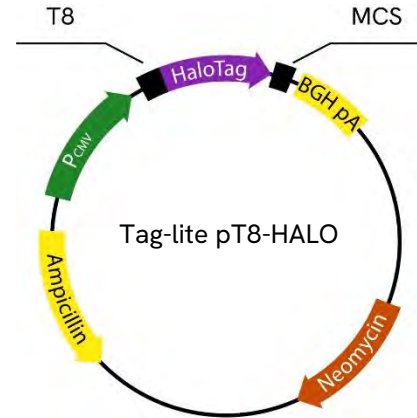
Tag name: HaloTag

Promoter: CMV

Host: Mammalian

Antibiotic resistance for E. Coli: Ampicillin

Mammalian selection marker: Neomycin



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

1	AAGCTTTGAG	ACATGGCCTT	ACCAGTGACC	GCCTTGCTCC	TGCCGCTGGC	CTTGCTGCTC	CACGCCGCCA	GGCCGGCCGC	CGCTAGCGGC	ATCGAGCAGA
	TTCGAAACTC	TGTACCGGAA	TGGTCACTGG	CGGAACGAGG	ACGGCGACCC	GAACGACGAG	GTGCGGCCGT	CCGGCCGGCG	GCGATCGCCG	TAGCTCGTCT
101	AGCTGATCAG	CGAGGAGGAC	CTGGCCGGCA	TCGATGCAGA	AATCGTACT	GGCTTCCAT	TCGACCCCA	TTATGTGAA	GTCTGGGCG	AGCGCATGCA
	TCGACTAGTC	GCTCCTCCTG	GACCGCCGCT	AGCTACGTCT	TTAGCCATGA	CCGAAAGGTA	AGCTGGGGT	AATACACCTT	CAGGACCCGC	TCGCGTACGT
201	CTACGTCGAT	GTTGGTCCGC	GCGATGGCAC	CCCTGTGCTG	TTCTGCACG	GTAACCCGAC	CTCCTCTAC	GTGTGGCGCA	ACATCATCCC	GCATGTTGCA
	GATGCAGCTA	CAACCAGGCG	CGCTACCGTG	GGGACACGAC	AAGGACGTGC	CATTGGGCTG	GAGGAGATG	CACACGCGTG	TGTAGTAGGG	CGTACAACGT
301	CCGACCCATC	GCTGCATTGC	TCCAGACCTG	ATCGGTATGG	GCAAATCCGA	CAAACCAGAC	CTGGGTATT	TCCTCGACGA	CCACGTCCGC	TTCATGGATG
	GGCTGGGTAG	CGACGTAACG	AGGTCTGGAC	TAGCCATACC	CGTTTAGGCT	GTTTGGTCTG	GACCAATAA	AGAAGTGTCT	GGTGCAGGCG	AAGTACCTAC
401	CCTTCATCGA	AGCCCTGGGT	CTGGAAGAGG	TCGTCCTGGT	CATTACGAC	TGGGGCTCCG	CTCTGGGTTT	CCACTGGGCC	AAGCGCAATC	CAGAGCCGCT
	GGAAGTAGCT	TCGGGACCCA	GACCTTCTCC	AGCAGGACCA	GTAAGTGTCT	ACCCCGAGGC	GAGACCCAAA	GGTGACCCGG	TTCGCTTAG	GTCTCGGCCA
501	CAAAGTATT	GCATTATGG	AGTTCATCCG	CCCTATCCCG	ACCTGGGACG	AATGGCCAGA	ATTTGCCCGC	GAGACCTTCC	AGGCCTTCCG	CACCACCGAC
	GTTTCCATAA	CGTAAATACC	TCAAGTAGGC	GGGATAGGGC	TGGACCCTGC	TTACCGGTCT	TAAACGGGCG	CTCTGGAAGG	TCCGGAAGGC	GTGGTGGCTG
601	GTCGGCCGCA	AGCTGATCAT	CGATCAGAAC	GTTTTTATCG	AGGGTACGCT	GCCGATGGGT	GTCGTCCGCC	CGCTGACTGA	AGTCGAGATG	GACCATTACC
	CAGCCGGCGT	TCGACTAGTA	GCTAGTCTTG	CAAAAATAGC	TCCCAGCGAC	GGCTACCCA	CAGCAGGCCG	GCGACTGACT	TCAGCTCTAC	CTGGTAATGG
701	GCGAGCCGTT	CCTGAATCCT	GTTGACCGCG	AGCCACTGTG	GCGCTTCCCA	AACGAGCTGC	CAATCGCCGG	TGAGCCAGCG	AACATCGTCG	CGCTGGTCTA
	CGCTCGGCAA	GGACTTAGGA	CAACGGCGCT	CGGTGACAC	CGCGAAGGGT	TTGCTCGACG	GTTAGCGGCC	ACTCGGTCCG	TTGTAGCAGC	GCGACCAGCT
801	AGAATACATG	GACTGGCTGC	ACCAGTCCCC	TGTCCTCAAG	CTGCTGTCT	GGGGACCCCG	AGGCGTCTG	ATCCCACCGG	CCGAAGCCCG	TCGCCTGGCC
	TCTTATGTAC	CTGACCGACG	TGGTGAGGGG	ACAGGGCTTC	GACGACAAGA	CCCCGTGGGG	TCCGCAAGAC	TAGGTGGCC	GGCTTCGGCG	AGCGGACCCG
901	AAAAGCCTGC	CTAAGTCAA	GGCTGTGGAC	ATCGGCCCGG	GTCTGAATCT	GCTGCAAGAA	GACAACCCGG	ACCTGATCGG	CAGCGAGATC	GCGCGCTGGC
	TTTTCGGACG	GATTGACGTT	CCGACACCTG	TAGCCGGGCC	CAGACTTAGA	CGACGTTCT	CTGTTGGGCC	TGACTAGCC	GTCGCTTAG	CGCGCACCCG
1001	GTGCGACGCT	GGAGATTCC	GGCGATTGCA	TATCCAGCGC	AGTGGCGGCC	GCTCGAGTC	TAGAGGGCCC	GTTTAAAC		
	CACGCTGCGA	CCTCTAAAGG	CCGCTAAGCT	ATAGTCTGTG	TCACCGCCGG	CGAGCTCAG	ATCTCCGGG	CAAAATTTG		

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