

R&D, Administration and Europe Office

Cisbio Bioassays
 Phone: +33 (0)4 66 79 67 05
 Fax: +33 (0)4 66 79 19 20
 E-mail: bioassays@cisbio.com

Japan Office

Sceti Medical Labo K.K.
 Phone: +81 (0)3 5510 2932
 Fax: +81 (0)3 5510 0130
 E-mail: reagent@scetimedilabo.co.jp

USA Office

Cisbio US, Inc.
 Phone : 888 963 4567
 Fax : 781 687 1500
 E-mail : htrfinfo@cisbio.us



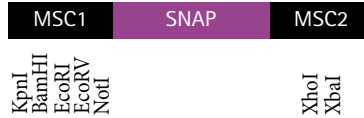
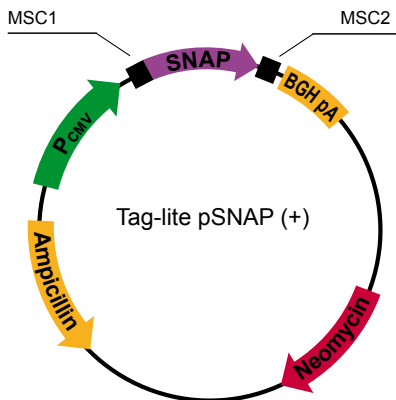
www.htrf.com

Plasmid Information

For in vitro research use only

Plasmid code: PSNPPOS rev 00 (April 2009)

Storage temperature : -20°C
 Quantity: 10 µg/10 µL H₂O



MSC1/MSC2: multiple cloning sites allowing N/C terminal fusion with the gene of interest

Plasmid name:	Tag-lite pSNAP (+)
Plasmid size:	5983bp
Tag name:	SNAP
SNAP-ORF size:	546bp

Promoter	CMV
Host cell	Mammalian
Antibiotic resistance for E. Coli	Ampicillin
Mammalian selection marker	Neomycin

SNAP-ORF sequence:

```

                                                    DpnI
                                                    ~~~~~
1  ATGGACAAG  ACTGCGAAAT  GAAGCGCACC  ACCCTGGATA  GCCCTCTGGG  CAAGCTGGAA  CTGTCTGGGT  GCGAACAGGG  CCTGCACGAG  ATCAAGCTGC
   TACCTGTTTC  TGACGCTTTA  CTTGCGGTGG  TGGGACCTAT  CGGGAGACCC  GTTCGACCTT  GACAGACCCA  CGCTTGTCCT  GGAGGTGCTC  TAGTTCGACG
101 TGGGCAAAGG  AACATCTGCC  GCCGACGCCG  TGGAAAGTGC  TGCCCCAGCC  GCCGTGCTGG  GCGGACCAGA  GCCACTGATG  CAGGCCACCG  CCTGGCTCAA
   ACCCGTTTCC  TTGTAGACGG  CGGCTGCGGC  ACCTTCACGG  ACGGGGTCCG  CGGCAGACCC  CGCTGTGCTT  CCGTGACTAC  GTCCGGTGCC  GGACCGAGTT
201 CGCCTACTTT  CACCAGCCTG  AGGCCATCGA  GGAGTTCCTC  GTGCCAGCCC  TGCACCACCC  AGTGTCCAG  CAGGAGAGCT  TTACCCGCCA  GGTGCTGTGG
   GCGGATGAAA  GTGGTCGGAC  TCCGGTAGCT  CCTCAAGGGA  CACGGTCGGG  ACGTGGTGGG  TCACAAGGTC  GTCTCTCGA  AATGGCGCGT  CCACGACACC
301 AAACCTGCTG  AAGTGGTGAA  GTTCGGAGAG  GTCATCAGCT  ACCAGCAGCT  GGCCGCCCTG  GCCGGCAATC  CCGCCGCCAC  GCCCGCCCTG  AAAACCCGCC
   TTTGACGACT  TTCACCATT  CAAGCCTCTC  CAGTAGTCGA  TGGTCTGCGA  CCGGCCGGAC  CGGCCGTTAG  GCGCGCGGTG  GCGCGCGCAC  TTTTGGCGGG
   ~~~~~
   DpnI
   ~~~~~
401 TGAGCGGAAA  TCCCGTGCCC  ATTCTGATCC  CCTGCCACCG  GGTGGTGTCT  AGCTCTGGCG  CCGTGGGGGG  CTACGAGGGC  GGGCTCGCCG  TGAAAGAGTG
   ACTCGCTTTT  AGGCACGGG  TAAGACTAGG  GGACGGTGGC  CCACCACAGA  TCGAGACCGC  GGCACCCCCC  GATGCTCCCG  CCCGAGCGCG  ACTTCTCAC
501 GCTGTGGGCC  CACGAGGGCC  ACAGACTGGG  CAAGCTTGGG  CTGGGC
   CGACGACCGG  GTGCTCCCG  TGTCTGACCC  GTTCGGACCC  GACCCG
    
```

Full sequence (ORF + plasmid):

Information available on: www.htrf.com

File under a txt format can be downloaded from the Tag-lite section of the website.

MSDS:

Information available on: www.htrf.com

Copyright © 2009 CIS bio international, France

HTRF®, TRACE®, Tag-lite, and the HTRF® logos are trademarks belonging to CIS bio international.
 HTRF® products are manufactured under one or more of the following patents and foreign equivalent : EP 0 180 492 / US 4,927,923 / US 5,220,012 / US 5,432,101 – EP 0 321 353 / US 5,457,185 / US 5,534,622 / US 5,346,996 / US 5,162,508 – EP 0 539 477 / US 5,512,493 – EP 0 539 435 / US 5,627,074 – EP 0 569 496 / US 5,527,684.

CIS bio international hereby grants to those buying HTRF® products from CIS bio international or its affiliates / distributors, a worldwide, non-exclusive, royalty-free, limited license to use HTRF® technology with said products for in-house life science research only. Signal amplification and correction using HTRF® technology are covered by the following US, patents or patent applications and foreign equivalents : US 5,512,493 – US 5,527,684 – US 6,352,672.

SNAP-tag and CLIP-tag are trademarks exclusively licensed to New England Biolabs. Plasmids encoding SNAP-tag and CLIP-tag and their substrates are made under license from New England Biolabs and may be covered by one or more of the following published patent applications filed internationally in US, EP and JP and also in additional jurisdictions : WO2002/083937 WO2004/031405 WO2004/104588 WO2004/031404 WO2005/085470 WO2008/012296 US, WO2008/122604 and EP patents 1410023 and 1696234.

The CMV promoter is covered under US. Patents 5,168,062 and 5,385,839 and its use is permitted for research purposes only. Any other use of the CMV promoter requires a license from the University of Iowa Research Foundation, 214 Technology Innovation Center, Iowa City, IA 52242.

