

**Japan Office**

**USA Office**

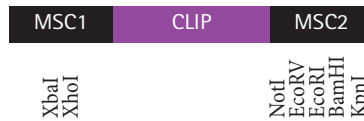
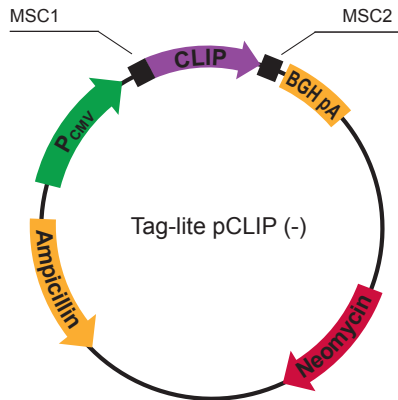
www.htrf.com

## Plasmid Information

For in vitro research use only

Plasmid code: PCLPNEG rev 00 (April 2009)

Storage temperature : -20°C  
 Quantity: 10 µg/10 µL H<sub>2</sub>O



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

Plasmid name:	Tag-lite pCLIP (-)
Plasmid size:	5982bp
Tag name:	CLIP
CLIP-ORF size:	549bp

<b>Promoter</b>	CMV
<b>Host cell</b>	Mammalian
<b>Antibiotic resistance for E. Coli</b>	Ampicillin
<b>Mammalian selection marker</b>	Neomycin

### CLIP-ORF sequence:

```

          NcoI
          ~~~~~
KpnI
~~~~~
1  ACCATGGACA AAGACTGCGA AATGAAGCGC ACCACCCCTGG ATAGCCCTCT GGGCAAGCTG GAAGTGTCTG GGTGCGAACA GGCCTGCAC GAGATCATCT
   TGGTACCTGT TTCTGACGCT TTACTTCGCG TGGTGGGACC TATCGGGAGA CCCGTTCGAC CTTGACAGAC CCACGCTTGT CCCGACGCTG CTCTAGTAGA

101 TCCTGGGCAA AGGAACATCT GCCGCCGACG CGTGGAAGT GCCTGCCCA GCGCCGTGC TGGGCGGACC AGAGCCACTG ATCCAGGCCA CCGCTGGCT
   AGGACCCGTT TCCTTGTAGA CGGCGGCTGC GGCACCTTCA CGGACGGGGT CGGCGGCACG ACCCGCCTGG TCTCGGTGAC TAGGTCCGGT GCGGGACCGA

201 CAACGCCTAC TTTACCAGC CTGAGGCCAT CGAGGAGTTC CCTGTGCCAG CCCTGCACCA CCCAGTGTTT CAGCAGGAGA GCTTTACCCG CCAGGTGCTG
   GTTGCGGATG AAAGTGGTGC GACTCCGGTA GCTCCTCAAG GGACACGGTC GGGACGTGGT GGGTACACAAG GTCGTCTCT CGAAATGGGC GGTCCACGAC

301 TGAAACTGC TGAAAGTGGT GAAGTTCGGA GAGGTACATCA GCGAGAGCCA CTTGGCCGCC CTGGTGGGCA ATCCCGCCGC CACCGCCGCC GTGAACACCG
   ACCTTTGACG ACTTTCACCA CTTCAGCCTC CTCCAGTAGT CGTCTCCGTT GGACCGGCGG GACCACCCGT TAGGGCGGCG GTGGCGGCGG CACTTGTGGC

          DpnI
          ~~~~~
401 CCCTGGACGG AAATCCCCTG CCCATTCTGA TCCCTGCCA CCGGGTGGTG CAGGGGCGACA GCGACGTGGG GCCTACCTG GCGGGGCTCG CCGTGAAAGA
   GGGACCTGCC TTAGGGGAC GGTAAGACT AGGGACCGGT GGCCACCAC GTCCCGCTGT CGTGCACCC CGGGATGGAC CCGCCGAGC GGCACCTTCT

          KpnI
          ~~~~~
501 GTGGTGTCTG GCCACGAGG GCCACAGACT GGGCAAGCCT GGGCTGGGT
   CACCGACGAC CGGGTCTCT CGGTGTCTGA CCCGTTCCGA CCCGACCCA
    
```

### Full sequence (ORF + plasmid):

Information available on: [www.htrf.com](http://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](http://www.htrf.com).