



Post Office Box 2746 Pinetop, Arizona 85935
Tel (928) 367-1200 Fax (928) 367-1205
www.nanolight.com

pUC19-Rm-GFP

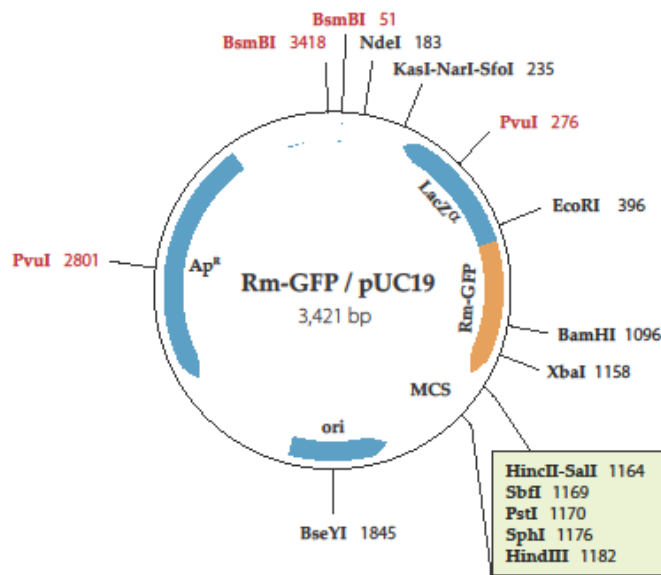
Renilla mullerie GFP - coding sequence in green, ATG start codon in **bold**
Key restriction enzymes are underlined
PUC19 sequences in blue

```
GAGCGGATAACAATTTACACAGGAAACAGCTATGACCATGATTACGCCAAGCTTGCATGCCTGCAGGTCGACTCTAGAT
TAGTGATGGTGATGGTGATGAACCCATTTCGTGTAAGGATCCTAGTGGTTTTCTATAGATGTCATTTGAGCAATAGCAGT
CTCATGCTGTTCAACGAACCCCGTCTTCTACGTAAGTCTTTTCCAAACGATGTTGAATAAAATGATACGAAGGAAACT
CCTTTACTACACCTTTTCGACTTCATTAATGTTTTTCATGTGACATGAATAATATTTCCAGAGTTTGTATATAGACAAGA
ATTACTTCGCCGACCAAGACGCCATTATTCATGTACATGGCTTCAAATGAAGGCTCTATTCCCTAAGATAGTCTTCTGCAT
GACGGGACCATCATCCGGGAAGTTACTACCTTTGTATTCCACTCTGTAGACGAACCTTGTCTTCTATTAATTTATATCTG
AACGAATTTCAACAAGTCCGCCATCTTCGTAACGTAATGTTTCGTTTCATACATAAATCCTGCTGGAAATGATTGTATAAAA
TAATCTGATATATCATTTCGGATATTTTCGTGAAAGTACGGTTGCCATATTGAAAAGCTGGTGACACAATATCAAATGCAA
AGGCAGTGGGGCCCTTTTCGTGACACGAATCTGAACCAGTTGATTGCCGAATAAAATATTCCCTTTGCCGCAACCCTCCA
TTGTAAAAACATGTTTGTAAACAATTCCTTCCAGATTTACTTTTATACGACATTACTTCTTGTAAACAAGTGTTCCTCAAT
ATTTGTTTACTCATTTTGAATTCACTGGCCGTCGTTTTACAACGTCGTGACTGGGAAAACCTGGCGTTACCCAATTAA
TCGCCTTGCAGCACATCCCCCTTTTCGCCAGCTGGCGTAATAGCG
```

Coding Sequence:

```
814 atgagtaaacaatatgaagaacacttggtttacaagaagtaatg
    M S K Q I L K N T C L Q E V M
769 tcgtataaagtaaactctggaaggaattgtaacaaccatgTTTT
    S Y K V N L E G I V N N H V F
724 acaatggagggttgccggcaaaggggaatTTTTattcggcaatcaa
    T M E G C G K G N I L F G N Q
679 ctggttcagattcgtgtcacgaaaggggccccactgccttttgca
    L V Q I R V T K G A P L P F A
634 tttgatattgtgtcaccagcttttcaatatggcaaccgtactttc
    F D I V S P A F Q Y G N R T F
589 acgaaatatccgaatgatatatcagattTTTTatacaatcattt
    T K Y P N D I S D Y F I Q S F
544 ccagcaggatttatgtatgaacgaacattacgttacgaagatggc
    P A G F M Y E R T L R Y E D G
499 ggacttgttgaaattcgttcagatataaatttaatagaagacaag
    G L V E I R S D I N L I E D K
454 ttcgtctacagagtggaatacaaaggtagtaacttcccggatgat
    F V Y R V E Y K G S N F P D D
409 ggtcccgtcatgcagaagactatcttaggaatagagccttcattt
    G P V M Q K T I L G I E P S F
364 gaagccatgtacatgaataatggcgtcttggtcggcgaagtaatt
    E A M Y M N N G V L V G E V I
319 cttgtctataaactaaactctgggaaatattattcatgtcacatg
    L V Y K L N S G K Y Y S C H M
```

274 aaaacattaatgaagtcgaaaggtgtagtaaaggagtttccttcg
 K T L M K S K G V V K E F P S
 229 tatcattttattcaacatcgtttggaaaagacttacgtagaagac
 Y H F I Q H R L E K T Y V E D
 184 gggggggttcggtgaacagcatgagactgctattgctcaaagaca
 G G F V E Q H E T A I A Q M T
 139 tctataggaaaaccactaggatccttacacgaatgggttcacac
 S I G K P L G S L H E W V H H
 94 catcaccatcactaa 80
 H H H H *



Red colored restriction sites: Multiple cutters
 Black colored restriction sites: Single cutters