



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

pGLUC Basic-1

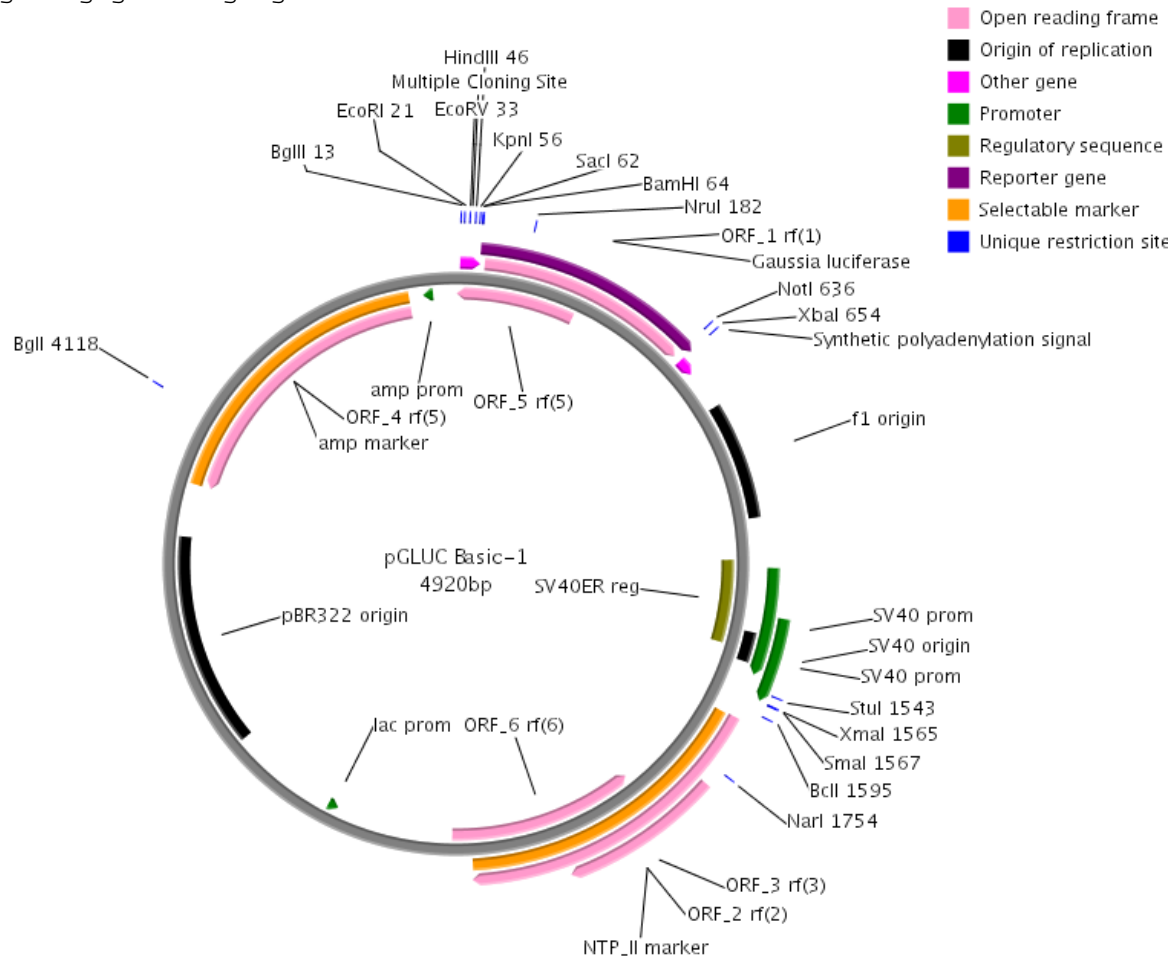
A promoterless vector with a MCS site upstream of the humanized Gaussia luciferase coding sequence (with secretion signal).

Size- 4920 bases

```
gacggatcgggagatcttgggaattctgcagatatcctcgagcccaagctt
ggtaccgagctcggatccagccaccatgggaggtcaaagttctgtttgcc
tgatctgcatcgtgtggccgaggccaagcccaccgagaacaacgaagac
ttcaacatcgtggcgtggccagcaacttcgacgaccacggatctcgatgc
tgaccgcggaagtgtgccggcaagaagctgccgctggaggtgctcaaag
agatggaagccaatgcccggaaagctggctgcaccaggggctgtctgatc
tgctgtcccacatcaagtgcacgcccagatgaagaagttcatcccagg
acgctgccacacctacgaaggcgacaaagagtcgacacagggcggcatag
gagggcgatcgtcgacattcctgagattcctgggttcaaggacttggag
cccatggagcagttcatcgacaggtcgatctgtgtgtggactgcacaac
tggctgcctcaaagggcttgccaacgtgcagtgctctgacctgctcaaga
agtggctgccgcaacgctgtgacaccttgccagcaagatccagggccag
gtggacaagatcaagggggcgggtggtgactaagcggccgctcgagcatg
catctagaaataattcttactgtcatgccaaagtaagatgcttttctgtgc
tgcaatagcaggcatgctggggatgctgggtgggtctatggcttctgaggc
ggaaagaaccagctggggctctagggggatccccacgcgccctgtagcg
gcgcattaagcgcggcgggtgtgggtggttacgcgcagcgtgaccgctaca
cttgccagcgccttagcgcggcctcttctcgcttcttcccttcttctct
cgccacgttcgcccggcttccccgtcaagctctaaatcggggcatccctt
tagggttccgatttagtgctttacggcacctcgacccccaaaaaacttgat
tagggtaggggttacgtagtgggcatcgccctgatagacgggttttctg
ccctttgacgttggagtcacgcttcttaatagtggaactctgttccaaa
ctggaacaacactcaaccctatctcggctctattcttttgatttataaggg
attttggggatttcggcctattggttaaaaaatgagctgatttaacaaaa
atthaacgcgaattaattctgtggaatgtgtgtcagttaggggtgtggaaa
gtccccaggctccccaggcaggcagaagtatgcaaagcatgcatctcaat
tagtcagcaaccaggtgtggaaagtccccaggctccccagcaggcagaag
tatgcaaagcatgcatctcaattagtcagcaaccatagtcgccccctaa
ctccgcccctcccccccctaactccgcccagttccgcccattctccgccc
catggctgactaattttttttatgatgagaggccgaggccgctctgc
ctctgagctattccagaagtagtgaggaggcttttttggaggcctaggct
tttgcaaaaagctcccgggagcttgtatatccattttcggatctgatcaa
gagacaggatgaggatcgtttcgcgatgattgaacaagatggattgcacgc
aggttctccggccgcttgggtggagaggctattcggctatgactgggcac
aacagacaatcggctgctctgatgccgcccgtgttccggctgtcagcgcag
gggcgcccgggttctttttgtcaagaccgacctgtccgggtgccctgaatga
actgcaggacgaggcagcgcggctatcgtggctggccacgacgggcttc
cttgccgagctgtgctcgacgttgtcactgaagcgggaagggactggctg
ctattgggcgaagtgcggggcaggatctcctgtcatctcaccttgctcc
tgccgagaaaagtatccatcatggctgatgcaatgcccggctgcatacgc
```

ttgatccggctacctgccattcgaccaccaagcgaacatcgcatcgag
cgagcacgtactcggatggaagccggctctgtcgatcaggatgatctgga
cgaagagcatcaggggctcgcgccagccgaactgttcgccaggctcaagg
cgcgcatgcccgcagcggcaggatctcgtcgtgacccatggcgatgcctgc
ttgccgaatatcatgggtggaaaatggccgcttttctggattcatcgactg
tggccggctgggtgtggcggaccgctatcaggacatagcgttggctaccc
gtgatattgctgaagagcttggcggcgaatgggctgaccgcttctcgtg
ctttacggatcgcgcctcccgattcgcagcgcacgccttctatcgct
tcttgacgagttcttctgagcgggactctgggggttcgaaatgaccgacca
agcgacgcccacctgccatcacgagatctcgattccaccgccccttct
atgaaaggttgggcttcggaatcgtttccgggacgcggctggatgatc
ctccagcgcggggatctcatgctggagttcttcgccaccccaacttgtt
tattgcagcttataatgggtacaaaataagcaatagcatcacaatttca
caataaagcattttttcactgcattctagttgtgggttgtccaaactc
atcaatgtatcttatcatgtctgtataccgtcgacctctagctagagctt
ggcgtaatcatggcatagctgttctcgtgtgaaattgttatccgctca
caattccacacaacatacagagccggaagcataaagtgtaaagcctgggg
gcctaagttagtgagctaaactcacattaattgcttgcgctcactgccgc
tttccagtcgggaaacctgtcgtgccagctgcattaatgaatcggccaac
gcgcggggagagggcgggttgcgctattgggcgctcttcgcttctcgtc
actgactcgtcgcctcggctcgttcggctgcccgcagcggatcagctca
ctcaaaggcggtaatacgggtatccacagaatcaggggataacgcaggaa
agaacatgtgagcaaaaggccagcaaaaggccaggaaccgtaaaaaggcc
gcgttgcctggcgtttttccataggctccgccccctgacgagcatcacia
aaatcgacgctcaagtcaaggtggcgaaacccgacaggactataaagat
accaggcgtttccccctggaagctccctcgtgcgctctcctgttccgacc
ctgccgcttaccggatacctgtccgcctttctcccttcgggaagcgtggc
gctttctcaatgctcacgctgttaggtatctcagttcgggtgtaggctcgttc
gctccaagctgggctgtgtgcacgaacccccgcttcagcccgaccgctgc
gccttatccggtaactatcgtcttgagtcacaaccggtaagacacgactt
atcgccactggcagcagcactggtaacaggattagcagagcagaggtatg
taggcggtctacagagttcttgaagtgggtggcctaactacggctacact
agaaggacagtatctgggtatctgcgctctgctgaagccagttaccttcgg
aaaaagagttggtagctcttgatccggcaaacaaaccaccgctggtagcg
gtgggttttttggtttgcaagcagcagattacgcgcagaaaaaaaggatct
caagaagatcctttgatctttctacgggggtctgacgctcagtggaacga
aaactcacgttaagggatcttgggtcatgagattatcaaaaaggatcttca
cctagatccttttaaattaaaaatgaagttttaaataatctaaagtata
tatgagtaaacttggctcagcagttaccaatgcttaatcagtgaggcacc
tatctcagcagatctgtctatctcgttcatccatagttgcctgactccccg
tcgtgtagataactacgatacgggagggcttaccatctggccccagtgct
gcaatgataccgcgagacccacgctcaccggctccagatttatcagcaat
aaaccagccagccggaagggccgagcgcagaagtggctcctgcaactttat
ccgcctccatccagctctattaattgttgcgggaagctagagtaagtagt
tcgccagttaatagtttgcgcaacgcttgttgccattgctacaggcatcgt
gggtgcacgctcgtcgtttgggtatggcttcatcagctccgggttcccaac
gatcaaggcgagttacatgatccccatgttgtgcaaaaaagcggtagc
tcttccggctcctccgactcgttgtcagaagtaagtggccgcagtggtatc
actcatgggtatggcagcactgcataattctcttactgtcatgccatccg
taagatgctttctgtgactgggtgagtaactcaaccaagtcattctgagaa
tagtgatgcggcgaccgagttgctcttgcccggcgtcaatacgggataa
taccgcgccacatagcagaactttaaagtgtcatcattggaaaacgctt
cttcggggcgaaaactctcaaggatcttaccgctgttgagatccagttcg
atgtaaccactcgtgcacccaactgatcttcagcatcttttactttcac
cagcgtttctgggtgagcaaaaacaggaaggcaaaatgccgcaaaaaagg
gaataagggcgacacggaaatgttgaatactcatactcttctttttcaa
tattattgaagcatttatcaggggtattgtctcatgagcggatacatatt

tgaatgtatttagaaaaataaacaataggggttccgcgcacatttcccc
 gaaaagtgccacctgacgtc



Features of pGLUC Basic-1 (4920 bp)

- T7 promoter bases: 20-38
- Polylinker bases: 12-63
- Gaussia luciferase reporter gene: 63-653
- SP6 promoter: 669-686
- Synthetic polyadenylation site: 653-697
- SV40 promoter bases: 1253-1573
- SV40 origin of replication: bases 1352-1437
- Neomycin ORF : bases 1609-2403
- SV40 PolyA: bases 2458-2830
- ColE1 origin: bases 3090-3763, Ampicillin ORF: bases 3908-4768

The plasmid map shown above was constructed with the help of the following link:
http://wishart.biology.ualberta.ca/PlasMapper/jsp/displayPlasmidMap.jsp?fileName=plasMap115_1101977745436.jpg&fileFormat=jpg
[Xiaoli Dong, Paul Stothard, Ian J. Forsythe, and David S. Wishart "PlasMapper: a web server for drawing and auto-annotating plasmid maps" Nucleic Acids Res. 2004 Jul 1;32\(Web Server issue\):W660-4.](#)

The CMV promoter is covered by U.S. 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

Gaussia luciferase is covered by US Patent # 6,232, 107 and IPO patents issued to Prolume Inc. Pittsburgh, PA.

This plasmid is being sold for research purposes only. Commercial use will require licenses from Prolume for the Gaussia gene and from the Univ. of Iowa Research foundation for use of the CMV promoter.