



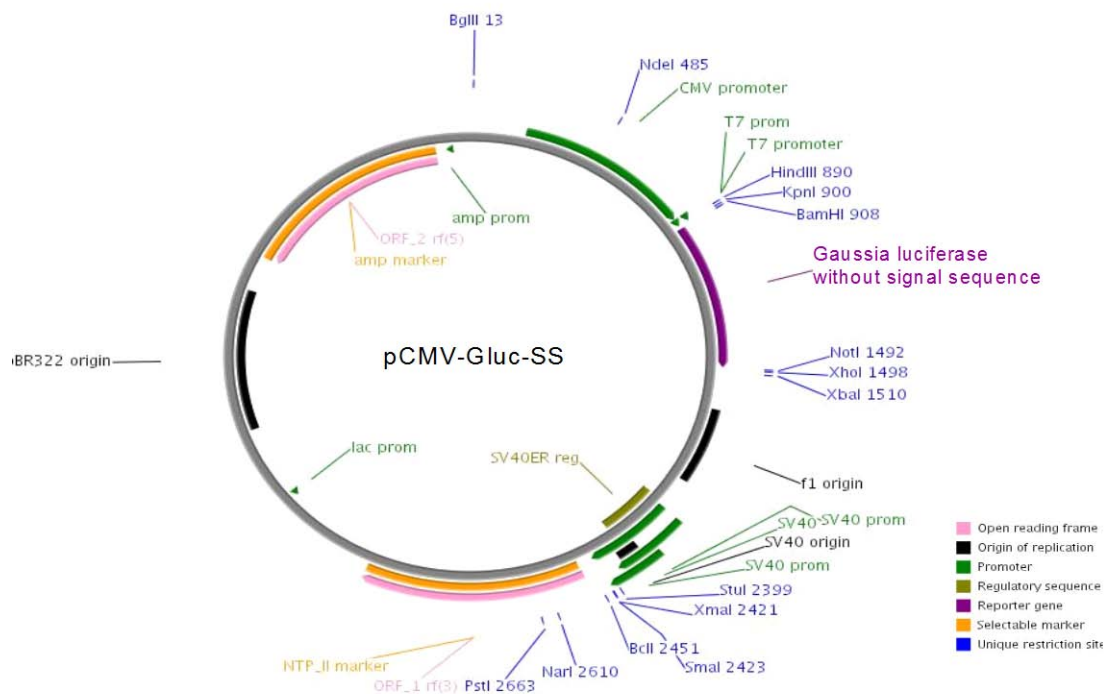
Post Office Box 2746 Pinetop, Arizona 85935  
Tel (928) 367-1200 Fax (928) 367-1205  
[www.nanolight.com](http://www.nanolight.com)

**Sequence, plasmid map and features of pCMV-GLUC without signal sequence vector**  
**Size 5710 bases,**

gacggatcgggagatctcccgatcccctatggtcgactctcagtacaatctgctctgatgccgcatagttaagccagtatct  
gctccctgcttgtgtgttggaggctcgctgagtagtgcgcgagcaaaatttaagctacaacaaggcaaggcttgaccgacaat  
tgcatgaagaatctgcttagggttaggcgttttgcgctgcttcgcgatgtacgggcccagataacgcgcttgacattgattat  
tgactagttattaatagtaatcaattacggggctcattagttcatagcccataataggagtccgcgcttacataacttacggt  
aaatggccccgctggctgaccgcccacgacccccgcccattgacgtcaataatgacgtatgttcccataagtaacgccaata  
gggactttccattgacgtcaatgggtggactatttacggtaaaactgccacttggcagtacatcaagtgtatcatatgccaa  
gtacgccccctattgacgtcaatgacggtaaatggccccgctggcattatgccagtacatgaccttatgggactttcctac  
ttggcagtacatctacgtatttagtcatcgctattaccatggtgatgcggttttggcagtacatcaatggcggtggatagcgg  
tttgactcacggggatttccaagtctccacccattgacgtcaatgggagtttgttttggcaccaaaatcaacgggactttc  
caaaatgtcgtaaactccgccccattgacgcaaatggcggttaggcgtgtacggtgggaggtctataaagcagagctct  
ctggctaactagagaaccactgcttactggcttatcgaatataatagcactcactataggagaccgaagcttggtaaccga  
gctcggatccaaatgaagcccaccgagaacaacgaagacttcaacatcgtggcgtggccagcaacttcgacgaccacggatc  
tcgatgctgaccgcggaagtggccggcaagaagctgccgctggaggtgctcaaagagatggaagccaatgcccgaaagc  
tggctgcaccaggggctgtctgatctgctgtcccacatcaagtgcacgcccgaagatgaagaagttcatcccaggacgctgc  
cacacctacgaaggcgacaaaagagtcgacacagggcgcataggcgaggcgatcgtcgacattcctgagattcctgggttca  
aggacttggagccatggagcagttcatcgcacaggtcgatctgtgtgtggactgcacaactggctcctcaagggttgc  
caactgtcagtgcttctgacctgctcaagaagtggctgcccgaacgctgtgagcactttgcccagcaagatccaggccaggtg  
gacaagatcaagggggcgggtggtagttagtaatgctcagcagcactctagaaataatttcttactgctcatgccaagtaag  
atgcttttctgtgctgcaatagcaggtatgctggggtgctgggtgctctatggcttctgaggcggaaaagaaccagctgggg  
ctctagggggtatccccacgcccctgtagcggcgcataaagcggcggggtgtggtggttacgcgcagcgtgaccgctaca  
cttggcagcgccttagcgcgcctcttctgctttcttcccttcttctctgcccagcttgcgggctttccccgtcaagctc  
taaactcggggcatcccttaggggtccgatttagtgctttacggcacctcgacccccaaaaaacttgattaggggtgatggttc  
acgtagtgggcatcgccctgatagacgggttttgcctttgacgttggagtcacgcttctttaaagtgaggactctgttc  
caactggaacaactcaaccctatctcggctctattcttttgatttataaagggttttggggatttccgctattggttaa  
aaaatgagctgatttaaaaaaatttaacgcgaattaatctgtggaatgtgtgtcagttagggtgtggaaagtcccaggc  
tcccaggcaggcagaagatgcaaagcatgcatctcaattagtcagcaaccaggtgtggaagtccccaggctcccagca  
ggcagaagatgcaaagcatgcatctcaattagtcagcaaccatagtcggccctaaactccgcccataactc  
cgcccagttccgcccattctccgccccatggctgactaatttttttatgagagggcggagggcgcctctgctctga  
gctattccagaagtagtgaggaggcttttggaggcctaggctttgcaaaaagctcccgggagcttgatataccattttc  
ggatctgatcaagagacaggatgaggatcgtttcgcatgattgaacaagatggattgacgcaggttctccggccgcttggg  
tggagaggctattcggctatgactgggcaacaacagacaatcggctgctctgatgccgcccgtgtccggctgtcagcgcaggg  
gcgcccgggttcttttgtcaagaccgacctgtccgggtgcctgaatgaactgcaggacgaggcagcgcggctatcgtggctg  
gccacgacgggcttcttgcgcagctgtgctcgacgttgtcactgaagcgggaaggactggctgctattgggcaagtg  
cggggcaggatctcctgtcatctcaccttgtcctgcccagaaaagatccatcatggctgatgcaatgcggcggctgcatac  
gcttgatccggctacctgcccattcgaccaccaagcgaaacatcgcatcgagcagcagcactcggatggaagccggtctt  
gtcgatcaggatgatctggacgaagagcatcaggggtcgcgcagccgaactgttcgcccaggctcaaggcgcgcagccccg  
acggcagggatctcgtcgtgacctatggcgatgctgcttgcgcaatcatggtggaaaatggcgcctttctggttcat  
cgactgtggcgggtgggtgtggcggaccgctatcaggacatagcgttggctaccggtgatattgctgaagagcttggcggc  
gaatgggctgaccgcttctcgtgctttacgggtatcgccgctccgattcgcagcgcacgcttctatcgcttcttgacg  
agttcttctgagcgggactctgggggtcgaaatgaccgaccaagcgcagcgcaccaactgccatcacgagatttcgattccacc  
gccgcttctatgaaaggttgggcttcggaatcgttttccgggacgcggctggatgatcctccagcgcggggatctcatgc  
tggagttcttgcgccaccccaactgtttattgcagcttataatggttacaaaataaagcaatagcatcacaatttcacaaa  
taaagcattttttcactgacttctagttgtggtttgtccaaactcatcaatgtatcttatcatgtctgtataaccgctgacc  
tctagctagagcttggcgtaatcatggtcatagctgttctgctgtgaaattgttatccgctcacaattccacacaacatac  
gagccggaagcataaagtgtaaagcctggggtgcctaataagtgagtaactcacattaattgctgtgcgctcactgcccgc  
ttccagtcgggaaacctgtcgtgccagctgcatatgaatcggccaacgcgcggggagaggcgtttgctatggggcgc  
tcttccgcttctcgtcactgactcgtcgcctcggctcgttccggctgcggcagcgggtatcagctcactcaaggcgggtaa  
tacggttatccacagaatcaggggataacgcaggaagaacatgtgagcaaaaggccagcaaaaggccaggaaccgtaaaaa

ggccgcggttgctggcggtttttccataggctccgccccctgacgagcatcacaaaaatcgacgctcaagtcaaggtggcga  
aaccgcagaggactataaagataaccaggcggtttccccctggaagctccctcgtgcgctctcctgttccgaccctgccgctta  
ccggatacctgtccgcctttctcccttcgggaagcgtggcgctttctcaatgctcacgctgtaggtatctcagttcgggtga  
ggcgttcgctccaagctgggctgtgtgcacgaaccccccggtcagcccgaccgctgcgcttatccggtaactatcgtctt  
gagccaacccggtgaagacacgacttatcgccactggcagcagccactggtaacaggattagcagagcgaggtatgtaggcg  
gtgctacagagttcttgaagtgggtggcctaactacggctacactagaaggacagtatttggatctgctgctctgctgaagcc  
agttaccttcggaaaaagagtggtagctcttgatccggcaaaacaaaccaccgctggtagcgggtggttttttggtttgcaag  
cagcagattacgcgcagaaaaaaggatctcaagaagatccttggatcttttctacgggctgacgctcagtggaacgaaa  
actcacgttaagggattttggatcatgagattatcaaaaaggatcttcacctagatccttttaaattaaaaatgaagttttaa  
atcaatctaaagtatatatgagtaaaacttggctgacagttaccaatgcttaatcagtgaggcacctatctcagcgatctgt  
ctatttcgttcatccatagttgcctgactccccgctcgtgtagataactacgatacgggagggcttaccatctggccccagtg  
ctgcaatgataccgcgagaccacgctcaccggctccagatttatcagcaataaaccagccagccggaagggccgagcgcag  
aagtggctcctgcaactttatccgcctccatccagtctattaattggtgcccgggaagctagagtaagtagttcggcagttaat  
agtttgcgcaacggtgttgccattgctacaggcatcgtgggtgcacgctcgtcgtttggatggcttcattcagctccgggtt  
ccaacgatcaaggcgagttacatgatccccatggttgcaaaaaagcgggttagctccttcggctcctccgatcgttgcag  
aagtaagtggccgcagtggtatcactcatggttatggcagcactgcataattctcttactgtcatgccatccgtaagatgc  
tttctgtgactggtgagtactcaaccaagtcattctgagaatagtgtatgcggcgaccgagtgctcttgcccggcgtcaa  
tacgggataataccgcgccacatagcagaactttaaagtgtcatcattggaaaacgttcttcggggcgaactctcaag  
gatcttaccgctgttgagatccagttcgatgtaaccactcgtgcacccaactgatcttcagcatctttactttcaccagc  
gtttctgggtgagcaaaaacaggaaggcaaaatgccgcaaaaaaggggaataagggcgacacggaaatgttgaatactcatac  
tcttctttttcaatattattgaagcatttatcaggggtattgtctcatgagcggatacatatttgaatgtatttagaaaaa  
taaacaaataggggttccgcgcacatttccccgaaaagtgccacctgacgctc

## Plasmid Map of pCMV-Gluc minus SS (5710 bp)



### Features:

pCMV-Gluc without the signal sequence (pCMV-Gluc minus SS; 5710 bp)

CMV promoter bases: 209-863

Gaussia luciferase gene without the signal peptide: 915-1424

T7 promoter bases: 864-882

Polylinker bases: 889-907

Synthetic polyadenylation site: 1443-1487

SV40 promoter bases: 2028-2363

SV40 origin of replication: bases 2142-2227

Neomycin ORF : bases 2399-3193

SV40 PolyA: bases 3248-3620

ColE1 origin: bases 3880-4553

Ampicillin ORF: bases 4698-5558

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](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.**

**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.**