



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

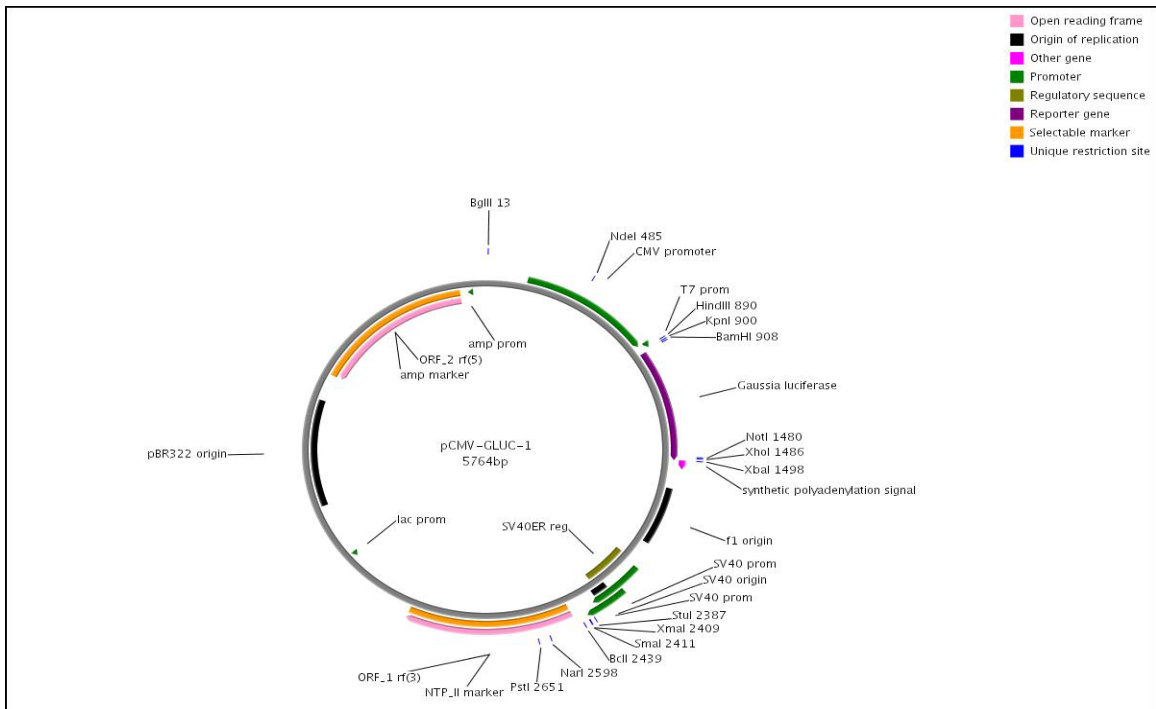
pCMV-GLUC-1

Humanized Gaussia luciferase with secretion signal

gacggatcgggagatctcccgatcccctatgggtcgactctcagtacaatc
tgctctgatgcccgcataagcagatctgctccctgcttgtgtgtt
ggaggctcgtgagtagtgccgagcaaaatctaagctacaacaaggcaag
gcttgaccgacaattgcatgaagaatctgcttaggggttaggcgctttgcg
ctgcttcgcatgtacgggcccagatatacgcgcttgacattgattattgac
tagttattaatagtaatacaattacgggggtcattagttcatagcccatata
tggagttccgcttacataacttacggtaaatggcccgcctggctgaccg
cccaacgacccccgcccattgacgtcaataatgacgtatgtcccattag
aacgccaatagggactttccattgacgtcaatgggtggactatttacgg
aaactgcccacttggcagtagcatcaagtgtatcatatgccaagtacgccc
cctattgacgtcaatgacggtaaatggcccgcctggcattatgcccagta
catgaccttatgggactttcctacttggcagtagcatctacgtattagtca
tcgctattaccatgggtgatgcccgttttggcagtagcatcaatgggctgga
tagcgggttgactcacggggatttccaagtctccacccccattgacgtcaa
tgggagttgttttggcaccaaaatcaacgggactttccaaaatgtcgta
acaactccgccccattgacgcaaatgggcccgtaggcgtgtacgggtgggag
gtctatataagcagagctctctggctaactagagaaccactgcttactg
gcttatcgaaattaatacagactcactataggagaccgaagcttggtagc
gagctcggatccagccaccatgggagtagcaagttctgtttgccctgatct
gcatcgtctgtggccgaggccaagcccaccgagaacaacgaagacttcaac
atcgtggccgtggccagcaacttcgacgaccacggatctcgatgctgaccg
cgggaagtggccggcaagaagctgcccgtggagggtgctcaaagagatgg
aagccaatgcccggaaagctggctgcaccaggggctgtctgatctgcctg
tcccacatcaagtgcacgcccagaatgaagaagttcatcccaggacgctg
ccacacctacgaaggcagacaagaagctccgcacagggcggcataggcaggg
cgatcgtcgacattcctgagattcctgggttcaaggacttggagcccatg
gagcagttcatcgacaggtcgatctgtgtgtggactgcacaactggctg
cctcaaagggttggcaactgacgtgttctgacctgctcaagaagtggc
tgccgcaacgctgtgacgacctttgccagcaagatccagggccaggtggac
aagatcaagggggcgggtggtagtaagcggccgctcgagcatgcatcta
gaaataattcttactgtcatgccaagtaagatgcttttctgtgctgcaat
agcaggcatgctggggatgcccgtgggtctatggcttctgagggcggaaag
aaccagctggggctctagggggatccccacgcgcccgttagcggcgcac
taagcgcggcgggtgtgggttacgcgcagcgtgaccgctacacttgcc
agcgccttagcgcggcctccttctgctttcttcccttcccttctcgccac
gttcgcccggctttcccgtcaagctctaaatcggggcatccctttagggt
tccgatttagtgctttacggcacctcgacccccaaaaaacttgattagggt
gatggttcacgtagtgggccatcgccctgatagacgggtttttcgccctt
gacgttggagtagcaggttctttaaagtggactcttgttccaaactggaa
caacactcaaccctatctcggtctattcttttgatttataagggattttg
gggatttcggcctattgggttaaaaaatgagctgatttaacaaaaattaa
cgcgaattaattctgtggaatgtgtgtcagttagggtgtggaaagtccc
aggctcccaggcaggcagaagtagcaaaagcatgcatctcaattagtca
gcaaccaggtgtggaaagtcccaggctcccagcaggcagaagtagca
aagcatgcatctcaattagtcagcaaccatagctcccggccctaactccgc

ccatcccgccttaactccgcccagttccgcccatttctccgccccatggc
tgactaattttttttatttatgcagaggccgaggccgcctctgcctctga
gctattccagaagtagtgaggaggcttttttggaggcctaggcttttgca
aaaagctcccgggagcttgtatatccattttcggatctgatcaagagaca
ggatgaggatcgtttcgcatgattgaacaagatggattgcacgcaggttc
tccggccgcttgggtggagaggctattcggctatgactgggcacaacaga
caatcggctgctctgatgccgccgtgttccggctgtcagcgcagggggc
ccggttctttttgtcaagaccgacctgtccgggtgccctgaatgaactgca
ggacgaggcagcgcggctatcgtggctggccacgacggggcgttccctgcg
cagctgtgctcgacgttgtcactgaagcgggaagggactggctgctattg
ggcgaagtgcggggcaggatctcctgtcatctcaccttgctcctgcccga
gaaagtatccatcatggctgatgcaatgcggcggctgcatacgttgatc
cggctacctgcccattcgaccaccaagcgaacatcgcatcgagcgagca
cgtactcggatcgaagcgggtcttgtcgatcaggatgatctggacgaaga
gcatcaggggctcgcccgaccgaactgttccgaggctcaaggcgcgca
tgcccgacggcgaggatctcgtcgtgacccatggcgatgctgcttgccg
aatacatggtggaaaatggccgcttttctggattcatcgactgtggccg
gctgggtgtggcggaccgctatcaggacatagcgttggctaccctgata
ttgctgaagagcttggcggcgaatgggctgaccgcttccctcgtgctttac
ggatcgccgctcccgatcgcagcgcacgccttctatcgcttcttga
cgagttcttctgagcgggactctgggggttcgaaatgaccgaccaagcgac
gccaacctgcccacacgagatctcgattccaccgccccttctatgaaa
ggttgggcttcggaatcgttttccgggacgcccggctggatgatcctccag
cgcggggatctcatgctggagttcttcgcccaccccaacttgtttattgc
agcttataatggttacaataaagcaatagcatcacaatttcacaaata
aagcattttttcactgcattctagttgtggtttgtccaaactcatcaat
gtatcttatcatgtctgtataccgctcgacctctagctagagcttggcgta
atcatggtcatagctgttctcctgtgtgaaattgttatccgctcacaattc
cacacaacatacagcgcgggaagcataaaagtgtaaagcctgggggtgcctaa
tgagtgaactaactcacattaattgcgttgcgctcactgcccgccttcca
gtcgggaaacctgtcgtgcccagctgcattaatgaatcggccaacgcgcgg
ggagaggcgggttgcgtattgggcgctcttccgcttccctcgtcactgac
tcgctgcgctcggctcgttccgctgcccgcgagcggatcagctcactcaaa
ggcggtaatacggttatccacagaatcaggggataacgcaggaaagaaca
tgtgagcaaaaggccagcaaaaggccaggaaccgtaaaaaggccgcgcttg
ctggcggtttttccataggctccgccccctgacgagcatcacaataatcg
acgctcaagtcaaggtggcgaacccgacaggactataaagataaccagg
cgtttccccctggaagctccctcgtgcgctctcctgttccgaccctgccg
cttaccggatacctgtccgcctttctcccttcgggaagcgtggcgctttc
tcaatgctcacgctgtaggtatctcagttcgggtgtaggtcgttcgctcca
agctgggctgtgtgcacgaacccccgcttcagcccgaccgctgcccctta
tccggtaactatcgtccttgagttcaacccggtaagacacgacttatcgcc
actggcagcagccactggtaacaggattagcagagcggatgttagggc
gtgctacagagttcttgaagtggcctaactacggctacactagaagg
acagtatctggatctgcgctctgctgaagccagttaccttcggaaaaag
agttggtagctcttgatccggcaaacaaaccaccgctggtagcgggtggtt
ttttggttgcaagcagcagattacgcgcagaaaaaaaggatctcaagaa
gatcctttgatctttctacggggtctgacgctcagtggaacgaaaactc
acgttaagggatcttggatcatgagattatcaaaaaggatcttcacctaga
tccttttaattaaaaatgaagttttaaatacaatctaaagtatatatgag
taaacttggctgacagttaccaatgcttaatacagtgaggcacctatctc
agcgatctgtctatctcgttccatccatagttgcctgactccccgctcgtg
agataactacgatacgggagggttaccatctggccccagtgctgcaatg
ataccgagacccacgctcaccggctccagatttatcagcaataaacca
gccagccggaagggccgagcgcagaagtggctcctgcaactttatccgct
ccatccagctctatcaattggttgcgggaagctagagtaagtagttcgcca
ggttaatagtttgcgcaacggtgttgccattgctacaggcatcgtgggtgc

acgctcgtcgtttggtatggcttcattcagctccgggtcccaacgatcaa
ggcgagttacatgatccccatggtgtgcaaaaaagcggtagctccttc
ggtcctccgatcgttgtcagaagtaagttggccgcagtggtatcactcat
ggttatggcagcactgcataattctcttactgtcatgccatccgtaagat
gcttttctgtgactgggtgagtactcaaccaagtcattctgagaatagtgt
atgcggcgaccgagttgctcttgccggcgtcaatacgggataataccgc
gccacatagcagaactttaaaagtgtcatcattggaaaacgttcttcgg
ggcgaaaactctcaaggatcttaccgctggtgagatccagttcgatgtaa
cccactcgtgcacccaactgatcttcagcatcttttactttcaccagcgt
ttctgggtgagcaaaaacaggaaggcaaaatgccgcaaaaaaggaataa
ggcgacacggaaatggtgaatactcatactcttcctttttcaatattat
tgaagcatttatcagggttattgtctcatgagcggatacatatttgaatg
tatttagaaaaataaacaatataggggtccgcgcacatttccccgaaaag
tgccacctgacgtc



FEATURES:

pCMV-GLUC-1 (5764 bp)

CMV promoter bases: 209-863

Gaussia luciferase gene: 920-1477

T7 promoter bases: 864-882

Polylinker bases: 889-907

SP6 promoter: 1513-1530

Synthetic polyadenylation site: 1497-1541

SV40 promoter bases: 2082-2417

SV40 origin of replication: bases 2196-2281

Neomycin ORF : bases 2453-3247

SV40 PolyA: bases 3302-3674

ColE1 origin: bases 3934-4607

Ampicillin ORF: bases 4752-5612

The plasmid map shown above was constructed with the help of the following link
[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.