



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

pCMV-KDEL-GLUC-1

Humanized Gaussia luciferase with C-terminal KDEL endoplasmic reticulum retention signal for intracellular expression

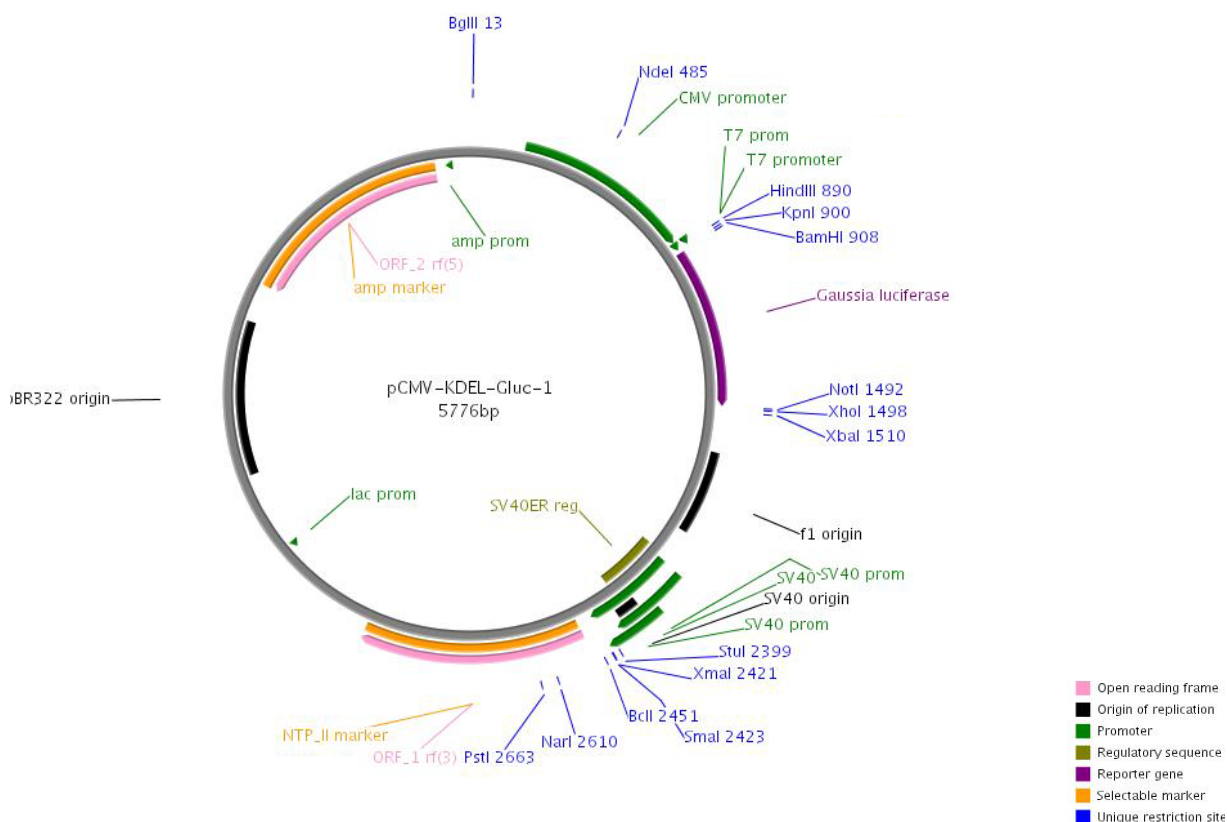
Size 5776 bases,

```
gacggatcgggagatctcccgatcccctatggctcgcactctcagtacaatc
tgctctgatgccgcatagttaagccagtatctgctccctgcttgtgtgtt
ggaggtcgcctgagtagtgcgcgagcaaaatttaagctacaacaaggcaag
gcttgaccgacaattgcatgaagaatctgcttaggggttaggcgttttgcg
ctgcttcgcgatgtacgggcccagatatacgcgcttgacattgattattgac
tagttattaatagtaatcaattacgggggtcattagttcatagcccata
tggagttccgcgttacataaacttacggtaaatggcccgcctggctgaccg
cccaacgacccccgccattgacgtcaataatgacgtatgttcccatagt
aacgccaatagggactttccattgacgtcaatgggtggactatttacggg
aaactgcccacttggcagtagacatcaagtgtatcatatgccaagtacgccc
cctattgacgtcaatgacggtaaatggcccgcctggcattatgcccagta
catgaccttatgggactttcctacttggcagtagacatctacgtattagtca
tcgctattaccatgggtgatgcggttttggcagtagacatcaatggcgctgga
tagcgggttgactcacggggatttccaagtctccacccattgacgtcaa
tgggagtttgttttggcaccaaaatcaacgggactttccaaaatgtcgtg
acaactccgccccattgacgcaaatgggcggtaggcgtgtacgggtgggag
gtctatataagcagagctctctggcctaactagagaacccactgcttactg
gcttatcgaaattaatagcactcactatagggagaccaagcttgggtacc
gagctcggatccagccaccatgggagtagcaagttctgtttgcccctgatct
gcatcgcctgtggccgagggcaagcccaccgagaacaacgaagacttcaac
atcgtggccgtggccagcaacttcgcgaccacggatctcgtatgctgaccg
cgggaagtgtcccggcaagaagctgcccgtggaggtgctcaaagagatgg
aagccaatgcccggaaagctggctgcaccaggggctgtctgatctgcctg
tcccacatcaagtgcacgcccagaatgaagaagttcatcccaggagcgtg
ccacacctacgaaggcgacaaagagtcgacacagggcgccataggcgagg
cgatcgtcgacattcctgagattcctgggttcaaggacttggagcccattg
gagcagttcatcgacaggtcgatctgtgtgtggactgcacaactggctg
cctcaaagggttgcacaacgtgcaggttctgacctgctcaagaagtggc
tgccgcaacgctgtgacgactttgcccagcaagatccagggccaggtggac
aagatcaaggggccggtgggtgacaaagacgaactgagcggccgctcgagc
atgcatctagaaataattcttactgtcatgccaagtaagatgcttttctg
tgctgcaatagcaggcatgctggggatgcggtgggctctatggcttctga
ggcggaaagaaccagctggggctctagggggtatccccacgcccctgta
gcggcgcatgaagcgcggcggtgtgggtgttacgcgcagcgtgaccgct
acacttgccagcgccttagcgcggcctcctttcgttttcttcccttctt
tctcgcacgcttcgcccgtttccccgtcaagctctaaatcggggcatcc
ctttagggttccgatttagtgctttacggcacctcgacccccaaaaaactt
gattagggtgatggttacgtagtgggcatcgccctgatagacggtttt
tcgccccttgacgttggagtccacgttcttaaatagtggtactctgttcc
aaactggaacaacactcaaccctatctcggctctattcttttgattataa
gggattttggggatttcggcctattggttaaaaaatgagctgatttaaca
aaaatttaacgcgaatttaattctgtggaatgtgtgtcagttagggtgtgg
aaagtccccaggctccccaggcaggcagaagtagcgaagcatgcatctc
aattagtcagcaaccaggtgtggaaagtccccaggctccccagcaggcag
aagtagcgaagcatgcatctcaattagtcagcaaccatagtcggcccc
taactccgcccattcccggccctaactccgcccagttccgcccattctccg
ccccatggctgactaatttttttatttatgacagaggccgaggccgcctc
```

tgctctgagctattccagaagtagtgaggaggcttttttggaggcctag
gcttttgcaaaaagctcccgggagcttgtatatccattttcggatctgat
caagagacaggatgaggatcgtttcgcatgattgaacaagatggattgca
cgcaggttctccggccgcttgggtggagaggctattcggctatgactggg
cacaacagacaatcggctgctctgatgccgccgtgtccggctgtcagcg
cagggcgcccggttctttttgtcaagaccgacctgtccggtgccctgaa
tgaactgcaggacgaggcagcgcggctatcgtggctggccacgacgggcg
ttccttgcgcagctgtgctcgcagcttgtcactgaagcgggaagggactgg
ctgctattgggcaagtgccggggcaggatctcctgtcatctcaccttgc
tcctgcccagaaaagtatccatcatggctgatgcaatgcccggctgcata
cgcttgatccggctacctgccattcgaccaccaagcgaacaatcgcatc
gagcgcagcagctactcggatggaagccggctcttgtcgatcaggatgatct
ggacgaagagcatcaggggctcgcgccagccgaactgttcgccaggctca
aggcgcgcgatgcccgacggcgaggatctcgtcgtgacccatggcgatgcc
tgcttgccgaatatcatggtggaanaatggccgcttttctggattcatcga
ctgtggccggctgggtgtggcggaccgctatcaggacatagcgttggcta
cccgtgatattgtcgaagagcttggcggcgaatgggctgaccgcttctc
gtgctttacggtatcgcgctcccgatcgcagcgcacgccttctatcg
ccttcttgacgagttcttctgagcgggactctggggttcgaaatgaccga
ccaagcgcgcccacctgccatcacgagatttcgattccaccgcccgcct
tctatgaaaggttgggcttcggaatcgttttccgggacgcccggctggatg
atcctccagcgcggggatctcatgctggagtcttctgcccccactt
gtttattgcagcttataatgggtacaaataaagcaatagcatcacaatt
tcacaaataaagcatttttttactgcattctagtgtggtttgtccaaa
ctcatcaatgtatcttatcatgtctgtataaccgtcgacctctagctagag
cttggcgtaatcatggtcatagctgtttcctgtgtgaaatgttatccgc
tcacaattccacacaacatacagagccggaagcataaaagtgtaaagcctgg
ggtgcctaagtgagtgagctaaactcacattaattgctgtgcgctactgcc
cgctttccagtcgggaaacctgtcgtgccagctgcattaatgaatcggcc
aacgcgcggggagaggcgggttgcgtattgggcgctcttccgcttctc
ctcactgactcgtcgcgctcggctcgttcggctgcggcgagcggatcagc
tactcaaaggcggtaatacgggttatccacagaatcaggggataaacgcag
gaaagaacatgtgagcaaaaggccagcaaaaggccaggaacctgaaaaag
gccgcgttgctggcgttttccataggtccgccccctgacgagcatca
caaaaatcgacgctcaagtgcagaggtggcgaaccgcagagactataaa
gataaccagcgtttccccctggaagctccctcgtgcgctctcctgttccg
acctgcccgttaccggatacctgtccgccttctcccttcgggaagcgt
ggcgttttctcaatgctcacgctgtaggtatctcagttcgggtgtaggtcg
ttcgtccaagctgggctgtgtgcacgaaccccccttccagcccgaccgc
tgcgccttatccggtaactatcgtcttgagtccaacccggtaagacacga
cttatcgccactggcagcagccactggtaacaggattagcagagcggagt
atgtaggcggtgctacagagttcttgaagtgggtggcctaactacggctac
actagaaggacagtatttggtatctgcgctctgctgaagccagttacctt
cggaaaaagagttggtagctcttgatccggcaaacaaaccaccgctggta
gcggtggttttttgtttgcaagcagcagattacgcgcagaaaaaaagga
tctcaagaagatcctttgatcttttctacggggtctgacgctcagtgga
cgaaaactcacgttaagggattttgggtcatgagattatcaaaaaggatct
tcacctagatccttttaaatataaaatgaagttttaaatcaatctaaagt
atataatgagtaaaacttggctcgcagcttaccaatgcttaatcagtgaggc
acctatctcagcgatctgtctatctcgttcatccatagttgcctgactcc
ccgtcgtgtagataaactacgatacgggagggcttaccatctggccccagt
gctgcaatgataccgcgagaccacgctcaccggctccagatttatcagc
ataaaaccagccagccggaaggccgagcgcagaagtggctcctgcaactt
tatccgcctccatccagctctaatgttgcggggaagctagagtaagt
agttcggcagttaatagtttgcgcaacggtgttgccattgctacaggcat
cgtgggtgtcacgctcgtcgtttgggtatggcttcattcagctccggttccc
aacgatcaaggcgcagttacatgatccccatgttgtgcaaaaaagcgggtt
agctccttcggctcctccgatcgttgtcagaagtaagttggccgcagtggt
atcactcatgggttatggcagcactgcataattctcttactgtcatgccat
ccgtaagatgcttttctgtgactgggtgagtactcaaccaagtcattctga
gaatagtgatgcggcaccgagttgctcttgcccggcgtcaatacggga

taataccgcgccacatagcagaactttaaaagtgctcatcattggaaaaac
 gttcttcggggcgaaaactctcaaggatcttaccgctgttgagatccagt
 tcgatgtaaccactcgtgcacccaactgatcttcagcatcttttacttt
 caccagcgtttctgggtgagcaaaaacaggaaggcaaaaatgccgcaaaaa
 agggaataagggcgacacgaaaatggtgaatactcatactcttcttttt
 caatattattgaagcatttatcagggttattgtctcatgagcggatacat
 atttgaatgtatttagaaaaataaacaataggggttccgcgcacatttc
 ccgaaaagtgccacctgacgtc

Plasmid Map of pCMV-KDEL-Gluc-1 (5776 bp)



Features:

pCMV-KDEL-Gluc-1 (5776 bp)
 CMV promoter bases: 209-863
 Gaussia luciferase gene: 920-1486
 T7 promoter bases: 864-882
 Polylinker bases: 889-907
 Synthetic polyadenylation site: 1509-1553
 SV40 promoter bases: 2094-2429
 SV40 origin of replication: bases 2208-2293
 Neomycin ORF : bases 2465-3259
 SV40 PolyA: bases 3314-3686
 ColE1 origin: bases 3946-4619
 Ampicillin ORF: bases 4764-5624

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.