



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

## **PGLUC-KDEL-Basic-1**

Humanized Gaussia Luciferase, promoterless  
with C-terminal KDEL endoplasmic retention signal

### **Size- 4932 bases**

```
gacggatcgggagatccttgggaattctgcagatatcctcgagcccaagctt
ggtaccgagctcggatccagccacc atgggagtc aaagtctgtttgccc
tgatctgcatcgctgtggccgagggccaagcccaccgagaacaacgaagac
ttcaacatcgtggccgtggccagcaacttcgcgaccacggatctcgatgc
tgaccgcggaagtggcccggaagaagctgcccgtggaggtgctcaaag
agatggaagccaatgcccggaaagctggctgcaccaggggctgtctgatc
tgcctgtcccacatcaagtgcacgcccgaagatgaagaagtcatcccagg
acgctgccacacctacgaaggcgacaaagagtcgacacagggcggcatag
gcgaggcgatcgtcgacattcctgagattcctgggttcaaggacttgag
cccatggagcagttcatcgacaggtcgatctgtgtgtggactgcacaac
tggctgacctcaaagggcttgccaacgtgcagtgttctgacctgctcaaga
agtggctgcccgaacgctgtgacgaccttgccagcaagatccagggccag
gtggacaagatcaagggggcgggtggtgacaaagcgaacttaggcggccgc
tcgagcatgcatctagaaataattcttactgtcatgccaagtaagatgct
ttctgtgctgcaatagcaggcatgctggggatgcccgtgggctctatggc
ttctgaggcggaaagaaccagctggggctctagggggtatccccacgccc
cctgtagcggcgcattaaagcggcgggtgtggtggttacgcgcagcgtg
accgctacacttgccagcgccttagcggccgctcctttcgctttcttccc
ttcctttctcgccaggttcgcccggctttccccgtcaagctctaaatcggg
gcatccctttaggggtccgatttagtgctttacggcacctcgaccccaaa
aaacttgataggggtgatggttcacgtagtgggcatcgccctgatagac
ggtttttcgccctttgacgcttgaggtccacgttctttaatagtggactct
tgttccaaactggaacaacactcaacctatctcggctctattcttttgat
ttataagggattttggggatttcggcctattgggttaaaaaatgagctgat
ttaacaaaaatttaacgcgaattaattctgtggaatgtgtgtcagttagg
gtgtggaagtccccaggctccccaggcaggcagaagtatgcaaagcatg
catctcaattagttagcaaccagggtgtggaagtccccaggctccccagc
aggcagaagtatgcaaagcatgcatctcaattagttagcaaccatagtcc
cgccccctaactccgcccataccgcccctaactccgcccagttccgcccata
tctccgccccatggctgactaattttttttatattatgcagaggccgaggc
cgctctgcctctgagctattccagaagtagtgaggaggcttttttgag
gcctaggcttttgcaaaaagctcccgggagcttgtatattccatcttcgga
tctgatcaagagacaggatgaggatcgtttcgcagatgattgaacaagatgg
attgcacgcagggttctccggccgcttgggtggagaggctattcggctatg
actgggcacaacagacaatcggtgctctgatgccgcccgtgttccggctg
tcagcgcaggggcccgggttctttttgtcaagaccgacctgtccggctgc
cctgaatgaactgcaggacgaggcagcgggctatcgtggctggccacga
cgggcttccctgcccagctgtgctcgacgttgtcactgaagcgggaagg
gactggctgctattgggcgaagtgcggggcaggatctcctgtcatctca
```

ccttgctcctgccgagaaagtatccatcatggctgatgcaatgccggcggc  
tgcatacgccttgatccggctacctgcccatccgaccaccaagcgaaacat  
cgcacgagcgcgacgactcggatggaagccggctctgtcgatcagga  
tgatctggacgaagagcatcaggggctcgcgccagccgaactgttcgcca  
ggctcaaggcgcgcatgcccagggcgaggatctcgtcgtgacccatggc  
gatgcctgcttgccgaatatcatgggtggaaaatggccgcttttctggatt  
catcgactgtggccggctgggtgtggcggaccgctatcaggacatagcgt  
tggctaccgctgatattgctgaagagcttggcggcgaatgggctgaccgc  
ttcctcgtgctttacggatcgcgcgctcccgatccgacgcgcatcgcctt  
ctatcgccttcttgacgagttcttctgagcgggactctggggttcgaaat  
gaccgaccaagcgcagcgcceaacctgccatcacgagatttcgattccaccg  
cgccttctatgaaaggttgggcttcggaatcgttttcggggacgcggc  
tggatgatcctccagcggggatctcatgctggagtcttcgcccacc  
caacttgtttattgcagcttataatgggttacaataaagcaatagcatca  
caaatttcacaaataaagcatttttttactgcattctagttgtggttg  
tccaaactcatcaatgtatcttatcatgtctgtataaccgtcgacctag  
ctagagcttggcgtaatcatgggtcatagctgtttcctgtgtgaaattgtt  
atccgctcacaattccacacaacatacagagccggaagcataaagtgtaaa  
gcctgggggtgcctaataagtgagtgagtaactcacattaattgctgtgcgctc  
actgcccgctttccagtcgggaaacctgtcgtgccagctgcattaatgaa  
tcggccaacgcgcggggagaggcgggttgcgctatggggcgtcttcgct  
tctcgtcactgactcgtcgcgctcggctcgttcggctgcggcgagcgg  
atcagctcactcaaaggcggtaatacgggttatccacagaatcaggggata  
acgcaggaagaacatgtgagcaaaaggccagcaaaaggccaggaaccgt  
aaaaaggcgcgcttgctggcgtttttccataggctccgccccctgacga  
gcatcacaaaaatcgacgctcaagttagaggtggcgaaaccgcagggac  
tataaagataaccaggcgtttccccctggaagctcctcgtgcgctctcct  
gttccgacctgcccgttaccggataacctgtccgctttctcccttcggg  
aagcgtggcgtcttctcaatgctcacgctgtaggtatctcagttcgggtg  
aggctcgttcgctccaagctgggctgtgtgcacgaacccccgttcagccc  
gaccgctcgccttatccggtaactatcgtcttgagttcaaccggtaag  
acacgacttatcgccactggcagcagccactggtaacaggattagcagag  
cgaggatgtaggcgggtgctacagagttcttgaagtgggtggcctaactac  
ggctacactagaaggacagtatcttggtatctgcgctctgctgaagccagt  
taccttcggaaaaagagttggtagctcttgatccggcaaaacaaaccg  
ctggtagcgggtgggttttttggtttgcaagcagcagattacgcgcagaaaa  
aaaggatctcaagaagatcctttgatcttttctacggggctgacgctca  
gtggaacgaaaactcacgtaagggatcttggtcatgagattatcaaaaa  
ggatcttcacctagatccttttaaatataaaatgaagtcttaaatcaatc  
taaagtatatatgagtaaaccttggctgacagttaccaatgcttaacag  
tgaggcacctatctcagcagatctgtctatctcgttcacatagttgcct  
gactccccgtcgtgtagataactacgatacgggagggcttaccatctggc  
cccagtgctgcaatgataccgcgagaccacgctcaccggctccagattt  
atcagcaataaaccagccagccggaaggccgagcgcagaagtggctctg  
caactttatccgctccatccagctctattaattggttgcggggaagctaga  
gtaagtagttcgccagttaatagtttgcgcaacgttgttgccattgctac  
aggcatcgtgggtgtcacgctcgtcgtttgggtatggcttcattcagctcgc  
gttcccaacgatcaaggcaggttacatgatccccatgttgtgcaaaaaa  
gcggttagctcctcggctcctccgatcgttgtcagaagtaagtggccgc  
agtgttatcactcatggttatggcagcactgcataattctcttactgtca  
tgccatccgtaagatgcttttctgtgactgggtgagtaactcaaccaagtca  
ttctgagaatagtgatgcccgcagcagttgctcttgcccggcgtcaat  
acgggataataccgcgccacatagcagaactttaaagtgtcatcattg  
gaaaacgcttcttcggggcgaaaactctcaaggatcttaccgctgttgaga  
tccagttcgatgtaaccactcgtgcacccaactgatcttcagcatcttt  
tactttcaccagcgtttctgggtgagcaaaaaacaggaaggcaaaatgccg  
caaaaaaggggaataagggcgacacggaaatgttgaatactcatactctc



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. 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.