



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

## **PUC18-Gluc\without\SS**

expressing native *Gussia* luciferase without the secretion signal

gcgccaatcgcgcaaacgcctctccccgcgcgttggccgattcattaatgcagctggcagcagaggtt  
cccgactggaaagcgggcagtgagcgcgaacgcaattaatgtgagtttagctcactcatttaggcaccccagg  
ctttacactttatgcttccggctcgtatgttgtgtggaattgtgagcggataacaatttcacacagggaa  
cagctatgaccatgattacgaattcgagctcggtagccggggatccatgaaaccaactgaaaacaatgaa  
gatttcaacattgtagctgtagctagcaactttgctacaacggatctcgatgctgaccgtggtaaatg  
ccggaaaaaaattaccacttgaggtactcaagaaatggaagccaatgctaggaagctggctgcactag  
gggatgtctgatgtcctgtcacacatcaagtgtacacccaaaatgaagaagttatccaggaagatgc  
cacacctatgaaggagacaaagaaagtgcacagggaggaataggagaggctattgttgacattcctgaaa  
ttcctgggtttaaggatttggaaacctggaacaattcattgcacaagttgacctatgtgtagactgcac  
aactggatgcctcaaaggtcttgccaatgtgcaatgttctgatttactcaagaaatggctgccacaaaga  
tgtgcaacttttgctagcaaaattcaaggccaagtggacaaaataaagggtgccgggtggatcatcacc  
atcaccatcactaatctagagtcgacctgcagcagatgcaagcttggcactggcctgcttttacaacgtc  
gtgactgggaaaaccctggcgttaccacaacttaatgccttgcagcacatcccccttccgacgctggcg  
taatagcgaagaggcccgaccgatcgcccttccaacagttgcgagcctgaatggcgaatggcgctg  
atgcggtattttctccttacgcactctgtgcggtatttcacaccgcataatggtgcactctcagtacaatct  
gctctgatgccgcatagttaagccagccccgacacccgccaacacccgctgacgcgcccctgacgggctg  
tctgctcccggcatccgcttacagacaagctgtgaccgtctccgggagctgcatgtgtcagaggttttca  
ccgtcatcaccgaaacgcgcgagacgaaagggcctcgtgatagcctatttttataggttaatgtcatga  
taataatgggtttcttagacgtcaggtggcacttttccgggaaatgtgcgcggaaccctatttgtttatt  
tttctaaatacattcaaataatgtatccgctcatgagacaataaccctgataaatgcttcaataatattga  
aaaaggaagagatagattcaacatttccgctgctgccccttatcccttttttgcggcattttgccttc  
ctgtttttgctcaccagaaacgctgggtgaaagttaaagatgctgaagatcagttgggtgcagagatggg  
ttacatcgaactggatctcaacagcggtaagatccttgagagttttcggcccgaagaacgttttccaatg  
atgagcacttttaagttctgctatgtggcgcgggtattatccgctattgacgcgggcaagagcaactcg  
gtcgccgcatacactattctcagaatgacttgggtgagtagtaccagtcacagaaaagcactctacgga  
tggcatcagagtaagagaattatgcagtgctgccataaccatgagtgataaacactgcccgaacttactt  
ctgacaacgatcggagaccgaaggagctaaccgcttttttgcacaacatgggggatcatgtaactcgcc  
ttgatcgttgggaaccggagctgaatgaagccatacacaacgacgagcgtgacaccacgatgcctgtagc  
aatggcaacaacgttgcgcaaacatattaactggcgaactacttactctagcttcccggcaacaattaata  
gactggatggaggcggataaagttgcaggaccacttctgcgctcggcccttccggctggctggtttattg  
ctgataaatctggagccggtagcgtgggtctcgcggtatcattgcagcactggggccagatggtaagcc  
ctcccgtatcgtagttatctacacgacggggagtcaggcaactatggatgaacgaaatagacagatcgct  
gagataggtgcctcactgattaagcattggtaactgtcagaccaagtttactcatatatacttttagattg  
atttaaaacttcattttttaatttaaaaggatctaggtgaagatcctttttgataatctcatgaccaa  
cccttaacgtgagttttcgttccactgagcgtcagaccccgtagaaaagatcaaaggatcttcttgagat  
ccttttttctgcgcgtaactctgctgcttgcaaacaaaaaacaccgctaccagcgggtgggtttgtttgc  
cggatcaagagctaccaactctttttccgaaggtaactggcttcagcagagcgcagataccaaaactg  
ccttctagtgtagccgtagttaggccaccacttcaagaactctgtagcaccgcctacatacctcgctctg  
ctaactcctgttaccagtggtgctgcccagtgggcgaatagtcgtgtcttaccgggttggactcaagacgat  
agttaccggataaaggcgcagcgggtcgggctgaacgggggggttcgtgcacacagcccagcttggagcgaac  
gacctacaccgaactgagatacctacagcgtgagctatgagaaagcgcacgcttcccgaaggggagaaag  
gcggacaggtatccggtaagcggcaggggtcggaaacaggagagcgcacgagggagcttccaggggaaacg  
cctggatctttatagtcctgtcgggtttcggccactctgacttgagcgtcgatttttgtgatgctcgtc  
agggggggcggagcctatggaaaaacgcccagcaacgcggccttttacgggttccctggccttttgcctg

tttgctcacatggttctttcctgcggttatcccctgattctgtggataaccgtattaccgcctttgagtgag  
ctgataccgctcgccgcagccgaacgaccgagcgcagcgagtcagtgagcgaggaagcggaaga

Gaussia luciferase coding sequence (without signal sequence) cloned in  
BamHI and XbaI sites

257 atgaaaccaactgaaaacaatgaagatttcaacattgtagctgta  
M K P T E N N E D F N I V A V  
302 gctagcaactttgctacaacggatctcgatgctgaccgtggtaaa  
A S N F A T T D L D A D R G K  
347 ttgcccggaaaaaaattaccacttgaggactcaaagaaatggaa  
L P G K K L P L E V L K E M E  
392 gccaatgctaggaaagctggctgcactaggggatgtctgatatgc  
A N A R K A G C T R G C L I C  
437 ctgtcacacatcaagtgtacacccaaaatgaagaagtttatccca  
L S H I K C T P K M K K F I P  
482 ggaagatgccacacctatgaaggagacaaagaaagtgcacagggga  
G R C H T Y E G D K E S A Q G  
527 ggaataggagaggctattggtgacattcctgaaattcctggggtt  
G I G E A I V D I P E I P G F  
572 aaggatttggaaacccatggaacaattcattgcacaagttgaccta  
K D L E P M E Q F I A Q V D L  
617 tgtgtagactgcacaactggatgcctcaaaggctcttgccaatgtg  
C V D C T T G C L K G L A N V  
662 caatggtctgatttactcaagaaatggctgccacaaagatgtgca  
Q C S D L L K K W L P Q R C A  
707 acttttgctagcaaaattcaaggccaagtggacaaaataaagggt  
T F A S K I Q G Q V D K I K G  
752 gccgggtggtgatcatcaccatcaccatcactaa-784  
A G G D H H H H H H \*