



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

pUC18-Gluc\with\SS
expressing native Gaussia luciferase with the N-terminal secretion signal

gcgccaatacgcacaaccgcctctccccgcgcgttggccgattcattaatgcagctggcagcagaggtt
cccgactggaaagcgggcagtgagcgcacaacgcaattaatgtgagttagctcactcatttaggcaccccagg
ctttacactttatgcttccggctcgtatggtgtgtggaattgtgagcggataacaatttcacacaggaaa
cagctatgacctgattacgaattcgagctcggtagcccgggatccatgggagtgaaagtctttttgccc
cttatttgtattgctgtggccgagggcacaacaaactgaaaacaatgaagatttcaacattgtagctgtag
ctagcaactttgctacaacggatctcgatgctgaccgtggttaattgcccggaaaaaaattaccacttga
ggtagtcaaagaaatggaagcgaatgctaggaagctggctgcactaggggatgtctgatatgcctgtca
cacatcaagtgtacacccaaatgaagaagtttatcccaggaagatgccacacctatgaaggagacaaag
aaagtgcacagggaggaataggagaggctattgttgacattcctgaaattcctgggtttaaggatttggga
accatggaacaattcattgcacaagttgacctatgtgtagactgcacaactggatgcctcaaaggtctt
gccaatgtgcaatggtctgatttactcaagaaatggctgccacaaagatgtgcaacttttgctagcaaaa
ttcaaggccaagtggacaaaataaagggtgccgggtggtgatcatcaccatcaccatcactaatctagagt
cgacctgcaggcatgcaagcttggcactggccgtcgttttacaacgtcgtgactgggaaaaccctggcgt
taccacaacttaatcgcttgcagcacatcccccttccgcagctggcgtaatagcgaagaggcccgacc
gatcgcccttcccaacagttgcgagcctgaatggcgaatggcgcctgatgaggattttctccttacgc
atctgtgaggatttcacaccgcataatggtgcactctcagtagcaatctgctctgatgccgcatagttaag
ccagccccgacaccgcacaacaccgcctgacgcgcctgacgggcttgtctgctcccggcatccgcttac
agacaagctgtgaccgtctccgggagctgcatgtgtcagagggttttaccgctcatcccgaaacgcgcga
gacgaaagggcctcgtgatacgcctatttttatagggttaatgtcatgataataatgggtttcttagagctc
aggtggcacttttcggggaaatgtgagcgggaaccctatttgtttattttctaaatacattcaaataatg
tatccgctcatgagacaataaccctgataaatgcttcaataatattgaaaaaggaagagatagagattc
aacatttccgctgtcgcccttattcccttttttgcggcattttgccttccctgtttttgctcaccagaaac
gctggtgaaagtaaaagatgctgaagatcagttgggtgcagagtggttacatcgaactggatctcaac
agcggtaagatccttgagagttttcgccccgaagaacgttttccaatgatgagcacttttaagttctgc
tatgtggcgcgggtattatccgctattgacgcggggcaagagcaactcggtagccgcatacactattctca
gaatgacttgggttagtactcaccagtcacagaaaagcattctacggatggcatgacagtaagagaatta
tgagctgctgccataaccatgagtgataaacactgcccgaacttacttctgacaacgatcggaggaccga
aggagctaacgccttttttgcacaacatgggggatcatgtaactcgccttgatcgttgggaaccggagct
gaatgaagccataccaaacgacgagcgtgacaccacgatgcctgtagcaatggcaacaacgcttgcgcaaa
ctattaactggcgaactacttactctagcttcccggaacaattaatagactggatggaggcggataaag
ttgcaggaccacttctgcgctcggcccttccggctggctgggtttattgctgataaatctggagccgggtga
gcgtgggtctcgcgggtatcattgagcactggggccagatggtaagccctcccgatcgtagttatctac
acgacggggagtcaggcaactatggatgaacgaaatagacagatcgtgagataggtgcctcactgatta
agcattggtaactgtcagaccaagtttactcatatatacttttagattgatttaaaacttcatttttaatt
taaaaggatctaggtgaagatcctttttgataatctcatgaccaaaaatcccttaacgtgagtttctgctc
cactgagcgtcagaccccgtagaaaagatcaaaggatcttcttgagatccttttttctgagcgtaatct
gctgcttgcacaacaaaaaacaccgcctaccagcgggtggtttgtttgcccggatcaagagctaccaactct
ttttccgaaggtaactggcttcagcagagcgcagataccaaatactgtccttctagtgtagccgtagttta
ggccaccacttcaagaactctgtagcaccgcctacatacctcgtctgctaatcctgttaccagtggtctg
ctgccagtgggcgaatagtcgtgcttaccgggttggactcaagacgatagttaccggataaaggcgcagcg
gtcgggctgaacgggggggttctgtgcacacagcccagcttggagcgaacgacctacaccgaactgagatac
ctacagcgtgagctatgagaaagcgcacgccttcccgaaggagaaaggcggacaggtatccggtaagcg
gcaggggtcggaaacaggagagcgcacgagggagcttccaggggaaacgcctgggtatctttatagtcctgt
cgggttccgccacctctgacttgagcgtcgatttttgtgatgctcgtcaggggggcccggagcctatggaaa

aacgccagcaacgcggcctttttacggttcctggccttttgctggccttttgctcacatgttctttcctg
cgttatcccctgattctgtggataaccgtattaccgcctttgagtgagctgataccgctcgccgcagccg
aacgaccgagcgcagcgagtcagtgagcgaggaagcggaaga

Gluc coding sequence containing the signal sequence (bases 257 - 832)
clones in BamHI and XbaI sites

atgggagtgaaagttctttttgccccttatttgtattgctgtggccgaggccaaaccaact
·M·G·V·K·V·L·F·A·L·I·C·I·A·V·A·E·A·K·P·T·
gaaaacaatgaagatttcaacattgtagctgtagctagcaactttgctacaacggatctc
·E·N·N·E·D·F·N·I·V·A·V·A·S·N·F·A·T·T·D·L·
gatgctgaccgtggtaaattgcccgaaaaaattaccacttgaggactcaaagaaatg
·D·A·D·R·G·K·L·P·G·K·K·L·P·L·E·V·L·K·E·M·
gaagccaatgctaggaaagctggctgcactaggggatgtctgatatgcctgtcacacatc
·E·A·N·A·R·K·A·G·C·T·R·G·C·L·I·C·L·S·H·I·
aagtgtacacccaaaatgaagaagtttatcccaggaagatgccacacctatgaaggagac
·K·C·T·P·K·M·K·K·F·I·P·G·R·C·H·T·Y·E·G·D·
aaagaaagtgcacagggaggaataggagaggctattgttgacattcctgaaattcctggg
·K·E·S·A·Q·G·G·I·G·E·A·I·V·D·I·P·E·I·P·G·
ttaaaggatttggaaacccatggaacaattcattgcacaagttgacctatgtgtagactgc
·F·K·D·L·E·P·M·E·Q·F·I·A·Q·V·D·L·C·V·D·C·
acaactggatgcctcaaaggtcttgccaatgtgcaatgttctgatttactcaagaaatgg
·T·T·G·C·L·K·G·L·A·N·V·Q·C·S·D·L·L·K·K·W·
ctgccacaaagatgtgcaacttttgctagcaaaattcaaggccaagtggaacaaaataaag
·L·P·Q·R·C·A·T·F·A·S·K·I·Q·G·Q·V·D·K·I·K·
ggtgccggtggtgatcatcaccatcaccatcactaa
·G·A·G·G·D·H·H·H·H·H·H·*·