



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

pUC18-human-PtGFP

Blue is the pUC18 vector sequence
Underlined are the key restriction sites that were inserted with the PCR product
Green is the humanized PtGFP sequence

AAGCTTGCATGCCTGCAGGTCGACTCTAGAGGATCCCCCCACAGATCTATGAACCGCAACGTGCTGAAGA
ACACCGGCCTGAAGGAGATCATGAGCGCCAAGGCCAGCGTGGAGGGCATCGTTAACAACCACGTGTTTCAGC
ATGGAGGGCTTCGGCAAGGGCAACGTGCTGTTTCGGCAACCAATTGATGCAGATCCGCGTGACCAAGGGCGG
CCCCCTGCCCTTCGCCTTCGACATCGTGAGCATCGCCTTCCAGTACGGCAACCGTACGTTACCAAGTACC
CCGACGACATCGCCGACTACTTTCGTGCAGAGCTTCCCCGCCGGCTTCTTCTACGAGCGCAACCTGCGCTTC
GAGGACGGCGCCATCGTGGACATCCGCAGCGACATCAGCCTGGAGGACGACAAGTTCCACTACAAGGTGGA
GTACCGCGGCAACGGCTTCCCCAGCAACGGGCCCGTGATGCAGAAGGCCATCCTGGGCATGGAGCCCAGCT
TCGAGGTGGTGTACATGAACAGCGGCGTGTGGTGGGCGAGGTGGACCTGGTGTACAAGCTTGAGAGCGGC
AACTACTACAGCTGCCACATGAAGACCTTCTACCGTTCTGAAGGGCGGCGTGAAGGAGTTCCCCGAGTACCA
CTTCATCCACCACCGCCTCGAGAAGACCTACGTGGAGGAGGGCAGCTTCGTGGAGCAGCACGAGACCGCCA
TCGCCCAGCTGACCACCATCGGCAAACCCCTGATACTAGTGGGGGGATCCTCTAGAGTCGACCTGCANGC
ATGCAAGCTTGGCACTGGCCGTCGTTTTACAACGTCGTGACTGGGAAAACCTGGCGTTACCCAACCTTAAT
CGCCTTGCAGCACATCCCCCTTTCGCCAGCTGGCGTAATAGCGAAGAGGCCCGCACCGATCGCCCTTCCCA
ACAGNTGCGCAGCCTGAATGGNGAA

7 atgaaccgcaacgtgctgaagaacaccggcctgaaggagatcatg
M N R N V L K N T G L K E I M
52 agcgccaagggcagcgtggagggcatcgtaacaaccacgtgttc
S A K A S V E G I V N N H V F
97 agcatggagggcttcggcaagggcaacgtgctgttcggcaaccaa
S M E G F G K G N V L F G N Q
142 ttgatgcagatccgctgaccaagggcgccccctgcccttcgcc
L M Q I R V T K G G P L P F A
187 ttcgacatcgtgagcatcgcccttcagtacggcaaccgtacgttc
F D I V S I A F Q Y G N R T F
232 accaagtaccccgacgacatcgccgactacttcgtgcagagcttc
T K Y P D D I A D Y F V Q S F
277 cccgagcgttcttctacgagcgcaacctgcgcttcgaggacggc
P A G G F F Y E R N L R F E D G
322 gccatcgtggacatccgagcgacatcagcctggaggacgacaag
A I V D I R S D I S L E D D K
367 ttccactacaaggtggagtaccgcggaacggcttccccagcaac
F H Y K V E Y R G N G F P S N
412 gggcccgatgcagaaggccatcctgggcatggagcccagcttc
G P V M Q K A I L G M E P S F
457 gaggtggtgtacatgaacagcggcgtgctggtggcgaggtggac

E V V Y M N S G V L V G E V D
502 ctggtgtacaagcttgagagcggcaactactacagctgccacatg
L V Y K L E S G N Y Y S C H M
547 aagaccttctaccgttcgaagggcggcgtgaaggagttccccgag
K T F Y R S K G G V K E F P E
592 taccacttcatccaccaccgcctcgagaagacctacgtggaggag
Y H F I H H R L E K T Y V E E
637 ggcagcttcgtggagcagcagagaccgccatcgcccagctgacc
G S F V E Q H E T A I A Q L T
682 accatcggcaaaccctgatactagtgggggggatcctctag
T I G K P L I L V G G I L *