



TGACCAAGTCAGCTTGGCACTGGCCGTCGTTTTACAACGTCGTGACTGGGAAAACCCCTGGCGTTACCCAACTTAA
 TCGCCTTGCAGCACATCCCCCTTTCGCCAGCTGGCGTAATAGCGAAGAGGCCCGCACCGATCGCCCTTCCCAACA
 GTTGCAGCAGCCTGAATGGCGAATGGGAAATTGTAACGTTAATATTTTTGTTAATATTTTTGTTAAAAATTCGCGTTA
 AATTTTTGTTAAATCAGCTCATTTTTTTAACCAATAGGCCGAAAATCGGCAAAATCCCTTATAAAATCAAAAAGATAG
 ACCGAGATAGGGTTGAGTGTGTTCCAGTTTGGAAACAAGAGTCCACTATTAAAGAACGTGGACTCCAACGTCAAA
 GGGCGAAAACCGTCTATCAGGGCGATGGCCACTACGTGAACCATCACCTAATCAAGTTTTTTGGGGTCGAGG
 TGCCGTAAAGCACTAAATCGGAACCCTAAAGGGATGCCCGATTTAGAGCTTGACGGGGAAAAGCCGGCGAACGTG
 GCGAGAAAGGAAGGGAAGAAAGCGAAAGGAGCGGGCGCTAGGGCGCTGGCAAGTGTAGCGGTCACGCTGCGCGTA
 ACCACCACACCCGCGCGCTTAATGCGCCGCTACAGGGCGCGTCAGGTGGCACTTTTCGGGGAAATGTGCGCGGA
 ACCCTATTTGTTTTATTTTTCTAAATACATTCAAATATGTATCCGCTCATGAGACAATAACCTGATAAATGCTT
 CAATAATATTGAAAAGGAAGAGTATGAGTATTCAACATTTCCGTGTGCGCCTTATTCCCTTTTTTTCGGCATT
 TGCCTTCTGTTTTTGTCTCACCCAGAAACGCTGGTGAAAGTAAAAGATGCTGAAGATCAGTTGGGTGCACGAGTG
 GGTACATCGAAGTGGATCTCAACAGCGGTAAGATCCTTGAGAGTTTTTCGCCCCGAAGAACGTTTTTCCAATGATG
 AGCACTTTTAAAGTTCTGCTATGTGGCGCGGTATTATCCCGTATTGACGCCGGGCAAGAGCAACTCGGTGCGCCG
 ATACACTATTCTCAGAATGACTTGGTTGAGTACTACCAGTCACAGAAAAGCATCTTACGGATGGCATGACAGTA
 AGAGAATTATGCAGTGCTGCCATAACCATGAGTGATAAACAAGTGGCCAACTTACTTCTGACAACGATCGGAGGA
 CCGAAGGAGCTAACCGCTTTTTTGCACAACATGGGGGATCATGTAACCTCGCCTTGATCGTTGGGAACCGAGCTG
 AATGAAGCCATACCAAACGACGAGCGTGACACCAGATGCCTGTAGCAATGGCAACAACGTTGCGCAAACTATTA
 ACTGGCGAACTACTTACTCTAGCTTCCCGGCAACAATTAATAGACTGGATGGAGGCGGATAAAAGTTGCAGGACCA
 CTTCTGCGCTCGGCCCTTCCGGCTGGCTGGTTTTATTGCTGATAAATCTGGAGCCGGTGAGCGTGGGTCTCGCGGT
 ATCATTGCAGCACTGGGGCCAGATGGTAAGCCCTCCCGTATCGTAGTTATCTACAGCAGGGGAGTCAGGCAACT
 ATGGATGAACGAAATAGACAGATCGCTGAGATAGGTGCCTCACTGATTAAGCATTGGTAACTGTCAGACCAAGTT
 TACTCATATATACTTTAGATTGATTTAAAACCTTCATTTTTAATTTAAAAGGATCTAGGTGAAGATCCTTTTTTGAT
 AATCTCATGACCAAAATCCCTTAACGTGAGTTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGA
 TCTTCTTGAGATCCTTTTTTCTGCGCGTAATCTGCTGCTTGCAACAAAAAACCCAGCTACCAGCGGTGGTT
 TGTTTGCAGGATCAAGAGCTACCAACTCTTTTTCCGAAGGTAACCTGGCTTCAGCAGAGCGCAGATACCAAATACT
 GTCCTTCTAGTGTAGCCGTAGTTAGGCCACCACTTCAAGAACTCTGTAGCACCGCTACATACCTCGCTCTGCTA
 ATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCGTGTCTTACCAGGTTGGACTCAAGACGATAGTTACCG
 GATAAGGCGCAGCGGTGGGCTGAACGGGGGTTCTGTCACACAGCCAGCTTGGAGCGAACGACCTACACCGAA
 CTGAGATACCTACAGCGTGAGCTATGAGAAAAGCGCCACGCTTCCCGAAGGGAGAAAAGCGGACAGGTATCCGGTA
 AGCGGCAGGGTCCGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGGAAAACGCTGGTATCTTTATAGTCTGTG
 GGGTTTTGCCACCTCTGACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCC
 AGCAACGCGGCCTTTTTACGGTTCTGGCCTTTTGTGCTGACATGTTCTTCTCGGTTATCCCT
 GATTCTGTGGATAACCGTATTACCGCCTTTGAGTGAGCTGATAACCGCTCGCCGAGCCGAACGACCGAGCGCAGC

GAGTCAGTGAGCGAGGAAGCGGAAGAGCGCCCAATACGCAAAACCGCCTCTCCCCGCGCGTTGGCCGATTTCATTAA
TGCAGCTGGCAGCAGAGGTTTCCCAGACTGGAAAGCGGGCAGTGAGCGCAACGCAATTAATGTGAGTTAGCTCACT
CATTAGGCACCCAGGCTTTACACTTTATGCTTCCGGCTCGTATGTTGTGTGGAATTGTGAGCGGATAACAATTT
CACACAGGAAACAGCTATGACCATGATTACGAATTTGGCCAAGTCGGCCTCTAATACGACTCACTATAGGGAGCT
CGTCGAGCGGGCCGCTCGACGAATTAATTCCAATCCCACAAAAATCTGAGCTTAAACAGCACAGTTGCTCCTCTCAG
AGCAGAATCGGGTATTCAACACCCTCATATCAACTACTACGTTGTGTATAACGGTCCACATGCCGGTATATACGA
TGACTGGGGTTGTACAAAGGCGGCAACAAACGGCGTTCCCAGGAGTTGCACACAAGAAATTTGCCACTATTACAGA
GGCAAGAGCAGCAGCTGACGCGTACACAACAAGTCAGCAAAACAGACAGGTTGAACTTCATCCCCAAAGGAGAAGC
TCAACTCAAGCCCAAGAGCTTTGTAAAGGCCCTAACAAAGCCCACAAAGCAAAAAGCCCACTGGCTCACGCTAGG
AACCAAAAGGCCCAGCAGTGATCCAGCCCCAAAGAGATCTCCTTTGCCCCGGAGATTACAATGGACGATTTCTCT
CTATCTTTACGATCTAGGAAGGAAGTTTGAAGGTGAAGGTGACGACACTATGTTCACTACTGATAATGAGAAGGT
TAGCCTCTTCAATTTAGAAAGAATGCTGACCCACAGATGGTTAGAGAGGCCTACGCAGCAGGTTCTCATCAAGAC
GATCTACCCGAGTAACAATCTCCAGGAGATCAAATACCTTCCCAAGAAAGGTTAAAGATGCAGTCAAAAAGATTGAG
GACTAATTGCATCAAGAACACAGAGAAAGACATATTTCTCAAGATCAGAAGTACTATTCCAGTATGGACGATTCAG
AGGCTTGCTTCATAAACCAAGGCAAGTAATAGAGATTGGAGTCTCTAAAAAGGTAGTTCTACTGAATCTAAGGC
CATGCATGGAGTCTAAGATTCAAATCGAGGATCTAACAGAACTCGCCGTGAAGACTGGCGAACAGTTTCATACAGA
GTCTTTTACGACTCAATGACAAGAAGAAAATCTTCGTCAACATGGTGGAGCACGACACTCTGGTCTACTCCAAAA
ATGTCAAAGATACAGTCTCAGAAGACCAAGGGCTATTGAGACTTTTCAACAAAGGATAATTTTCGGGAAACCTCC
TCGGATTCCATTGCCAGCTATCTGTCACTTCATCGAAAGGACAGTAGAAAAGGAAGGTGGCTCCTACAAATGCC
ATCATTGCGATAAAGGAAAGGCTATCATTCAAGATCTCTCTGCCGACAGTGGTCCCAAGATGGACCCCCACCCA
CGAGGAGCATCGTGAAAAAGAAGACGTTCCAACCACGTTTCAAAAGCAAGTGGATTGATGTGACATCTCCACTG
ACGTAAGGGATGACGCACAATCCCCTACTATCCTTCGCAAGACCCTTCTCTATATAAGGAAGTTCATTTCAATTTGG
AGAGGACACGCTCGAGAACATGaacgcaacgtgctgaaaaacaccggtctgaaagaaattatgagcgcgaaaag
cgagcgtggaaggcattgtgaacaaccatgtgttagcatggaaggctttggcaaaggcaacgtgctgtttggca
accagctgatgcagattcgcgctgaccaaaggcggtcgctgccggttgctgctggtggcggaagtggatctggtgata
agtatggcaaccgcacctttaccaaatatccggatgatattgaggattatgtgtgcagagctttccggcaggct
tttttatgaacgcaacctgcgctttgaagatggtgcgattgtggatattcgagcagatattagcctggaagatg
ataaatttcattataaagtggaatatcgcggtaacggctttccgagcaacggctccgggtgatgcagaaagcgattc
tgggcatggaaccgagctttgaagtgggtgatatgaacagcggtgtgctggtggcggaagtggatctggtgata
aactggaagcggaactattatagctgccatataaaaacctttatcgagcaagggcggtgtgaaagaatttc
cggaatatcattttatcatcatcgctggaaaaaacctatgtggaagaaggcagctttgtggaacagcatgaaa
ccgagattgagcagctgaccaccattggcaaacgctgggagcctgcatgaatgggtgTAACTGACTGACTGAA
AGCTTGGATCCTCTAGAGTCTCTGCTTTAATGAGATATGCGAGACGCCTATGATCGCATGATATTTGCTTTCAATT
CTGTTGTGCACGTTGTAAAAACCTGAGCATGTGTAGCTCAGATCCTTACCGCCGTTTCGGTTCATTCTAATGA
ATATATCACCCGTTACTATCGTATTTTTATGAATAATATTCTCCGTTCAATTTACTGATTGTACCCTACTACTTA
TATGTACAATATTAATAAATAAACAATATATTGTGCTGAATAGGTTTATAGCGACATCTATGATAGAGCGCCACA
ATAACAACAATTGCGTTTTATTATTACAAATCCAATTTTAAAAAAGCGGCAGAACCAGTCAAACCTAAAAGAC
TGATTACATAAATCTTATTCAAATTTCAAAGGCCCCAGGGGCTAGTATCTACGACACACCGAGCGGCGAACTAA
TAACGTTCACTGAAGGGAACCTCCGTTCCCCGCCGCGCATGGGTGAGATTCTTTGAAGTTGAGTATTGGCCG
TCCGCTCTACCGAAAGTTACGGGCACCATTCAACCCGTTCCAGCACGGCGGCCGGGTAACCGACTTGCTGCCCCG
AGAATTATGCAGCATTTTTTTGGTGTATGTGGGCCCCAAATGAAGTGCAGGTCAAACCTTGACAGTGACGACAAA
TCGTTGGGCGGTTCCAGGGCGAATTTTGCACAACATGTGAGGCTCAGCAGGACCTGCAGGCATGCAAGCTAGC
TTACTAGTGATGCATATTCTATAGTGTACCTAAATCTGCGGCCGC

blue = start codon
red = stop codon
underlined = restriction sites
underlined bold = NotI restriction site
yellow = CAMV 35S promoter
green = ORF of PtGFP
pink = initiation sequence
cyan = terminator