

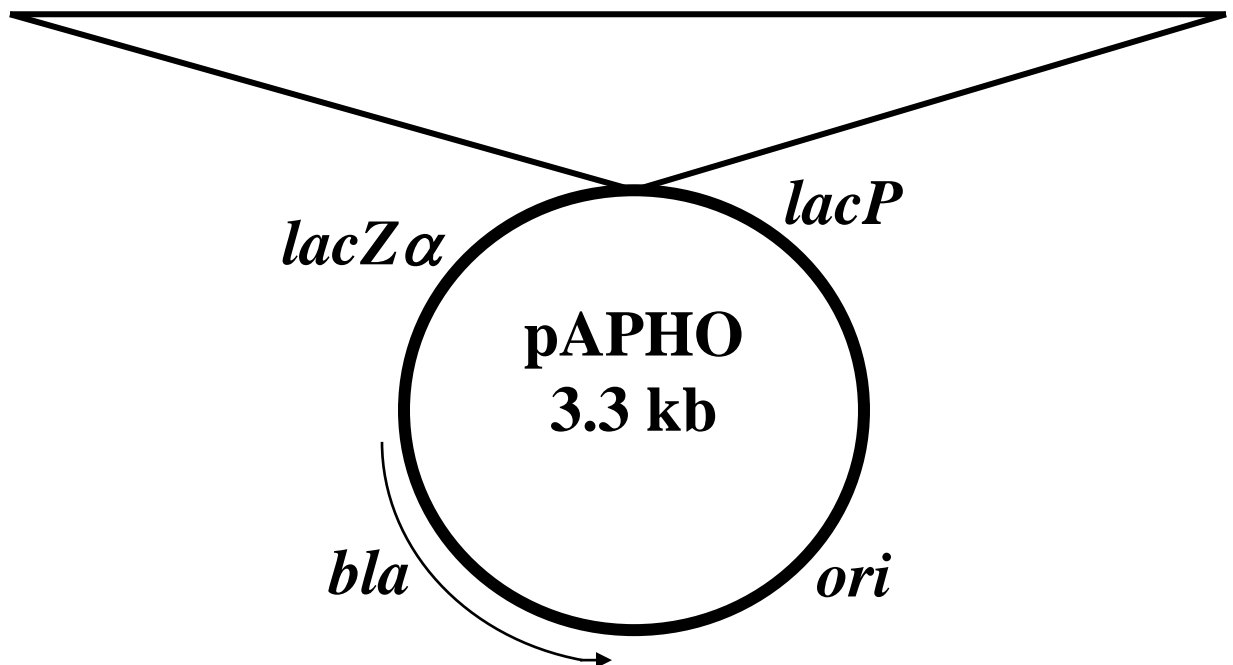


Post Office Box 2746 Pinetop, Arizona 85935
 Telephone (928) 367-1200 Fax (928) 367-1205

www.nanolight.com

pUC19 Aequorin Photoprotein

G AAT TCC AAG CTT ACA TCA GAC TTC GAC AAC CCA AGA TGG ATT GGA
 CGA CAC AAG CAT ATG TTC AAT TTC CTT GAT GTC AAC CAC AAT GGA AAA
 ATC TCT CTT GAC GAG ATG GTC TAC AAG GCA TCT GAT ATT GTC ATC AAT
 AAC CTT GGA GCA ACA CCT GAG CAA GCC AAA CGA CAC AAA GAT GCT GTA
 GAA GCC TTC TTC GGA GGA GCT GGA ATG AAA TAT GGT GTG GAA ACT GAT
 TGG CCT GCA TAT ATT GAA GGA TGG AAA AAA TTG GCT ACT GAT GAA TTG
 GAG AAA TAC GCC AAA AAC GAA CCA ACG CTC ATC CGT ATA TGG GGT GAT
 GCT TTG TTT GAT ATC GTT GAC AAA GAT CAA AAT GGA GCC ATT ACA CTG
 GAT GAA TGG AAA GCA TAC ACC AAA GCT GCT GGT ATC ATC CAA TCA TCA
 GAA GAT TGC GAG GAA ACA TTC AGA GTG TGC GAT ATT GAT GAA AGT GGA
 CAA CTC GAT GTT GAT GAG ATG ACA AGA CAA CAT TTA GGA TTT TGG TAC
 ACC ATG GAT CCT GCT TGC GAA AAG CTC TAC GGT GGA GCT GTC CCC TAA
 GAA GCT CTA CGG TGG TGA TGC ACC CTG GGA AGA TGA TGT GAT TTT GAA
TAA AAC ACT GAT GAA *TTC*



pAPHO - Aequorin photoprotein coding sequence for the mature apoprotein cloned into plasmid pUC19 from pB-IE on an *Eco*RI fragment (shown in *italics* on the sequence). Stop codon is