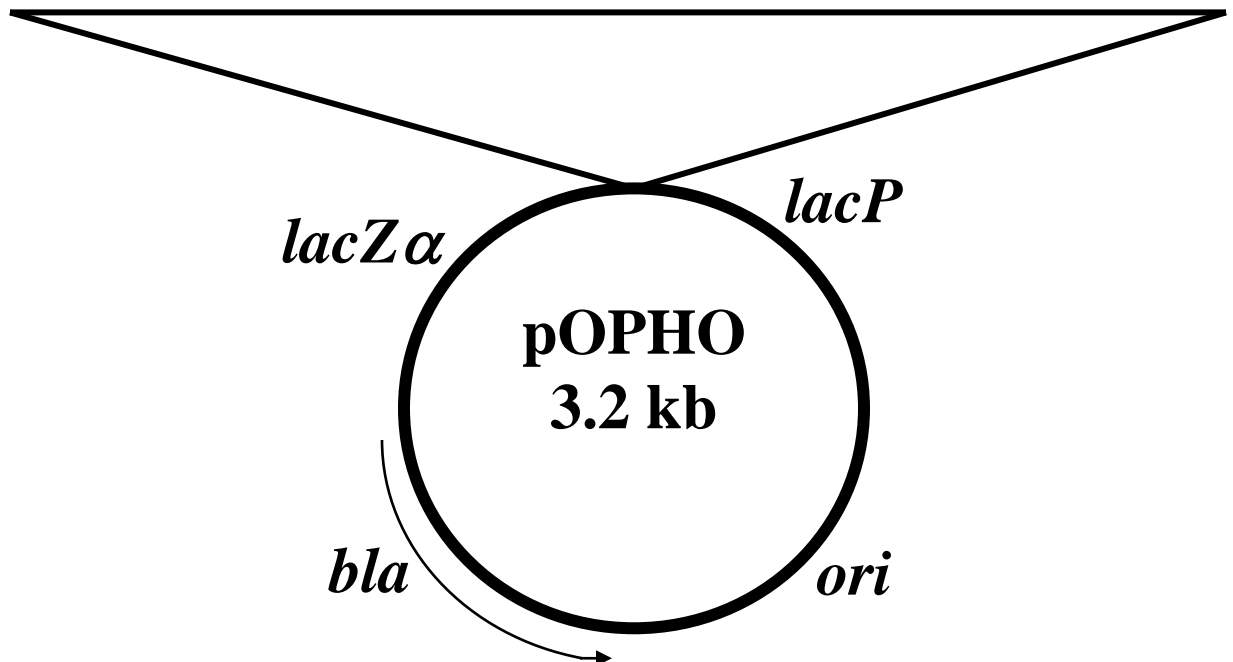




Post Office Box 2746 Pinetop, Arizona 85935  
 Tel: (928) 367-1200 Fax: (928) 367-1205  
[www.nanolight.com](http://www.nanolight.com)

### PUC 19 Obelin Photoprotein

GAA *TTC* ATG AGC AGC AAA TAT GCG GTG AAA CTG AAA ACC GAT TTT GAT  
 AAT CCG CGT TGG ATT AAA CGC CAT AAA CAT ATG TTT GAT TTT CTG GAT  
 ATT AAC GGC AAC GGC AAA ATT ACC CTG GAT GAA ATT GTG AGC AAA GCG  
 AGC GAT GAT ATT TGC GCG AAA CTG GAA GCG ACC CCA GAA CAG ACC AAA  
 CGC CAT CAG GTG TGC GTG GAA GCG TTT TTT CGC GGT TGC GGC ATG GAA  
 TAT GGC AAA GAA ATT GCG TTT CCG CAG TTT CTG GAT GGC TGG AAA CAG  
 CTG GCG ACC AGC GAA CTG AAA AAA TGG GCA CGC AAC GAA CCG ACC CTG  
 ATT CGC GAA TGG GGT GAT GCG GTG TTT GAT ATT TTT GAT AAA GAT GGC  
 AGC GGC ACC ATT ACC CTG GAT GAA TGG AAA GCG TAT GGC AAA ATT AGC  
 GGC ATT AGC CCG AGC CAG GAA GAT TGC GAA GCG ACC TTT CGC CAT TGC  
 GAT CTG GAT AAC AGC GGT GAT CTG GAT GTG GAT GAA ATG ACC CGT CAG  
 CAT CTG GGC TTT TGG TAT ACC CTG GAT CCG GAA GCG GAT GGC CTG TAT  
 GGC AAC GGC GTG CCG CTC GAG AAG GGC GAA *TTC*



pOPHO - Obelin photoprotein coding sequence cloned into plasmid pUC19 on an *EcoRI* fragment (shown in *italic* on the sequence) in the reverse orientation relative to the *lac* promoter. The