



Post Office Box 2746 Pinetop, Arizona 85935  
Telephone (928) 367-1200 Fax (928) 367-1205  
[www.nanolight.com](http://www.nanolight.com) E-mail: [info@prolume.com](mailto:info@prolume.com)

### pCMV-h-Pt GFP

gacggatcgggagatctcccgatcccctatgggtcgactctcagtacaatctgctctgatgccgcatagtta  
agccagtatctgctccctgcttgtgtgttggaggtcgctgagtagtgcgcgagcaaaatttaagctacaac  
aaggcaaggcttgaccgacaattgcatgaagaatctgcttaggggttaggcgcttttgccgctgcttcgcgatg  
tacggggccagatatacgcgcttgacattgattattgactagttattaatagtaatcaattacggggtcatta  
gttcatagcccataatggagttccgcgcttacataacttacggtaaatggcccgcctggctgaccgccc  
cgacccccgcccattgacgtcaataatgacgtatggtcccatagtaacgccaatagggactttccattgac  
gtcaatgggtggactatctacggtaaactgcccacttggcagtagatcaagtgtatcatatgccaagtacg  
ccccctattgacgtcaatgacggtaaatggcccgcctggcattatgccagtagatgaccttatgggactt  
tccctacttggcagtagatctacgtattagtcacgctattaccatgggtgatgcggttttggcagtagatca  
atggggcgtggatagcgggttgactcacggggatttccaaaatgctcgaacaactccgcccattgacgtcaatgggagttg  
tttggcaccaaaatcaacgggactttccaaaatgctcgaacaactccgcccattgacgtcaatgggagttg  
taggcggtgtagcgtgggaggtctatataagcagagctctctggctaactagagaaccactgcttactggc  
ttatcgaaaattaatacgcactcactatagggagaccgaagcttggtagcagctcgGATCTATGAACCGCAA  
CGTGCTGAAGAACACCGGCCTGAAGGAGATCATGAGCGCCAAGGCCAGCGTGGAGGGCATCGTTAACAACC  
ACGTGTTTCAGCATGGAGGGCTTCGGCAAGGGCAACGTGCTGTTTCGGCAACCAATTGATGCAGATCCGCGTG  
ACCAAGGGCGGCCCTTCCCTTCGCCTTCGACATCGTGAGCATCGCCTTCCAGTACGGCAACCGTACGTT  
CACCAAGTACCCCGACGACATCGCCGACTACTTCGTGCAGAGCTTCCCGCCGGCTTCTTCTACGAGCGCA  
ACCTGCGCTTCGAGGACGGCGCCATCGTGGACATCCGCAGCGACATCAGCCTGGAGGACGACAAGTTCAC  
TACAAGGTGGAGTACCGCGCAACGGCTTCCCAGCAACGGGCCCGTGATGCAGAAGGCCATCCTGGGCAT  
GGAGCCCAGCTTCGAGGTGGTGTACATGAACAGCGCGTGTGGTGGGCGAGGTGGACCTGGTGTACAAGC  
TTGAGAGCGGCAACTACTACAGCTGCCACATGAAGACCTTCTACCGTTTGAAGGGCGGCGTGAAGGAGTTC  
CCCAGTACCCTTCCATCCACCACCGCCTCGAGAAGACCTACGTGGAGGAGGGCAGCTTCGTGGAGCAGCA  
CGAGGCCGCCATCGCCAGCTGACCACCATCGGCAAGCCCCCTGGGATCCCTGCACGAGTGGGTGTAGGGGA  
TCCTctagaataaattcttactgtcatgccaagtaagatgcttttctgtgctgcaatagcagggcatgctgg  
ggatgcggtgggctctatggcttctgagggcggaaagaaccagctggggctctagggggatccccacgcgc  
cctgtagcggcgcattaaagcgcggcgggtgtggtggttacgcgcagcgtgaccgctacacttgccagcgc  
ctagcggccgctcctttcgctttcttcccttcttctcgccacgttcgcccggctttccccgctcaagctct  
aaatcggggcatccttttagggttccgatttagtgctttacggcacctcgacccccaaaaaacttgattag  
gtgatgggtcacgtagtgggccatcgccctgatagacgggttttccgcccattgacgttggagtcacgctt  
tttaatagtgtagctctgttccaaaactggaacaacactcaaccctatctcggctctattcttttgattata  
agggattttggggatttcggcctattgggttaaaaaatgagctgatttaacaaaaatttaacgcgaattaat  
tctgtggaatgtgtgctcagttaggggtgtggaagtcccaggctcccaggcaggcagaagtagcaaacg  
atgcatctcaattagtcagcaaccaggtgtggaagtcccaggctcccagcaggcagaagtagcaaacg  
catgcatctcaattagtcagcaaccatagtcggcccttaactccgcccattccgccccttaactccgccc  
gttccgcccattctccgcccattggctgactaattttttttatattatgacagaggccgaggccgcctctgcc  
tctgagctattccagaagtagtgaggaggcttttttggaggcctaggcttttgcaaaaagctcccgggagc  
ttgtatatccatcttccggatctgatcaagagacaggatgaggatcgtttcgcatgatgcaacaagatggat  
tgcacgcaggttctccggccgcttgggtggagaggctattcggctatgactgggcacaacagacaatcggc  
tgctctgatgccgctgttccggctgtcagcgcaggggcccggcttctttttgtcaagaccgacctgtc  
cgggtgccctgaatgaactgcaggacgaggcagcgcggctatcgtggctggccacgacgggcttcttgcg  
cagctgtgctcgacgttgtcactgaagcgggaagggactggctgctattgggccaagtgcggggcaggat  
ctcctgtcatctcaccttgcctcctgcccagaaagtatccatcatggctgatgcaatgcggcggctgcatac  
gcttgatccggctacctgcccattcgaccaccaagcgaacaatcgcatcgagcgcagcgtactcggatgg  
aagccggcttctgtagcaggtgatctggacgaagagcatcaggggctcgcgccagccgaactgttcgccc

aggctcaaggcgcgcatgccccgacggcgaggatctcgtcgtgacccatggcgatgcctgcttgccgaatat  
catgggtggaaaatggcgcgttttctggattcatcgactgtggccggctgggtgtggcggaccgctatcagg  
acatagcgttggctaccctgatattgctgaagagcttggcggcgaatgggctgaccgcttctcgtgctt  
tacggatcgccgctcccgatcgcagcgcacgccttctatcgccttcttgacgagttctctgagcggg  
actctgggggttcgaaatgaccgaccaagcgcgcaccaacctgccatcacgagatctcgattccaccgccc  
cttctatgaaagggttgggcttcggaatcgtttccgggacgcggctggatgatcctccagcgcggggatc  
tcatgctggagttcttcgcccacccaacttgtttattgcaacttataatgggtacaaaataagcaatagc  
atcacaaatcacaataaagcatttttttctactgcattctagttgtgggttgtcctcctcctcctcctc  
atcttatcatgtctgtataaccgtcgacctctagctagagcttggcgtaatcatgggtcatagctgttctc  
tgtgaaattggttatccgctcacaattccacacaacatacgcagccggaagcataaagtgtaaagcctgggg  
gcctaagtgagtgagctaaactcacattaattgctgtgctcactgcccgttccagtcgggaaacctgtc  
gtgccagctgcattaatgaatcggccaacgcgcggggagagggcttgcgtattgggctcctccgctt  
cctcgtcactgactcgtcgcctcggctcgtcggctgcggcgagcggatcagctcactcaaaggcgta  
atacgggttatccacagaatcaggggataacgcaggaagaacatgtgagcaaaaggccagcaaaaggccag  
gaaccgtaaaaaggccgcgcttgcggcttttccataggctccgccccctgacgagcatcacaaaaatc  
gacgctcaagtcaagggtggcgaaaccgcagaggactataaagataaccaggcgttccccctggaagctcc  
ctcgtgctcctcctgttccgacctgcccgttaccggatacctgtccgccttctccttccggaagcgt  
ggcgttctcaatgctcacgctgtaggatctcagttcgggtgtaggtcgttcgctccaagctgggctgtg  
tgcacgaacccccctcagccccgaccgctgcgccttaccggtaactatcgtcttgagccaacccggta  
agacacgacttatcgccactggcagcagccactggtaacaggattagcagagcgaggtatgtaggcgggtgc  
tacagagttcttgaagtgggtggcctaactacggctacactagaaggacagtatctgggtatctgcgctctgc  
tgaagccagttaccttcggaaaaagagttggtagctcttgatccggcaaaacaaaccaccgctggtagcgg  
ggttttttgtttgcaagcagcagattacgcgcagaaaaaaggatctcaagaagatcctttgatctttc  
tacggggtctgacgctcagtggaacgaaaactcacgtaagggttttgggtcatgagattatcaaaaagga  
tcttcacctagatccttttaaatataaaatgaagttttaaatacaatctaaagtataatgagtaaaacttg  
tctgacagttaccaatgcttaatcagtgaggcacctatctcagcgatctgtctatctcgttcatccatagt  
tgctgactccccgtcgtgtagataactacgatacgggagggcttaccatctggccccagtgctgcaatga  
taccgcgagaccacgctcaccggctccagatttatcagcaataaaccagccagccggaagggccgagcgc  
agaagtggctcctgcaactttatccgcctccatccagctctaatattgttgccgggaagctagagtaagtag  
ttcgccagttaatagttgcgcaacgctgttgccattgctacaggcatcgtgggtgtcacgctcgtcgtttg  
gtatggcttcatcagctccggctcccaacgatcaaggcgagttacatgatccccatgtgtgcaaaaaa  
gcggttagctccttcggctcctccgatcgttgtcagaagtaagttggccgcagtggttatcactcatgggtat  
ggcagcactgcataattctcttactgtcatgccatccgtaagatgcttttctgtgactgggtgagtaactcaa  
ccaagtcattctgagaatagtgatgctggcgaccgagttgctcttgcccggcgtcaatacgggataatacc  
gcgccacatagcagaactttaaaagtgtcatcattggaaaacgcttcttcggggcgaaaactctcaaggat  
cttaccgctgttgagatccagttcgatgtaaccactcgtgcacccaactgatcttcagcatcttttactt  
tcaccagcgtttctgggtgagcaaaaacaggaaggcaaaatgccgcaaaaagggaataagggcgacacgg  
aaatggtgaatactcactcttcttttcaatattattgaagcatttatcaggggttatgtctcatgag  
cggatacatatttgaaatgatttagaaaaataaacaataaggggttccgcgcacatttccccgaaaagtg  
cacctgacgtc

