

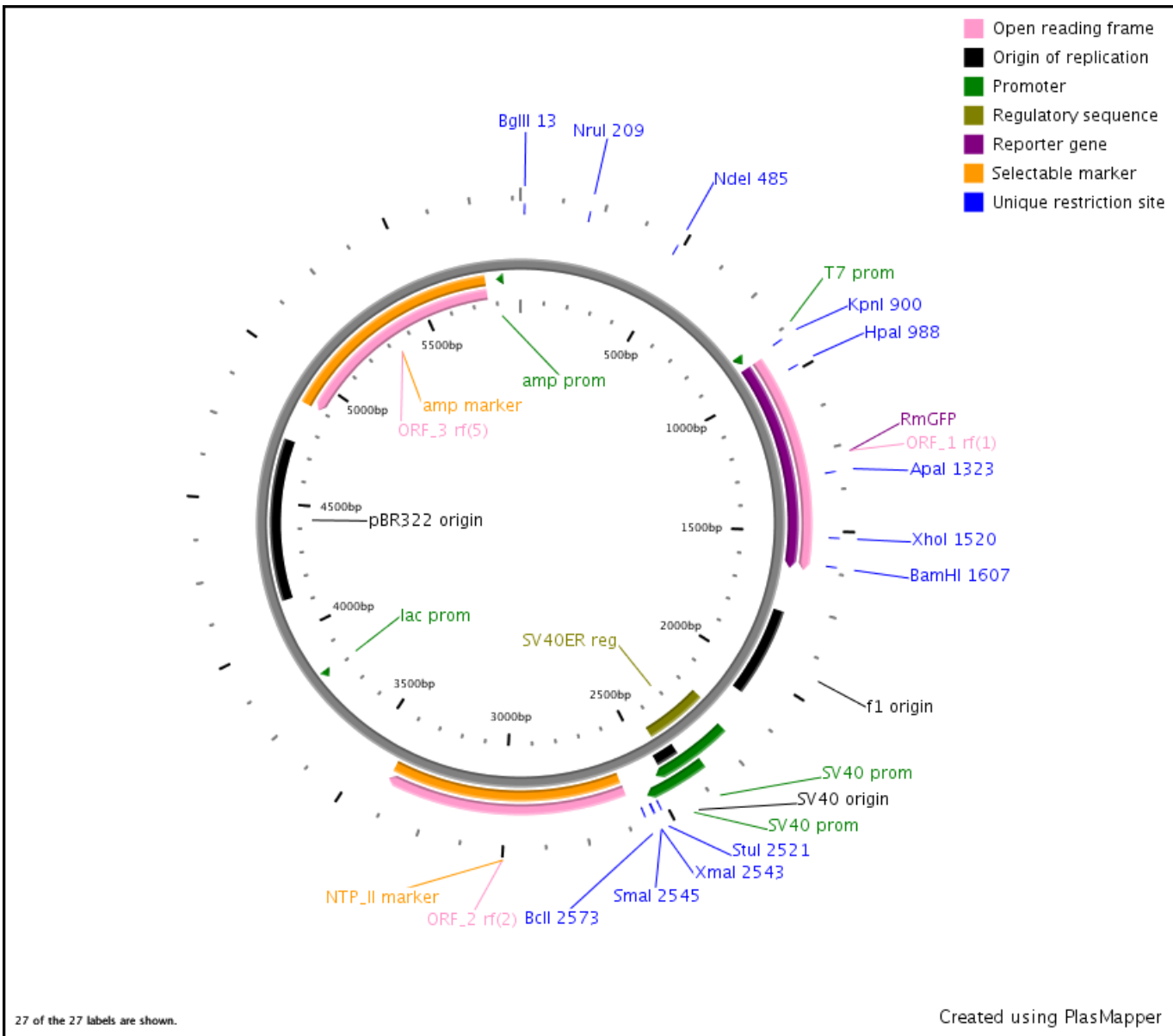


PO Box 2746, Pinetop, AZ 85935
Telephone (928) 367-1200 Fax (928) 367-1205
www.nanolight.com E-mail: info@prolume.com

pCMV-human-R.Mullerei GFP

gacggatcgggagatctcccgatcccctatgggtcgactctcagtacaatctgctctgatgccgcatagtta
agccagtatctgctccctgcttgtgtgttggaggtcgctgagtagtgcgcgagcaaaatttaagctacaac
aaggcaaggcttgaccgacaattgcatgaagaatctgcttaggggttaggcggttttgcgctgcttcgctgatg
tacgggcccagatatacgcgcttgacattgattattgactagttattaatagtaatacaattacgggggtcatta
gttcatagcccataatggagttccgcgttacataacttacggtaaatggcccgcctggctgaccgccc
cgacccccgcccattgacgtcaataatgacgtatgttcccatagtaacgccaatagggactttccattgac
gtcaatgggtggactatttacggtaaacgcccacttggcagtagatcaagtgtatcatatgccaaagtacg
ccccctattgacgtcaatgacggtaaatggcccgcctggcattatgccagtagatgaccttatgggactt
tccacttggcagtagatctacgtattagtcacgctattaccatgggtgatgcccgttttggcagtagatca
atgggctgtagtagcgggtttgactcacggggatttccaagtctccacccttagcgtcaatgggagtttg
tttggcaccaaaatcaacgggactttccaaaatgtcgtacaactccgcccattgacgcaaatgggctg
taggcgtgtacgggtgggaggtctatataagcagagctctctggctaactagagaaccactgcttactggc
ttatcgaaattaatacagactcactataggagacccaagcttggtagcagctcgGATCTATGAGCAAGCA
GATCCTGAAGAACACCTGCCTGCAGGAGGTGATGAGCTACAAGGTGAACCTGGAGGGCATCGTTAACAACC
ACGTGTTACCCATGGAGGGCTGCGGCAAGGGCAACATCCTGTTTCGGCAACCAATTGGTGCAGATCCGCGTG
ACCAAGGGCGCCCCCTGCCCTTCGCCTTCGACATCGTGAGCCCCGCCTTCCAGTACGGCAACCGTACGTT
CACCAAGTACCCCAACGACATCAGCGACTACTTCATCCAGAGCTTCCCCGCCGGCTTCATGTACGAGCGCA
CCCTGCGCTACGAGGACGGCGGCCTGGTGGAGATCCGCGAGCATCAACCTGATCGAGGACAAGTTCGTG
TACCGCGTGGAGTACAAGGGCAGCAACTTCCCGACGACGGGCCCGTGTATGCAGAAGACCATCCTGGGCAT
CGAGCCCAGCTTCGAGGCCATGTACATGAACAACGGCGTGTCTGGTGGGCGAGGTGATCCTGGTGTACAAGC
TTAACAGCGGCAAGTACTACAGCTGCCACATGAAGACCCTGATGAAGAGCAAGGGCGTGGTGAAGGAGTTC
CCCAGCTACCACTTCATCCAGCACCGCCTCGAGAAGACCTACGTGGAGGACGGCGGCTTCGTGGAGCAGCA
CGAGACCGCCATCGCCAGATGACCAGCATCGGCAAGCCCCTGGGATCCCTGCACGAGTGGGTGTGATAct
agaaataatttctactgtcatgccaagtaagatgcttttctgtgctgcaatagcaggcatgctggggatgc
gggtgggctctatggcttctgagggcggaaagaaccagctggggctctagggggatcccacgcgcctgta
gcggcgcattaagcgcggcgggtgtggtggttacgcgcagcgtgaccgctacacttgccagcgcctagcg
cccgtcctttcgttttcttcccttcccttctcgcacagttcgcggctttcccctcaagctctaaatcg
gggcatccctttagggttccgatttagtgctttacggcacctcgacccccaaaaaacttgattaggggtgatg
gttcacgtagtgggcccacgcccctgatagacgggttttccgccccttgacgttggagttccaggttctttaat
agtggactcttgttccaaactggaacaacactcaaccctatctcggctctattcttttgatttataagggat
tttgggatttccggcctattggttaaaaaatgagctgatttaacaaaaatttaacgcgaatattctgtg
gaatgtgtgtcagttagggtgtggaagtccccaggctccccaggcaggcagaagtagcaaaagcatgcat
ctcaattagtcagcaaccaggtgtggaagtccccaggctccccaggcaggcagaagtagcaaaagcatgca
tctcaattagtcagcaaccatagtcggcccctaaactccgcccctccgcccctaaactccgcccagttccg
cccattctccgcccctggctgactaattttttttatattatgacagaggccgaggccgcctctgcctctgag
ctattccagaagtagtgaggaggttttttggaggcctaggcttttgcaaaaagctcccgggagcttgtat
atccattttcggatctgatcaagagacaggatgaggatcgtttccgatgattgaacaagatggattgcacg
cagggttctccggccgcttgggtggagaggctattcggctatgactgggcacaacagacaatcggctgctct
gatgccgcccgtgttccggctgtcagcgcaggggcccgggttctttttgtcaagaccgacctgtccgggtgc
cctgaatgaactgcaggacgaggcagcgcggctatcgtggctggccacgacgggcttccctgcgagctg
tgctcgacgttgtcactgaagcgggaaggactggctgctattgggccaagtgcggggcaggatctcctg
tcatctcaccttgctcctgccgagaaagtagtccatcatggctgatgcaatgcggcggctgcatacgttga

tccggctacctgccattcgaccaccaagcgaacatcgcatcgagcgcgacgactcggatggaagccg
gtcttgtcgatcaggatgatctggacgaagagcatcaggggctcgcgccagccgaactgttcgccaggctc
aaggcgcgcgatgcccgcggcgaggatctcgtcgtgaccatggcgatgcctgcttgccgaatatcatggg
ggaaaatggccgcttttctggattcatcgactgtggccggctgggtgtggcggaccgctatcaggacatag
cgttggctacctgctgatattgctgaagagcttggcggcgaatgggctgaccgcttcctcgtgctttacgg
atcgccgctcccgatcgcagcgcacgccttctatcgccttcttgacgagttcttctgagcgggactctg
gggttcgaaatgaccgaccaagcgcgcccacactgccatcacgagatttcgattccaccgccccttcta
tgaaagggtgggcttcggaatcgttttccgggacgcgggctggatgatcctccagcgcggggatctcatgc
tggagtcttccgccaccccaacttgtttattgacgcttataatgggtatacaataaagcaatagcatcaca
aatctcacaataaagcatttttttcaactgacttctagttgtgggttggccaaactcatcaatgtatctta
tcatgtctgtataaccgctcgacctctagctagagcttggcgtaaatcatgggtcatagctgtttcctgtgtgaa
attgttaccgctcacaattccacacaacatacagcgggaagcataaagtgtaaagcctgggggtgcctaa
tgagtgcgtaactcacattaattgctgctcactgcccgtttccagtcgggaaacctgctgctgcca
gctgcattaatgaatcggccaacgcgcggggagaggcgggttgcgtattgggcgctcttcgcttcctcgc
tcaactgactcgcctcggctcgttcggctcggcgagcgggtatcagctcactcaaaggcggtaatacgg
ttatccacagaatcaggggataacgcaggaaagaacatgtgagcaaaaggccagcaaaaggccaggaaccg
taaaaaggccgcgcttgcctggcgttttccataggctccgccccctgacgagcatcacaaaaatcgacgct
caagtgcagagggtggcgaaacccgacaggactataaagataaccaggcgtttccccctggaagctccctcgtg
cgctctcctgttccgacctgcccgttaccggatacctgtccgctttctcccttcgggaagcgtggcgct
ttctcaatgctcacgctgttaggtatctcagttcgggtgtaggtcgttcgctccaagctgggctgtgtgcacg
aaccctccgcttcagcccgcagcctgccccttaccggtaactatcgtcttgagccaaccggtaagacac
gacttatcgccactggcagcagccactggtaacaggattagcagagcagaggtatgtaggcgggtgctacaga
gttcttgaagtggtggcctaactacggctacactagaaggacagtatgtggatctgcgctctgctgaagc
cagttaccttcggaaaaagagttggtagctcttgatccggcaaaacaaaccaccgctggtagcgggtgggttt
ttgtttgcaagcagcagattacgcgcagaaaaaaggatctcaagaagatcctttgatctttctacggg
gtctgacgctcagtggaacgaaaactcacgtaagggattttgggtcatgagattatcaaaaaggatcttca
cctagatccttttaaattaaaaatgaagttttaaatacaatctaaagtataatgagtaaaacttggctgac
agttaccaatgcttaatcagtgaggcacctatctcagcgcgactgtctatcttctgctccatagttgctg
actccccgctcgtgtagataaactacgatacgggagggcttaccatctggccccagtgctgcaatgataccgc
gagaccacgctcaccggctccagatttatcagcaataaaccagccagccggaagggccgagcagagaagt
ggctcctgcaactttatccgctccatccagctctattaattgttgcgggaagctagagtaagtagttcggc
agttaatagtttgcgcaacgcttgttgcattgctacaggcatcgtgggtgtcacgctcgtcgtttggatgg
cttcatcagctccgggttcccaacgatcaaggcaggttacatgatccccatggtgtgcaaaaaagcgggt
agctccttcggctcctccgatcgttgtcagaagtaagttggccgcagtggttatcactcatgggttatggcagc
actgcataattctcttactgtcatgccatccgtaagatgcttttctgtgactgggtgagtagtcaaccaagt
cattctgagaatagtgatgccccgaccgagttgctcttgcggcgctcaatacgggataataccgcgcca
catagcagaactttaaaagtgctcatcattggaaaacgcttcttcggggcgaaaactctcaaggatcttacc
gctgttgagatccagttcgatgtaacccactcgtgcacccaactgatcttcagcatcttttactttacca
gcgtttctgggtgagcaaaaacaggaaggcaaaatgccgcaaaaagggaataagggcgacacggaaatgt
tgaatactcatactcttcccttttcaatattatgaagcatttatcaggggtattgtctcatgagcggata
catatttgaatgtatttagaaaaataaacaataaggggttccgcgcacatttccccgaaaagtgccacctg
acgtc



- █ Open reading frame
- █ Origin of replication
- █ Promoter
- █ Regulatory sequence
- █ Reporter gene
- █ Selectable marker
- █ Unique restriction site

