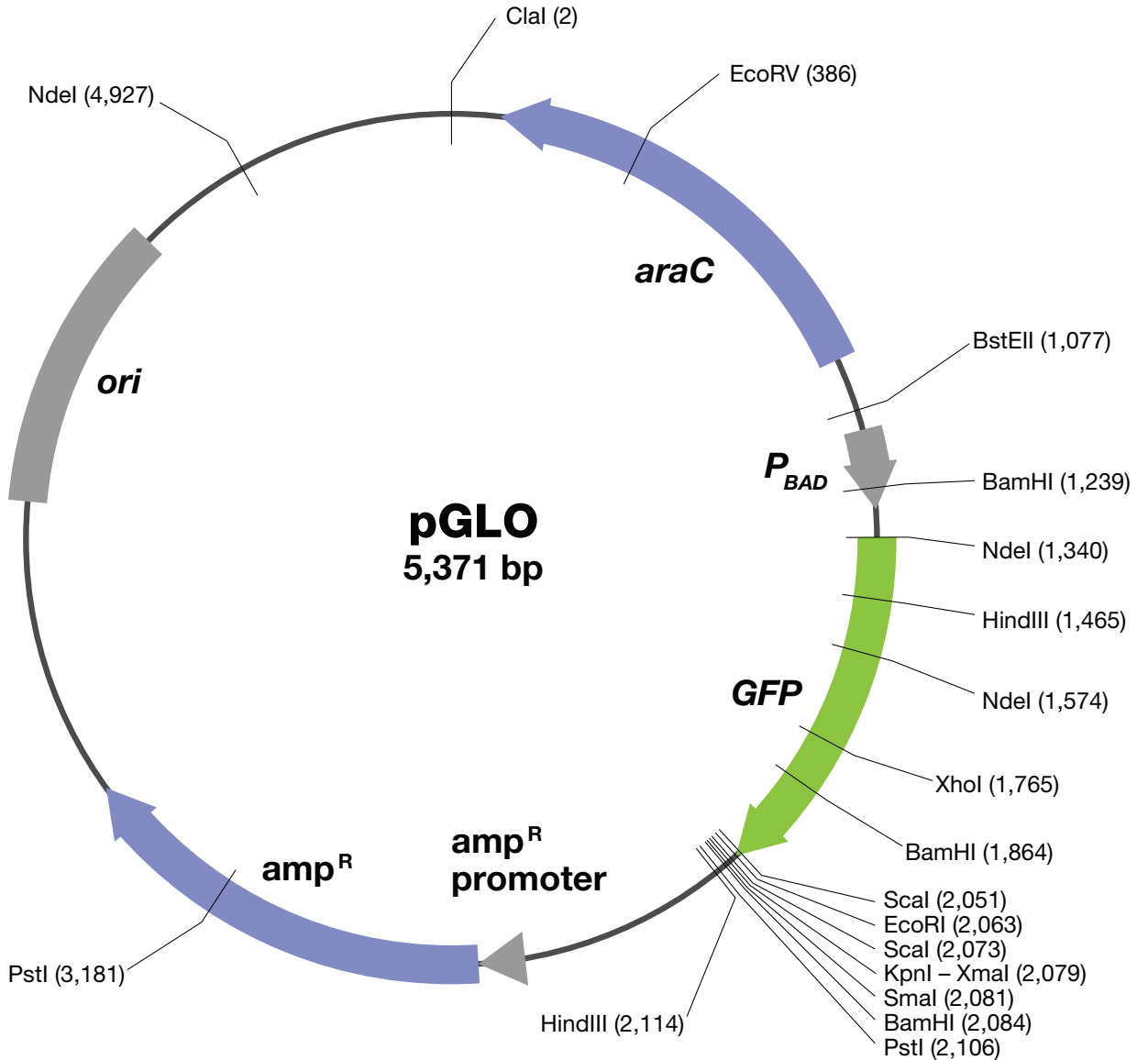


Plasmid Map

The features of the pGLO plasmid are illustrated in the plasmid map below. Specific sequences, including gene and promoter sequences, are shown as thick lines along the circle with arrowheads that point in the direction of operation or transcription. Short lines perpendicular to the circle indicate restriction sites and are labeled with the names of the associated restriction enzymes and the number of the nucleotide just before the cut site.



pGLO Plasmid Sequence

ATCGATGCATAATGTGCCTGTCAAATGGACGAAGCAGGGATTCTGCAAACCCATGCTACTCCGCTCAAGCCGTCAAATTTGCTGATTGCTTACCAATATGA
CAACTTGACGGCTACATCATTCACTTTTTCTTACAACCGGCACGGAACCTCGCTCGGGCTGGCCCCGGTGCATTTTTAAATACCCGCGAGAAATAGAGT
TGATCGTCAAACCAACATTGCGACCCGACGGTGGCGATAGGCATCCGGGTGGTCTCAAAGCAGCTTCGCCTGGCTGATACGTTGGTCTCGCGCC
AGCTTAAAGCCTAATCCCTAACTGCTGGCGGAAAAGATGTACAGACGCGACGGCGACAAGCAAACATGCTGTGCGACGCTGGCGATATCAAATTTG
CTGTCTGCCAGGTGATCGCTGATGACTGACAAGCCTCGCTACCCGATTCCATCCGTTGGATGGAGCGACTGTTAATCGCTTCCATGCGCCGAC
TAACAATTGCTCAAGCAGATTATCGCCAGCAGCTCCGAATAGCGCCCTTCCCTTGGCCGGCTTAATGATTTGCCAAACAGGTGCTGAAATGCGG
CTGGTGGCTTCAATCCGGGCGAAAGAACCCCGTATTGGCAAATATTGACGGCCAGTTAAGCCATTCAATGCCAGTAGGCGCGGACGAAAGTAAACCC
ACTGGTATACCAATTCGCGAGCCTCCGGATGACGACCGTAGTGATGAATCTCTCTGGCGGGAACAGCAAAATATCACCCGGTCCGCAACAAATTTCTC
GTCCCTGATTTTTACCACCCCTGACCCGCAATGGTGAATTTAGAATATAACCCCTTTCATCCAGCGGCTCGGTGATAAAAAATCGAGATAACCGTT
GGCCTCAATCGCGCTTAAACCCGCCACCAGATGGGCATTAACGAGTATCCCGGCAGCAGGGGATCATTTCGCTTACGCCATATTTTCACTACTCC
GCCATTAGAGAAAGAACCAATTGTCATATTGCATCAGACATTGCCGTCACTGCGCTTTTACTGGCTTCTCGCTAACCAACCCGGTAAACCCCGCTT
ATTAAGCATTCTGTAAACAAAGCGGGACCAAGCCATGACAAAAACCGGTAAACAAAGTGTCTATAATCACGGCAGAAAAGTCCACATTGATTATTTGCA
CGGGCTCACATTTGCTATGCCATAGCATTTTTATCCATAAGATTTTACCGGATCCAGGCTTTCATCGCTTTTATCGCAACTCTTACCGACCTGTT
TTTTGGGCTAGAAATAATTTGTTAACTTTAAGAAGGAGATACATATGGCTAGCAAAGGAGAAGAACTTTTCACTGGAGTTGCCCAATCTTGTGAT
TAGATGGTGTATTAATGGGCACAAATTTCTGTCACTGGAGAGGGTGAAGGTGATGCTACATACGGAAGCTTACCCCTTAAATTTATTTGCACTACTGGAA
AATACCTGTTCCATGGCCAACTTGTCACTACTTTCTTATGGTGTCAATGCTTTTCCGTTATCCGGATCATATGAAACGGCATGACTTTTTCAAGA
GTGCCATGCCCGAAGTTATGTACAGGAACGACTATATCTTTCAAAGATGACGGAACTACAAGACGCTGAGTGAAGTTTGAAGTGCATACCC
TTGTTAATCGTATCGAGTTAAAAGGTATTGATTTAAGAAGATGAAACATTCTCGGACACAACTCGAGTACAATACTCACACAATGTATACATCAC
GGCAGACAACAAAAGAAATGGAATCAAAGCTAACTTCAAATTCGCCACAACATTGAAGATGGATCCGTTCACTAGCAGACCATTATCAACAAAATACTC
CAATTGGCGATGGCCCTGTCTTTTACCAGACAACCTTACCTGTGACACAATCTGCCCTTTGAAAGATCCCAACGAAAAGCGTGACCACATGGTCC
TTCTTGAGTTTTTGAACCTGCTGGGATTACACATGGCATGGATGAGCTCTACAATAATGAAATTCGAGCTCGGTCCGGGACGACCCCGCAAACTTCC
TGCAGGCATGCAAGCTTGGCTGTTTTGGCGGATGAGAGAAGATTTTACGCTGATACAGATTAATCAGAACGCAAGCGGTCTGATAAAACAGAAATTT
GCCTGGCGGCGATAGCGCGGTGGTCCCACCTGACCCCATGCCAACTCAGAAGTGAACGCGCGTAGCGCGGATGGTAGTGTGGGGTCCCCCATGC
GAGAGTAGGGAATGCCAGGCATCAAATAAAACGAAAGGCTCAGTGAAGACTGGCCCTTTCGTTTTATCTGTTGTTTGTGGTGAACGCTCTCCTGA
GTAGCAAAATCCGCGGGGAGCGGATTTGAACGTTGCGAAGCAACGCGCCGAGGGTGGCGGGCAGGACCCCGCAAACTTCCGATGAGTAACTAAA
TTAAGCAGAAGGCCATCTGACGGATGGCTTTTTGCGTTTCTACAACCTCTTTGTTATTTTCTAAATACATTCAAATATGATCCGCTCATGAGACAATA
ACCTGATAAATGCTTCAATAATATGAAAAGGAAGAGATGAGTATTAACATTTCCGCTGTCGCCCTTATCCCTTTTTTGGCGCATTTTGCCTTCTGTT
TTTGCTCACCCAGAACGCTGGTGAAGTAAAAGATGCTGAAGACTGATGGGTGCACGAGTGGGTACACTGAACTGGATCTCAACAGCGGTAAAGATC
AAGGATCTAGGTGAAGATCCTTTTTGATAATCTCATGACCAAACTCCCTTAACTGAGTTTTTCCGTTCCACTCAGACCCCGTGAAGAAAGATCAAAG
AATCGGTCGCGCATACACTATTCTCAGAATGACTTGGTTGAGTACTACCAGTACAGAAAAGCATCTTACGGATGGCATGACAGTAAGAGAATTATG
CAGTGTGCCATAACCATGAGTGATAACACTGCGGCCAATTTACTTCTGACAACGATCGGAGGACCGAAGGAGCTAACCGCTTTTTTGCACAACATGGG
GGATCATGTAACCTCGCTTGTGTTGGGAACCGGAGCTGAATGAAGCCATACAAAACGACGAGCGTGACACCACGATGCCTGCAGCAATGGCAACAA
CGTTGCGCAAACTATAACTGGCGAATCTTACTCTAGCTTCCCGGCAACATTAATAGACTGGATGGAGCGGATAAAGTTGCAGGACCATTCTCGCG
CTCGGCCCTTCCGGCTGGCTGGTTATTGCTGATAAATCTGGAGCCGGTGGAGCTGGGCTCGCGGTATCATTGCAGCACTGGGGCCAGATGGTAAGC
CCTCCCGTATCGTAGTTATCTACACGACGGGAGTCAAGGCAACTATGGATGAACGAAATAGACAGATCGCTGAGATAGGTGCCTCACTGATTAAGCATTG
GTAACCTGTCAGACCAAGTTTACTCATATATACTTTAGATTGATTTACGCGCCCTGTAGCGGGCATTAAAGCGCGGGTGTGGGTGGTTACCGCGCAGCT
GACCGATAAGCCTTGCAGCGCCCTAGCGCCCGCTCTTCCCTTCCCTTTTCCCTTTTCCGCGCTTCCGCGGTTCCCGGTAACCTTCAAACT
GGGGGCTCCCTTTAGGGTCCGATTTAGTGTCTTACGGCACCTCGACCCCAAAAACCTTGAATTTGGGTGATGGTTACGTTAGTGGGCCATCGCCCTGAT
AGACGGTTTTTCCGCCCTTGGAGTCCAGCTTTAATAGTGGACTCTTGTCCAAACTGGAACAACACTCAACCCATCTCGGGCTATTCTTTT
GATTTATAAGGGATTTTGGCGATTTCCGGCTATTGGTTAAAAAATGAGCTGATTTAAACAAAAATTAACGCGAATTTAAACAAAATTTAACGTTTACAATTTAA
AAGGATCTAGGTGAAGATCCTTTTTGATAATCTCATGACCAAACTCCCTTAACTGAGTTTTTCCGTTCCACTCAGACCCCGTGAAGAAAGATCAAAG
GATCTTCTGAGATCCTTTTTTCTGCGCGTAACTGCTGCTTGCACAACAAAAAACCCCGCTACCAGCGGTGGTTTTGTTGCCGGATCAAGAGCTACC
AATCTTTTTCCGAAGGTAACCTGGCTTACAGCAGAGCGCAGATACCAAACTACTGCTTCTAGTGTAGCCGATGTTAGGCCACCATTCAAGAATCTGTGA
GCACCGCTACATACCTCGCTCTGCTAATCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTGTCTTACCGGGTGGACTCAAGACGATAGTTA
CCGGATAAGCGCAGCGGTGCGGCTGAACGCGGGGTTCTGTCACACAGCCAGCTTGGAGCAGCAACCTGAGTGGAGCAGCAACCTGAGATACCTACAGCT
GAGCATTGAGAAAGCGCCACGCTTCCGAAGGGAGAAAGGCGGACAGGTATCCGGTAAGCGGCAGGGTCCGAACAGGAGAGCGCACAGGGGAGCT
TCCAGGGGAAACGCTTGTATCTTATAGTCTGTCGGGTTTCCGCCACTCTGACTTGGAGCTGATTTTTGTGATGCTCGTACAGGGGGCGGAGCCCT
ATGAAAAACGCCAGCAACGCGGCCCTTTTACGGTCTCTGGCCTTTTGTGGCCTTTTGTCTCACATGTTCTTCTCGCTTATCCCTGATTCTGTGGATA
ACCGTATTACCGCCTTTGAGTGAAGTGTGATACCGCTCGCCGACGCGAAGCAGCCAGCGCAGCGAGTCACTGAGCGAGGAAGCGGAAGAGCGCTGA
TGCGGTATTTTCTCCTTACGCATCTGTGCGGTATTTACACCCGCATATGGTGCCTCTCAGTACAATCTGCTCTGATGCCGCATAGTTAAGCCAGTATACAC
TCCGCTATCGCTACGTGACTGGGTGATGGCTGCGCCCGGACACCCGCAACACCCGCTGACGCGCCCTGACGGGCTTGTCTGCTCCGGCATCCGC
TTACAGACAAGCTGTGACCGTCTCCGGGAGCTGCATGTGTGACAGGTTTCCACCTCATACCCGAAACGCGCGAGGCGAGCAAGGATGGCGCCCAA
CAGTCCCGCGCCAGGGGCTGCCACCATACCCAGCCGAAACAAAGCTCATGAGCCCGAAGTGGCGAGCCGATCTCCCTTCGGTGTGATGTC
GGCGATATAGCGCCAGCAACCGCACCTGTGGCGCCGGTGTGCGCGCCAGTGCCTCGCGCTAGAGGATCTAATTTCTCATGTTTACAGCTTATC

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23
France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050
Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670
The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

