



100

1 **CTCGAGC**GGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAAGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCGCTGGTGACCCGCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCAGTACATGACCTTATGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCGAGAAAGTTGGGGGAGGGGTCGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAAACTGGGAAAGTGAT
701 GTCGTGTAAGTGGCTCCGCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTT**BstEII (860)**
AgeI (857)NcoI (865) **BamHI (883)** **Acc65I (895)**
901 CTCTAGACTCGAGATGCATGATATCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGC
1001 TTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAAACAAGTTAAACAACAACAAATGCATTCAATTTTATGTTTCAGG
1101 TTCAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTCTAAAATACAGCATAGCAAAACTTAACTCCAAT
1201 CAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTTCTTTCA
1301 TGGAGTTTAAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTGCTTTCATTTCTTTATGTTTTAAATGCAGTACCTCCACATTCCCTTTTATGAAAA
1401 TATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTA
1501 GTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTG
1601 CACGCAGTTGCCGCGGGTTCGCGCAGGGCGAACTCCCGCCCGCAGGCTGCTCGCGGATCTCGGTATGGCCGGCCGGAGGCGTCCCGAAAGTTCGTG
116 V C N G A P D R L A F E R G W P Q E G I E T M A P G S A D R F N T
1701 GACACGACCTCCGACCACTCGCGTACAGTCTGTCAGGCGCGCACCCACAGCCAGGGTGTGTCCGGCACCACCTGGTCTGGACCGCGCTGA
82 S V V E S W E A Y L E D L G R V W V W A L T N D P V V Q D Q V A S I
1801 TGAACAGGGTACGTCTGTCGGGACACACCGCGAAGTCTGCTCCACGAAGTCCCGGGAGAACCAGCCGGTCCGAGTCCAGAACTGACCGCTCCGGC
49 F L T V D D R V V G A F D D E V F D R S F G L R D T W F E V A G A
1901 GACGTCGCGCGGGTGGACCCGGAACGGCACTGGTCAACTTGGCCATGATGGCCCTCTATAGTGAGTTCGTATTATACTATGCCATATACTATGCCGA
16 V D R A T L V P V A S T L K A M
2001 TGATTAATTGTCAATCCGGTTGCTTTGAATTAGCGGTGTTTTACAAACACCTAAAAAAGGGTTTAAAAGATACCTTTGAACCCTAAGAAGCCCGAGAA
2101 TTAGTCCGCTCAAACTCAAGGGGACAAATTCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCACCCACCAATGTGAACAAACTCCAA
2201 CGCCATTACATCCCTCCCCCGCCGACTAGCCGTCTCAAAGCCGAGGTGACTATTGCGGCCGATAGGACCACGGGTACAGGAAGCAGCAGCC
2301 GGTGAGGGACCAGGCCCTCTTCTTTGTGGGTGACTCACCCGCCGCTCCACCGGGTCCCGCTCCTCATTGAGCTCCTTGAACAGGGCCCGG
2401 GAGCGCCATCTTCCACGCACGCAACTGGTGCCGACGGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACGCGGCCGCCAAGCCAGATCGTGCC
2501 GGGTGTGGGGCCACATGGCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCCTCGCCCATCCAGGTGGCGTCCGACATGTGCT
2601 CCGAAGCGGGCGGGCCAGCCGCACTCCTGTCCCTCATTCTCCCAACCATGACTCTCCGGCTCCGGGCGAGCAAGCCCGCACCTCCCTTT
2701 GTTAGCCCTATTGCTGAACGGCAATCGAAGGAGCAGGGCAACAACAACAAAAAAGACCAGAGTGGCGCGGAGTAGCACGGCGGGCGGCGC
2801 GGACACCAGCTAGGCCTCAAGCCGACACGAGGCGAGGCTACGGGTTGCCGCTAGGCTCGCACTCTGCTCCCGCGCCGCCCAACTCGAAGCGGG
2901 AATGCTCGAGCTAATCCCCGCGACGACAGCGGGGCCCGGCTCGGAGCAGGACTCCAGCTCGGGCGCCGGGAAGCCACACCCGCCCTCAC
3001 TGCGTTCTGACGGCAAGCTTCGGCGAAGAAGCTCTGGCCCTGCGGGTCTGTTGGTCTTTATAGCCGCTCGGGTCCAGGCCCGCCCGGCAATCAG

3101 CGCCGCCCGCCCCGAGCCGCTCCTCCGGTGGGCGCGGGACCCGCCCTGCTGTGGGGAGGGGCGGCGCTGGAGGCCCTCGCGCGCTCTGGCGGAA
3201 CTAGTCGACGGACTGGGCTACGGGCGCCCCGAGAGGCGCAGCCAGAGGCCGCGCTAGGAAGGGGCGGCGCCGAGAACACGATCCCTCCCACCCCC
3301 CTCGGACGTGACTCGGACCACATCCCGGGGTCGCTAGGGCCCTCCCTTCTGCTCCTTCCCCAGCCTGGCGCGCTCTGGGGCGCCGTGACTCAGCCAG
3401 AATGTTGGCAATGGGGAGGGCGGAACGGGAAGTGGAGGACGCGGATGGAAAAGTCGAAACGAAGGAAGCTGAGTTTCGCCTGCAGGTTAATTAAGAA
3501 CATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAT
3601 CGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGC
3701 CGCTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTC
3801 CAAGCTGGGCTGTGTGCACGAACCCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCG
3901 CCACTGGCAGCAGCCACTGGTAAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAA
4001 GAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGG
4101 TTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTCTGACGCTCAGTGAACGAAAAC
4201 TCACGTTAAGGGATTTTGGTCATGGCTAGT