



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG  
101 AAACAAAACAAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA  
201 AATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTC  
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC  
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG  
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC  
601 AGAGCGCACATCGCCACAGTCCCGAGAAAGTTGGGGGAGGGGTCCGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGGTAACTGGGAAAGTGAT  
701 GTCGTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC  
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCCGGTACCATGGGAATTC AAGCTTGGATCCAGATCTGGTAC  
901 CTCTAGACTCGAGATGCATGATATCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGC  
1001 TTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAAACAAGTTAAACAACAACAAATGCATTCATTTTATGTTTCAGG  
1101 TTCAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTAACTCCAAT  
1201 CAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTGCCAATGTGCATTAGCTGTTGACGCTCACCTCTTTCA  
1301 TGGAGTTTAAGATATAGTGATTTTTCCAAGGTTTGAAGTACTCTTTCATTTCTTTATGTTTTAAATGCAGTACCTCCACATTCCCTTTTTAGTAAAA  
1401 TATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTA  
1501 GTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTCAGGTTAAGCTCCAGGCTTCTTGTGATGCA  
1601 CCAAGTCTTGGGCTTCTGGAACCTCAACATCAGCTGTACAGTGAATCCAGTCTTTTCATAAAAAAGGCAGGTTTCTGGGAGCAGAAGTTCCAGAAGG  
191 W T R P G E P V E V D A T V T F G L R E Y F P L N R P A S T E L F  
1701 GCAGGAACTCCAGCCCTTTCAGCAGTTC AACTCCAGGCAGAACAACAGCAGATCCCAGACCTTTCCCTGGTGGT CAGGGCTCACTCCAACAGTTGCCA  
157 A P V G A R E A A E V G P L V V A S G L G K G Q H D P S V G V T A L  
1801 GAAACCAAGCTGGCTCTTTGGCCTGTGTGGTCCAGCAGACCTCCATTTGTTGTTGTGCTCCAGAGCTGCTCCAGAGAGCTCAGCCATTCTTGGTCC  
124 F W A P E K P R H P A L L G E M Q Q Q A A L R S G S L E A M R P G  
1901 AATTTAGCAAAAACAGCACCAGCTTCAACAGACTCAGGTGTTGTCCAACAGTCAACAGCAGCTCCATCATCTGCAACCCAACTTTTCCAATGTCCAGT  
91 I E A F V A G A E V S E P T T W V A V A G D D A V W V K G I D L  
2001 CCCACTGGTGAGGAAGAGTCTTGCAGTTCCTCACCTCTCAATGTGCCTGT CAGGTCAAGTGTGCCTTGTG CAGGGTAGCTGCAAAAGCAG  
57 G V R T L F L E Q L E T V R E I H R D P D V T H R T A P Y D A F A A  
2101 CAGCCAGTGTCTCACAGCTCTTGAACATCATCTCTGGTTGCCAGCCTCACTGTGGGTTTGTACTCAGTCATGGTGGCCCTCTATAGTGAGTCGTATT  
24 A L T R V A R P V D D R T A L R V T P K Y E T M  
2201 ATACTATGCCGATATACTATGCCGATGATTAATGTCAATCCGTTGCTTTGAATTAGCGGTGGTTTTCAACACCTAAAAAAGGGTTAAAAGATACC  
2301 TTTGAACCGTAAGAAGCCCGAGAATTAGTCCGCTCAAACTCAAGGGGACAAATCCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCA  
2401 CCCACCAATGTGAACAACTCCAACGCCATTACATCCCCTCCCCCGCGGACTAGCCGTGCTCAAAAGCCCGAGGTGACTATTGCGGCCGATAGGAC  
2501 CACGGGGTACAGGAAGCAGCAGCCGGT GAGGGACCAGGCCCTTCTCTTTGTGTGGTGACTCACCCGCCGCTCCACCGGGCTGCCGCTCTCCATT  
2601 TTGAGTCTCTTGAACAGGGCCCGGAGCGCCATCTTTCCACGCACGCAACTGGTGCCGGACGGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACG  
2701 CGGCGCCCAAGCCAGATCGTGCCGGTGTGGGGCCACATGGCCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCCCTCGCCCC  
2801 ATCCAGTGGCGTCGACATGTGCTCCGAAGCGGGCGGGCCAGCCGCCACTCCTGTCCCTCCATTCTCCCAACCATGACCTCTCCGGCTCCGGG  
2901 CGAGCAAGCCCGCACCTCCCTTTGTTAGCCCTATTGCTGAACGGCAATCGAAGGCAGCAGGGCAACAACAACAAAAAAGACCAGAGTGCCG  
3001 CCGGAGTAGCACGCGCGGCGGCGGACACCACGCTAGGCCCTCAAGCCGACACGAGGCGAGGCTACGGGTTGCCGCTAGGCTCGCACTCTGCCTCC

3101 CGCGCCGCCGCAACTCGAAGCGGGAATGCTCGCAGCTAATCCCCGCCGACGACAGCGGGGCCGCGCTCGGAGCAGGACCTCCAGCTCGGGGCC  
3201 GCGGAAGCCACACCCGCCCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAGAACGTCTGGCCCTGCGGGTCGTTGTGGTCTTTATAGCCGCTCGG  
3301 CGTCAGGCCCGCCCGGCAATCAGCGCCGCCCGAGCCCGCTCTCCGGTGGGCGCGGGACCCGCCCTGCTGTGGGGAGGGCGCGCTG  
3401 GAGGCCCTCGCGCTCTGGCGGAAGTACTAGTACGACGGACTGGGCTACGGGCCGCCCGAGAGGCGCAGCCAGAGGCCGCGCTAGGAAGGGCGGGCGCC  
3501 GAGAACACGATCCCTCCCACCCCTCGGACGTGACTCGGACCACATCCCGGGTCGCTAGGGCCCTCCCTTCTGCTCCTTCCCCAGCTGGCGCG  
3601 CTCTGGGGCGCCGTGACTCAGCCAGAATGTTGGCAATGGGGAGGGCGGAACGGGGAAGTGGAGGACGCGGATGGAAAAGTCGAAACGAAGGAAGCTGA  
3701 GTTTCGCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCG  
3801 CCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTC  
3901 GTGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCATAGCTCACGCTGTAGGTATC  
4001 TCAGTTCGGTGTAGTTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTC  
4101 CAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGG  
4201 TGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCA  
4301 AACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGG  
4401 GTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGT