



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCCGCTGGTGACCCGCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGTACTGGCTCCGCTTTTTCCGAGGTTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCAGGTCACCATGGTGAAGGCGAGGAGCTGTTACCAGGG
120 M V S K G E E L F T G
901 GTGGTGCCATCTGGTTCGAGCTGGACGGCGACGTAACGGCCACAAGTTCAGCGTGTCCGGCAGGGCGAGGGCGATGCCACCTACGGCAAGCTGACCC
120 V V P I L V E L D G D V N G H K F S V S G E G E G D A T Y G K L T
1001 TGAAGTTCATCTGCACCACCGCAAGCTGCCGTGCCCTGGCCACCCTCGTACCACCTGACCTACGGCGTGCAGTGTTCAGCCGCTACCCCGACCA
145 L K F I C T T G K L P V P W P T L V T T L T Y G V Q C F S R Y P D H
1101 CATGAAGCAGCAGACTTCTTCAAGTCCGCCATGCCCGAAGGTACGTCCAGGAGCGACCATCTTCTTCAAGGACGACGGCAACTACAAGCCCGCGCC
178 M K Q H D F F K S A M P E G Y V Q E R T I F F K D D G N Y K T R A
1201 GAGGTGAAGTTCGAGGGCGACACCCTGGTGAACCGCATCGAGCTGAAGGGCATCGACTTCAAGGAGGACGGCAACATCCTGGGCGACAAGCTGGAGTACA
112 E V K F E G D T L V N R I E L K G I D F K E D G N I L G H K L E Y
1301 ACTACAACAGCCACAACGTCTATATCATGCGCCACAAGCAGAAGACGGCATCAAGTGAACCTCAAGATCCGCCACAACATCGAGGACGGCAGCGTGCA
145 N Y N S H N V Y I M A D K Q K N G I K V N F K I R H N I E D G S V Q
1401 GCTCGCCGACCACTACCAGCAGAACACCCCATCGGCGACGGCCCGTGTGCTGCCGACAACCACTACCTGAGCACCCAGTCCGCCCTGAGCAAAGAC
178 L A D H Y Q Q N T P I G D G P V L L P D N H Y L S T Q S A L S K D
1501 CCCAACGAGAAGCGGATCACATGGTCTGCTGGAGTTCGTGACCGCCCGGGATCACTCTCGGCATGGACGAGCTGTACAAGTAAAGCTAGCTGGCCA
212 P N E K R D H M V L L E F V T A A G I T L G M D E L Y K
1601 GACATGATAAGATACATTGATGAGTTTGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTG
1701 TAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTA
1801 AAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATA
1901 AGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAAGATATAGTGTATTTTCCAAGGTTTGA
2001 ACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAA
2101 ATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACCTTTAATAGA
2201 AATTGGACAGCAAGAAAGCGAGCTTCTAGCTCAGGTTTAAAGCTCCAGGCTTCTTGTGCATGCACCAAGTTCTTGGGCCTTCTGGAACCTCAACATCAGCT
200 A G P K R T M C W T R P G E P V E V D A
2301 GTCACAGTGAATCCAGTCTTTTATAAAAAGCAGGTTTCTGGGAGCAGAAGTTTCCAGAAAGGCAGGAATCCAGCCCTTTCAGCAGCTTCAACTCCAG
178 T V T F G L R E Y F P L N R P A S T E L F A P V G A R E A A E V G P
2401 GCAGAACAAAGCAGATCCAGACCCTTCCCTGGTGGTCAGGGCTCACTCCAACAGTTGCCAGAAACCAAGCTGGCTTTTTGGCCTGTGGTGCCAG
145 L V V A S G L G K G Q H D P S V G V T A L F W A P E K P R H P A L
2501 CAGACCTTCCATTTGTTGTTGTGCTGCCAGCCTGCTTCCAGAGAGCTCAGCCATTCTGGTCCAATTTCCAGCAAAAACAGCACCAGCTTCAACAGACTCA
112 L G E M Q Q Q A A L R S G S L E A M R P G I E A F V A G A E V S E
2601 GGTGTTGTCAAACCTGCAACAGCAGCTCCATCATCTGCAACCCAAACTTTTCCAATGTCCAGTCCCCTCTGGTGAGGAAGAGTTCTTGCAGTTCTGTCA
78 P T T W V A V A A G D D A V W V K G I D L G V R T L F L E Q L E T V
2701 CCCTCTCAATGTGCTGTGAGGGTCACTGTGTGCTTGTGTCAGGGTGTGCTGCAAAAAGCAGCAGCCAGTGTCTCACAGCTCTTGGAAATCATCTCT
45 R E I H R D P D V T H R T A P Y D A F A A A L T R V A R P V D D R
2801 GGTGCCAGCCTCAGTGGGTTTGTACTAGTCTGTTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTC
12 T A L R V T P K Y E T M
2901 AATCCGGTTGCTTTGAATTAGCGGTGGTTTTCAACAACCTAAAAAGGGTTAAAAGATACCTTTGAACCGTAAGAAGCCCGAGAATTAGCTCCGCTC
3001 AAAACTCAAGGGGACAAATTCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAATGTGAACAACTCCAACGCCATTACATC
3101 CCCTCCCCCGCCGCGACTAGCCGTGCTCAAAAAGCCGAGGTGACTATTGCGCCGATAGGACCACGGGGTACAGGAAGCAGCAGCCGGTGAGGGACCA
3201 GGCCCTCTTCTTTGTGTGGGTGACTCACCCGCCGCTCCACCGGGTGCCTGCTCCATTTTGTGCTCTTGAACAGGGCCCGGGAGCGCCATCT

3301 TTCCACGCACGCAACTGGTGCCGGACGGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACGCGGCCGCCAAGCCAGATCGTGCCGGGTGCTGGGGC
3401 CACATGGCCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCCTCGCCCATCCAGGTGGCGTCGGACATGTGCTCCGAAGCGGGC
3501 GGGCCCCAGCCGCACTCCTGTCCCTCATTCTCCCAACCATGACTCTCCGGGCTCCGGGCGAGCAAGCCCCGCACCCTCCCTTTGTTAGCCCTAT
3601 TGCTGAACGGCAATCGAAGGCAGCAGGGCAACAACAACAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGCGCGGCGGACACCAGCT
3701 AGGCCTCAAGCCGGACACGAGGCGAGGCTACGGGGTTGCCGTAGGCCTCGACTCTGCCTCCCGCGCCGCCCAACTCGAAGCGGGAATGCTCGCAGC
3801 TAATCCCCGCGACGACAGCGGGGCCGCGCTCGGAGCAGGACCTCCAGCTCGGCGGCCGCGGAAGCCACACCCGCCCTCACCTGCGTTCTGACG
3901 GCAAGCTTCGGCGAAGAAGCTCTGGCCCTGCGGGTCGCTTGTGGTCTCTTATAGCCGCTCGGCGTCAGGCCCGCCCGCCAATCAGCGCCGCCGCC
4001 CGAGCCGCTCCTTCGGTGGGCGGGACCCGCCCTGCTGTGGGGAGGGGCGGCGCTGGAGGCCCTCGCGCTCTGGCGAACTAGT**CGACGGA**
4101 CTGGGTACGGCCGCCCCGAGAGGCGAGCCAGAGGCCGCTAGGAAGGGGCGGCGCCGAGAACACGATCCCTCCCCACCCCTCGGACGTGAC
4201 TCGGACCACATCCCGGGTTCGCTAGGGCCCTCCCTTCTGCTCTTTCCCCAGCCTGGCGCTCTGGGGCGCCGTGACTCAGCCAGAATGTTGGCAAT
4301 GGGGGAGGGCGGAACGGGGAAGTGGAGGACCGGATGGAAAAGTCGAAACGAAGGAAGCTGAGTTTCGCCTGCAGGTTAATTAAGAACATGTGAGCAA
4401 AGGCCAGCAAAGGCCAGGAACCGTAAAAAGCCGCTTGGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGT
4501 CAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGAT
4601 ACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTG
4701 TGTGCACGAACCCCGTTCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCA
4801 GCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTG
4901 GTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAACCCGCTGGTAGCGGTGTTTTTTGTTTG
5001 CAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTACGGGTCTGACGCTCAGTGAACGAAAACTCACGTTAAGGG
5101 ATTTTGGTCATGGCTAGT