



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGTAAGTGGCTCCGCTTTTCCGAGGTTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCAGGTCACCATGGTGAAGGCGAGGAGCTGTTACCAGGG
125 M V S K G E E L F T G
901 GTGGTGCCATCTGGTTCGAGCTGGACGGCGACGTAACGGCCACAAGTTACGCGTGTCCGGCAGGGCGAGGGCGATGCCACCTACGGCAAGCTGACCC
125 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
45 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 CATGAAGCAGCAGCTTCTTCAAGTCCGCCATGCCGAAAGGTACGTCCAGGAGCGACCATCTTCTTCAAGGACGACGGCAACTACAAGCCCGCGCC
78 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 ACTACAACAGCCACAACGTCTATATCATGCGCCACAAGCAGAAGAACGGCATCAAGTGAACCTCAAGATCCGCCACAACATCGAGGACGGCAGCGTGA
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 GCTCGCCGACCACTACCAGCAGAACACCCCATCGGCGACGGCCCGTGTGCTGCCGACAACCCTACCTGAGCACCAGTCCGCCCTGAGCAAAGAC
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 TAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTA
1801 AAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATA
1901 AGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTCCCAAGGTTTGA
2001 ACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATGAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAA
2101 ATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACCTTTAATAGA
2201 AATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTGCACGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCC
125 D Q E E A V F H V C N G A P D R L A F E
2301 CGCCCCACGGTGTCTCGCCGATCTCGGTATGCGCCGGCCGGAGGCGTCCCGGAAGTTCTGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCA
103 R G W P Q E G I E T M A P G S A D R F N T S V V E S W E A Y L E D L
2401 GGCCGCGACCCACCCAGGCGGTTGTCGGCACCACCTGGTCTGGACCGCGCTGATGAACAGGGTACAGTCCGTCGTCGCGGACCAACCGGCGAA
70 G R V W V W A L T N D P V V Q D Q V A S I F L T V D D R V V G A F
2501 GTCGTCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGATCAGAAGTCCGACCGTCCGGCGACGTCGCGCGGGTGGACCCGGAACGGCACTGGTC
37 D D E V F D R S F G L R D T W F E V A G A V D R A T L V P V A S T
2601 AACTTGGCCATGATGGCCCTCTATAGTGAAGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAATCCGGTTGCTTTGAATTAGCGGT
3 L K A M
2701 GGTTCCTACAACCTAAAAAGGGTTAAAAGATACCTTTGAACCGTAAGAAGCCCGAGAATTAGCTCCGCTCAAAACTCAAGGGGACAAATTCAAA
2801 AATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAATGTGAACAACTCCAACGCCATTACATCCCTCCCGCCGCGACTAGCCGT
2901 GCTCAAAAGCCGAGGTGACTATTGCGGCCGATAGGACCACGGGTACAGGAAGCAGCAGCCGGTGGAGGACCAGGCCCTCTTCTTTGTGGTGGTAC
3001 TCACCCGCGCTCCACCGGGTCCGCGTCTCCATTTTGTGCTCTTGAACAGGGCCCGGAGCGGCATCTTCCACGACGCAACTGGTGGCGGA
3101 CGGGATGGCTCACCTAGTTAGGAGGAGGCAACGCGGCCGCAAGCCAGATCGTCCGGTGTGGGGCCACATGGCTCGGACGCTAACCCC
3201 AGCCTGTTGCTTCGGGAAAAACCCAGGCTCGCCCATCCAGTGGCGTGGACATGTGCTCCGAAGGCGGGGCGCCAGCCGCACTCTGTCCC

3301 TCCATTCTCCCAACCATGACCTCTCCGGGCTCCGGGCGAGCAAGCCCCGACCCCTCCCTTTGTTAGCCCTATTGCTGAACGGCAATCGAAGGCAGCA
3401 GGGCAACAACAACAAAAAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGGCGGGCGGACACCACGCTAGGCCTCAAGCCGGACACGAGGCGA
3501 GGCTACGGGGTTGCCGCTAGGCCTCGCACTCTGCCTCCCGCGCCGCCCAACTCGAAGCGGGAATGCTCGCAGCTAATCCCCGCCGACGACAGCGGGGC
3601 CCGGCCGCTCGGAGCAGGACCTCCAGCTCGGCGGCCCGGGAAGCCACCCGCCCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAGAAGCTCTGG
3701 CCCTGCGGGTCGCTTGTGGTCTTTATAGCCGCTCGGCGTCAGGCCCGCCCGGCAATCAGCGCCCGCCCGAGCCCGCTCCTTCGGTGGGCGC
3801 GGGACCCCGCCCTGCTGTGGGGAGGGGCGGCCGCTGGAGGCCCTCGCGCTCTGGCGAACTAGT**CGACGGACTGGGCTACGGGCCGCCCGAGAG**
3901 **GCGCAGCCAGAGGCCGCTAGGAAGGGGCGGGCGCCGAGAACACGATCCCTCCCCACCCCTCGGACGTGACTCGGACCACATCCCGGGTTCGCTA**
4001 **GGGCCCTCCCTTCTGCTCCTTTCCCCAGCCTGGCGGCTCTGGGGCGCCGTGACTCAGCCAGAATGTTGGCAATGGGGAGGGCGGAACGGGAAGTGG**
4101 **AGGACCGGATGAAAAAGTCGGAACGAAGGAAGCTGAGTTTCGCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGCCAGCAAAAGCCAGGAACCGT**
4201 **AAAAAGCCCGTGTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGAC**
4301 **TATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGG**
4401 **AAGCGTGGCGTTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCCC**
4501 **GACCGCTGCGCCTTATCCGTAACCTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAG**
4601 **CGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGT**
4701 **TACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGCAAGCAGCAGATTACGGCAGAAAA**
4801 **AAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGT**