



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAACTAGCAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCCGCTGGTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGGAGGGTTCGGCAATTGAACCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTGCCCGCAGAACACAGCTGAAGCTTACCAGTGCACCATGGTCTGGGGCCCTGCATGCTGCTGCTGCT
1 M V L G P C M L L L L
901 GCTGCTGCTGGGCTGAGGCTACAGCTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCAGGACTTCTGGAACCGCAGGCGAGCCGAGGCCCTGGGT
11 L L L G L R L Q L S L G I I P V E E E N P D F W N R E A A E A L G
1001 GCCCCAAGAAGTGCAGCCTGCACAGACAGCCCAAGAACCTCATCATCTTCTGGGCGATGGGATGGGGGTGTCTACGGTGCAGCTGCCAGGATCC
45 A A K K L Q P A Q T A A K N L I I F L G D G M G V S T V T A A R I
1101 TAAAAGGCAGAAAGGACAACTGGGCGCTGAGATACCCCTGGCTATGAGCCGCTTCCCATATGTGGCTCTGTCCAAGACATACAATGTAGACAAACA
78 L K G Q K K D K L G P E I P L A M D R F P Y V A L S K T Y N V D K H
1201 TGTGCCAGACAGTGGAGCCACAGCCACGGCCTACCTGTGCGGGTCAAGGGCAACTTCCAGACCATTGGCTTGTGAGTGCAGCCGCCGCTTAAACCAGTGC
111 V P D S G A T A T A Y L C G V K G N F Q T I G L S A A A R F N Q C
1301 AACACGACACGCGCAACGAGGTCTATCCGTGATGAATCGGGCAAGAAAGCAGGGAAGTCAAGTGGGAGTGGTAACCACCACAGTGCAGCACGCT
145 N T T R G N E V I S V M N R A K K A G K S V G V V T T T R V Q H A
1401 CGCCAGCCGGCACCTACGCCACACGGTGAACCGCAACTGGTACTCGGACGCCGACGTGCCTGCCTCGGCCCGCAGGAGGGTGCAGGACATCGCTAC
178 S P A G T Y A H T V N R N W Y S D A D V P A S A R Q E G C Q D I A T
1501 GCAGCTCATCTCCAACATGGACATTGATGTGATCCTGGGTGGAGGCCGAAAGTACATGTTCCGATGGGAACCCAGACCCTGAGTACCCAGATGACTAC
211 Q L I S N M D I D V I L G G R K Y M F R M G T P D P E Y P D D Y
1601 AGCCAAGTGGGACAGGCTGGACGGGAAGAATCTGGTGCAGGAATGGCTGGCGAAGCGCCAGGGTGGCCGGTATGTGTGGAACCGCACTGAGCTCATGC
245 S Q G G T R L D G K N L V Q E W L A K R Q G A R Y V W N R T E L M
1701 AGGCTTCCCTGGACCCGCTGTGACCCATCTCATGGGTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACACTGGACCCCTCCT
278 Q A S L D P S V T H L M G L F E P G D M K Y E I H R D S T L D P S L
1801 GATGGAGATGACAGAGGCTGCCCTGCGCTGCTGAGCAGGAACCCCGCGCTTCTTCTCTCGTGGAGGGTGGTCGCATCGACCACGGTATCACGAA
311 M E M T E A A L R L L S R N P R G F F L F V E G G R I D H G H H E
1901 AGCAGGGCTTACCGGCACTGACTGAGACGATCATGTTGACGACGCCATTGAGAGGGCGGGCCAGCTCACCAGCGAGGAGGACAGCTGAGCCTCGTCA
345 S R A Y R A L T E T T I M F D D A I E R A G Q L T S E E D T L S L V
2001 CTGCCGACACTCCACGCTTCTCCTCGAGGCTACCCCTGCGAGGGAGCTCCATCTCGGGCTGGCCCTGGCAAGCCCGGGACAGGAGGCTCA
378 T A D H S H V F S F G G Y P L R G S S I F G L A P G K A R D R K A Y
2101 CACGGTCTCCTATACGGAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCCGCGGATGTTACCAGAGCGAGAGCGGGAGCCCGAGTATCGGCAG
411 T V L L Y G N G P G Y V L K D G A R P D V T E S E S G S P E Y R Q
2201 CAGTCAGCAGTCCCCTGGACGAAGAGACCCACGACGGCAGGACGTGGCGGTGTTCCGCGCGGCCCGCAGGCGACCTGGTTCACGGCGTGCAGGAGC
445 Q S A V P L D E E T H A G E D V A V F A R G P Q A H L V H G V Q E
2301 AGACCTTCATAGCGCAGTCAATGGCCTTCCGCGCTGCTGGAGCCCTACACCGCCTGCGACCTGGCGCCCCCGCGGCACCACCGACGCGCGCACCC
478 Q T F I A H V M A F A A C L E P Y T A C D L A P P A G T T D A A H P
2401 GGGCGGTCCCGTCCAAGCTCTGGATTGAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACATAAGTGCAGTGAAA
511 G R S R S K R L D •
2501 AAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAAATGCATTATTTTATG
2601 TTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACCTTAAACC
2701 TCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCT
2801 TCTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAACAGTCTTCATTCTTTATGTTTAAATGCACTGACCTCCACATTCCTTTTTT
2901 AGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTCATAATATCCCC
3001 CAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTTGCCA
125 D Q E E A V
3101 CAAAGTGCACGAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGTGTCTCGCCGATCTCGGTCATGGCCGGCCGGAGGCGTCCCGGAA
118 F H 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
3201 GTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACACCCAGGCCAGGGTGTGTCGGGACCACTGGTCTGGACC
85 N T 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

3301 GCGCTGATGAACAGGGTCACGTCGTCCCGACCACACCGGGAAGTCGTCCTCCACGAAGTCCCGGGAGAACCCGAGCCGGTCGGTCCAGAACTCGACCG
51 A S I 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
3401 CTCGGGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGGCCCTCCTATAGTGAGTGTATTATACTATGCCGATATACT
18 G A V D R A T L V P V A S T L K A M
3501 ATGCCGATGATTAATTGTCAATCCGGTTGCTTGAATTAGCGGTGGTTTTCAACACCTAAAAAGGGTTTAAAAGATACCTTTGAACCGCTAAGAAGC
3601 CCGAGAATTAGCTCCGCTCAAACTCAAGGGGACAAATTCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAAATGTGAACAA
3701 ACTCCAACGCCATTACATCCCCTCCCCCGCCGACTAGCCGTGCTCAAAGCCCGAGGTGACTATTGCGGCCGATAGGACCACGGGTACAGGAAGC
3801 AGCAGCCGGTGAGGGACCAGGCCCTTCTCTTTGTGTGGTGACTCACCCGCCGCTCCACCGGGTCCCGCGTCTCCATTTGAGCTCCTTGCAACAG
3901 GGCCCGGAGCGGCCATCTTTCACGCACGCAACTGGTGCCGGACGGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACGCGGCCGCCAAGCCAGA
4001 TCGTGCCGGGTGCTGGGGCCACATGGCCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCTCGCCCATCCAGGTGGCGTGGAC
4101 ATGTGCTCCGAAGCGGGCGGGCCAGCCGCCACTCTGTCCCTCATTCTCCCAACCATGACCTCTCCGGGCTCCGGGCAGCAAGCCCGCACCC
4201 TCCCTTTGTTAGCCCCTATTGCTGAACGGCAATCGAAGGCAGCAGGGCAACAACAACAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGGC
4301 GCGGCGCGACACCACGCTAGGCCTCAAGCCGGACACGAGCGAGGCTACGGGTTGCCGCTAGGCCTCGACTCTGCCTCCCGCGCCGCCGCAACTCG
4401 AAGCGGAATGCTCGCAGCTAATCCCCGCCGACGACAGCGGGGCCGGCCGCTCGGAGCAGACCTCCAGCTCGGCGGCCGCGGAAGCCACACCCGCC
4501 CCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAGAAGCTCTGCCCTGCGGGTGCCTGTGGTCTCTTATAGCCGCTCGGCGTCAGGCCGCCCGGC
4601 CAATCAGCGCCGCCGCCGAGCCCGCTCTTCCGGTGGCGCGGGACCCCGCCCTGCTGTGGGGAGGGCGGCCGCTGGAGGCCCTCGCGCGCTCT
4701 GCGGAACTAGTCGACGGACTGGGTACGGGCCGCCCGAGAGGCGCAGCCAGAGCGCGCTAGGAAGGGCGGGCGCCGAGAACACGATCCCTCCC
4801 CACCCCTCGGACGTGACTCGGACCACATCCCGGGTGCCTAGGGCCCTCCTTCTGCTCTTTCCCCAGCTGGCGGCTCTGGGGCGCGGTGACT
4901 CAGCCAGAATGTTGGCAATGGGGGAGGGCGGAACGGGGAAGTGGAGGACGCGGATGGAAAAGTCGAAAACGAAGGAAGCTGAGTTTCGCTGCAGTTAA
5001 TTAAGAACATGTGAGCAAAAGCCAGCAAAAGCCAGGAACCGTAAAAAGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCA
5101 CAAAAATCGACGCTCAAGTCAGAGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCG
5201 ACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGTTTTCTCATAGCTACGCTGTAGGTATCTCAGTTCGGTGTAGGTCC
5301 TTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTACGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTCCAACCCGGTAAGACACGA
5401 CTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTAC
5501 ACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACACCCGCTGGTA
5601 GCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAA
5701 CGAAAACCTACGTTAAGGGATTTGGTCATGGCTAGT