



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCCGCTGGTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGGAGGGTTCGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGTAAGTGGCTCCGCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCAGGTCACCATGGTTCTGGGCGCCTGCATGCTGCTGCTGCTG
120 M V L G P C M L L L L
901 CTGCTGCTGGCCTGAGGCTACAGCTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCAGGACTTCTGGAACCGCGAGGCAGCCGAGGCCCTGGGTG
120 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 CCGCAAGAAGCTGCAGCCTGCACAGACAGCCGCAAGAACCTCATCATCTTCTGGGCGATGGGATGGGGGTGTCTACGGTACAGCTGCCAGGATCCT
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 L
1101 AAAAGGGCAGAAGAAGGCAAACTGGGGCTGAGATACCCCTGGCTATGGACCGTTCATATGTGGCTCTGTCCAAGACATACAATGTAGACAAACAT
78 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 GTGCCAGACAGTGGAGCCACAGCCACGGCTACCTGTGCGGGTCAAGGGCAACTCCAGACCATTGGCTTGTGAGTGCAGCCGCGCTTAAACAGTGA
112 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 A R F N Q C
1301 ACACGACACGCGCAACGAGGTATCTCCGTGATGAATCGGGCCAAGAAAGCAGGGAAGTCAAGTGGGAGTGGTAACCACCACAGTGCAGCACGCTC
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 S
1401 GCCAGCCGGCACCTACGCCACAGGTGAACCGCAACTGGTACTCGGACGCCAGCTGCCTGCCTCGGCCCGCAGGAGGGTGCAGGACATCGCTACG
178 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 CAGCTCATCTCCAACATGGACATTGATGTGATCCTGGTGGAGGCCGAAAGTACATGTTTCGCATGGGAACCCAGACCCTGAGTACCAGATGACTACA
212 Q L I S N M D I D V I L G G G R K Y M F R M G T P D P E Y P D D Y
1601 GCCAAGGTGGGACAGGCTGGACGGGAAGAATCTGGTGCAGGAATGGTGGCGAAGCGCCAGGGTGCCTGGTGTGGAACCGCACTGAGCTCATGCA
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 Q
1701 GGCTTCCCTGGACCCGTCTGTGACCCATCTCATGGTCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACACTGGACCCCTCCCTG
278 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 ATGGAGATGACAGAGGCTGCCCTGCGCCTGCTGAGCAGGAACCCCGGGCTTCTTCTCTCTGTTGGAGGGTGGTGCATCGACCACGGTATCACGAAA
312 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 GCAGGGCTTACCGGCACTGACTGAGACGATCATGTTGACGACGCCATTGAGAGGGCGGGCCAGCTCACCAGCGAGGAGGACACGCTGAGCCTCGTAC
345 S R A Y R A L T E T I M F D D A I E R A G Q L T S E E D T L S L V T
2001 TGCCGACCCTCCACGCTTCTTCTCGGAGGCTACCCCTGCGAGGAGCTCCATCTTCCGGCTGGCCCTGGCAAGGCCGGACAGGAGGCGCTAC
378 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 R A K A Y
2101 ACGGCTCTCTATACGAAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCCGGGCGGATGTTACCGAGAGCGAGAGCGGGAGCCCGAGTATCGGCAGC
412 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 AGTCAGCAGTGCCTGGACGAAGAGACCCACGAGGCGAGGAGCTGGCGGTGTTGCGCGCGCCCGCAGGCGACCTGGTTCACGGCGTGCAGGAGCA
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 Q
2301 GACCTTATAGCGCAGTGCATGGCTTCCGCGCTGCTGGAGCCCTACACCGCTGCGACCTGGCGCCCCCGCGGACACCACGCGCCGCGACCCG
478 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 GGGCGTCCCGTCCAAGCGTCTGGATTGAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACACTAGAATGCAGTAAAA
512 G R S R S K R L D •
2501 AAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTATTTTATGT
2601 TTCAGGTTTCAAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCT
2701 CCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTT
2801 CTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGTGTGAACTAGCTCTTCATTTCTTTATGTTTAAATGCACTGACCTCCACATTCCCTTTTAA
2901 GTAAAATATTAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTACATAATATCCCC
3001 AGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTCAGGTTTAAAGCTCCAGGCTTCTTGT
200 A G P K R T
3101 CATGCACCAAGTCTTGGGCTTCTGGAACCTCAACATCAGCTGTACAGTGAATCCAGTCTTTCATAAAAAGGAGGTTTCTGGGAGCAGAAGTTCC
193 M C 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
3201 AGAAAGGCAAGAACTCCAGCCCTTTCAGCAGCTTCACTCCAGGCAAGCAACAGCAGATCCAGACCCTTCCCTGGTGGTCAAGGCTCACTCCAACAG
159 L F 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

3301 TTGCCAGAAACCAAGCTGGCTCTTTTGGCTGTGTGGTCCAGCAGACCTTCCATTTGTTGTTGCTGCCAGCCTGCTTCCAGAGAGCTCAGCCATTCT
126 A L 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
3401 TGGTCCAATTTTCAGCAAAAACAGCACCAGCTTCAACAGACTCAGGTGTTGTCCAAACTGCAACAGCAGCTCCATCATCTGCAACCCAAACTTTTCCAATG
93 P G I E A F V A G A E V S E P T T W V A V A A G D D A V W V K G I
3501 TCCAGTCCCACCTCTGGTGAGGAAGATTCTTGCACTTCTGTCACCCTCTCAATGTGCTGTCCAGGGTCAACTGTGCCTTGTTCAGGGTAGTCTGCAA
59 D L 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
3601 AAGCAGCAGCCAGTGTCTCACAGCTCTTGGAAACATCATCTCTGGTTGCCAGCCTCACTGTGGGTTTGTACTCAGTCATGGTGGCCCTCTATAGTGAGT
26 A A A L T R V A R P V D D R T A L R V T P K Y E T M
3701 CGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAATCCGGTTGCTTTGAATTAGCGGTGGTTTTCACAAACCTAAAAAGGGTTAAAA
GATACCTTTGAACCGCTAAGAAGCCCGAGAATTAGCTCCGCTCAAAACTCAAGGGGACAAATTCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAG
3901 TCTCCACCACCAAATGTGAACAAACTCCAACGCCATTACATCCCTCCCCCGCCGCGACTAGCCGTGCTCAAAGCCCGAGGTGACTATTGCGGCCGA
4001 TAGGACCACGGGGTACAGGAAGCAGCAGCCGGTGAGGGACCAGGCCCTTCTCTTTGTGTGGTGACTACCCGCCCGCTCCACCGGGTCCCGCGTCC
4101 TCCATTTTGTAGCTCCTTGAACAGGGCCCGGAGCGGCCATCTTCCACGCACGCAACTGGTGCCGGACGGGATGGCCTCACCTAGTTAGGGAGGACGG
4201 GCAACGCGGGCCGCAAGCCAGATCGTGCCGGGTGCTGGGGCCACATGGCCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCCT
4301 CGCCCATCCAGGTGGCTCGGACATGTCTCCGAAGCGGGCGGGCCCGCCAGCCGCACTCCTGTCCCTCCATTCTCCCAACCATGACCTCTCCGGGC
4401 TCCGGGCGAGCAAGCCCGCACCTCCCTTTGTAGCCCTATTGCTGAACGGCAATCGAAGGCAGCAGGGCAACAACAACAAAAAAGACCAGA
4501 GTGCGCCCGAGTAGCACGCGCGGGCGGGACACCACGCTAGGCCCAAGCCGGACAGGGCGAGGCTACGGGGTTGCCGCTAGGCCTCGCACTCT
4601 GCCTCCCGCGCCCGCAACTCGAAGCGGAATGCTCGCAGCTAATCCCGCCGACGACAGCGGGGCCCGCCGCTCGGAGCAGGACCTCCAGCTCGG
4701 CGGCCCGGAAGCCACACCCGCCCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAGAAGCTTGCCCTGCGGGTCTGTGGTCTTTATAGCCG
4801 CGTCGGGTCAGGCCCGCCCGCAATCAGCGCCGCCCGCCGAGCCGCTCTTCCGGTGGGCGGGACCCCGCCCTGCTGTGGGGAGGGGGCGG
4901 CGCTGGAGGCCCTCGCGGCTCTGGCGGAAGTCTGACGGACTGGGCTACGGGCCCGCCCGAGAGGGCGAGCCAGAGGCCGCGCTAGGAAGGGGGC
5001 GGGCGCGAGAACACGATCCCTCCCCACCCCTCGGACGTGACTCGGACCACATCCCGGGTCTGCTAGGGCCCTCCCTTCTGCTCTTTCCCCAGCCT
5101 GGCAGCTCTGGGGCGCGTACTCAGCCAGAATGTTGCAATGGGGGAGGGCGGAACGGGGAAGTGGAGGACCGGATGGAAAAGTCGGAACGAAGGA
5201 AGCTGAGTTTCGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGGTTTTTCCATAG
5301 GCTCCGCCCCCTGACGAGCATCAAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGC
5401 TCCCTCGTGCCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTTCTCCCTTCGGGAAGCGTGGCGTTTCTCATAGCTCAGCTGTA
5501 GGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGTGGGCTGTGTGCACGAACCCCGTTAGCCCGACCGCTGCGCTTATCCGGTAACTATCGTCT
5601 TGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTG
5701 AAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGAT
5801 CCGGCAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTC
5901 TACGGGGTCTGACGCTCAGTGGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGT