



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACG
101 AAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGGC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGACTGGCTCCGCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCCGGTACCATTGAAATCAAGGTGCTGTTTGCCTCATCTGT
120 I A V A E A K P T E I N E D L N I A A V A S N F A T T D L E T D L
1001 TCACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGTGCCTGGCAAAAACCTCCCCCAGA
45 F T N W E T M N V I S T D T E Q V N T D A D R G K L P G K K L P P D
1101 TGTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTTGCCCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAA
78 V L R E L E A N A R R A G C T R G C L I C L S H I K C T P K M K K
1201 TTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTTCA
112 F I P G R C H T Y E G E K E S A Q G G I G E A I V D I P E I P G F
1301 AGGATAAGGAGCCACTGGACAGTTTATTGCTCAAGTGGACCTCTGTCTGATTGCACCACTGGCTGTCTGAAGGGCTTCCAATGTCCAGTCTGTA
145 K D K E P L D Q F I A Q V D L C A D C T T G C L K G L A N V Q C S D
1401 CCTCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTTTCCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGACAGATGAGCT
178 L L K K W L P Q R C T T F A S K I Q G R V D K I K G L A G D R •
1501 AGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTG
1601 CTTTATTTGTAACCATTATAAGCTGCAATAAACAGTTAACAACAACAATTGCATTATTTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGGTTTTTTA
1701 AAGCAAGTAAACCTCTACAATGTGGTATGGAATTCTAAATACAGCATAGCAAACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAG
1801 GGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGATTTTCCC
1901 AAGGTTTGAAGTACTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTATGTAATAATTAATACATCATT
2001 GCAATGAAAATAAATGTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAACC
2101 TTTAATAGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCTCAGGTTAAAGCTCCAGGCTTCTTGTGCATGCACCAAGTTCTTGGGCTTCTGGAACCTCA
200 • A G P K R T M C W T R P G E P V E
2201 ACATCAGCTGTACAGTGAATCCAGTCTTTTCAAAAAGGCAGGTTTCTGGGAGCAGAAGTTTCCAGAAAGGCAGGAAGTCCAGCCCTTTCAGCAGCTT
181 V D A 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
2301 CAACTCCAGGCAGAAACAAGCAGATCCAGACCTTTCCCTGGTGGTCAGGGCTCACTCCAACAGTTGCCAGAAACCAAGCTGGCTCTTTTGGCCTGTG
148 V G P 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
2401 TGGTGCCAGCAGACTTCCATTTGTTGTGCTGCCAGCCTGTTCCAGAGAGCTCAGCCATTCTTGGTCCAATTTCCAGAAAACAGCACCAGCTTCA
115 P A L 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
2501 ACAGACTCAGGTGTTGTCAAACTGCAACAGCAGCTCCATCATCTGCAACCCAACTTTTCCAATGTCAGTCCCCTCTGGTGGGAAGAGTTCTTGCA
81 V S E 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
2601 GTTCTGTACCCTCTCAATGTGCTGTGAGGTTCAACTGTGCTGTTTGCAGGGTAGTCTGAAAAGCAGCAGCCAGTGTCTCACAGCTCTTGGAAC
48 E T V 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
2701 ATCATCTCTGGTTGCCAGCCTCACTGTGGGTTTGTACTCAGTCATGGTGGCCCTCTATAGTGAGTCGATTATACTATGCCGATATACTATGCCGATGA
15 D D R T A L R V T P K Y E T M
2801 TTAATTGCAATCCGGTTGCTTTGAATTAGCGGTGGTTTTCCAAACCTAAAAAGGGTTAAAAGATACCTTTGAACCGCTAAGAAGCCCGAGAATTA
2901 GCTCCGCTCAAACTCAAGGGGACAAATCCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAATGTGAACAACTCCAACGC
3001 CATTACATCCCCTCCCCCGCCGACTAGCCGTGCTCAAAAGCCGAGGTGACTATTGCGGCCGATAGGACCACGGGTTCACAGGAAGCAGCAGCCGGT
3101 GAGGGACCAGGCCCTTCTCTTTGTGTGGTGACTCACCCGCCGTCACCCGGGTGCCCGTCTCCATTTTGTGCTCTTGAACAGGGCCCGGGAG
3201 CGGCCATCTTCCACGCACGCAACTGGTGGCGACGGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACGCGCGCCGCAAGCCAGATCGTGCCGGG

3301 TGCTGGGGCCACATGGCCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCCTGCCCATCCAGGTGGCGTCGGACATGTGCTCCG
3401 AAGGCGGGCGGGCCCCAGCCGCACTCTGTCCCTCCATTCTCCCAACCATGACCTCTCCGGGCTCCGGGCGAGCAAGCCCCGACCCCTCCCTTTGTT
3501 AGCCCCTATTGCTGAACGGCAATCGAAGGCAGCAGGGCAACAACAACAAAAAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGCGGGCGGGGA
3601 CACCACGCTAGGCCTCAAGCCGACACGAGGCGAGGCTACGGGGTTGCCGCTAGGCCTCGACTCTGCCTCCCGCGCCGCCGCAACTCGAAGCGGGAAT
3701 GCTCGCAGCTAATCCCCGCCGACGACAGCGGGCCCGCCGCTCGGAGCAGGACCTCAGCTCGGCGGCCGGAAGCCACACCCGCCCTCACCTGC
3801 GTTCTGACGGCAAGCTTCGGCGAAGAACGTCTGGCCCTGCGGGTCGCTTGTGGTCTTTATAGCCGCTCGGCGTCAGGCCCGCCCCGGCCAATCAGCGC
3901 CGCCCGCCCGAGCCCGCTCCTCCGGTGGGCGGGACCCCGCCCTGCTGTGGGGAGGGCGGCCGCTGGAGGCCCTCGCGCTCTGGCGGAACTA
4001 GTCGACGGACTGGGTACGGGCCCCCCGAGAGGCGCAGCCAGAGGCCGCTAGGAAAGGGCGGGCGCGAGAACACGATCCCTCCCCACCCCTC
4101 GGACGTGACTCGGACCACATCCCGGGTTCGCTAGGGCCCTCCCTTCTGCTCCTTTCCCCAGCCTGGCGGCTCTGGGGCGCCGTACTCAGCCAGAAT
4201 GTTGGCAATGGGGAGGGCGGAACGGGGAAGTGGAGGACCGGATGGAAAAGTCGAAACGAAGGAAGCTGAGTTTCGCTGCAGGTTAATTAAGAACAT
4301 GTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGA
4401 CGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCCTGCCG
4501 TTACCGGATACCTGTCCGCTTTCTCCCTTCGGAAGCGTGGCGCTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAA
4601 GCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCA
4701 CTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAA
4801 CAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGATTGGTAGCTCTTGATCCGGCAAACAACCACCGTGGTAGCGGTGGTTT
4901 TTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCA
5001 CGTTAAGGGATTTTGGTCATGGCTAGT