



1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATGCCACAGTCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGGTAAACTGGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCCCGCCCTACCTGAGGCC
301 GCCATCCACGCCGTTGAGTTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTGCCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCGCTACCTGAGATCACCGGTACCCTGCTACCCTGCGTGTCCACTCCTGGTGGCTCTCGT
601 CCTTCTTGTGTGGAATTGAGACCTCTGATGCAGGTCCTATGGTGCAATGTGGAAGACAGTATCTGCTGCCAGGACTACATCCGTACCCTCTGCCA
13▶ L L A V A I Q T S D A G P Y G A N V E D S I C C Q D Y I R H P L P
701 TCACGTTTGTGAGGAGTTCTTCTGGACCTCAAATCCTGCCGAAGCCTGGCGTTGTTTTGATAACCGTCAAGAACCAGATATCTGTGCCGATCCCA
47▶ S R L V K E F F W T S K S C R K P G V V L I T V K N R D I C A D P
801 GGCAGGTCTGGTGAAGAAGCTACTCCATAAAGTCTAGGGAGGAGGACCTGATGACCATGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTT
80▶ R Q V W V K K L L H K L S •
901 TGGACAAACCACAAGTGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAA
1001 GTTAAACAACAATTGCATTATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAAT
1101 TCTAAAATACAGCATAGCAAACTTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTGGCC
1201 AATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTCTTCAATTTCTTTATGTTTT
1301 AAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATC
1401 CAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTC
1501 TAGCTTTAGTTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGTTCGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCAT
141▶ • N R T Y K L P I L E E I T T K V L K G N M E I L V F C D P A Y
1601 AGTCAGAGATGAGCTCTGACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTG
109▶ D S I L E R C M G C P S V V R I S R D V E D S Y P H R V A V I T D
1701 AAAGTCTTCTGCCGTTGCTCAGCAGACCAATGGCAATGGCTCAGCACAGACAGTACCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATG
76▶ F D K Q G N S V A S G I A I A E A C V T V R G I Y A E I H V A S I
1801 ATCTCCAGTCTTGGTCTGATGGCCGCCGACATGGTGTGTTGCTCCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGAT
42▶ I E G T K T R I A A G V H H K N D E Y L M T I K E T A V E V L E L D
1901 CCTGCTGAGAGATGTTGAAGTCTTCTGTTGGCCCTCTATAGTGTGCTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGC
9▶ Q Q S I N F T K M
2001 GTGGATGGCGTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCTACCGCCATTGCGTCAATGGGGC
2101 GGAGTTGTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCA
2201 AACCGCTATCCACGCCATTGATGACTGCCAAAACCGCATCATGGAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATA
2301 AGGTCATGACTGGGCATAATGCCAGCGGGCCATTTACCGTATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCCAAGTG
2401 GGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTACTATGGGAACATACGTATTATTGACGTCAATGGGGGGGG
2501 TCGTTGGCGGTGAGCCAGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAA
2601 CCGTAAAAAGGCCGCTTGTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAGAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACA
2701 GGACTATAAAGATACCAGGCTTTCCCGTGAAGCTCCCTCGTGCCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTT
2801 CGGAAGCGTGGCGTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCTGCTCAAGCTGGGCTGTGTGCAGAACCCCGTTCA
2901 GCCCAGCCGCTGCGCTTATCCGGTAACTATCGTCTTGTAGTCAACCCGGTAAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAAACAGGATTAGC
3001 AGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGC
3101 CAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGTATCCGGCAAAACAAACCCGCTGGTAGCGGTGGTTTTTTTGTGGAAGCAGCAGATTACGCGCAG
3201 AAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTATGGCTAGTTAA
3301 TTAACATTTAAATCAGCGGCCGAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTTGTGGAATCGTAACTAACATACGCTCTCCATCA
3401 AAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCGAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA