



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAAACAAACG
101 AAACAAAACAACTAGCAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCCGCTGGCTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGTAAGTCCGCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTGCCGCCAGAACACAGCTGAAGCTTCA CCGGTCACCATGGAATCAAGGTGCTGTTTGCCTCATCTGT
12 M E I K V L F A L I C
901 ATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCAACTTTGCCACCACAGATCTTGAGACTGACCTGT
12 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 TTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATCTGGCTTCA
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 AGGATAAGGAGCCACTGGACAGTTTATTGCTCAAGTGGACCTCTGTCTGATTGCACCACTGGCTGTCTGAAGGGCTTCCAATGTCCAGTGTCTGA
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 CCTCCTGAAGAAGTGGCTTCCCAGAGGTGTACCACCTTTGCCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGACAGATGAGCT
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 AGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTG
1601 CTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAACAACAACAATTGCATTATTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGTTTTTTA
1701 AAGCAAGTAAACCTCTACAATGTGGTATGGAATTCTAAATACAGCATAGCAAACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAG
1801 GGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGATTTTCCC
1901 AAGGTTTGAAGTACTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCCACATCCCTTTTATGTAATAATTCAGAAATAATTTAAATACATCATT
2001 GCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACC
2101 TTTAATAGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTTGCACAAAAGTGCACGCAGTTGCCGGCCGGTGCAGCAGG
125 D Q E E A V F H V C N G A P D R L
2201 GCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTATGGCCGGCCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACA
106 A F E 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
2301 GCTCGTCCAGGCCGCGCACCCACCCAGGCCAGGGTGTGTCGGCACCACTGGTCTGGACCGGCTGATGAACAGGGTACGTCGTCCCGGACCAC
73 E D L 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
2401 ACCGGCAAGTGTCTCCACGAAGTCCCGGGAACCCGAGCCGGTCCGAGAACTCGACCGTCCGGCGACGTCGCGCGGGTGAACCCGGAACG
40 G A F 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
2501 GCACTGGTCAACTTGGCCATGATGGCCCTCTATAGTGTGATGATTACTATGCGGATATACTATGCGGATGATTAATTGTCAATCCGGTTGCTTTGA
6 A S T L K A M
2601 ATTAGCGTGGTTTTCACAACTTAAAAAGGGTTAAAAAGATACCTTTGAACCGCTAAGAAGCCGAGAAATTAGCTCCGCTCAAACTCAAGGGGACA
2701 AATTCAAAAATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAATGTGAACAAAATCCAACGCCATTACATCCCCTCCCCCGCCGG
2801 ACTAGCCGTGCTCAAAAGCCGAGGTGACTATTGCGCCGATAGGACCACGGGTACAGGAAGCAGCAGCCGGTGAAGGACCAGGCCCTTCTCTTTGT
2901 GTGGGTGACTCACCCGCCGCTCCACCGGGTCCCGCTCCTCATTGAGTCTTGAACAGGGCCGGAGCGGCATTTTCCACGCACGCAACT
3001 GGTGCCGACGGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACGCGGCGCCCAAGCCAGATCGTGCCGGGTGCTGGGGCCACATGGCCTCGGCAC
3101 GCTAACCCAGCCTGGTTGCTTCGGGAAAACCCAGGCTCGCCCATCCAGGTGGCGTCCGACATGTGCTCCGAAGGGGGGGCCAGCCGCCAC
3201 TCCTGTCCCTCCATTCTCCCAACCATGACCTCTCCGGGCTCCGGGCGAGCAAGCCCGCACCTCCCTTTGTTAGCCCTATTGCTGAACGGCAATCG

3301 AAGGCAGCAGGGCAACAACAACAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGGCGGGCGGGACACCACGCTAGGCCTCAAGCCGGAC
3401 ACGAGGCGAGGCTACGGGGTTGCCGCTAGGCCTCGCACTCTGCCTCCCAGCGCCGCCGCAACTCGAAGCGGAATGCTCGCAGCTAATCCCCGCCGACGA
3501 CAGCGGGCCCGCCCGCTCGGAGCAGGACCTCCAGCTCGGCGGCCCGGAAGCCACACCCGCCCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAG
3601 AACGTCTGGCCCTGCGGGTCTGTGGTCTTATAGCCGCTCGGCGTCAGGCCGCCCGGCAATCAGCGCCGCCCGAGCCCGCTCCTTCC
3701 GGTGGGCGGGACCCGCCCTGCTGTGGGGAGGGCGGCCGCTGGAGGCCCTCGCGGCTCTGGCGGAAGTAGT**CGACGGACTGGGCTACGGCCGC**
3801 **CCCCGAGAGGCGCAGCCAGAGCCGCGCTAGGAAGGGCGGGCGCCGAGAACACGATCCCTCCCCACCCCTCGGACGTGACTCGGACCACATCCCGC**
3901 **GGTTCGCTAGGGCCCTCCTTCTGCTCCTTCCCCAGCCTGGCGGCTCTGGGGCGCGTGACTCAGCCAGAATGTTGGCAATGGGGAGGGCGGAACG**
4001 **GGGAAGTGAGGACGCGGATGGAAAAGTCGAAACGAAGGAAGCTGAGTTTCGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGCCAGCAAAAGGCC**
4101 **AGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACC**
4201 **CGACAGGACTATAAGATACCAGGCGTTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTCT**
4301 **CCCTTCGGGAAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCC**
4401 **G TTCAGCCCGACCCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGA**
4501 **TTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCT**
4601 **GAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACG**
4701 **CGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTA**
4801 **GT**
-