



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAAACAAACG
101 AAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA
201 AATGGCCCCGCTGGTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCATTGACGTC
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGGC
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGGAGGGTTCGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT
701 GTCGTGACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCCGGTACCATTGAAATCAAGGTGCTGTTTGCCTCATCTGT
12 M E I K V L F A L I C
901 ATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGT
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 TGTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAA
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 AGGATAAGGAGCCACTGGACAGTTTATTGCTCAAGTGGACCTCTGTCTGATTGCACCACTGGCTGTCTGAAGGCCCTGCCAATGCCAGTCTGTA
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 CCTCTGAAGAAGTGGCTTCCAGAGGTGTACCACCTTTGCCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGACAGATGAGCT
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 AAGCAAGTAAACCTCTACAATGTGGTATGGAATTCTAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAG
1801 GGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGATTTTCCC
1901 AAGGTTTGAACACTAGCTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCACATCCCTTTTATGTAATAATCAGAAATAATTAATACATCATT
2001 GCAATGAAAATAAATGTTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACC
2101 TTTAATAGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCGAATTCTCGACTCATTCTTTGCCCTCGGACGAGTGTGGGGCGTGGTTTTCCACTATCGG
342 E K A R P R T S P R R N G S D A
2201 CGAGTACTTCTACACAGCCATCGGTCCAGACGGCCGCGCTTCTGCGGGCGATTGTGTACGCCGACAGTCCCGGCTCCGGATCGGACGATTGCGTCGCA
325 L V E V C G D T W V A A S R R A I Q T R G V T G A G S R V I A D C
2301 TCGACCTGCGCCAAGTGCATCATCGAAATGCGGTCAACCAAGCTCTGATAGAGTTGGTCAAGACCAATGCGGAGCATATACGCCGGAGCCGCGGC
292 R G Q A W A A D D F N G D V L S Q Y L Q D L G I R L M Y A R L R P
2401 GATCTGCAAGCTCCGGATCCCTCCGCTCGAAGTAGCGGCTGTCTGCTCATAAGCAACCAAGCCCTCCAGAAGAAGATGTTGGCAGCTCGTATT
258 S G A L E P H R R E F Y R T Q Q E M C A L W P R W F F I N A V E Y Q
2501 GGGAAATCCCCGAACATCGCCTCGCTCCAGTCAATGACCGCTGTTATGCGGCCATTGTCCTGAGGACATTGTTGGAGCCGAAATCCGCGTGCACGAGGTG
225 S D G F M A E S W D I V A T I R G N D T L V N N S G F D A H V L H
2601 CCGGACTTCGGGGCAGTCTCGGCCAAAGCATCAGCTCATCGAGAGCTGCGCGACGGACGCACTGACGGTGTCTGCATCACAGTTTGCCAGTGATAC
192 R V E P C D E A W L M L E D L A Q A V S A S V T D D M V T Q W H Y
2701 ACATGGGGATCAGCAATCGGCATATGAAATCACGCCATGTAGTGTATTGACCGATTCTTGGCGTCCGAATGGGCGAACCCTCGTCTGGCTAAGAT
158 V H P D A I A C I F D R W T T Y Q G I G Q P G F P G S T Q S L D
2801 CGGCCGACGAGTTCGATCAGCTCCGCGACGGGTTGCAGAACAGCGGGGAGTTCCGGTTTCAGGCAGGCTTGAACGTCGACACCTGTGACCGGGC
125 A A A I A D M L E A V P Q L V A P L E T E P L D Q L T V G Q A R R
2901 GGAGATGCAATAGGTGAGGCTCTCGTGAATTTCCCAATGTCAAGCACTTCCGGAATCGGGAGCGGGCCGATGCAAAGTCCGATAAACATAACGATCT
92 S I C Y T L S E S F E G I D L V E P I P L A A S A F H R Y V Y R D
3001 TTGTAGAAACATCGGCGAGCTATTTACCCGACGACATATCCACGCCCTCTACATCGAAGCTGAAAGCAGGAGATTCTTCCGCTCCGAGAGCTGCA
58 K Y F G D A C S N V R L V Y G R G G V D F S F A R S E E G E S L Q M
3101 TCAGGTCCGAGACGCTGCAACTTTTCGATCAGAAACTTCGACAGACGTCGCGGTGAGTTTCAGGCTTTTCATGATGGCCTCTATAGTGAGTCTG
25 L D S V S D F K E I L F K A V S T A T A L E P K K M
3201 ATTATACTATGCCGATATACTATGCCGATGATTAATTGCAATCCGGTTGCTTTGAATTAGCGGTGGTTTTTCAACACCTAAAAAGGGTTTAAAGAT

3301 ACCTTTGAACCGCTAAGAAGCCCGAGAATTAGCTCCGCTCAAACTCAAGGGGACAAATTCCAAAATGACTCCAGCGCCAGGCTGGCCTGACTAGTCT
3401 CCACCCACCAAATGTGAACAACTCCAACGCCATTACATCCCCTCCCCCGCCGACTAGCCGTGCTCAAAAGCCGAGGTGACTATTGCGGCCGATAG
3501 GACCACGGGGTCACAGGAAGCAGCAGCCGGTGAGGGACCAGGCCCTCTTCTTTGTGTGGTGACTCACCCGCCCGCTCCACCGGGCTGCCGCTCTCC
3601 ATTTTGAAGCTCTTGCAACAGGGCCCGGAGCGGCATCTTTCCACGCACGCAACTGGTGCCGGACGGGATGGCCTACCCTAGTTAGGGAGGCAGGGCA
3701 ACGCGGCGCCGCAAGCCAGATCGTGCCGGTGTGGGGCCACATGGCCTCGGCACGCTAACCCAGCCTGGTTGCTTCGGGAAAAACCCAGGCCTCGC
3801 CCCATCCAGGTGGCGTCGGACATGTGCTCCGAAGCGGGCGGGCCCCAGCCGCACTCCTGTCCCTCATTCTCCCAACCATGACCTCTCCGGGCTCC
3901 GGGCGAGCAAGCCCCGACCCTCCCTTTGTAGCCCTATTGCTGAACGGCAATCGAAGGCAGCAGGGCAACAACAACAAAAAAAAAAGACCAGAGTG
4001 CGGCCGGAGTAGCAGCGGGCGGGCGGACACCACGCTAGGCCTCAAGCCGGACACGAGGCAGGCTACGGGGTTGCCGTAGGCCTCGCACTCTGCC
4101 TCCCGCGCCGCCGCAACTCGAAGCGGGAATGCTCGCAGCTAATCCCCGCCGACGACAGCGGGGCCCGCCGCTCGGAGCAGGACCTCCAGCTCGGCGG
4201 CCCGCGGAAGCCACACCCGCCCTCACCTGCGTTCTGACGGCAAGCTTCGGCGAAGAAGCTTGCCCTGCGGGTTCGCTTGTGGTCTCTTATAGCCGCT
4301 CGGCGTCAGGCCCGCCCGCCAATCAGCGCCCGCCCGAGCCGCTCTTCGGTGGCGGGACCCCGCCCTGCTGTGGGGAGGGCGGGCCG
4401 CTGGAGGCCCTCGCGCTCTGGCGGAAGTACTAGTCGACGGACTGGGCTACGGGCCGCCCGAGAGGCGCAGCCAGAGGCCGCTAGGAAGGGCGGGC
4501 GCCGAGAACAGATCCCTCCCCACCCCTCGGACGTGACTCGGACCACATCCCGGGTTCGCTAGGGCCCTCCCTTCTGCTCTTTCCCCAGCCTGGC
4601 GCGCTCTGGGGCGCCGTGACTCAGCCAGAATGTTGGCAATGGGGAGGGCGGAACGGGGAAGTGGAGGACGCGGATGGAAAAGTCGGAACGAAGGAAGC
4701 TGAGTTTCGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGCCAGCAAAAGGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCT
4801 CCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCC
4901 CTCGTGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGTTTCTCATAGCTCACGCTGTAGGT
5001 ATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGA
5101 GTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAG
5201 TGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCG
5301 GCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTAC
5401 GGGGTCTGACGCTCAGTGAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTAGT