



1 CTCGAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAAACAAACG  
101 AAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAAGGACCTGCAGGCGTTACATAACTTACGGTA  
201 AATGGCCCCGCTGGTGACCGCCCAACGACCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAACGCCAATAGGGACTTTCATTGACGTC  
301 AATGGGTGGAGTATTTACGGTAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCCCTATTGACGTCAATGACGGTAAATGGCC  
401 CGCCTGGCATTATGCCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATCGCTATTACCATGATGATGCGGTTTTGG  
501 CAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGATTTCCAAGTCTCCACCCATTGACGTCAATGGGAGTTTGTGTTTACTAGTCAGTGGCC  
601 AGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGGAGGGTTCGGCAATTGATCCGGTGCCTAGAGAAGTGGCGCGGGTAACTGGGAAAGTGAT  
701 GTCGTGTAAGTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTTGCCGTGAACGTTTCCCTATCAGTGATAGAGATCTCCC  
801 TATCAGTGATAGAGATCTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTACCAGGTCACCATGTTCTGGGGCCCTGCATGCTGCTGCTGCTG  
12 M V L G P C M L L L L  
901 CTGCTGCTGGCCTGAGGCTACAGTCTCCCTGGGCATCATCCAGTTGAGGAGGAGAACCAGGACTTCTGGAACCGCGAGGCAGCCGAGGCCCTGGGTG  
12 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 CCGCAAGAAGCTGCAGCCTGCACAGACAGCCGCAAGAACCTCATCATCTTCTGGGCGATGGGATGGGGGTGTCTACGGTACAGTGCAGGATCCT  
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 AAAAGGGCAGAAGAAGGACAACTGGGGCTGAGATACCCCTGGCTATGGACCGTTCATATGTGGCTCTGTCCAAGACATACAATGTAGACAAACAT  
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 R F N Q C  
1301 ACACGACACGCGCAACGAGGTATCTCCGTGATGAATCGGGCCAAGAAAGCAGGGAAGTCAAGTGGGAGTGGTAACCACCACACGAGTGCAGCACGCTC  
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 GCCAGCCGACCTACGCCACACGGTGAACCGCAACTGGTACTCGGACGCCGACGTGCTGCCTCGGCCCGCAGGAGGGTGCAGGACATCGCTACG  
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 GCCAAGTGGGACAGGCTGGACGGGAAGAATCTGGTGCAGGAATGGTGGCGAAGCGCCAGGGTGCCTGATGTGGAACCGCACTGAGTGCATGCA  
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 ATGGAGATGACAGAGGCTGCCCTGCGCCTGCTGAGCAGGAACCCCGGGCTTCTTCTCTCTGAGGGTGGTGCATCGACCACGGTATCACGAAA  
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 TGCCGACCCTCCACGCTTCTTCTCGGAGGCTACCCCTGCGAGGAGCTCCATCTCGGGCTGGCCCTGGCAAGGCGGGACAGGAGGCGCTAC  
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 A R D R K A Y  
2101 ACGGTCTCTATACGAAACGGTCCAGGCTATGTGCTCAAGGACGGCGCCGCGGATGTTACCGAGAGCGAGAGCGGGAGCCCGAGTATCGGCAGC  
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 AGTCAGCAGTGCCTGGACGAAGAGACCCACGAGGCGAGGAGTGGCGGTGTTGCGCGCGCCCGCAGGCGACCTGTTTACGGCGTGCAGGAGCA  
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 TTCAGTTTCAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCAAAATACAGCATAGCAAACTTTAACCT  
2701 CCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCCTCACCTT  
2801 CTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGTTTGAAGTACTCTTCATTTCTTTATGTTTAAATGCACTGACCTCCACATTCCCTTTTAA  
2901 GTAAAATATTAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCC  
3001 AGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCGAATTCGACTCATTCCCTTGGCCCT  
3101 GGACGAGTGTGGGGCGTGGTTTCCACTATCGGCGAGTACTTCTACACAGCCATCGTTCAGACGGCGGCTTCTGCGGGCGATTGTGTACGCCGA  
336 P R T S P R R N G S D A L V E V C G D T W V A A S R R A I Q T R G V  
3201 CAGTCCCGCTCCGGATCGGACGATTGCGTGCATCGACCCTGCGCCCAAGCTGCATCATCGAAATGCGCTCAACCAAGCTCTGATAGAGTTGGTCAAG  
303 T G A G S R V I A D C R G Q A W A A D D F N G D V L S Q Y L Q D L

3301 ACCAATGCGGAGCATATACGCCCGAGCCGCGGCGATCTGCAAGTCCGGATGCCTCCGCTCGAAGTAGCGCGTCTGCTGCTCCATACAAGCCAACCAC  
270 G I R L M Y A R L R P S G A L E P H R R E F Y R T Q Q E M C A L W  
3401 GGCTCCAGAAGAAGATGTTGGCGACCTCGTATTGGGAATCCCCGAACATCGCTCGCTCCAGTCAATGACCCTGTTATGCGGCCATTGTCCGTCAGGA  
236 P R W F F I N A V E Y Q S D G F M A E S W D I V A T I R G N D T L V  
3501 CATTGTTGGAGCCGAAATCCGCGTGCACGAGGTCCGGACTTCGGGGCAGTCTCGGCCAAAGCATCAGTCAATCGAGAGCTGCGCGACGGACGCACT  
203 N N S G F D A H V L H R V E P C D E A W L M L E D L A Q A V S A S  
3601 GACGGTGTCTCCATCACAGTTTCCAGTGATACATGGGGATCAGCAATCGCGCATATGAAATCACGCCATGTAGTGTATTGACCATTCTTGCGGT  
170 V T D D M V T Q W H Y V H P D A I A C I F D R W T T Y Q G I G Q P  
3701 CCGAATGGCCGAACCCGCTCGTCTGGTAAGATCGGCCGAGCGATCGATCCATGAGTCCGCGACGGTTGCAGAACAGCGGGCAGTTCCGGTTTCAG  
136 G F P G F G S T Q S L D A A A I A D M L E A V P Q L V A P L E T E P  
3801 GCAGGTCTTCAACGTGACACCTGTGCACGGCGGGAGATGCAATAGGTGAGGCTCTCGCTGAATCCCAATGTCAAGCACTTCCGGAATCGGGAGCGC  
103 L D Q L T V G Q A R R S I C Y T L S E S F E G I D L V E P I P L A  
3901 GGCCGATGCAAAGTGCCGATAAACATACGATCTTTGTAGAAACATCGCGCAGTATTTACCCGAGGACATATCCACGCCCTCTACATCGAAGCTG  
70 A S A F H R Y V Y R D K Y F G D A C S N V R L V Y G R G G V D F S  
4001 AAAGCACGAGATTCTCGCCCTCGAGAGCTGCATCAGGTGCGAGACGCTGTGCAACTTTTCGATCAGAACTTCGCGACAGACGTCGCGGTGAGTTCAG  
36 F A R S E E G E S L Q M L D S V S D F K E I L F K A V S T A T L E P  
4101 GCTTTTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAATCCGGTTGCTTTGAATTAGCGGTG  
3 K K M  
4201 GTTTTCACAACACCTAAAAAAGGGTTTAAAAGATACCTTTGAACCGCTAAGAAGCCCGAGAATTAGTCCGCTCAAACTCAAGGGGACAAATTCCAAAA  
4301 ATGACTTCCAGCGCCAGGCTGGCCTGACTAGTCTCCACCCACCAAATGTGAACAAACTCCAACGCCATTACATCCCCTCCCCCGCCGCGACTAGCCGTG  
4401 CTCAAAGCCCGAGGTGACTATTGCGGCCGATAGGACCACGGGGTACAGGAAGCAGCAGCCGGTGAGGGACCAGGCCCTCTTCTTTGTGTGGGTGACT  
4501 CACCCGCCGCTCCACCGGGTCCCGCTCCTCATTTGAGTCTTGAACAGGGCCCGGAGCGGCCATCTTCCACGCACGCAACTGGTGCCGGAC  
4601 GGGATGGCCTCACCTAGTTAGGGAGGCAGGGCAACGCGCGCCGCAAGCCAGATCGTGCCGGGTGCTGGGGCCACATGGCCTCGGCACGCTAACCCCA  
4701 GCCTGTTGCTTCGGGAAAAACCCAGGCTCGCCCATCCAGGTGGCGTCGGACATGTGCTCCGAAGGCGGGCGGGCCAGCCGCACTCCTGTCCCT  
4801 CCATTCTCCCAACCATGACCTCTCCGGGCTCGGGGAGCAAGCCCGCACCTCCCTTTGTTAGCCCTATTGCTGAACGGAATCGAAGGCAGCAG  
4901 GGCAACAACAACAAAAAAGACCAGAGTGCGGCCGGAGTAGCACGCGGGCGGGCGGACACCACGCTAGGCCTCAAGCCGACACGAGGGCAG  
5001 GCTACGGGGTTGCCGCTAGGCCTCGCACTCTGCCTCCCGCGCCGCCGCAACTCGAAGCGGAATGCTCGCAGTAATCCCCGCCGACGACAGCGGGCC  
5101 CGGCCGCTCGGAGCAGGACCTCCAGCTCGGGGCGCCGGAAGCCACACCCGCCCTCACCTGCGTTCTGACGGCAAGTTCGGGAAGAAGCTGTGGC  
5201 CCTGCGGTGCTTGTGGTCTCTTATAGCCGCTCGGGTCCAGGCCGCCCGCCAATCAGCGCCGCCCGCCGAGCCGCTCTTCCGGTGGGCGCG  
5301 GGACCCCGCCCTGCTGTGGGGAGGGGCGCCGCTGGAGGCCCTCGCGCTCTGGCGAACTAGT CGACGGACTGGGCTACGGGCCGCCCGAGAGG  
5401 CGCAGCCAGAGGCCGCTAGGAAGGGGCGGGCGCCGAGAACACGATCCCTCCCCACCCCTCGGACGTGACTCGGACCACATCCC GCGGTGCTAG  
5501 GGCCCTCCCTTCTGCTCTTTCCCCAGCTGGCGCTCTGGGGCGCGTACTCAGCCAGAATGTTGGCAATGGGGAGGGCGGAACGGGAAAGTGA  
5601 GGACGCGGATGAAAAAGTCGAAACGAAGGAAGCTGAGTTTCGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTA  
5701 AAAAGGCCGCTTGTGGCTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACT  
5801 ATAAAGATACCAGGCTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGA  
5901 AGCGTGGCGTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTGCTCCAAGTGGGCTGTGTGCAGAACCCCGTTAGCCG  
6001 ACCGCTGCGCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAC  
6101 GAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTT  
6201 ACCTTCGAAAAAGATTGGTAGCTCTTGTATCCGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAA  
6301 AAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGT