



1 CCTGCAGGCGTTACATAAATTACGGTAAATGGCCCGCTGGCTGACCGCCCAACGACCCCCGCCATTGACGTCAATAATGACGTATGTTCCCATAGTAA
101 CGCCAATAGGGACTTTCCATTGACGTCAATGGGTGGAGTATTTACGGTAAACTGCCCACTTGGCAGTACATCAAGTGTATCATATGCCAAGTACGCCCC
201 TATTGACGTCAATGACGGTAAATGGCCCGCTGGCATTATGCCAGTACATGACCTTATGGGACTTTCCTACTTGGCAGTACATCTACGTATTAGTCATC
301 GCTATTACCATGATGATGCGGTTTTGGCAGTACATCAATGGGCGTGGATAGCGGTTTACTCACGGGGATTTCGAAGTCTCCACCCATTGACGTCAATG
401 GGAGTTTGTTTGACTAGTCAGGGCCCCAACCCCCAAGCCCCATTTCACAACACGCTGGCGCTACAGGCGCGTGACTTCCCTTGTCTTGGGGCGGG
501 GGGCTGAGACTCCTATGTGCTCCGGATTGGTCAGGCACGGCCTTGGCCCCGCTCCTGCCACCGCAGATTGGCCGCTAGGCTCCCGAGCGCCCTGGC
601 TCCGAGGGCCGGCACCATAAAAGAAGCCCTAGCCACGTCCCTCGCAGTTCGGCGGTCCCAGGGTCTGTCTCAAGCTTCCGCCAGAACACAGg
701 taagtgccgtgtgtggttcccggggctggcctctttacgggttatggccttgcgtgccttgaattacttccatgcccttgctgcagtacgtgattc
801 ttgatcccagactcgggttgaagtgggtgggagagttcgaggccttgcgcttaaggagccccttcgctcgtgcttgagttgaggcctggcttggggc
901 ctggggccgcccgtgtaactctggtggcaccttcgcccctgctcgtgctttcgtctagcattttaaatttttgataaccagctgcgacg
1001 cttttttctggcgagatagtctttaaagtcggggcaagatctgcacactggtatttcggtttttggggccgggggcgagggggcccgctgcgtccc
1101 agcgcacatgctcggcgaggcggggcctgcgagcgcggccacgagaatcgagcggggtagtctcaaactggccggcctgctctggtgcctgcctcgc
1201 gcccgctgtatcggccgctggggcggcaaggctggccggctcggcaccagttcgtgagcggaaagatggccgcttcccggcctgctgcaggggagc
1301 tcaaatggaggacgcggcggcgggagagcggggcgggtgagtcacccacacaaaggaaaaggcctttccttctcatccgtcgttcatgtgactcca
1401 cggagtaccgggcccgtccaggcacctcgattagtctcgcagcttttgagtagctgctcttttaggttggggggaggggttttatgcatggagtttcc
1501 ccacactgagtggggtggagactgaagagttaggccagcttggcacttgatgtaattctccttgaatttgcctttttgagtttgatcttgcctcattc

Agel (1672)

1601 tcaagcctcagacagtggttcaaagtttttttcttccatttcagGTGTCGTGAAAACACCCTAAAAGCCACCGTCAACATGTCAAGCTCTTCTGGC
1701 TCCTTCTCAGCCTTGTGCTGTAAGTGTGCTCAGTCCACCATTGAGGAACAGGCCAAGACATTTTTGGACAAGTTAACCACGAAGCCGAAGACCTGTT
7 L L L S L V A V T A A Q S T I E E Q A K T F L D K F N H E A E D L F
1801 CTATCAAAGTTCATTGCTTCTTGAATTATAACACCAATATTACTGAAGAGAATGTCCAAAACATGAATAATGCTGGGGACAAATGGTCTGCCTTTTTA
40 Y Q S S L A S W N Y N T N I T E E N V Q N M N N A G D K W S A F L
1901 AAGGAACAGTCCACACTTGCCCAAATGTATCCACTACAAGAAATCAGAATCTCACAGTCAAGCTTCAAGTCTCAGCTGCAGGCTCTTCAAGAAATGGGTCTTCA
74 K E Q S T L A Q M Y P L Q E I Q N L T V K L Q L Q A L Q Q N G S S
2001 TGCTCTCAGAAGACAAGACAAACGGTTGAACACAATTCTAAATACAATGAGCACCATCTACAGTACTGGAAAAGTTTGAACCCAGATAATCCACAAGA
107 V L S E D K S K R L N T I L N T M S T I Y S T G K V C N P D N P Q E
2101 ATGCTTACTTGAACAGGTTTGAATGAAATAATGGCAAACAGTTTACTACAATGAGAGGCTCTGGGCTTGGGAAAGCTGGAGATCTGAGGTCCGC
140 C L L G L E P G L N E I M A N S L D Y N E R L W A W E S W R S E V G
2201 AACGAGCTGAGCCATTATATGAAGAGTATGTGGTCTTGAAGAAATGAGATGGCAAAGCAATCATTATGAGGACTATGGGATTATTGGAGAGGAGACT
174 K Q L R P L Y E E Y V L K N E M A R A N H Y E D Y G D A W R G D
2301 ATGAAGTAAATGGGTAGATGGCTATGACTACAGCCGGCCAGTTGATTGAAGATGTGGAACATACCTTTGAAGAGATTAACCATTATATGAACATCT
207 Y E V N G V D G Y D Y S R G Q L I E D V E H T F E E I K P L Y E H L
2401 TCATGCCTATGTGAGGGCAAAGTTGATGAATGCCTATCCTTCTATATCAGTCCAATTGGATGCCTCCCTGCTCATTTGCTTGGTGATATGTGGGGTAGA
240 H A Y V R A K L M N A Y P S Y I S P I G C L P A H L L G D M W G R
2501 TTTTGGACAAATCTGACTCTTTGACAGTTCCTTTGGACAGAAACCAACATAGATGTTACTGATGCAATGGTGGACCAGGCCTGGGATGCACAGAGAA
274 F W T N L Y S L T V P F G Q K P N I D V T D A M V D Q A W D A Q R
2601 TATTCAAGGAGCCGAGAAGTTCTTTGTATCTGTTGGTCTTCTAATATGACTCAAGGATTCTGGGAAAATCCATGCTAACGGACCCAGGAAATGTTCA
307 I F K E A E K F F V S V G L P N M T Q G F W E N S M L T D P G N V Q
2701 GAAAGCAGTCTGCCATCCCACAGCTTGGGACCTGGGGAAGGGCGACTTCAAGATCCTTATGTGCACAAAGGTGACAATGGACGACTTCTGACAGCTCAT
340 K A V C H P T A W D L G K G D F R I L M C T K V T M D D F L T A H
2801 CATGAGATGGGCATATCCAGTATGATATGGCATATGCTGCACAACCTTTTCTGCTAAGAAATGGAGCTAATGAAGGATCCATGAAGCTGTTGGGGAAA
374 H E M G H I Q Y D M A Y A A Q P F L L R N G A N E G F H E A V G E
2901 TCATGTCATTTCTGACGACACACCTAAGCATTTAAATCCATTGGTCTTCTGTACCCGATTTTCAAGAAGACAATGAAACAGAAATAAACTTCTGCT
407 I M S L S A A T P K H L K S I G L L S P D F Q E D N E T E I N F L L
3001 CAAACAAGCACTCACGATTGTTGGGACTCTGCCATTACTTACATGTTAGAGAAGTGGAGGTGGATGGTCTTAAAGGGGAAATCCCAAAGACCAGTGG
440 K Q A L T I V G T L P F T Y M L E K W R W M V F K G E I P K D Q W
3101 ATGAAAAAGTGGTGGGAGATGAAGCGAGAGATAGTTGGGGTGGTGAACCTGTGCCCATGATGAAACATACTGTACCCCGCATCTCTGTTCCATGTTT
474 M K K W W E M K R E I V G V V E P V P H D E T Y C D P A S L F H V
3201 CTAATGATTACTATTTCATTCGATATTACACAAGGACCCTTTACCAATTCAGTTTCAAGAAGCACTTTGTCAAGCAGCTAAACATGAAGGCCCTCTGCA
507 S N D Y S F I R Y Y T R T L Y Q F Q F Q E A L C Q A A K H E G P L H

S N D Y S F I R Y Y T R T L Y Q F Q F Q E A L C Q A A K H E G P L H
3301 CAAATGTGACATCTCAAACCTACAGAAGCTGGACAGAACTGTTCAATATGCTGAGGCTTGAAAATCAGAACCTGGACCCTAGCATTGAAAAATGTT
540▶ K C D I S N S T E A G Q K L F N M L R L G K S E P W T L A L E N V
3401 GTAGGAGCAAAGAACATGAATGTAAGGCCACTGCTCAACTACTTTGAGCCCTTATTTACCTGGCTGAAAGACCAGAACCAAGAAATCTTTTGTGGGATGGA
574▶ V G A K N M N V R P L L N Y F E P L F T W L K D Q N K N S F V G W
3501 GTACCGACTGGAGTCCATATGCAGACCAAAGCATCAAAGTGAGGATAAGCCTAAAATCAGCTCTTGAGATAAAGCATATGAATGGAACGACAATGAAAT
607▶ S T D W S P Y A D Q S I K V R I S L K S A L G D K A Y E W N D N E M
3601 GTACCTGTTCCGATCATCTGTTGCATATGCTATGAGGCAGTACTTTTTAAAAAGTAAAAAATCAGATGATTCTTTTTGGGGAGGAGGATGTGCGAGTGGCT
640▶ Y L F R S S V A Y A M R Q Y F L K V K N Q M I L F G E E D V R V A
3701 AATTTGAAACCAAGAATCTCCTTTAATTTCTTTGCTACTGCACCTAAAAATGTGTCTGATATCATTCTAGAACTGAAGTTGAAAAGGCCATCAGGATGT
674▶ N L K P R I S F N F F V T A P K N V S D I I P R T E V E K A I R M
3801 CCCGGAGCCGTATCAATGATGCTTTCCGTCTGAATGACAACAGCCTAGAGTTTCTGGGGATACAGCCAACACTTGGACCTCCTAACAGCCCCCTGTTTC
707▶ S R S R I N D A F R L N D N S L E F L G I Q P T L G P P N Q P P V S
3901 CATATGGCTGATTGTTTTGGAGTTGTGATGGGAGTGATAGTGGTTGGCATTGTATCTGATCTTCACTGGGATCAGAGATCGGAAGAAGAAAAATAAA
740▶ I W L I V F G V V M G V I V V G I V I L I F T G I R D R K K K N K

NheI (4099)

GCAAGAAGTGGAGAAAATCCTTATGCCTCCATCGATATTAGCAAAGGAGAAAATAATCCAGGATCCAAAACTGATGATGTTGACAGCTCCTTTTAGC
774▶ A R S G E N P Y A S I D I S K G E N N P G F Q N T D D V Q T S F •
4101 TAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATT
GCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAATTCATTTCATTTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTT
4201
4301 AAAGCAAGTAAACCTCTACAATGTGGTATGGAATGTTAATTAAGTACCATGACCAAAATCCCTAACGTGAGTTTTCGTTCCACTGAGCGTCAGAC
4401 CCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTTCTGCGGTAATCTGCTGCTTGAACAACAAAAAACCACCCTACCAGCGGTGTTTGT
4501 TGCCGATCAAGAGCTACCAACTCTTTTTCCGAAGTAACTGGCTTACAGAGCGCAGATACCAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCA
4601 CCACTTCAAGAACTCTGTAGCACCCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCGTGTCTTACCGGTTG
4701 GACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTGGGCTGAACGGGGGTTCTGTGCACACAGCCAGCTTGAGGCAACGACCTACACCGAACTGA
4801 GATACCTACAGCGTGAGCTATGAGAAAGCCACGCTTCCGAAGGAGAAAGGCGGACAGGTATCCGGTAAGCGGCAGGGTCGGAACAGGAGAGCGCAC
4901 GAGGGAGCTTCCAGGGGAAACGCCTGGTATCTTTATAGTCTGTCTGGGTTTCCGCACCTCTGACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGG
5001 CGGAGCTATGAAAAACGCCAGCAACGCGGCTTTTTACGGTTCCTGGCCTTTTGTGCTGCTTGTGCTCACATGTTCTTAATTAACCTGCAGGGCCTGA
5101 AATAACCTCTGAAAGAGGAACTTGGTTAGGTACCTTCTGAGGCTGAAAGAACCAGCTGTGGAATGTGTGTCAGTTAGGGTGTGGAAGTCCCCAGGCTCC
5201 CCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCAGGTGTGGAAGTCCCAGGCTCCCAGCAGGCAGAAGTATGCAAAGCATGC
5301 ATCTCAATTAGTCAGCAACCATAGTCCCACTAGTTCGCCAGAGCGCGGAGGGCTCCAGCGGCCCCCTCCCCACAGCAGGGGCGGGTCCCGCGC
5401 CCACCGAAGGAGCGGGCTCGGGCGGGCGGCTGATTGGCGGGGCGGCTGACGCCGACGGCTATAAGAGACCACAAGCGACCCGAGGGCCAG
5501 ACGTCTTCGCCGAAGCTTCCGTCAGAACGAGGTGAGGGGCGGTGGCTTCCGGGGCCCGAGCTGGAGGTCTGCTCCGAGCGGGCGGGCC
5601 CGTGTCTGCGGGGATTAGTGCAGCATTCCCGTTCGAGTTGCGGGCGGCGGGAGGAGAGTGCAGGCTAGCGGCAACCCGTAGCCTCGC
5701 CTCGTGTCCGGCTTGAAGCCTAGCGTGGTGTCCGCGCCGCCGCGTCTACTCCGGCCACTCTGGTCTTTTTTTTTTTTGTGTTGTTGCCCTGCT
5801 GCCTTCGATTGCCGTTACGAATAGGGGCTAACAAAGGAGGGTGCGGGCTTGTCTGCCCGAGCCCGGAGAGTTCATGGTTGGGGAGGAATGGAGGA
5901 CAGGAGTGGCGCTGGGGCCCGCCCTTCGGAGCACATGTCGACGCCACTGGATGGGGCAGGCCTGGGTTTTTCCGAAGCAACCAGGCTGGG
6001 TTAGCGTCCGAGGCCATGTGCCCCAGCACCCGGCACGATCTGGCTTGGCGGCGCGGTTGCCCTGCCTCCCTAACTAGGGTGAAGCCATCCCGTCCG
6101 GCACCAGTTGCGTGCCTGAAAGATGGCCGCTCCGGGCCCTGTTGCAAGGAGCTCAAATGGAGGACGCGGCAGCCGGTGGAGCGGGCGGGTGAAGTCA
6201 CCCACACAAAGGAAGAGGGCTGTCCCTCACCGGCTGCTGCTTCTGTGACCCGTTGCTCTATCGGCCAATAGTCACCTCGGGCTTTTGTAGCACGG
6301 CTAGTCGCGCGGGGGAGGGATGTAATGGCTTGGAGTTTGTTCACATTTGGTGGTGGAGACTAGTCAGGCCAGCCTGGCGCTGGAAGTCATTTTTG
6401 GAATTTGCTCCCTTGTGTTTTGAGCGGAGCTAATCTCGGGCTTCTAGCGGTTCAAAGGTATCTTTTAAACCCTTTTTAGGTGTTGTGAAAACCACCG

AgeI (6514)

CTAATTCAAAGCAACCGGTACAGTACCCCTGCCCGCCCGAGGTGAAAGCGGGTGTGAGGAGCGCGCGGCGGAGGTGATTTGAACATTCAGATA
▶ 1▶ M P P A P P G G E S G C E E R G A A G H I E H S R Y

6601 CCTATCATTACTCGATGCTGTTGATAACAGCAAGATGGCTTTGAACTCAGGGTACCACCAGCTATTGGACCTTACTATGAAAACCATGGATACCAACCG
26▶ L S L L D A V D N S K M A L N S G S P P A I G P Y Y E N H G Y Q P
6701 GAAAACCCCTATCCCGCACAGCCCACTGTGGTCCCCACTGTCTACGAGGTGCATCCGGCTCAGTACTACCCGTCCTCCCGTGCCTCCAGTACGCCCGGAGG
60▶ E N P Y P A Q P T V V P T V Y E V H P A Q Y Y P S P V P Q Y A P R
6801 TCCTGACGCAGGCTTCCAACCCCGTGTGCACGCAGCCAAAATCCCCACTCCGGGACAGTGTGCACCTCAAAGACTAAGAAAGCAGCTGTGCATCCTT
93▶ V L T Q A S N P V V C T Q P K S P S G T V C T S K T K K A L C I T L
6901 GACCCTGGGACCTTCTCGTGGGAGCTGCCTGGCCGCTGGCCTACTCTGGAAGTTCATGGGACGCAAGTGTCCAACTCTGGGATAGAGTGCGACTCC
126▶ T L G T F L V G A A L A A G L L W K F M G S K C S N S G I E C D S
7001 TCAGGTACCTGCATCAACCCCTAACTGGTGTGATGGCGTGCACACTGCCCCGGGGGAGGACGAGAATCGGTGTGTTGCGCTCTACGGACCAAACT
160▶ S G T C I N P S N W C D G V S H C P G G E D E N R C V R L Y G P N
7101 TCATCTTCAGGTGTACTCATCTCAGAGGAAGTCTGGCACCCTGTGTGCCAAGACGACTGGAACGAGAAGTACGGGCGGGCGCTGCAGGGACATGGG
193▶ F I L Q V Y S S Q R K S W H P V C Q D D W N E N Y G R A A C R D M G
7201 CTATAAGAATAATTTTACTTAGCAGGAATAGTGGATGACAGCGGATCCACCAGCTTTATGAAACTGAACACAAGTCCCGGCAATGTGCATATCTAT
226▶ Y K N N F Y S S Q G I V D D S G S T S F M K L N T S A G N V D I Y
7301 AAAAACTGTACCACAGTGTGCCTGTTCTTCAAAGCAGTGGTTTCTTACGCTGTATAGCCTGCGGGGTCAACTTGAACCAAGCCGCAGAGCAGGA
260▶ K K L Y H S D A C S S K A V V S L R C I A C G V N L N S S R Q S R
7401 TTGTGGGCGGAGAGCGCGCTCCCGGGGCGCTGGCCCTGGCAGTGCAGCTGCAGCTCCAGAAGTCCACGTGTGCGGAGGCTCCATCATCACCCCGA
293▶ I V G G E S A L P G A W P W Q V S L H V Q N V H V C G G S I I T P E
7501 GTGGATCGTGACAGCCGCACTGCGTGGAAAACTCTTAAACATCCATGGCATTGGACGGCATTGGCGGGATTTGAGACAATCTTTCATGTTCTAT
326▶ W I V T A A H C V E K T C P L N N P W H W T A F A G I L R Q S F M F Y
7601 GGAGCCGATACCAAGTAGAAAAAGTATTCTCATCAAATTATGACTCCAAGACCAAGAACAATGACATTGCGCTGATGAAGCTGCAGAGCCTCTGA
360▶ G A G Y Q V E K V I S H P N Y D S K T K N N D I A L M K L Q K P L
7701 CTTTCAACGACCTAGTGAACCCAGTGTGCTGCCAACCCAGGCATGATGCTGCAGCCAGAACAGCTCTGCTGGATTTCCGGGTGGGGGCCACCGAGGA
393▶ T F N D L V K P V C L P N P G M M L Q P E Q L C W I S G W G A T E E
7801 GAAAGGAAGACTCAGAAGTGTGAACGCTGCCAAGGTGCTTCTATTGAGACACAGAGATGCAACAGCAGATATGTCTATGACAACCTGATCACACCA
426▶ K G K T S E V L N A A K V L L I E T Q R C N S R Y V Y D N L I T P
7901 GCCATGATCTGTGCCGGCTTCTGCAGGGGAACGTCGATTCTTGCAGGGTGCAGTGGAGGGCCTTGGTCACTTGAAGAACAATATCTGGTGGCTGA
460▶ A M I C A G G F Q G N V D S C Q G G D S G G P L V T S K N N I W W L
8001 TAGGGATACAAGCTGGTCTGTGCCAAAAGCTTACAGACCAGGAGTGTACGGGAATGTGATGGTATTACGCGATGGATTTATCGACAATAG
493▶ I G D T S W G S G C A K A Y R P G V Y G N V M V F T D W I Y R Q M R

XbaI (8182)

8101 GGCAGACGGCTAATCCGCTAGGAGCAGGTTTCCCAATGACACAAAACGTGCAACTTGAACCTCCGCCTGGTCTTCCAGGTCTAGAGGGGTAACACTTT
526▶ A D G •
8201 GTACTGCGTTTGGCTCCACGCTCGATCCACTGGCGAGTGTTAGTAACAGCACTGTTGCTTCGTAGCGGAGCATGACGGCGTGGAACCTCTCCTTGTA
8301 ACAAGGACCCACGGGGCCAAAAGCCACGCCACAGGGCCCGTCACTGTGTGCAACCCAGCAGCGGACTTTACTGCGAAACCCACTTTAAAGTGACATT
8401 GAAACTGGTACCCACACACTGGTGACAGGCTAAGGATGCCCTTCAGGTACCCCGAGGTAACACGCGACACTCGGGATCTGAGAAGGGGACTGGGGCTTCT
8501 ATAAAAGCGCTCGGTTTAAAAGCTTCTATGCCTGAATAGGTGACCGGAGTGGCACCTTTCCTTTGCAATTACTGACCTATGAATACAACCTGACTGT
8601 TTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGGCCACCATGAAGAACTGAACTGACAGCAACTTCTGTTG
8701 AGAAGTTTCTCATTGAAAAATTTGATTCTGTTTCTGATCTCATGCAGTGTCTGAAGGTGAAGAAAGCAGAGCCTTTTCTTTTGTGTTGGAGGAAGG
12▶ E K F L I E K F D S V S D L M Q L S E G E E S R A F S F D V G G R G
8801 TTATGTTCTGAGGGTCAATCTTGTGCTGATGGTTTTTACAAGACAGATATGTTTACAGACACTTTGCCCTGCTGCTGCTGCCAATCCAGAAGTTCTG
45▶ Y V L R V N S C A D G F Y K D R Y V Y R H F A S A A L P I P E V L
8901 GACATTGGAGAATTTTCTGAATCTCTCACCTACTGCATCAGCAGAAGAGCACAAAGGAGTCACTCTCCAGGATCTCCCTGAAACTGAGTCCAGCTGTT
79▶ D I G E F S E S L T Y C I S R R A Q G V T L Q D L P E T E L P A V
9001 TGCAACCTGTTGCTGAAGCAATGGATGCCATTGCAGCAGTGTCTGAGCCAACTCTGGATTTGGTCCCTTTGGTCCCCAAGGCATTGGTCACTAC
112▶ L Q P V A E A M D A I A A A D L S Q T S G F G P F G P Q G I G Q Y T
9101 CACTTGGAGGATTTTCAATTTGTGCCATTGCTGATCCTCATCTACTGGCAGACTGTGATGGATGACACAGTTTCTGTTCTGTTGCTCAGGCACTG
145▶ T W R D F I C A I A D P H V Y H W Q T V M D D T V S A S V A Q A L
9201 GATGAACCTCATGCTGTGGCAGAAGATTGCTGAAGTGCAGACCTGGTCCATGCTGATTTTGAAGCAACAATGTTCTGACAGACAATGGCAGAATCA
179▶ D E L M L W A E D C P E V R H L V H A D F G S N N V L T D N G R I
9301 CTGCACTATTGACTGGTCTGAAGCCATGTTGGAGATTCTCAATATGAGGTGGCAACATTTTTTTTTGGAGACCTTGGCTGGCTGCATGGAACAACA
212▶ T A V I D W S E A M F G D S Q Y E V A N I F F W R P W L A C M E Q Q
9401 AACAAGATTTTTGAAGAAGACCCAGAAGTGGTGGTCCCCAGACTGAGAGCTTACATGCTCAGAATGGCTGGACCACTGTATCAATCTCTG
245▶ T R Y F E R R H P E L A G A G S P R L R A Y M L R I G L D Q G L Y Q S L
9501 GTTGATGGAACCTTTGATGATGCTGCTGGGCACAAGGAAGATGTGATGCCATTGTGAGGTCTGGTCTGGAACCTGTTGAAGAACTCAAATGGAAGAA
279▶ V D G N F D D A A W A Q G R C D A I V R S G A G T V G R T Q I A R
9601 GGTCTGCTGCTTTGGACTGATGGATGTTGAAGTCTGGCTGACTCTGGAACAGGAGACCCTCCACAAGACCCAGAGCCAAGGAATGAATATTAGC
312▶ R S A A V W T D G C V E V L A D S G N R R P S T R P R A K E •
9701 TAGATTATCCCTAATACCTGCCACCCCACTTAACTCAGTGGTGAAGAAGCGTCTCAGAAGTGTGTTTCAATTGGCCATTTAAGTTTAGTAGTAAAA
9801 GACTGGTAAATGATAACAATGCATCGTAAAACCTCAGAAGGAAAGGAGAATGTTTTGTGACCCTTTGGTTTTCTTTTTGCGTGTGGCAGTTTAAAG
9901 TTATTAGTTTTTAAAATCAGTACTTTTTAATGAAAACAACCTTGACCAAAAATTTGTACAGAATTTTGGAGCCATTAAAAAGTTAAATGAGAAACCTG

10001 TGTGTTCTTTGGTCAACACCGAGACATTTAGGTGAAAGACATCTAATTCTGGTTTTACGAATCTGGAACTTCTTGAAAATGTAATTCTTGAGTTAACA
10101 CTTCTGGGTGGAGAATAGGGTTGTTTTCCCCCACATAATTGGAAGGGGAAGGAATATCATTTAAAGCTATGGGAGGGTTGCTTTGATTACAACACTGGA
10201 GAGAAATGCAGCATGTTGCTGATTGCCTGTCACTAAAACAGGCCAAAACTGAGTCCTTGGGTTGCATAGAAAAGCTG

