



1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA  
101 GAGAAGGTGGCGCGGGGTAAACTGGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC  
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC  
301 GCCATCCACGCCGTTGAGTCCGCTTCTGCCGCTCCCGCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

**NcoI (560)**  
**AgeI (552)**

501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGGGCCTACCTGAGATCACCGGTACCCATGGCTCGCTACAGACTGCACTCCTGGTTGCCTCGT  
601 CCTCCTTGTGTGGCGTTCAAGCAACTGAGGCAGGCCCTACGGCGCAACATGGAAGACAGCGTCTGCTGCCGTGATTACGTCCGTACCGTCTGCC  
13▶ L L A V A L Q A T E A G P Y G A N M E D S V C C R D Y V R Y R L P  
701 CTGCGCGTGGTGAACACTTCTACTGGACCTCAGACTCCTGCCGAGGCTGGCGTGGTGTGCTAACCTTCAGGGATAAGGAGATCTGTGCCGATCCCA  
47▶ L R V V K H F Y W T S D S C P R P G V V L L T F R D K E I C A D P

**NheI (866)**

801 GAGTGCCTGGGTGAAGATGATTCTCAATAAGCTGAGCCAATGAAGAGCCTACTCTGATGACCGTGGCTAGCTGGCCAGACATGATAAGATACATTGATG  
80▶ R V P W V K M I L N K L S Q •  
901 AGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAA

1001 ACAAGTTAAACAACAACAAATTGCATTCATTTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATG  
1101 GAATTCTAAAATACAGCATAGCAAACTTAACTCCAATCAAGCCTCTACTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGT  
1201 TGCCAATGTGCATTAGCTGTTTGAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTACTCTTCATTTCTTTATG  
1301 TTTTAAATGCACTGACCTCCACATTCCCTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAG  
1401 AATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAGCGAG

1501 CTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAGCAGTCAGGA  
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  
1601 GCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGTGACCACCTGATGGATCTGCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGG  
110▶ A Y 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  
1701 TGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTACGACAGACAGTGACCTGCCAATGTAGGCCTCAATGTGGACAGCAGA  
77▶ D 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  
1801 GATGATCTCCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGTCTGTTGCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCC  
44▶ I 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  
1901 AGATCCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAAA  
10▶ L D Q Q S I N F T K M  
2001 CAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCTCAATAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATG

2101 GGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCCATTTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGA  
2201 GTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCC  
2301 CATAAGGTCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTCAATTGACGTCAATAGGGGGGCTACTTGGCATATGATACACTTGTACTGCCA  
2401 AGTGGGAGTTTACCCTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGGGGC  
2501 GGGGTCGTTGGCGGTGACGACAGGCGGGCCATTTACCCTAAGTTATGTAACGCTGAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCA  
2601 GGAACCGTAAAAGGCCGCTTGTGGCGTTTTTCATAGGCTCCGCCCCCTGACGAGCATCAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCC

2701 GACAGGACTATAAAGATACCAGGCTTTCCCTGGAAGTCCCTCGTGGCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTC  
2801 CCTTCGGGAAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCCG  
2901 TTCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGAT  
3001 TAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTG  
3101 AAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTGATCCGGCAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGC

3201 GCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAACGAAAACACGTTAAGGGATTTTGGTCATGGCTAG

---

3301 TTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCC

3401 ATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCCAGTGAAGTGCAAGTGCCAGGTGCCAGAACATTTCTCTATCGAA