



EcoRI (23)

NotI (2) **XbaI (19)** **SdaI (38)** **SpeI (45)** **EcoNI (67)** **PvuII (76)**

1 **GCGGCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGGCCACTAGTCCACTCCCAGAGCCTTGCCAGGCAGCTGCAATCACCAACCAGCATCC**

101 **TTTGGGTTTGACCCACTGAGCACATGACCCCAATTAGTCTCTGGCAGCATCCCCTGCTCTCTGTTACATCAGAGAGCACAGAGTAGCCGATATAAAT**

201 **GCTACTGGATGCTGGAGGTGCGAAGACAGACAAGTCCCACACAGCAGCTTGGTGACACCTAGCAGACACC** **NeoI (270)**
ATGGAAATCAAGGTGCTGTTTGCCCTCAT
▶ M E I K V L F A L I

BglIII (385)

301 **CTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGTGTGTGGCCTCCAACCTTGCCACCACAGATCTTGAGACTGAC**
10▶ C 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

401 **CTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCC**
44▶ L 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

501 **CAGATGCTCTGAGGGAGCTGGAGGCAATGCCAGAAGGGCTGGTGCACAAGAGGCTGCCTCATTTGCCTCTCCACATTAAGTGCACCCCTAAGATGAA**
77▶ P D 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

601 **GAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCCTGGC**
110▶ K 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

701 **TTCAAGGATAAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGCTTGCCTGCAATGTCCAGTGCT**
144▶ F 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

801 **CTGACCTCCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTTGCCAGCAAGATTGAGGTAGGGTGGACAAAATCAAGGCTTGGCTGGGGACAGATG**
177▶ S D 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 •

NheI (904)

901 **ATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAACCAACTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATG**
210▶

1001 **CTATTGCTTATTTGTAACCATTATAAGCTGCAATAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTCAAGGGGAGGTGTTGGAGGT**

1101 **TTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATC**

1201 **CTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGT**

SspI (1383)

1301 **GTATTTTCCAAGTTTTGAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAA**

1401 **ATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAA**

1501 **CAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTGACGCGAGTTGCCGGCCGG**
125▶ • D Q E E A V F H V C N G A P

1601 **GTGCGCAGGGCGAACTCCCGCCCCACGGTGTCTCGCGATCTCGGTATGGCCGGCCCGGAGCGTCCCGGAAGTTCGTGGACACGACCTCCGACCAC**
110▶ D R L 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 E S W

1701 **TCGGCGTACAGCTCGTCCAGGCGCGCACCCACCCAGGCCAGGGTGTGTCGGCACCTGCTGCTGGACCGCGCTGATGAACAGGGTCACGTCGT**
76▶ E A Y L 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

SgrAI (1811)

1801 **CCCGGACCACCCGGCGAAGTCTCCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGTCAGAACTCGACCGCTCCGGCGACGTCGCGCGGGTGAG**
43▶ R V V 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

1901 **CACCGAACGGCACTGTTCAACTTGGCCATGATGGCTCCTCCTGTCAGGAGAGAAAGAGAAGAAGGTTAGTACAATTGCTATAGTGAGTTGATTATAC**
10▶ V P V A S T L K A M

2001 **TATGCAGATATACTATGCCAATGATTAATTGTCAA** **ACTAGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAA**

2101 **AAGGCCGCTTGTGGCTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTAT**

2201 **AAAGATAACAGCGTTTTCCCTGGAAGCTCCTCTGTCGCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTCGGGAAG**

2301 **CGTGGCGTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCCCGAC**

2401 **CGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGA**

2501 **GGTATGTAGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTAC**

2601 **CTTCGAAAAAGAGTTGGTAGCTCTTATCCGGCAAAACCAACCCTGCTGCTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAA**

2701 **GGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACAT**

2801 **TTAAATCA**