



```

                    PvuI (7)
                    SgfI (6)
1  GGATCTGCGATCGTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
EcoNI (96)
101  GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

                    HindIII (245)
                    Psp1406I (203)
                    PvuII (239)
                    Bsu36I (291)
                    EcoNI (287)
201  GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTCACGCGCCCGGCCCTACCTGAGGCC
301  GCCATCCACGCGGTGAGTGCAGTTCGCCGCTCCCGCCTGTGGTGCCTCCTGAACTGCGTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

                    NgoMIV (441)
401  GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

                    SphI (568)
                    BspLU11I (559)
                    KasI (535)
                    AgeI (552)
                    BsrGI (562)
501  TCTGTTCTGCGCGCTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTAACATGTACAGCATGCGACTCGCATCCTGTGCACATTGACA
                    M Y S M Q L A S C V T L T
601  CTTGTGCTCCTGTCAACAGCGCACCCACTTCAAGCTCCACTTCAAGCTCTACAGCGGAAGCAGCAGCAGCAGCAGCAGCAGCAGCAGCAGCAGCAGCAGC
14▶ L V L L V N S A P T S S S T S S S T A E A Q Q Q Q Q Q Q Q Q Q Q Q
                    PvuII (708)
701  ACCTGGAGCAGCTGTTGATGGACCTACAGGAGCTCCTGAGCAGGATGGAGAATTACAGGAACCTGAAACTCCCCAGGATGCTCACCTTCAAATTTTACTT
47▶ H L E Q L L M D L Q E L L S R M E N Y R N L K L P R M L T F K F Y L
                    BglIII (824)
                    HindIII (890)
801  GCCAAGCAGGCCACAGAATTGAAAGATCTTCAGTGCCTAGAAGTGAACCTTGGACCTCTGCGGCATGTTCTGGATTTGACTCAAAGCAAAAGCTTTCAA
80▶ P K Q A T E L K D L Q C L E D E L G P L R H V L D L T Q S K S F Q
901  TTGGAAGATGCTGAGAATTCATCAGCAATATCAGAGTAACTGTTGAAAACTAAAGGGCTCTGACAACACATTTGAGTGCCAATTCGATGATGAGTCA
114▶ L E D A E N F I S N I R V T V V K L K G S D N T F E C Q F D D E S
                    NheI (1095)
1001  CAACTGTGGTGGACTTTCTGAGGAGATGGATAGCCTTCTGTCAAAGCATCATCTCAACAAGCCCTCAATAACTATGTACCTCCTGCTTACAACACCGTAG
147▶ A T V V D F L R R W I A F C Q S I I S T S P Q •
                    MscI (1101)
1101  CTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACACTAGAATGCAGTGAIAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCT

                    HpaI (1233)
1201  TTATTTGTAACCATTATAAGCTGCAATAAAACAGTTAAACAACAACAATTGCATTCAATTTATGTTTCAGGTTCAAGGGGAGGTGTGGAGGTTTTTAA

                    EcoRI (1329)
1301  GCAAGTAAACCTCTACAAATGTGGTATGGAATTTCTAAAATACAGCATAGCAAACTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGG
1401  ATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAGATATAGTGATTTTCCAA

                    SspI (1568)
                    SwaI (1582)
1501  GGTTTGAAGTGTCTCTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATAGTAAATATTCAGAAATAATTTAAATACATCATTGC

                    EcoO109I (1643)
1601  AATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAACCTT
1701  TAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGAACCTTGAAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCAT
                    N R T Y K L P I L E E I T T K V L K G N
                    BstXI (1872)
1801  TCATCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACTGATGGATCTGTCCACCTCATC
120▶ M E I L V F C D P A Y D S I L E R C M G C P S V V R I S R D V E D
1901  AGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTACGACAGACAGTACCCTG
87▶ S Y P H R V A V I T D F D K Q G N S V A S G I A I A E A C V T V R
                    StuI (2007)
2001  CCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCGCTTGGCTGCTGATGGCCGCCGACATGGTGTGTTGTCTCATAGAGCATGGTGA
53▶ G I Y A E I H V A S I I E G T K T R I A A G V H H K N D E Y L M T I
                    BspHI (2157)
                    BbsI (2153)
                    XmnI (2149)
2101  TCTTCTCAGTGGCAGCTCCACAGCTCCAGATCTGCTGAGAGATGTTGAAGTCTTTCATGATGGCCCTCTATAGTGAAGTCTATTATACTATGCCGA
20▶ K E T A V E V L E L D Q Q S I N F T K M
                    AseI (2215)
2201  TATACTATGCCGATGATTAATTGCAAACAGCGTGGATGGGCTCCAGC TATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCCACCG

                    SpeI (2370)
2301  TACACGCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAAGTCCCGTGGATTTAAGTAGTCAAAAACAACTCCCATGACGTC

                    SnaBI (2498)
2401  ATGGGGTGGAGACTTGAAGTCCCGTGGATGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAAAGCGATGACTAATA

```

2501 CGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACT

NdeI (2603)

2601 TGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGA

PacI (2789)

PstI (2782)

SdaI (2781)

BspLU11

2701 ACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAA TTAAGAA

2801 CATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGCGCTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAAT

2901 CGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCTGC

3001 CGCTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCCGGTGTAGGTGCTTCGCTC

ApaLI (3113)

3101 CAAGCTGGGCTGTGTGCACGAACCCCGTTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTCCAAACCGGTAAGACACGACTTATCG

3201 CCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAA

3301 GAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAAACACCGCTGGTAGCGGTGG

3401 TTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAC

EagI (3549)

PacI (3529) SwaI (3538)

NotI (3548)

3501 TCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTT

3601 GTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCAAGTGAAGTGCAGGTGCCAGAACA

3701 TTTCTCTATCGAA