



PvuI (7)
SgfI (6)
1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **HindIII (245)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTCACGCGCCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

AgeI (552) **BspLU11I (560)** **EcoO109I (599)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCAACATGTCTGAAGAAATGTTTATGCAAATCTCAAATCCA
1 M S E E I V Y A N L K I Q
601 GGACCCTGATAAAAAAGAAGAAACCCAGAAGTCTGACAAATGTGGGGAAAAGTATCCGCCGATGCTCCCATTCACAGCAAAAAACAGCTTTGATTCTG
13 D P D K K E E T Q K S D K C G G K V S A D A S H S Q Q K T V L I L
701 ATTCTTCTATGCCTTCTGCTGTTCAATGGAATGGGGTCTTAGGAGGCATCTTTTATAACAATTTGGCAACAGAAATGATAAAATCGAATCAATTGCAAA
47 I L L C L L L F I G M G V L G G I F Y T T L A T E M I K S N Q L Q
801 GGGCAAGGAAGAACTTCAAGAAAATGTTTCCCTACAGCTGAAGCACAATCTCAACAGCTCCAAGAAAATCAAGAACCTTTCTGCCATGCTGCAAAGCAC
80 R A K E E L Q E N V S L Q L K H N L N S S K K I K N L S A M L Q S T
901 AGCCACCCAGCTGTGCCGAGAGCTGTATAGCAAAGAACCAGAGCAAAATGTAACCATGTCCAAGGGTTGAGAATGGTACAAGGACAGCTGTTATCT
113 A T Q L C R E L Y S K E P E H K C K P C P K G S E W Y K D S C Y S
1001 CAACTCAATCAGTATGGAACATGGCAAGAGGTGTCATGGCCTGCTCTGCTCGGAATGCCAGCTCCTGAAGGTTAAGAACAAGGATGTGCTGGAATTTA
147 Q L N Q Y G T W Q E S V M A C S A R N A S L L K V K N K D V L E F

DraIII (1166)
1101 TAAAGTACAAGAAGCTACGCTATTTTTGGCTTGCATTGTTGCCAGAAAAGATCGCACACAATATCCACTAAGTGAGAAGATGTTCTCTCTGAAGAGTC
180 I K Y K K L R Y F W L A L L P R K D R T Q Y P L S E K M F L S E E S

ScaI (1229) **Tth111I (1246)** **NdeI (1293)**
1201 TGAAAGAAGCACAGATGACATAGATAAGAAGTACTGCCGATATAGACAGGTCAATGTTTATTATACATACTGCACCTGATGAGAACAATATCATATG
213 E R S T D D I D K K Y C G Y I D R V N V Y Y T Y C T D E N N I I C

NheI (1397)
1301 GAAGAGACAGCCAGCAAGGTGCAAGTGGAAAAGTGTGTTGAATGGCCTCCAGAGGATAGCAGGTAGCTACTCTGAGTTGCTTGTCCCTAGCTGCTCTGCT
247 E E T A S K V Q L E S V L N G L P E D S R •

MscI (1403)
1401 AGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTG

HpaI (1535)
1501 CTTTATTTGTAACCATTATAAGCTGCAATAAACAGTTAAACAACAACATTCATTCTTTATGTTTCAGGTTCAAGGGGAGGTGTGGGAGGTTTTTTA

EcoRI (1631)
1601 AAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTAAATAACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTCTGAG
1701 GGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGTATTTCC

SapI (1813) **SspI (1870)** **SwaI (1884)**
1801 AAGTTTTGAACAGCTCTTCATTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCTTTTATGTAATAATTCAGAAATAATTTAAATACATCATT

EcoO109I (1945)
1901 GCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGGAACC

2001 TTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGGC
141 • N R T Y K L P I L E E I T T K V L K G

SacI (2145) **BstXI (2174)**
2101 ATTCACTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCA
121 N 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
2201 TCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCAGCAGACCCAATGGCAATGGCTTCAAGCAGACAGTACCC
87 D 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 (2309)
2301 TGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCACTTGGTCTGATGGCCGCCGACATGGTGTCTGTTGCTCCTCATAGAGCATGGT
54 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

BbsI (2455)
2401 GATCTTCTCAGTGGCAGCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTATAGTGTGATTATACTATGCC
21 I K E T A V E V L E L D Q Q S I N F T K M

AseI (2517) **SacI (2574)**
2501 GATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCAC

