



PvuI (7)
SgfI (6)
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTT

AgeI (552) **BspHI [m] (560)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCATCATGATCTTCCACACAGGAACAACGAAGCCTACCTGGT

MscI (626) **BbsI (667)**
601 GCTGCTTTGCTGTATAGGAACCTGGCTGGCCACCTGCAGTTGTCCTTCGGTGCCCAATATCGAAGGAAGACTTAAGAACTACAATTGACCTCTTGAAA
13▶ L L C C I G T W L A T C S L S F G A P I S K E D L R T T I D L L K
701 CAAGAGTCTCAGGATCTTTATAACAATATAGCATAAAGCAGGCATCTGGGATGTCAGCAGACGAATCAATACAGTCCCGTGTTCAGCCTGGACCGGG
47▶ Q E S Q D L Y N N Y S I K Q A S G M S A D E S I Q L P C F S L D R
801 AAGCATTAAACCAATCTCGGTATCATAGCACATCTGGAGAAAGTCAAAGTGTGAGCGAGAACACAGTAGATACTTCTGGGTGATAAGATGGCTAAC
80▶ E A L T N I S V I A H L E K V K V L S E N T V D T S W V I R W L T
901 AAACATCAGCTGTTCAACCCACTGAATTTAAACATTTCTGTCCCTGGAAACTGATGAATCCTATGATTGTAAGTGTTCGCTTACGGTTTTAAAG
113▶ N I S C F N P L N L N I S V P G N T D E S Y D C K V F V L T V L K

NheI (1094)
1001 CAGTTCTCAAAGTGCATGGCAGAAGTGCAGGCTAAGGACAATACTACATGCTGAGTGTGGGGGGGGTGCAGTGTCTCAGCAGTGCCTGTCTAGC
147▶ Q F S N C M A E L Q A K D N T T C •

MscI (1100)
1101 TGGCCAGACATGATAAGATACATTGATGAGTTTGACAAACCACAAGTGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTT

HpaI (1232)
1201 TATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAATTGCATTATTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAG

EcoRI (1328)
1301 CAAGTAAACCTCTACAAATGTGGTATGGAATCTAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTGAACTCTTTCTGAGGGA
1401 TGAATAAGGCATAGGCATCAGGGGCTGTGCCAATGTGCATTAGCTGTTGACGCTCACCTCTTTCATGGAGTTTAAAGATATAGTATTTTCCCAAG

SapI (1510) **SspI (1567)** **SwaI (1581)**
1501 GTTTGAAGTAGCTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCA

EcoO109I (1642)
1601 ATGAAAATAAATGTTTTTATTAGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTT
1701 AATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTCTCGGTGACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATT
141▶ • N R T Y K L P I L E E I T T K V L K G N

SacI (1842) **BstXI (1871)**
1801 CATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCA
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 R I S R D V E D
1901 GAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGCCCAATGGCAATGGCTTACGACAGACAGTACCCTGC
86▶ 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 G

StuI (2006)
2001 CAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGTCTGTGCTCATAGAGCATGGTGTAT
53▶ 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

BbsI (2152) **XmnI (2148)**
2101 CTTCTCAGTGGCGACCTCCACCAGTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGCTGGCCCTCTATAGTGTGATTATACTATGCCGAT
20▶ K E T A V E V L E L D Q Q S I N F T K M

AseI (2214) **SacI (2271)**
2201 ATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCCT

SpeI (2369)
2301 ACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAA

SnaBI (2497)
2401 TGGGGTGGAGACTTGAAAATCCCGTGAAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATGTAATAGCGATGACTAATAC

