



PvuI (7)
SgfI (6) **MfeI (82)**
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAAGTGGCGCGGGGTAAACTGGAAAAGTGTGCTGTACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTT

NcoI (575)
AgeI (552) **BspHI (560)** **BstXI (569)**
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCATCATGAAACCCACTGCCATGGCACTGATGTGCCTGCTGCT
SphI (635) 1▶ M K P T A M A L M C L L L

601 GGCTGCCGTGTGGATACAGGATGTTGACAGCAAGAGCATGCTTACGGTCTCCAATAGCTGCTGCTTGAACACCTTGAAGAAAGAGCTTCCCCTGAAGTTT
13▶ A A V W I Q D V D S K S M L T V S N S C C L N T L K K E L P L K F
PvuII (743)
701 ATCCAGTGTACAGAAAGATGGGCTCCTCCTGTCTGATCCCCCAGCTGTGGTATTACAGGCTGAACAAAGGTAGAGAAAGCTGCGCCTCAACTAACAAAA
47▶ I Q C Y R K M G S S C P D P P A V V F R L N K G R E S C A S T N K
MscI (849)
NheI (843)
801 CGTGGTTCAAAATCACCTGAAGAAGGTGAACCCCTGCTAACCCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGA
80▶ T W V Q N H L K K V N P C •

HpaI (981) **MfeI (992)**
901 ATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAAACAATTGCA

EcoRI (1077)
1001 TTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCAAAATACAGCATAGCA
1101 AAACTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTG

SapI (1259)
1201 CAGCCTCACCTTCTTTTATGAGTAAAGATATAGTGTATTTCCCAAGTGTGAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCAC

SspI (1316) **SwaI (1330)** **EcoO109I (1391)**
1301 ATTCCCTTTTATGATAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTC
1401 ATAATATCCCCAGTTTATGAGTGGACTTAGGGAACAAAGAACCTTAAATAGAAATTTGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGT
141▶ • N R T Y

SacI (1591)
1501 ACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCT
136▶ 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 A Y D S I L E R
BstXI (1620)
1601 GCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTG
103▶ 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 D F D K Q G N
StuI (1755)
1701 CTCACAGCAGACCAATGGCAATGGCTTCAGCAGACAGTACCCTGCCAATGTAGGCCTCAATGTGACAGCAGAGATGATCTCCCAGTCTTGGTCC
69▶ 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 I I E G T K T R
XmnI (1897)
1801 TGATGGCCGCCCCGACATGGTGTCTGTTGTCCTCATAGAGCATGGTGTCTTCTCAGTGGCGACTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAA
36▶ 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 L D Q Q S I N F
BbsI (1901) **AseI (1963)**
1901 GGTCTTCACTGGTGGCCCTCTATAGTGAAGTATTACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCT
3▶ T K M

SacI (2020)
2001 IATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTT

SpeI (2118)
2101 GGAAAGTCCCGTTGATTTACTAGTCAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCAT

SnaBI (2246)
2201 TGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCCATAGGTCATGTACTGGGCATA

