



PvuI (7)
SgfI (6) **MfeI (82)**
1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCCGCAATTGAACGGGTGCCTA
101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGTGCTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **HindIII (245)**
PvuII (239) **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCGCCCTACCTGAGGGC
301 GCCATCCACGCCGGTTGAGTCGGCTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCAGCTTTGCCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT

SgrAI (551) **BstXI (579)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACCGCGAAGGAGGGCCACCATGAACGCCAAGGTCGTGGTCTGCTGG
1 M N A K V V V V L
601 TCCTCGTGTGACCGGCTCTGCCTCAGCGACGGGAAGCCCGTCAAGCTGAGCTACAGATGCCCATGCCGATCTTCGAAAGCCATGTTGCCAGAGCCAA
10 V L V L T A L C L S D G K P V S L S Y R C P C R F F E S H V A R A N
701 CGTCAAGCATCTCAAATTTCTCAACTCCAAAGTGTGCCCTTCAAGTGTAGCCCGCTGAAGAACAACAACAGACAAGTGTGATTGACCCGAAGCTA
43 V K H L K I L N T P N C A L Q I V A R L K N N N R Q V C I D P K L
MscI (848)

HindIII (825) **NheI (842)**
801 AAGTGGATTACAGGAGTACCTGGAGAAAGCTTTAAACAAGTAACTAGCTAGCCGAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAGTAGAA
77 K W I Q E Y L E K A L N K •

HpaI (980) **MfeI (991)**
901 TGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAAATTGCAT

EcoRI (1076)
1001 TCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAA
1101 AACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGC

SapI (1258)
1201 AGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTTCCCAAGGTTGAAGTAGCTCTTCATTCTTTATGTTTTAAATGCACTGACCTCCACA

SspI (1315) **SwaI (1329)** **EcoO109I (1390)**
1301 TTCCCTTTTATGATAAAATATTAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTCA
1401 TAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTA
141 • N R T Y

SacI (1590)
1501 CTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTG
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 (1619)
1601 CACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGC
102 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 S
StuI (1754)
1701 TCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGTCT
69 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 (1896)
1801 GATGGCCGCCCCGACATGGTGTGTTGTCCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAG
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
BspHI (1904)
BbsI (1900) **AseI (1962)**
1901 GTCTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTT
2 T K M

SacI (2019)
2001 ATCTGACGGTTCATAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTG

SpeI (2117)
2101 GAAAGTCCGTTGATTTACTAGTCAAACAACCTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCCAAT

SnaBI (2245)
2201 GATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCCAAGTAGGAAAGTCCCATAGGTCATGTAAGTGGGCATAA

