



PvuI (7)
SgfI (6)
MfeI (82)

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCCGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTGTGCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)

Psp1406I (203)
PvuII (239)
Bsu36I (291)

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCGCCCTACCTGAGGGCC

301 GCCATCCACGCCGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

AgeI (552)
BspHI (560)
BglIII (585)
BstBI (598)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACCGGTATCATGAACTCGCTGGAACAAGCGGAAGATCTCAAGGCTTT

▶
1▶ M N S L E Q A E D L K A F

BbsI (605)
XmnI (616)
BsrGI (682)

601 CGAAAGAAGACTTACTGAATATATTCATTGCTTGAACCTGCCACTGGACGTTGGAGAATGCTTCTTATAGTGGTGTCTGTCTGTACAGCTACTGGTGCC

701 TGGAACTGGCTAATAGATCCTGAGACACAAAAGGTGTCCTTCTTACGTCATTATGGAACCATCCGTTTTTACTATTAGCTGATAACTCTAATAGGCT

801 TGTTCTTTGCTGGAATACACAAAAGAGTAGTCGACCATCAATTATAGCTGCTCGATGTCGAACGGTATTAGCAGAATACAATATGCTGTGATGATAC

80▶ L F F A G I H K R V V A P S I I A A R C R T V L A E Y N M S C D D T

MscI (952)

NheI (946)

901 AGGAAAGCTAATTTTGAACCTAGACCTCATGTTCAATGACAGTCCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAC

113▶ G K L I L K P R P H V Q •

HpaI (1084)
MfeI (1095)

1001 AGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACCAAGTTAAACAACAACAATT

EcoRI (1180)

1101 GCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATA

1201 GCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGT

SapI (1362)

1301 TTGCAGCCTCACCTTCTTTTCATGGAGTTTAAAGATAGTGTATTTTCCCAAGGTTTGAAGTACTCTTCATTCTTTATGTTTTAAATGCACTGACCTCC

SspI (1419)
Swal (1433)
EcoO109I (1494)

1401 CACATTCCTTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCC

1501 TTCATAATATCCCCAGTTTAGTGTGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGG

141▶ • N R T

SacI (1694)

1601 TGTAATTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTC

137▶ Y 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

BstXI (1723)

1701 TCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCG

104▶ R 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

StuI (1858)

1801 TTGCTCACAGCAGACCAATGGCAATGGCTTCAACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGG

70▶ N 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

1901 TCCTGATGGCCGCCCCGACATGGTGCTTGTGCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTT

37▶ R 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

BbsI (2004)
XmnI (2000)
AseI (2066)

2001 GAAGTCTTCATGGTGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCA

4▶ F T K M ▶

SacI (2123)

2101 GCTTATCTGACGGTTCACATAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACAT

SpeI (2221)

2201 TTTGAAAGTCCCCTGATTTTACTAGTCAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCC

SnaBI (2349)

2301 CATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCCAAGTAGGAAAGTCCCATAAAGTTCATGTAAGTGGG

