



PvuI (7)
SgfI (6)
MfeI (82) **EcoNI (96)**

1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCGCCCGCCGCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCCTGTTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441) **NaeI (441)**

401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560)

KasI (535)
BstEII (555) **AgeI (552)**

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCCTAGCTGAGATCACCGGTACCATTGGAACATAGCACGTTTCTCTCTGGCCTGGTACTGGC
 601 TACCTTCTCTCGCAAGTGAAGTCCCTTCAAGATACCTATAGAGGAAGTGGAGACAGAGTGTGTGTAATTGCAATACCAGCATCACATGGGTAGAGGGA
 701 ACGGTGGGAACTGCTCTCAGACATTACAAGACTGGACCTGGGAAAACGCATCCTGGACCCACGAGGAATATATAGTGTAAATGGGACAGATATATACA
 801 AGGACAAAGAATCTACCGTCAAGTTCATTATCGAATGTCCAGAGCTGTGTGGAGCTGGATCCAGCCACCGTGGCTGGCATCATTGTCAGTGTGCAT

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

BamHI (858)

901 TGCCACTCTGCTCCTTGTGTTGGAGTCTTCTGCTTGGACATGAGACTGGAAGGCTGTCTGGGGCTGCCGACACACAAGCTCTGTTGAGGAATGAC
 113▶ A T L L L A L G V F C F A G H E T G R L S G A A D T Q A L L R N D

Tth111I (997)

1001 CAGGTCTATCAGCCCTCCGAGATCGAGATGATGCTCAGTACAGCCACCTTGGAGGAACTGGGCTCGGAACAAGTGAACCTGAGACTGGTGGCTTCTAG
 147▶ Q V Y Q P L R D R D D A Q Y S H L G G N W A R N K •

NheI (1099)

MscI (1105)

1101 CTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTAT

HpaI (1237) **MfeI (1248)**

1201 TGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAATTGCATTCTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTT

EcoRI (1333)

1301 TAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTG
 1401 AGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTC

SspI (1572) **SwaI (1586)**

1501 CCAAGGTTTGAAGTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAAATATTAGAAAATAATTTAAATACATCA

EcoO109I (1647)

1601 TTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAA
 1701 CCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTCTCTGGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTG

141▶ • N R T Y K L P I L E I T T K V L K

SacI (1847) **BstXI (1876)**

1801 CCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCT
 121▶ G 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
 1901 CATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCCTTGTCTACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTGC
 88▶ 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

StuI (2011)

2001 CCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGTCTTGTTCCTCATAGAGCATG
 55▶ R 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

BspHI (2161)

BbsI (2157)

XmnI (2153)

2101 GTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGCTTCTCATGATGGCCCTCCTATAGTGAGTCGTATTATACTATG
 21▶ T I K E T A V E V L E L D Q Q S I N F T K M

AseI (2219)
SacI (2276)

2201 CCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACATAACAGAGCTCTGCTTATATAGACCTCC

SpeI (2374)

2300 CACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGAGTTGTTACGACATTTTGGAAAGTCCCCTGATTACTAGTCAAAAACAACTCCCATTG
 2399 ACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCAGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGA

