



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

1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAAGTGGCGGGGTAAGTGGAAAGTGTGCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245) **Bsu36I (291)**
EcoNI (287)

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCCTTCTGCCGCTCCCGCTGTGGTGCCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC
 401 GGGCCTTTGTCCGGGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

NcoI (560)
BstEII (555)

501 TCTGTTCTGGCGGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTACCATTGGAGTCGGCCGACTTCTACGAGGTGGAGCCGGGCC
 601 CCCGATGAGCAGTCACTCCAGAGCCCCCGCACGCGCCAGCAACGCCCTTTGGCTTTCCCGGGGCGGGGCCCCGCGCCCGCCACAGCCACCT
 13▶ P M S S H L Q S P P H A P S N A A F G F P R G A G P A P P A P P
 701 GCCGCCCCGAGCGCTGGGCGGCATCTGCGAGCACGAGAGCTTATAGACATCAGCGCTACATCGACCCGGCCGCTTCAACGACGAGTTCCTGGCCG
 47▶ A A P E P L G G I C E H E T S I D I S A Y I D P A A F N D E F L A
 801 ACCTTTCCAGCAGCCGACAGCAGGAGAAGGCAAGGCGGCGGGCCCCGCGGTGGCGCGGTGACTTTGACTACCCGGAGCCCGCGGGGCC
 80▶ D L F Q H S R Q Q E K A K A A A G P A G G G G D F D Y P G A P A G P
 901 CGCGGCGCGGTGATGTCGCGGGGGCGCACGGCCCCCTCCCGGCTACGGCTGTGCGGCGGGCGGCTACCTGGACGGCAGGCTGGAGCCCTGTACGAG
 113▶ G G A V M S A G A H G P P P G Y G C A A A G Y L D G R L E P L Y E

KasI (1008)

1001 CGCGTCGGGGCGCCCGCTACGCGCGTGGTGTATCAACAAGAGCCCCGCGAGGAGGACGAGGCGAAGCAGCTGGCGCTGGCCGGCCTTCCCTACC
 147▶ R V G A P A L R P L V I K Q E P R E E D E A K Q L A L A G L F P Y

MluI (1133)

1101 AGCCACCGCCGCCACCGCCACCGCCGACCGCTCCTCCGCGCACCTGGCCGCCCCCACTTGCAGTTCAGATCGCGCACTGCGCCAGACCAC
 180▶ Q P P P P P P P P H P H A S P A H L A A P H L Q F Q I A H C G Q T T

MscI (1215)
SfiI (1282)

1201 CATGCACCTACAGCTGGCACCACACCGCCCGCCAGCCCGTGGCCAGCCGACGCTGCGCCCGCCTTGGGTGCTGCGGGCTGCCTGGCCCCGGG
 213▶ M H L Q P G H P T P P P T P V P S P H A A P A L G A A G L P G P G

BssHII (1301)
XcmI (1385)

1301 AGCGCGCTCAAGGCTTGGCCGGTGGCACCACCGCTCCGCACGGGAGGCGGGCGGTGGCAGCGGTGCGGGCAAAGCCAAGAAGTCGGTGG
 247▶ S A L K G L A G A H P D L R T G G G G G G S G A G A G K A K K S V
 1401 ACAAGAACAGCAACGAGTACCGGTACGGCGGGAACGCAACACATCGCGGTGCGCAAGAGCGAGATAAAGCCAACACGCAACCGTGGAGACGCAACA
 280▶ D K N S N E Y R V R R E R N N I A V R K S R D K A K Q R N V E T Q Q
 1501 GAAGTGTGGAGTTGACAGTGAACATGACCGCTGCGCAAGCGGTGGAACAGCTGAGCCGTAACGGACACGCTGCGGGGCATCTCCGCCAGCTG
 313▶ K V L E L T S D N D R L R K R V E Q L S R E L D T L R G I F R Q L

NcoI (1622)
NheI (1645)
MscI (1651)

1601 CCTGAGAGCTCCTGGTCAAGGCCATGGGCACTGCGCGTGAAGCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAATA
 347▶ P E S S L V K A M G N C A •

HpaI (1783) **MfeI (1794)**

1701 GAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACATTTG

EcoRI (1879)

1801 CATTCAATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAAATCTAAAATACAGCATAG
 1901 CAAAACCTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTT
 2001 TGCAGCCTCACCTCTTTTATGAGTAAAGATATAGTATTTTCCCAAGGTTTGAAGTCTTCTTCAATTTCTTTATGTTTTAAATGCACTGACCTCCC

SspI (2118) **Swal (2132)**

2101 ACATTCCTTTTTTAGTAAATATTTCAGAAATAATTTAAATACATCATTGCAATGAAAAATAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCT
 2201 TCATAATATCCCCAGTTTGTAGTGTGGACTTAGGGAACAAAGGAACCTTAAATGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGT
 2301 GTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAAGGATAGTCAAGATGAGCTCT
 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 (2422)

2401 CTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCGTACAGCCACAATGGTGTCAAAGTCTTCTGCCCGT
 103▶ 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 N

StuI (2557)

2501 TGCTCACAGCAGACCCAATGGCAATGGCTTTCAGCACAGACAGTACCCTGCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGT
 70▶ 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

XmnI (2699)

2601 CCTGATGGCCCGCCGACATGGTGTGTTGTCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCTGCTGAGAGATGTTG
 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

