



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

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCTCCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMI (441)
NaeI (441)

401 GGGCCTTTGTCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560)
BstEII (555)
KasI (535) **AgeI (552)** **Neol (582)**

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGGCTACCTGAGATCACCGGTACCATTGGGCTCACAGACCCTGCCCATGGTCACATGCAGAC
 601 CCTCATCTTCTTAGACCTGGAAGCCACTGGCCTGCCTTCTGCTCGGCCGAAGTCACAGAGCTGTGCCTGCTGGCTGTCCACAGAGCTGCTCTGGAGAAC
 13▶ L I F L D L E A T G L P S S R P E V T E L C L L A V H R R A L E N
 701 ACTTCCATTTCTCAGGGACATCCACCTCAGTGCACAGACCGCCCGTGTGGTGGACAAGCTCTCTGTGCATTGCTCCAGGAAAGCTGTAGCCCTG
 47▶ T S I S Q G H P P P V P R P P R V V D K L S L C I A P G K A C S P

MscI (872)
BalI (872)

801 GGGCCAGTGAGATCACAGTCTGAGCAAAGCTGAGCTGGAAGTACAGGGCGTCAACGCTTCGATGACAACTGGCCATCCTGCTCCGAGCCTTCTGCA
 80▶ G A S E I T G L S K A E L E V Q G R Q R F D D N L A I L L R A F L Q

BstEII (937)

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

Asp718I (1006)
Acc65I (1006)
StuI (1040)
Eco147I (1040) **NheI (1054)**

1001 CTAGATGGTACCTTCTGTGGACAGCATCGCTGCCCTAAAGGCTTGGAAACAAGCTAGCAGCCCTCAGGGAATGGTTCGAGGAAAAGCTACAGCCTGG
 147▶ L D G T F C V D S I A A L K A L E Q A S S P S G N G S R K S Y S L
 1101 GCAGCATCTACACCCGCTGTACTGGCAAGCACCGACAGACTACATACTGCTGAAGGTGATGTTCTAACCTGCTCAGCATCTGTGAGTGAAGCCACA
 180▶ G S I Y T R L Y W Q A P T D S H T A E G D V L T L L S I C Q W K P Q
 1201 GGCCCTACTGCAGTGGGTGGACGAACATGCCCGCCTTTAGCACCGTCAAGCCATGTACGGCACTCCGGCTACCCTGGAACAACCACTAAGGCCA
 213▶ A L L Q W V D E H A R P F S T V K P M Y G T P A T T G A T T N L R P

MscI (1325)
BalI (1325)

1301 CATGCTGCCACAGCTACTACACCCTGGCCACAGCCAATGGAAGTCCCAGCAATGGCAGGAGCAGGGCAGCCTAAGAGTCTCCTCCAGAGAAGGTCCCAG
 247▶ H A A T A T T P L A T A N G S P S N G R S R R P K S P P P E K V P
 1401 AAGCCCATCACAGGAGGGGCTGCTGGCCCACTGAGCCTGCTGACCCTCCTGACCTTGGCAATAGCCACTCTGTATGGACTTCTCCTGGCCTCACCTGG
 280▶ E A P S Q E G L L A P L S L L T L L T L A I A T L Y G L F L A S P G

NheI (1533)
EcoRI (1527)

1501 GCAGTAAGTCAAGAGGGGAAATATGATGAATTCGCTAGCTCGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAA
 313▶ Q •

HpaI (1696)

1601 ATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTA

MfeI (1707)
EcoRI (1792)

1701 ACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATCTA
 1801 AAATACAGCATAGCAAACTTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATG
 1901 TGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTCTCTCATTCTTTATGTTTAAAT

SspI (2031) **SwaI (2045)**

2001 GCACTGACCTCCACATTCCCTTTTTAGTAAAAATTCAGAAAATATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGA
 2101 TGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTGTGGACTTAGGAACAAGGAACCTTTAATAGAAAATGGACAGCAAGAAAGCGAGCTTCTAGC

SacI (2306)
BstXI (2335)

2201 TTTAGTTCCTGGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTC
 141▶ • N R T 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
 2301 AGAGATGAGTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAG
 108▶ S I L E 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

StuI (2470)
Eco147I (2470)

2401 TCCTTCTGCCGTTGCTCACAGCAGACCAATGGAATGGCTTACGACAGACAGTACCCTGCAATGTAGGCTCAATGTGGACAGCAGAGATGATCT
 74▶ D K Q G 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

