



PvuI (7)
SgfI (6) 1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
MfeI (82)
101 GAGAAGGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTCACGCGCCCGCCCTACCTGAGGCC
HindIII (245)
Bsu36I (291)
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTT

BstEII (555)
AgeI (552) **NcoI (560)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTACCATTGGGTGAAACTCTGGGAGATTCTCCTGTTGACCCAGA
1 M G E T L G D S P V D P E

601 GCATGGTGCCCTTCGCTGATGCACTGCCTATGAGCACTTCAAGAAATACCATGGTTGACACAGAGATGCCATTCTGGCCCACTTCGGAATCAGC
13 H G A F A D A L P M S T S Q E I T M V D T E M P F W P T N F G I S
XmnI (796)
701 TCTGTGGACCTCTCCGTGATGGAAGACCACTCGCATTCTTTGACATCAAGCCCTTTACCACAGTTGATTCTCCAGCATTCTGCTCCACTATGAAG
47 S V D L S V M E D H S H S F D I K P F T T V D F S S I S A P H Y E
801 ACATTCCATTCAAGAGCTGACCAATGGTTGCTGATTACAATATGACCTGAAGCTCAAGAATACCAAAGTGCATCAAAGTAGAACCTGCATCTCC
80 D I P F T R A D P M V A D Y K Y D L K L Q E Y Q S A I K V E P A S P

BspHI (939)
StuI (934)
901 ACCTTATTATTCTGAAAAGACCCAGCTCTACAACAGGCCTCATGAAGAACCTTCTAACTCCCTCATGGCCATTGAGTGCCGAGTCTGTGGGGATAAGCA
113 P Y Y S E K T Q L Y N R P H E E P S N S L M A I E C R V C G D K A

1001 TCAGGCTTCCACTATGGAGTTCATGCTTGTGAAGGATGCAAGGGTTTTTCCGAAGAACCATCCGATTGAAGCTTATTATGATAGGTGTGATCTTA
147 S G F H Y G V H A C E G C K G F F R R T I R L K L I Y D R C D L N
HindIII (1069)

BamHI (1103) 1101 GCCGGATCCAAAAAAGTAGAAATAAATGTCAGTACTGTGCTTTTTCAGAAAGTGCCTTGTGGGGATGTCTCACAATGCCATCAGGTTTGGGCGGAT
180 C R I H K K S R N K C Q Y C R F Q K C L A V G M S H N A I R F G R M
ScaI (1133)
BglII (1232) **EcoRV (1243)**
1201 GCCACAGCGCCGAGAAGGAGAAGCTGTTGGCGGAGACTCAGTGATATCGACCAGCTGAACCCAGAGTCTGCTGATCTGCGAGCCCTGGCAAAGCATTG
213 P Q A E K E K L L A E I S S D I D Q L N P E S A D L R A L A K H L
1301 TATGACTCATAATAAAGTCTTCCCGCTGACCAAAGCAAGCGAGGCGGATCTTGACAGGAAAGACAACGGACAAATCACCATTGTCATCTACGACA
247 Y D S Y I K S F P L T K A K A R A I L T G K T T D K S P F V I Y D

EcoRI (1401) 1401 TGAATTCCTTAATGATGGGAGAAGATAAATCAAGTTCAAACATATCACCCCTGCAGGAGCAGAGCAAAGAGGTGGCCATCCGAATTTTTCAAGGGTG
280 M N S L M M G E D K I K F K H I T P L Q E Q S K E V A I R I F Q G C
1501 CCAGTTTCGATCCGTAGAAGCGGTGCAAGAGATCACAGAGTATGCCAAAAATCCCTGGTTTCATTAACCTTGATTTGAATGACCAAGTACTCTGCTC
313 Q F R S V E A V Q E I T E Y A K N I P G F I N L D L N D Q V T L L

1601 AAGTATGGTGTCCATGAGATCATCTACAGATGCTGGCCTCCCTGATGAATAAAGATGGAGTCTCATCTCAGAGGGCCAAGGATTGATGACCCAGGGAGT
347 K Y G V H E I I Y T M L A S L M N K D G V L I S E G Q G F M T R E
1701 TCCTCAAAGCCTGCGGAAGCCCTTTGGTGACTTTATGGAGCCTAAGTTTGTGTTGCTGTGAAGTTCATGCACTGGAATTAGATGACAGTACTGCTGGC
380 F L K S L R K P F G D F M E P K F E F A V K F N A L E L D D S L A
BspHI (1685)

1801 TATATTTATAGCTGTATTATTCTCAGTGGAGACCGCCAGGCTTGTGACGTGAAGCCCATCGAGGACATCAAGACAACCTGCTGACGGCCCTGGAA
413 I F I A V I I L S G D R P G L L N V K P I E D I Q D N L L Q A L E
EcoO109I (1889)

1901 CTGACGCTCAAGCTGAATCACCCAGAGTCTCTCAGCTGTTCCGAAGGTGCTCCAGAAGATGACAGACCTCAGGAGATCGTACAGAGACGTCGACG
447 L Q L K L N H P E S S Q L F A K V L Q K M T D L R Q I V T E H V Q
Bsu36I (1968) **BbrPI (1990)**

2001 TACTGCATGTGATCAAGAAGACAGAGACAGACATGAGCCTTACCCCTGCTCCAGGAGATCTACAAGACTTGTATTAGCAGGAAAGTCCGCTAGCTGG
480 L L H V I K K T E T D M S L H P L L Q E I Y K D L Y •
BglII (2057) **NheI (2091)**
2101 CCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTAT

2201 TTGTAACCATTATAAGCTGCAATAAACAAAGTTAAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAA
HpaI (2229) **MfeI (2240)**

2301 GTAAAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACCTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGA
EcoRI (2325)

