



125

PvuI (7)
SgfI (6)
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
NgoMI (441)
NaeI (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGCTCAACTCTACGCTTTTGTTCGTTT

KasI (535) **AgeI (552)** **SphI (560)** **SfiI (576)**
Bsp120I (575) **DraIII (588)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCCTAGCTGAGATCACCGGTGACAGTGCAGCCGAGGTGGGCCAAGGGCCACGATGTGGCT
601 TGGAGTCTGCTGACCCCTTCTGCTCTGTTCAGCCTTGAGGGTCAAGAAAACCTTTTCAACATCAACAGTGTGACATGAAGAGCCTGCCGGACTGGACG
13> G V L L T L L L C S S L E G Q E N S F T I N S V D M K S L P D W T
701 GTGCAAAATGGGAAGAACCTGACCTGCAGTGTCTCGCGGATGTCAGCACCACCTCTCAGCTCAAGCCTCAGCACCAGATGCTGTTCTATAAGGATGACG
47> V Q N G K N L T L Q C F A D V S T T S H V K P Q H Q M L F Y K D D
801 TGCTGTTTTACAACATCTCTCCATGAAGAGCACAGAGATTTTTATTCTGAAGTCCGGATCTATGACTCAGGGACATATAAATGTACTGTGATTGT
80> V L F Y N I S S M K S T E S Y F I P E V R I Y D S G T Y K C T V I V
901 GAACAACAAAGAGAAAACCTGCGAGTACCAGGTGTTGGTGGAGGAGTGCCAGTCCAGGGTGACTGGACAAGAAAGAGGCCATCCAAGGTGGG
113> N N K E K T T A E Y Q V L V E G V P S P R V T L D K K E A I Q G G

EcoO109I (1034)
1001 ATCGTGAGGGTCAACTGTTCTGTCCAGAGGAAAAGGCCCAATACACTTCACAATTGAAAACTTGAACATAATGAAAAATGGTCAAGCTGAAAAGAG
147> I V R V N C S V P E E K A P I H F T I E K L E L N E K M V K L K R

XhoI (1109) **BamHI**
1101 AGAAGAATTCGAGACCAGAATTTGTGATACTGGAATTCGCCGTTGAGGAACAGGACCGGTTTTATCCTTCCGATGTCAAGCTAGGATCATTCTGG
180> E K N S R D Q N F V I L E F P V E E Q D R V L S F R C Q A R I I S G

NdeI (1204) **BstEII (1239)**
1201 GATCCATATGCAGACCTCAGAATCTACCAAGAGTGAAGTGGTACCGTACCGGAACTCTTCTACACCAAGTTCACATCAGCCCCACCGGAATGATC
213> I H M Q T S E S T K S E L V T V T E S F S T P K F H I S P T G M I

ApaI (1326)
1301 ATGGAAGGAGCTCAGCTCCACATTAAGTGCACCATTCAAGTGACTCACCTGGCCAGGAGTTCCAGAAATCATAATTGAGAAGGACAAGGGCATTGTGG
247> M E G A Q L H I K C T I Q V T H L A Q E F P E I I I Q K D K A I V

NcoI (1442) **Eco72I (1467)**
MscI (1439) **BbrPI (1467)**
1401 CCCACAACAGACATGGCAACAAGGCTGTGTAAGTGCAGTGTGATGGCCATGGTGGAGCACAGTGGCAACTACACGTGCAAAGTGGAGTCCAGCCGATATCCAA
280> A H N R H G N K A V Y S V M A M V E H S G N Y T C K V E S S R I S K
1501 GGTGAGCAGCATCGTGGTCAACATAACAGAACTATTTTCCAGCCGAAGTGAATCTTCTTACACATCTGGACCAAGTGAAGCTGAACCTGTCC
313> V S S I V V N I T E L F S K P E L E S S F T H L D Q G E R L N S
1601 TGCTCCATCCCAGAGCACCTCCAGCCAACCTCACCATCCAGAAGGAAGATACGATTGTGTACAGACTCAAGATTTCCAGAGATAGCTCAAAGTCCG
347> C S I P G A P P A N F T I Q K E D T I V S Q T Q D F T K I A S K S
1701 ACAGTGGGACGTATATCTGCACTGCAGGTATTGACAAAGTGGTCAAGAAAAGCAACACAGTCCAGATAGTCGATGTGAAATGCTCTCCCAACCAGGAT
380> D S G T Y I C T A G I D K V V K K S N T V Q I V V C E M L S Q P R I

ClaI (1857)
1801 TTCTTATGATGCCAGTTTGGAGTCATAAAAGGACAGACCATCGAAGTCCGTTGCGAATCGATCAGTGGAACTTTGCCTATTTCTTACCAACTTTAAAA
413> S Y D A Q F E V I K G Q T I E V R C E S I S G T L P I S Y Q L L K
1901 ACAAGTAAAGTTTTGGAGAATAGTACCAAGAATCAAAATGATCTGCGGTATTCAAAGACAACCCACTGAAGACGTGCAATACCAAGTGTGTTGCAGATA
447> T S K V L E N S T K N S N D P A V F K D N P T E D V E Y Q C V A D

BsiBI (2073)
BsaBI (2073)
2001 ATTGCCATCCCACGCCAAAATGTTAAGTGAAGTCTGAGGGTGAAGGTGATAGCCCCGTGGATGAGGTCCAGATTTCTATCCTGTCAAGTAAGGTGGT
480> N C H S H A K M L S E V L R V K V I A P V D E V Q I S I L S S K V V
2101 GGAGTCTGGAGGACATTGTGCTGCAATGTGCTGTGAATGAAGTATGCTGCTCCATCACCTATAAGTTTTACAGAGAAAAGAGGGCAACCCCTCTAT
513> E S G E D I V L Q C A V N E G S G P I T Y K F Y R E K E G K P F Y
2201 CAAATGACCTCAAATGCCACCCAGGCATTTTGGACCAAGCAGAAGGCTAA CAAGGAACAGGAGGGAGAGTATTACTGCACAGCCTTCAACAGAGCCAACC
547> Q M T S N A T Q A F W T K Q K A N K E Q E G E Y Y C T A F N R A N

NcoI (2353)
2301 ACGCTCCAGTGTCCCAGAAAGCAAATACTGACAGTCAAGTATTCTGCCCCATGGAAAGAAAGACTTATTGCAAGTGGTATCATCGGAGTATCAT
580> H A S S V P R S K I L T V R V I L A P W K K G L I A V V I I G V I I
2401 TGCTCTCTGATCATTGCGGCAAAATGTTATTTCTGAGGAAAGCCAAGGCAAGCAGATGCCAGTGGAAATGTCAGGCCAGCAGTACCCTTCTGAAC
613> A L L I I A A K C Y F L R K A K A K Q M P V E M S R P A V P L L N

Tth111I (2560)
2501 TCCAACAACGAGAAAATGTCAGATCCCAATATGGAAGCTAACAGTCAATTACGGTCAACAATGACGATGTCGAAACCATGCAATGAAACCAATAAATGATA
647> S N N E K M S D P N M E A N S H Y G H N D D V G N H A M K P I N D

