



PvuI (7)
SgfI (6) 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA **MfeI (82)**

101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC

Psp1406I (203) 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC **HindIII (245)**

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCCTCCCGCCTGTGGTGCCTCCTGAAGTCGCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTCTCAACTCTACGTCTTTGTTTCGTTT **NgoMIV (441)**

501 TCTGTTTGTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTGACCTGAGTGTGGAGGGGACAGGACCGGAGGAAAC **KasI (535)** **BstEII (555)** **AgeI (552)**
1 M A A V L E G Q E P E E T

601 TGCAGCTGCTGCAGAGGATGCTGCAACATCCACCTTGGAGGCTGTGGATGCCGGGCTGGAGCACCTTTCTCCCGGCTGGGAACAGTTGAACTTGGAC
13▶ A A A A E D A A T S T L E A V D A G P G A P F L P A G N Q L N L D
701 CTACGCCAGGGGCTGCCATCGCTACAGTACCTGTGTTCCAGCAGCCCCACAGTGTGCAAGTGGAGTTTTTGGCGCTCAGCACCCATGAGGACC
47▶ L R P G G G C H R L Q Y L C S Q P P Q L L Q V E F L R L S T H E D
801 CTCAGCTGCTGGATGACACCCTAGCCAAGGTTCCGTGGAGTCTGCTGCGCCTTCGCTCCCTGGTCTCAAAGTGGCCAGAGCCGGGTGCCCTGGGTGC
80▶ P Q L L D D T L A K V P W S L L R L R S L V L K G G Q S R G A L G A
901 CTGCTCCATGGAACCTGACCACTCTGCTGCTGGCCTGAGTGATTTGGCCTGCTGGCTCACCTAGACCTGAGCTTCAACAGGCTGGAGACGCTGCCT
113▶ C L H G G T L T T L P A G L S D L A C L A H L D L S F N R L E T L P
1001 ACCTGTGTCGGGAAGTGCAGGCTGGATGCGTTGCTGCTTTCTACAACACTCTCAGAGCTGCCCGAGGCGCTGGGGACCTGCCCGCCCTTACCT
147▶ T C V P E L H G L D A L L L S H N H L S E L P E A L G A L P A L T

1101 TCCTCACCGTGACACACAACCGCTAGAAAGACTACCCCTAACACTCGGGTCCCTGTCCACCCTTACGCGTCTCGATCTCTGAGAACCTTCTAGACAC **SmaI (1147)** **XbaI (1191)**
180▶ F L T V T H N R L E R L P L T L G S L S T L Q R L D L S E N L L D T

1201 CATAACCTCTGAGATTGAAACCTGCGCAGCCTCAGTGAAGTCAATCTGGCCTCAACCGGCTGCAGAGTCTCCAGCTTCCCTCGCGGGGCTGCGGTCC **FspI (1223)**
213▶ I P S E I G N L R S L S E L N L A S N R L Q S L P A S L A G L R S
1301 CTGCGGCTCCTTGTCTGCACAGCAACCTCCTGACCTCAGTGCCTGCTGGCCTGGTCCACTGCCACTGATCACTCGGCTTGAAGGACAAACCGG
247▶ L R L L V L H S N L L T S V P T G L V H L P L I T R L D L R D N R

1401 TCCGAGACCTGCCTGCTGAGTACTAGACGCTCCCTTTGTGCGCCTGCAGGGGAAACCTTGGGCGAGGCTCTCCAGCACCCCAAGTCCCCAGACAT **SdaI (1443)**
280▶ L R D L P A E L L D A P F V R L Q G N P L G E A S P A P P S P P D I
1501 ATCCAGGTTCCAGAAATGCCAGGCTATTGCTGACCTCAGATTTGGACAGCTTCTTGTGACTCCCCATGGCTGTTCTGTGACCCTGGCCTGTGGTGC
313▶ S Q V P E M P R L L L T S D L D S F L V T P H G C S V T L A C G V

1601 CGCCTACAGTTTCCAGCAGGAGCCACCACCGCTGTACCATCCACTATCGACTCTGGTTGCCGAGCCAGGCTGGTCTCTGCGGACCTCATGACT **BspHI (1692)**
347▶ R L Q F P A G A T T T P V T I H Y R L W L P E P G L V S L G P H D
1701 TCCTGCTTAGTAGTGTCTGGAGCTGCAGCCCCATGGGTGGCTTCCAGCAAGATGTGAGCTTATGGCTGCTTGTCCCTCCAGAGTCCGTCGCTG
380▶ F L L S S V L E L Q P H G V F Q Q D V S L W L L F V P P R V R R C
1801 TCGTGAGGTAGTTGTGAGGACACGGAGTAACAACCTGGAATGACCTAGAGACCAACTGGAGGAAGAGGCACCCAAGAGGCTCTGGCTCGCTGCCAG
413▶ R E V V V R T R S N N T W N D L E T Q L E E E A P K R L W A R C Q
1901 GTGCCCACTTCTCTGGTTCTGGTCTTTTACGCCAGTATCCAACACTTGCCTTTTGCACCAGAGGGGAGCACTTCTGCTCCTCAGGTCATCCTG
447▶ V P H F S W F L V V L R P V S N T C L L P P E G A L L C S S G H P

2001 GGGTCAGAGTCACTTTTCCCTGGGTTACAGAAGAGCCTCGGCAAGTCTCCATGCAGGTAGTGCATATGGCTGGACTAGAGCTGAGAACCCTTCTGGA **NdeI (2065)**
480▶ G V R V T F P P G V T E E P R Q V S M Q V V H M A G L E L R T L L E
2101 AGAATCAGAGGCATCAGTAAGCCCTGCTGTGCTTTTACAGAGCGGCCCCCAGCTTCTACAACCTGTCACTGTGAGTTGCCCTGCCTCCTGGC
513▶ E S E A S V S P L L C L S Q S G P P S F L Q P V T V Q L P L P P G

2201 GTTACAGGCTCAGTCTAGACCGCTCCACCTGCATCTGCTTACCGAACCCCTGACAACCCTGGGATGATATCACCCTCAGGTTGGCACTGGAAT **XbaI (2214)** **EcoRV (2272)**
547▶ V T G F S L D R S H L H L L Y R T P L T T T W D D I T T Q V A L E

2301 TCACCCACTGTATGCACGTTCCAGGTACGCACCTTCTCCTGGTACTGGCTCTGGTATACCACAAAACCTGTGTGGGGGGCTGGCCCGAAGGCCTG **XcmI (2363)** **Bst1107I (2355)** **SfiI (2385)**
580▶ F T H L Y A R F Q V T H F S W Y W L W Y T T K T C V G G L A R K A W

2401 GGAACGGTGCAGCTGCACCGTGTGAATCTCATTGCACTGCAGAGGCGCCGAGACCTGAGCAGGTTCTGCTGCAGTGCCTGCCACGGAACAAGTAGAT **DraIII (2416)** **KasI (2444)**
613▶ E R L R L H R V N L I A L Q R R R D P E Q V L L Q C L P R N K V D
2501 GCCACCCTGAGTCGGCTGCTGGTTCGCTACCGTGGCCCTGAACCTCTGAGACTGTGGAGATGTTTGGGGTGAAGTCTTTGACGCTTTGAGCGGG
647▶ A T L S R L L V R Y R G P E P S E T V E M F E G E K F F A A F E R

2601 GCATTGATGATAGTGCAGACCGTCCAGACTGTGTGGATGGCAGAATCTGCTTTTGTCTTATTACACCTGAAGAAGCTAAAGGAGGTATACATCACACC **BbsI (2653)** **Bst1107I (2686)**
680▶ G I D V D A D R P D C V D G R I C F V F Y S H L K N V K E V Y I T T

2701 AGCCTTGGACCGGGAAGCTCAGGATGTTGAGGACAGGTGTCTTTTATCGGGGTTCACTTCTGTGGAGTACCTGCAGAGGCAAGGCTGCCCGCAG **PshAI (2732)** **Acc65I (2769)**
713▶ A L D R E A Q D V R G Q V S F Y R G S L P V E V P A E A E A A R Q
2801 AGGAAGGTACAGATGCACTGTGGATGGCCACCTTGCCTCAAGCTACCGAGACTCCGGGTGCCAGGGGAGTGGGACGGGACTGACTTCTCCTTGA
747▶ R K G T D A L W M A T L P I K L P R L R G A O G S G O G T D F S L

