



**PvuI (7)**  
**SgfI (6)**  
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA  
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

**Psp1406I (203)** **HindIII (245)** **Bsu36I (291)**  
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCGCTACCTGAGGCC  
301 GCCATCCACGCGGGTTGAGTCGCGTTTCTGCCGCCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

**NgoMIV (441)** **NaeI (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

**KasI (535)** **AgeI (552)** **BspLU11I (560)**  
501 TCTGTTTCTGCGCGCTTACAGATCCAAGCTGTGACCGCGGCTACCTGAGATCACCGGTCAACATGTCCCTGCTTTTTCTCGATGCAACTCCATCGTCAC  
1▶ M S L L F S R C N S I V T  
601 CGTCAAGAAGGATAAGCGACACATGGCTGAAGTGAATGCTTCCCTCTCAAGCACTTTGTCACTGCCAAGAAAAGATCAATGGAATCTTTGAGCAGCTG  
13▶ V K K D K R H M A E V N A S P L K H F V T A K K K I N G I F E Q L

**BbsI (733)** **BstEII (765)**  
701 GGGCCTACATCCAAGAGAGCGCCAGCTTCTTGAAGACCCACAGGAACACAGAAGTGGACCCGGTTACCACGGAAGAGCAGGTCCTGGACGTCAAAG  
47▶ G A Y I Q E S A S F L E D T H R N T E L D P V T T E E Q V L D V K

**Asp718I (800)** **Acc65I (800)**  
801 GGTACCTGTCCAAGGTCAGGGGTATCAGCGAAGTGTGGCCAGGCGGCACATGAAGTGGCTTTTTTTGGCCGGACGAGCAATGGGAAGAGCACCGTGAT  
80▶ G Y L S K V R G I S E V L A R R H M K V A F F G R T S N G K S T V I

**StuI (989)**  
901 CAATGCCATGCTCTGGGACAAAGTTCTGCCATCTGGGATTGGTATACCACCAATTGCTTCTCGGGTGGGGGACAGATGGCCATGAGGCCCTTCTCCT  
113▶ N A M L W D K V L P S G I G H T T N C F L R V G G T D G H E A F L  
1001 CTCACAGAGGGCTCAGAAGAGAAGAGTGTCAAGACTGTGAACCAACTGGCCATGCCCTCCATCAGGACGAGCAGTTGCATGCAGGCAGCATGGTGA  
147▶ L T E G S E E K K S V K T V N Q L A H A L H Q D E Q L H A G S M V

**BsrBI (1126)** **ClaI (1168)**  
1101 GTGTGATGTGGCCAACTCCAAGTGTCCGCTCCTGAAGATGACCTCGTGTGATGGACAGCCCTGGGATCGATGTTACCACGAGCTGGACAGCTGGAT  
180▶ S V M W P N S K C P L L K D D L V L M D S P G I D V T T E L D S W I

**DraIII (1285)**  
1201 TGATAAGTTTTGCCTGGATGCTGATGTGTTTGTGCTGGTGGCCAACTCAGAGTCCACGCTGATGCAGACGGAGAAGCAGTTCTTCCACAAAGTGAAGTGA  
213▶ D K F C L D A D V F V L V A N S E S T L M Q T E K Q F F H K V S E

**Eco47III** **AfeI (1397)**  
1301 CGTCTCTCCCGCCCAACATCTTCATCCTGAACAACCGCTGGGATGCGTCTGCCTCGGAGCCTGAGTACATGGAGGAGGTGCGGCGGCAGCACATGGAGC  
247▶ R L S R P N I F I L N N R W D A S A S E P E Y M E E V R R Q H M E

**XcmI (1406)**  
1401 GCTGCACCAGCTTTCTGGTGGATGAGCTGGGCGTGGTGGATCGAGCTCAGGCTGGGGACCGGATCTTCTCGTGTCTGCCAAGGAGGTTCTCAGCGCCAG  
280▶ R C T S F L V D E L G V V D R A Q A G D R I F F V S A K E V L S A R  
1501 GGTCCAGAAAGCCAGGGCATGCCAGAAGGAGGCGGCTCTCGCAGAAAGTTTCAAGTGVAGGATGTTTGAAGTTTTCAGAAATTCAGAGGAGGAGGTTGAG  
313▶ V Q K A Q G M P E G G G A L A E G F V R M F E F Q N F E R Q F E  
1601 GAGTGCATTTCCAGTCTGCAGTAAAGACCAAATTTGAGCAGCACACAGTCCGGGCAAGCAGATTGCAGAGGCGGTTCTGTCATCATGGATCCCTGC  
347▶ E C I S Q S A V K T K F E Q H T V R A K Q I A E A V R L I M D S L

**BsrBI (1747)**  
1701 ACATCGCAGCTCAGGAGCAGCGGGTTATTGCCTAGAAATGCGGGAAGAGCGGCAAGACCGGCTGAGGTTTATTGACAAGCAGCTGGAGCTCCTGGCTCA  
380▶ H I A A Q E Q R V Y C L E M R E E R Q D R L R F I D K Q L E L L A Q

**KasI (1877)**  
1801 AGACTACAAGCTGCGAATTAAGCAGATTACGGAGGAAGTGAAAGGCAGGTGCCACAGCCATGGCTGAAGAGATCAGGCGCCTCTGTGCTAGTTGAC  
413▶ D Y K L R I K Q I T E E V E R Q V S T A M A E E I R R L S V L V D  
1901 GAGTACCAGATGGACTTCCACCCATCCCGAGTTGTCTCAAGTTTATAAGAAGCAGCTGCACCGCCATATAGAGGAAGGCTGGGCCGGAACCTGTCTG  
447▶ E Y Q M D F H P S P V V L K V Y K N E L H R H I E E G L G R N L S  
2001 ACCGCTGTCCACTGCCATTGCCAGTTCAGTGCAGACTATGCAGCAGGACATGATAGACGGCTTGAAGCCCCTTCTCTGTATCTATGCGGAATCAGAT  
480▶ D R C S T A I A S S L Q T M Q Q D M I D G L K P L L P V S M R N Q I

**XcmI (2190)**  
2101 AGACATGCTGGTCCCTCGACAGTGTCTCCCTCAGCTATGACCTGAATTGTGACAAGCTGTGTGCTGACTTTCAGGAGGACATCGAGTTCCACTTCTCC  
513▶ D M L V P R Q C F S L S Y D L N C D K L C A D F Q E D I E F H F S

**Bsp120I (2232)**  
2201 CTTGGATGACTATGCTAGTGAACAGGTTCTGGGCCCCAAGAATAGCCGCGGGCCTTGTAGGCTACAGTGATCAGGTTTACAGGCTCCTCCTCCTCTGA  
547▶ L G W T M L V N R F L G P K N S R R A L L G Y S D Q V Q R P L P L  
2301 CACGTCCAACCCAGCATGCCCCCTTGCACAGAGCTCCCTCACAGAGGAGTCAATGGTCTCCATGGTTACTGGCTGGCCTCTCTGAGCTCTAG  
580▶ T P A N P S M P P L P Q S S L T Q E E L M V S M V T G L A S L T S R  
2401 GACCTCCATGGGCATTCTTGTGGTGGAGGAGTGGTGTGGAAGGCAGTGGGCTGGAGACTCATCGCCCTCTCCTTTGGACTGTATGGCCTCCTGTACGTC  
613▶ T S M G I L V V G G V W K A V G W R L I A L S F G L Y G L L Y V  
2501 TATGAGCGACTGACCTGGACCACAAAGCCAAAGAGAGGGCTTCAAGCGCCAGTTTGTGGAATACGCCAGTGAAGAAGTACAGCTCATCAGTTCAGTTACA  
647▶ Y E R L T W T T K A K E R A F K R Q F V E Y A S E K L Q L I S Y  
2601 CCGGCTCTAACTGCAGCCACCAAGTCCAGCAGGAATTGTCTGGACATTTGCTCATCTGTGCCAGCAAGTTGACATCACCCGAGATAATCTGGAGCAGGA  
680▶ T G S N C S H Q V Q Q E L S G T F A H L C Q Q V D I T R D N L E Q E

