



**PvuI (7)**  
**SgfI (6)** 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
**MfeI (82)**  
101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGTGCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**HindIII (245)**  
**Psp1406I (203)** 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCGCCCTACCTGAGGCC  
**PvuII (239)**  
**Bsu36I (291)**  
301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

---

**NgoMIV (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT

---

**AgeI (552)** 501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACCGTAGGAGGGCCATCATGAGACTCATGAGAAACATTTACATATTT  
**BspHI (568)**  
601 TGTAGTATTGTTATGACAGCAGAGGGTGTGCTCCAGAGTCCAGAGAAAGGGAAGTGTGACCAACTGCTCCAACATGTCTCTAAGAAAGGTTCCCG  
11▶ C S I V M T A E G D A P E L P E E R E L M T N C S N M S L R K V P  
1▶ M R L I R N I Y I F

**XcmI (715)**  
701 CAGACTTGACCCAGCCACAACGACTGGATTTATCTATAACCTCCTTTTCAACTCCAGAGTTCAGATTTTATTCTGTCTCCAACTGAGAGTTTT  
44▶ A D L T P A T T T L D L S Y N L L F Q L Q S S D F H S V S K L R V L

**PvuII (825)**  
801 GATTCTATGCCATAACAGAATTCAACAGCTGGATCTCAAAACCTTTGAATTCAACAAGGAGTTAAGATATTTAGATTTGTCTAATAACAGACTGAAGAGT  
77▶ I L C H N R I Q Q L D L K T F E F N K E L R Y L D L S N N R L K S

**BglII (935)**  
901 GTAACCTGGTATTACTGGCAGGTCTCAGGTATTTAGATCTTTCTTTAATGACTTTGACACCATGCCTATCTGTGAGGAAGCTGGCAACATGTCACACC  
111▶ V T W Y L L A G L R Y L D L S F N D F D T M P I C E E A G N M S H

**AvrII (1007)**  
1001 TGGAAATCCTAGTTTGTAGTGGGGCAAAAATACAAAATCAGATTTCCAGAAAATGCTCATCTGCATCTAAATACTGCTTCTTAGGATTGAGAAGTCT  
144▶ L E I L G L S G A K I Q K S D F Q K I A H L H L N T V F L G F R T L  
1101 TCCTCATTATGAAGAAGTAGCCTGCCATCTTAAACACAACAAAAGTGCACATTGTTTTACCAATGGACACAAAATTTCTGGGTTCTTTTTCGTTGATGGA  
177▶ P H Y E E G S L P I L N T T K L H I V L P M D T N F W V L L R D G  
1201 ATCAAGACTTCAAAAATATTAGAAATGACAAATATAGATGGCAAAAGCCAATTTGTAAGTTATGAAATGCAACGAAATCTTAGTTTAGAAAATGCTAAGA  
211▶ I K T S K I L E M T N I D G K S Q F V S Y E M Q R N L S L E N A K

**XmnI (1383)**  
1301 CATCGTTCTATTGCTTAATAAAGTTGATTTACTCTGGGACGACCTTTTCTTATCTTACAATTTGTTGGCATAATCATGAGTGAACACTTTTCAGATCCG  
244▶ T S V L L L N K V D L L W D D L F L I L Q F V W H T S V E H F Q I R

**BsrGI (1487)**  
1401 AAATGTGACTTTTGGTGGTAAGGCTTATCTTGACCACAATTCATTTGACTACTCAAATACTGTAATGAGAAGTATAAAAATGGAGCATGTACATTTTCAGA  
277▶ N V T F T G G K A Y L D H N S F D Y S N T V M R T I K L E H V H F R  
1501 GTGTTTACATTCACAGGATAAAATCTATTTGCTTTTGACCAAAAATGGACATAGAAAACCTGACAATATCAAATGCACAAATGCCACACATGCTTTTCC  
311▶ V F Y I Q Q D K I Y L L T K M D I E N L T I S N A Q M P H M L F  
1601 CGAATTATCTACGAAATCCAATATTTAAATTTGCAATAATATCTAACAGACGAGTTTTTTAAAAGAAGTATCCAAGGCTCCTCACTGAAAACCTCT  
344▶ P N Y P T K F Q Y L N F A N N I L T D E F F K R T I Q L P H L K T L  
1701 CATTTTGAATGGCAATAAACTGGAGACACTTTCTTTAGTAAGTTGCTTTGCTAACAACACACCCTTGAACACTTGGATCTGAGTCAAATCTATTACAA  
377▶ I L N G N K L E T L S L V S C F A N N T P L E H L D L S Q N L L Q

**MscI (1824)**  
1801 CATAAAAATGATGAAAATGCTCATGGCCAGAACTGTGGTCAATATGAATCTGTACATAAATTTGCTGATTCTGTCTCAGGTGCTTGCCAAAA  
411▶ H K N D E N C S W P E T V V N M N L S Y N K L S D S V F R C L P K  
1901 GTATTCAAAATCTTGACCTAAAATAACCAAAATCCAAAATGTACCTAAAGAGACTATTCATCTGATGGCCTTACGAGAAGTAAATATTGCATTTAATTT  
444▶ S I Q I L D L N N N Q I Q T V P K E T I H L M A L R E L N I A F N F  
2001 TCTAACTGATCTCCCTGGATGCAGTCATTTAGTAGACTTTAGTCTGAACATTGAAATGAACTTATTCTCAGCCATCTCTGGATTTTGTTCAGAGC  
477▶ L T D L P G C S H F S R L S V L N I E M N F I L S P S L D F V Q S

**BspHI (2194)**  
2101 TGCCAGGAAGTTAAACTCTAAATGCGGGAAGAAATCCATTCCGGTGTACCTGTGAATTAATAAATTTTATTTCAGCTTGAACATATTTCAGAGGTCATGA  
511▶ C Q E V K T L N A G R N P F R C T C E L K N F I Q L E T Y S E V M

**Bsu36I (2243)**  
2201 TGGTTGGATGGTCAGATTCATACACCTGTGAATACCTTTAAACCTAAGGGGAACTAGGTTAAAAGACGTTTCTCCACGAATTATCTTGAACACAGC  
544▶ M V G W S D S Y T C E Y P L N L R G T R L K D V H L H E L S C N T A  
2301 TCTGTTGATTGTACCATTGTGGTTATTATGCTAGTCTGGGTTGGCTGTGGCCTTCTGCTGTCTCCACTTTGATCTGCCCTGGTATCTCAGGATGCTA  
577▶ L L I V T I V V I M L V L G L A V A F C C L H F D L P W Y L R M L

**MscI (2475)**  
**NheI (2469)**  
2401 GGTCAATGCACACAAACATGGCACAGGGTTAGGAAAACAACCAAGAACAACCTCAAGAGAAATGTCTAAGCTAGCTGGCCAGACATGATAAGATACATTG  
611▶ G Q C T Q T W H R V R K T T Q E Q L K R N V •  
2501 ATGAGTTTGGACAAACCAACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAA

**HpaI (2607)** MfeI (2618)  
 2601 TAAACAAGTTAAACAACAACAAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGT  
 2701 ATGGAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGC  
**SapI (2885)**  
 2801 TGTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTGAAGCTAGCTCTTCATTTCTTT  
 2901 ATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGG  
**EcoO109I (3017)**  
 3001 CAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAAGC  
 3101 GAGCTTCTAGCTTTAGTTCCTGGTGTACTTGGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCA  
 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  
**SacI (3217)** **BstXI (3246)**  
 3201 GGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAA  
 111 P A Y D 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  
**StuI (3381)**  
 3301 TGGTGTCAAAGTCCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGC  
 78 T D F 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  
 3401 AGAGATGATCTCCCAGTCTGGTCTGATGGCCGCCGACATGGTCTTGTGTCCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGC  
 45 S I I E G T K T R I A A G V H H K N D E Y L M T I K E T A V E V L  
**XmnI (3523)** **AseI (3589)**  
 3501 TCCAGATCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCA  
 11 E L D Q Q S I N F T K M  
**SacI (3646)**  
 3601 AAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTGCGTCA  
**SpeI (3744)**  
 3701 ATGGGGCGGAGTTGTTACGACATTTGGAAAGTCCCGTTGATTTACTAGTCAAACAAACTCCCATGACGTCAATGGGGTGGAGACTTGGAAATCCCCG  
**SnaBI (3872)**  
 3801 TGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAG  
**NdeI (3977)**  
 3901 TCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTG  
 4001 CCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGG  
**PaeI (4163)**  
**PstI (4156)**  
**SdaI (4155)**  
 4101 GCGGGGGTGTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAAITAAAGAACATGTGAGCAAAGGCCAGCAAAGG  
 4201 CCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAA  
 4301 CCCGACAGGACTATAAAGATACCAGCGCTTTCCCCTGGAAGCTCCCTCGTGCCTCTCTGTCCGACCCTGCCGCTTACCGGATACCTGTCCGCCCTT  
**ApaLI (4487)**  
 4401 CTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGTTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCC  
 4501 CCGTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTGAGTCAACCCGGTAAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAG  
 4601 GATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTG  
 4701 CTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTA  
 4801 CGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGC  
**EagI (4923)**  
**NotI (4922)**  
 4901 TAGTTAATTAACATTTAAATCAGCGGCCGAATAAAATATCTTTATTTTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTC  
**PaeI (4903)**  
 5001 TCCATCAAAACAAAACGAAACAAAACAAACTAGCAAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA