



**PvuI (7)**  
**SgfI (6)**
**MfeI (82)** **EcoNI (96)**

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**Psp1406I (203)**
**HindIII (245)** **Bsu36I (291)**

201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCCCTACCTGAGGCC

301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

---

**NgoMIV (441)**

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

---

**KasI (535)**
**NcoI (560)**

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTCCACCATGGTGAATTACTGCAAGCAAATTGTTCTTCTCTCTGG

601 ATTAGAATATATGAATGACTATAACTTTAGAGCACTTAAGTCCCTTACTGAACCATGACCTAAAAGTACTGAAACATGCAAGTGAAGTATTGACGAATT

13▶ L E Y M N D Y N F R A L K S L L N H D L K L T K N M Q D D Y D R I

**XmnI (769)**
**BbsI (769)**

701 AACATTGCTGACTTAATGGAAGAAAAGTCCAGAGGATGCTGGATTGAGCAAACGATAGAAGTATGTAAGACATTCCAGAACTTGCAGCTCGTGTGG

47▶ N I A D L M E E K F P E D A G L S K L I E V C E D I P E L A A R V

**BstXI (878)**
**HindIII (890)**

801 ATATACTTAGAAAAGAGATGGAGAAAAGTAAAAATAAAACAAAAATAAAATCAGAATCCAGCCCTCCTCCACTTACCTCCAGCCTGATGGAAGCTTGGGA

80▶ D I L R K E M E K V K N K T K I K S E S S P P P L T S S L M E A W E

**MscI (983)**

901 AGTAGAACAGCTATGGTACAGCTTCATCAGAGGTACACCGTCTAAGATCTCACTGAAAAGGCATCCACTAAGCTCGCTAGCTGGCCAGACATGATAAG

113▶ V E P A M V T A S S E V H R L R S H •

1001 ATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATA

---

**HpaI (1115)**
**MfeI (1126)**

1101 AGCTGCAATAAACAGTTAAACAACAACATTGCATTCAATTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACA

---

**EcoRI (1211)**

1201 AATGTGGTATGGAATCTAAAATACAGCATAGCAAACCTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCA

---

**SapI (1393)**

1301 TCAGGGGCTGTTGCCAATGTCATTAGCTGTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTAGCTCTTC

---

**SspI (1450)**
**SwaI (1464)**

1401 ATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATGAAAATATTGAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTT

---

**EcoO109I (1525)**

1501 TTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGC

---

1601 AAGAAAGCGAGCTTCTAGCTTTTAGTTCTGCTGACTTGGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTTCATCTCAATGAGCACAA

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

**SacI (1725)**
**BstXI (1754)**

1701 AGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCGACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGAC

114▶ C D 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

**StuI (1889)**

1801 AGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTACGACAGACAGTGCCTGCCAATGTAGGCTCAATG

81▶ A V I 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

1901 TGGACAGCAGAGATGATCTCCCGTCTGGTCTGATGGCCGCCGACATGGTCTGTTGCTCCTCATAGAGCATGGTGTCTCTCAGTGGCGACCT

47▶ H V A 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

**BbsI (2035)**
**AseI (2097)**

2001 CCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATT

14▶ V L E L D Q Q S I N F T K M

**SacI (2154)**

2101 AATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCCAT

---

**SpeI (2252)**

2201 TTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCCATTTGACGTCAATGGGGTGGAGACTTGGGA

---

**SnaBI (2380)**

2301 AATCCCCGTGAGTCAAACCCTATCCACGCCATTGATGTAAGTGCACAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTGCACAA

**NdeI (2485)**

2401 TAGGAAAGTCCCATAAGGTCATGTA  
2501 ATGTA

**PacI (2671)**

**PstI (2664)**

**SdaI (2663)**

**BspLU11I (2681)**

2601 GTCAATGGGCGGGGTCGTTGGGCGT  
2701 GCAAAAGGCCAGGAACCGTAAAAAG  
2801 TGGCGAAACCCGACAGGACTATAAAG

**ApaLI (2995)**

2901 CCGCCTTTCTCCCTTCGGGAAGCGT  
3001 CGAACCCCGTTAGCCCGACCGCTGCG  
3101 GGTAACAGGATTAGCAGAGCGAGGTA  
3201 GCGCTCTGCTGAAGCCAGTTACCTTC  
3301 GCAGATTACGCGCAGAAAAAGGATCT

**EagI (3431)**

**PacI (3411)**

**SwaI (3420)**

**NotI (3430)**

3401 GTCATGGCTAGTTAATTAACATTTAAATC  
3501 ATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTG