



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

1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**Psp1406I (203)**
**HindIII (245)**
**EcoNI (287)**

201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCCGCCCTACCTGAGGCC  
 301 GCCATCCAGCGCGGTTGAGTCGCGTTTCTGCCGCCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGTCAAGTGCAGACC

---

**NgoMIV (441)**  
**NgoMI (441)**  
**NaeI (441)**

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

---

**PciI (560)**  
**BspLU11I (560)**

501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTCAACATGTGGCAGCTGCTCCTCCCACTGCTCTGCTACTTCT  
1▶ M W Q L L L P T A L L L L

---

**SphI (611)**
**BglII (624)**
**XhoI (673)**

601 AGTTTCAGCTGGCATGCGGACTGAAGATCTCCCAAAGGCTGTGGTGTCTGGAGCCTCAATGGTACAGGGTGTCTGAGAAGGACAGTGTGACTCTGAAG  
 13▶ V S A G M R T E D L P K A V V F L E P Q W Y R V L E K D S V T L K

---

**XhoI (774)**  
**StuI (770)**  
**Eco147I (770)**

701 TGCCAGGGAGCCTACTCCCTGAGGACAATCCACACAGTGGTTTCACAATGAGAGCCTCATCTCAAGCCAGGCTCGAGCTACTTCATTGACGCTGCCA  
 47▶ C Q G A Y S P E D N S T Q W F H N E S L I S S Q A S S Y F I D A A

---

**SalI (802)**

801 CAGTCGACGACAGTGGAGAGTACAGGTGCCAGACAAACCTCTCCACCCTCAGTGACCCGGTGCAGCTAGAAGTCCATATCGGCTGGCTGTTGCTCCAGGC  
 80▶ T V D D S G E Y R C Q T N L S T L S D P V Q L E V H I G W L L L Q A  
 901 CCCTCGGTGGGTGTTCAAGGAGGAAGACCTATTACCTGAGGTGTACAGCTGGAAGAACTGCTCTGCATAAGGTACATATTTACAGAATGGCAAA  
 113▶ P R W V F K E E D P I H L R C H S W K N T A L H K V T Y L Q N G K

---

**PstI (1074)**

1001 GGCAGGAAGTATTTTCATCATAATTCTGACTTCTACATTCAAAAGCCACACTCAAAGACAGCGGCTCCTACTTCTGCAGGGGCTTTGTTGGGAGTAAAA  
 147▶ G R K Y F H H N S D F Y I P K A T L K D S G S Y F C R G L V G S K

---

**Asp718I (1178)**  
**Acc65I (1178)**

1101 ATGTGCTTTCAGAGACTGTGAACATCACCATCACTCAAGGTTTGGCAGTGTCAACCATCTCATATTCTTTCCACCTGGGTACCAAGTCTCTTTCTGCTT  
 180▶ N V S S E T V N I T I T Q G L A V S T I S S F F P P G Y Q V S F C L

---

**BstBI (1259)**  
**AsuII (1259)**

1201 GGTGATGGTACTCCTTTTTGCGAGTGGACACAGGACTATATTTCTGTGGAAGACAAACATTCGAAGCTCAACAAGAGACTGGAAGGACCATAAATTTAAA  
 213▶ V M V L L F A V D T G L Y F S V K T N I R S S T R D W K D H K F K

---

**MscI (1345)**  
**NheI (1339)**

1301 TGGAGAAAGGACCCTCAAGACAAATGACCCCCATCCCATGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGC  
 247▶ W R K D P Q D K •

---

**HpaI (1477)**
**MfeI (1488)**

1401 AGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAAATTGCATTCA

---

**EcoRI (1573)**

1501 TTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATTCTAAAATACAGCATAGCAAAC  
 1601 TTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGC  
 1701 CTCACCTTCTTTCATGGAGTTTAAAGATATAGTGTATTTTCCAAAGTTTGAAGTACTCTTCATTCTTTATGTTTTAAATGCACTGACCTCCACATTC

---

**SspI (1812)**

1801 CCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAA  
 1901 TATCCCCAGTTTAGTAGTTGACTTAGGAACAAAGGAACCTTTAATAGAAATTTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTT  
141▶ • N R T Y K

---

**SacI (2087)**

2001 GAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCAC  
 135▶ L P I L E E I T T K V L K G N M E I L V F C D P A Y D S I L E R C

---

**BstXI (2116)**

2101 ATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCA  
 101▶ M G C P S V V R I S R D V E D S Y P H R V A V I T D F D K Q G N S V

---

**StuI (2251)**  
**Eco147I (2251)**

2201 CAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCTGAT  
 68▶ A S G I A I A E A C V T V R G I Y A E I H V A S I I E G T K T R I

