



**PvuI (7)**  
**SgfI (6)**

1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGTGCTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**HindIII (245)**  
**Psp1406I (203)**

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

---

**PvuII (239)**  
**Bsu36I (291)**

401 GGGCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT

---

**NgoMIV (441)**

501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCACCATGGTGTCTTTGCCTCGGCTGTGCGCCTATGGGGCTG

1 M V S L P R L C A L W G C

---

**BstEII (555)**  
**AgeI (552)**

601 CTTGTTGACAGCGGTCCACTAGGGCAGTGTGTACTGTCAGTGACAAACAGTACCTCCACGATGGCCAGTGTGTATTTGTCCAGCCAGGAAGCCGA

13 L L T A V H L G Q C V T C S D K Q Y L H D G Q C C D L C Q P G S R

---

**BssHII (584)**  
**BbsI (726)**

701 CTGACAAGCCACTGCACAGCTCTTGAGAAGACCAATGCCACCATGTGACTCAGCGGAATTCTCAGCCAGTGAACAGGGAGATTCGCTGTCAACAGC

47 L T S H C T A L E K T Q C H P C D S G E F S A Q W N R E I R C H Q

---

**EcoRI (757)**  
**Tth11I (859)**

801 ACAGACACTGTGAACCAATCAAGGGCTTCGGGTTAAGAAGGAGGGCACCGCAGAATCAGACACTGTCTGTACCTGTAAGGAAGGACAACACTGCACCAAG

80 H R H C E P N Q G L R V K K E G T A E S D T V C T C K E G Q H C T S

901 CAAGGATTGCGAGGCATGTGCTCAGCACACGCCCTGTATCCCTGGCTTTGGAGTTATGGAGATGGCCACTGAGACCACTGATACCGTCTGTATCCCTGC

113 K D C E A C A Q H T P C I P G F G V M E M A T E T T D T V C H P C

---

**BstAPI (899)**  
**BstBI (1033)**

1001 CCAGTCGGCTTCTTCCAATCAGTCATCACTTTTCGAAAAGTGTATCCCTGGACAAGCTGTGAGGATAAGAAGTGGAGGCTACAGAAAGGAACGA

147 P V G F F S N Q S S L F E K C Y P W T S C E D K N L E V L Q K G T

---

**PshAI (1104)**  
**BsaBI (1164)**

1101 GTCAGACTAATGTCACTGTGGTTAAAGTCCCGGATGCGAGCCCTGCTGGTCATTCTGTGATGGGCATCCTCATCACCATTTTCGGGGTGTCTTCT

180 S Q T N V I C G L K S R M R A L L V I P V V M G I L I T I F G V F L

1201 CTATATCAAAAAGTGGTCAAGAAACCAAGGATAATGAGATGTTACCCCTGCGGCTCGACGGCAAGATCCCCAGGAGATGGAAGATTATCCCGGTCAT

213 Y I K K V V K K P K D N E M L P P A A R R Q D P Q E M E D Y P G H

---

**BsrBI (1387)**

1301 AACACCGTGTCTCCAGTGCAGGAGACTGCACGGGTGTCAGCCTGTACACAGGAGGATGGTAAAGAGAGTCGCATCTCAGTGCAGGAGCGGAGGTTGA

247 N T A A P V Q E T L H G C Q P V T Q E D G K E S R I S V Q E R Q V

---

**NheI (1448)**

1401 CAGACAGCATAGCCTTGAGGCCCTGGTCTGAACCTGGAAGTCTTTGCTAGCTGCCAGACATGATAAGATACATTGATGAGTTTGACAAACCACAA

280 T D S I A L R P L V •

---

**HpaI (1586)**

1501 CTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACA

---

**MfeI (1597)**  
**EcoRI (1682)**

1601 TTGCATTCAATTTATGTTTCAGTTTCAGGGGAGGTGTGGGAGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTCAAAAATACAGCA

1701 TAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCT

---

**SapI (1864)**

1801 GTTTGCAGCCTCACCTTCTTTTATGAGTAAAGATATAGTGTATTTTCCAAGGTTTGAAGTCTTCTTCTTTATGTTTTAAATGACTGACCT

---

**SspI (1921)**  
**SwaI (1935)**

1901 CCCACATTCCCTTTTAGTAAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGC

2001 CCTTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAACCTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTC

141 • N R

---

**SacI (2196)**

2101 GGTGACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGC

138 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 P A Y D S I L

---

**BstXI (2225)**

2201 TCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATGTGCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCC

104 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 T D F D K Q G

---

**StuI (2360)**

2301 CGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTT

71 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 S I I E G T K

