



**PvuI (7)**  
**SgfI (6)** **MfeI (82)**  
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**Psp1406I (203)** **HindIII (245)**  
**PvuII (239)** **Bsu36I (291)**  
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCAGAGGGCTCGCATCTCTCTTACCGCGCCCGCCGCCCTACCTGAGGCC  
301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

---

**NgoMIV (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

---

**KasI (535)** **AgeI (552)** **BspHI (560)**  
501 TCTGTTTGGCGCGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTCATCATGAAACTAGACATGACTGGGGACTGCACGCCAGTGC  
601 GGTGCTGATGGCCGAGTGTGACCGTACTGGAGCAGTTCCTGTCCAGGCTCCACGGGGCTCTCCCGGATGCAAGGGGCTGCCACATAGCCAGTTC  
13▶ V L M A A V L T V T G A V P V A R L H G A L P D A R G C H I A Q F  
701 AAGTCCCTGTCTCCACAGGAGCTGCAGGCCTTAAAGAGGGCCAAAGATGCCTTAGAAGAGTGCCTTCTGCTGAAGGACTGCAGGTGCCACTCCCGCCTC  
47▶ K S L S P Q E L Q A F K R A K D A L E E S L L L K D C R C H S R L

---

**Bsu36I (816)** **NcoI (845)**  
**BspXI (845)**  
801 TCCCCAGGACCTGGGACCTGAGGCACTGCAGGTGAGGGAGCGCCCATGGCTTTGGAGGCTGAGCTGGCCCTGACGCTGAAGTTTGGAGGCCACCGC  
80▶ F P R T W D L R Q L Q V R E R P M A L E A E L A L T L K V L E A T A  
901 TGACACTGACCCAGCCCTGGTGGACGTCTGGACCAGCCCTTACACCTGCACCATATCCTCTCCAGTTCGGGGCTGTATCCAGCCTCAGCCACG  
113▶ D T D P A L V D V L D Q P L H T L H H I L S Q F R A C I Q P Q P T

---

**Bsp120I (1003)** **XhoI (1081)**  
1001 GCAGGGCCAGGACCCGGGGCCCTCCACCATTGGCTGTACCGGCTCCAGGAGGCCAAAAAGGAGTCCCTGGCTGCCTCAGGCGCTCTGTACCT  
147▶ A G P R T R G R L H H W L Y R L Q E A P K K E S P G C L E A S V T

---

**NheI (1198)**  
1101 TCAACCTCTCCGCTCCTCACGCGAGACCTGAATTGTGTGGCAGTGGGACCTGTGTGTCTGACCCTCCACCAGTGCATGCAACCTGAGATTTTATGC  
180▶ F N L F R L L T R D L N C V A S G D L C V •

---

**MscI (1204)**  
1201 TAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATT

---

**HpaI (1336)** **MfeI (1347)**  
1301 GCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGGAGGTGGGAGGTTTTTT

---

**EcoRI (1432)**  
1401 AAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCATAAATACAGCATAGCAAACCTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGA  
1501 GGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCC

---

**SspI (1671)** **SwaI (1685)**  
1601 CAAGGTTTGAAGTACTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTATGTAATAATTCAGAAATAATTTAAATACATCAT  
1701 TGCAATGAAAATAAATGTTTTTATTAGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTGTAGTTGGACTTAGGGAACAAGGAAC  
1801 CTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGC  
141▶ • N R T Y K L P I L E E I T T K V L K G

---

**SacI (1946)** **BstXI (1975)**  
1901 CATTCACTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTC  
121▶ N M E I L V F 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  
2001 ATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGCAAAATGCTTCTGCCGTTGCTCAGCAGCAGCCCAATGGCAATGGCTTCAGCAGACAGACTGACC  
88▶ D S Y P H R V 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  
2101 CTGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCAGTCTGGTCTGATGGCCGCCCCGACATGGTGTCTGTTGCTCATAGAGCATGG  
54▶ R G I Y A E I 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

---

**BspHI (2260)**  
**BbsI (2256)** **XmnI (2252)**  
2201 TGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTTCTCATGATGGCCCTCTATAGTGAGTCTGATTATACTATGC  
21▶ I K E T A V E V L E L D Q Q S I N F T K M

---

**SacI (2375)**  
2301 CGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCCA

---

**SpeI (2473)**  
2401 CCGTACAGCCTACCGCCATTTGCGTCAATGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGAC  
2500 GTCAATGGGGTGGAGACTTGAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACT

**SnaBI (2601)**  
2600 AATACGTAGATGTA CTGCAAGTAGGAAAGTCCATAAGGTCATGTA CTGGGCATAATGCCAGGCGGGCCATTACC GTCAATTGACGTCAATAGGGGGCG

---

**NdeI (2706)**  
2700 TACTTGGCATATGATACACTTGATGTA CTGCAAGTGGGCAGTTTACC GTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTAT

---

**SdaI (2884) PacI (2892)**  
2800 GGGAAACATACGTCAATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGT CAGCCAGGCGGGCCATTACC GTAAGTTATGTAACGCCTGCAGGTTAA T

---

**BspLU11I (2902)**  
2898 TAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAC  
2998 AAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGA  
3098 CCCTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGGTG TAGGTCGT

---

**ApaLI (3216)**  
3198 TCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCCGTT CAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGAC  
3298 TTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGCCTAACTACGCTACA  
3398 CTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAG  
3498 CGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAC

---

**EagI (3652)**  
3598 GAAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGGCCGAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGG  
PacI (3632) SwaI (3641) NotI (3651)

3698 TTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCCAGTGCAAGTGCAGGTGCC  
3798 AGAACATTTCTCTATCGAA