



**PvuI (7)**  
**SgfI (6)** 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
MfeI (82) EcoNI (96)

101 GAGAAGTGGCGCGGGTAAACTGGAAAGTGTGTCGTGACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

**HindIII (245)** **Bsu36I (291)**  
**Psp1406I (203)** 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCGCCGCCGCCCTACCTGAGGCC  
PvuII (239) EcoNI (287)

301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

**NgoMIV (441)** 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

**AgeI (552)** **BspLU11I (560)** **SspI (596)**  
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTCAACATGTTGCAGATGGCTGGGCGAGTCTCCAAAATGAATA  
1▶ M L Q M A G Q C S Q N E Y

**SphI (615)** **BspLU11I (667)**  
601 TTTTGACAGTTTGTTCATGCTTGCATACCTTGTCAACTTCGATGTTCTTCTAATACTCCTCCTAAACATGTCAGCGTTATTGTAATGCAAGTGTGACC  
13▶ F D S L L H A C I P C Q L R C S S N T P P L T C Q R Y C N A S V T  
701 AATTCAGTAAAGGAACGAATGCGATTCTCTGGACCTGTTGGACTGAGCTTAATAATTTCTTGGCAGTTTCTGCTAATGTTTTGCTAAGGAAGA  
47▶ N S V K G T N A I L W T C L G L S L I I S L A V F V L M F L L R K

**XmnI (895)**  
801 TAAGCTCTGAACCATTAAAGGACGAGTTTAAAAACACAGGATCAGGTCTCCTGGGCATGGCTAACATTGACCTGGAAAAGAGCAGGACTGGTGAATGAAAT  
80▶ I S S E P L K D E F K N T G S G L L G M A N I D L E K S R T G D E I

**XhoI (916)** **StuI (912)** **BbsI (946)** **SalI (973)**  
901 TATTCTCCGAGAGGCTCGAGTACACGGTGAAGAATGCACCTGTGAAGACTCATCAAGAGCAAACCGAAGTTCGACTCTGACCATTGCTTTCCACTC  
113▶ I L P R G L E Y T V E E C T C E D C I K S K P K V D S D H C F P L

**PvuII (1067)**  
1001 CCAGCTATGGAGGAAGGCGCAACCATTCTGTACCACGAAAACGAATGACTATTGCAAGAGCCTGCCAGCTGCTTTGAGTGCTACGGAGATAGAGAAAT  
147▶ P A M E E G A T I L V T T K T N D Y C K S L P A A L S A T E I E K

**MscI (1149)** **XhoI (1130)** **NheI (1143)**  
1101 CAATTTCTGCTAGGTAATTAACCATTTGCACTCGAGCAGTGCCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGA  
180▶ S I S A R •

**HpaI (1281)** **MfeI (1292)**  
1201 ATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAAACAACAACATTGCA

**EcoRI (1377)**  
1301 TTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTAAAATACAGCATAGCA  
1401 AAACTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTG

**SapI (1559)**  
1501 CAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTGCTCTTCATTTCTTTATGTTTAAATGCACTGACCTCCAC

**Spl (1616)** **SwaI (1630)** **EcoO109I (1691)**  
1601 ATTCCCTTTTATGAAAAATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTC  
1701 ATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGT  
141▶ • N R T Y

**SacI (1891)**  
1801 ACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCT  
136▶ 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 E R

**BstXI (1920)**  
1901 GCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTG  
103▶ 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 N

**StuI (2055)**  
2001 CTCACAGCAGACCAATGGCAATGGCTTTCAGCAGACAGTACCCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGTCC  
69▶ 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 T R

**XmnI (2197)**  
2101 TGATGGCCGCCGACATGGTGTGTTGTCCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACAGCTCCAGATCCTGCTGAGAGATGTTGAA  
36▶ I A A G V H H K N D E Y L M T I K E T A V E V L E L D Q Q S I N F

**BbsI (2201)** **AseI (2263)**  
2201 GGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCT  
3▶ T K M

**SacI (2320)**  
2301 TATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCCACCGTACACGCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTT

**SpeI (2418)**  
2401 GGAAAGTCCCCTGTTGATTTACTAGTCAAAAACAACTCCCATTCAGTCAATGGGGTGGAGACTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCCAT

2501 TGATGTA CTG CCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTA CTG CCAAGTAGGAAAGTCCCATAAGGTCATGTA CTG GGCATA  
**SnaBI (2546)**

2601 ATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTA CTG CCAAGTGGGCAGTTTACCGTAAATAC  
**NdeI (2651)**

2701 TCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGGTCTGTTGGGCGGT CAGCCAG

2801 GCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTG  
**PstI (2830)** **SdaI (2829)** **BspLUIII (2847)** **PacI (2837)**

2901 CTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGG

3001 CGTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCTGTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGCGCTTTC

3101 TCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTT CAGCCCGACCGCTGCGCTTA  
**ApaLI (3161)**

3201 TCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCG

3301 GTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAG

3401 AGTTGGTAGCTCTTGATCCGGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAA

3501 GATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGG  
**EagI (3597)** **PacI (3577)** **SwaI (3586)** **NotI (3596)**

3601 CCGCAATAAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAAC

3701 AAAC TAGCAAAAATAGGCTGTCCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA