



100

**PvuI (7)**  
**SgfI (6)** MfeI (82) EcoNI (96)  
1 GGATCTGGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

**HindIII (245)** **Bsu36I (291)**  
**Psp1406I (203)** **PvuII (239)** EcoNI (287)  
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCGCCCGCCCTACCTGAGGCC  
301 GCCATCCACGCGGTTGAGTGCCTTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

**NgoMI (441)** **NaeI (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT

**NcoI (560)** **BstEII (555)**  
**KasI (535)** **AgeI (552)**  
501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGGCTACTCTGAGATCACCGGTACCATGGCGGCTCTGAGAAGTTGGTGACTCGGAGCGTCTG  
1▶ M A A L R S W V T R S V C

**NcoI (676)**  
601 CTCTCTGTTCAAGTACAGACAGCGTTTTCTGTCTTGGCTAACTCTAAGAAACGCTGTTTCTCAGAATTGATAAAACCATGGCACAAAACCTGTGTTGACT  
13▶ S L F R Y R Q R F P V L A N S K K R C F S E L I K P W H K T V L T

**DraIII (713)**  
701 GGATTTGGCATGACACTGTGTGGGTTCTATTGCTCAGAAATCGGAGCCTCAITCTCTCAGTAACGAAGCATTGATGAGGAGGGCTGTGCTTTGGTAA  
47▶ G F G M T L C A V P I A Q K S E P H S L S N E A L M R R A V S L V

**BstZ17I (861)** **Bst1107I (861)**  
801 CAGATAGCACCTCTACCTTTCTGTCTCAAACACCTACGCGCTGATTGAAGCAATCACCGAGTATACTAAGGCTGTTTACACGTTAGTGTCTCTGTACCG  
80▶ T D S T S T F L S Q T T Y A L I E A I T E Y T K A V Y T L V S L Y R

**EcoRI (927)**  
901 ACAATATAACAAGTTTACTTGGGAAGATGAATCCCAGGAGGAAGATGAGGTGTGGCAGGTAATTATAGGAGCCAGAGTTGAGATGACTTCAAACAGCAG  
113▶ Q Y T S L L G K M N S Q E E D E V W Q V I I G A R V E M T S K Q Q

**SspI (1001)**  
1001 GAATATTTGAAGTTGGAGACCCTTGGATGACAGCGGTTGGCCTTTCAGAGATGGCTGCTGAGGCTGCCTATCAAACCTGGAGCAGATCAGGCCTCCATAA  
147▶ E Y L K L E T T W M T A V G L S E M A A E A A Y Q T G A D Q A S I  
1101 CTGCCAGGAATCACATCCAGTTGGTGAAGTACAGGTACAGGAGGTGCCCGAGCTCTCCAGAAAGGCAAAACCAAGTTGGCTGAAGCACAGACAAAAGA  
180▶ T A R N H I Q L V K S Q V Q E V R Q L S Q K A E T K L A E A Q T K E

**EcoO109I (1275)** **XcmI (1298)**  
1201 GCTGCACCAGAAAGCACAGGAAGTGAGTGATGAGGGGGCTGACCAGGAAGAGGAGGCCTACCTGCGTGAAGATTGAGGGCCTGAGCCAGTGCCTGTCC  
213▶ L H Q K A Q E V S D E G A D Q E E E A Y L R E D •

**MscI (1320)**  
**DraIII (1303)** **NheI (1314)**  
1301 ATCCACTGTGGGGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTAAAAAATGCTTTATTGTGA

**HpaI (1452)** MfeI (1463)  
1401 AATTTGTGATGCTATTGCTTTATTGTAACCATTATAAGCTGCAATAACAAGTTAACAACAACAATTGCATTCATTTATGTTTCAGGTTTCAGGGGGAG

**EcoRI (1548)**  
1501 GTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTTCAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTAC  
1601 TTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTTAA

**SspI (1787)**  
1701 ATATAGTGATTTTCCCAAGGTTGAACTAGCTCTTCATTTCTTTATGTTTTAAATGCACCTGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAT

**SwaI (1801)** **EcoO109I (1862)**  
1801 AATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACT  
1901 TAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGTTACTTGAGGGGGATGAGTTCCTCAATGGTG  
141▶ • N R T Y K L P I L E E I T

**SacI (2062)** **BstXI (2091)**  
2001 GTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGACATGCCACAGGGGCTGACCACCTCGA  
126▶ 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 M G C P S V V R I  
2101 TGGATCTGTCCACCTCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCAGCAGACCAATGGCAATGGCTTC  
93▶ 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 A S G I A I A E  
2201 AGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCGAGTCTGGTCTGATGGCGCCCGACATGGTCTTGTG  
60▶ 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 A A G V H H K N

**BspHI (2376)** **XmnI (2368)**  
2301 TCCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGCTTCATGATGGCCCTCTATAGTAG  
26▶ D E Y L M T I K E T A V E V L E L D Q Q S I N F T K M ◀

