



100

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

Psp1406I (203) 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCGCTACCTGAGGCC **HindIII (245)** **Bsu36I (291)**
301 GCCATCCACGCGGTTGAGTCGCGTTTCTGCCGCCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT
NgoMIV (441)
NgoMI (441)
NaeI (441)

501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGGCTACTCTGAGATCACCGGTACCATGGGGCGGGCAGAGATGCCATCCTGGACGCTTTGA **NcoI (560)**
BstEII (555)
KasI (535) **AgeI (552)** 1▶ M G R A R D A I L D A L E
601 AAACCTGTGAGGGATGAACCTAAAAAGTTCAAGTGAAGCTGCTGACAGTGAAGTGCAGAGAGGCTATGGGCGCATCCACGCGGGGCCCTGCTGCAG **Bsp120I (686)**
13▶ N L S G D E L K K F K M K L L T V Q L R E G Y G R I P R G A L L Q
701 ATGACGCGCATAGATCTCACTGACAACTTGTGAGTACTATCTGGAGTGCATGGCTGGAGCTCACAATGACTGTGCTTAGAGACATGGGCTTACAGG **BglIII (711)**
47▶ M D A I D L T D K L V S Y Y L E S Y G L E L T M T V L R D M G L Q
801 AGCTGGCTGAGCAGCTGCAACGACTAAAGAAGAGTCTGGAGCTGTGGCAGCTGCAGCCAGTGTCCCTGCTCAGAGTACAGCCAGAAGCAGGACTTTGT **DraIII (892)**
80▶ E L A E Q L Q T T K E E S G A V A A A A S V P A Q S T A R T G H F V
901 GGACCAGCACAGGCAAGCACTCATTGCCAGGGTACAGAAAGTGGACGGAGTGTGGATGCTTTGCATGGCAGTGTGCTGACTGAAGGACAGTACCAGGCA **StuI (1091)**
113▶ D Q H R Q A L I A R V T E V D G V L D A L H G S V L T E G Q Y Q A **Eco147I (1091)**
1001 GTTCGTGCAGAGACCACCAGCCAAGACAAGATGAGGAAGCTCTTCAGCTTTGTTCCATCCTGGAACCTGACCTGCAAGGACTCCCTCCTCCAGGCTTTGA **XmnI (1035)**
147▶ V R A E T T S Q D K M R K L F S F V P S W N L T C K D S L L Q A L
MscI (1152)
BalI (1152)
1101 AGGAAATACATCCCTACTTGGTGTGAGCTGGAGCAGAGCTGAGGGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAC **NheI (1146)**
180▶ K E I H P Y L V M D L E Q S •

1201 AGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACAAAT **HpaI (1284)** **MfeI (1295)**

1301 GCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTAAAATACAGCATA **EcoRI (1380)**
1401 GCAAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCCAATGTGCATTAGCTGT
1501 TTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTACCTCTTCATTTCTTTATGTTTAAATGCACTGACCTCC

1601 CACATTCCCTTTTATGATAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATAGGCAGAATCCAGATGCTCAAGGCC **SspI (1619)** **SwaI (1633)**
1701 TTCATAATATCCCCAGTTTGTAGTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTTGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTCTGG
1801 TGTACTTGGGGGATGAGTTCCCTCAATGGTGGTTTTGACCAGCTTGCATTCATCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTC **BstXI (1923)**
137▶ 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 E
1901 TCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCC **StuI (2058)**
104▶ 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 **Eco147I (2058)**
2001 TTGCTCACAGCAGACCAATGGCAATGGCTTCCAGCAGACAGTGAACCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTTGG **XmnI (2200)**
70▶ 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 T
2101 TCCTGATGGCCGCCCCGACATGGTGTCTCATAGAGCATGGTGTGATCTCTCAGTGGCGACCTCCACAGCTCCAGATCCTGCTGAGAGATGTT **BspHI (2208)**
37▶ R 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 **VspI (2266)**
AseI (2266)
2201 GAAGTCTTCATGATGGCCCTCCTATAGTGTGATGATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCA **XmnI (2200)**
4▶ F T K M
2301 GCTTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACAT

