



100

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

Psp1406I (203) **HindIII (245)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCCTTACGCGCCCGCCGCTACCTGAGGGCC
301 GCCATCCACGCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCCGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

KasI (535) **NcoI (560)** **BstEII (555)**
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTACCATTGGCAGCTGGGCAAATGGGCACGAAGAGTGGGTGGG
1▶ M A A G Q N G H E E W V G

BsrGI (683)
601 CAGCGCATACCTGTTTGTGAGTCTCGCTGGACAAGGTGGTCTGTGCGATGCCTACGCGCACCCAGCAGAAGGTGGCAGTGTACAGGGCTCTGCAG
13▶ S A Y L F V E S S L D K V V L S D A Y A H P Q Q K V A V Y R A L Q

BsrBI (714)
701 GCTGCCTTGGCAGAGAGCGCGGAGCCGGACGTGCTGCAGATGCTGAAGATCCACCGCAGCGACCCGACGCTGATCGTGCAGCTGCGATTCTGCGGGC
47▶ A A L A E S G G S P D V L Q M L K I H R S D P Q L I V Q L R F C G

NotI (871) **DraIII (889)**
801 GGCAGCCCTGTGGCGCTTCTCCGCGCTACCGCAGGGGGCGCTGCGCGCCGCGCTGCAGAGGAGCCTGGCGGCGCGCTCGCCAGCACTCGGTGCC
80▶ R Q P C G R F L R A Y R E G A L R A A L Q R S L A A A L A Q H S V P

KasI (922) **NgoMIV (919)**
901 GCTGCAACTGGAGCTGCGCGCCGGCGCCGAGCGCTGGACGCTTTGCTGGCGGACGAGGAGCGCTGTTTGAGTTGCATCCTAGCCAGCAGCCCGACCGG
113▶ L Q L E L R A G A E R L D A L L A D E E R C L S C I L A Q Q P D R

Eco47III (959)
1001 CTCGGGATGAAGAACTGGCTGAGCTGGAGGATGCGCTGCGAAATCTGAAGTGCGGCTCGGGGCGCGGGTGGCGACGGGGAGGTGCGTTCGGCCCCCT
147▶ L R D E E L A E L E D A L R N L K C G S G A R G G D G E V A S A P

Bsp120I (1061)
1101 TGCAGCCCCGGTGCCTCTCTGTCGGAGGTGAAGCCGCCGCCGCCGCTTCCAGCTGCTGCTGCGGCGCTTCTGTCAGGGTCCAGCTGTAGTGAATCGCC
180▶ L Q P P V P S L S E V K P P P P P P A Q T F L F Q G Q P V V N R P

SandI (1295)
1201 GCTGAGCCTGAAGACCAACAGACGTTCCGCGCTCTGTGGTCTCAAATGGCGCAAGGTGGGGCGCTCACTGCAGCGAGGCTGCCGGCGCTCGGGAC
213▶ L S L K D Q Q T F A R S V G L K W R K V G R S L Q R G C R A L R D

StuI (1352)
1301 CCGGCGTGGACTCGCTGGCCTACGAGTACGAGCGGAGGACTGTACGAGCAGGCTTCCAGCTGCTGCGGCGCTTCTGTCAGGGTCCAGGGCCGCGCCG
247▶ P A L D S L A Y E Y E R E G L Y E Q A F Q L L R R F V Q A E G R R

XhoI (1426) **XcmI (1481)**
1401 CCACGCTGCAGCGCTGGTGGAGCACTCGAGGAGAACGAGCTCACCAGCCTGGCAGAGGACTTGTGGGCTGACCGATCCCAATGGCGGCTGGCCTA
280▶ A T L Q R L V E A L E E N E L T S L A E D L L G L T D P N G G L A •

MscI (1559)
1501 GACCAGGGTGCAGCCAGCTTTTGGAGAACCCTGGATGGCTTAGGGTCTTCCTCGTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAAC
313▶

HpaI (1691)
1601 CACAAC TAGAATGCAGTGA AAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAAC

MfeI (1702) **EcoRI (1787)**
1701 AACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATCTAAAATA
1801 CAGCATAGCAAACTTTAACTCCTCAAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCAT
1901 TAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTGCTCTTCATTTCTTTATGTTTAAATGCAT

SspI (2026) **SwaI (2040)**
2001 GACCTCCACATTCCTTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTC
2101 AAGGCCCTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAG
141▶

BstXI (2330)
2201 TTCCTGGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGA
139▶ N R 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

StuI (2465)
2301 TGAGCTCTGACATGCCACAGGGGCTGACCCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTT
106▶ L 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

2401 CTGCCCCTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCCTGCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCA
73▶ Q G 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

