



PvuI (7)
SgfI (6) 1 GGATCTGCGATCGTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA MfeI (82) EcoNI (96)

101 GAGAAGGTGGCGGGGTAACCTGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMI (441)
NaeI (441) 401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT

KasI (535) 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTGAGTGCAGGATGAAGATGATACATCACCTTAAATATTA SphI (560) AgeI (552) SspI (592)

601 AACTCGAAACCAGCTCTCGTCTCCGTTGGCTCTGCATCTCCTCCTGGTGGCGTGTGATGGCTTTGATTCTGCTGATCCTGTCCGTTGGGATGGTTGTC 1▶ M Q D E D G Y I T L N I K

13▶ T R K P A L V S V G S A S S S W W R V M A L I L L I L C V G M V V

Eco47III (794)
AfeI (794) 701 GGGCTGGTGGCTCTGGGATTTGGTCTGTGCATGCAGCGCAATTACCTACAAGTGTAGAATGAAAATCGCACAGGAACCTGCAACAATTAGCAAAGCGCT 47▶ G L V A L G I W S V M Q R N Y L Q D E N E N R T G T L Q Q L A K R

801 TCTGTCAATATGTGGTAAACAATCAGAACTAAAGGCCACTTTCAAAGGTCATAAATGCAGCCCTGTGACACAACTGGAGATATTAGGAGATAGCTG 80▶ F C Q Y V V K Q S E L K G T F K G H K C S P C D T N W R Y Y G D S C

ScaI (944) 901 CTATGGGTTCTTACAGCACAACCTAACATGGGAAGAGAGTAAGCAGTACTGCATGCAATGCTACTCTCCTGAAGATTGACAACCGGAACATTGTG 113▶ Y G F F R H N L T W E E S K Q Y C T D M N A T L L K I D N R N I V

1001 GAGTACATCAAAGCCAGGACTCATTAAATTCGTTGGTGGGATTTATCTGCCAGAAAGTCAATGAGGCTGGAAGTGGGAGGATGGCTCGTTATCTCAG 147▶ E Y I K A R T H L I R W V G L S R Q K S N E V W K W E D G S V I S

1101 AAAATATGTTTGAAGTTTGGAAAGTGGAAAAGGAAATATGAATTGTGCTTATTTTATAATGGGAAAATGCACCCTACCTTCTGTGAGAACAAACATTA 180▶ E N M F E F L E D G K G N M N C A Y F H N G K M H P T F C E N K H Y

MscI (1278)
NheI (1272) 1201 TTTAATGTGTGAGAGGAAGGCTGGCATGACCAAGGTGGACCAACTACCTTAATGCAAAGAGGTGGACAGGATGCTAGCTGGCCAGACATGATAAGATACA 213▶ L M C E R K A G M T K V D Q L P •

1301 TTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTG

HpaI (1410) MfeI (1421) 1401 CAATAAACAAGTTAAACAACAACATTCATTCTTTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGT

EcoRI (1506) 1501 GGTATGGAATTCTAAAATACAGCATAGCAAACCTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGG 1601 GGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTGCTTTCATTT

SspI (1745) **SwaI (1759)** 1701 TTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTGAGAAATAATTTAAATACATCATTGCAATGAAAAATAATGTTTTTTATT

EcoO109I (1820) 1801 AGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAA

1901 AGCGAGCTTCTAGCTTTAGTTTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAG 141▶ 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

SacI (2020) **BstXI (2049)** 2001 TCAGGAGCATAGTCAGAGATGAGCTCTGTCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCA 112▶ 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 D S Y P H R V A V

StuI (2184) **Eco147I (2184)** 2101 CAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTGCACCTGCCAATGTAGGCCCTCAATGTGGAC 79▶ I T D F D K 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

2201 AGCAGAGATGATCTCCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGCTTGTGCTCCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACC 46▶ 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 I K E T A V E V

XmnI (2326) **BspHI (2334)** **VspI (2392)** **AseI (2392)** 2301 AGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGATGGCCCTCTATAGTGTGATTATACTATGCCGATATACTATGCCGATGATTAATTG 12▶ L E L D Q Q S I N F T K M

SacI (2449) 2401 TCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTGCG

