



PvuI (7)
SgfI (6) 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

MfeI (82) EcoNI (96)

101 GAGAAGGTGGCGGGGTAAGTGGAAAGTGTGTCGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
PvuII (239) 201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC

Bsu36I (291) EcoNI (287)

301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441) 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

BspLU11I (560)
AgeI (552) 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCCTACCTGAGATCACCGGTCAACATGTTGCGCTGTACGTGTTGGTAATGGGAGTTTCTGC

1► M L R L Y V L V M G V S A

601 CTTACCCCTTCAGCTGCGGCACACAGGGGCTGCCAGAAGCTGCCGGTTCGTTGGGAGGCATTACAAGCGGAGTTTCCAGGCTGGAAGGGGAGCCTGTA

13► F T L Q P A A H T G A A R S C R F R G R H Y K R E F R L E G E P V

Bsu36I (702) 701 GCCCTGAGGTGCCCCAGGTGCCCTACTGGTTGTGGCCCTCTGTGAGCCCCGCATCAACCTGACATGGCATAAAAATGACTCTGCTAGGACGGTCCCAG

47► A L R C P Q V P Y W L W A S V S P R I N L T W H K N D S A R T V P

Bsp120I (821) 801 GAGAAGAAGACACGGATGTGGCCAGGACGGTCTGTGGCTTCTGCCAGCCTTGCAGGAGGACTCTGGCACCTACGTCTGCACACTAGAAATGC

80► G E E E T R M W A Q D G A L W L L P A L Q E D S G T Y V C T T R N A

901 TTCTTACTGTGACAAAATGTCCATTGAGCTCAGAGTTTTGAGAATACAGATGCTTTCCTGCCGTTTCATCTCATACCCGAAAATTTAACCTTGTCAACC

113► S Y C D K M S I E L R V F E N T D A F L P F I S Y P Q I L T L S T

EcoRI (1030) 1001 TCTGGGGTATTAGTATGCCTGACCTGAGTGAATTCACCCGTGACAAAATGACGTGAAGATTCAATGGTACAAGATTCTCTTCTTTGGATAAAGACA

147► S G V L V C P D L S E F T R D K T D V K I Q W Y K D S L L L D K D

DraIII (1144) 1101 ATGAGAAATTTCTAAGTGTGAGGGGACCCTCACTTACTCGTACAGATGTGGCCCTGGAAGATGCTGGCTATTACCGCTGTGCTGACATTTGCCCA

180► N E K F L S V R G T T H L L V H D V A L E D A G Y Y R C V L T F A H

1201 TGAAGCCAGCAATAACAATCACTAGGAGTATTGAGCTACGCATCAAGAAAAAAGAGAGACCATTCTGTGATCATTTCCCCCTCAAGACCATA

213► E G Q Q Y N I T R S I E L R I K K K K E E T I P V I I S P L K T I

1301 TCAGCTTCTGCGGTCAAGACTGACAATCCCGTGAAGGTGTTTCTGGGAACCGGCACACCCTTAACCACCATGCTGTGGTGGACGGCAATGACACCC

247► S A S L G S R L T I P C K V F L G T G T P L T T M L W W T A N D T

XmaI (1416) 1401 ACATAGAGAGCGCTACCCGGGAGGCGCGTGACCGAGGGGCCACGCCAGGAATATTAGAAAATAATGAGAATACATTGAAGTCCATTGATTTTTGA

280► H I E S A Y P G G R V T E G P R Q E Y S E N N E N Y I E V P L I F D

1501 TCCTGTACAAGAGAGGATTTGCACATGGATTTAAATGTGTTGCCATAATACCCTGAGTTTTTCAGACACTACGCACACAGTCAAGGAAGCCTCCTCC

313► P V T R E D L H M D F K C V V H N T L S F Q T L R T T V K E A S S

1601 ACGTTCTCCTGGGCATTGTGCTGGCCCACTTCACTGGCCCTTCTGGTTTTGGGGGAATATGGATGCACAGCGGTGAAAACAGAACTGGAAAAG

347► T F S W G I V L A P L S L A F L V L G G I W M H R R C K H R T G K

MscI (1775)
NheI (1769) 1701 CAGATGGTCTGACTGTGCTATGGCTCATCATCAAGACTTTCAATCCTATCCCAAGTGAATAAATGGAGCTAGCTGGCCAGACATGATAAGATACATTG

380► A D G L T V L W P H H Q D F Q S Y P K •

1801 ATGAGTTTGGACAAACCACTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAA

HpaI (1907) MfeI (1918) 1901 TAAACAAGTTAAACAACAATTGCATTCTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGT

EcoRI (2003) 2001 ATGGAACTCTAAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGC

SapI (2185) 2101 TGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTGAACCTAGCTCTTCATTCTTT

SspI (2242) SmaI (2256) 2201 ATGTTTTAAATGCACTGACCTCCACATTCCTTTTGTAGTAAAATATTAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAG

EcoO109I (2317) 2301 CAGAAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTGTAGTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTTGGACAGCAAGAAAGC

2401 GAGCTTCTAGCTTTAGTTCTGGTGTACTTGGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCA

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 D

BstXI (2546) 2501 GGAGCATAGTCAGAGATGAGCTCTGACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAA

111► 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 I

StuI (2681) 2601 TGGTGTCAAAGTCTTCTGCCCCTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGC

78► 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 A

