



PvuI (7)
SgfI (6) **MfeI (82)**
1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCCGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGTCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGACGCGTCAGCATGCGGTGGAACACTTTCTGGGCATCCTGTG

▶ ▶ M R W N T F W G I L C

601 CCTCAGCCTCCTAGCTGTTGGCACTTGCAGGACGATGCCGAGAACATTGAATACAAAGTCTCCATCTCAGGAACAGTGTAGAGTTGACGTGCCCTCTA

11▶ L S L L A V G T C Q D D A E N I E Y K V S I S G T S V E L T C P L **XbaI (696)**

701 GACAGTGACGAGAACTTAAAATGGGAAAAAATGGCCAAGAGCTGCCTCAGAAGCATGATAAGCACCTGGTGTCTCCAGGATTTCTCGAAGTCGAGGACA

45▶ D S D E N L K W E K N G Q E L P Q K H D K H L V L Q D F S E V E D **DraIII (763)**

801 GTGGCTACTACGTCGTCTACACACCAGCCTCAAATAAAAAACAGTACTTGTACTGAAAGCTCGAGTGTGTGAGTACTGTGTGGAGGTGGACCTGACAGC

78▶ S G Y Y V C Y T P A S N K N T Y L Y L K A R V C E Y C V E V D L T A **XhoI (860)** **ScaI (872)**

901 AGTAGCCATAATCATCATTGTTGACATCTGTATCACTCTGGGCTTGCTGATGGTCATTATTACTGGAGCAAGAATAGGAAGGCCAAGCCAAAGCCTGTG

111▶ V A I I I I V D I C I T L G L L M V I Y Y W S K N R K A K A K P V

1001 ACCCGAGGAACCGGTGCTGTAGCAGGCCAGAGGGCAAACAAAGGACCGGCCACCACTGTTCCCAACCCAGACTATGAGCCATCCGCAAAGGCCAGC

145▶ T R G T G A G S R P R G Q N K E R P P P V P N P D Y E P I R K G Q **BsrBI (1045)**

AgeI (1009) **MscI (1174)**
1101 GGGACCTGTATTCTGGCCTGAATCAGAGAGCAGTCTGACAGATAGGAGAGACATCGCCTTCTGTAGACCTAGCTGGCCAGACATGATAAGATAACATTGA

178▶ R D L Y S G L N Q R A V • **NheI (1168)**

1201 TGAGTTTGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAAT

HpaI (1306) **MfeI (1317)**
1301 AAACAAGTTAAACAACAACAAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGAGGTGTGGAGGTTTTTAAAGCAAGTAAAACTCTACAAATGTGGTA

EcoRI (1402)
1401 TGGAAATCTAAAAACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCT

▶ ◀

1501 GTTGCCAAATGTCATTAGCTGTTTGCAGCCTCACCTTCTTTTATGAGTAAAGATATAGTGTATTTCCCAAGTTTGAAGTACTGCTTTCATTTCTTTA

SspI (1641) **SwaI (1655)**
1601 TGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAAATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGC

EcoO109I (1716)
1701 AGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACCTTTAATAGAAATGGACAGCAAGAAAGCC

1801 AGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAGCAGTCAG

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 P

1901 GAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAAT

111▶ 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 **BstXI (1945)**

2001 GGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCTCAATGTGACAGCA

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 **StuI (2080)**

2101 GAGATGATCTCCCGAGTCTGGTCTGATGGCCGCCGACATGGTGTCTGTTGCTCATAGAGCATGGTGTCTCTCAGTGGCGACCTCCACCAGCT

44▶ 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 L E

BbsI (2226) **XmnI (2222)** **AseI (2288)**
2201 CCAGATCTGCTGAGAGATGTTGAAGGCTTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAA

11▶ L D Q Q S I N F T K M ◀

2301 AACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAA
←
2401 TGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCCTTGATTTACTAGTCAAAAACAACTCCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGT
←
2501 GAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGT
←
2601 CCCATAAGGTCATGTACTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGTACTGC
←
2701 CAAGTGGGCGAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGG
←
2801 CGGGGTCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAAGAACATGTGAGCAAAAGGCCAGCAAAGGC
←
2901 CAGGAACCGTAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAAC
←
3001 CCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTC
←
3101 TCCCTTCGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTGCTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCC
←
3201 CGTTCAGCCCAGCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGG
←
3301 ATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGC
←
3401 TGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTAC
←
3501 GCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTACGTAAAGGGATTTTGGTCATGGCT
←
3601 AGTTAATTAACATTTAAATCAGCGGCCCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCT
←
3701 CCATCAAAAACAAAACGAAAACAAAACAACTAGCAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA