



PvuI (7)
SgfI (6)
 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACCGCGCCGCCCTACCTGAGGGC
 301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

AgeI (552) **BspLU11I (560)**
 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCCTACCTGAGATCACCGGTCAACATGTGTGTAATTTTTCCGGTAGAAATCGACGTGCCCA
 1▶ M C V I F P V E I D V S Q

BstXI (635) **BsrGI (647)** **BspHI (692)**
 601 GACGATTATTCGAGATTGTCAGGTGGACAAAACAAACCAGAGAGTTGGTGTACATTAACAAGATTATGAACACGCAATTGACAAAACCGTTTCTCATGATG
 13▶ T I I R D C Q V D K Q T R E L V Y I N K I M N T Q L T K P V L M M

EcoO109I (711)
 701 TTTAACATTTCCGGTCTATACGAAGCGTTACGCGCAAGAACAACAATTTGCGCGACAGAATAAAATCAAAGTTCGATGAACAATTTGATCAACTAGAAC
 47▶ F N I S G P I R S V T R K N N N L R D R I K S K V D E Q F D Q L E
 801 GCGATTACAGCGATCAAATGGATGGATTCCACGATAGCATCAAGTATTTAAAGATGAACACTATTCGGTAAGTTGCCAAAATGGCAGCGTGTGAAAAG
 80▶ R D Y S D Q M D G F H D S I K Y F K D E H Y S V S C Q N G S V L K S

BspHI (924) **HindIII (953)** **SalI (985)**
 901 CAAGTTTGCTAAAATTTAAAGAGTCATGATTATACCGATAAAAAAGTCTATTGAAGCTTACGAGAAATACTGTTTGCCCAAATTTGGTCGACGAACGCAAC
 113▶ K F A K I L K S H D Y T D K K S I E A Y E K Y C L P K L V D E R N
 1001 GACTACTACGTGGCGGTATGCGTGTGAAGCCGGGATTTGAGAACCGCAGCAACCAAGTGTCTTTTCGAGTACAACCCGATTGGTAACAAAGTTATTG
 147▶ D Y Y V A V C V L K P G F E N G S N Q V L S F E Y N P I G N K V I
 1101 TGCCGTTTGCTCACGAAATTAACGACACGGGACTTTACGAGTACGACGTCGTAGCTTACGTGGACAGTGTGCAGTTTGTGCGAACAATTTGAAGAGTT
 180▶ V P F A H E I N D T G L Y E Y D V V A Y V D S V Q F D G E Q F E E F
 1201 TGTGCAGAGTTTAAATTTGCCGTCGTCGTTCAAAAATTCGAAAAGGTTTTATATTACAACGAAGCGTCGAAAACAAAAGCATGATCTACAAGGCTTTA
 213▶ V Q S L I L P S S F K N S E K V L Y Y N E A S K N K S M I Y K A L
 1301 GAGTTTACTACAGAATCGAGCTGGGGCAAATCCGAAAAGTATAAATGGAAAATTTTTGTAACGGTTTTATTATGATAAAAAATCAAAGTGTGTATG
 247▶ E F T T E S S W G K S E K Y N W K I F C N G F I Y D K K S K V L Y

ApaLI (1421) **MscI (1468)**
SpeI (1417) **NheI (1462)**
 1401 TTAAATTCACAATGTAAGTACTAGTCACTCAACAAAATGTAATATTAACACAATTAATAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTT
 280▶ V K L H N V T S A L N K N V I L N T I K •

1501 TGGACAAACCACAACCTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAA

HpaI (1600) **EcoRI (1696)**
 1601 GTTAACAACAACAATTGCATTATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGGGTATGGAAT
 1701 TCTAAAATACAGCATAGCAAACCTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCC

SapI (1878)
 1801 AATGTGCATTAGCTGTTGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTCCCAAGGTTGAACTAGCTCTTCAATTTCTTATGTTTT

SwaI (1949)
 1901 AAATGCACTGACCTCCACATTCCTTTTTAGTAAATATTAGAAAATATTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATC

EcoO109I (2010)
 2001 CAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTC

SacI (2210) **BstXI (2239)**
 2101 TAGCTTTAGTTCTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCAT
 -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 A Y
 2201 AGTCAGAGATGAGCTCTCTGCATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTG
 109▶ 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 T D

StuI (2374)
 2301 AAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTACGACAGACAGTACCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATG
 76▶ 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 S I
 2401 ATCTCCCAGTCTTGGTCTGATGGCCGCCGACATGGTCTTGTGCTCATAGAGCATGGTATCTTCTCAGTGGCAGCTCCACCAGCTCCAGAT
 42▶ 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 L D

BbsI (2520)
XmnI (2516) **AseI (2582)**
 2501 CCTGCTGAGAGATGTTGAAGGCTTTCATGGTGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCC
 9 Q Q S I N F T K M

2601 GTGGATGGCGTCTCCAGCTTATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTCGCTCAATGGGGC
SacI (2639)

2701 GGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCA
SpeI (2737)

2801 AACCGCTATCCACGCCCATTTGATGTAAGTCCAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCAAAGTAGGAAAGTCCCAT
SnaBI (2865)

2901 AGGTCATGTAAGTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGTAAGTCCAAAGT
NdeI (2970)

3001 GGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATATTGACGTCAATGGGCGGGG

PacI (3156)
PstI (3149)
SdaI (3148) **BspLU11I (3166)**
 3101 TC GTTGGGCGGT CAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAA
 3201 CCGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACA
 3301 GGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTT

ApaLI (3480)
 3401 CGGGAAGCGTGGCGCTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCA
 3501 GCCCGACCCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGC
 3601 AGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGC
 3701 CAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGCAAGCAGCAGATTACGCGCAG

PacI (3896)
 3801 AAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAA

EagI (3916)
SwaI (3905) **NotI (3915)**
 3901 TTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCA
 4001 AAACAAACGAAACAAACAACTAGCAAAATAGGCTGTCCCGAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA