



PvuI (7)
SgfI (6)
MfeI (82) **EcoNI (96)**

1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGTCTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACC GTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245) **Bsu36I (291)**

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTCACGCGCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

BstEII (555)

501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCACCATGGCCCTCCACCTGCCGGCTCCTCAGTGTGCACT

1 ▶ M A P P T C R L L S A A L

AgeI (552) **NcoI (560)** **NgoMIV (578)**

601 GGTCTGTGCTGCTGCTGCTGGCCACCAACCACCAGGCTACAGGGGCTGTTGTGGCCAGTGAAGTGCCTGTCAATGCCTGAAGACCCTGCCAAGGGTTGAC

13▶ V L L L L L A T N H Q A T G A V V A S E L R C Q C L K T L P R V D

BbsI (678)

EcoO109I (735) **PshAI (753)**

701 TTCAAGAACATCCAGAGCTTGAGTGTGACGCCCCAGGACCCACTGCGCCAGAGGATCATAGCCACTCTCAAGGGCGGTCAAAAAATTTGCTTTG

47▶ F K N I Q S L S V T P P G P H C A Q T E V I A T L K G G Q K V C L

NheI (883)

801 ACCTGAAGCCCCCTGGTTAGAAAATCATCAAAAGATACTGAACAAAGGCAAGGCTAACTGACCTGAAAAGGAGGAGCCTGCTAGCTGGCCAGACAT

80▶ D P E A P L V Q K I I Q K I L N K G K A N •

901 GATAAGATACATTGATGATTTGGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACC

HpaI (1021) **MfeI (1032)**

1001 ATTATAAGCTGCAATAAACAAGTTAACAACAACAAATTGCATTCAATTTATGTTTTAGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACC

EcoRI (1117)

1101 TCTACAAATGTGGTATGGAATTTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCA

SapI (1299)

1201 TAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGAGCCTCACCTCTTTTCATGAGTAAAGATATAGTGTATTTCCCAAGGTTTGAAGTAA

SspI (1356) **SwaI (1370)**

1301 CTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAA

EcoO109I (1431)

1401 TGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTG

1501 GACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCTCAATGA

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

SacI (1631) **BstXI (1660)**

1601 GCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTG

116▶ V F C 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

StuI (1795)

1701 CCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCAGACAGTACCCTGCCAATGTAGGCC

83▶ R V A V 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

1801 TCAATGTGGACAGCAGAGATGATCTCCCAAGTCTGGTCTGATGGCCGCCGACATGGTGTCTGTTGTCCTCATAGAGCATGGTGTCTTCTCAGTGG

49▶ E I H V 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

BbsI (1941)

1901 CGACCTCCACCAGCTCCAGATCTGCTGAGAGATGTTGAAGGCTTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCG

16▶ V E V L E L D Q Q S I N F T K M

XmnI (1937)

2001 ATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACATAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACC

SacI (2060)

2101 GCCATTTGCGTCAATGGGGCGGAGTTGTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGGTGGAGA

SpeI (2158)

2201 CTTGAAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACT

SnaBI (2286)

2301 **NdeI (2391)**
 GCCAAGTAGGAAAGTCCCATAAGGTCATGTA
 2401 **NdeI (2391)**
 CACTTGATGTA
 2501 **PstI (2570)** **SdaI (2569)** **PacI (2577)** **BspLU11I (2587)**
 ATTGACGTC
 2601 **PstI (2570)** **SdaI (2569)** **PacI (2577)** **BspLU11I (2587)**
 AGGCCAGCAAAGGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGT
 2701 **PstI (2570)** **SdaI (2569)** **PacI (2577)** **BspLU11I (2587)**
 CAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTCCCCCTGGAAAGCTCCCTCGTGGCTCTCTGTTCGACCTGCCGCTTACCGGAT
 2801 **PstI (2570)** **SdaI (2569)** **PacI (2577)** **BspLU11I (2587)**
 ACCTGTCCGCCTTCTCCCTTCGGAAGCGTGGCGCTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTGCTTCGCTCCAAGCTGGGCTG
 2901 **ApaLI (2901)**
 TGTGCACGAACCCCGTTCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCGGTAAGACACGACTTATCGCCACTGGCAGCA
 3001 **ApaLI (2901)**
 GCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGTACACTAGAAGAACAGTATTTG
 3101 **ApaLI (2901)**
 GTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGTTTTTTGTTTG
 3201 **ApaLI (2901)**
 CAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAACGAAAACACGTTAAGGG
 3301 **EagI (3337)** **PacI (3317)** **SwaI (3326)** **NotI (3336)**
 ATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTAATCGTA
 3401 **EagI (3337)** **PacI (3317)** **SwaI (3326)** **NotI (3336)**
 ACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGA
 3501 **EagI (3337)** **PacI (3317)** **SwaI (3326)** **NotI (3336)**
 A