



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

1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACACGCGCCCGCCCTACCTGAGGCC

301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

KasI (535)
AgeI (552) **BspLU11I (560)**

501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGGCTACCTGAGATCACCGGTCAACATGTTTCCATTGTGCTTTTTCCACCTGTTTCTTC

1▶ M F P F V L F S T L F S S

AvrII (613)

601 CATATTTACTGAACCTAGGAGAAATCGTTGGGTCTGCAACTCCTCTGACGCCACCGTTTGGTACGACTACTGTGATAACTGAAATTTCTGTTTCAATA

13▶ I F T E P R E N R W V C N S S D A T V W Y D Y C D N L K F P V S I

701 CGTCTGAGCCCTGTATAACATTGAAGGAAGCGTGAATACTCTATCTACTACATTCCAAGAAGAGATATAAGGAGCTTGATTTCAATATCTATC

47▶ R S E P C I T L K G S R G I L Y L Y Y I P R R D I R S L Y F N I Y

FspI (827)

801 TATCTACTAAGTCCATGAATTTCCATTGCGCAAGAAGTTATTTGCCGAGGATATGATGATGATTTTTCTTTTGCAGAGCTCTGAAGGGAGAGACTGT

80▶ L S T K S M N F P L R K E V I C R G Y D D D F S F C R A L K G E T V

901 GAACACAACAATAACAATTTCTCTTCAGGGGAATACGATTTTCTAAGGGACAGTACAACGTATTACAGAAGCCATCGAGGGGAACACTGAAGAAAAGCTC

113▶ N T T I Q F S F R G I R F S K G Q Y N C I T E A I E G N T E E K L

MscI (1054)

1001 TTTTGCTTGAACCTTACTATCATACTACCTGATTTCAATTAGAATGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAA

147▶ F C L N F T I I H Y P D F N •

HpaI (1186) **MfeI (1197)**

1101 CTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAAACAACAACAA

EcoRI (1282)

1201 TTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCAAAATACAGCA

1301 TAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCT

SapI (1464)

1401 GTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTCTCTTCAATTTCTTTATGTTTTAAATGCACTGACCT

SspI (1521) **SwaI (1535)**
EcoO109I

1501 CCCACATTCCTTTTTAGTAAAAATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGC

1601 CCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCT

1701 GGTGACTTGAGGGGATGAGTTCTCAATGGTGGTTTTGACCAGCTTGCATTATCTCAATGAGCACAAGCAGTCAAGGAGCATAGTCAGAGATGAGC

138▶ 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 D S I L

BstXI (1825)

1801 TCTCTGCACATGCCACAGGGGTGACCCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCC

104▶ 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 F D K Q G

StuI (1960)

1901 CGTTGCTCACAGCAGACCAATGGCAATGGCTTCCAGCAGACAGTACCCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTT

71▶ 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 I E G T K

2001 GGTCTGATGGCCGCCCCGACATGGTCTTGTCTCATAGAGATGGTGTGATCTTCTCAGTGGCGACCTCCACCTCCAGATCCCTGCTGAGAGATG

38▶ T R I A A G V H K N D E Y L M T I K E T A V E V L E L D Q S I

BbsI (2106)
XmnI (2102) **AseI (2168)**

2101 TTGAAGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGATGGCGTCTC

4▶ N F T K M

2201 CAGCTTATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTGCGTCAATGGGGCGGAGTTGTTACGAC

SpeI (2323)

2301 ATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCAGC

SnaBI (2451)

2401 CCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCAAAGTAGGAAAGTCCCATAAAGTTCATGTAAGT

NdeI (2556)

2501 GCATAATGCCAGGCGGGCCATTTACCGTCAATGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGTACTGCAAGTGGGCGAGTTTACCGTA

2601 AATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGGTCGTTGGGCGGTCA

PacI (2742)

PstI (2735)

SdaI (2734)

BspLU11I (2752)

2701 GCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGCCAGGAACCGTAAAAAGGCCG

2801 CGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATA

2901 CCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGCG

ApaLI (3066)

3001 CTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCCGACCGCTGCG

3101 CTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGT

3201 AGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGA

3301 AAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTC

PacI (3482) SmaI (3491)

3401 AAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC

EagI (3502)

NotI (3501)

3501 AGCGGCGCAATAAAATATCTTTATTTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAAC

3601 AAAACAACTAGCAAAATAGGCTGTCCCAGTGCAAGTGCAAGTGCCAGAACATTTCTCTATCGAA