



PvuI (7)
SgfI (6)
MfeI (82)

1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTATGTCGTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)

Psp1406I (203)
PvuII (239)
Bsu36I (291)

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCTGCTTCTCAACTCTACGCTTTTGTTCGTTT

BstEII (555)

AgeI (552)
NcoI (560)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTACCATGGGAGCATTTTAGACAAGCCAAAGATGGAGAAGCA

601 TAATGCCAGGGCAGGGGAATGGTTACGATACGGCCTAAGCAGCATGCAAGTTGGCGAGTTGAAATGGAGGACGCACACACGGCTGTGATCGGTTTG

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

Tth111I (714)

701 CCAAGTGGACTTGAGACATGGTCATTCTTGTGTATATGATGGGCATGCTGGTTCTCAGGTTGCCAATACTGCTGTGAGCACTTGTAGATCACATCA

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

PstI (824)

801 CCAATAACAGGATTTTCAGAGGATCTGCAGGAGCACCTTCTGTGGAGAACGTAAGAATGGAATCAGAACAGGGTTTCTGGAGATTGATGAACACATGAG

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

PvuII (944)
XhoI (999)

901 AGTTATGTCAGAGAAGAAACATGGTGCAGATAGAAGCGGGTCAACAGCTGTGGCGTCTTAATCTCTCCCAACATACTTATTTCACTAAGTGGAGAC

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

EcoRI (1082)

1001 TCGAGAGTTTACTTTGTAGGAATAGAAAAGTTCACTTCTTACACAAGACCATAAACCAAGTAACCCGCTGGAAAAAGAACGAATTGAGAATGCAGGGG

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

Bsp120I (1145)
NcoI (1176)

1101 GCTCGGTGATGATTACAGCGTGTCAATGGCTCTCTGGCTGTATCGAGGCGCCTTGGGATTTTCGATTACAAATGTGCCATGGAAAAGTCCCACAGAGCA

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

1201 GCTTGTCTCCCCAGAGCCGAAGTCCATGATATTGAAAGGCTGAAGAAGATGACCAGTTCATCATCCTTGCATGCGATGGCATCTGGGACGTCATGGGG

213▶ L V S P E P E V H D I E R S E E D D Q F I I L A C D G I W D V M G

SapI (1303)
SalI (1375)

1301 AACGAAGAGCTCTGTGACTTTGTGAGATCCAGACTTGAAGTCACTGATGACCTTGAGAAAAGTTGCAATGAAGTAGTCGACACCTGCTGTATAAGGGAA

247▶ N E E L C D F V R S R L E V T D D L E K V C N E V V D T C L Y K G

BsaBI (1419)
ApaLI (1439)

1401 GTCGAGACAACATGAGTGTGATTTTGTCTGTTTTCCAAGTGCACCCAAAGTCTCGGAGAGGCGGTGAAGAAGGAGGCGGAGCTGGACAAGTACCTGGA

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

BbrPI (1555)

Tth111I (1546)
FspI (1562)

1501 GAGCAGAGTAGAAGAAATCATAAGAAGCAGGTGGAAGGCGTCCCTGACTTAGTCCACGTGATGCGCACGTTAGCCAGTGAAGCACTCCCCAGCTCCCA

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

1601 CCAGGGGTGAATTGGCAAGCAAGCGGAATGTAATTGAAGCCGTTTACAATAGACTGAACCTTACAAAAATGACGACACTGATTCTGCGTCAACCGATP

347▶ P G G E L A S K R N V I E A V Y N R L N P Y K N D D T D S A S T D

NheI (1721)

BsrBI (1713)
MscI (1727)

1701 ATATGGTAAAGCCGCTCACGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTA

380▶ D M W •

HpaI (1859)
MfeI (1870)

1801 TTTGTAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAAACAACAATTGCATTCTTTTATGTTTCAGGTTCA

EcoRI (1955)

1901 GGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACCTTAACTCCAATCAAG

2001 CCTCTACTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGA

SapI (2137)
SspI (2194)

2101 GTTTAAGATATAGTGTATTTTCCAAGTTTGAAGTCTCTCATTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATGAAAAATTT

Swal (2208)

2201 CAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGTTTAGTAG

2301 TTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAGCGAGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTC
141 • N R T Y K L P I L E E
BstXI (2498)

2401 AATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGTGACC
129 I T T K V L K G N M E I L V F C D P A Y D S I L E R C M G C P S V

2501 ACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCCTTCTGCCCGTTGCTCACAGCAGACCCAATGGCAA
95 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 N S V A S G I A I

StuI (2633)

2601 TGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCCTGATGGCCGCCCGACATGGTG
62 A E A C V T V R G I Y A E I H V A S I I E G T K T R I A A G V H H
BbsI (2779)

2701 CTTGTTGCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCCTA
29 K N D E Y L M T I K E T A V E V L E L D Q Q S I N F T K M
XmnI (2775)

AseI (2841)

2801 TAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCCTAAACGA

SpeI (2996)

2901 GCTCTGTTATATAGACCTCCACCGTACACGCTACCGCCATTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTA
3001 GICAAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCAT
SnaBI (3124)

3101 CATCATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCG

NdeI (3229)

3201 TCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCCTAAATACTCCACCCATTGACGTCAATGGA
3301 AAGTCCCTATTGGCGTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTGCTTGGGCGGTACGCCAGGCGGGCCATTTACCGTAAGTTA

PacI (3415)

PstI (3408)

SdaI (3407)

BspLU11I (3425)

3401 TGTAACGCTGCAGGTTAATAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCCTTTTCCATAGGCTCC
3501 GCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCTTTCCCTGGAAGCTCCCT
3601 CGTGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCATAGCTCAGCTGTAGGTAT

ApaLI (3739)

3701 CTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTACGCCGACCCTGCGCTTATCCGGTAACTATCGTCTTGAGT
3801 CCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGTACAGAGTCTTGAAGTG
3901 GTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGC
4001 AAACAACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGG

EagI (4175)

PacI (4155) SwaI (4164) NotI (4174)

4101 GGTCTGACGCTCAGTGGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTT
4201 TCATTACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCC
4301 CCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA