



PvuI (7) SgfI (6) MfeI (82) EcoNI (96)
 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGACTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) HindIII (245) Bsu36I (291)
 PvuII (239) EcoNI (287)
 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCGCCCGCCCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441) NgoMI (441) NaeI (441)
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCCTGCTTGCCTCAACTCTACGCTTTTGTTCGTTT

KasI (535) NcoI (560) BstEII (555) AgeI (552) XcmI (566)
 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCCCTACCTGAGATCACCGGTACCATGGCCACTGGCCACGGAGTATGTGTTCTC
 1 M A I A H L A T E Y V F S

BspLUII (676)
 601 GGACTTCTTGCTGAAGGAGCCACCCAGCCCAAGTTCAAGGGGCTGCGACTGGAGCTGGCGGTGGACAAGATGGTCACATGTATTGCCGTGGGTCTACCT
 13 D F L L K E P T E P K F K G L R L E L A V D K M V T C I A V G L P

BsiBI (732) BsaBI (732) Asp718I (742) BglII (731) Acc65I (742)
 701 CTGCTGCTCATCTCGCTGGCCTTCGCTCAGGAGATCTCCATCGGTACCCAGATAAAGTCTTCTCCCCGAGTTCTTCTCCTGGCAGAGGCTGCCTTTG
 47 L L L I S L A F A Q E I S I G T Q I S C F S P S S F S W R Q A A F

BsrGI (822)
 801 TGGATTCACTACTGCTGGGCTGCTGTACAGCAGAAGAGCTCCCTGCAGAGCGAGTCTGAAACCTCCCACTGTGGCTGCACAAGTCTTCCCTACATCCT
 80 V D S Y C W A A V Q Q K S S L Q S E S G N L P L W L H K F F P Y I L
 901 ACTGCTGTTTGCATACCTCTGTACCTGCCGCACTCTTCTGGCGCTTCTGCGAGCTCCACACCTCTGCTCAGACCTGAAAGTTATCATGGAGAACTT
 113 L L F A I L L Y L P A L F W R F S A A P H L C S D L K F I M E E L

Tth111I (1000) XhoI (1040)
 1001 GACAAAGTCTACAACCGCCATCAAGGCTGCCAAGAGTGTCTGAGATTTGGACCTAAGAGACGGACCTGGACCCCGAGGAGTACTGAGAATGTGGGGC
 147 D K V Y N R A I K A A K S A R D L D L R D G P G P P G V T E N V G

EcoRV (1113) Eco32I (1113) ScaI (1151) BspHI (1187)
 1101 AGAGTCTGTGGGAGATATCTGAAAGCCACTTCAAGTACCCAATCGTGGAGCAGTACTTGAAGACAAAAAAGAACTCTAGTCATTTAATCATGAAATACAT
 180 Q S L W E I S E S H F K Y P I V E Q Y L K T K K N S S H L I M K Y I

NgoMIV (1205) NgoMI (1205) NaeI (1205)
 1201 TAGCTGCCGGCTGGTGACATTTGTGGTTACTGTTGGCATGTATCTACTTGAGCTATTACTTACGCTCTCTTCACTCTCGGACGAGTTTCTGTGCAGC
 213 S C R L V T F V V I L L A C I Y L S Y Y F S L S S L S D E F L C S

PvuI (1338) PvuII (1378)
 1301 ATCAAATCAGGCGTCTGAAAAATGACAGCACCATCCCGATCGCTTCCAGTGAAGCTCATCGCCGTGGGCATCTTCCAGCTGCTCAGCCTCATTAACC
 247 I K S G V L K N D S T I P D R F Q C K L I A V G I F Q L L S L I N
 1401 TCATTGTGTATGCTCTGTGATTTCCCGTGGTCTACAGTCTTCCATCCCATCCCGCAGAAAAACGGACATTTCAAAGTGTATGAAATCTGCCAC
 280 L I V Y A L L I P V V V Y T F F I P F R Q K T D I L K V Y E I L P T
 1501 CTTGATGTTCTACATTTCAAGTCTGAAGGCTACAATGACTTGAGCCTCTACAACCTTTTCTGGAAGAGAACATAAGTGAGCTCAAATCGTACAAGTGT
 313 F D V L H F K S E G Y N D L S L Y N L F L E E N I S E L K S Y K C

BsiBI (1680) BsaBI (1680)
 1601 CTGAAGGTGCTGGAGAACATTAAGAATGGGAGGGCATTGACCCCATGCTACTCCTGACAAAACCTGGGCATGATTAAGATGGACATCATTGATGGAA
 347 L K V L E N I K S N G Q G I D P M L L L T N L G M I K M D I I D G
 1701 AAATTTCCACGTCCCTACAGACCAAGGGAGAGGACAGGGCAGCCAGAGAGTGGAGTTCAAAGATTTGGACCTGAGCAGCGAGGCTGCAGCAACAATGG
 380 K I P T S L Q T K G E D Q G S Q R V E F K D L D L S S E A A A N N G

NheI (1859)
 1801 GGAGAAGAACTCTGCCAGAGGCTTCTGAATCCGTCTGTAATGGTTTCTTCTTGAAGCTAGTGGCCAGACATGATAAGATACATTGATGAGTTTGG
 413 E K N S R Q R L L N P S C

HpaI (1997)
 1901 ACAAACCACAAGTGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAACAAGTT

MfeI (2008) EcoRI (2093)
 2001 ACAAACAACAATTGCATTCATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTC
 2101 AAAATACAGCATAGCAAACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAAT
 2201 GTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTCTTCATTTCTTTATGTTTTAAA

SspI (2332) SmaI (2346)
 2301 TGCACTGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAAATCCAG

EcoO109I (2407)
 2401 ATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAG

2501 CTTTAGTTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGT
 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 D

BstXI (2636)
 2601 CAGAGATGAGCTCTGTCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGCAAA
 108 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 F

StuI (2771)
Eco147I (2771)
 2701 GTCCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATC
 75 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 I

2801 TCCCCAGTCTGGTCTGATGGCCGCCGACATGGTCTTGTGTCCCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCT
 41 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 Q

BspHI (2921) **VspI (2979)**
XmnI (2913) **AseI (2979)**
 2901 GCTGAGAGATGTTGAAGGCTTTCATGATGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAAACAGCGTG
 8 Q S I N F T K M

3001 GATGGCGTCTCCAGCTTATCTGACGGTTCCTAAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCTACCGCCATTTGCGTCAATGGGGCGGA

SpeI (3134)
 3101 GTTGTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAA

SnaBI (3262)
Eco105I (3262)
 3200 CCGCTATCCACGCCATTGATGTACTGCCAAAACGCATCATCATGGTAATAGCGATGACTAATCGTAGATGTACTGCCAAGTAGGAAAGTCCCATAG

NdeI (3367)
 3300 GTCATGTACTGGGCATAATGCCAGGCGGGCATTACCCTGATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGTACTGCCAAGTGGG
 3400 CAGTTTACCGTAAATACTCCACCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGGGCGGGGTC

SdaI (3545) **PacI (3553)** **BspLU11I (3563)**
 3500 GTTGGGCGGTCAGCCAGGCGGGCATTACCCTAAGTTATGTAACGCCCTG C A G G T T A A T T A A G A A C A T G T G A G C A A A A G G C C A G A A A A G G C C A G G A A
 3598 CCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAATAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACA
 3698 GGACTATAAAGATACCAGGCGTTTCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTT

ApaLI (3877)
 3798 CGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCCGTTCA
 3898 GCCCGACCCTGCGCCTTATCCGTAACATATCGTCTTGTAGTCCAAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGC
 3998 AGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGC
 4098 CAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAG

PacI (4293)
 4198 AAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCAATGGCTAGTTAA

EagI (4313)
NotI (4312)
SwaI (4302)
 4298 TTAACATTTAAATC AGCGGCGCAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCA
 4398 AAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCCAGTGCAAGTGCAAGTGCAGGTCAGAACATTTCTCTATCGAA