



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

1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCAGTTCGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

KasI (535)
NcoI (560)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTACCATGGGAAGGTGCTATCCAAGATCTTCGGGAACAAGGA
 1 M G K V L S K I F G N K E

BamHI (605)
NgoMIV (631)
BsrGI (650)
BstEII (675)
XcmI (683)

601 AATGCGGATCCTCATGCTGGCCTGGACGCAGCCGCAAGACAACGATCCTGTACAAGTTGAAGCTGGCCAATCGGTGACCACCATCCCCACGGTGGGC
 13 M R I L M L G L D A A G K T T I L Y K L K L G Q S V T T I P T V G

BsrBI (775)
EagI (772)
SmaI (799)

701 TTCAACGTGGAGACGGTGACTTACAAAAACGTCAGTTCAACGTGTTGGATGTGGGCGCCAGGACAAGATCCGGCCGCTCTGGCGGCATTACTACACCG
 47 F N V E T V T Y K N V K F N V W D V G G Q D K I R P L W R H Y Y T
 801 GGACCCAGGTCTGATCTTCGTGGTAGACTGCGCCGACCGCAGCCGATCGACGAGGCCCGCCAGGAGCTGCACCGCATTATCAATGACCGGGAGATGAG
 80 G T Q G L I F V V D C A D R D R I D E A R Q E L H R I I N D R E M R

BstAPI (930)

901 GGACGCCATCATCCTCATCTTCGCCAACAGCAGGACCTGCCGATGCCATGAAACCCATGAGATCCAGGAGAACTGGGCCTGACCCGGATTCCGGAC
 113 D A I I L I F A N K Q D L P D A M K P H E I Q E K L G L T R I R D

HpaI (1065)

1001 AGAACTGGTATGTGCAGCCCTCCTGTGCCACCTCCGGGACGGACTCTATGAGGGGCTCACATGGTTAACCTTAACCTAACAATCCTAATGAGCGTCT
 147 R N W Y V Q P S C A T S G D G L Y E G L T W L T S N Y K S •

MscI (1111)
NheI (1105)

1101 CCACCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGA

HpaI (1243)
MfeI (1254)

1201 TGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACAATTGCATTATTTATGTTTCAGGTTTCAGGGGAGGTGTGGAG

EcoRI (1339)

1301 GTTTTTTAAAGCAAGTAAACCTCTACAAATGGTATGGAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCT
 1401 TTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGT

SspI (1578)
SwaI (1592)

1501 ATTTTCCAAGGTTTGAAGTCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATAGTAAATATTCAGAAATAATTTAAAT
 1601 ACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTGTAGTGTGGACTTAGGGAACA
 1701 AAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACC
 141 • N R T Y K L P I L E E I T T K V

SacI (1853)
BstXI (1882)

1801 AGCTTGCCATTCATCTCAATGAGCACAAGCAGTCCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCTGTGATGATCTGT
 123 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 R I S R D
 1901 CCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTCTGCCCCTGCTCACAGCAGACCCAATGGCAATGGCTTCCAGCAGAC
 90 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 A E A C V

StuI (2017)

2001 AGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCGAGTCTTGGTCTGTGATGGCCGCCCGACATGGTGTGTTGTCCTCATAG
 57 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 K N D E Y

BbsI (2163)
XmnI (2159)

2101 AGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGCTTCTCATGGTGGCCCTCTATAGTGAGTCGTATTAT
 23 L M T I K E T A V E V L E L D Q Q S I N F T K M

AseI (2225)
SacI (2282)

2201 ACTATGCCGATATACTATGCCGATGATTAATTGCAAAACAGCGTGGATGGCGTCTCCAGCTIATCTGACGGTCACTAAACGAGCTCTGCTTATATAGA

SpeI (2380)

2301 CCTCCCACGTACACGCCTACCGCCATTTGCGTCAATGGGGCGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCC
 2401 ATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCC

