



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

1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245) **EcoNI (287)**

201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACCGCGCCCGCCGCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCAGTTCGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560)

KasI (535)
BstEII (555) **AgeI (552)**

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTACCATGGCAGGTGGGGATGGAATGCTACGCCGTCTGCTGAG
1▶ M A G G D G M L R R L L R

MscI (623)
SphI (651) **DraIII (667)**

601 GCTGCACCGCACCAGATCGCGGTGGCCATAGACAGTGCCTTTCCGCTGCTGCATGCTCTAGCCGACCACGACGTGGTCCCTGAGGACAAGTTCCAGGAG
 13▶ L H R T E I A V A I D S A F P L L H A L A D H D V V P E D K F Q E

StuI (734)
XmaI (767)

701 ACGCTCCGCTGAAGGAGAAGGAAGGCTGCCCCAGGCCTTCCACGCCCTGCTGCTGGCTCTGACCCGGGACAGTGGGGCCATCCTGGATTTCTGGA
 47▶ T L R L K E K E G C P Q A F H A L L S W L L T R D S G A I L D F W

BsrBI (826)

801 GGATTCTCTTAAAGACTACAATCTGGAGCGGTACAGCCGCTGCATAGCATCCTGGACGGCTTCCCAAAGATGTGGACCTAAACCAGTCCCGGAAAGG
 80▶ R I L F K D Y N L E R Y S R L H S I L D G F P K D V D L N Q S R K G

SacII (926)
XhoI (977)

901 GAGAAAGCCCTTGTGGTCCCAAGGCCGCGGTACTGCCACCCAGACCCCAAGAGAAAAGCACTGGAGGAGCCTCGAGCCACCCACCAAGCAACT
 113▶ R K P L A G P K A A V L P P R P P T K R K A L E E P R A T P P A T
 1001 CTGGCCTCAAAGAGCGTCTCCAGCCAGGCTCCACCTGAAGACTAAGCCCTAAGAAGCCAGATGGCAACTGGAGTCACAGCATTCTCTTTGGAA
 147▶ L A S K S V S S P G S H L K T K P P K K P D G N L E S Q H L P L G

BsaBI (1188)

EcoRI (1103) **Neol (1112)**
BamHI (1187)

1101 ACGGAATTCAGACCATGGCAGCTTCTGTCCAGAGAGCTGTGACCGTGGCCTCTGGGGATGTTCCAGGAACCCGAGGGGCGTGGAAAGGATCCTTATCCA
 180▶ N G I Q T M A A S V Q R A V T V A S G D V P G T R G A V E G I L I Q

BstBI (1270)

1201 GCAGGTGTTTGTAGTCAGGAAGATCCAAGAAGTGCAATCAGGTTGGGGAGAGTTTTATACACCAACAAGTTCGAAGACCCAGTGGCAATTTGAAGAAC
 213▶ Q V F E S G R S K K C I Q V G G E F Y T P N K F E D P S G N L K N

BstXI (1389)

1301 AAGCCCCGGAGTGGTAGCAGCCTAAAGCCAGTGGTCCGAGCCAAGGGAGCCAGGTCACTATACCTGGTAGAGATGAGCAGAAAAGTGGGCCAGCAGTGTG
 247▶ K A R S G S S L K P V V R A K G A Q V T I P G R D E Q K V G Q Q C

HpaI (1438)
PshAI (1498)

1401 GGGTTCCTCCCTTCCATCCCTCCCAAGTGGAGCCAGGTTAACCAGACAGGATGAGTGTGCCGTGTGCCACGACGGAGGTGAGCTCATCTGTTGTGA
 280▶ G V P P L P S L P S E P Q V N Q N E D E C A V C H D G G E L I C C D

XmaI (1508)

1501 CGGCTGTCCCGGCGCTTCCACCTGGCTTGCCTGTCCCACTCTGCAGGAGATCCCAAGTGGCTCTGGAGATGCTCCTGCTGCCTCCAGGGCAGAGTC
 313▶ G C P R A F H L A C L S P P L Q E I P S G L W R C S C C L Q G R V

PshAI (1670)
Bsp120I (1698)

1601 CAACAGAACCTGTCCAGCCTGAGGTGTCCAGGCCCGGAGCTACCTGCAGAGACCCGATCCTCGTGGGACTGAGGTTCAGGTTTCAGAGAAAACCGGG
 347▶ Q Q N L S Q P E V S R P P E L P A E T P I L V G L R S A S E K T R

NdeI (1742)

1701 GCCATCCAGGGAGCTCAAAGCCAGCTCTGATGCTGTGCACATATGTGAACCTGCTGGCCCCGACCCCTGCAGTCCCTGCTGGAGCCTTCAGCACT
 380▶ G P S R E L K A S S D A A V T Y V N L L A P H P A A P L L E P S A L

DraIII (1898)

BstAPI (1897)

BssHII (1850)
ApaLI (1895)

1801 GTGCCCTCTACTGAGTGTGGGAATGAGGGGCGCCAGGTCCAGCACCAAGCGCGCATGTCAGTGTGTGGCGATGGCACCGAGGTGTTGGCGTGTGCA
 413▶ C P L L S A G N E G R P G P A P S A R C S V C G D G T E V L R C A

BstAPI (1973)

1901 CACTGTGCCGCTGCTTCCACTGGCGCTGCCACTTCCCGACGGCCGCCCGGGGACCAATCTCCGCTGCAAATCCTGCTGCTGAGACTCGACTC
 447▶ H C A A A F H W R C H F P T A A A R P G T N L R C K S C S A D S T

Bsp120I (2038)

2001 CCACGCCAGGCACACCGGGCGAAGCTGTACCCACCTCTGGGCCCGTCCAGCACCTGGGCTTGGCAAGGTAGGGGACGACTCTGCTAGTCACGACCCTGT
 480▶ P T P G T P G E A V P T S G P R P A P G L A K V G D D S A S H D P V

2101 TCTACATAGGGACGACCTGGAGTCCCTCCTCAATGAGCACTCATTGACGGCATCCTGCAGTGGGCCATCCAGAGCATGTCACGCCCGCTGGCCGAGACA
 513▶ L H R D D L E S L N E H S F D G I L Q W A I Q S M S R P L A E T

MscI (2224)

NheI (2218)

2201 CCACCTTCTCTTCTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTT
 547▶ P P F S S •

