



PvuI (7)
SgfI (6) **MfeI (82)**
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **HindIII (245)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTTCGAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

BstEII (555)
AgeI (552) **NcoI (560)**
501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTCAACCTGGCTCTCTTCAGTGCCAGTCTCCATACATTAACC
13▶ I I P F T G P I Q G G L Q E G L Q V T L Q G T T K S F A Q R F V V

PstI (634) **BstEII (651)**
601 GATCATCCCCTTTACTGGACCAATCCAAGGAGGGCTGCAGGAGGGACTTCAGTGACCCTCCAGGGGACTACCAAGAGTTTGCACAAAGGTTTGTGGTG
13▶ I I P F T G P I Q G G L Q E G L Q V T L Q G T T K S F A Q R F V V

XmnI (708)
701 AACTTTCAGAACGCTTCAATGAAATGACATTGCCTTCCACTTCAACCCCCGGTTTGAGGAAGGAGGTATGTGGTTTGAACACGAAGCAGAACGGAC
47▶ N F Q N S F N G N D I A F H F N P R F E E G G Y V V C N T K Q N G

Bsu36I (808)
EcoO109I (804)
801 AGTGGGGTCTGAGGAGAGAAAGATGCAGATGCCCTTCCAGAAGGGGATGCCCTTTGAGCTTTGCTTCTCGTGCAGAGGTCAGAGTTCAAGTGATGGT
80▶ Q W G P E E R K M Q M P F Q K G M P F E L C F L V Q R S E F K V M V

DraIII (943)
901 GAACAAGAAATCTTTGTGCAGTACCAACACCGGTACCTACCTCGTGGACACCATCGTGTCTCCGGTCTTGAAGCTGTCTTTATCACCTTC
113▶ N K K F F V Q Y Q H R V P Y H L V D T I A V S G C L K L S F I T F

NcoI (1034) **BstXI (1049)**
1001 CAGACTCAGAATTTCTGCTGCCACCAGGCACCCATGGCTCAAACCTACCATCCATATGGTTCACAGCACCCCTGGACAGATGTTCTTACTCTGGAA
147▶ Q T Q N F R P A H Q A P M A Q T T I H M V H S T P G Q M F S T P G

BspHI (1184)
1101 TCCCTCTGTGGTGTACCCACCCAGCCTATACCATACCTTTCTACACCCCATCCAAATGGGCTTACCCTGCAAGTCCATCATGATCAGGCCAA
180▶ I P P V V Y P T P A Y T I P F Y T P I P N G L Y P S K S I M I S G N

1201 TGTCTTGCAGATGCTACGAGTTCCATATCAACCTTCCGCTGTGGAGGTGACATTGCTTCCACCTGAACCCCGTTTCAATGAGAATGCTGTTGCCGA
213▶ V L P D A T R F H I N L R C G G D I A F H L N P R F N E N A V V R

SapI (1330)
1301 AACACTCAGATCAAACTCCTGGGGGAGGAAAGAGCGAAGTCTGCTTGGGAGGATGCCCTTCAGTCGAGGCCAGAGCTTCTCGGTGGATCATATGTG
247▶ N T Q I N N S W G Q E E R S L L G R M P F S R G Q S F S V W I I C

BspLU11I (1436) **XbaI (1485)**
1401 AAGGTCACTGCTTCAAGGTAGCTGTGAATGGTCAACACATGTGTGAATATTACCACCGCTGAAGAACTTGCAGGATATCAACACTCTAGAAGTGGCGGG
280▶ E G H C F K V A V N G Q H M C E Y Y H R L K N L Q D I N T L E V A G

BbrPI (1516) **AvrII (1544)** **NheI (1557)**
1501 TGATATCCAGCTGACCCACGTGCAGACATAGGCAAGTCTCTGGCCTAGGGATAAGGGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGAC
313▶ D I Q L T H V Q T •

HpaI (1695)
1601 AAACCACAAGTGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAA

MfeI (1706) **EcoRI (1791)**
1701 CAACAACAATTGCATTATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAA
1801 AATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCCAATGT

SapI (1973)
1901 GCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAACAGCTCTTCATTCTTTATGTTTAAATG

SwaI (2044)
2001 CACTGACCTCCACATTCCTTTTATGTAATAATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGAT

EcoO109I (2105)
2101 GCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTCTAGCT

2201 TTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAGCAGTCAGGAGCATAGTCA
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

