



PvuI (7)
SgfI (6) MfeI (82) EcoNI (96)
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245) **Bsu36I (291)**
Psp1406I (203) PvuII (239) EcoNI (287)
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACAGCGCCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTGGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

KasI (535) **AgeI (552)** **BspHI (560)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTCATCATGACACCACTGAACGTCCTTCTCCCAAGGGTGTG
1▶ M T P P E R L F L P R V C

XcmI (657) **Bsp120I (653)** PvuII (689)
EcoO109I (652)
601 TGGCACCACCTACACCTCCTCCTTCTGGGGCTGCTGCTGGTTCTGCTGCCTGGGGCCAGGGGCTCCCTGGTGTGGCCTCACACCTTCAGCTGCCAG
13▶ G T T L H L L L L G L L L V L L P G A Q G L P G V G L T P S A A Q

NsiI (721)
701 ACTGCCCCTCAGCACCCCAAGATGCATCTTCCACAGCACCTCAAACCTGCTGCTCACCTCATTGGAGACCCAGCAAGCAGAAGTCACTGCTCTGGA
47▶ T A R Q H P K M H L A H S T L K P A A H L I G D P S K Q N S L L W

RsrII (809)
801 GAGCAAAACACGGACCGTCCCTCCAGGATGGTTTCTCCTTGAGCAACAATTCTCCTGGTCCCCACAGTGGCATCTACTTCTGCTACTCCAGGT
80▶ R A N T D R A F L Q D G F S L S N N S L L V P T S G I Y F V Y S Q V
901 GGTCTTCTGGGAAAGCCTACTCTCCAAAGGCCACTCTCCCACTCTACCTGGCCATGAGGTCCAGCTTCTCCTCCAGTACCCCTTCCATGTG
113▶ V F S G K A Y S P K A T S S P L Y L A H E V Q L F S S Q Y P F H V

PstI (1036)
1001 CCTCCTCAGTCCAGAAAGTGGTGTATCCAGGGCTGCAGGAACCTGGCTGCACTCGATGTACCACGGGGCTGCGTTCAGCTCACCCAGGGAGACC
147▶ P L L S S Q K M V Y P G L Q E P W L H S M Y H G A A F Q L T Q G D

BstXI (1110) **ScaI (1147)** **NheI (1184)** **MscI (1190)**
1101 AGCTATCCACCCACAGATGGCATCCCCACCTAGTCCCTCAGCCCTAGTACTGTCTTTGGAGCCTTCGCTCTGTAGAAGTCTAGCTGGCCAGACA
180▶ Q L S T H T D G I P H L V L S P S T V F F G A F A L •

1201 TGATAAGATACATTGATGAGTTTGGACAAACCACAAGTGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAAC

HpaI (1322) MfeI (1333)
1301 CATTATAAGCTGCAATAAACAAGTTAAACAACAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAC

EcoRI (1418)
1401 CTCTACAAATGTGGTATGGAAATCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGC
1501 ATAGGCATCAGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAAC

SspI (1657) **SwaI (1671)**
1601 GCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAAATATTAGAAATAATTTAAATACATCATTGCAATGAAAATAA

EcoO109I (1732)
1701 ATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATT
1801 GGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCTCAATG
141▶ • N R T Y K L P I L E E I T T K V L K G N M E I

SacI (1932) **BstXI (1961)**
1901 AGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCACCTCATCAGAGTAGGGGT
116▶ L V F C D P A Y D 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

StuI (2096)
2001 GCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTGTCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCAATGATGAGC
83▶ 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 T V R G I Y A
2101 CTCAATGTGGACAGCAGAGATGATCTCCAGCTTGGTCTGATGGCCGCCGACATGGTCTGTTGCTCATAGAGCATGGTATCTTCTCAGTG
50▶ 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 L M T I K E T

XmnI (2238)
2201 GCGACCTCCACAGCTCCAGATCCTGCTGAGAGATGTTGAAGGCTTTCATGATGGCCCTCTATAGTGTGATTATACTATGCGGATATACTATGCC
16▶ A V E V L E L D Q Q S I N F T K M

AseI (2304) **SacI (2361)**
2301 GATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTAC

SpeI (2459)
2401 CGCCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAACAAACTCCCATGACGTCAATGGGGTGGAG

