



PvuI (7) SgfI (6) MfeI (82)
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
101 GAGAAAGTGGCGGGGTAACGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
201 Psp1406I (203) PvuII (239) HindIII (245) Bsu36I (291)
GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACACGGCCCGCCCTACCTGAGGGC
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTCCTGACCCGCTGCTTGTCAACTCTACGCTTTTGTTCGTTT

KasI (535) AgeI (552) SphI (567)
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTAG AGGGCCAGCATGCGCCGGCCAGCCGAGACTACGGCAAGT
1 M R R A S R D Y G K

FspI (603) KasI (685)
601 ACCTGCGCAGTCCGAGGAGATGGGCGAGCGGCCCGGGCTCCACACGAGGGTCCGCTGCACCCCGCGCTTCTGCACCGGCTCCGGCGCCGCCACCCGC
11 Y L R S S E E M G S G P G V P H E G P L H P A P S A P A P A P P P A

BstXI (711)
701 CGCCTCCCGCTCCATGTTCTGGCCCTCCTGGGCTGGGACTGGGCGAGGTGGTCTGCAGCATCGCTCTGTTCTGTACTTTTCGAGCGCAGATGGATCCT
44 A S R S M F L A L L G L G L G Q V V C S I A L L F L Y F R A Q M D C P
801 AACAGAATATCAGAAGACAGCACTACTGCTTTTATAGAACTCCATGAGACTCCAGAAAACGCAGATTGTCAGGACTCGACTCTGGAGAGTGAAGACAC
78 N R I S E D S T H C F Y R I L R L H E N A D L Q D S T L E S E D T

SdaI (909) DraIII (962) Eco47III (978)
901 TACCTGACTCCTGCGAGGAGATGAAACAAGCCTTTAGGGGGCGTGCAGAAGGAACTGCAACACATTGTGGGGCCACAGCGCTTCTCAGGAGCTCCAGC
111 L P D S C R R M K Q A F Q G A V Q K E L Q H I V G P Q R F S G A P A

Bsu36I (1043)
1001 TATGATGGAAGGCTCATGGTTGGATGTGGCCAGCGAGGAAGCTGAGGCCAGCCATTTGCACACCTCACCATCAATGCTGCCAGCATCCCATCGGGT
144 M M E G S W L D V A Q R G K P E A Q P F A H L T I N A A S I P S G

PvuI (1132) Acc65I (1125) BglIII (1149) HpaI (1187)
1101 TCCATAAAGTCACTCTGTCTCTTGGTACCACGATCGAGGCTGGGCCAAGATCTCTAACATGACGTTAAGCAATGGAAAATAAGGGTTAACCAAGATG
178 S H K V T L S S W Y H D R G W A K I S N M T L S N G K L R V N Q D

BspHI (1237) PvuII (1274)
1201 GCTTCTATTACCTGTACGCCAACATTTGCTTTCCGTCATGAAACATCGGGAAGCGTACCTACAGACTATCTTCAGCTGATGGTGTATGTCGTTAAAC
211 G F Y Y L Y A N I C F R H E T S G S V P T D Y L Q L M V Y V V K T

EcoRI (1370)
1301 CAGCATCAAAATCCCAAGTTCTCATAACCTGATGAAAGGAGGGAGCAGAAAACTGGTCGGGCAATTCTGAATTCACCTTTTATTCCATAAATGTTGGG
244 S I K I P S S H N L M K G G S T K N W S G N S E F H F Y S I N V G

BspEI (1465) BsaBI (1463) BamHI (1462)
1401 GGATTTTTCAAGCTCCGAGCTGGTGAAGAAATTAGCATTAGGTGTCCAACCTTCCCTGCTGGATCCGGATCAAGATGCGACGTACTTTGGGGCTTTC
278 G F F K L R A G E E I S I Q V S N P S L L D P D Q D A T Y F G A F

MscI (1559) NheI (1553)
1501 AAGTTCAGGACATAGACTGAGACTCATTTCGTGGAACATTAGCATGGATGTCCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAC
311 K V Q D I D •

HpaI (1691)
1601 CACAAC TAGAATGCAGTGA AAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAACAAAC

MfeI (1702) EcoRI (1787)
1701 AACAAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAAATTCATAAATA
1801 CAGCATAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCAT
1901 TAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTCTTCATTTCTTTATGTTTAAATGCAT

SspI (2026) SmaI (2040)
2001 GACCTCCACATTCCTTTTTAGTAAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTC

EcoO109I (2101)
2101 AAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAG
141 •

BstXI (2330)
2201 TTCCTGGTGTACTTGAGGGGATGAGTTCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGA
139 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 S I
2301 TGAGCTCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTT
106 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 D K

