



PvuI (7)
SgfI (6) 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
MfeI (82) EcoNI (96)

101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCAGAGGGCTCGCATCTCTCTTACCGCGCCCGCCCTACCTGAGGCC
PvuII (239) EcoNI (287) **Bsu36I (291)**

301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
NaeI (441) 401 GGGCCTTTGTCCGGCGCTCCCTTGGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT

SphI (560)
KasI (535) 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTGAGCATGCCAAGAAGCCGACGCCCATCCAGCTGAACCC
AgeI (552) PvuII (589)
1 M P K K K P T P I Q L N P

HpaI (619) 601 GGCCCCGACGGCTCTGCAGTTAACGGGACCAGCTCTGCGGAGACCAACTTGGAGGCCTTGCAAGAAGCTGGAGGAGCTAGAGCTTGTGAGCAGCAG
13 A P D G S A V N G T S S A E T N L E A L Q K K L E E L E L D E Q Q
701 CGAAAGCGCCTTGAGGCCTTTCTTACCCAGAAGCAGAAGGTGGGAGAAGCTGAAGGATGACGACTTTGAGAAGATCAGTGAAGCTGGGGGCTGGCAATGGCG
47 R K R L E A F L T Q K Q K V G E L K D D D F E K I S E L G A G N G

XcmI (819) 801 GTGTGGTGTTCAGGCTCTCCACAAGCCTTCTGGCCTGGTCATGGCCAGAAGCTAATTCATCTGGAGATCAAACCCGCAATCCGGAACCAGATCATAAG
80 G V V F K V S H K P S G L V M A R K L I H L E I K P A I R N Q I I R

BsaBI (975) 901 GGAGCTGCAGTTCTGCATGAGTGAACCTCTCCGTACATCGTGGGCTTCTATGGTGGCTTCTACAGCATGGCGAGATCAGTATCGATGGAGCACATG
113 E L Q V L H E C N S P Y I V G F Y G A F Y S D G E I S I C M E H M

EcoRI (1040) 1001 GATGGAGTTCTCTGGATCAAGTCTGAAGAAAGCTGGAAGAATTCTGAACAAATTTAGGAAAAGTTAGCATTGCTGTAATAAAAAGCCCTGACATATC
147 D G G S L D Q V L K K A G R I P E Q I L G K V S I A V I K G L T Y
1101 TGAGGGAGAAGCAAGATCATGCACAGAGATGTCAAGCCCTCAACATCTAGTCAACTCCCGTGGGAGATCAAGCTCTGTGACTTTGGGTCAGCGG
180 L R E K H K I M H R D V K P S N I L V N S R G E I K L C D F G V S G

XcmI (1220) 1201 GCAGCTCATCGACTCCATGGCCAACTCCTTCTGTTGGCACAAGGTCTACATGTCCGCAAGAAAGCTCCAGGGGACTCATTACTCTGTGCAGTCAGACATC
213 Q L I D S M A N S F V G T R S Y M S P E R L Q G T H Y S V Q S D I
1301 TGGAGCATGGGACTGTCTCTGGTAGAGATGGCGGTTGGGAGGTATCCCATCCCTCCAGATGCCAAGGAGCTGGAGCTGATGTTGGGTGCCAGGTGG
247 W S M G L S L V E M A V G R Y P I P P P D A K E L E L M F G C Q V

XmaI (1438) 1401 AAGGAGATGCGGTGAGACCCACCCAGGCCAAGGACCCCGGAGGCCCTTAGCTCATACGGAATGGACAGCCGACCTCCCATGGCAATTTTTGAGTT
280 E G D A A E T P P R P R T P G R P L S S Y G M D S R P P M A I F E L
1501 GTTGATTACATAGTCAACGAGCCTCCTCAAAACTGCCAGTGGAGTTCAGTCTGGAATTTCAAGATTTGTGAATAAATGCTTAATAAAAAACCC
313 L D Y I V N E P P P K L P S G V F S L E F Q D F V N K C L I K N P

NeoI (1481)

BglII (1648) 1601 GCAGAGAGCAGATTTGAAGCAACTCATGTTTATCAAGAGATCTGATGCTGAGGAAAGTGGATTTTGCAGGTTGGCTCTGCTCCACCATCG
347 A E R A D L K Q L M V H A F I K R S D A E E V D F A G W L C S T I

NheI (1765) 1701 GCCTAACAGCCAGCACACCAACCCATGCTGCTGGCGTCTAAGTGTGGGAAGCAACAAAGAGCTAGCTGGCCAGACATGATAAGATACATTGATGA
380 G L N Q P S T P T H A A G V •

1801 GTTTGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAA

HpaI (1903) **MfeI (1914)** 1901 CAAGTTAAACAACA AATTGCATTCAATTTATGTTTCAGGTT CAGGGGAGGTGTGGGAGT TTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGG
EcoRI (1999)

2001 AATTCTAAAATACAGCATAGCAAACCTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTT
2101 GCCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTTAAAGATATAGTGTATTTCCCAAGGTTTGAAGTCTCTTCAATTTCTTTATGT

SspI (2238) **SwaI (2252)** 2201 TTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGA
2301 ATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATGGACAGCAAGAAAGCGAGC
2401 TTCTAGCTTTAGTTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCTCAATGAGCACAAAGCAGTCAGGAG
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

SacI (2513) 2501 CATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGT
110 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 R V A V I T
2601 GTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTACGACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAG
77 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 E I H V A S

