



PvuI (7)
SgfI (6)
MfeI (82)

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGCGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTGTCTGTACTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)

Psp1406I (203)
PvuII (239)
Bsu36I (291)

201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACCGCGCCGCCCTACCTGAGGGC

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

AgeI (552)
BspLUII (568)
SapI (581)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCTACCTGAGATCACCGGTAGGAGGGCCAACATGTCCAAGCTGCTCTTCTGCTCACTTGGC

1► M S K L L F L S L A

BsrBI (611)

601 CTCTGGGCGACCGCTCCCTGGTTACTGAAACAGCTCTCGTGGCTCTAAGCCAGCCAGAGTGCATGCCATGCTTCTCGGTATCCCGTGGCCGTGG

11► L W A S R S P G Y T E T A L V A L S Q P R V Q C H A S R Y P V A V

701 ACTGCTCTGGACTCCTCTCCAGGCTCCCACTCCACCAGATCCACGTCTTCTCATTGCCACTTACAGGCTCGGTGGCCACCAGCAGCAGAGCCAGCC

44► D C S W T P L Q A P N S T R S T S F I A T Y R L G V A T Q Q Q S Q P

801 CTGCCTACAACGGAGCCCCAGGCTCCCGATGCACCATCCCCGACGTGCACCTGTTCTCCACGGTGCCTACATGCTAAATGTCAGTGCAGTGCACCA

77► C L Q R S P Q A S R C T I P D V H L F S T V P Y M L N V T A V H P

BspEI (959)
FspI (976)
SdaI (995)

901 GCGGGCGCCAGCAGCAGCTCCTAGCCTTTGTGGCTGAGCGAATCATCAAGCCGACCTCCGGAAGCGTGGCCTGCGCACAGCGGGACAGCGCCTGC

111► G G A S S S L L A F V A E R I I K P D P P E G V R L R T A G Q R L

1001 AGGTGCTCTGGCATCCCCCTGCTTCTGGCCCTTCCGGACATCTTCTCTCAAGTACCGACTCCGCTACCGGCGCCGAGGAGCCTCTCACTTCCGCCA

144► Q V L W H P P A S W P F P D I F S L K Y R L R Y R R R G A S H F R Q

SandI (1103)
Bsu36I (1130)

1101 GGTGGGACCCATTGAAGCCAGACTTTACCCTCAGGAACCTGAAACCCCATGCCAAGTATTGCATCCAGGTGTCAGCTCAGGACCTCACAGATTATGGG

177► V G P I E A T T F T L R N S K P H A K Y C I Q V S A Q D L T D Y G

BstXI (1203)
Acc65I (1268)
AvrII (1293)

1201 AAACCAAGTGACTGGAGCCTCCCTGGGCAAGTAGAAAGTGCACCCATAAGCCC**GTTCTGGAGTAGGGGTACCTGGGGTGGGCCTGGCTGTGCTTAGGA**

211► K P S D W S L P G Q V E S A P H K P 1► V P G V G V P G V G 1► L A V P R

NheI (1352)

SphI (1348)
NdeI (1383)

1301 GTAGCAGTCTGACTGGGCTCAGTGCCAGCAGCTCTCTCGAATCTCTGCATGCTAGCTGGAACGCACATGCACCAGCGGCACATATGAATCTACTAAG

6► S S S P D W A Q C Q Q L S R N L C M L A W N A H A P A G H M N L L R

XcmI (1442)

1401 AGAAGAAGAGGATGAAGAGACTAAAAATAATGTGCCCCGTATCCAGTGTGAAGATGGTTGTGACCCACAAGGACTCAAGGACAACAGCCAGTTCTGCTTG

39► E E E D E E T K N N V P R I Q C E D G C D P Q G L K D N S Q F C L

BamHI (1504)
NcoI (1588)

1501 CAAAGGATCCGCCAAGGCTGGCTTTTTATAAGCACCTGCTTACTCTGACATCTTCAAAGGGGAGCCTGCTCTACTCCCTGATAGCCCATGGAGCAAC

73► Q R I R Q G L A F Y K H L L D S D I F K G E P A L L P D S P M E Q

1601 TTCACACCTCCCTACTAGGACTCAGCAACTCTCCAGCCAGAGGATCACCCCGGAGACCAACAGATGCCAGCCTGAGTTCTAGTCAGCAGTGGCA

106► L H T S L L G L S Q L L Q P E D H P R E T Q Q M P S L S S S Q Q W Q

BstBI (1726)
SfiI (1739)

1701 GCGCCCCCTTCTCCGTTCCAAGATCCTTCGAAGCCTCCAGGCTTTTTGGCCATAGCTGCCCGGGTCTTTGCCACGGAGCAGCAACTCTGACTGAGCCC

139► R P L L R S K I L R S L Q A F L A I A A R V F A H G A A T L T E P

NheI (1842)

EcoRI (1836)

1801 TTAGTGCCAACAGCTTAAGGATGCCAGGTTCCCATGAATCG**CTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAA**

173► L V P T A •

HpaI (1980)
MfeI (1991)

1901 TGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAAATTGCAT

EcoRI (2076)

2001 TCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGA**AATTCTAAATACAGCATAGCAA**

2101 AACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCCAATGTGCATTAGCTGTTTGC

SapI (2258)

2201 AGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGTTTGAACAGCTCTTCATTCTTTATGTTTTAAATGCACTGACCTCCACA

