

PvuI (7)
SgfI (6) 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
MfeI (82)
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGTCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACC GTATATAAGTGCAGTAGTCGCC

Psp1406I (203) 201 GTGAACGTTCTTTTTTCGAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACCGCGCCCGCCCTACCTGAGGCC
HindIII (245)
Bsu36I (291)
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTT
NgoMIV (441)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGCTCCGGACACCATGGAGTTCCTGCCAGACCCAGCATCCCC
BspEI (558)
1 M E F L P G P Q H P P
Bsu36I (680)
601 AGGGCCCCAACCATGGACTTGGAGGAGCCAAAGGCCCGAGGTGCCCTCTGAGAACCACCCTCTAACACCAGTCGGCCTTAGGGCTGGAGGCCCT
11 G P P T M D L E E P K G P E V P S E N H P S N T Q S A L G P G G P
701 GTCACCCTGAGTGAGATGGAGCTGGACACATCCAGTGTGAGGAGCTGGTTCAGCAGCTAGAAGCTTACCCAGCGACCTGGTGGCCATTCCAGATG
45 V T L S E M E L D T S S V Q E L V Q Q L E A L P S D L G G P F P D
801 GAGCCCCCTGTCCCTGCACATTGCTACCGCCAAGGACTTGCACCCAGGAGAACCAGATGCTGGTGGGCTTGTGCGCCGAGGCAGGCGGGGACGA
78 G A P C P L H I A T G Q G L A T Q E N P D A G G L L S A E A G G D D
BstAPI (961)
901 CCTTCTGGGTCTGCTGAGGGATGAGGCCTCTTCGCTGCCAATCTGTCCCCAAGATCTGCACAGACTGCACCCCGCCTTCTGCAGCCGCCAGAGGAC
111 L L G L L R D E A S S P A Q S V P Q D P A Q T A P R L L Q P P E D
XmaI (1011)
SandI (1007)
1001 CCTGATGGGACCCGGATGGATGGAAGGGCATCAGCAGAGCCAGCAGACAGCAGGAGCTCCAGCAGTTCGCCGGAGCCCTGGCTGGAGACAGCACCCC
145 P D G D P G W M E G A S A E P A D S R S S S S S P E P W L E T A P
Acc65I (1126)
1101 TGGTGACACAGCAGGAACACCTGTGGGTACCCAGAGCCGGGAGACCTGGCCTCGTGCCCTGCCGTACCCAGGTCACAGGTCACAGGTCCTGTGGGCGGAGGA
178 L V T Q Q E P P V G T Q S R E T L A S C P A V T E V P G P C G P E E
1201 GCTCATGGATGGTGTCTCTCCGGGCGAAATACCTGGGCTCCACCCAGCTGTGTCAGAGCGCAGCCCGGCGCCAGCACACGCATGGGGCAGGCGCAG
211 L M D G V I F G A K Y L G S T Q L L S E R S P A P S T R M G Q A Q
1301 GAGGCCATGGACCGCTCAAGGCCCTGAAGGCCGAGACCCAGCCATGGTGAAGTGGACATCTTTATCTCCACCAAGCGTCAAAGTGTGGCTGCTG
245 E A M D R V K A P E G E T Q P M V E V D I F I S T K R V K V L A A
MscI (1493)
1401 ACTCCAGGATGCCCTGATGGACCATGCCCTGCAGACCATCTCTACATCGCAGACATTGGGCTGTACTGGTCTTGTATGGCCCGGAGGGCTGGCCAG
278 D S Q D A L M D H A L Q T I S Y I A D I G P V L V L M A R R R L A R
BbsI (1552)
1501 GAGGACAACCTCCCAGGACCCAGCGGCGACTCTACAAGATGCTGTGCCATGCTTCCACTCAGAGGATGCCAGCTCATAGCACAAGCCATCGGCCAG
311 R T T P Q D R Q R R L Y K M L C H V F H S E D A Q L I A Q A I G Q
1601 GCCTTCAGCATCGCTACAGCCAGTTCTACAGGAGAACAGGATCGACCCAGCCAGGTGGGACGACGCCCTCAACCCTGCGAGCCACCCGCACAACG
345 A F S I A Y S Q F L Q E N R I D P S Q V G T Q P S T A A S H P H N
1701 GCGACCTGGACCACTTCTGCAACAGCCAGAAGTGTAGGGAGTTTGCATCCAGAAGCGCCAGGTGAAGGCCTGGGCGTCCGCTAGTTGAATCAGGCTG
378 G D L D H F C N S Q N C R E V C I Q K R P G E G L G V A L V E S G W
Eco47III (1856)
1801 GGGCTCGCTACTGCCACTGCGGTGCTTCCAACTGCTCCATGGGGCCCTGCCAGCGCTGCGGGCCCTGAGCATAGGGACCGCTCACTGCCATC
411 G S L L P T A V I A N L L H G G P A E R C G A L S I G D R V T A I
1901 AATGGAACCAGCCTGGTGGGCTGTGCTGGCTGCCTGCCAGGCTGCGGTGCGGAGGTGCGCGACTTTCATCCGTGACGCTGAGTATTATCCACTGCC
445 N G T S L V G L S L A A C Q A A V R E V R R L S S V T L S I I H C
BstEII (2004) **BbrPI (2032)**
AgeI (2001) **NgoMIV (2024)**
2001 CACCGGTACCACGGCCGTCATCCGCCGCCACAGTGCAGGAGCAGTGGGCTTCTGTGTGGAGGATGGAATCATTGTAGCCTGCTGCGTGGGGCGC
478 P P V T T A V I R R P H V R E Q L G F C V E D G I I C S L L R G G A
BssHII (2176)
2101 TGCTGAGCGTGGTGGAGTCCGCGTTGGACACCGCATCATTGAGGTCAACGGGAGAGCGTGGTGGCTATGCCCATGCGCGCATCATCCAGCTGCTACA
511 A E R G G V R V G H R I I E V N G Q S V V A M P H A R I I Q L L T
MscI (2297)
Bst1107I (2272) **NheI (2291)**
2201 GAGACAAGAGAAATCCATCAAGACCATGCCAGCTGCCACCTACCGACTGCTCACGGGACAGGAACAACAGTATACCTATGACTGCCCCGTAGCTGG
545 E T R E I H I K T M P A A T Y R L L T G Q E Q P V Y L •
2301 CCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAAGTGAATGCAAGTGAATAATGCTTTATTGTAATTTGTGATGCTATTGCTTTAT

2401 TTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCTTTTATGTTTCAGGTTGAGGGGAGGTGTGGAGGTTTTTAAAGCAA
HpaI (2429) **MfeI (2440)**

