



PvuI (7)
SgfI (6) 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
MfeI (82) **EcoNI (96)**
101 GAGAAGTGGCGCGGGTAAACTGGAAAGTGATGCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **HindIII (245)** **EcoNI (287)**
201 GTGAACGTTCTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC
301 GCCATCCAGCGCGGTTGAGTCGCGTTCGCCGCCCTCCCGCTGTGGTGCCTCCTGAAGCTCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560) **BstEII (555)** **Bsp120I**
KasI (535) **AgeI (552)** **XmaI (595)**
501 TCTGTTCTGCGCGGTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTACCATTGGCGATTGGCGCTGCTGGCCACGGGTGTCCTCCCGG
1▶ M A I R R C W P R V V P G

Psp1406I (631) **XmaI (642)** **BssHIII (689)**
601 GCCCGCGCTGGGATGGCTGTTCTGCTGCTGCTGAACGTTCTGGCCCGGGCCGCGCTCCCGCGCTCCTCGACTTCCCGGCTCCGGTCTGCGCGCAGGAG
13▶ P A L G W L L L L L N V L A P G R A S P R L L D F P A P V C A Q E

ScaI (724) **Tth111I (765)**
701 GGGCTGAGCTGCAGAGTCAAGAATAGTACTTGTCTGGATGACAGCTGGATCCACCCCAAAAACCTGACCCCGTCTTCCCAAAAAACATCTATATCAATC
47▶ G L S C R V K N S T C L D D S W I H P K N L T P S S P K N I Y I N
801 TTAGTGTTCCTCTACCCAGCAGGAGAATTAGTCCCTGTGTTGATGTTGAGTGGACCTGCAGACAGATGCCAGCATCTGTACNITCAGGTTGCAGA
80▶ L S V S S T Q H G E L V P V L H V E W T L Q T D A S I L Y L E G A E

BsrBI (928)
901 GCTGTCCGTCTGCAGCTGAACCAATGAGCGGCTGTGTGTCAGTTCCAGTTTCTGTCCATGCTGCAGCATCACCGTAAGCGGTGGCGGTTTTCTTC
113▶ L S V L Q L N T N E R L C V K F Q F L S M L Q H H R K R W R F S F

SandI (1073)
1001 AGCCACTTTGTGGTAGATCCTGGCCAGGAGTATGAAGTACTGTTACCACCTGCCGAAGCCCATCCCTGATGGGGACCCAAACCACAAATCCAAGATCA
147▶ S H F V V D P G Q E Y E V T V H H L P K P I P D G D P N H K S K I

XcmI (1175)
1101 TCTTTGTGCTGACTGTGAGGACAGCAAGATGAAGTACTACCTCATGCGTGTGAGCTCAGGCAGCCTTTGGGATCCCAACATCACTGTGGAGACCTTGA
180▶ I F V P D C E D S K M K M T T S C V S S G S L W D P N I T V E T L D
1201 CACACAGCATCTGCGAGTGGACTTACCCTGTGGAATGAATCCACCCCTACCAGTCTGCTGGAAAGTTTCTCGACTCAGAGAACCACAGCTGCTTT
213▶ T Q H L R V D F T L W N E S T P Y Q V L L E S F S D S E N H S C F

SspI (1314) **EcoRI (1336)** **BbrPI**
1301 GATGTCGTTAAACAAATTTGGCGCCAGGCAAGAAGAATCCATCAGCGAGCTAATGTCACATTCACCTAAGCAAGTTTCACTGGTGTGCCATCACCC
247▶ D V V K Q I F A P R Q E E F H Q R A N V T F T L S K F H W C C H H
1401 ACCTGTCAGTCCAGCCCTTCTTACGAGCTGCCTAAATGACTGTTGAGACAGCTGTGACTGTGCCCTGCCAGTAATCTCAAATACACAGTTCCCAA
280▶ H V Q V Q P F F S S C L N D C L R H A V T V P C P V I S N T T V P K

XcmI (1553) **BstXI (1553)**
1501 GCCAGTTGCAGACTACATTCCTCCCTGTGGGTGTATGGCCTCATCACTCATCGCCATTCTGCTGGTGGATCTGTCTATGCTGATCATCTGTATGACC
313▶ P V A D Y I P L W V Y G L I T L I A I L L V G S V I V L I I C M T

KasI (1612)
1601 TGGAGGCTTTCTGGCGCCATCAAGAGAAACATGGTGTGACTCCAAAATCAATGGCATCTTGCCCGTAGCAGACCTGACTCCCCACCCTGAGGCCCA
347▶ W R L S G A D Q E K H G D D S K I N G I L P V A D L T P P P L R P
1701 GGAAGTCTGGATCGTCTACTCGGCCGACCCCTCTATGTGGAGTGGTCTAAAGTTTCCGCCAGTTCCTGATCACTGCCTGTGGCACTGAAGTAGC
380▶ R K V W I V Y S A D H P L Y V E V V L K F A Q F L I T A C G T E V A

BspHI (1841)
1801 CCTTGACCTCTGGAAGAGCAGGTTATCTCTGAGGTGGGGTCACTGACCTGGGTGAGCCGACAGAAGCAGGAGATGGTGGAGAGCAACTCCAAAATCATC
413▶ L D L L E E Q V I S E V G V M T W V S R Q K Q E M V E S N S K I I
1901 ATCCTGTGTTCCCGAGGCACCAAGCAAAGTGGAAAGTATCTTGGTGGGGTGTGAGCCTGCTGTCCAGCTACGGTGTGACCACTGGAAGCCTGCTGGGG
447▶ I L C S R G T Q A K W K A I L G W A E P A V Q L R C D H W K P A G
2001 ACCTTTCACTGCAGCCATGAACATGATCCTGCCAGACTTCAAGAGGCCAGCTGCTTCCGACACTACGTTGTTGCTACTTCACTGAGGATCTGTAGTGA
480▶ D L F T A A M N M I L P D F K R P A C F G T Y V V C Y F S G I C S E

Acc65I (2135) **BspEI (2171)**
2101 GAGGGATGTCCCCGACCTCTTCAACATCACCTCAGGTTACCCACTCATGGACAGATTTGAGGAGGTTTACTTCCGGATCCAGGACCTGGAGATGTTTGA
513▶ R D V P D L F N I T S R Y P L M D R F E E V Y F R I Q D L E M F E

NgoMIV (2261)
2201 CCCGGCCGGATGCACCATGTGAGAGAGCTCACAGGGGACAATTACCTGCAGAGCCCTAGTGGCCGGCAGCTCAAGGAGGCTGTGCTTAGGTTCCAGGAGT
547▶ P G R M H H V R E L T G D N Y L Q S P S G R Q L K E A V L R F Q E

BglIII (2358)
2301 GGCAAACCCAGTGCCCCGACTGGTTTCGAGCGTGAGAACCCTCTGCTTAGCTGATGGCCAAAGATCTCCCTCCCTGGATGAAGAAGTGTGTAAGACCCACT
580▶ W Q T Q C P D W F E R E N L C L A D G Q D L P S L D E E V F E D P L
2401 GCTGCCACAGGGGGAGGAATTGTCAAACAGCAGCCCTGGTGGCGGAACTCCATCTGACGGCTGCCTGTGGTAGATGTCTGTGTCAGTGAGGAAGAA
613▶ L P P G G G I V K Q Q P L V R E L P S D G C L V V D V C V S E E E
2501 AGTAGAATGGCAAAGCTGACCTCAGCTATGGCCACAGAGAGCTAGTGGPTCACACCTCCAAAGCATGGTGTGCCAGCAGGACAGGTCCTGCAG
647▶ S R M A K L D P Q L W P Q R E L V A H T L Q S M V L P A E Q V P A
2601 CTCATGTGGTGGAGCCTCTCATCTCCAGACGGCAGTGGAGCAGCTGCCAGCTGCCATGACAGAGGACAGCGAGGCTTGGCCGCTGCTGGGGTCCA
680▶ A H V V E P L H L P D G S G A A A Q L P M T E D S E A C P L L G V Q

