



PvuI (7)
SgfI (6)
1 GGATCTGCATCGCTCCGGTGCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTCACGCGCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

AgeI (552) **BspHI (568)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTAGGAGGGCCATCATGATGCCTCCCTGGCTCCTGGCTAGGACT
1► M M P P W L L A R T
SspI (669) **NsiI (683)**

601 CTGATCATGGCACTGTTCTTCTCCTGCCTGACACCAGGAAGCTTGAATCCCTGCATAGAGGTAGTTCCTAATATTACCTACCAATGCATGGATCAGAAAC
11► L I M A L F F S C L T P G S L N P C I E V V P N I T Y Q C M D Q K
701 TCAGCAAAGTCCCTGATGACATTCCTTCTCAACCAAGAACATAGATCTGAGCTTCAACCCCTTGAAGATCTTAAAAAGCTATAGCTTCTCCAATTTTTTC
44► L S K V P D D I P S S T K N I D L S F N P L K I L K S Y S F S N F S

BbsI (846)
801 AGAACTTCAGTGGCTGGATTATCCAGGTGTGAAATTGAAACAATTGAAGACAAGGCATGGCATGGCTTACACCACCTCTCAAACCTGATACTGACAGGA
77► E L Q W L D L S R C E I E T I E D K A W H G L H H L S N L I L T G
901 AACCTATCCAGAGTTTTTCCCGAGGAAGTTTCTGGACTAAACAAGTTTAGAGAATCTGGTGGCTGTGGAGACAAAATTGGCCTCTCTAGAAAGCTTCC
111► N P I Q S F S P G S F S G L T S L E N L V A V E T K L A S L E S F
1001 CTATTGGACAGCTTATAACCTTAAAGAACTCAATGTGGCTCACAAATTTTATACATTCCTGTAAGTTACCTGCATATTTTTCCAATCTGACGACCTAGT
144► P I G Q L I T L K K L N V A H N F I H S C K L P A Y F S N L T N L V
1101 ACATGTGGATCTTTTATAACTATATTCAAACCTTACTGTCAACGACTTACAGTTTCTACGTGAAAATCCACAAGTCAATCTCTTTAGACATATCT
177► H V D L S Y N Y I Q T I T V N D L Q F L R E N P Q V N L S L D I S
1201 TTGAACCAATTGACTTCATTCAAGACCAAGCCTTTGAGGAAATTAAGCTCCATGAACTGACTGAACTGAAAGGTAATTTAATAGCTCAAATATAATGAAAA
211► L N P I D F I Q D Q A F Q G I K L H E L T L R G N F N S S N I M K
1301 CTTGCCTTCAAACCTGGCTGGTTTACACATCCATCGGTTGATCTTGGGAGAAATTTAAAGTAAAGGAATCTGAAATTTTTGAACCCTCTATCATGGA
244► T C L Q N L A G L H I H R L I L G E F K D E R N L E I F E P S I M E

HpaI (1433)
1401 AGGACTATGTGATGTGACCATTGATGAGTTTCAAGTTAACAATGATTTTTTCAAGTATGATTTGTTAAGTTCCATTGCTTGGCGAATGTTTCTGCA
277► G L C D V T I D E F R L T Y T N D F S D D I V K F H C L A N V S A
1501 ATGTCTCTGGCAGGTGTATCTATAAAATCTAGAAGATGTTCTAAACATTTCAAATGGCAATCCTTATCAATCATTAGATGCAACTTAAGCAGTTTC
311► M S L A G V S I K Y L E D V P K H F K W Q S L S I I R C Q L K Q F
1601 CAACTCTGGATCTACCCTTTCTAAAAGTTTGACTTTAACTATGAACAAAGGGTCTATCAGTTTTAAAAAGTGGCCCTACCAAGTCTCAGCTATCTAGA
344► P T L D L P F L K S L T L T M N K G S I S F K K V A L P S L S Y L D
1701 TCTTAGTAGAAAATGCACTGAGCTTTAGTGGTTGCTGTTCTTATTCTGATTTGGGAACAACAGCCTGAGACACTTAGACCTCAGCTTCAATGGTGCCATC
377► L S R N A L S F S G C C S Y S D L G T N S L R H L D L S F N G A I

BstXI (1810) **SapI (1827)** **EcoRI (1878)**
1801 ATTATGAGTGCCAAATTTCAAGGCTAGAAAGCTGCAGCACCTGGATTTTTCAGCACTCTACTTTAAAAAGGGTACAGAAATTTCTCAGCGTTCTTATCCC
411► I M S A N F M G L E E L Q H L D F Q H S T L K R V T E F S A F L S
1901 TTGAAAAGCTACTTTACCTTGACATCTTATACTAACACCAAAATGACTTCGATGGTATATTTCTTGGCTTACCAGTCTCAACACATTAATAATGGC
444► L E K L L Y L D I S Y T N T K I D F D G I F L G L T S L N T L K M A
2001 TGGCAATTTCTTCAAAGACAACACCCTTTCAAATGTCTTTGCAAAACAACAACTTACATTCCTGGATCTTTCTAAATGTCAATTTGGAACAATATCT
477► G N S F K D N T L S N V F A N T T N L T F L D L S K C Q L E Q I S

PvuII (2187)
2101 TGGGGGTATTTGACACCCTCCATAGACTTCAATTATTAATATGAGTCAACAATCTATTGTTTTGGATTATCCATTATAACCGCTGATTATCCC
511► W G V F D T L H R L Q L L N M S H N N L L F L D S S H Y N Q L Y S
2201 TCAGCACTCTTGATTGCAGTTTCAATCGCATAGAGACATCTAAAGGAATACTGCAACATTTTCAAAGAGTCTAGCCTTCTTCAATCTTACTAATAATTC
544► L S T L D C S F N R I E T S K G I L Q H F P K S L A F F N L T N N S
2301 TGTGCTTGTATATGTGAACATCAGAAATTCCTGCAGTGGGTCAAGGACCAGAAGCAGTTCTTGGTGAATGTTGAACAAATGACATGTGCAACACCTGTA
577► V A C I C E H Q K F L Q W V K D Q K Q F L V N V E Q M T C A T P V

BsrGI (2449)
2401 GAGATGAATACCTCCTTAGTGTGGATTTAATAATTCTACCTGTTATATGTACAAGACAATCATCAGTGTGTGAGTGGTGTGATTGTGGTATCCA
611► E M N T S L V L D F N N S T C Y M Y K T I I S V S V V S V I V V S

BamHI (2580)
2501 CTGTAGCATTCTGATATACCACTTCTATTTTACCTGATACTTATTGCTGGCTGTAAGGTAACAGCAGAGGAGAAAAGCGGATCCTACCCTATGATGT
644► T V A F L I Y H F Y F H L I L I A G C K K Y S R G E S G S Y P Y D V

MscI (2623)
NheI (2617)
2601 GCCAGACTACGCCATAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTG
677► P D Y A •

2701 **HpaI (2755)**
TGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAACAAACAACAATTGCATTATTTTATGTTTCAGGTTTCAGGGG

2801 **EcoRI (2851)**
GAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAAATTCTAAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTC

2901
TACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTT

3001 **SapI (3033)** **SspI (3090)**
AAGATATAGTGTATTTTCCCAAGGTTTGAACAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCTTTTATGTAATAATATTGCA

3101 **SwaI (3104)** **EcoO109I (3165)**
AATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGG

3201
ACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGAGGGGGATGAGTTCCTCAATG

3301 **SacI (3365)** **BstXI (3394)**
GTGTTTTGACCAGTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGACATGCCACAGGGGCTGACCACCC

3401
127 T T K V L K G N M E I L V F C D P A Y D S I L E R C M G C P S V V R
94 I S R D V E D S Y P H R V A V I T D F D K Q G N S V A S G I A I A

3501 **StuI (3529)**
TTCAGCACAGACAGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGCTTG

3601 **BbsI (3675)**
61 E A C V T V R G I Y A E I H V A S I I E G T K T R I A A G V H H K
3601 TTGTCCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCTGCTGAGAGATGTTGAAGGTCTTCATGGTGGCCCTCTATAGT

3701 **AseI (3737)** **SacI (3794)**
27 N D E Y L M T I K E T A V E V L E L D Q Q S I N F T K M
GAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTACTAAACGAGCTC

3801 **SpeI (3892)**
TGCTTATATAGACCTCCACCGTACACGCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTACGACATTTTGAAAAGTCCCGTTGATTTACTAGTCA

3901
AAACAACTCCCATTGACGTCAATGGGGTGAGACTTGAAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTAAGTCCAAACCGCATCATC

4001 **SnaBI (4020)**
ATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCATTACCCTCAT

4101 **NdeI (4125)**
TGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGACTGCCAAGTGGGCAGTTTACCCTAAATACTCCACCATTGACGTCAATGAAAAGT

4201
CCCTATTGGCGTTACTATGGGAACATACGTATTATTGACGTCAATGGGCGGGGCTGTTGGCGGTCAGCCAGGCGGGCATTACCCTAAGTTATGTA

4301 **PaeI (4311)** **SdaI (4303)**
ACGCTGCAGGTTAATAAGAACATGTGAGCAAAAAGCCAGCAAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGCTCCGCC

4401
CCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTG

4501
CGCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCA

4601 **ApaLI (4635)**
GTTGCGGTGATAGTTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCCAA

4701
CCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGTACAGAGTCTTGAAGTGGTGG

4801
CCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGTATCCGGCAAAC

4901
AAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTC

5001 **EagI (5071)** **PaeI (5051)** **SwaI (5060)** **NotI (5070)**
TGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCAT

5101
TACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCCAG

5201
TGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA