



PvuI (7)
SgfI (6)
MfeI (82) **EcoNI (96)**

1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245) **Bsu36I (291)**

201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCCCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCAGTTCGCCGCTCCCGCCTGTGGTGCCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441) **NaeI (441)**

401 GGGCCTTTGTCCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT

KasI (535)
AgeI (552)
BspHI (568)

501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTAGGAGGGCCATCATGAGACTCATCAGAAACATTTACATATTT
 601 TGTAGTATTGTTATGACAGCAGAGGGTGTGCTCCAGAGCTGCCAGAAGAAAGGGAAGTATGACCAACTGCTCCAACATGTCTCTAAGAAAGGTTCCCG
 11▶ C S I V M T A E G D A P E L P E E R E L M T N C S N M S L R K V P

XcmI (715)

701 CAGACTTGACCCAGCCACAACGACACTGGATTTATCCTATAACCTCCTTTTTCAACTCCAGAGTTGAGATTTTCTGCTCCAACTGAGAGTTTT
 44▶ A D L T P A T T T L D L S Y N L L F Q L Q S S D F H S V S K L R V L

PvuII (825)

801 GATTCTATGCCATAACAGAATTCACAGCTGACTCTAAAACCTTTGAATTCACAAGGAGTTAAGATATTTAGATTTGTCTAATAACAGACTGAAGAGT
 77▶ I L C H N R I Q Q L D L K T F E F N K E L R Y L D L S N N R L K S

BglII (935)

901 GAACTTGGTATTTACTGGCAGGTCTCAGGTATTTAGATCTTTCTTTAATGACTTTGACACCATGCCTATCTGTGAGGAAGCTGGCAACATGTCACACC
 111▶ V T W Y L L A G L R Y L D L S F N D F D T M P I C E E A G N M S H

AvrII (1007)

1001 TGGAAATCCTAGGTTTGTGAGTGGGCAAAAATACAAAATCAGATTTCCAGAAAATGCTCATCTGCATCTAAATACTGTCTTCTTAGGATTCCAGAACTCT
 144▶ L E I L G L S G A K I Q K S D F Q K I A H L H L N T V F L G F R T L
 1101 TCCTCATTATGAAGAAGGTAGCCTGCCATCTTAAACACAACAAAAGTGCACATTTGTTTACCAATGGACACAAAATTTCTGGGTTCTTTTGCCTGATGGA
 177▶ P H Y E E G S L P I L N T T K L H I V L P M D T N F W V L L R D G
 1201 ATCAAGACTTCAAAAATATTAGAAATGACAAAATATAGATGGCAAAAAGCAATTTGTAAGTTATGAAATGCAACGAAATCTTAGTTTAAAAATGCTAAGA
 211▶ I K T S K I L E M T N I D G K S Q F V S Y E M Q R N L S L E N A K
 1301 CATCGTTCTATTGCTTAATAAAGTTGATTTACTCTGGGACGACCTTTTCTTATCTTACAATTTGTTTGGCATAATCAGTGGAAACATTTCCAGATCCG
 244▶ T S V L L L N K V D L L W D D L F L I L Q F V W H T S V E H F Q I R

BsrGI (1487)

1401 AAATGTGACTTTTGGTGGTAAGGCTTATCTTGACCACAATTCATTTGACTACTCAAATACTGTAATGAGAAGTATAAAATGGAGCATGTACATTTTCCAGA
 277▶ N V T F G G K A Y L D H N S F D Y S N T V M R T I K L E H V H F R
 1501 GTGTTTTACATTCACAGGATAAAAATCTATTTGCTTTTACCAAAAATGGACATAGAAAACCTGACAATATCAAATGCACAAAATGCCACACATGCTTTTCC
 311▶ V F Y I Q Q D K I Y L L L L T K M D I E N L T I S N A Q M P H M L F
 1601 CGAATTATCCTACGAAATTCCTAATTTAAATTTTGCCTAATAATCTTAAACAGAGAGTTGTTTAAAAGAAGTATCCAACTGCCTCAGTTGAAAACCTCT
 344▶ P N Y P T K F Q Y L N F A N N I L T D E L F K R T I Q L P H L K T L
 1701 CATTTTGAATGGCAATAAAGTGGAGACACTTTCTTTAGTAAGTTGCTTTGCTAACAACACACCTTGGAAACACTTGGATCTGAGTCAAATCTATTACAA
 377▶ I L N G N K L E T L S L V S C F A N N T P L E H L D L S Q N L L Q

MscI (1824)

1801 CATAAAAATGATGAAAATGCTCATGGCCAGAAAAGTGGTCAATATGAATCTGTCATACAATAAATGTCTGATTCTGTCTTCCAGGTGCTTGCCAAAA
 411▶ H K N D E N C S W P E T V V N M N L S Y N K L S D S V F R C L P K
 1901 GTATTCAAAATACTTGACCTAAATAATAACCAAAATCCAACTGTACCTAAGAGACTATTCATCTGATGGCCTTACGAGAACTAAAATTTGCATTAATTT
 444▶ S I Q I L D L N N N Q I Q T V P K E T I H L M A L R E L N I A F N F
 2001 TCTAACTGATCTCCCTGGATGCAGTCATTTAGTAGACTTTGAGTCTGAACATTGAAATGAACCTTCTCAGCCCATCTCTGGATTTTGTTCAGAGC
 477▶ L T D L P G C S H F S R L S V L N I E M N F I L S P S L D F V S

BspHI (2194)

2101 TGCCAGGAAGTAAAACTCTAAATGCGGGAAGAAATCCATTCCGGTGTACCTGTGAATAAAAAATTTTCAATCAGCTTCAAACATATTCCAGAGGTCATGA
 511▶ C Q E V K T L N A G R N P F R C T C E L K N F I Q L E T Y S E V M

Bsu36I (2243)

2201 TGGTTGGATGGTCAGATTCATACACCTGTGAATACCCTTTAAACCTAAGGGGAACTAGGTTAAAAAGACGTTTCATCTCCACGAATTATCTTGAACACAGC
 544▶ M V G W S D S Y T C E Y P L N L R G T R L K D V H L H E L S C N T A
 2301 TCTGTTGATTGTACCATTTGTGGTTATTATGCTAGTCTGGGGTGGCTGTGGCCCTTCTGCTGTCTCCACTTTGATCTGCCCTGGTATCTCAGGATGCTA
 577▶ L L I V T I V V I M L V L G L A V A F C C L H F D L P W Y L R M L
 2401 GGTCAATGCACACAAATCGCACAGGGTTAGGAAAAACCAAGAACTCAAGAGAAATGTCCGATTCCACGCATTTATTTCATACAGTGAACATG
 611▶ G Q C T Q T W H R V R K T T Q E Q L K R N V R F H A F I S Y S E H

XbaI (2534)

2501 ATTCTCTGTGGGTGAAGAATGAATGATCCCAATCTAGAGAAGGAAGATGGTCTATCTTGTATTTGCCTTTATGAAAGCTACTTTGACCCTGGCAAAAG
 644▶ D S L W V K N E L I P N L E K E D G S I L I C L Y E S Y F D P G K S

HindIII (2618)

2601 CATTAGTAAAAATATTGTAAGCTTCATTGAGAAAAGCTATAAGTCCATCTTTGTTTTGTCTCCAACTTTGTCCAGAATGAGTGGTGCCATTATGAATTC
 677▶ I S E N I V S F I E K S Y K S I F V L S P N F V Q N E W C H Y E F
 2701 TACTTTGCCCAACCAATCTCTTCCATGAAAATCTGATCATATAATCTTATCTTACTGGAACCCATTCCATTCTATTGCATTTCCACCAGGTATCATA
 711▶ Y F A H H N L F H E N S D H I I L I L L E P I P F Y C I P T R Y H

2801 AACTGAAAGCTCTCTGGAAAAAAGCATACTTGAATGGCCCAAGGATAGGCGTAAATGTGGGCTTTCTGGGCAAACCTTCGAGCTGCTATTAATGT
744▶ K L K A L L E K K A Y L E W P K D R R K C G L F W A N L R A A I N V

PstI (2931) XhoI (2962)

2901 TAATGTATTAGCCACCAGAGAAATGTATGAACTGCAGACATTACAGAGTTAAATGAAGAGTCTCGAGGTTCTACAATCTCTGATGAGAACAGATTGT
777▶ N V L A T R E M Y E L Q T F T E L N E E S R G S T I S L M R T D C

MscI (3049) NheI (3043)

3001 CTATAAAATCCCACAGCTCTTGGGAAGTTGGGGACCACATACAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGA
811▶ L •

HpaI (3181) MfeI (3192)

3101 ATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAAACAAGTTAAACAACAACATTGCA
3201 TTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCA
3301 AAACCTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTG
3401 CAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTTCCAAGGTTTGAAGTCTCTTCATTCTTTATGTTTTAAATGCACTGACCTCCCAC

EcoO109I (3591)

3501 ATTCCCTTTTATGAAAAATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTC
3601 ATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTAAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGT
141▶ • N R T Y

SacI (3791)

3701 ACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCT
136▶ 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 Y D S I L E R

BstXI (3820)

3801 GCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTG
103▶ 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 D F D K Q G N

StuI (3955)

3901 CTCACAGCAGACCAATGGCAATGGCTTCAGCAGACAGTACCCTGCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCC
69▶ 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 I I E G T K T R

4001 TGATGGCCGCCCGACATGGTGTCTGTTGCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCTGCTGAGAGATGTTGAA
36▶ I A A G V H H K 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

BspHI (4105) AseI (4163)

4101 GGTCTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGC T
3▶ T K M

SacI (4220)

4201 TATCTGACGGTTCACTAACGAGCTCTGTTATATAGACCTCCACCCTACACGCCTACCGCCATTGCGTCAATGGGGCGGAGTTGTTACGACATTTT

SpeI (4318)

4301 GGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCCA

SnaBI (4446)

4400 TTGATGTAAGTCCCAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCCAAGTGGAAAGTCCCATAGGTCATGTAAGTGGGCAT

NdeI (4551)

4500 AATGCCAGGGGGCCATTTACCGTCATTGACGTCAATAGGGGGTACTTGGCATATGATACACTTGATGTAAGTGGGCGAGTTTACCGTAAATA
4600 CTCCACCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATATTGACGTCAATGGGCGGGGCTGTTGGGCGGTACGCCA

PstI (4730) SdaI (4729) PacI (4737)

4700 GCGGGCCATTTACCGTAAGTTATGTAACGCCCTG C A G G T T A A T T A A G A A C A T G T G A G C A A A G G C C A G A A A G G C C A G A A C C G T A A A A G G C C G C G
4798 TTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACC
4898 AAGCGTTTTCCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCT

ApaLI (5061)

4998 TTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCC
5098 TTATCCGGTAACTATCGTCTTGAAGTGGTGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAA
5198 GCGGTGTACAGAGTTCTTGAAGTGGTGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAA
5298 AAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAA

PacI (5477) NotI (5496)

5398 GAAGATCCTTTGATCTTTTCTACGGGCTCTGACGCTCAGTGGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAAATTAACATTTAAATC AG

EagI (5497)

5498 CGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAA
5598 AACAAACTAGCAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA