



PvuI (7)
1 GGATCTGCATCGCTCCGGTGCCTGTCAGTGGGCGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGTGCCTA
101 GAGAAGGTGGCGGGGTAACCTGGGAAAGTGTGCTGTACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

SgfI (6) **MfeI (82)**

HindIII (245)
201 GTGAACGTTCTTTTTGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCAGAGGGGCTCGCATCTCTCTTTCACGGCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGGTTGAGTCCGGTTCTGCCGCTCCCGCTTGGTGGCTCTGAAGTGGTCCGCCGCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGGCTGACCCTGCTTGTCAACTCTACGCTTTTGTTCGTTT

KasI (535) **NgoMIV (557)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACTCTGAGATCACCGGGCCGCCACCATTGGAAATCAGCTTGTGGCATTGCTCCCTCC
1▶ M G N Q L A G I A P S

Scal (666) **XhoI (688)**
601 CAGATCCTTTCTGTAGAGAGTTATTTTTCAGATATTCATGACTTTGAATATGATAAAAGCCTGGGGAGTACTCGGTTTTTAAAGTTGCTCGAGCCAAGC
12▶ Q I L S V E S Y F S D I H D F E Y D K S L G S T R F F K V A R A K

StuI (707) **BamHI (740)**
701 ACCGAGAAGCGCTGGTCTGTGAAGGTTTTGCAATTCAGGATCCCATGCTTTAACACAGCTATAAACAAGAGCTGGAGGAACTGAAAATCAGGCT
45▶ H R E G L V V V K V F A I Q D P T L P L T S Y K Q E L E E L K I R L
801 TAATTCTGCACAGAAATTGCTACCTTTCCAGAAAGCATCAGAAAAAGCATCTGAGAAAGCAGCTATGCTCTTTAGGCAGTATGTGCGAGACAATCTCTAT
78▶ N S A Q N C L P F Q K A S E K A S E K A A M L F R Q Y V R D N L Y

Eco4VII (940)
901 GATCGCATCAGTACCCTGCTTCTGAATAACATTGAGAAGGCTGGATTGCTTTCCAGATCCTGACAGCTGTGGACCAAGCACACAAATCTGGAGTTC
112▶ D R I S T R P F L N N I E K R W I A F Q I L T A V D Q A H K S G V

BstEII (1031)
1001 GTCATGGGACATCAAGACTGAGAATGTGATCACCAGTTGGAATTGGGTTCTTCTAACTGATTTTCCAGTTTTAAGCCCACTTATCTTCCAGAAGA
145▶ R H G D I K T E N V M V T S W N W V L L T D F A S F K P T Y L P E D
1101 CAACCCGGCAGATTTCAATTATTTCTTTGACACATCACGGAGGAGAATTGCTATATTGCTCTGAACGTTTTGTTGATGGTGGGATGTTTCCACTGAG
178▶ N P A D F N Y F F D T S R R R T C Y I A P E R F V D G G M F A T E
1201 TTAGAATATATGAGAGATCCTTCAACTCCGCTTGTAGACTTAAATAGCAATCAGAGAACAAGAGGAGAGTTGAAGAGAGCAATGGACATCTTTTCAGCAG
212▶ L E Y M R D P S T P L V D L N S N Q R T R G E L K R A M D I F S A
1301 GTTGTGTGATAGCTGAGCTTTTACAGAAGGTGACCATTATTTGATCTCTCAACTTTTGGCTTATAGAAATGGACATTTTTTCCCTGAACAAGTGCCT
245▶ G C V I A E L F T E G V P L F D L S Q L L A Y R N G H F F P E Q V L
1401 AAATAAAATGAAGATCACAGTATCAGAGAATTGTAACCTCAGATGATTCACCGTGGCCAGATAAAGCTTTAGAGGCAGAAGATTACTTAAAAACAGCAG
278▶ N K I E D H S I R E L V T Q M I H R E P D K R L E A E D Y L K Q Q
1501 CGTGGCAATGCCTTCTGAAATATTTTACACTTTTCCAGCCCTACATGGCCCAAGTTGCAAGGAAACGTTTCTTCTGCAGATGAGCGTATTCTG
312▶ R G N A F P E I F Y T F L Q C P Y M A Q T F A K E T F L S A D E R I L
1601 TTATACGGAAGGATTTGGGCAACATTATTCACAATCTCTGTGGACATGATCTGCCAGAAAAAGCCGAAGGAGAGCCTAAGGAAAAATGGCTGTTATCTT
345▶ V I R K D L G N I I H N L C G H D L P E K A E G E P K E N G L V I L
1701 GGTATCTGTTATAACATCTGCCTACAGACCCTTAAATACTGTGATTCAAAACAGCTGCTTTGGAACGATTCTTCAATTTGGCTCCAAGATTAAGTGT
378▶ V S V I T S C L Q T L K Y C D S K L A A L E L I L H L A P R L S V

AvrII (1857)
1801 GAAATCCTTTGGATCGTATTACTCCATATCTTTGCATTTCCAGCAATGACTCTGTTCTAGGGTGGGGCTGAAGCCTTGAGGACGTTGACCAAAGTTC
412▶ E I L L D R I T P Y L L H F S N D S V P R V R A E A L R T L T K V

EcoRV (1929) **BspEI (1943)**
1901 TTGCTCTCGTCAAAGAGTTCTCGTAATGATCAATATTTTCCGGAATACATTTCTGCCAGGCATAGCCCACTTAGCCCAAGATGATGCTACTATCGT
445▶ L A L V K E V P R N D I N I Y P E Y I L P G I A H L A Q D D A T I V
2001 TAGACTAGCCTATGCTGAAAACATAGCTCTGCTGGCAGAAACAGCTCTGAGATTCTGGAATTAGTACAGTAAAAAACTTAATATGAAAAATGACCCC
478▶ R L A Y A E N I A L L A E T A L R F L E L V Q L K N L N M E N D P
2101 AATAATGAAGAAATAGATGAGGTTACACATCCAATGGAATTAACACAGAGCTCCAAGCCTTACATGAAATGGTCCAGCAGAAAGTTGTTACTTTGC
512▶ N N E E I D E V T H P N G N Y D T E L Q A L H E M V Q K V V T L
2201 TAAGTGACCTGAAAATATTGTAACAAACCTTGATGAAAAATGGAATAACACGCGTGTGTGATTCTTTGGACGTCAGAAAGCCAACGATGTTTTGTT
545▶ L S D P E N I V K Q T L M E N G I T R L C V F F G R Q K A N D V L L
2301 GTCCACATGATTACTTTCTAAATGATAAGAATGATTGGCATCTACGTGGAGCATTTTTTGATAGTATAGTTGGTGTGCTGCCTATGTTGGCTGGCAA
578▶ S H M I T F L N D K N D W H L R G A F F D S I V G V A A Y V G W Q
2401 AGCTCCTCAATTCAAGCCTCTGCTGCAACAAGGTCTTAGTGATGCTGAGGAAATTTGTCATTGTGAAAGCTTTTATGCCCTTACTTGTATGTGCGAGT
612▶ S S S I L K P L L Q Q G L S D A E E F V I V K A L Y A L T C M C Q
2501 TAGGACTGCTACAAAACCCCATGTTTACGAATTTGCCAGTGATATTGCCCTTCTGTGTCATCCCAATTTATGGATACGTTATGGTCCGCTGGGATT
645▶ L G L L Q K P H V Y E F A S D I A P F L C H P N L W I R Y G A V G F
2601 TATCACAGTGGTACGCTGTCAAAATAGTACAGCTGATGCTACTGTAACCTGATGCCTTATCTTGACCCATATATTCCCAACCAATAACAGATGAA
678▶ I T V V A R Q I S T A D V Y C A K L M P Y L D P Y I T Q P I I Q I E
2701 AGAAAACTGTTCTGCTCAGTGTTTTAAAGGAACAGTAAGTCGTTCTATATTTGATTATGCTTTGAGGTCTAAAGATATTACTAGCTTGTTCAGACATC
712▶ R K L V L L S V L K E P V S R S I F D Y A L R S K D I T S L F R H

BamHI (2855)
2801 TTCACATGCGTCAGAAGAAACGAAATGGTTCTCTTCCGACTGCCTCCGCCAGAGGATCCTGCCATAGCACAGCTTCTGAAGAAGTTGCTCTCACAGGG
745▶ L H M R Q K K R N G S L P D C P P P E D P A I A Q L L K K L L S Q G
2901 AATGACAGAGGAAGAGGAAGACAACTTCTGGCACTGAAAGACTTTCATGATGAAATCTAATAAAGCAAAGGCCAATATAGTGACAGAGCCATCTTCAT
778▶ M T E E E E D K L L A L K D F M M K S N K A K A N I V D Q S H L H
3001 GATAGTAGTCAGAAAGGTGTAATTGACTTGGCAGCTTTAGGCATAACTGGGAGACAAGTTGATCTTGTAAAACCAAACAAGAACCAGATGACAAACGGG
812▶ D S S S Q K G V I D L A A L G I T G R Q V D L V K T K Q E P D D K R
3101 CCAGAAAACATGAAAACAGACTCAAATGAAATGAAGAATGAAAAGCATGTTGGTCTACTGGACCCAAACATGCCACAGCCCTACTTAAAGG
845▶ A R K H V K Q D S N V N E E W K S M F G S L D P P N M P Q A L P K G

PshAI
3201 GAGTGATCAGGAGGTGATTACAGACTGGGAAACCTCCTCGTTCGAGTCTCTGCTGGCATTGTTGTCCTTTGTCAACTTCTTACAGGTTCCAGAAGTG
878▶ S D Q E V I Q T G K P P R S E S S A G I C V P L S T S S Q V P E V

3301 **AgeI (3329)** ACAACTGTCCAAAATAAAAAACAGTAATACCGGTTTTAAGTAGTACAATCTTACCATCCACCTATCAGATTCGAATTACAACCTGTAACCTGAACTTC
912▶ T T V Q N K K P V I P V L S S T I L P S T Y Q I R I T T C K T E L **XcmI (3489)**

3401 AGCAACTCATCCAGCAAAGCGGGAGCAGTGCAATGCTGAGAGAATAGCTAAGCAGATGATGGAAAATGCTGAATGGGAGAGTAAACCACCACCACCTGG
945▶ Q Q L I Q Q K R E Q C N A E R I A K Q M M E N A E W E S K P P P P G

3501 ATGGCGTCTAAAGGGCTGTTAGTTGCCATCTTCATGAGCATAAATCTGCTGTGAATCGAATTAGAGTCTCTGATGAACACTCACTTTTTGCAACATGT
978▶ W R P K G L L V A H L H E H K S A V N R I R V S D E H S L F A T C **BglIII (3663)**

3601 TCAATGATGGCACAGTGAAGTCTGGAACAGTCAAAAGATGGAGGGGAGACCACCCTACCAGATCTATTCTTACATACAGCCGAATTGGAGGACGAG
1012▶ S N D G T V K I W N S Q K M E G K T T T T R S I L T Y S R I G G R

3701 TCAAGACGCTCACATTCTGCAAGGCTCCCACTATTTAGCCATAGCATCTGATAATGGTGTCTGCCAGCTTCTTGAATTGAGGCTTCTAAGCTGCCAA
1045▶ V K T L T F C Q G S H Y L A I A S D N G A V Q L L G I E A S K L P K **XbaI (3833)**

3801 **EcoRI (3829)** GTCTCTAAAATCCATCCTCTACAAAGCAGAATTCTAGATCAGAAGGAGGACGGTTGTGTGGATATGCATCACTTCAACTCTGGAGCACAGTCTGTT
1078▶ S P K I H P L Q S R I L D Q K E D G C V V D M H H F N S G A Q S V **NsiI (3867)**

3901 CTTGCCTATGCCACTGTGAATGGCTCTCTGGTTGGCTGGGACCTTAGGTCTTCAAGCAATGCGTGGACTTTAAAGCATGATTTAAAGTCGGGCCTCATCA
1112▶ L A Y A T V N G S L V G W D L R S S S N A W T L K H D L K S G L I **Acc65I (4053)**

4001 CTTCTTTGCTGTGGACATCCACCAATGCTGGCTCTGCATTGGTACAAGCAGTGGTACCATGGCTTGTGGGACATGAGGTTCCAGTTGCCAATTTCAAG
1145▶ T S F A V D I H Q C W L C I G T S S G T M A C W D M R F Q L P I S S **XcmI (4198)**

4101 TCACTGTATCCTTCCAGGGCTCGAATCAGACGCTCTCAATGCACCCTCTGTATCAGTCTGGGTGATTGCAGCTGTTCCAGGCAACAACGAAGTGTCC
1178▶ H C H P S R A R I R R L S M H P L Y Q S W V I A A V Q G N N E V S **ApaLI (4249)**

4201 ATGTGGGACATGGAGACTGGTGACAGAAGATTTACTCTCTGGGCCAGCAGTGCACCACCTTCTGAATTACAGCCTTCTCCTCATAGCGTCCATGGTA
1212▶ M W D M E T G D R R F T L W A S S A P P L S E L Q P S P H S V H G

4301 TCTACTGTAGTCCGAGATGAAAATCCCTACTAACAGCTGGCTCAGATATGAAAATAAGGTTTTGGGACTTGGCTTACCAGAAAGGTCCTATGT
1245▶ I Y C S P A D G N P I L L T A G S D M K I R F W D L A Y P E R S Y V **SpeI (4413)**

4401 TGTTCAGGAAGTACTAGTTCCCCATCTGTGTCCTACTACAGGAAAATAATTGAAGGCACTGAAGTTGTCAGGAAATTCAGAATAAGCAGAAAGTAGGA
1278▶ V A G S T S S P S V S Y Y R K I I E G T E V V Q E I Q N K Q K V G **ScaI (4410)**

4501 CCAAGTGTGACACCCCTCGAAGGGGCCAGAGTCCCTGCCCGTGGGACATCATGACATCACTGATGTCGCCACATTCCAGACCACACAGGGCTTCA
1312▶ P S D D T P R R G P E S L P V G H H D I I T D V A T F Q T T Q G F **Bsp120I (4523)**

4601 TCGTAACTGCTTCTAGAGATGGGATTGTGAAGGTGTGAAATAAAACCTGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACA
1345▶ I V T A S R D G I V K V W K • **MscI (4655)**

4701 **NheI (4649)** ACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAAACA
HpaI (4787) MfeI (4798)

4801 **EcoRI (4883)** ATTGCAATCATTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTCTAAAATACAGC
4901 ATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGC
5001 TGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAAGATATAGTGTATTTTCCAAAGTTTGAACATAGCTCTTCAATTTCTTTATGTTTTAAATGCACTGACC

5101 **SwaI (5136)** TCCCATTCCCTTTTAGTAAAATATTAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAAATCCAGATGCTCAAGG
5201 CCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCC
1411▶ • N R

5301 TGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAAGGAGCATAGTCAGAGATGAG
138▶ T Y 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

5401 CTCTCTGCACATGCCACAGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTCCCTGACAGCCACAATGGTGTCAAAGTCCCTCTGC
105▶ E R 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 **StuI (5561)**

5501 CCGTTGCTCACAGCAGACCAATGGCAATGGCTTCCAGCAGACAGTACCCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAGTCT
71▶ G N 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

5601 TGGTCTGATGGCCGCCGACATGGTCTTGTGCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCTGCTGAGAGAT
38▶ T R 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 **AseI (5769)**

5701 **AseI (5769)** GTTGAAGTCTTCTCATGATGGCCCTCTATAGTGAAGTCTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCT
5▶ N F T K M

5801 CCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCTACCGCCATTTGCGTCAATGGGCGGAGTTGTTACGA

5901 **SpeI (5924)** CATTGTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCCATGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCAC

6001 **SnaBI (6052)** GCCCATTGATGACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCCATAGGTCATGACTG

6101 GGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACCTTGATGTACTGCCAAGTGGGCAGTTACCGT
6201 AAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTC

PacI (6343)
6301 AGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCC
6401 GCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGAT
6501 ACCAGGCGTTTTCCCTGGAAAGCTCCCTCGTGCGCTCTCTGTTCCGACCCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGC

ApaLI (6667)
6601 GCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCCCGACCGCTGC
6701 GCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATG
6801 TAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGG
6901 AAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCT

PacI (7083) SmaI (7092)
7001 CAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAAT

EagI (7103)
NotI (7102)
7101 CAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGGTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAA
7201 CAAAACAACTAGCAAAATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTATCGAA