



**PvuI (7)**  
**SgfI (6)**  
1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACC GTATATAAGTGCAGTAGTCGCC

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

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

**NgoMIV (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGCTCAACTCTACGCTTTTGTTCGTTT

**AgeI (552)** **BspHI (560)**  
501 TCTGTTCTGCGCGCTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCATCATGAACAGCAGCAGCGCAACATCACCTACGCCAGTCG

601 CAAGCGCGGAAGCCGGTGCAGAAAACAGTAAAGCCAATCCCAGTGAAGGAATCAAGTCAAATCCTTCAAGCGGCATAGAGACCGACTTAATACAGAG  
13▶ K R R K P V Q K T V K P I P A E G I K S N P S K R H R D R L N T E

**NheI (712)** **AseI (742)**  
701 TTGACCGTTTGGCTAGCCTGCTGCCTTCCCAACAAGTGTATTAATAAGTTGGACAAACTTTCAGTTCTTAGGCTCAGCGTCAGTTACCTGAGAGCCA

47▶ L D R L A S L L P F P Q D V I N K L D K L S V L R L S V S Y L R A

**StuI (885)**  
801 AGAGCTTCTTTGATGTTGCATTAATACTCCCTACTGAAAGAAACGGAGGCCAGGATAACTGTAGAGCAGCAAATTCAGAGAAGGCCGTAACCTACA

80▶ K S F F D V A L K S S P T E R N G G Q D N C R A A N F R E G L N L Q

**EcoRI (907)**  
901 AGAAGGAGAATTCTTATTACAGGCTCTGAATGGCTTTGTATTAGTTGTCACTACAGATGCTTTGGTCTTTTATGCTTCTTCTACTATACAAGATTATCTA

113▶ E G E F L L Q A L N G F V L V V T T D A L V F Y A S S T I Q D Y L

1001 GGGTTTCAGCAGTCTGATGTCATACATCAGAGTGTATATGAACCTATCCATACCGAAGACCGAGCTGAATTCAGCGTCAGCTACCTGGGCATTAATC

147▶ G F Q Q S D V I H Q S V Y E L I H T E D R A E F Q R Q L H W A L N

1101 CTTCTCAGTGTACAGAGTCTGGACAAGGAATTGAAGAAGCCACTGGTCTCCCCAGACAGTAGTCTGTTATAACCCAGACCAGATTCTCCAGAAAAC

180▶ P S Q C T E S G Q G I E E A T G L P Q T V V C Y N P D Q I P P E N S

**NdeI (1221)**  
1201 TCCTTAAATGGAGAGGTGCTTCATATGTCGTCTAAGGTGTCTGCTGGATAATTCATCTGGTTTTCTGGCAATGAATTTCCAAGGGAAGTTAAAGTATCTT

213▶ P L M E R C F I C R L R C L L D N S S G F L A M N F Q G K L K Y L

1301 CATGGACAGAAAAAGAAAGGAAAGATGGATCAATACTCCACCTCAGTTGGCTTTGTTTGGCATAGTACTCCACTTCAGCCACCATCCATACTTAAAA

247▶ H G Q K K K G K D G S I L P P Q L A L F A I A T P L Q P P S I L E

**BspEI (1400)**  
1401 TCCGGACCAAAAATTTTATCTTTAGAACCAACACAACTAGACTTCACACCTATTGGTTGTGATGCCAAAGGAAGAATTGTTTTAGGATATACTGAAGC

280▶ I R T K N F I F R T K H K L D F T P I G C D A K G R I V L G Y T E A

**ApaLI (1506)**  
1501 AGAGCTGTGCACGAGAGGCTCAGGTTATCAGTTTATTCATGCAGCTGATATGCTTTATTGTGCCGAGTCCCATATCCGAATGATTAAGACTGGAGAAAAGT

313▶ E L C T R G S G Y Q F I H A A D M L Y C A E S H I R M I K T G E S

1601 GGCATGATGTTTTCCGGCTTTCACAAAAACAACCGACTGGGTCAGTCTAATGCACGCCTGCTTTATAAAAAAGAACGACGATTATATCA

347▶ G M I V F R L L T K N N R W T W V Q S N A R L L Y K N G R P D Y I

1701 TTGTAACCTCAGAGACCACTAACAGATGAGGAAGGAACAGAGCATTTACGAAAACGAAATACGAAGTTGCCTTTTATGTTTACCACTGGAGAAGCTGTGTT

380▶ I V T Q R P L T D E E G T E H L R K R N T K L P F M F T T G E A V L

**BamHI (1833)** **SpeI (1864)**  
1801 GTATGAGCAACCAACCTTTTCTGCCATAATGGATCCCTTACCCTAAGGACTAAAAATGGCACTAGTGGAAAAGACTCTGCTACCACATCCACTCTA

413▶ Y E A T N P F P A I M D P L P L R T K N G T S G K D S A T T S T L

1901 AGCAAGGACTCTCTCAATCCTAGTCCCTCCTGGCTGCCATGATGCAACAAGATGAGTCTATTTATCTCTATCCTGCTCAAGTACTTCAAGTACTGCAC

447▶ S K D S L N P S S L L A A M M Q Q D E S I Y L Y P A S S T S S T A

2001 CTTTTGAAAACAACCTTTTCAACGAATCTATGAATGAATGCAGAAAATGGCAAGATAATACTGCACCGATGGGAAATGATACTATCCTGAAACATGAGCA

480▶ P F E N N F F N E S M N E C R N W Q D N T A P M G N D T I L K H E Q

**Bsu36I (2110)** **BstEII (2137)** **AvrII (2196)**  
2101 AATTGACCAGCCTCAGGATGGAACCTATTTGCTGGAGGTCACCCAGGGCTCTTTCAAGATAGTAAAAACAGTGACTTGACAGCATAATGAAAAACCTA

513▶ I D Q P Q D V N S F A G G H P G L F Q D S K N S D L Y S I M K N L

2201 GGCATTGATTTTGAAGACTCAGACACATGCAGAATGAAAAATTTTCAGAAATGATTTTTCTGGTGGGTTGACTTCAGAGCATTGACTTAACGGATG

547▶ G I D F E D I R H M Q N E K F R N D F S G E V D F R D I D L T D

2301 AAATCCTGACGTATGTCCAAGATTCTTTAAGTAAGTCTCCCTTCATACCTTACAGATTCAACAGCAACAGTCTTGGCTCTGAACTCAAGCTGTATGGT

580▶ E I L T Y V Q D S L S K S P F I P S D Y Q Q Q Q S L A L N S S C M V

**XbaI (2415)**  
2401 ACAGGAACACCTACATCTAGAACAGCAACAGCAACATCACAAAAGCAAGTAGTAGTGGAGCCACAGCAACAGCTGTGTCAGAAGATGAAGCACATGCAA

613▶ Q E H L H L E Q Q Q Q H H Q K Q V V V E P Q Q Q L C Q K M K H M Q

2501 GTTAATGGCATGTTTAAAAATTTGAACTTAACCAATTCGTGCTTTCAATTTGCCACAGCAAGACCCACAACAATATGCTTTACAGACTTACATG

647▶ V N G M F E D I R W N S N Q F V P F N C P Q Q Q D P Q Q Y N V F T D L H

2601 GGATCAGTCAAGAGTTCCCTACAAATCTGAAATGATTCTATGCCTTATACACAGAACTTTATTTCTGTAATCAGCCTGTATTACCACAACATTCCAA

680▶ G I S Q E F P Y K S E M D S M P Y T Q N F I S C N Q P V L P Q H S K

2701 ATGTACAGAGCTGGACTACCCTATGGGGAGTTTGAACCATCCCCATACCCCACTACTTCTAGTTTGAAGATTTTGTCACTTGTTTACAACCTCCTGAA  
713▶ C T E L D Y P M G S F E P S P Y P T T S S L E D F V T C L Q L P E  
BstXI (2802) BspLU11I (2848)  
2801 AACCAAAAGCATGGATTAATCCACAGTCAGCCATAAATACTCCTCAGACATGTTATGCTGGGGCCGTGTCGATGTATCAGTGCCAGCCAGAACCTCAGC  
747▶ N Q K H G L N P Q S A I I T P Q T C Y A G A V S M Y Q C Q P E P Q  
BbrPI (2905)  
2901 ACACCCACGTGGGTGAGTGCAGTACAATCCAGTACTGCCAGGCCAACAGGCATTTTTAAACAAGTTTCAAGTGGAGTTTTAAATGAAACATATCCAGC  
780▶ H T H V G Q M Q Y N P V L P G Q Q A F L N K F Q N G V L N E T Y P A  
3001 TGAATTAATAACATAAATAACACTCAGACTACCACACATCTTCAGCCACTTCATCATCCGTGAGAAGCCAGACCTTTTCTGATTTGACATCCAGTGA  
813▶ E L N N I N N T Q T T T H L Q P L H H P S E A R P F P D L T S S G  
MluI (3132)  
3101 TTCCTGTAATCCAAGCCAATTTTACCCCTGACGCGTTCTAGCTCGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGA  
847▶ F L •  
3201 AAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACA  
HpaI (3301) EcoRI (3397)  
3301 AGTTAAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAA  
3401 TTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGC  
SapI (3579)  
3501 CAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTTATGAGTAAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTGCTCTTCATTTCTTTATGTTT  
SspI (3636) SmaI (3650)  
3601 TAAATGCACTGACCTCCACATTCCTTTTTAGTAAAAATTCAGAAAAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAAT  
EcoO109I (3711)  
3701 CCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTT  
3801 CTAGCTTTAGTTCTGGTGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCA  
141▶ N R 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  
SacI (3911) BstXI (3940)  
3901 TAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGT  
109▶ Y D S I L 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  
StuI (4075)  
4001 CAAAGTCCTTCTGCCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCCTCAATGTGGACAGCAGAGAT  
76▶ F D K Q 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  
4101 GATCTCCCAGTCTTGGTCTGATGGCCGCCCCGACATGGTGTCTGTGCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGA  
43▶ I E G T K 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  
BspHI (4225)  
XmnI (4217) AseI (4283)  
4201 TCCTGCTGAGAGATGTTGAAGTCTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAG  
9▶ D Q Q S I N F T K M  
SacI (4340)  
4301 CGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGG  
SpeI (4438)  
4401 CGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTC  
SnaBI (4566)  
4501 AAACCGCTATCCACGCCATTGATGTACTGCCAAAACGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCCAT  
NdeI (4671)  
4601 AAGGTCATGTAAGTGGCATAATGCCAGGCGGGCATTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCTCAAGT  
4701 GGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTACTATGGAAACATACGTCAATTATTGACGTCAATGGGCGGGG  
PacI (4857)  
PstI (4850)  
SdaI (4849) BspLU11I (4867)  
4801 GTCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGACGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAGGCCAGGA  
4901 ACCGTA AAAAGGCCGCTTGTGGCGTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGAC  
5001 AGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGGCTCTCTGTCCGACCTGCGGCTTACCGGATACCTGTCCGCTTTCTCCCT  
ApaLI (5181)  
5101 TCGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTC

5201 AGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAG

5301 CAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAG

5401 CCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGGTTTTTTTGGTTGCAAGCAGCAGATTACGCGCA

5501 GAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTA PacI (5597)

**EagI (5617)**

SwaI (5606) **NotI (5616)**

5601 ATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGGTTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATC

5701 AAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA