



**PvuI (7)**  
**SgfI (6)** 1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA **EcoNI (96)**

101 GAGAAAGTGGCGCGGGTAAACTGGAAAAGTATGTCGTGTACTGGCTCCGCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

**HindIII (245)** **Bsu36I (291)**  
**Psp1406I (203)** **PvuII (239)** **EcoNI (287)**  
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTCACGCGCCCGCCGCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

**NgoMIV (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTGCCTGACCCTGTTGCTCAACTCTACGCTTTTGTTCGTTT

**AgeI (552)** **NcoI (568)**  
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCGCTACCTGAGATCACCGGTAGGAGGGCCACCATGGGCTGTGACCGGAACTGTGGGCTCATC  
1► M G C D R N C G L I

**BbsI (674)**  
601 GCTGGGGCTGTCATTGGTGTCTCTGGCTGTGTTGGAGGTATTCTAATGCCAGTTGAGACCTGCTTATCCAGAAGACAATAAAAAGCAAGTTGTCC  
11► A G A V I G A V L A V F G G I L M P V G D L L I Q K T I K K Q V V

701 TCGAAGAAGTACAATTGCTTTTAAAAATTGGGTTAAACAGGCACAGAAGTTTACAGACAGTTTTGGATCTTTGATGTGCAAAATCCACAGGAAGTAT  
44► L E E G T I A F K N W V K T G T E V Y R Q F W I F D V Q N P Q E V M

**SnaBI (842)**  
801 GATGAACAGCAGCAACATTCAAGTTAAGCAAAGAGGTCCTTACGTACAGAGTTCGTTTTCTAGCCAAGGAAATGTAACCCAGGACGCTGAGGACAAC  
77► M N S S N I Q V K Q R G P Y T Y R V R F L A K E N V T Q D A E D N

**BstBI (933)** **Bsp119I (933)** **PvuII (997)**  
901 ACAGTCTCTTTCCTGCGAGCCCAATGGTGCCATCTTCGAACTTCACTATCAGTTGGAACAGAGGCTGACAACCTCACAGTTCTCAATCTGGCTGTGGCAG  
111► T V S F L Q P N G A I F E P S L S V G T E A D N F T V L N L A V A

1001 CTGCATCCCATATCTATCAAAATCAATTTGTTCAAATGATCCTCAATCATTATTAACAAGTCAAAATCTTCTATGTTCCAAGTCAGAACTTTGAGAGA  
144► A A S H I Y Q N Q F V Q M I L N S L I N K S K S S M F Q V R T L R E

**BamHI (1118)**  
1101 ACTGTTATGGGGCTATAGGGATCCATTTTTGAGTTTGGTTCGTAACCTGTTACTACCACAGTTGGTCTGTTTTATCCTTACAACAATACTGCAGATGGA  
177► L L W G Y R D P F L S L V P Y P V T T T V G L F Y P Y N N T A D G

1201 GTTTATAAAGTTTTCAATGGAAAAGATAACATAAGTAAAGTTGCCATAATCGACACATATAAAGGTAAGGAATCTGTCCTATTGGGAAAGTCACTGGC  
211► V Y K V F N G K D N I S K V A I I D T Y K G K R N L S Y W E S H C

**AseI (1305)**  
1301 ACATGATTAATGGTACAGATGCAGCCTCATTCCACCTTTTGTGAGAAAAGCCAGGTATTGCAGTCTTTTTCTTCTGATATTGCAGGTCAATCTATGC  
244► D M I N G T D A A S F P P F V E K S Q V L Q F F S S D I C R S I Y A

**StuI (1462)**  
1401 TGTATTTGAATCCGACGTTAATCTGAAAGGAATCCCTGTGTATAGATTTGTTCTTCCATCCAAGGCCCTTGCCTCTCCAGTTGAAAACCCAGACAATAT  
277► V F E S D V N L K G I P V Y R F V L P S K A F A S P V E N P D N Y

**NdeI (1540)**  
1501 TGTTTCTGCACAGAAAAAATTATCTCAAAAAATTGTACATCATATGGTGTGCTAGACATCAGCAAATGCAAAGAAGGGAGACCTGTGTACATTTCACTTC  
311► C F C T E K I I S K N C T S Y G V L D I S K C K E G R P V Y I S L

1601 CTCATTTCTGTATGCAAGTCCCTGATGTTTCAAGAACCTATTGATGGATTAACCCAAATGAAGAAGACATAGGACATACTGGATATTGAACCTATAAC  
344► P H F L Y A S P D V S E P I D G L N P N E E H R T Y L D I E P I T

**XmnI (1758)**  
1701 TGGATTCACCTTACAATTTGCAAAACGGCTGCAGGTCAACCTATTGGTCAAGCCATCAGAAAAAATCAAGTATTAAGAATCTGAAGAGGAATATATT  
377► G F T L Q F A K R L Q V N L L V K P S E K I Q V L K N L K R N Y I

**BspLU11I (1852)**  
1801 GTGCCTATTCTTTGGCTTAATGAGACTGGGACCATTGGTGTGAGAAGGCAACATGTTTCAAGTCAAGTAACTGGAAAAATAAACCTCCTTGGCCTGA  
411► V P I L W L N E T G T I G D E K A N M F R S Q V T G K I N L L G L

**MscI (1995)**  
1901 TAGAAATGATCTTACTCAGTGTGGTGTGGTGTGTTGTTGCTTTTATGATTTTCAATTTGTGCATGCAGATGAAAAACAATAAAATAAGCTAGCTGGCC  
444► I E M I L L S V G V V M F V A F M I S Y C A C R S K T I K • —

2001 AGACATGATAAGATACATTGATGAGTTTGGACAACCACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTT

**HpaI (2127)**  
2101 GTAACATTATAAGTGAATAAACAAGTTAAACAACAATTCGATTCATTTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGT

**EcoRI (2223)**  
2201 AAAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAAT

2301 AAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTCTTTCATGGAGTTTAAAGATATAGTGATTTTCCCAAGGTTTG

2401 AACTAGCTCTTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCACATTCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAA

2501 AATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAG

2601 AAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCT  
 141 • N R T Y K L P I L E E I T T K V L K G N M E

2701 CAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGTCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTA  
 118 I L V F C D P A 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

2801 GGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATG  
 85 P H R V A V I T D F D K Q G N S V A S G I A I A E A C V T V R G I

2901 TAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGTCTGTTGCTCATAGAGCATGGTGTCTTCT  
 51 Y A E I H V A S I 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

3001 CAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACT  
 18 T A V E V L E L D Q Q S I N F T K M

3101 ATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCCTACACG

3201 CCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAACAAACTCCCATTGACGTCAATGGGG

3301 TGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGA

3401 TGTACTGCCAAGTAGGAAAGTCCATAAGGTATGTAAGTGGGCATAATGCCAGGCGGGCCATTTACCCTGATTGACGTCAATAGGGGGCTACTTGGCAT

3501 ATGATACACTTGTACTGCAAGTGGGCAGTTTACCCTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATAC

3601 GTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCCTAAGTTATGTAACGCTGCAGGTTAAITAAAGAACATGTG  
 SdaI (3675) PacI (3683) BspLU11I (3693)

3701 AGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGC

3801 TCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGCTTA

3901 CCGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCT

4001 GGGCTGTGTGCACGAACCCCGTTAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGGTAAGACACGACTTATCGCCACTG

4101 GCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAG

4201 TATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGTTTTTT

4301 TGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAACTCACGT

4401 TAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAAATATCTTTATTTTTCATTACATCTGTGTGTTGTTTTTTGTGTGA  
 EagI (4443) PacI (4423) SmaI (4432) NotI (4442)

4501 ATCGTAACTAACATACGCTCTCCATCAAACAAACGAAACAAACAAACTAGCAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTC

4601 TATCGAA