



**PvuI (7)**  
**Sgfl (6)** **EcoNI (96)**  
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGTCCGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGGGGTAACCTGGGAAAGTATGTCGTGTACTGGCTCCGCCCTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

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

301 GCCATCCAGCGGGTTGAGTCGCGTTCGCCCGCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCGCCGCTAGGTAAGTTTAAAGTCCAGGTCGAGACC

**NgoMI (441)**  
**NaeI (441)**  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGAGCTACCTAGACTCAGCCGGCTTCCACGCTTTGCTGACCTGCTTGTCAACTCTACGCTTTGTTTCGTTT

**SphI (560)**  
**AgeI (552)**  
501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGGCGCTACCTGAGATCACCGTTCAGCATGCAGAGCACTTCTAATCATCTGTGGCTTTTATCTGA  
1 M Q S T S N H L W L L S D

**HindIII (690)**  
601 TATTTTAGCCAAGGAGCTACTGCAATGTCTTTCGTGGAAGACATAAGAAAAGTGGTATTTATTTGCTATCAAAGTATTTAATAACATAAGCTTCTCT  
13 I L G Q G A T A N V F R G R H K K T G D L F A I K V F N N I S F L  
701 CGTCCAGTGGATGTTCAAATGAGAGAATTTGAAGTGTGAAAAAAGTCAATCACAAAATATTGTCAAATTTTGTCTATTGAAGAGGAGACAACAACA  
47 R P V D V Q M R E F E V L K K L N H K N I V K L F A I E E E T T T

**ScaI (807)** **EcoRI (892)**  
801 GACATAAAGTACTTATTATGGAATTTTGTCCATGTGGGAGTTTATACACTGTTTAGAAGAACCTTCTAATGCCTATGGACTACCAGAATCTGAATTCCT  
80 R H K V L I M E F C P C G S L Y T V L E E P S N A Y G L P E S E F L

**EcoRV (964)**  
901 AATTGTTTTGCGAGATGTGGTGGTGGGAATGAATCATCTACGAGAGAATGGTATAGTGCACCGTGATATCAAGCCAGGAAATATCATGCGTGTATAGGG  
113 I V L R D V V G G M N H L R E N G I V H R D I K P G N I M R V I G

**BsrGI (1016)**  
1001 GAAGATGGACAGTCTGTGTACAAACTCAGAGTTTTGGTGCAGCTAGAGAATTAGAAGATGATGAGCAGTTTGTCTCTGTATGGCACAGAAGAATATT  
147 E D G Q S V Y K L T D F G A A R E L E D D E Q F V S L Y G T E E Y

**XcmI (1197)**  
1101 TGCACCCTGATATGATGAGAGAGCAGTCTAAGAAAAGATCATCAGAAGAAATATGGAGCAACAGTTGATCTTTGGAGCATTGGGGTAACATTTTACCA  
180 L H P D M Y E R A V L R K D H Q K K Y G A T V D L W S I G V T F Y H

**EcoO109I (1236)**  
1201 TGCAGTACTGGATCACTGCCATTTAGACCTTTGAAGGCTCGTAGGAATAAAGAAGTGTATATAAAAATATTACAGGAAAGCCTTCTGGTGCATA  
213 A A T G S L P F R P F E G P R R N K E V M Y K I I T G K P S G A I

**BstXI (1327)**  
1301 TCTGGAGTACAGAAAAGCAGAAAATGGACCAATTTGACTGGAGTGGAGACATGCCTGTTTCTGAGTCTTCTCGGGTCTTCTAGTTCTACTTACCCTG  
247 S G V Q K A E N G P I D W S G D M P V S C S L S R G L Q V L L T P

**SpeI (1465)**  
1401 TTCTTGCAAACATCCTTGAAGCAGATCAGGAAAAGTGTGGGGTTTTGACCAAGTTTTTTCAGAAAAGTGTATATCTTACCGAATGGTAATTCATGT  
280 V L A N I L E A D Q E K C W G F D Q F F A E T S D I L H R M V I H V

**BspHI (1566)**  
1501 TTTTTCGCTACAACAAATGACAGCTCATAAGATTTATATACATAGCTATAATACTGCTACTATATTTTCATGAACTGGTATATAAACAAACAAAATATT  
313 F S L Q Q M T A H K I Y I H S Y N T A T I F H E L V Y K Q T K I I  
1601 TCTTCAAATCAAGAACTTATCTACGAAGGGCAGCCTTGTCTTGAAGCTGGAAGGCTGGCACAACATTTCCCTAAAAGTACTGAGGAAAACCTATAT  
347 S S N Q E L I Y E G R R L V L E P G R L A Q H F P K T T E E N P I

**NheI (1795)**  
1701 TTGTAGTAAAGCGGGAACCTCTGAATACCATAGGATTAATATATGAAAAATTTCCCTCCCTAAAGTACATCCAGTTATGATTTAGACGGGGATGCTAG  
380 F V V S R E P L N T I G L I Y E K I S L P K V H P R Y D L D G D A S

**PstI (1836)**  
1801 CATGGCTAAGCAATAACAGGGTGTGTGTTATGCCTGCAGAATTTGCCAGTACTTACTGCTTTATCAGGAATTAATGCGAAAAGGGATACGATGGCTG  
413 M A K A I T G V V C Y A C R I A S T L L L Y Q E L M R K G I R W L  
1901 ATTGAATTAATTAAGATGATTACAATGAAACTGTTCAAAAAAGACAGAAGTTGTGATCACATTGGATTTCTGTATCAGAAACATTGAAAAACTGTGA  
447 I E L I K D D Y N E T V H K K T E V V I T L D F C I R N I E K T V  
2001 AAGTATATGAAAAGTTGATGAAGTCAACCTGGAAGCGCAGAGTTAGGTGAAATTTAGACATACACCAAAATTTGAGACTTTCCAGTTCTCAGGG  
480 K V Y E K L M K I N L E A A E L G E I S D I H T K L L R L S S S Q G

**EcoRV (2122)**  
2101 AACAAATAGAAACAGTCTTCCAGGATTCGACAGCAGATTATCTCCAGTGGATCACTGGCAGACGCATGGGCACATCAAGAAGGCACTATCCGAAAGAC  
513 T I E T S L Q D I D S R L S P G G S L A D A W A H Q E G T H P K D  
2201 AGAAATGTAGAAAACTACAAGTCTGTTAAATTCATGACAGAGATTTACTATCAGTTCAAAAAAGACAAAGCAGAACGTAGATTAGCTTATAATGAAG  
547 R N V E K L Q V L L N C M T E I Y Y Q F K K D K A E R R L A Y N E  
2301 AACAAATCCAAAATTTGATAAGCAAAAAGTATTACCATGCCACAAAAGCTATGACGCCTTTACAGATGAATGTGTTAAAAAGTATGAGGCAATTTTT  
580 E Q I H K F D K Q K L Y Y H A T K A M T H F T D E C V K K Y E A F L  
2401 GAATAAGTCAGAAGATGGATAAGAAAGTCTTATCTTAGGAAACAGTATTATCTGCTGACTAATCAGTGTTTGATATTGAAGAAGAAGTATCAAAA  
613 N K S E E W I R K M L H R K K L L S L T N Q C F D I E E E V S K  
2501 TATCAAGAATATACTAATGAGTTACAAGAACTCTGCCAGAAAATGTTTACAGCTTCCAGTGAATCAAAACATACCATGACCCCAATTTATCCAAGTT  
647 Y Q E Y T N E L Q E T L P Q K M F T A S S G I K H T M T P I Y P S  
2601 CTAACACATTAGTAGAAATGACTCTTGGTATGAAGAAATTAAGGAAGAGATGGAAGGGTGGTTAAAGAACTGCTGAAAATAACACATTTTGAAGAA  
680 S N T L V E M T L G M K K L K E E M E G V V K E L A E N N H I L E R

**NeoI (2714)** **Psp1406I (2734)** **BamHI (2749)** **NdeI (2789)**  
2701 GTTTGGCTCTTAAACATGGATGGTGGCCTTCGCAACGTTGACTGCTTGGATCCTATCCCTATGATGTGCCAGACTATGCTGGCTATCCATATGATGTT  
713 F G S L T M D G G L R N V D C L G S Y P Y D V P D Y A G Y P Y D V

**MscI (2852)**

NheI (2846)

2801 CCTGATTATGCTGGATACCCCTTATGATGTGCCAGACATATGCCTAAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAAC  
747 ▶ P D Y A G Y P Y D V P D Y A •

**HpaI (2984)**

2901 AGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACAAAT

**EcoRI (3080)**

3001 GCATTCAATTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAAATCTAAAATACAGCATA

3101 GCAAACTTTAACCTCCAATCAAGCTCTACTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGT

3201 TTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTCCCAAGTTTGAAGTACTCTTCATTCTTTATGTTTTAAATGCACTGACCTCC

**SwaI (3333)** **EcoO109I (3394)**

3301 CACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCC

3401 TTCATAATATCCCCAGTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTCTGG  
141 ◀ • N R T

**SacI (3594)**  
**Ecl136II (3594)**

3501 TGTAAGTGGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTC  
137 ◀ 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 E

**BstXI (3623)**

3601 TCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCG  
104 ◀ 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 G

**StuI (3758)**

3701 TTGCTCACAGCAGCCCAATGGCAATGGCTTCAGCAGACAGTACCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGG  
70 ◀ 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 T

3801 TCCTGATGGCCGCCCGACATGGTCTTGTCTCCTCATAGAGCATGGTGTCTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTT  
37 ◀ 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 N

**BspHI (3908)**

3901 GAAGTCTTCATGATGGCCCTCTATAGTGTGATGATTATACTATGCGGATATACTATGCCGATGATTAATTTGTCAAACAGCGTGGATGGCGTCTCCA  
4 ◀ F T K M

**SacI (4023)**  
**Ecl136II (4023)**

4001 GCTTATCTGACGGTTCCTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACAT

**SpeI (4121)**

4101 TTTGGAAAGTCCCGTTGATTTACTAGTCAAAACAACTCCCATGACGTCAATGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCC

**SnaBI (4249)**

4201 CATTGATGACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCATGACTGGGC

**NdeI (4354)**

4301 ATAATGCCAGGCGGCCATTTACCGTCAATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGACTGCCAAGTGGGCGAGTTACCGTAAA

4401 TACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTGACGTCAATGGGCGGGGTCGTTGGCGGTGACG

**PstI (4533)**  
**SdaI (4532)**  
**SbfI (4532)** **PciI (4550)**  
**BspLU11I (4550)**

4501 CAGGCGGGCCATTTACCGTAAGTTATGTAACGCCCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCCGC  
4601 TTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAATAATCGACGCTCAAGTCAGAGTGGCGAAACCGACAGGACTATAAAGATACC

4701 AGGCGTTTTCCCTGGAAGTCCCTCGTGCCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCT

**ApaLI (4864)**

4801 TTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGTTGAGTGTGAGTTCGCTCCAAGTGGGCTGTGTGCACGAACCCCGTTCAGCCCGACCGCTGCGCC

4901 TTATCCGGTAACTATCGTCTTGAGTCAACCCGGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAG

5001 GCGGTGCTACAGAGTTCTTGAAGTGGTGCCTAACTACGGCTACACTAGAAGACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAA

5101 AAGAGTTGGTAGCTCTTGATCCGGCAAAACAAACCCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAA

**SwaI (5289)** **NotI (5299)**

5201 GAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGAACGAAACTCAGCTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAG

**EagI (5300)**

5301 CGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAACGAAACAA

5401 AACAACTAGCAAAATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA