



**EcoRI (19)**

**NotI (2)**

**XbaI (15)**

**SdaI (29)**

**SspI (49)**

1 GCGGCCGCTATGCATCTAGAATTCCTGCAGGAGCTTGAATAAAATGAATATTAGAAGCTGTTAGAATAAGAGAA

75 AATGACAGAGGAAAAGTGAAGGGGAGAACTGAAAGTGGGAAATTCCTCTGAGGCAGAAAGGACCATCCCTTATA

149 AATAGCACAGGCCATGAAGGAAGATCATTCTCACTGCAGCCTTTGACAGCCTTTGCCTCATCTTGCAGGTAGCA

**BspHI (245)**

223 GCCGACACCAGCCTGGCTTCCATCATGATGGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGA  
M M E I K V L F A L I C I A V A E

**BglIII (363)**

297 GGCAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTTGCCACCACAGATCTTG  
17▶ A K P T E I N E D L N I A A V A S N F A T T D L

371 AGACTGACCTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCT  
42▶ E T D L F T N W E T M N V I S T D T E Q V N T D A

445 GACAGGGGCAAGCTGCCTGGCAAAAAGTCCCCCAGATGTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGC  
67▶ D R G K L P G K K L P P D V L R E L E A N A R R A

519 TGGTTGCACAAGAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTG  
91▶ G C T R G C L I C L S H I K C T P K M K K F I P

**EcoRV (660)**

593 GCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCCA  
116▶ G R C H T Y E G E K E S A Q G G I G E A I V D I P

667 GAGATTCCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCAC  
141▶ E I P G F K D K E P L D Q F I A Q V D L C A D C T

741 CACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCCAGAGGTGTA  
165▶ T G C L K G L A N V Q C S D L L K K W L P Q R C

**NheI (882)**

815 CCACTTTTGCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCT  
190▶ T T F A S K I Q G R V D K I K G L A G D R •

889 GGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATT

963 TGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTG

1037 CATTCAATTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTG

1111 GTATGGAATTAATTCTAAAATACAGCATAGCAAAAACCTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTC  
→

1185 TGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTT  
1259 CATGGAGTTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTGCTCTTCATTTCTTTATGTTTTAAATGCACT

**SspI (1361)**

1333 GACCTCCCACATTCCCTTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTT  
1407 TTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTGTAGTTGGACTTAGGGAACA

1481 AAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCAC  
1254 • D Q E E A V

1555 AAAGTGCACGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCA  
1184 F H V C N G A P D R L A F E R G W P Q E G I E T M

1629 TGGCCGGCCCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCG  
934 A P G S A D R F N T S V V E S W E A Y L E D L G

1703 CGCACCCACACCCAGGCCAGGGTGTGTCCGGCACCACTGGTCTGACCGCGCTGATGAACAGGGTCACGTC  
684 R V W V W A L T N D P V V Q D Q V A S I F L T V D

**SgrAI (1789)**

1777 GTCCCGGACCACACCGGCGAAGTCTGCTCCACGAAGTCCCGGGAGAACCAGCGGTCGGTCCAGAACTCGA  
444 D R V V G A F D D E V F D R S F G L R D T W F E V

1851 CCGCTCCGGCGACGTCGCGCGGGTGGACACCGGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCCTGTC  
194 A G A V D R A T L V P V A S T L K A M ←

1925 AGGAGAGGAAAGAGAAGAAGTTAGTACAATTGCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCA

1999 ATGATTAATTGTCAAAGTGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAAC  
←

2073 CGTAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTC

2147 AAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCGCT  
2221 CTCCTGTTCCGACCCTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCAT  
2295 AGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCCCGT  
2369 TCAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCAC  
2443 TGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGG  
2517 CCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAG  
2591 AGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTA  
2665 CGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAACGAAAAAC  
2739 TCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA