



EcoRI (23)

NotI (2) XbaI (19) SdaI (38) SpeI (45) SspI (59)

1 GCGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGGCCCACTAGTCGCTACAATATTTTCTGAACGGA

75 AGAATAAATAAACTTGTCTGTAAAGAAAACCCAGGTAAGGAAAGTGGCAGTCCAGACTGCCCGGAAGTTCC

149 TGGAGGCTAAGGCCTCACCCCGTCGCTTGATAGGACCTGCTGAGCCACATGACTAAGGCACGATCGCTCCGG

AflII (272)

223 ACGTGTAAAGGTGCTGGGTTCCAAGATGGCTGCCCGCGGAGGCCCGACTTAAGTATGTCACTTCCGCACCA

297 GCGAGAAAAGCGGACCCCTCAGCCAATGAGGCCATAGGGCGGGGCTAGGCCATGATGGGCTTCAAACCTACCCA

371 ATAGGGCGTCCGAACTAAAGCGCCTACAAAGTAACGTCACGTCGAGTTGCAGAGCGCCGGCAGCGGGGCAGAG

445 GTGGCCAAGCCAATGCGATGGCTGGGGCGGGGTCGGACGCTCTATAAGTTGTCGATAGCGGGCACTCCGCCCT

NcoI (531)

519 AGATTCTAAGGACCATGGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAACCCACT

1 M E I K V L F A L I C I A V A E A K P T

BglII (646)

593 GAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCAAACCTTGGCCACCACAGATCTTGAGACTGACCTGTT

21 E I N E D L N I A A V A S N F A T T D L E T D L F

667 CACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGC

45 T N W E T M N V I S T D T E Q V N T D A D R G K

741 TGCCTGGCAAAAACTCCCCCAGATGTCCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGTTGCACAAGA

70 L P G K K L P P D V L R E L E A N A R R A G C T R

815 GGCTGCCTCATTTGCCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACAC

95 G C L I C L S H I K C T P K M K K F I P G R C H T

889 TTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCT

119 Y E G E K E S A Q G G I G E A I V D I P E I P G

963 TCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGCTG

144 F K D K E P L D Q F I A Q V D L C A D C T T G C L

1037 AAGGGCCTTGCAATGTCCAGTGTCTGACCTCTGAAGAAAGTGGCTTCCCAGAGGTGTACCACTTTTGCCAG

169 K G L A N V Q C S D L L K K W L P Q R C T T F A S

NheI (1165)

1111 CAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGAT

193 K I Q G R V D K I K G L A G D R

1185 AAGATACATTGATGAGTTTGACAAAACCACAACCTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTG

1259 ATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCAATTTATG

1333 TTTGAGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATTAAT

1407 TCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATA

1481 AGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAG

1555 ATATAGTGTATTTTCCAAAGTTTGAAGTACTAGCTCTTCATTTCTTATGTTTTAAATGCACTGACCTCCACATT

SspI (1644)

1629 CCCTTTTGTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAG

1703 AATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTGTAGTGGACTTAGGGAACAAAGGAACCTTTAA

1777 TAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTGCACGCAG

125 D Q E E A V F H V C

1851 TTGCCGCGGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTTCATGGCCGGCCCGGA

113 N G A P D R L A F E R G W P Q E G I E T M A P G S

1925 GGCCTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACACC

89 A D R F N T S V V E S W E A Y L E D L G R V W V W

SgrAI(2072)

1999 AGGCCAGGGTGTGTCGGCACCACCTGGTCCTGGACCGCGCTGATGAACAGGGTACGTCGTCGGGACCACA

64 A L T N D P V V Q D Q V A S I F L T V D D R V V

2073 CCGGCAAGTCGCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGAGTCCAGAACTCGACCGCTCCGGCGAC

39 G A F D D E V F D R S F G L R D T W F E V A G A V

2147 GTCGCGCGGGTGGACCCGAAACGGCACTGGTCAACTGGCCATGATGGCTCCTCCTGTCCAGGAGAGGAAAGA

15 D R A T L V P V A S T L K A M

2221 GAAGAAGGTTAGTACAATTGCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCAATGATTAATTGTC
2295 AAACTAGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCG
2369 CGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGG
2443 CGAAACCCGACAGGACTATAAGATACCAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGAC
2517 CCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTA
2591 GGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGC
2665 TGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCAC
2739 TGTTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCT
2813 AACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCT
2887 TGATCCGGCAAACAACCACCGCTGGTAGCGGTGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAA
2961 AGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGA
3035 TTTTGGTCATGGCTAGTTAATTAACATTTAAATCA