



EcoRI (23) SdaI (38)
NotI (2) XbaI (19) BamHI (29) SpeI (45) NotI (78)
1 CGCGCCGCGTCGACGATATCTAGAATTCGGATCCTGACAGGGCCACTAGTTCGCCAGAGCGCGGAGGGCCTCCAGCGGCCGCCCTCCCCACAGCAG
101 GGGCGGGGTCGCCGCCACCGAAGGAGCGGGCTCGGGGCGGGCGGCGCTGATTGGCCGGGCGGGCCTGACGCCAGCGGCTATAAGAGACCACAAG
HindIII (230) EcoNI (276)
201 CGACCCGACAGGGCCAGACGTTCTTCGCCGAAGCTTCGAGGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCCGCCATCCACGCCGGTT
301 GAGTCGCGTTCGCGCCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCGCCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACCGGGCCTTTGTCCGGCG
401 CTCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGCTTTGTTCGTTTTCTGTTCTGCGCCGTT
NeoI (515)
501 ACAGATCCAAGCCACCATGGAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATA
1 M E I K V L F A L I C I A V A E A K P T E I N E D L N I
BglIII (630)
601 GCTGCTGTGGCCTCAACTTTGCCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGA
29 A A V A S N F A T T D L E T D L F T N W E T M N V I S T D T E Q V
701 ACACAGATGCTGACAGGGGAAGCTGCCTGGCAAAAACTCCCCAGATGCTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAAGAGG
62 N T D A 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 G C T R G
801 CTGCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTAAAAGGAGTCTGCTCAG
95 C L I C L S H I K C T P K M K K F I P G R C H T Y E G E K E S A Q
901 GGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTG
129 G G I G E A I V D I P E I P G F K D K E P L D Q F I A Q V D L C A
1001 ATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTCTCTGACCTCCTGAAGAAGTGGCTTCCCCAGAGGTGTACCACTTTTGCCAGCAAGAT
162 D C T T G C L 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 K I
NheI (1149)
1101 TCAGGGTAGGGTGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGACAAAACACAA
195 Q G R V D K I K G L A G D R
1201 CTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACA
1301 TTGCATTCATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATAC
1401 AGCATAGCAAACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATT
1501 AGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGTATTTTCCAAGTTTGAACCTAGCTCTTCATTTCTTTATGTTTAAATGCACTG
SspI (1628)
1601 ACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAAATAATGTTTTTTATTAGGCAGAATCCAGATGCTCA
1701 AGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCC
1801 TCAGTCTGCTCCTCTGCCACAAAGTGCACGCAAGTGCAGGCTGCGCGCCGGGTGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTGATGGC
125 D Q E E A V 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 A
1901 GGCCCGGAGGCTCCCGAAGTTCTGTTGACACGCTCCGACCACTCGGCGTACAGCTGCTCCAGGCGCGCACCCACCCAGGCCAGGGTGTGTCGG
91 P G S 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 A L T N D P
SgrAI (2056)
2001 GCACCACCTGGTCTGGACCGGCTGATGAACAGGGTACGTCGTCGCCGACACACCGGCGAAGTGCCTCCACGAAGTCCCGGGAGAACCCGAGCCG
58 V V Q D Q V A S I F L T V D D R V V G A F D D E V F D R S F G L R
2101 GTCGGTCCAGAACTGACCGCTCCGGCAGCTGCGCGCGGTGAGCACCAGGACCGGCACTGGTCAACTTGGCCATGATGGCTCCTCCTGTCAGGAGAGGA
25 D T W F E V A G A V D R A T L V P V A S T L K A M
2201 AAGAGAAGAAGTTAGTACAATTGCTATAGTGTGATTACTATGCAGATATACTATGCCAATGATTAATTGTCAAAGTGGGCTGCAGGTTAATT
2301 AAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTGTGCGGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACA
2401 AAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGAC
2501 CCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTCGGGAAGCGTGGCGTTTTCTCATAGCTCACGCTGATAGGTATCTCAGTTCGGTGTAGGTCGTT
2601 CGCTCAAAGTGGGCTGTGTGCACGAACCCCGTTACGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGTAAGACAGACT
2701 TATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCTAACTACGGCTACAC
2801 TAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGTATCCGGCAACAAACCACCGCTGGTAGC
2901 GGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGAACG
3001 AAAACTCACGTTAAGGGATTTGGTCATGGCTAGTTAATTAACATTTAAATCA