



EcoRI (23)
NotI (2) XbaI (19) SdaI (38) SpeI (45)
1 CGCGCCGCGTCGACGATATCTAGAATTCGGATCTGCAGGGCCACTAGTTACCGACCACCCGCAAACAGCAGGGTCCCTGGGCTTCCCAAGCCGCGCA

SphI (188)
101 CCTCTCCGCCCCGCCCTGCGCCCTCTTCTCGCGTCTGCCCTCTCCCCACCCGCTTCTCCCTCCCCGCCCAGCGGCGCATGCGCCGCGCTCGG
201 AGCGTGTTTTTATAAAGTCCGGCCGCGCCAGAACTTCAGTTTGTGGCTGCGGCAGCAGGTAGCAAAGTGACGCCGAGGGCTGAGTGCTCCAGTAG

NcoI (328)
301 CCACCGCATCTGGAGAACCAGCGGTTACCATGGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGA
M E I K V L F A L I C I A V A E A K P T E I N E

BglII (443)
401 AGACCTCAATATAGCTGCTGTGGCCTCCAACCTTTGCCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGAC
24 D L N I 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
501 ACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTG
58 T E Q V 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
601 GTTGACAAGAGGCTGCCTCATTGGCTCTCCACATTAAGTGACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGTGAAAA
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 G R C H T Y E G E K
701 GGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTG
124 E S A Q 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
801 GACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCTGAAGAAAGTGGCTTCCCCAGAGGTGTACCACTT
158 D L C A 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

NheI (962)
901 TTGCCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTT
191 F A S K I Q G R V D K I K G L A G D R •
1001 GGCAAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAACAAG
1101 TTAACAACAACAATTGCATTCAATTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATT

SspI (1441)
1201 AATTCTAAAAACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTT
1301 GCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTACTCTTCAATTTCTTTATGT
1401 TTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGA
1501 ATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTTGGACAGCAAGAAAGCGAGC
1601 TTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAAGTGACGACAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGAT
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
1701 CTCGGTCAATGGCCGCGCCGAGGCGTCCCGGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCGCGCACCCACCCAGGCG
96 E T M A 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

SgrAI (1869)
1801 AGGGTGTGTCCGGCACCACCTGGTCTGGACCGCGCTGATGAACAGGGTACGTCGTCCCGACCACACCGGCGAAGTCTGCTCCACGAAGTCCCGGG
62 L T N D P 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
1901 AGAACCCGAGCCGGTCCGAGAACTCGACCCGCTCCGGCGACGTCGCGCGGTTGAGCACCAGGAAACGCACTGGTCAACTTGGCCATGATGGCTCCTCC
29 F G L R 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
2001 TGTCAGGAGAGGAAAGAGAAGGTTAGTACAATTGCTATAGTGTGATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAA ACTAGGG
2101 CTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGCGTTTTTCCATAGGCTCCGCCCCCT
2201 GACGAGCATCAGAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCCTGGAAGCTCCCTCGTGCCT
2301 CTCCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTC
2401 GGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTACGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCG
2501 GTAAGACACGACTTATGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGTACAGAGTTCTTGAAGTGGTGGCTA
2601 ACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAACAAC
2701 CACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGGTCTGAC
2801 GCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCTAGGTTAATTAACATTTAAATCA