



EcoRI (23) SdaI (38)
NotI (2) XbaI (19) BamHI (29) SpeI (45) NotI (78)
1 CGCGCCGCGTTCGACGATATCTAGAATTCGGATCCTGCAGGGCCACTAGTTCGCCAGAGCGCGAGGGCTCCAGCGGCCGCCCTCCCCACAGCAG
101 GGGCGGGTCCCAGCCACCGAAGGAGCGGGCTCGGGGCGGGCGGCTGATTGGCCGGGGCGGGCTGACGCCGACGCGGCTATAAGAGACCACAAG
201 CGACCCGAGGGCCAGACGTTCTTCGCCGAGAGTCGTGGGGTTTCTGCTTCAACAGTGCTTGACGGAACCCGGCGCTGTTCCCCACCCCGCGCGC
301 CGCCATAGCCAGCCTCCGTACCTCTTACCACCTCGGACTGCCCAAGGCCCGCGCGCTCCAGCGCCGCGAGCCACCGCGCGCGCGC
401 GCCTCTCCTTAGTCGCCGCATGGAAATCAAGGTGCTGTTGCCCTCATCTGTATTGCTGTTGCTGAGGCAAACCCACTGAAATCAATGAAGACCTCAA
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
BglII (534)
501 TATAGTCTGTGGCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTCAACCACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAG
27 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 T E Q
601 GTGAACACAGATGCTGACAGGGCAAGCTGCCTGGCAAAAACTCCCCAGATGCTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAA
61 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 G C T
701 GAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGC
94 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 E S A
801 TCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCCTGGCTTCAAGGATAAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGT
127 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 D L C
901 GCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCATGTCCAGTGTCTGACCTCCTGAAGAAAGTGGCTTCCCCAGAGGTGACCCTTTTGCAGCA
161 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 F A S
NheI (1053)
1001 AGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACC
194 K I Q G R V D K I K G L A G D R •
1101 ACAACTAGAATGCAGTAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAACAAGTTAACAACA
1201 ACAATTGCATTATTATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTAATTCTAAA
1301 ATACAGCATAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTG
1401 CATTAGCTGTTTCGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTCTCTCATTCTTTATGTTTAAATGC
SspI (1532)
1501 ACTGACCTCCACATTCCCTTTTATGTAATAATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAATGTTTTTTATTAGGCAGAATCCAGATG
1601 CTCAAGGCCCTTCATAATATCCCCAGTTTAGTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTT
1701 ATCCTCAGTCTGCTCCTCTGCCACAAAGTGCACGAGTTGCCGGCGGGTGCAGCGGGGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCA
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
1801 GCGCGGCCGGAGGCGTCCCGAAGTTCGTGGACAGCCTCCGACCACTCGGCGTACAGCTCGTCCAGGCGCGCACCCACCCAGCCAGGGTGTG
93 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 L T N
SgrAI (1960)
1901 TCCGGCACCACTGGTCTGACCGCGCTGATGAACAGGGTCACGTCGTCCGGACCACACCGCGGAAGTCTCTCCACGAAGTCCCGGGAGAACC
59 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 F G L
2001 GCCGGTGGTCCAGAAGTCCGACCGTCCGGCGAGCTCGCGCGGGTGGACCGAACGGCACTGGTCAACTGGCCATGATGGCTCCTCTGTCCAGGAG
26 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
2101 AGGAAAGAGAAGAAGGTTAGTACAATTGCTATAGTGTGATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAACTAGGGCTGCAGGTT
2201 AATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGCCGCTTGTGCGGTTTTCCATAGGCTCCGCCCCCTGACGAGCAT
2301 CACAAAAATCGACGCTCAAGTCAAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCGTTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTT
2401 CGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGT
2501 CGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCAACCCGGTAAGACAC
2601 GACTTATCGCCACTGGCAGCAGCCTGTTAAGCAGGATTAGCAGAGCGAGGTATGTAGGCGGTGTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCT
2701 AACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGTCCGGCAACAAACCACCGCTGG
2801 TAGCGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGG
2901 AACGAAAACCTCACGTTAAGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA