



EcoRI (19)
NotI (2) **XbaI (15)** **SdaI (29)** **SpeI (36)** **NcoI (70)**
1 CGGCCCGCTATGCATCTAGAATTCTCTGAGGGCCCACTAGTCTCCAGGCATGACTCCAACAATGCATCCCATGGGATTGGGGTCCCCAGATCTGGGG
101 CTTGTAGGCTGACTCTCCCTGTGCACACGCTCTACACGCATGCGTGCACCATTGCTGCCCGCCCTTGACAGGGAGTCAGCAGGGAGGACTG
201 GGTATGCCCTGCTTATCAGCAGCTTCCAGCTTCTCTGCTGATTCTTAGAGGCTGGGGTCTAGAACGAGCTGGTGCACGTGGCTTCCCAAAGT
301 CTCTCAGATAATGAGAGGAAATGCAGTCATCAGTTTGAGAAGGCTAGGGATTCTGGCCATAGCTCAGACCTGCGCCACCATCTCCCTCAGGCAGCC
BspHI (437)
401 CTTGGCTGGTCCCTGCGAGCCGTGGAGACTGCCAGTcATGATGAAATCAAGGTGCTGTTGCCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCAC
501 TGAATCAATGAAGACCTCAATATAGCTGTGGCCTCCAATTTGCCACCACAGATCTGAGACTGACCTGTTCACTCACTGGGAGACCATGAATGTG
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 T N W E T M N V
601 ATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGCTCTGAGGAGCTGGAGGCCAATG
55▶ I S T D 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
701 CCAGAAGGCTGGTTGACAAGAGGCTGCCTCATTGCTCTCCACATTAAAGTGCACCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACCTTA
88▶ A R R A 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
EcoRV (852)
801 TGAAGTGAAAAGGAGTCTGCTCAGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTCAAGGATAAGGAGCCACTGGACCAGTTT
121▶ E G E K 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
901 ATTGCTCAAGTGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCCAGA
155▶ I A Q V 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
NheI (1074)
1001 GGTGTACCCTTTTCCAGCAAGATTGAGGTAGGGTGGACAAAATCAAGGTCTGGCTGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATA
188▶ R C T T F A S K I Q G R V D K I K G L A G D R •
1101 ATTGATGAGTTTGGACAAACCACAAC TAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGT
1201 GCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAACCTCTACAATG
1301 TGGTATGGAATTAATCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCA
1401 TCAGGGGCTGTTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAGTTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTGCTTTC
SspI (1553)
1501 ATTTCTTATGTTTTAAATGACTGACCTCCCACATCCCTTTTTAGTAAAAATTTCAGAAATAATTTAAATACATCATTGCAATGAAAATAATGTTTT
1601 TTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAACCTTTAATAGAAATGGACAGC
1701 AAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCTCTGCCACAAAGTGACGACGAGTTGCCGGCCGGTCCGCGAGGGCGAACTCCCGCCCCACGG
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
1801 CTGCTCGCGATCTCGGTTCATGGCCGGCCCGAGGCGTCCCGAAAGTTCTGGACACGACCTCCGACCACTCGGGGTACAGCTGCTCAGGCGCGCAC
100◀ Q E G I 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
SgrAI (1981)
1901 CACACCCAGGCCAGGTTGTTCCGGCACCCTGGTCTGGACCGGCTGATGAACAGGGTACGTCGTCGCCGACACACCGGCGAAGTCTGCTCCCA
66◀ W V W A 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
2001 CGAAGTCCCGGAGAACCCGAGCCGCTCCGAGAACTGACCTCCGCGGAGTCCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCCAT
33◀ F D R S 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
2101 GATGGCTCTCTGTGAGGAGGAAAGAGAAGGTTAGTACAATTGCTATAGTGAAGTTGATTATACTATGCAGATATACTATGCAATGATTAAT
2201 GTCAACTAGGGCTGCAGTTAATTAAGAACATGTGAGCAAAAGCCAGCAAAAGCCAGGAACCGTAAAAAGCCGCTTGTGGCGTTTTTCCATAGG
2301 CTCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCT
2401 CCTCGTGGCTCTCTGTTCCGACCTGCGCTTACCGGATACCTGTCCGCTTTCTCCCTCGGGAAGCGTGGCGCTTCTCATAGCTCAGCTGTAG
2501 GTATCTCAGTTCGGTGTAGTCTGTTCCGCTCAAGCTGGCTGTGTGCACGAACCCCGTTCCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTT
2601 GAGTCAACCCGGTAAGACACGACTTATGCCACTGGCAGCAGCCACTGTTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTCTACAGATCTTGA
2701 AGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATC
2801 CGGCAAAACAACCCTGTTAGCGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCTTTTGTCTTTTCT
2901 ACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA