



EcoRI (23)

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

1 CGCGCCGCTGCACGATATCTAGAATTCGGATCCTGACAGGCCACTAGTTTCCAGAGCTTTTCGAGGAAGTTTCTTCAACTCAAATTCATCCGCTGAT
101 AATTTTCTTATATTTTCTAAAGAAGGAAGAGAAGCGCATAGAGGAGAAGGAAATAATTTTTAGGAGCTTTCTTACGGCTATGAGGAATTTGGGGCT

Tth111I (271)

201 CAGTTGAAAAGCTAAACTGCCTCTCGGAGGTTGGGCGGGCAACTACTTTTCAGCGGCACGGAGACGGCGTCTACGTGAGGGGTGATAAGTGACGC
301 AACACTCGTTGCATAAAATTTGCGCTCCGCCAGCCCCGGAGCATTAGGGGCGTTGGCTTTGTTGGGTGAGCTTGTGTCCCTGTGGGTGACGTGGT
401 TGGTGATTGGCAGGATCCTGGTATCCGCTAACAGGtactggcccacagccgtaaacacccgctgagaggggggaatgggtgaggtcaagctgg
501 aggccttctggggttgggtgggcccgtgaggggaggggagggcgaggtgacgcgacaccggccttctgggagagtgggccttgttgacctaagggggg
601 cgagggcagttggcagcgcgacgcccacagaaactaacagacattaaccaacagcgattccgtcgctttacttgggaggaaggcggaaaagaggtag

NruI (786)

701 tttgtgtggcttctggaaccctaaatttggaaatcccagtatgagaatggtgtcccttctgtgtttcaatgggatttttacttcgagagcttctgtgggt

BspEI (877)

801 ttggttttgtttcagtttgcctaaccctgcttaggtttgaggcagattgagttcggtcggggagtttgaatatccggaacagtttagtgggaaag
901 ctgtggacgcttgtaagagagcgtctgattttccgctgttgacgttgaacctgaaatgacgaatttcgtattaagtacttagccttgtaaaattg
1001 aggggaggcttgcggaatattaactgatttaagcattttgaaggaatagttgctaattttgaagaatattaggtgtaaaagcaagaataacaatgatcc

NcoI (1140)

1101 tgaggtgacacgcttatgttttacttttaactagGTCACCATGGAAATCAAGGTGCTGTTTGCCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCAC
1 M E I K V L F A L I C I A V A E A K P T

BglII (1255)

1201 TGAAATCAATGAAGACCTCAATATAGTGTGTGGCCTCCAACCTTGCCACCACAGATCTTGAGACTGACCTGTTCACTCACTGGGAGACCATGAATGTG
20 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
1301 ATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACCTCCCCAGATGCTCCTGAGGGAGCTGGAGGCCAATG
54 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
1401 CCAGAAGGGCTGGTGCACAAGAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTA
87 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
1501 TGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCCCTGGCTTCAAGGATAAGGAGCCACTGGACAGTTT
120 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
1601 ATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCTGAAGAAGTGGCTTCCCGAGA
154 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 (1774)

1701 GGTGTACCACCTTTTCCAGCAAGATTGAGGTGAGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGTGGCCAGACATGATAAGATAC
187 R C T T F A S K I Q G R V D K I K G L A G D R
1801 ATTGATGAGTTTGGACAAACCACAACACTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCT

1901 GCAATAACAAGTTAACAACAACAATTGCATTCATTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATG

2001 TGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCA

2101 TCAGGGGCTGTGCAATGTGCATTAGTGTGTTGACGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTGTCTCTTC
2201 ATTTCTTTATGTTTAAATGCACTGACCTCCACATTCCTTTTATGATAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTT
2301 TTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTGTAGTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTTGGACAGC
2401 AAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCTCTGCCACAAAGTGCACGAGTTGCCGGCCGGTGCAGCAGGGCAACTCCCGCCCCACGG
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
2501 CTGCTCGCCGATCTCGGTCATGGCCGGCCGGAGGCGTCCCGAAGTTCGTGGACACGACTCCGACCACTCGGCTGACAGCTCGTCCAGGCCGCGCAC
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 (2681)

2601 CACACCCAGGCCAGGGTGTGTCGGCACCACTGGTCTGGACCGCTGATGAACAGGGTCACGTCGTCGCCGACACACCGGCGAAGTCGTCCTCCA
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
2701 CGAAGTCCCGGAGAACCCGAGCCGGTCCGTCAGAACTCGACCGCTCCGGCGACGTCGCGCGCGGTGAGCACCAGAACGGCACTGTTCAACTGGCCAT
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
2801 GATGGCTCCTCTGTCAGGAGAGAAAGAGAAGGTTAGTACAATTGCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCAATGATTAAT

2901 GTCAACTAGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAGGCCGCGTGTGGCGTTTTTCCATAGG

3001 CTCCGCCCTCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTCGGAAGCT

3101 CCCTCGTGGCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAG

3201 GTATCTCAGTTCGGTGTAGGCTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTT

3301 GAGTCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGTAACAGGATTAGCAGAGCGAGGTATGTAGGGCGTGTACAGAGTCTTGA

3401 AGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATC

3501 CGGCAACAACCACCGCTGGTAGCGGTGTTTTTTGTTTGAAGCAGCAGATTACGCCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCT

3601 ACGGGGTCTGACGCTCAGTGAACGAAAACACGTTAAGGGATTTGGTCATGGCTAGTTAATTAACATTTAATCA