



EcoRI (19)

NotI (2) XbaI (15) SdaI (29) SpeI (36)

1 CCGCCGCTATGCATCTAGAATTCCTGCAGGGCCACTAGTTCCTTTGAAAGCAGTCGAGGGGTGCTAGGTGTGGCAGGGACGAGCTGGCCGGCGCTCG
101 CTGGGTGCACCGCACCAGCGGACAGCCACGGCGGGAGGACTACAACCTCCCGGCACACCCCGCCGCCCTCTACTCCAGAAGGCCGGGG

AscI (244)

201 GGTGGACCGCCTAAGAGGGCTGCGCTCCCGACATGCCCGCGCGCCATTAACCGCCAGATTTGAATCGCCGGACCCGTTGGCAGAGGTGGCGGG

BspHI (306)

301 CCGCATCATGATGGAATCAAGGTGCTGTTGCCCTCATCTGATTGCTGTTGCTGAGGCAAAACCACTGAAATCAATGAAGACCTCAATATAGCTGCT
M 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 A A

BglII (424)

401 GTGGCTCCAACCTTGGCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAG
32 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 N T
501 ATGCTGACAGGGGCAAGCTGCTGGCAAAAACTCCCCCAGATGCTCTGAGGGAGCTGGAGGCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCT
65 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 C L
601 CATTTCCTCTCCACATTAAGTGACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGG
98 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 G G

EcoRV (721)

701 ATTGGAGAGCAATTGTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCA
132 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 D C
801 CCACTGGCTGTCTGAAGGGCTTGCATGTCAGTGTCTGACCTCTGAAGAAGTGGCTTCCCGAGAGGTGTACCACCTTTGCCAGCAAGATTCAGGG
165 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 Q G

NheI (943)

901 TAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAC TAGAA
198 R V D K I K G L A G D R
1001 TGCAGTGAIAAATGCTTTATTGTGAAATTTGTGATGCTATTGCTTTATTGTAACCATTATAAGCTGCAATAAACAAGTTAACAAACAATTCAT
1101 TCATTTTATGTTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTAATCTAAAATACAGCATA

1201 GCAAAATTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGT
1301 TTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTCCCAAGTTTGAAC TAGCTCTTCATTCTTTATGTTTTAAATGCACTGACCTCC

SspI (1422)

1401 CACATTCCTTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCC
1501 TTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGAAACAAAGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTC
125 D
1601 CTGCTCCTCTGCCACAAAGTGACCGCAGTGTCCCGCCGGGTGCGCGAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTTCATGGCCGGCCCC
123 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 P G
1701 GAGGCGTCCCGGAAGTTCGTGGACACGACCTCCGACCACTCGGGCTACAGCTCGTCCAGGCCGCGCACCCACACCAGGCCAGGTGTTGTCCGGCACCA
89 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 V V

SgrAI (1850)

1801 CCTGGTCTGGACCGCGTGTGATGAACAGGGTCACTGCTCCCGGACCACACCGGCAAGTCTGCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGT
56 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 D T
1901 CCAGAACTCGACCCGCTCCGGCGACGTCGCGCGCGTGGACACCGGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCCTGTCAGGAGAGGAAAGAGA
23 W F E V A G A V D R A T L V P V A S T L K A M
2001 AGAAGGTAGTACAATTGCTATAGTGAAGTGTATTACTATGCAGATATACTATGCCAATGATTAATTGTCAAAGTGGGCTGCAGGTTAATTAAGAAC

2101 ATGTGAGCAAAAGCCAGCAAAAGGCCAGGAACCGTAAAAGGCCGCGTGTGCTGGCTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATC

2201 GACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCC

2301 GCTTACCGGATACCTGTCCGCTTTCTCCCTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGGTGAGGTGCTCGCTCC

2401 AAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTCCAACCCGTAAGACACGACTTATCGC

2501 CACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAG

2601 AACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACACCACCCTGGTAGCGGTGGT

2701 TTTTTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGTCTGACGCTCAGTGAACGAAAATC

2801 CACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA