



EcoRI (23)

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

1 CGCGCCGCGTCGACGATATCTAGAATTCGGATCTCGAGGGCCaCTAGtGTTTCATCGGAGCCAGGTTTACTCCCTTAAGTGGAAATTTCTTCCCCAC
101 TCCCTCTTGCTTTCTCCAAGGAGGAACCCAGGCTACTGAAAGTCCGGCTGGGGCGGGGACTGTGGTTTCAGGGTAGAACTGCGTGTGGAACGGGA
201 CAGGGAGCGGTTAGAAGGGTGGGGCTATTCCGGGAAGTGGTGGGGGAGGGAGCCAAAAGTACACCTAGTCCACTCATTATCCAGCCCTCTTATTCT

BsrGI (380)

301 CGGCCCCGCTCTGCTTCAGTGGACCCGGGAGGGCGGGGAAGTGGAGTGGGAGACCTAGGGTGGGCTTCCCGACCTTGTGTACAGGACCTCGACCTAG
401 CTGGCTTTGTTCCCATCCCCACGTTAGTTGTTGCCCTGAGGCTAAAAGTACAGCCAGGGGCCCAAGTTCAGACTGCCCTCCCCCTCCCCGGAG
501 CCAGGGAGTGGTTGGTGAAGGGGAGGCCAGCTGGAGAACAACGGGTAGTCAGGGGTTGAGCGATTAGACCTTGTACCTACCCAGGAATGGTTG
601 GGGAGGAGGAGGAAGAGGTAGGAGGTAGGGGAGGGGGCGGGGTTTTGTACCTGTACCTGCTCCGGCTGTGCTAGGGCGGGCGGGGAGTGGGG

AgeI (702)

701 GACCGGTATAAAGCGGTAGGCGCTGTGCCGCTCCACCTCTCAAGCAGCCAGCGCTGCCTGAATCTGTTCTGCCCTCCCCACCATTTACCACCA

NcoI (801)

801 CCTATGAAATCAAGGTGCTGTTTCCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGCCTC
▶ 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 V A S

BglII (916)

901 CAACTTTGCCACCACAGATCTTGAGACTGACCTGTTCACTCACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGAC
33▶ 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 D A D
1001 AGGGGCAAGCTGCCTGGCAAAAACTCCCCAGATGCTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTGGC
67▶ 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 I C
1101 TCTCCACATTAAGTGACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGA
100▶ 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 I G E
1201 GGCAATTGTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCACTGGACCAAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGC
133▶ 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 T T G
1301 TGTCTGAAGGGCTTGCATGTCCAGTGTCTGACCTCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTTTCCAGCAAGATTGAGGGTAGGGTG
167▶ 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 R V

NheI (1435)

1401 ACAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTGA
200▶ D K I K G L A G D R •
1501 AAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCAATTTA
1601 TGTTTCAGGTTCCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTAATCTAAAATACAGCATAGCAAACCT
1701 TTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCC
1801 TCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTACTCTTCATTTCTTATGTTTTAAATGCACTGACCTCCACATTCC

SspI (1914)

1901 CTTTTTAGTAAAATATTAGAAATAATTTAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTCATAAT
2001 ATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCAGGCTTCTAGCTTATCCAGTCTGCTCCT
125▶ • D Q E E
2101 CTGCCACAAAGTGCACGAGTTGCCGCGGGGTCGCGCAGGGCGAACTCCC GCCCCACGGCTGCTCGCGATCTCGGTCATGGCCGCGCCGGAGGGGTC
120▶ 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 S A D
2201 CCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCGCGCACCCACACCAGCCAGGGTGTGTCCGGCACCCTGGTCC
87▶ 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 Q D

SgrAI (2342)

2301 TGGACCGCGTATGAACAGGGTACGTCGTCGCCGACACACCGCGAAGTCTCTCCACGAAGTCCCGGGAGAACCAGCGGCTCGGTCGAGAACT
53▶ 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 W F E
2401 CGACCGTCCGGCAGCTCGCGCGGGTGAACCGGAACGGCACTGGTCAACTGGCCATGATGGCTCCTCTGTCAGGAGAGGAAAGAGAAGAGTT
20▶ V A G A V D R A T L V P V A S T L K A M
2501 AGTACAATTGCTATAGTGTGATTATTAATACTATGCAGATATACTATGCAATGATTAATTGTCAAAGTAGGGCTGCAGGTTAATTAAGAACATGTGAGC
2601 AAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAAATCGACGCTCA
2701 AGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGCGTTTTCCCCTGGAAGCTCCCTCGTGGCTCTCCTGTTCCGACCTGCCGTTACCG
2801 GATACCTGTCCGCTTTTCTCCCTTCGGGAAGCGTGGCGTTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGGTAGGTGTTGCTCCAAGCTGGG
2901 CTGTGTGCACGAACCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCAACCCGGTAAGACACGACTTATCGCACTGGCA
3001 GCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTAT
3101 TTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACACCCTGGTAGCGGTGGTTTTTTTTGT
3201 TTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGCTGACGCTCAGTGGAAACGAAAACCTCACGTTAA
3301 GGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA