



EcoRI (19)

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

1 CGCGCCGCTATGCATCTAGAATTCCTGCAGGGCCCACTAGTGTCCACGCCTCCAGCAGGGGACGCCCGGGCTGGGGGCGGGAGTCAGACCGCGCCTG

BspEI (106)

101 GTACCATCCGGACAAAGCCTGCGCGGCCCGCCCGCCATTGGCCGTACCGCCCCGCGCCGCCCCATCCCGCCCTCGCGCCGGTCCGGCGCT

201 TAAAGCAATAGGAACCGCCGCTTGTCCCGTCACGGCCGGGGCAGCAATTGTGGCGGCTCGGGGCTCGTGCTCTTTTCGCGCAAAAAGGATT

NcoI (382)

301 TGGCGGTAAAGTGGCCGGACTTTGAGGCAGCGCGGCCGGGGCGGAGCGGGATCGAGCCCTCGCCGAGGCTGCCGCATGGGCCCGCGCCCG

BspHI (438)

SphI (460)

401 CCGCCGCTGTACCCGGGCGCGCGGGCCGTGAGCGTCATGATTCTGGGCCCTGCATGCTGCTGCTGCTGCTGCTGGGCTGAGGCTACAGCTCT

M I L G P C M L L L L L L L L L G L R L Q L

501 CCCTGGGCATCATCCAGTTGAGGAGGAGAACCCGGACTTCTGGAACCGCGAGGCAGCCGAGGCCCTGGGTGCCCAAGAAGTGCAGCCTGCACAGAC

21 S L G I I P V E E E N P D F W N R E A A E A L G A A K K L Q P A Q T

601 AGCCGCCAAGAACCTCATCATCTTCTGGCGATGGGATGGGGTGTCTAGCGGTGACAGCTGCCAGGATCTAAAAGGGCAGAAGAAGGACAAACTGGGG

54 A A K N L I I F L G D G M G V S T V T A A R I L K G Q K K D K L G

701 CCTGAGATACCCCTGGCTATGGACCGCTTCCCATATGTGGCTCTGCAAGACATACAATGTAGACAAACATGTGCCAGACAGTGGAGCCACAGCCACGG

88 P E I P L A M D R F P Y V A L S K T Y N V D K H V P D S G A T A T

801 CCTACCTGTGCGGGTCAAGGGCAACTTCCAGACCATTGGCTTGTAGTGCAGCCGCCGCTTTAACAGTGAACACGACACGCGGCAACGAGGTTCATCTC

121 A Y L C G V K G N F Q T I G L S A A A R F N Q C N T T R G N E V I S

901 CGTGATGAATCGGGCCAAGAAGCAGGGAAGTCAAGTGGGVTAAACCACACAGAGTGCAGCACGCCTGCCAGCCGGCACCTACGCCACACGGTG

154 V M N R A K K A G K S V G V T T T R V Q H A S P A G T Y A H T V

1001 AACCGCAACTGGTACTCGGACGCCGACGTGCCTCGCCCGCCAGGAGGGTGCAGGACATCGCTACGCGCTCATCTCCAACATGGACATTGATG

188 N R N W Y S D A D V P A S A R Q E G C Q D I A T Q L I S N M D I D

1101 TGATCCTGGTGGAGGCCGAAAGTACATGTTTCGCATGGGAACCCAGACCTGAGTACCAGATGACTACAGCCAAGTGGGACCAGGCTGGACGGGAA

221 V I L G G G R K Y M F R M G T P D P E Y P D D Y S Q G G T R L D G K

1201 GAATCTGGTGCAGGAATGCCTGGCGAAGCCAGGCTGCCGCTGTGTGGAAACCGACTGAGCTCATGAGGCTTCCCTGGACCCGCTGTGACCCAT

254 N L V Q E W L A K R Q G A R Y V W N R T E L M Q A S L D P S V T H

1301 CTCATGGGCTCTTTGAGCCTGGAGACATGAAATACGAGATCCACCGAGACTCCACTGGACCCCTCCCTGATGGAGATGACAGAGGCTGCCCTGCGCC

288 L M G L F E P G D M K Y E I H R D S T L D P S L M E M T E A A L R

SacII (1420)

1401 TGCTGAGCAGGAACCCCGCGCTTCTTCTTCTCGTGGAGGGTGGTGCATCGACCACGGTTCATCACGAAAGCAGGGCTTACCGGGCACTGACTGAGAC

321 L L S R N P R G F F L F V E G G R I D H G H H E S R A Y R A L T E T

1501 GATCATGTTGACGACGCCATTGAGAGGGCGGCCAGCTACCAGCGAGGAGGACACGCTGAGCCTGCTACTGCCGACCACTCCCAGTCTTCTCCTTC

354 I M F D D A I E R A G Q L T S E E D T L S L V T A D H S H V F S F

1601 GGAGGCTACCCCTGCGAGGGAGCTCCATCTTGGGCTGGCCCTGGCAAGGCCGGGACAGGAAGGCTACACGGTCTCCTATACGAAACGGTCCAG

388 G G Y P L R K G S S I F G L A P G K A R D R K A Y T V L L Y G N G P

1701 GCTATGTCTCAAGGACGGCGCCCGCGGATGTTACCGAGAGCGAGAGCGGGAGCCCGAGTATCGGCAGCAGTGCAGGCTGCCCTGGACGAAGAGAC

421 G Y V L K D G A R P D V T E S E S G S P E Y R Q Q S A V P L D E E T

1801 CCACGAGGCGAGGACGTGGCGGTGTTCCGCGCGGCCCGCAGGCGCACCTGGTTCACGGCGTGCAGGAGCAGACCTTCATAGCGCACGTATGGCCTTC

454 H A G T G E D V A V F A V C A H L V H G V Q E Q T F I A H V M A F

1901 GCCGCTGCTGGAGCCCTACACCGCTGCGACCTGCGCCCGCCCGCCGACGCGCCGCGGCGGCTTCCAGTCCAAAGGCTCTGATG

488 A A C L E P Y T A C D L A P P A G T T D A A H P G R S R S K R L D

NheI (2004)

2001 GAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATG

521 TATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAAATTGCATTCATTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGT

2201 TTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATC

2301 CTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGT

SspI (2483)

2401 GTATTTTCCCAAGTTTGAAGTACTAGCTTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTCAGAAATAATTTAA

2501 ATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTGTAGTGTGGACTTAGGGAA

2601 CAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCTCAGTCTGCTCCTCTGCCACAAAGTGCAGCAGTTGCCGGCCGG

125 D Q E E A V F H V C N G A P

2701 GTCGCGCAGGGCAACTCCCGCCCCACGGCTGCTCGCGGATCTCGGTCATGCGCCGCGCCGAGGCGTCCCGGAAGTTCGTGGACACGACCTCCGACCAC

110 D R L A F E R G W P Q E G I E T M A P G S A D R F N T S V V E S W

2801 TCGGCGTACAGCTGTCAGGCGCGCACCCACACCCAGGCGAGGCTGTTGTCGGCACCACCTGGTCTGGACCCGCTGATGAACAGGGTACGCTGCT

76 E A Y L E D L G R V W V W A L T N D P V V Q D Q V A S I F L T V D D

SgrAI (2911)

2901 CCCGGACCACCCGGCAAGTCTCTCCACGAAGTCCCGGAGAACCCGAGCCGGTCCGGTCCAGAACTCGACCGCTCCGGCGACGTCGCGCGGGTGGAG

43 R V V G A F D D E V F D R S F G L R D T W F E V A G A V D R A T L

3001 CACCGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCTGTGAGGAGGAAAGAGAAGGTTAGTACAATTGCTATAGTGTGATTATTATAC

10 V P V A S T L K A M

3101 TATGCAGATATACTATGCCAATGATTAATTGTCAAAGTGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAA

3201 AAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTAT

3301 AAAGATACAGGCGTTTTCCCTGGAAAGTCCCTCGTGCCTCTCCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCTTCGGGAA

3401 CGTGGCGCTTTCTCATAGCTACGCTGTAGGTATCTCAGTTCGGTGTAGGCTGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGAC

3501 CGCTGCGCTTATCCGGTAACTATCGTCTTGTGATCAACCCGGTAAGACACGACTTATCGCCACTGGCAGAGCCACTGGTAACAGGATTAGCAGAGCGA

3601 GGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTAC
3701 CTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGGTTGCAAGCAGCAGATTACGCGCAGAAAAAAA
3801 GGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACAT
3901 TTAAATCA