



EcoRI (23)

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

1 CCGCGCCGCTCGACGATATCTAGAAATTCGGATCCTGCAGGGCCACTAGTATAGGGCTGTCTGGAGCCACTCCAGGGCCACAGAAATCTTGTCTCTGAC

101 TCAGGGTATTTTGTCTTTTGTGTAATGCTCTTCTGACTAATGCAAACCATGTGTCCATAGAACCAGAAGATTTTCCAGGGGAAAAGGTAAGGA

201 GGTGGTGAAGTGTCTGGTCTGCCCTCCAGGGCTTCCCTGGGTTAAGAGCCAGGCAGGAAGCTCTCAAGAGCATTGCTCAAGAGTAGAGGGGGCCT

301 GGGAGGCCAGGGAGGGATGGGAGGGAAACCCAGGCTGCCCCCAACCATGTCCTCCACCCTCTCAACCTCCCTCCACGGCCTGGAGAGTGGG

BstXI (455)

401 ACCAGGTATGGAGGCTTGAGAGCCCTGTTGGAGGAAGCCACAAGTCCAGGAACATGGGAGTCTGGGAGGGGGCAAAGGAGGCAGGAACAGGCCATCA

501 GCCAGGACAGGTGTAAGGCAGGCAGGAGTGTCTGCTGGGAAAAGTGGGATCAAGCACCTGGAGGGCTCTCAGAGCAAAGACAAACTGAGGTCC

Acc65I (661) **NcoI (682)**

601 CTGCCACTCTACAGAGCCCCACGCCGCCAGCTATAAGGGCCATGCACCAAGCAGGGTACCCAGGCTGCAGAGGTGCCATGGAAATCAAGGTGCT
 1► M E I K V L

BglII (797)

701 GTTGGCCCTCATCTGTATTGCTGTGTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGTGTGCCCTCCAACCTTTGCCACCACAGAT
 6► 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 N F A T T D

801 CTTGAGACTGACCTGTTCCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCA
 40► 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 R G K L P G

901 AAAAACTCCCCCAGATGTCTGAGGGAGCTGGAGGCCAATGCCAGAAAGGCTGGTTGCACAAGAGGCTGCCTCATTTGCCTCTCCACATTAAGTGCAC
 73► 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 L S H I K C T

1001 CCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGTGAAGAGGAGTCTGCTCAGGGAGGATTGGAGAGGCAATTTGTTGATATCCCA
 106► 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 A I V D I P

1101 GAGATTCCTGGCTCAAGGATAAGGACCACTGGACAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCA
 140► 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 C L K G L A

1201 ATGTCCAGTGTCTGACCTCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTTTCCAGCAAGATTGAGGGTAGGGTGGACAAAATCAAGGGTCTGGC
 173► 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 D K I K G L A

NheI (1316)

1301 TGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTG
 206► G D R •

1401 AAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGA

1501 GGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAAATTAATCTAAAATACAGCATAGCAAACTTTAACCTCCAATCAAGCC

1601 TCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTTATGGAGT
 1701 TTAAGATATAGTGTATTTTCCCAAGTTTGAAGTCTTCTTATTTTATGTTTAAATGCACTGACCTCCACATTCCTTTTATGATAAATATTCA
 1801 GAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTATAAATATCCCCAGTTTATGATGTT
 1901 GGACTTAGGGAACAAGGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCTGCCACAAAGTGCACGCA
 125► D Q E E A V F H V C

2001 GTTGGCGCCGGTCCGCGAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGTGCATGGCCGGCCGGAGGCTCCCGAAGTCTGTGGACAGC
 114► 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 R F N T S V

2101 ACCTCCGACCACTCGGGTACAGCTGCTCCAGGCCGCGCACCCAGGCCAGGGTGTGTCCGGCACCACTGGTCTGGACCGCGCTGATGAACA
 80► 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 Q V A S I F L

SgrAI (2223)

2201 GGGTACGTCGTCGGGACACACCGGCAAGTCTCTCCACGAAGTCCCGGAGAACCAGGCGGTGCGTCCAGAACTCGACCGCTCCGGCGACGTC
 47► 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 V A G A V D

2301 CGCGCGGTGAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCTGTGAGGAGGAAAGAGAAGGTTAGTACAATTGCTATAGTGA
 14► R A T L V P V A S T L K A M

2401 GTTGATATACTATGCAGATATACTATGCAATGATTAATTGTCAAAGTGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCC

2501 AGGAACCGTAAAAAGCCGCTGTTGCTGGCTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACC

2601 CGACAGGACTATAAGATACCAGGCTTTCCCTGGAAGTCCCTCGTGCCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCT

2701 CCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCC

2801 GTTCAGCCCAGCGCTGCCTTATCCGGTAACTATCGTCTTGTGTCACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGA

2901 TTAGCAGAGCGAGGTATGAGGGGTGCTACAGAGTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCT

3001 GAAGCCAGTTACCTTCGAAAAAGAGTGTGAGTCTTGTATCCGCAAAACACCCGCTGGTAGCGGTGTTTTTTGTTGCAAGCAGCAGATTACG

3101 CGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTA

3201 GTTAATTAACATTTAAATCA