



EcoRI (23)

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

1 GCGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGGCCCACTAGTTGATTTCCTTCATCCCTGGCACAC

75 GTCCAGGCAGTGTCAATCCATCTCTGCTACAGGGGAAAAACAAATAACATTTGAGTCCAGTGGAGACCGGGAGC

149 AGAAGTAAAGGGAAGTGATAACCCCCAGAGCCCGGAAGCCTCTGGAGGCTGAGACCTCGCCCCCTTGC GTGAT

BbrPI (266)

223 AGGGCTACTCGGAGCCACATGACCAAGGCACTGTCGCCTCCGCACGTGTGAGAGTGCAGGGCCCAAGATGGCTG

297 CCAGGCCTCGAGGCTGACTCTTCTATGCTACTTCCGTACCGGCAGAAAGGGGGCCCTCCAGCCAATGAGGC

371 TGGGGGCGGGCCTTACCTTGATAGGCACTCGAGTTATCCAATGGTGCCTGCGGGCCGGAGCGACTAGGAAT

SfiI (479)

445 AACGTCATGCCGAGTTGCTGAGCGCCGGCAGCGGGGGCCGAAACCAATGCGATGGCCGGGGCGGAG

NcoI (573)

519 TCGGGCGCTCTATAAGTTGCTGATAGGCGGGCACTCCGCCCTAGTTTCTAAGGACCATGGAAATCAAGGTGCTG

1 M E I K V L

593 TTTGCCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCACTGAAATCAATGAAGACCTCAATATAGCTGCTGT

7 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

BglII (688)

667 GGCCTCCAACCTTTGCCACCACAGATCTTGAGACTGACCTGTTCCACCACTGGGAGACCATGAATGTGATTAGCA

31 A S N F A T T D L E T D L F T N W E T M N V I S

741 CTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGCTCTG

56 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

815 AGGGAGCTGGAGGCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTTGCCCTCCACATTAAGTG

81 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

889 CACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAG

105 T P K M K K F I P G R C H T Y E G E K E S A Q G

963 GGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTATT

130 G I G E A I V D I P E I P G F K D K E P L D Q F I

1037 GCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCT

155 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

1111 CCTGAAGAAGTGGCTTCCCCAGAGGTGTACCACTTTTCCAGCAAGATTGAGGGTAGGGTGGACAAAATCAAGG

179 L K K W L P Q R C T T F A S K I Q G R V D K I K

NheI (1207)

1185 GTCTGGCTGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAA

204 G L A G D R •

1259 CTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGC

1333 TGCAATAAACAAGTTAACAACAACAATTGCATTCAATTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTT

1407 TTAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAAACTTTAACCTC

1481 CAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTG

1555 CATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAACCTAGC

SspI (1686)

1629 TCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAAATATTCAGAAATAATTTA

1703 AATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCC

1777 CCCAGTTTGTAGTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTA

1851 GCTTATCTCAGTCTGCTCCTCTGCCACAAAGTGCACGAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCG

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

1925 CCCCCACGGCTGCTCGCGATCTCGGTCTGCGCCGGCCGGAGGCGTCCCGGAAGTTCGTGGACACGACCTCCG

103 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

1999 ACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACCCAGGCCAGGGTGTGTCGGCACCACTGTGCTC

78 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 (2114)

2073 TGGACCGCGTGTGAACAGGGTACGTCGTCGCCGACCACACCGCGAAGTGCCTCCACGAAGTCCCGGGA

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

2147 GAACCCGAGCCGGTCGGTCCAGAACTCGACCGCTCCGGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGG
29 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
2221 TCAACTTGGCCATGATGGCTCCTCCTGTCAGGAGAGAAAGAGAAGAAGGTTAGTACAATTGCTATAGTGAGTT
4 L K A M
2295 GTATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAACTAGGGCTGCAGGTTAATTAAGAACATGT
2369 GAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCC
2443 CCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGC
2517 GTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTC
2591 TCCCTTCGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTCCGCTC
2665 AAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTC
2739 CAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAG
2813 GCGGTGCTACAGAGTTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCT
2887 CTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGG
2961 TGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTA
3035 CGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAA
3109 TCA