



EcoRI (23)

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

1 GCGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGGCCCACTAGTGACCTTTATTTCATAAGAGATGATG
75 TATTCTTGATACTACTTCTCATTGCAAATTCGAATTATTATTAATTTTCATATCAATTAGAATAATATATCTTC

BsrGI (177)

149 CTTC AATTTAGTTACCTCACTATGGGCTATGTACAACTCCAAGAAAAAGTTAGTCATGTGCTTTGCAGAAGAT

HindIII (225)

223 AAAAGCTTAGTGTA AACAGGCTGAGAGTATTTGATGTAAGAAGGGGAGTGGTTATATAGGTCTTAGCCAAAAC
297 ATGTGATAGTCACTCCAGGGTTGCTGGAAAAGAAGTCTGTGACACTCATTAACCTATTGGTGCAGAATTTGAA

NcoI (385)

371 TGATCTAAAGGAGACCATGGAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCA
M E I K V L F A L I C I A V A E A K P

BglII (500)

445 CTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTTGCCACCACAGATCTTGAGACTGACCTG
20 T E I N E D L N I A A V A S N F A T T D L E T D L
519 TTCACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAA
45 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
593 GCTGCCTGGCAAAAACTCCCCCAGATGCTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAA
69 L P G K K L P P D V L R E L E A N A R R A G C T
667 GAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCAC
94 R G C L I C L S H I K C T P K M K K F I P G R C H
741 ACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCCAGAGATTCCTGG
119 T Y E G E K E S A Q G G I G E A I V D I P E I P G
815 CTTC AAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTC
143 F K D K E P L D Q F I A Q V D L C A D C T T G C
889 TGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCCAGAGGTGTACCACTTTTGGC
168 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

NheI (1019)

963 AGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATG
193 S K I Q G R V D K I K G L A G D R
1037 ATAAGATACATTGATGAGTTTGGACAAACCACA ACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTG
1111 TGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACA CAATTGCATTCAATTTTA
1185 TGTTTCAGGTT CAGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTA

1259 ATTCTAAAATACAGCATAGCAAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAA
1333 TAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTTCATGGAGTTTA
1407 AGATATAGTGTATTTTCCAAGGTTTGAAGTACTGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACA

SspI (1498)

1481 TTCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGC
1555 AGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTGTAGTGTGGACTTAGGGAAACAAAGGAACCTTT
1629 AATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCCTGCTCCTCTGCCACAAAGTGCACGC

125 D Q E E A V F H V C

1703 AGTTGCCGGCCGGGTGCGCGAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGGCCCC
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
1777 GAGGCGTCCCGGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACAC
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
1851 CCAGGCCAGGGTGTGTCCGGCACCACCTGGTCTGACCGCGCTGATGAACAGGGTCACGTCGTCGCCGACCA
65 W A L T N D P V V Q D Q V A S I F L T V D D R V V

SgrAI (1926)

1925 CACCGGCGAAGTCGTCCTCCACGAAGTCCCGGGAGAACCCGAGCCGGTCCGGTCCAGA ACTCGACCGCTCCGGCG
40 G A F D D E V F D R S F G L R D T W F E V A G A
1999 ACGTCGCGCGGGT GAGCACCGGAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCCTGT CAGGAGAGGAAA
15 V D R A T L V P V A S T L K A M
2073 GAGAAGAAGGTTAGTACAATTGCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCAATGATTAATTG

2147 TCAAACTAGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGC
2221 CGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGT
2295 GGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCG
2369 ACCCTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTG
2443 TAGGTATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCCGACC
2517 GCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCC
2591 ACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGG
2665 CTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCT
2739 CTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAA
2813 AAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGG
2887 GATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA