



EcoRI (23) SdaI (38)
NotI (2) XbaI (19) BamHI (29) SpeI (45)
1 GCGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCGAGGGCCACTAGTCGGGAGCGCTACTTCTTCCGAGTGAGAGACAGAAAGAGAGGACCCGAGT
101 CTCACAGCCCTGAGGAACTGACACGCAGACCCCACTCCAGTCCCGGGGGCCCAACGTGAGGGGAGGACCTGGACGTTACCGGCGAAACGGTTTCCA
PvuII (228)
201 GGTGAGAGGTCACCCGAGGGACAGGCAGCTGCTCAACCAATAGGACCAGCTCTCAGGGCGGATGCTGCCTCTCATTGGCGCCGTTAAGAATGACCAGTA
BstBI (357) BstXI (379)
301 GCCAATGAGTCGGCCTGGGGGGCTACCAGTGACGTGAGTTGCGGAGGAGCCGCTTCGAATCGGCAGCGCCAGCTTGGTGGCATGAACCAACCAGCGG
401 CCTCCAACGAGTAGCGAGTTCACCAATCGGAGGCCTCCACGACGGGGTTCGGGGGAGGATATAAGCCGAGTCGGCGACCGCGCTCGATACTGGCT
501 GTGACTACACTGACTTGGACACTTGGCCTTTTGCGGGTTTGTAGAGGTAAGCGTCGCGGCTGCTTCCAGGCCTACCCTGATTTTGGTTCGTGGCTCCTCC
601 TGACCCTGAGACCTCTGTGCGCCTCAGATCAGAACCGTCGTCGCGTTTCGGGGCTACAGCCTGTTGCTGGACTCTGTGAGACACCTGACCGACCGCTGAG
NcoI (727)
701 CGACTGACTGGTCCACAGCGCCGGCACCATGGAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAA
M E I K V L F A L I C I A V A E A K P T E I N E
BglII (842)
801 GACCTCAATATAGCTGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACCATGAATGTATTAGCACTGACA
25▶ D L N I A A V A S N F A T T D L E T D L F T N W E T M N V I S T D
901 CAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCTGGCAAAAACTCCCCCAGATGCTCCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGG
58▶ T E Q V N T D A D 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
1001 TTGCACAAGAGGCTGCCTCATTTGCCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCTGGCAGGTGCCACACTTATGAAGGTGAAAAG
91▶ C T 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 T Y E G E K
1101 GAGTCTGCTCAGGGAGGATTGGAGAGCAATTGTTGATATCCAGAGATTCTGGCTCAAGGATAAGGAGCCACTGGACCAAGTTTATTGCTCAAGTGG
125▶ E S A Q G 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 A Q V
1201 ACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTCCAATGTCCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTT
158▶ D L C A D C T T G 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
NheI (1361)
1301 TGCCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTG
191▶ A S K I Q G R V D K I K G L A G D R •
1401 GACAAACCACAAC TAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAACAAGT
1501 TAACAACAACAATTGCATTCAATTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATTA
SspI (1840)
1601 ATTCTAAAATACAGCATAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTG
1701 CCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTATTTTCCCAAGGTTTGAAGTAGCTCTTCAATTTCTTTATGTT
1801 TTAAATGCACTGACCTCCACATTCCTTTTTAGTAAAATATTAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAA
1901 TCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCT
2001 TCTAGCTTATCCAGTCTGCTCCTCTGCCACAAAGTGCACGAGTTGCCGCGCGGTGCGCGAGGGCGAACTCCCCCCCCACGGCTGCTCGCCGATC
125◀ • D Q E E A V F H V C N G A P D R L A F E R G W P Q E G I
2101 TCGGTGATGGCCGCGCGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGCGGTACAGCTCGTCCAGGCGCGCACCCACACCCAGGCCA
95◀ E T M A P G 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 W A L
SgrAI (2268)
2201 GGGTGTGTCCGGCACCACCTGGTCTGGACCGCGCTGATGAACAGGGTACGTCGTCGCCGACCAACCGCGAAGTCTCCTCCACGAAGTCCCGGGA
62◀ T N D P V V Q D 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
2301 GAACCCGAGCCGGTCCGAGAACTGACCGCTCCGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGGTCAACTGGCCATGATGGCTCCTCCT
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 L K A M ◀
2401 GTCAGGAGAGGAAAGAGAAGGTTAGTACAATTGCTATAGTGAGTTGATTATACTATGCAGATATACTATGCAATGATTAATTGTCAA ACTAGGGC
2501 TGAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTG
2601 ACGAGCATCAAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCGTTTTCCCCTGGAAGCTCCCTCGTGGCTC
2701 TCCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGATAGTATCTCAGTTCG
2801 GTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGACGAACCCCGTTAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTCCAACCCGG
2901 TAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCTAA
3001 CTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAAACAAAC
3101 ACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGGTCTGACG
3201 CTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA