



**EcoRI (23) SdaI (38)**  
**NotI (2) XbaI (19) BamHI (29) Acc65I (75)**

1 GCGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGGCCCTGAAATAACCTCTGAAAGAGGAACTTGGTTAGGTACCTTCTGAGGCTGAAAGAACCA  
101 GCTGTGGAATGTGTGTCAGTTAGGGTGTGAAAGTCCCCAGGCTCCCCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCAGGTGT  
201 GAAAGTCCCCAGGCTCCCCAGCAGGCAGAAGTATGCAAAGCATGCATCTCAATTAGTCAGCAACCATAGTCCACTAGTTCCAGATGGTAAATATACAC  
301 AAGGATTTAGTCAAACAATTTTTTGGCAAGAATATTATGAATTTGTAATCGTTGGCAGCCAATGAAATACAAAGATGAGTCTAGTTAATAATCTACA  
401 ATTATTGGTTAAAGAAGTATATTAGTCTAATTTCCCTCGGTTTGTCTAGCTTTTCTCTTCTGTCAACCCACACGCTTTGGCACCATGGAAATCAAG  
501 GTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTGCACCA  
5 V L 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  
**BglIII (602)**  
601 CAGATCTTGAGACTGACCTGTTCCACCAACTGGGAGACCATGAATGTGATTAGCACTGCACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCC  
38 T D 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  
701 TGGCAAAAACCTCCCCCAGATGCTCTGAGGGAGCTGGAGGCCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTGCTCTCCACATTAAG  
71 G 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  
**MfeI (889)**  
801 TGCACCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATA  
105 C T 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  
901 TCCAGAGATTCTGGCTCAAGGATAAGGAGCCACTGGACAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCT  
138 I P 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  
1001 TGCCAATGTCAGTCTGCTGACCTCCTGAAGAAGTGGCTTCCCAGAGGTGTACCACCTTTTCCAGCAAGATTGAGGGTAGGGTGGACAAAATCAAGGGT  
171 A 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  
**NheI (1121)**  
1101 CTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTGGACAAACCACAACCTAGAATGCAGTAAAAAATGCTTTAT  
205 L A G D R •  
**MfeI (1270)**  
1201 TTGTGAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTCAATAAAACAGTTAACACAACAATTGCATTCTTTTATGTTTCAGGTTTCAG  
1301 GGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCCAAATC  
1401 AAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTTCAT  
1501 GGAGTTTAAAGATATAGTGTATTTCCCAAGGTTTGAAGTCTCTTCTTTATGTTTAAATGCACTGACCTCCCACATTCCTTTTATGTAATAAT  
1601 ATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATAGGCAGAATCCAGATGCTCAAGGCCCTTCAATAATCCCCAGTTTAG  
1701 TAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCTGCCACAAGTGC  
125 • D Q E E A V F H  
1801 ACGCAGTTGCCGCGGGTGCAGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTCATGGCCGCGCCGAGGCGTCCCGGAAAGTTCGTGG  
115 V C 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  
1901 ACACGACCTCCGACCACTCGGCTACAGCTCGTCCAGGCCGCGCACCCAGCCAGGGTGTGTTCCGGCACCACTGGTCTGGACCGCTGAT  
82 V 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  
**SgrAI (2028)**  
2001 GAACAGGGTACGTCGTCGCCGACACACCGGCGAAGTCTGCTCCACGAAGTCCCGGAGAACCAGCCGGTCCGAGAACCTGACCGCTCCGGCG  
49 F L 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  
**MfeI (2191)**  
2101 ACGTCGCGCGGTTGAGCACCAGGCAACGGCACTGGTCAACTTGGCCATGATGGCTCCTCCTGTCAGGAGAGGAAAGAGAAGAAGGTTAGTACAATTGCTAT  
15 V D R A T L V P V A S T L K A M  
2201 AGTGAGTTGATTATACTATGAGATATACTATGCCAATGATTAATTGTCAAAGTGGGCTGACAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAA  
2301 AGGCCAGGAACCGTAAAAAGCGCGTGTGCGGCTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCG  
2401 AAACCCGACAGGACTATAAAGATACAGGCGTTTCCCCTGGAAGCTCCCTCGTGGCTCTCCTGTCCGACCTGCCGTTACCGGATACCTGTCCGCC  
2501 TTTCTCCCTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTGCGGTAGGTGTTGCTCAGGCTGGGCTGTGTGCACGAAC  
2601 CCCCCTTACGCCGACCGCTGCGCCTTATCCGGTAATATCGTCTTGGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAA  
2701 CAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCT  
2801 CTGCTGAAGCAGTTACCTTCCGAAAAAGAGTTGGTAGCTCTTGTATCCGGCAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGA  
2901 TTACGGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTACGGGGTCTGACGCTCAGTGAACGAAACTCACGTTAAGGGATTTTGGTCAT  
3001 GGCTAGTTAATTAACATTTAATCA