



**SdaI (38)**

**NotI (2) Sall (9) XbaI (19) BamHI (29)**

1 GCGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAAGGCTTACATTTGAATGTTGGCTACATATGTATGAGTTTTCAACTCCAGGAGAAAACGTCTC  
101 TTTAAAAGAGAACAACAAAAGCTAACAGAAAATACAAGTGTGACATTGGCCTTAGTTCGACCAAGAAAGCAATTCATCTTGTTCCTTTGTGGTAT  
201 ACAGATAAAGAAAAATAAATCACTACAACGAAAGCAAAATCTCTTCAGCGTCTCTAATACATCTTCCAAATCAGTGTGTCTGACCTTTTCTTAAGACTT  
301 TAACCATCACAAAGAAACAGTGGGGAGGGAGTGTGCTGCCTAGTAGTTAAAGGGCAGGAGAATTCAGTGGTGTGAGAAGGGATTAGTGAGAGCTGG  
401 AAGAGAGGACCAGCCCCTCCAGTGTGAGGAATCTGGCTTGGGATTACTGTCTGGCAGAAAATCTCTTCGGGCAATTAACAGCTGGCATCAGGGGAAAA  
501 GCAGACATCCAACAACACTAGCTCTGAAGGAGATCAGCAGAGAAAACCTTCAGGGATTGACTGGTACTGGTGTGAGCAGCTCTGTGgtgggtaccctgtgacc

**NcoI (665)**

601 aaagctctaggaacatgaaggagatttgccttgctataaacctgtttcctattctcctttcatttccatgggtaactattactatggtagtccaactag  
701 tggatgcttttgtaaatgacatctatggaaagtcttttggatcagggtgatctttttatgtatgtatgtgatggatgggtgcacgagagcagg  
801 tgcccagattctcaaggaggcttcagttaacaaggagtgggagtgtatgtatgtggtgcaaggcactgaagtcagtctctctgtaagagcactctatg  
901 ctctttaccactgtgccttctccccagccaagaatagtattcttatgggtagaaatttaataagaaaactcaaagaccaggagagtgtgattctgtca  
1001 tctagctattatgcctgcagatatttaaggtgaataattgttttgactattgttttagaaatgttttcacatgaaagattccatttccggagtgggtt  
1101 gaaaaagtatgcaaaagaacttttgcaactctgtttttgcctttctgtttttcagCTGATTTTCATCTGAGCACCCCTGTCTTCTCCATGCAAAGAGCA

**NcoI (1239)**

1201 GCATAGGAGACCTGTGTTCTGAACTCTTGCTTCGAGAACCATGGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCCACT  
1301 GAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCAACTTTGCCACCACAGATCTTGAGACTGACCTGTTCCACCAACTGGGAGACCATGAATGTGA  
21▶ E I N E 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  
1401 TTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGTCCTGAGGGAGCTGGAGGCCAATGC  
54▶ I S 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 R E L E A N A  
1501 CAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTAT  
87▶ R R A G 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  
1601 GAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCCAGAGATTCCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTA  
121▶ E G E K 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  
1701 TTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAAATGTCCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCCAGAG  
154▶ I 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 L K K W L P Q R

**NheI (1873)**

1801 GTGTACCACCTTTTCCAGCAAGATTCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACA  
187▶ C T T F A S K I Q G R V D K I K G L A G D R •  
1901 TTGATGAGTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTG  
2001 CAATAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGT  
2101 GGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCAT  
2201 CAGGGGCTGTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTCTTCA  
2301 TTTCTTTATGTTTAAATGCACTGACCTCCACATTCCCTTTTATGATAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTT  
2401 TATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTGTGGACTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCA  
2501 AGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTGCTGCCACAAAGTGCACGAGTTGCCGGCCGGTGCAGGCGGAACTCCCGCCCCACGGC  
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  
2601 TGCTCGCCGATCTCGGTGATGGCCGGCCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCGCGCACCC  
99◀ Q E G I 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  
2701 ACACCCAGGCCAGGGTGTGTCGGCACCACTGGTCTGGACCGCTGATGAACAGGGTACGCTCGTCCGGACCAACCCGCGAAGTGTCTCCAC  
66◀ V 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 G A F D D E V  
2801 GAAGTCCCGGAGAACCAGCCGGTCCGAGTCCGAGTCCGACCGCTCCGGCAGCTCGCGCGGTTGAGCACCAGGACCGGACTGGTCAACTTGGCCATG  
33◀ F D R S 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  
2901 ATGCTCCTCTGTGAGGAGGAAAGAGAAGAAGTTAGTACAATTGCTATAGTGTATTATACTATGAGATATACTATGCCAATGATTAATTG  
3001 TCAAAGTACTAGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGCCGCGTGTGGCGTTTTTCCATAGGC  
3101 TCCGCCCCCTGACGAGCATCACAATAATCGACGCTCAAGTCAAGGTGGCGAAACCCGACAGGACTATAAGATAACCAGGCGTTTTCCCCTGGAAGCTC  
3201 CCTCGTGCCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGTTTTCTCATAGCTCAGCTGTAGG

3301 TATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTTCAGCCCGACCGCTGGCCTTATCCGGTAACTATCGTCTTG  
3401 AGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAA  
3501 GTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCC  
3601 GGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTA  
3701 CGGGGTCTGACGCTCAGTGAACGAAAACCTACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA