



SdaI (38)

1 CGGCGCGCTCGACGATATCTAGAATTCGGATCCTGCAGGGCCACTAGTGTGGATTACAGGTGTGAGCTATTGTCCCAGCCAAAAGGAAAAGTTT
101 ACTGTAGTAACCTTCCGGACTAGGGACCTCGGCCTCAGCCTCAGGTACCTAGGTGCTTTAGAAAGGAGGCCACCCAGGCCATGACTACTCTTGGC
201 ACAGGGAGCCCTGCACACAGATGTGCTAAGCTCTCGTCCAGCCAGAGGGAGGAGGCTGAGCCAGTCAGAAGGAGATGGGCCCCAGAGAGTAAGAAA
301 GGGGGAGGAGGCCAAAGCTGATCCAAAAGGTGGTCTAAGCAGTCAAGTGGAGGAGGTTCCAATCTGATGGCGGAGGGCCAAAGCTCAGCCTAACGAG
401 GAGGCCAGGCCACCAAGGGGCCCTGGAGACTTGTTCCCTTGCCCTTGTGGTTTTTGCATTTCTGTCCCTTGTCTCATTGCGGAAGTTCTT

NcoI (552)

501 CTTCTTACCTGCACCCAGAGCCTGCCAGAGAAGACAAGGGCAGAAAGCACCATGGAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGA
601 GGCAAAACCCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTGGCCACCACAGATCTTGAGACTGACCTGTTCAACCACTGGGAG
16▶ A K P 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 F T N W E
701 ACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCTGGCAAAAACTCCCCAGATGCTCCTGAGGGAGC
50▶ T M N V 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
801 TGGAGGCCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTGCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCTGGCAG
83▶ L E A N A 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
901 GTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTCTGGCTTCAAGGATAAGGAGCCA
116▶ C H 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 F K D K E P
1001 CTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTGAAGGGCCTTCCAATGTCCAGTGTCTGACCTCCTGAAGAAGT
150▶ L D Q F 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

NheI (1186)

1101 GGCTTCCCAGAGGTGTACCACTTTTGCAGCAAGATTGAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGTGGCCAGAC
183▶ W L P Q R C T T F A S K I Q G R V D K I K G L A G D R •
1201 ATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTGTGAAATTTGTGATGCTATTGCTTTATTTGTAA
1301 CCATTATAAGCTGCAATAACAAGTAAACAACAACAATTGCATTCATTTTATGTTTACAGTTCAGGGGAGGTGTGGGAGTTTTTTAAAGCAAGTAAAA
1401 CCTCTACAAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAAT
1501 AAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTCTTTTCATGGAGTTTAAAGATATAGTGATTTTTCCCAAGGTTTG
1601 AACTAGCTCTTCAATTTCTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAA
1701 AATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAG
1801 AAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCAGCTCTGCTCTGCCACAAAGTGACGCAGTTGCCGGCCGGTCCGCGAGGGCGAACTC
125▶ • D Q E E A V F H V C N G A P D R L A F E
1901 CGCCCCCACGGTGTCTCGCGATCTCGGTCTATGGCCGGCCGAGGCTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCC
104▶ R 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 W E A Y L E D
2001 AGGCCGCGACCCACACCAGGCCAGGGTGTGTCCGGCACCCTGGTCTGGACCGCGTGATGAACAGGGTCACGTCGTCGCGGACCACACCGGCGA
70▶ L G R V W 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
2101 AGTCGCTCCACGAAGTCCCGGGAACCCGAGCCGGTCCGAGTCCGAGAACTCGACCGCTCCGGCGACGTCGCGCGCGGTGAGCACCGGAACGGCACTGGT
37▶ D D E V 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
2201 CAACTGGCCATGATGGCTCCTCTGTGAGGAGGAAAGAGAAGAAGTTAGTACAATTGCTATAGTGAGTTGATTATACTATGCAGATATACTATGC
4▶ L K A M
2301 CAATGATTAATTGTCAAAGTGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCG
2401 TTTTCCATAGGCTCCGCCCTGACGAGCATCAAAAAATGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCGTTTC
2501 CCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTCTCCCTTCCGGAAGCGTGGCGCTTCTCATAG
2601 CTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCGACCGCTGCGCTTATCCGGT
2701 AACTATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTA
2801 CAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAGGAGTTGG
2901 TAGCTCTTGATCCGGCAAAACCAACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCT
3001 TTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA