



Sdal (38)

1 CGGGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGTACATGTTCTGCCAGCTGACTTTGTCAAGACAGTGATGTCTGTGTTCCAGCAGTTGTTC
101 TGAGTATCCTTTTCATTATCCACTGTCCTTTCTTCTTAAATTCACCCCCAACATTGTAATAGCTTCTTTCTTAAACTCTGTTCAAAGAACCAGCTTGA

SpeI (238)

201 GTGTGTCAGCTGCTTCCTGCTGGGGCTCGCAACCACTAGTGACCTTTATTATAAGAGATGATGATTCTTGATACTACTTCTCATTGCAAATTC
301 AATTATTATTAATTTTCATATCAATTAGAATAATATATCTTCTTCAATTTAGTTACCTCACTATGGGCTATGTACAACTCCAAGAAAAAGTTAGTCATG
401 TGCTTTGCAGAAGATAAAAGCTTAGTGTAACAGGCTGAGAGTATTTGATGTAAGAAGGGGAGTGTTATATAGGCTTAGCCAAAACATGTGATAGTC

NcoI (578)

501 ACTCCAGGGGTTGCTGGAAAAGAAGTCTGTGACACTCATTAACTATTGGTGAGAAATTTGAAATGATCTAAAGGAGACCATGGAAATCAAGGTGCTGTTT
601 GCCCTCATCTGTATTGCTGTTGCTGAGGCAAAACCACTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCTCCAACCTTTGCCACCACAGATCTTG
8) 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 T D L
701 AGACTGACCTGTTCCAACTGGGAGACCATGAATGTATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAA
41) 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 G K K
801 ACTCCCCAGATGCTCCTGAGGGAGCTGGAGGCAATGCCAGAAGGGCTGGTTGCACAAGAGGCTGCCTCATTGGCTCTCCACATTAAGTGCACCCCT
74) 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 C T P
901 AAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAGGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGA
108) 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 I P E
1001 TTCCTGGCTTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGTCTGAAGGGCTTGCCAATGT
141) 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 A N V
1101 CCAGTGTCTGACCTCCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTTTCCAGCAAGATTACAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGG
174) 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 L A G

NheI (1212)

1201 GACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAACCAACTAGAAATGCAGTGAAAAAATGCTTTATTTGTGAAAT
208) D R •
1301 TTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCAATTTATGTTTCAGGTTACAGGGGAGGTG
1401 TGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAAATTAATTCTAAAATACAGCATAGCAAAACTTTAACCTCAAATCAAGCTCTA
1501 CTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTTAA
1601 GATATAGTGTATTTTCCAAAGTTTGAAGTACTGCTCTTCATTTCTTTATGTTTTAAATGCAGTACCTCCACATTCCTTTTATAGTAAAATATTAGAAA
1701 TAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGAC
1801 TTAGGGAACAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGTCTGCTCCTCTGCCACAAGTGCACGCAGTTG
125) • D Q E E A V F H V C N
1901 CCGGCCGGTTCGCGCAGGGCGAACTCCCGCCCCACGGCTGCTCGCCGATCTCGGTATGGCCGGCCGGAGGCGTCCCGAAAGTTCGTGGACACGACCT
112) G A P D R L A F E R G W P C E G I E T M A P G S A D R F N T S V V E
2001 CCGACCACTCGGCGTACAGCTCGTCCAGGCGCGCACCCACCCAGGCGTGTTCGGCACCACCTGGTCTGGACCGCGTGTGAACAGGGT
79) S W E A Y L E D L G R V V W A L T N D P V V Q D Q V A S I F L T
2101 CACGTCTGCCGACACACCGCGAAGTCTGCTCCACGAAGTCCCGGGAGAACCCGAGCCGGTCCGTCAGAACTCGACCGCTCCGGCAGCTCGCGC
46) 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 V D R
2201 GCGGTGAGCACCAGGACGGCACTGGTCAACTGGCCATGATGGCTCCTCTGTGAGGAGGAAAGAGAAGGTTAGTACAATTGCTATAGTGAGTTG
12) A T L V P V A S T L K A M
2301 TATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAAGTGGGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGA
2401 ACCGTAAGGAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGAC
2501 AGGACTATAAAGATACCAGCGTTTTCCCTGGAAGCTCCCTCGTGGCTCTCCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCT
2601 TCGGAAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTGCTCCAAGTGGGCTGTGTGCACGAACCCCGTTC
2701 AGCCCCACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAG
2801 CAGAGCGAGGTATGTAGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAG
2901 CCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCA
3001 GAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTA
3101 ATTAACATTTAAATCA