



PstI (7) SpeI (14)

1 CTCTGAGGGCCACTAGTGC... 101 GCGGATGTTATCTACCAT... 201 CGAATCGGCGGCGCCAG... 301 GTATATAAGCCGAGTAG... 401 GCGGCTGTATTCTAGAC...

NeoI (548)

501 CTACGGCTGTGGCTGG... 601 GACAGCAAATGGGTCGG... 701 CCCTGGCGTTACCCA... 801 CGCAGCCTGAATGG... 901 TCGTCCCCTCAAAC... 1001 TCCGACGGGTTGT... 1101 CTGTGGTCAACGG... 1201 CGGTGATGGTGTG... 1301 GACTACACAAATC... 1401 GACTACCTACGGG... 1501 ATGCCGATCGCGT... 1601 CGACGGCAGCTG... 1701 CGAGGCGTTAAC... 1801 ACGCCGTGGCGT... 1901 GCGCATGGTGCCA... 2001 AGTGTGATCATCT... 2101 AGTATGAAGGG... 2201 GTCCATCAAAAAT... 2301 TACTGGCAGGCG... 2401 GGTCCGGCTTAC... 2501 GGAAGCAAACACC... 2601 CTGCACTGGATG... 2701 TACCGCAGCCG... 2801 GCAGTGGCGTCT... 2901 GGTAAATAAGCG... 3001 CCCGTGCACCGT... 3101 CGAAGCAGCGT... 3201 AGCCGAAAAAC... 3301 ACTGCCAGCTGG... 3401 GGATCTGCCATT... 3501 CGCGGCGACTTC... 985 ArgGlyAspPheGlnPheAsnIeSerArgTyrSerGlnGlnGlnLeuMetGluThrSerHisArgHisLeuLeuHisAlaGluGluGlyThrTrpLeuA

3601 ATATCGACGGTTTCCATATGGGGATTGGTGGCGACGACTCCTGGAGCCCGTCAGTATCGGGGAATTACAGCTGAGCGCCGGTTCGCTACCATTACCAGTT  
1018▶ snI leAspGlyPheHisMetGlyI leGlyGlyAspAspSerTrpSerProSerValSerAlaGluLeuGlnLeuSerAlaGlyArgTyrHisTyrGlnLe

EcoRI (3733)

3701 GGTCTGGTGTCAAAAAATAATAATCTAGTCGAGAATTCGCTAGCTCGACATGATAAGATAACATTGATGAGTTTGGACAAAACCACAAGTGAATGCAGTGAA  
1051▶ uValTrpCysGlnLys•••

3801 AAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAA

3901 GTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTAGATCCA

PacI (4014)

4001 TTTAAATGTTAATTAACCTAGCCATGACCAAAATCCCTTAACGTGAGTTTTTCGTTCCACTGAGCGTCAGACCCGTAGAAAAGATCAAAGGATCTTCTTGA

4101 GATCCTTTTTTCTGCGCGTAATCTGTGCTTGCACAAACAAAAAACACCGCTACCAGCGGTGGTTTGTTCGCGGATCAAGAGCTACCAACTCTTTTTTC

4201 CGAAGGTAAGTGGCTTCAGCAGAGCGCAGATACCAAATACTGTTCTTCTAGTGTAGCCGTAGTTAGGCCACCACTTCAAGAACTCTGTAGCACCGCCTAC

4301 ATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCGTGCTTACCAGGTTGGACTCAAGACGATAGTTACCGGATAAGGGC

4401 CAGCGGTCGGGCTGAACGGGGGTTCTGTGCACACAGCCAGCTTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGAGCTATGAGAAAGCG

4501 CCACGCTTCCCGAAGGGAGAAAGGCGGACAGGTATCCGGTAAGCGGCAGGGTCGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGAAAACGCCTGGTA

4601 TCTTTATAGTCCTGTGGGTTTTGCCACCTCTGACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGGC

PacI (4754)

4701 GCCTTTTTACGGTTCCTGGCCTTTTGCTGGCCTTTTGCTCACATGTTCTTAATTAATTTTTCAAAAAGTAGTTGACAATTAATCATCGGCATAGTATATC

4801 GGCATAGTATAATACGACTCACTATAGGAGGGCCATCATGGCCAAGTTGACCAGTGTGCTCCAGTGTCTCACAGCCAGGGATGTGGCTGGAGCTGTTGA

4900 GTTCTGGACTGACAGGTTGGGGTCTCCAGAGATTTTGTGGAGGATGACTTTGCAGGTGTGGTCAGAGATGATGTCACCCTGTTTCATCTCAGCAGTCCAG

5000 GACCAGGTGGTGCCTGACAACACCCTGGCTTGGGTGTGGGTGAGAGGACTGGATGAGCTGTATGCTGAGTGGAGTGAGGTGGTCTCCACCAACTTCAGGG

5100 ATGCCAGTGGCCCTGCCATGACAGAGATTGGAGAGCAGCCCTGGGGGAGAGAGTTTGCCTGAGAGACCCAGCAGGCAACTGTGTGCACTTTGTGGCAGA

PacI (5263)

5200 GGAGCAGGACTGAGGATAAGAATTGTAACAAAAACCCCGCCCGGGGGTTTTTTGTTAATTAA