



EcoRI (23) PstI (38)

NotI (2) XbaI (19) BamHI (29)

1 CGCGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGCCTGAGGCAAAGGGAGTGGCTACAGCCTGGCACGGTCGATTAAGCCCTGCTCTCCGGTCC
101 TGGGACACTTTCTTTTTCTCTTTTGTAGTCACAGGTCCTCTAACATGAGAATCAAGTATTTTACGCTGATTTCTTATAAAATTGTGAGAACTCCAT
201 AGGCGATGTACCGCTACTCCTACCTTAACCGTGTGTAAAGACAGCAAAACAAATGAACTATACTGCAAGATCTCTTCTATTTCCCTATTCAAACCTAA
301 AATGAAGAGGGAGGGGAGACATGGACAAGCAAGCATTCCACAGGCGCCCTGCCAACGCTGTCACTCAAACCAGGACCCAATCACAGACTTTTTAGCC
401 AAGCCTTATCCGCTCTCTTGAGAACTTTCTGCGTCCGCCATCCTGTAGGAAGGATTTGTACACTTTAACTCCCTCCCTGGTCTGAGTCCCACACTC
501 TCACCACCAGCACCTTCAGGAGCTGACCTTAACAGCTTACCCACAGGACCCCGAAGTTGCGTGCCTCCGCAACAGTGTCAATAGCAGCACCAGCA
601 CTTCCCCACACCTCCCCCTCAGGAATCCGTACTCTTAGCGAACCCAGAACTCTGGAGAGTTCTGGACAAGGGCGGAACCCACAACCTCCGATTACT
701 CAAGGGAGGCGGGGAAGCTCCACCAGACGCGAACTGCTGGAAGATTCTGGCCCAAGGCTCTCCGGCTCGCTGATTGGCCAGCGGAGAGTGGGGC
801 GGGCCGGTGAAGACTCCTTAAAGGCGCAGGGCGGCGAGCAGGGCACCAGACGCTGACAGCTACTCAGAATCAAATCTGGTTCATCCAGAGACAAGCGAA
901 GACAAGAGAAGCAGAGCAGCGGCGGTTCCCGATCCTCGGCCAGGACCAGCCTTCCCAGAGCATCCACGCGCGGAGCGCAACCTTCCCAGGAGCATC

BglII (270)

BsrGI (459)

NcoI (1074)

1001 CCTGCCGCGGAGCGCAACTTTCCCGGAGCATCCACGCCGCGGAGCGCAGCCTTCCAGAAGCAGAGCGCGGCCCATGGAAATCAAGGTGCTGTTTGCCC
M E I K V L F A
1101 TCATCTGTATTGCTGTTGCTGAGGCAAACCCTGAAATCAATGAAGACCTCAATATAGCTGCTGTGGCCTCCAACCTTTGCCACCACAGATCTTGAGAC
9L 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 E T
1201 TGACCTGTTCACTACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGAACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTC
42D 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 L
1301 CCCCAGATGTCCTGAGGGAGCTGGAGGCAATGCCAGAAGGGTGGTTGCACAAGAGGCTGCCTCATTTCCTCCACATTAAGTGCAACCCCTAAGA
76P 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 K
1401 TGAAGAAATTTATCCTGGCAGGTGCCACACTTATGAAGTGAAGAGGAGTCTGCTCAGGAGGGATTGGAGAGGCAATTGTTGATATCCAGAGATTTCC
109M 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 I P
1501 TGGCTTCAAGGATAAGGACCACTGGACCACTTATTGCTCAAGTGGACCTCTGTGCTGATTGCACCACTGGCTGCTGAAGGGCCTTCCCAATGTCCAG
142G 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 Q
1601 TGCTCTGACCTCTGAAGAAGTGGCTTCCCAGAGGTGTACCCTTTTCCAGCAAGATTGAGGGTAGGGTGACAAAATCAAGGGTCTGGCTGGGGACA
176C 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 D

BglII (1189)

NheI (1708)

1701 GATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGT
209R
1801 GATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAAACAATGCATTCAATTTATGTTTCAGGTTTCAGGGGGAGGTGTTGG
1901 AGGTTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATTAATTCTAAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTTG
2001 AATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATA

SspI (2187)

2101 TAGTGATTTTTCCCAAGGTTTGAAGTAGCTCTTCAATTTCTTATGTTTTAAATGCACTGACCTCCCACATTCCTTTTTAGTAAAAATATTCAGAAATAAT
2201 TTAATACATCATTGCAATGAAAAATAATGTTTTTATTAGGCAGAAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAG
2301 GGAACAAAGGAACCTTTAATAGAAATTTGACAGCAAGAAAGCGAGCTTCTAGCTTATCCTCAGCTGCTCCTCTGCCACAAAGTGACGCGAGTTGCCG
125D Q E E A V F H V C N G A
2401 CCGGGTCCGCGAGGGCGAACTCCCGCCCCACGGTCTCGCCGATCTCGGTGATGGCCGGCCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGA
111P 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 V V E S
2501 CCACTCGCGTACAGCTCGTCCAGGCCGCGCACCCACACCAGGCCAGGGTGTGTCGCGCACCACTGGTCTGGACCGCGCTGATGAACAGGGTCCAG
78W 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 F L T V

SgrAI (2615)

2601 TCGTCCCGGACCACCCGGCAAGTCGCTCCACGAAGTCCCGGAGAACCAGCGGGTCCGCTCCAGAACTGACCGCTCCGGCGACGTCCGGCGCGG
44D 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 A T
2701 TGAGCACCAGGACCGCACTGGTCAACTTGGCCATGATGGCTCCTCCTGTGAGGAGGAAAGAGAAGAAGGTTAGTACAATTGCTATAGTGAGTTGATT
11L V P V A S T L K A M

PstI (2851)

2801 ATACTATGCAGATATACTATGCAATGATTAATTGTCAAAGTGGCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCG
2901 TAAAAAGGCCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAATAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGA
3001 CTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTCTCCCTTCGG
3101 GAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAGCTGGGCTGTGTGCACGAATCCCCCGTTACGCC
3201 CGACCGCTGGCCTTATCCGGTAACTATCGTCTTGTAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGA
3301 GCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAG
3401 TTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCCCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAA
3501 AAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCTAGTTAATTA

3601 ACATTAAATCA