



2701 TTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAAATAATGTTTTTTATT
SspI (2745) SwaI (2759)

2801 AGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAA

2901 AGCGAGCTTCTAGCTTTAGTTCTCTGGTGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTATCTCAATGAGCACAAAGCAG
141 •••AsnArgThr TyrLysLeuProl | eLeuGlu | eThr Thr LysVal | LeuLysGlyAsnMetGlu | eLeuVal | PheCysA

BstXI (3049)

3001 TCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCA
 112 spProAlaTyrAspSer | eLeuGluArgCysMetGlyCysProSerValValArg | eSerArgAspValGluAspSerTyrProHisArgValAlaVal

3101 CAATGGTGTCAAAGTCCTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTGACCTGCCAATGTAGGCCCAATGTGGAC
 79 | | eThrAspPheAspLysGlnGlyAsnSerValAlaSerGly | eAla | eAlaGluAlaCysValThrValArgGly | eTyrAlaGlu | eHisVal

3201 AGCAGAGATGATCTCCAGTCTTGGTCTGATGGCCGCCGACATGGTGTCTTGTCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCAC
 46 AlaSer | eLeuGluGlyThrLysThrArg | eAlaAlaGlyValHisHisLysAsnAspGluTyrLeuMetThr | eLysGluThrAlaValGluValL

XmnI (3326)
AseI (3392)

3301 AGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTG
 12 eGluLeuAspGlnGlnSer | eAsnPheThrLysMet

3401 TCAAACACAGCGTGGATGGCGTCTCCAGCTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTGCG

SpeI (3547)

3501 TCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAACAAACTCCCATTCAGCTCAATGGGGTGGAGACTTGAAATC

SnaBI (3675)

3600 CCCGTGAGTCAAACCGCTATCCACGCCATTGATGTAAGTCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTAAGTCCAAAGTAGG

NdeI (3780)

3700 AAAGTCCATAAGGTCATGTAAGTCCATAATGCCAGGCGGGCCATTTACCGTATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGT

3800 ACTGCCAAGTGGGAGTTTACCATAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCATTATTGACGTCA

SdaI (3958) PacI (3966) BspLU11I (3976)

3900 ATGGGCGGGGTCGTTGGGCGGTGAGCCAGCGGGCCATTTACCGTAAGTTATGTAACGCCCTG C A G G T T A A T T A A G A A C A T G T G A G C A A A A G G C C A G C

3998 AAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTG

4098 GCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGCTTACCGGATACCTGTCC

ApaLI (4290)

4198 GCCTTTCTCCCTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCCGTTGTAAGTGCCTCGCTCAAGCTGGGCTGTGTGCAGC

4298 AACCCCCGTTGAGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAAGTCCAAACCGTAAGACAGACTTATCGCCACTGGCAGCAGCCACTGG

4398 TAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGC

4498 GCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGC

4598 AGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGT

EagI (4726)
PacI (4706) SwaI (4715) NotI (4725)

4698 CATGGCTAGTTAATTAACATTTAAATC AGCGGCCGCAATAAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACAT

4798 ACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCCGAGTGCAAGTGCCAGGTGCCAGAACATTTCTCTATCGAA