



PvuI (7)
SgfI (6) 1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGCGCAATTGAACGGGTGCCTA

101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTATGTCGTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) 201 GTGAACGTTCTTTTTTCGCAACGGGTTTCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTTCACGCGCCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGCTTTTGTTCGTTT

NgoMIV (441)

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTCATCATGAGCTCCAAACAGCACTGTGTCAAATAATGATGG

601 CCACTTAATTCCTGCCCTGGGCTTTGGCACCTATAAACCCAAGGAGGTTCCCAAGAGTAAGTCACTGGAGGCTGCATGCCTAGCTCTAGATGTTGGGTAC

701 CGCCATGTTGATACTGCTTATGCATACCAAGTAGAAGAGGAGATAGGACAGGCCATTCAAAGCAAGATTAAGCTGGGGTTGTAAGAGAGAAAGACCTGT

801 TCCTCACTACAAAGCTTTGGTGCACTTGTCTTCGACCAGAGCTGGTCAAGCCTGCCTTGAAAAAGTCACTGAAAAAGCTTCACTGGAATATGTTGATCT

901 TTACATTATGCATTACCCAGTGCCAATGAAAGTCAAGGGATAATGATTTCCAGTAAATGAGCAAGGGAAATCTCTGTTGGACACTGTGGATTCTGTGAC

1001 ACATGGGAGAGGTTGGAGGAGTAAAGGATGCAAGGATGGTCAAGTCCATTGGGGTGTCCAACCTTTAACACAGGCAGCTGGAGCGAATCCTCAATAAGC

1101 CAGGACTGAAGTACAAACCTGTCTGCAACCAGGTTGAATGTCTCTATTTGAACCAGCGTAAGCTACTGGATTAAGTGCATCAAAAGACATTGTTCT

1201 IGTTGCTTACGGTGTCTGCGGACCCAGCGATATAAAGAATGGGTGGACCAGAACTCCCAGTCTCTTGAATGATCCAGTCTTTGTGATGTGGCCAAA

1301 AAGAACAAGCGAAGTCCCTGCTTGTGACCTTCGATACCTGATTCAACGTGGGATTGTGCCCTGGCCAGAGTTTCAAAGAGAATGAGATGAGAGAGA

1401 ATTTGCAGTTTTTTGGATTTGAGTTGTCCTGAGGACATGAAAACACTAGATGGCCTGAACAAAACCTTTCGATACCTCCAGCAGAGTTCCTTGTGA

1501 CCACCCAGAGTATCCATTTGTGGAGGAATATTAACATGGGACCTAATCATGGCTTGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACA

1601 AACCACACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAAC

1701 AACACAATTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATCTAAA

1801 ATACAGCATAGCAAACTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTG

1901 CATTAGCTGTTTGACGCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTCCCAAGGTTTGAAGTACTCTTCTTCTTTATGTTTAAATGC

2001 ACTGACCTCCACATTCCCTTTTGTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATG

2101 CTAAGGCCCTTCATAATATCCCCAGTTTGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTTGACAGCAAGAAAGCGAGCTTCTAGCTT

2201 TAGTTCCTGGTGTACTTGAGGGGATGAGTTTCTCAATGGTGGTTTTGACCAGCTTGCATTATCTCAATGAGCACAAGCAGTCAAGGATAGTCAG

2301 AGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTC

2401 CTTCTGCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAAGCAGACAGTACCCTGCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCC

2501 CCAGTCTTGGTCCTGATGGCCGCCCGACATGGTGCTTGTTCCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCT
40 G T K T R I A A G V H H K N D E Y L M T I K E T A V E V L E L D Q Q

BbsI (2614)

XmnI (2610)

AseI (2676)

2601 GAGAGATGTTGAAGGTCTTCATGGTGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGAT
7 S I N F T K M

2701 GGCGTCTCCAGCTTATCTGACGGTTCACATAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTT

SpeI (2831)

2801 GTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGC

SnaBI (2959)

2901 TATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAAGTCA

NdeI (3064)

3001 TGTAAGTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGT

3101 TTACCGTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCATTATTGACGTCAATGGGGCGGGGTCTGTTG

PacI (3250)

PstI (3243)

SdaI (3242)

BspLU11I (3260)

3201 GGCGTCCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAA

3301 AAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTA

3401 TAAAGATACCAGGCTTTCCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAA

ApaLI (3574)

3501 GCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCCGA

3601 CCGCTGCCCTTATCCGGTAACTATCGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGTGAAGCCAGTTA

3701 AGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGTGAAGCCAGTTA

3801 CCTTCGGAAAAAGAGTTGGTAGCTCTTATCCGGCAAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAA

PacI (3990) SwaI (3999)

3901 AGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAACTCACGTTAAGGGATTTTGGTCAATGGCTAGTTAATTAACA

EagI (4010)

NotI (4009)

4001 TTTAAATCAGCGCGCAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAA

4101 AACGAAACAAACAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA