



PvuI (7)
SgfI (6) 1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
 MfeI (82) **EcoNI (96)**

101 GAGAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
HindIII (245)
Bsu36I (291)

201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACACGCGCCCGCCCTACCTGAGGCC
 EcoNI (287)

301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCTCCTGAAGTGCCTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTAACTCTACGTCTTTGTTTCGTTT

BspHI (560)

501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGGCTACCTGAGATCACCGGTCATCATGAGCCTCCCGTCCAGCCGCGGGCCCGTGTCCCGGG
 XmaI (594)

AgeI (552)
M S L P S S R A A R V P G

601 TCCTTCGGGCTCCTTGTGCGCGCTGCTCGCGTGTCTGCTCCTGCTGACGCCGCGGGCCCTCGCCAGCGCTGGTCCCTGTCTGTGTGTGACAGAG
 Bsp120I (655)
Eco47III (667)

701 CTGCGTTGCACTTGTACGCGTTACGCTGAGAGTAAACCCAAAACGATTGGTAAACTGCAGGTGTTCCCGCAGGCCCAGTGTCTCAAGGTGGAAG
 MluI (717)
PstI (757)
XcmI (788)

801 TGGTAGCTCCCTGAAGAACGGGAAGCAAGTTTGTCTGGACCCGGAAGCCCTTTTCTAAAGAAAGTCAATCCAGAAAATTTTGGACAGTGGAAACAAGAA
 MscI (913)

901 AAAGTCTAGCTAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGACAAAACCACAACCTAGAAATGCAGTGAATAAATGCTTTATTTGTAAATTTGT
 NheI (907)

113▶ N •

HpaI (1045)
MfeI (1056)

1001 GATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCATTTTATGTTTCAGGTTCAAGGGGAGGTGTGGG

EcoRI (1141)

1101 AGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACCTTAACTCCAATCAAGCCTCTACTTGAATC

1201 CTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTTTTCATGGAGTTAAGATATAGT

SapI (1323)
SspI (1380)
SwaI (1394)

1301 GTATTTTCCAAGGTTTGAAGTCTCTTCATTCTTTATGTTTTAAATGCAGTACCTCCACATTCCTTTTATGTAATAATTCAGAAATAATTTAA

1401 ATACATCATTGCAATGAAAATAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAA

1501 CAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGA
 SacI (1655)
BstXI (1684)

1601 CCAGCTTCCATTATCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCT
 SacI (1655)
BstXI (1684)

1701 GTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCAGCACAG
 StuI (1819)

1801 ACAGTGACCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCGAGTCTTGGTCTGATGGCGCCCGGACATGGTGTGTTGCTCCTCAT
 BbsI (1965)
XmnI (1961)

1901 AGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCTGTTGGCCCTCTATAGTGAGTCGTATT
 AseI (2027)
SacI (2084)

2001 ATACTATCGCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCCTAAACGAGCTCTGCTTATATA

AseI (2027)
SpeI (2182)

2101 GACCTCCCACCGTACACGCCCTACGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAAACTC

2201 CCATTGACGTCATGGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCCATTTGATGACTGCCAAAACCGCATCATCATGGTAATAG

SnaBI (2310)

2301 CGATGACTAATACGTAGATGACTGCCAAGTAGAAAGTCCCATAAAGTCAAGTCAAGTCAAGTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAAT

NdeI (2415)

2401 AGGGGGCGTACTTGGCATATGATACACTTGTACTGCAAGTGGGCGAGTTTACCGTAAATACTCCACCATTGACGTCAATGGAAAGTCCCTATTGGC

PstI (2594)
SdaI (2593)

2501 GTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAG

PacI (2601) BspLU11I (2611)

2601 GTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAG

2701 CATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCTTTCCCTGGAAGCTCCCTCGTGCGCTCTCCTG

2801 TTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTA

ApaLI (2925)

2901 GGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTCCAACCCGGTAAGA

3001 CACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACG

3101 GCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCCACCGC

3201 TGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAG

EagI (3361)

PacI (3341) SmaI (3350) NotI (3360)

3301 TGGAACGAAAACACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGCCGCAATAAAATATCTTTATTTTCATTACATCTGTG

3401 TGTGTTTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCAAACAACGAAACAACAACTAGCAAATAGGCTGTCCCCAGTGCAAGTGCA

3501 GGTGCCAGAACATTTCTCTATCGAA