



PvuI (7)
SgfI (6)
1 GGATCTGCATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATGCCACAGTCCCCGAGAAGTTGGGGGGAGGGGTCGGCAATTGAACCGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203)
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCCTCCCGCCTGTGGTGCTCTGAACTGCGTCCGCGCTTAGGTAAGTTTAAAGCTCAGGTCGAGACC
401 GGGCCTTTGTCCGGCGCTCCCTTGAGGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTCAACTCTACGTCTTTGTTTCGTTT
501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGGCGCCTACCTGAGATCACCGTAGGAGGCCATCATGAGACAGACTTTGCCTGTATCTACTTT

DraIII (644)
601 TGGGGGGCCTTTTGCCTTTGGGATGCTGTGTGCATCCTCCACCACCAAGTGCCTGTTAGCCATGAAGTTGCTGACTGCAGCCACCTGAAGTTGACTC
701 AGGTACCCGATGATCTACCCACAAACATAACAGTGTGAACCTTACCCATAATCAACTCAGAAGATTACCAGCCGCAACTTCACAAGGTATAGCCAGCT
801 AACTAGCTTGATGTAGGATTTAACACCATCTCAAACCTGGAGCCAGAATTGTCCAGAACTTCCCATGTTAAAAGTTTTGAACCTCCAGCACAAATGAG
901 CTATCTCAACTTTCTGATAAACCTTTGCCTTCTGCACGAATTTGACTGAACTCCATCTCATGTCCAACCTCAATCCAGAAAATAAAAATAATCCCTTTG

BglII (1025)
1001 TCAAGCAGAAGAATTTAATCACATTAGATCTGTCTCATAATGGCTTGTCATCTACAAAATTAGGAACTCAGGTTGAGTTGAAAAATCTCCAAGAGCTTCT

Eco47III (1121)
1101 ATTATCAAACAATAAAATCAAGCGCTAAAAAGTGAAGAACTGGATATCTTTGCCAATTCATCTTTAAAAAATTAGAGTTGTCATCGAATCAAATAAA

SandI (1268)
1201 GAGTTTTCTCCAGGGTGTTCACGCAATTGGAAGATTATTTGGCCTTTTCTGAACAATGTCCAGCTGGGTCCAGCCTTACAGAGAAGCTATGTTTGG
1301 AATTAGCAAAACACAGCATTGGAATCTGTCTCTGAGTAACAGCCAGCTGTCCACCACCAGCAATACAACCTTTCTTGGGACTAAAGTGACAAATCTCAC
1401 TATGCTCGATCTTCTCACAACAACCTAAATGTGGTTGGTAACGATTCTTTGCTTGGCTTCCACAAC TAGAATATTTCTTCTAGAGTATAATAATATA
1501 CAGCATTGTTTTCTCACTTTTGACGGGCTTTTCAATGTGAGGTACCTGAATTTGAAACGGTCTTTACTAAACAAAGTATTTCCCTTGCCTCACTCC
1601 CCAAGATTGATGATTTTTCTTTTCAGTGCTAAAATGTTTGGAGCACCTAACATGGAAGATAATGATATTCCAGGCATAAAAAGCAATATGTTACAGG
1701 ATTGATAAACCTGAAATACTTAAGTCTATCCAACCTCTTTACAAGTTTGCGAACCTTTGACAAATGAAACATTTGTATCACTTGCTCATTCTCCCTTACAC

ScaI (1871)
1801 ATACTCAACCTAACCAAGAATAAAATCTCAAAAATAGAGAGTGATGCTTTCTTTGGTTGGGCCACCTAGAAGTACTTGACCTGGGCCTTAATGAAATTG

XbaI (1931)
1901 GGCAAGAACTCACAGGCCAGGAATGGAGAGGTCTAGAAAATATTTTTGAAATCTATCTTCTCACAACAAGTACCTGCAGCTGACTAGGAACTCCTTTGC
2001 CTTGGTCCCAAGCCTTCAACGACTGATGCTCCGAAGGGTGGCCCTAAAAATGTGGATAGCTCTCCTTACCATTCCAGCCTCTTCGTAACCTGACCATT
2101 CTGGATCTAAGCAACAACAACATAGCCAACATAAATGATGACATGTTGGAGGGTCTTGAGAACTAGAAAATCTCGATTTGCAGCATAACAACCTTAGCAC
2201 GGCTCTGGAACACGCAAACCTGGTGGTCCCATTATTCTTCTAAAGGGTCTGTCTCACCTCCACATCCTTAACCTGGAGTCCAACGGCTTTGACGAGAT

ClaI (2340)
2301 CCCAGTTGAGGTCTTCAAGGATTTATTTGAACTAAAGATCATCGATTTAGGATTGAATAATTTAAACACACTTCCAGCATCTGTCTTTAATAATCAGGTG
2401 TCTCTAAAGTCATTGAACCTT CAGAAGAATCTCATAACATCCGTTGAGAAGAAGGTTTTCCGGCCAGCTTT CAGGAACCTGACTGAGTTAGATATGCGCT
2501 TTAATCCCTTTGATTGCACGTGTGAAAGTATTGCCTGGTTTGTAAATTGGATTAACGAGACCCATACCAACATCCCTGAGCTGTCAAGCCACTACCTTTG

BclI (2689)
2601 CAACACTCCACCTCACTATCATGGGTTCCAGTGAGACTTTTTGATACATCATCTTGCAAAGACAGTGCCCCCTTTGAACTCTTTTTCATGATCAATACC
2701 AGTATCCTGTTGATTTTTATCTTTATTGACTTCTCATCCACTTTGAGGGCTGGAGGATATCTTTTTATTGGAATGTTTCAGTACATCGAGTTCTTGGTT
2801 TCAAAGAAATAGACAGACAGACAGAACAGTTTGAATATGCAGCATATATAATTCATGCCTATAAAGATAAGGATTGGGTCTGGGAACATTTCTCTTCAAT
2901 GGAAAAGGAAGACCAATCTCTCAAATTTGTCTGGAAGAAAGGACTTTGAGGCGGGTGTTTTTGAACTAGAAGCAATTGTTAACAGCATCAAAAAGAC
3001 AGAAAAATATTTTTGTTATAACACACCATCTATTAAGACCCATTATGCAAAGATTCAAGGTACATCATGCACTTCAACAAGCTATTGAACAAAATC

3101 TGGATTCCATTATATTGGTTTTCTTGAGGAGATTCCAGATTATAAACTGAACCATGCACTCTGTTTGCAGAGGAATGTTTAAATCTCACTGCATCTT

3201 GAACTGGCCAGTTTCAGAAAGAACGGATAGGTGCCTTTCTGCATAAATTGCAAGTAGCACTTGGATCCTAAAACTCTGTGCATCTGTACAAGGgtaagtca

BamHI (3261) **BsrGI (3283)**

3301 ctgactgtctatgctctgggaagggtggcaggagatggggcagtgaggaaaagtggcactatgaaccACTAGTTTGACAATTAATCATAAGCATAGT

3401 ATAATACTCACTACTATAGcaattgtactaaccttcttctcttctcctgacag**GAGGAG**CCATCATGGCCGCTATGGAGATCGAGTGCCGCATCA

3500 CCGGCACCCTGAACGGCTGGAGTTCGAGCTGGTGGCGGCGGAGAGGGCACCCCGAGCAGGGCCGCATGACCAACAAGATGAAGAGCACCAAAGGCGC

3600 CCTGACCTTCAGCCCTACCTGCTGAGCCACGTGATGGCTACGGCTTCTACCACTTCGGCACCTACCCAGCGGCTACGAGAACCCCTTCTGCACGCC

3700 ATCAACAACGGCGGTACACCAACACCCGCATCGAGAAGTACGAGGACGGCGGTGCTGCACGTGAGCTTACGCTACCGCTACGAGCGCGCGGTGA

3800 TCGGCGACTTCAAGGTGATGGGCACCGGCTTCCCGAGGACAGCGTGATTTCCCGACAAGATCATCCGAGCAACGCCACCGTGGAGCACCTGCACCC

BsaBI (3906)

3900 TATGGGCGATAACGATCTGGATGGCAGCTTACCCGCACCTTACGCTGCGCGACGGCGCTACTACAGCTCCGTGGTGGACAGCCACATGCACTTCAAG

Bsp120I (4031)

4000 AGCGCCATCCACCCAGCATCTGCAGAACGGGGGCCCATGTTCCGCTTCCGCCGTGGAGGAGGATCACAGCAACACCGAGCTGGGCATCGTGGAGT

NheI (4138)

4100 ACCAGCACGCTTCAAGACCCCGGATGCAGATGCCTAAAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAGTAAATGC

HpaI (4276)

4200 AGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAAACAATTGCATTCA

4300 TTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAAAC

4400 TTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGC

4500 CTCACCTTCTTCATGGAGTTTAAAGATATAGTGTATTTCCCAAGTTTGAAGTACTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTC

4600 CCTTTTTAGTAAATATTAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTTATTAGGAGAATCCAGATGCTCAAGGCCCTTCATAA

4700 TATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTCTGGTGTACTT

141 ◆ ● ● ● AsnArgThr TyrLys

4800 GAGGGGATGAGTTCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCAC

135 ◆ LeuP roI l eLeuGl uGl u l eThr Thr LysVal l eLysGl yAsnMe tGl u l eLeuVal l PheCysAspP roAl aTyrAspSer l l eLeuGl uArgCysM

4900 ATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAAATGGTGTCAAAGTCTTCTGCCGTTGCTCA

101 ◆ e tGl yCysP roSer Val l eAl a l eAl aGl uAl aCysVal l Thr Val a rGl y l eTyrAl aGl u l eHi sValAl aSer l l e l eGl uGl yThr LysThr Arg l e

5000 CAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGTCTGTAT

68 ◆ l Al aSer Gl y l eAl a l eAl aGl uAl aCysVal l Thr Val a rGl y l eTyrAl aGl u l eHi sValAl aSer l l e l eGl uGl yThr LysThr Arg l e

5100 GGCGCCCGACATGGTGTCTTGTCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTC

35 ◆ Al aAl aGl yVal l Hi sHi sLysAsnAspGl uTyrLeuMe tThr l l eLysGl uThrAl aVal l Gl uVal l LeuGl uLeuAspGl nGl nSer l l eAsnPheThr L

5200 TTCATGATGGCCCTCTATAGTGTGATTACTATGCGATATACTATGCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATC

1 ◆ ysMe t

5300 TGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTGCGTCAATGGGGCGGAGTTGTACGACATTTGGAA

5400 AGTCCCGTTGATTTACTAGTCAAAAACAACTCCCAATTGACGTCAATGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGAT

SnaBI (5541)

5500 GTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTGATGACTGGGCATAATGC

NdeI (5646)

5600 CAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGACTGCCAAGTGGGAGTTTACCGTAAATACTCCA

5700 CCCATTGACGTCAATGAAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGTCGTTGGCGGTCAGCCAGGCGG

SdaI (5824)

5800 GCCATTTACCGTAAGTTATGTAACGCCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGCCGCTTGTGG

5900 CGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGCGTT

6000 TCCCCTGGAAGCTCCCTCGTGCCTCTCTGTCCGACCCTGCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCAT

ApaLI (6156)

6100 AGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTACGCCGACCGCTGCGCCTTATCCG

6200 GTAACATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGATGTAGGCGGTGC

6300 TACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTT

6400 GGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGGTTTTTTTGTGGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATC

NotI (6591)

6500 CTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGAAAACACAGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGC

6600 AATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAAC

6700 TAGCAAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTATCGAA