



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

101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

**HindIII (245)**  
**Psp1406I (203)** 201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCAGAGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC Bsu36I (291)

301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCTGTGGTGCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

**NgoMIV (441)**  
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

**BstEII (555)**  
**KasI (535)** 501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGGCTACCTGAGATCACCGGTACCATTGGCCCCAGCCTTCTGCTGCTGCTGCTGCTGCTGGCC AgeI (552)

601 ACAGGGTTGCGTCTCAGGCCCTCTGCTGG GAGCGCATAACAAAAGTGAGGCTCCTTGAAGGGGAGACTCTGCTGTGCGAGTGTCTCTATAAGGGCTAC Tth111I (666)

701 AAAAAACCGTGGAGGGCAAGTTTGGTGCAAAATCAGGAAGAAGAAGTGAGCCTGGCTTTGCCGAGTCTGGGTGAAAGGGCCCCGCTACTTGTCTC Bsp120I (781) PstI (796)

801 AGGACGATGCCAGGCCAAGGTGGTCAACATCACCATGGTGGCCCTCAAGCTCCAGGACTCAGGCCGATACTGGTGCATGCGCAACACCTCTGGGATCCT FspI (878) BamHI (893)

901 GTACCCCTTGATGGGCTTCCAGCTGGATGTGTCTCCAGCTCCCAAACTGAGAGGAACATTCTTTACACATCTGGACAACATCCTCAAGAGTGAAGT PvuII (919)

1001 GTCACAACCTGGCCAAGCCCTACCTCAGGCCCTGATGCCCTTTTACCCTGGTGTGATGGTGTTCACCCAGGACTCATCACCTTGCTAGGCTCTTAG Bsu36I (1022) XcmI (1046) AvrII (1087)

1101 CCTCCACCAGACCTGCCTCAAGACAGGCTACAGCTTCACTGCTACCAGCACCACCAGGCCAGGACCAGGACCATTGGGTCGCCAGACAGTACCAGC SandI (1161) SandI (1180) Tth111I (1194)

1201 GTCTCCAGCAATGCCAGGACTCCTCTGCTGGCCGAGAACATCCATCTCCACTAAGTCTGGGACTCAGCACCAGATCGCCACCACAGGGCTCTGCCTC SphI (875)

1301 ACCAGCAGATCTCTCAACAGACTACCCTCCATGCCCTCCATCAGGCACCAGGATGTTTACTCCACTGTGCTTGGGTTGGTGTGACCTCCTGGTGC BglII (1306)

1401 TGATGCTGATCATGGTCTATGGGTTTTGGAAGAAGAGACACATGGCAAGCTACAGCATGTGCAGCGATCCTTCTACACGTGACCCACCTGGAAGACCAGA BsaBI (1401) BbrPI (1475) BbsI (1490)

1501 GCCCTATGTGAAGTCTACTTGTCTGAGGCCACTTAAGCATGGGGTCTAGTCTGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAAC NheI (1547)

1601 TAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAATA HpaI (1685) MfeI (1696)

1701 TGCATTCATTTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATCTAAATACAGCAT EcoRI (1781)

1801 AGCAAACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTG

1901 TTTGAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGTTTGAAGTGTCTTCTTCTTATGTTTTAAATGCACTGACCTC SapI (1963)

2001 CCACATTCCTTTTATGATAAATATTGAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCC SspI (2020) SmaI (2034)

2101 CTTCAATAATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTCTAGCTTTAGTTTCCTG SacI (2295)

2201 GTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCT BstXI (2324)

2301 CTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCC 104 R C M G C P S V V R I S R D V E D S Y P H R V A V I T D F D K Q G

