



125

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

101 GAGAAGTGGCGCGGGTAAACTGGAAAGTGATGCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

Psp1406I (203) 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC EcoNI (287)

301 GCCATCCACGCGGTTGAGTCGCGTTTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCGGCTCTCCACGCTTGGCTGACCCTGCTTGTCTAACTCTACGCTTTGTTTCGTTT NgoMIV (441)

501 TCTGTTCTGCGCGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTACAGTGTGCCCTGCTGCTGCTGCCCTGCTGTGGGGGG SphI (560) SandI (598)

601 GTCCTGCAGGAGAAGCCAGTGTACGAGCTGCAAGTGCAGAAGTGGTACGGTGCAGGAGGGCTGTGCGTCTTGTGCCCTGCTCCTTCTCTTACCCC SdaI (603)

701 TGGAGATCCTGGTATTCCTCTCCCCACTCTACGTCTACTGGTCCGGGACGGGAGATCCATACTACGCTGAGGTTGTGCCACAACAACCCAGACA

801 GAAGAGTGAAGCCAGAGACCCAGGGCGATTCCGCCTCTTGGGGATGTCAGAAGAAGAACTGCTCCTGAGCATCGGAGATGCCAGAATGGAGGACAC

901 GGAAGCTATTTCTCCGCTGGAGAGAGGAAGGATGTAATAATAGCTACCAACAGAATAAGCTGAAGTGGAGGTGACAGCCTGATAGAAACCC

1001 GACATCCACTTTCTGGAGCCTCTGGAGTCCGGCCGCCCAAGGCTGAGCTGCAGCCTTCCAGGATCCTGTGAAGCGGGACCACCTCTCACATTCTCCT BamHI (1063)

1101 GGACGGGAATGCCTCAGCCCCCTGGACCCGAGACCACCGCTCCTCGGAGCTCACCTCACCCCAAGGCCGAGGACCATGGCACCAACCTCACCTG BsrBI (1140) NcoI (1179)

1201 TCAGATGAAACGCAAGGAGCTCAGGTGACCACGGAGAGAAGTGTCCAGCTCAATGTCTCCTATGCTCCACAGACCATCACCATCTTCAGGAACGGCATA BstEII (1224)

1301 GCCCTAGAGATCTGCAAAACACCTCATACCTTCCGGTCTGGAGGGCAGGCTCTGCGGCTGCTGTGATGCTCCAGCAACCCCTGCACACCTGA

1401 GCTGGTTCCAGGGTCCCCTGCCCTGAACGCCACCCCATCTCAATACCGGATCTTGAGCTTCGTCGAGTAAGGTGTCAGAAGAAGGAGGCTTCCAC

1501 CTGCCGCGCTCAGCACCCGCTGGGCTTCTGCAAAATTTTCTGAATCTCTCAGTTTACTCCCTCCCACAGTTGCTGGGCCCTCTGCTCCTGGGAGGCT Bsp120I (1575)

1601 GAGGGTCTGCACTGCAGATGCTCCTTTTCGAGCCCGCCGCCCCCTCCTGTGCTGGCGGCTTGGAGAGAAGCCGCTGGAGGGGAACAGCAGCCAGGGCT NgoMIV (1635)

1701 CATTCAAGGTCAACTCCAGCTCAGTGGGCCCTGGCCAAACAGCTCCCTGATCCTCCAGGGGGGCTCAGCTCCGACCTCAAAGTCAAGTGCAGGCTGCAAGGCCTG Bsp120I (1726) StuI (1793)

1801 GAACATCTATGGTCCCAGAGCGGCTGTGCTGCTGCAAGGGAGATCGAACCTCGGGACAGGAGTGGTCTGCAGCCCTTGGTGGTGTGGTGTG BsrBI (1818) SandI (1810)

1901 ATGGCCCTGCTCTGTATCTGTCTGTGCCTCATCTTTTAAATAGTAAAGCCCGCAGGAAGCAAGCAGCTGGGAGACCAGAGAAAATGGATGATGAAG BbsI (1996)

2001 ACCCCATTATGGGTACCATCACCTCGGTTCCAGGAAGAAGCCCTGGCCAGACAGCGCCGGAGATCAAGCATCTCCTCTGGGGATGCCCTCCCTTGGGA Acc65I (2011)

2101 AGAACAAAAGGAGCTCCATTATGCCTCCCTTAGTTTTTCTGAGATGAAGTCGAGGGAGCTAAGGACCAAGGCCCCAAGCACCACGGAGTACTCGGAG Bsu36I (2158) ScaI (2189)

2201 ATCAAGACAAGCAAGTGAGGATTTGCCAGAGTTCAGTCTGCTAGTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCAACTAGAAT NheI (2241)

2301 GCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAATTGCATT HpaI (2379) MfeI (2390)

2401 CATTTTATGTTTCAGGTTACAGGGGAGGTGTTGGAGGTTTTTAAAGCAAGTAAACCTCTACAATGTGGTATGGAATCTAAAATACAGCATAGCAA EcoRI (2475)

2501 ACTTTAACCTCCAATCAAGCCTCTACTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGCTGTTGCCAATGTGCATTAGCTGTTTGCA

2601 GCCTCACCTTCTTTCATGGAGTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACAT

2701 TCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCAT

2801 AATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGTAC

2901 TTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGC
 135 K L P I L E E I T T K V L K G N M E I L V F C D P A Y D S I L E R C

3001 ACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCCTTCTGCCGTTGCT
 102 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 N S

3101 CACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCCTG
 69 V A S G I A I A E A C V T V R G I Y A E I H V A S I I E G T K T R

3201 ATGCGCGCCCGACATGGTGCTTGTTCCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGG
 35 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 S I N F T

3301 TCTTCATGGTGGCCCTCCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTA
 24 K M

3401 TCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCTACCGCCATTGCGTCAATGGGCGGAGTTGTTACGACATTTTGG

3501 AAAGTCCCGTTGATTTACTAGTCAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCCATTTG

3601 ATGTAAGTCCAAACCGCATCATCATGGTAATAGCGGATGACTAATACGTAGATGACTGCAAGTAGGAAAGTCCATAAGGTCATGACTGGGCATAAT

3701 GCCAGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGACTGCAAGTGGGCAGTTTACCGTAAATACTC

3801 CACCCATTGACGTCAATGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTACGCCAGGC

3901 GGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAAGACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGTTGCT

4001 GGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCG

4101 TTTCCCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTCTCCCTTGGGAAGCGTGGCGCTTTCTC

4201 ATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCCTTATC

4301 CGGTAACATATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGT

4401 GCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAG

4501 TTGGTAGCTCTTGATCCGGCAAACAACCACCGCTGGTAGCGGTGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGA

4601 TCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGC

4700 CGCAATAAATATCTTTATTTTTCATTACATCTGTGTGGTTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACA

4800 AACTAGCAAATAGGCTGTCCCCAGTGCAAGTGACAGTGCCAGAACATTTCTCTATCGAA