



PvuI (7)
SgfI (6) 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA **MfeI (82)**
101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGTCCGCCTTTTTCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC

Psp1406I (203) 201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACGCGCCCGCCGCTACCTGAGGCC **HindIII (245)** **Bsu36I (291)**
301 GCCATCCACGCCGGTTGAGTCGCGTTCTGCGCCCTCCCGCTGTGGTGCCTCCTGAAGTGCCTCCGCTCTAGGTAAGTTAAAGCTCAGGTCGAGACC

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGCTTTGTTTCGTTT **NgoMIV (441)**

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCGCCTACCTGAGATCACCGGTATCATGAGTGGACCTTTCCAGCAGAGCCTACAAAAGGTCC **AgeI (552)**
1 M S G P F P A E P T K G P **BspHI (560)**
BspHI (674)

601 CCTCGCATGCAACCTGCTCCAAAAGTGAACCTCAAAGGACATCTTACTGGTGGCCCCACAAAAGCTTCTCATGAGGGAATCAAAGGCTTTGGG **Bsp120I (654)** **HindIII (666)**
13 L A M Q P A P K V N L K R T S S L V G P T Q S F F M R E S K A L G
BspHI (710) 701 GCTGTCAAATCATGAATGGCCTCTTCCATATTACCCTGGGGGACTGCTGATGATCCACAGGGGTCTTCGCACCCATCTGTTTGTGATGATGGTACC **Acc65I (794)**
47 A V Q I M N G L F H I T L G G L M I P T G V F A P I C L S V W Y

801 CTCTCTGGGAGGCATTATGTACATTATTTCCAGGATCACTCCTGGCAGCTGCAGCAGAAAAAACCCTCCAGGAAGAGTTTGGTCAAAGCAAAGTGATAAT **BsrGI (818)** **PstI (848)**
80 P L W G G I M Y I I S G S L L A A A A E K T S R K S L V K A K V I M
901 GAGCTCTAAGCCTCTTTGCTGCCATTTCTGGAATAATTTCTTCAATCATGGACATACTTAACATGACACTTCTCATTTTTTAAAAATGAGAAGACTG **ScaI (1085)**
113 S S L S L F A A I S G I I L S I M D I L N M T L S H F L K M R R L

1001 GAGCTTATCAAACCTTCAAAGCCGATGTTGATATCTACGACTGTGAACCATTAATTCCTCAGAGAAAACTCCCATCTACACAGTACTGTAAACAGCA **EcoRV (1030)** **ScaI (1085)**
147 E L I Q T S K P Y V D I Y D C E P S N S S E K N S P S T Q Y C N S

1101 TTCAGTCTGTGTTCTTGGCATTCTGTGCGCGATGCTGATCTCTGCTTCTTCCAGAACTTGTGACAGCTGGTATTGTGAGAAATGAGTGGAAAAGAAT **BsaBI (1131)**
180 I Q S V F L G I L S A M L I S A F F Q K L V T A G I V E N E W K R M

1201 GTGTACCAGATCCAAATCTAATGTGGTTCTGCTGTGAGCTGGAGAAAAAATGAGCAGACGATTAAAATGAAAGAAGAAATCATTGAGCTAAGTGGAGTA **BsaBI (1208)**
213 C T R S K S N V V L L S A G E K N E Q T I K M K E E I I E L S G V

1301 TCTTCCCAACCAAGAATGAAGAGGAAATTGAAATTATCCAGTGCAGGAGGAAGAAGAAGAAGCAGAAATAAATTTCCAGCACCTCCCAAGAGC **XmnI (1330)**
247 S S Q P K N E E E I E I I P V Q E E E E E A E I N F P A P P Q E

1401 AGGAATCCTTGCCAGTGGAAAATGAGATCGTCTTAACTCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAA **NheI (1442)** **MscI (1448)**
280 Q E S L P V E N E I A P •

1501 TGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACCAAGTTAACAAACAATTGCAT **HpaI (1580)** **MfeI (1591)**

1601 TCATTTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAA **EcoRI (1676)**
1701 AACTTTAACTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTGCCAATGTGCATTAGCTGTTTGC

1801 AGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTCTTCTATTCTTTTAAATGCACTGACCTCCACA **SapI (1858)**

1901 TTCCCTTTTATGATAAATATTGAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCA **SspI (1915)** **SwaI (1929)**
2001 TAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTA **XmnI (2496)**
141 • N R T Y

2101 CTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTG **BstXI (2219)**
136 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

2201 CACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGC **StuI (2354)**
102 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 N S

2301 TCACAGCAGACCAATGGCAATGGCTTACAGCAGACAGTACCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGTCTC **StuI (2354)**
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

2401 GATGGCCGCCGACATGGTCTTGTGCTCCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAAG **XmnI (2496)**
36 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

2501 GTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTT
2601 ATCTGACGGTTCACATAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTG

2701 GAAAGTCCCGTTGATTTACTAGTCAAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATT

2801 GATGTAAGTCCAAACCGCATCATCATGGTAATAGCGTAGTAATAACGTAGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAA

2901 TGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACT

3001 CCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGGTCTGTTGGGCGGTGAGCCAGG

3101 CGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGC
3201 TGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGC
3301 GTTCCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCT

3401 CATAGCTACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCCTGCGCCTTAT
3501 CCGGTAACATATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGG
3601 TGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGA
3701 GTTGGTAGCTCTTGATCCGGCAAACAACACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAG

3801 ATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAAACGAAAACCTACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGGC

3901 CGCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACA
4001 AACTAGCAAATAGGCTGTCCCCAGTGAAGTGCAGGTGCCAGAACATTTCTCTATCGAA

AseI (2562)
SpeI (2717)
SnaBI (2845)
NdeI (2950)
PacI (3136)
PstI (3129)
SdaI (3128)
BspLU11I (3146)
ApaLI (3460)
PacI (3876) SwaI (3885) EagI (3896)
NotI (3895)