



PvuI (7)
SgfI (6) **EcoNI (96)**
1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
101 GAGAAGGTGGCGCGGGTAAACTGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245) **Bsu36I (291)**
Psp1406I (203) **PvuII (239)** **EcoNI (287)**
201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCCCTACCTGAGGGCC
301 GCCATCCACGCGGTTGAGTCGCGTTCGCCGCTCCCGCCTGTGGTGCCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
NaeI (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCGCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCAACTCTACGTCTTTGTTTCGTTT

KasI (535) **AgeI (552)** **BspHI (560)**
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGCGCTACCTGAGATCACCGGTCATCATGACTAGCCGGAAACCAAGTTTCACTGTGTAATTG
1 M T S R E H Q V S L C N C

KasI (614) **BsrBI (678)** **XhoI (683)**
601 CGTCCCCCTACTCCGGCGCTCCTTTGCGACGCTCCCTGGAGAAAAGCAGCCCACTGCACGCGCTCAGTCGACTTCCGCTCTCGAGTGTCTCCAAGC
13 V P L L R R L L C D A P W R K A R P L H A L S R Y F R S R V S P S
701 AAGATGGCGGAGGAGCCGAGTCTGTGTGCGAGTTCCTACTTCAATTGCTGCTGGAGGGAAAGGACTTACGGATGTCTCCCAAGAAACAACCCCGCCG
47 K M A A E E P Q S V L Q L P T S I A A G G E G L T D V S P E T T T P

BbsI (807) **PstI (815)** **XmaI (825)**
801 AGCCCCCGTCTTCCGCTGCAATTTCCCGGGAACAGAGAACCTGCTGGCGACACCAAGAAAAAATTGACATTTTGCTAAAGGCTGTGGGAGACACTCC
80 E P P S S A A V S P G T E E P A G D T K K K I D I L L K A V G D T P
901 TATTATGAAAACAAAGAAGTGGCGAGTAGAGCGAACACCAACCATCCAAGGACTCATTGACTTCAAAAAGTTTCTTAAACTTGTGCCTCAGAACAG
113 I M K T K K W A V E R T R T I Q G L I D F I K K F L K L V A S E Q

BstXI (1041)
1001 TTGTTTATTTATGTAATCAGTCCTTTGCTCCTCCCAAGAAAGTGGAACTCTATGAGTGTGTTGGCAGTGTGTAAGCTGTTTACATT
147 L F I Y V N Q S F A P S P D Q E V G T L Y E C F G S D G K L V L H

MscI (1134) **NheI (1128)**
1101 ACTGCAAGTCTCAGCGTGGGGATGAACCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAACAAAA
180 Y C K S Q A W G •

HpaI (1266)
1201 TGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAACAAACAACATTGCATTCTTTTATGTTTC

EcoRI (1362)
1301 AGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTTAAAATACAGCATAGCAAACTTTAACCTCCA
1401 AATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGAGCCTCACCTTCTT
1501 TCATGGAGTTAAGATATAGTGTATTTCCCAAGGTTTGAAGTCTCTCATTCTTTATGTTTAAATGCACTGACCTCCACATTCCCTTTTAGTA

SspI (1601) **SwaI (1615)** **EcoO109I (1676)**
1601 AAATATTCAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTCATAATATCCCCAGT
1701 TTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTACTTGAGGGGGATGA
141 • N R T Y K L P I L

SacI (1876)
1801 GTTCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAGCAGCTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGG
131 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 M G C P

BstXI (1905)
1901 GCTGACCACCCTGATGGTCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTGCTCACAGCAGACCCA
98 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 V A S G

StuI (2040)
2001 ATGGCAATGGCTTACGACAGACAGTACCCCTGCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGTCTGATGGCCGCCCGA
64 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 I A A G V

BspHI (2190) **BbsI (2186)** **XmnI (2182)**
2101 CATGGTGCTTGTGCTCATAGAGCATGGTGTCTTCTCAGTGGCGACTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGATGGC
31 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 K M

AseI (2248)
2201 CCTCCTATAGTGAAGTCGATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTTCA

SacI (2305)
2301 TAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTGCGTCAATGGGCGGAGTTGTTACGACATTTGGAAAGTCCCGTTGA

