



PvuI (7)
SgfI (6) MfeI (82) EcoNI (96)
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
101 GAGAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245) **Bsu36I (291)**
Psp1406I (203) **PvuII (239)** EcoNI (287)
201 GTGAACGTTCTTTTTCGCAACGGGTTTGGCCGAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTCGCGTTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

KasI (535) **AgeI (552)** **BspLU11I (560)** **BstAPI (570)**
501 TCTGTTTGGCGCGTTACAGATCCAAGCTGTGACCGCGCGCTACCTGAGATCACCGGTCAACATGTGGCTGCAGAGCCTGCTGCTTGGGCACTGTGGC
1▶ M W L Q S L L L L G T V A

BsrBI (621)
601 CTGCAGCATCTCTGCACCCGCGCTCGCCAGCCAGCAGCAGCCCTGGGAGCATGTGAATGCCATCCAGGAGGCCCGGCGTCTCTGAACCTGAGT
13▶ C S I S A P A R S P S P S T Q P W E H V N A I Q E A R R L L N L S

BsrGI (794)
701 AGAGACACTGCTGCTGAGATGAATGAAACAGTAGAAGTCATCTCAGAAATGTTTGACCTCCAGGAGCCGACCTGCCTACAGACCCGCTGGAGCTGTACA
47▶ R D T A A E M N E T V E V I S E M F D L Q E P T C L Q T R L E L Y

BstAPI (801) **Bsp120I (834)** **MscI (851)**
801 AGCAGGGCCTGCGGGGAGCCTACCAAGCTCAAGGGCCCTTGACCATGATGCCAGCCACTACAAGCAGCACTGCCCTCAACCCCGAAACTTCTGT
80▶ K Q G L R G S L T K L K G P L T M M A S H Y K Q H C P P T P E T S C

NheI (997)
901 TGCAACCCAGACTATCACCTTTGAAAGTTTCAAAGAGAACCTGAAGACTTTCTGCTTGTATCCCTTTGACTGCTGGGAGCCAGTCCAGGAGTGAAGCT
113▶ A T Q T I T F E S F K E N L K D F L L V I P F D C W E P V Q E •

MscI (1003)
1001 AGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTG

HpaI (1135) **MfeI (1146)**
1101 CTTTATTTGTAACCATTATAAGCTGCAATAAACAGTTAAACAACAACATTGCATTATTTTTATGTTTCAGGTTACAGGGGAGGTGTGGGAGGTTTTTTA

EcoRI (1231)
1201 AAGCAAGTAAACCTCTACAAATGTGGTATGGAATCTAAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAG
1301 GGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTCTTTTCATGGAGTTAAGATATAGTGATTTTTCC

SapI (1413) **SspI (1470)** **SwaI (1484)**
1401 AAGGTTTGAAGTACTGCTCTTCATTTCTTTATGTTTTAAATGCAGTACCTCCACATTCCTTTTATGATAAAATATTCAGAAATAATTTAAATACATCATT
1501 GCAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACC
1601 TTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGGC
141▶ N R T Y K L P I L E E I T T K V L K G

SacI (1745) **BstXI (1774)**
1701 ATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCA
121▶ N M E I L V F C D P A Y D S I L E R C M G C P S V V R I S R D V E
1801 TCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCAGCAGACCCAATGGCAATGGCTTCAGCAGACAGCAGTGAACC
87▶ D S Y P H R V A V I T D F D K Q G N S V A S G I A I A E A C V T V R

StuI (1909)
1901 TGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCTGATGGCCGCCCGACATGGTGTCTTGTCTCATAGAGCATGGT
54▶ G I Y A E I H V A S I I E G T K T R I A A G V H H K N D E Y L M T

BbsI (2055) **XmnI (2051)**
2001 GATCTTCTCAGTGGGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCC
21▶ I K E T A V E V L E L D Q Q S I N F T K M

AseI (2117) **SacI (2174)**
2101 GATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCAC

SpeI (2272)
2201 CGTACACGCTACCGCCATTTGCGTCAATGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGT
2301 CAATGGGTTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAA

SnaBI (2400)
2401 TACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGGGGCCATTTACCGTCAATTGACGTCAATAGGGGGCGTA

NdeI (2505)
2501 CTTGGCATATGATACACTTGTACTGCCAAGTGGGAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGG

PacI (2691)

SdaI (2683)

2601 GAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAG

BspLU11I (2701)

2701 AACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAA

2801 ATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCCCT

2901 GCCGTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCG

ApaLI (3015)

3001 TCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAAGTCCCAACCCGGTAAGACACGACTTAT

3101 CGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAG

3201 AAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAACCCCGCTGGTAGCGGT

3301 GGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAAACGAAA

EagI (3451)

PacI (3431) SwaI (3440) **NotI (3450)**

3401 ACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTT

3501 TTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAAACTAGCAAATAGGCTGTCCCCAGTGCAAGTGCAGGTGCCAGAA

3601 CATTCTCTATCGAA