



PvuI (7)
SgfI (6)
 1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA
 101 GAGAAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGTAAGTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
 201 GTGAACGTTCTTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGGCTCGCATCTCTCTTACAGCGCCCGCCGCCCTACCTGAGGCC
 301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCTCCCGCCTGTGGTGCCTCTGAAGTGCCTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441) **NaeI (441)**
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTCTCAACTCTACGTCTTTGTTTCGTTT

NcoI (560) **BstEII (555)** **KasI (535)** **AgeI (552)** **BstXI (582)**
 501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGCGCCCTAGCTGAGATCACCGGTACCATGGAATCCAACACTCCATCCATCTGAATGGATCAGA
 1 M E S N Y S I H L N G S E

BamHI (638) **SpeI (696)**
 601 AGTGGTGGTTTATGATTCTACCATCTCCAGAGTCTGTGGATCCTCTCAATGGTGGTTGTCTCCATCACTTTCTCTTGGTGTGCTGGGCAATGGACTA
 13 V V V Y D S T I S R V L W I L S M V V V S I T F F L G V L G N G L

BspEI (719)
 701 GTGATTTGGGTAGCTGGATTCCGGATGCCACACTGTCAACTATCTGGTATCTGAATCTAGCATTGGCTGACTTTTTCTTTCACAGCAACTCTACCAT
 47 V I W V A G F R M P H T V T T I W Y L N L A L A D F S F T A T L P

DraIII (865) **BbsI (898)**
 801 TCCTTCTTGTGAAATGGCTATGAAAGAAAAATGGCCTTTTGGCTGGTCTGTGTAATAATAGTTCACATTGGTAGATGTAACCTGTTTGGAAAGTGT
 80 F L L V E M A M K E K W P F G W F L C K L V H I V V D V N L F G S V
 901 CTTCTTGATTGCTCTCATTGCCTTGACCGCTGCATTTGTGTTCTGCATCCAGTCTGGGCTCAGAACCACCGCACTGTGAGCCTGGCTAGGAAGTGGTT
 113 F L I A L I A L D R C I C V L H P V W A Q N H R T V S L A R K V V

Bsp120I (1003) **EcoRI (1064)** **BsrGI (1087)**
 1001 GTTGGCCCTGGATTTTGTCTGATTCTCACTTTGCCATTTTTATTTTCTGACTACTGTTAGAATTCCTGGAGGAGATGTGATTGTACATTCACT
 147 V G P W I F A L I L T L P I F I F L T T V R I P G G D V Y C T F N

BamHI (1102) **SphI (1193)**
 1101 TTGGATCCTGGGCTCAAACCTGATGAAGAAAAGTTGAACACAGCTATCACTTTTGTAACTAGAGGGATCATCAGGTTCTTTATTGGTTTCAGCATGCC
 180 F G S W A Q T D E E K L N T A I T F V T T R G I I R F L I G F S M P
 1201 CATGTCAATTGTTGCTGTTTGGATGGACTCATTGCTGCAAGATCAACAGAAGAACTTGTAAATCCAGCGCTCTTACGAGTCCCTTACAGCAGT
 213 M S I V A V C Y G L I A V K I N R R N L V N S S R P L R V L T A V
 1301 GTGGTTCCTTTTATCTGCTGGTTTCCCTTTCAGCTGTGGCCCTTTGGGACAGCTGGTTTAAAGAGACATTGCTTAGTGGTAGTTATAAAATTC
 247 V A S F F I C W F P F Q L V A L L G T V W F K E T L L S G S Y K I

HpaI (1411) **BspLU11I (1403)**
 1401 TTGACATGTTTGTAAACCAACAAGCTCATTGGCTTACTTCAATAGTTGTCTCAATCCGATGCTCTATGTTTTATGGCCAGGACTTTCGTGAGAGATT
 280 L D M F V N P T S S L A Y F N S C L N P M L Y V F M G Q D F R E R F

PstI (1589)
 1501 TATTCATTCCCTGCCTTATAGTCTTGAGAGAGCCCTGAGTGAGGATTCTGGTCAAACAGTATTCAAGCACCAGTCTACTTACCTCCTGCAGACATT
 313 I H S L P Y S L E R A L S E D S G Q T S D S S T S S T S P A D I

MscI (1649) **NheI (1643)**
 1601 GAGTTAAAGGCCCATGAAGCTGTGATAAGGGATGGTTTTGAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGA
 347 E L K A P •

HpaI (1781)
 1701 ATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAACAACAACAATTGCA

EcoRI (1877)
 1801 TTCATTTTATGTTTCAGGTTTCAGGGGAGGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTAAAATACAGCATAGCA
 1901 AAACCTTAACCTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTG
 2001 CAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTACTCTTCACTTTCTTTATGTTTAAATGCACTGACCTCCAC

SspI (2116) **SwaI (2130)**
 2101 ATTCCCTTTTATGTAATAATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTC
 2201 ATAATATCCCCAGTTTAGTAGTTGACTTAGGGAACAAGGAACCTTAAATAGAAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTGT
 141 • N R T Y

SacI (2391)
 2301 ACTTGAGGGGATGAGTTCTCAATGGTGGTTTTGACCAGCTTCCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCT
 136 K L P I L 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

BstXI (2420)
 2401 GCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTG
 103 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

StuI (2555)

2501 CTCACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCC
69 S 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

XmnI (2697)

2601 TGATGGCCGCCCGACATGGTGCTTGTGTCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAA
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

BspHI (2705)

BbsI (2701) **AseI (2763)**

2701 GGTCTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGC T
3 T K M

SacI (2820)

2801 TATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTT

SpeI (2918)

2901 GGAAAGTCCC GTT GATTACTAGTCAAAA CAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCCA

SnaBI (3046)

3000 TTGATGTA CTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTA CTGCCAAGTAGGAAAGTCCCATAAGGTCATGTA CTGGGCAT

NdeI (3151)

3100 AATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGTA CTGCCAAGTGGGCAGTTTACCGTAAATA

3200 CTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGTCTGTTGGCGGT CAGCCA

PstI (3330) **SdaI (3329)** PacI (3337) **BspLU11I (3347)**

3300 GCGGGCCATTTACCGTAAGTTATGTAACGCTG CAG G TT AA TTAAGAACATGTGAGCAA AAGGCCAGCAA AAGGCCAGGAACCGTAAA AAGGCCGCG

3398 TTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATC AAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACC

3498 AGGCGTTTTCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTTCTCCCTTCGGGAAGCGTGGCGCT

ApaLI (3661)

3598 TTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTT CAGCCGACCGCTGCGCC

3698 TTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAG

3798 GCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAA

3898 AAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAA

PacI (4077) **SwaI (4086)** **NotI (4096)**

3998 GAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAACGAAA ACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AG

EagI (4097)

4098 CGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAA

4198 AACAAACTAGCAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA