



EcoRI (23) **SdaI (38)**
NotI (2) **XbaI (19)** **BamHI (29)** **SpeI (45)**
1 **CGCGCCGCGTCGACGATATCTAGAATTCGGATCCTGCAGGGCCACTAGTGGAGCCGAGAGTAATTCATACAAAAGGACTCGCCCCTGCCTTGGGGAATC**
101 **CCAGGGACCGTCGTTAAACTCCCACTAACGTAGAACCAGAGATCGCTGCGTTCCCGCCCCCTCACCCGCCGCTCTCGTCATCACTGAGGTGGAGAAGA**
AgeI (289)
201 **GCATGCGTGAGGCTCCGGTTCCTCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGACAAGTTGGGGGAGGGGTCGGCAATTGAACCGTGCCTAG**
301 **AGAAGGTGGCGCGGGTAAACTGGGAAAGTGATGTCGTGTACTGGCTCCGCTTTTTCCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCCG**
401 **TGAACGTTCTTTTCGCAACGGGTTTGCCGCCAGAACACAGGtaagtgccgtgtgtggttcccggggcctggcctctttacgggttatggccttgcgt**
BstXI (522)
501 **gccttgaattacttccatgcccttggctgcagtagctgattcttgatcccagcctcgggttgaagtgggtgggagagttcgaggccttgcgcttaagg**
601 **agcccccttgcctcgtgcttgagttgaggcctggcttggcgctggggccgcccgtgctaactcgttggcaccttgcgcctgtctcgctgcttgcgt**
PvuII (734) **BglII (781)**
701 **aagtctctagccatttaaaatTTTTgataaccagctgcgacgctTTTTtctggcgagatagcttgtaaatcgggccaagatctgcacactgggtattt**
801 **cggtttttggggcgcggggcggcagcggggcccgtgctcccagcgcacatgttcggcgaggcggggcctgcgagcgcggccaccgagaatcggacgggg**
901 **gtagtctcaaaactggccggcctgctctgggtgcctggcctcgcgccgctgtatcggcccctgggaggcaaggctggcccggctggcaccagttgcg**
SacI (1043)
1001 **tgagcggaaagatggccgcttcccggcctgctgcaggagctcaaaaaggaggacgcggcgcccgggagagcgggggggtgagtcacccacacaaaagga**
XhoI (1180)
1101 **aaaggcctttccttctcatccgctcgttcatgtgactccacggagtaccgggcgctccaggcacctcgattagttctcagagcttttggagtacgtc**
1201 **gtcttttaggttgggggaggggttttatgcatggagtttcccacactgagtggtggagactgaagagttaggccagcttggcacttgatgtaattct**
1301 **ccttggaaattgcccTTTTgagttggatcttgctcattctcaagcctcagacagtggttcaaaagtttttcttccatttcagGTGTCGTGAAAAC**
NcoI (1415)
1401 **ACCCCTAAAAGCCACCATGGAAATCAAGGTGCTGTTTGCCTCATCTGTATTGCTGTTGCTGAGGCAAAAACCACTGAAATCAATGAAGACCTCAATATA**
M E I K V L F A L I C I A V A E A K P T E I N E D L N I
BglII (1530)
1501 **GCTGCTGTGGCCTCCAACCTTTGCCACCACAGATCTTGAGACTGACCTGTTACCAACTGGGAGACCATGAATGTGATTAGCACTGACACAGAGCAGGTGA**
29▶ **A A V A S N F A T T D L E T D L F T N W E T M N V I S T D T E Q V**
1601 **ACACAGATGCTGACAGGGGCAAGCTGCCTGGCAAAAACTCCCCCAGATGTCTGAGGGAGCTGGAGGCAATGCCAGAAGGGCTGGTTGCACAAGAGG**
62▶ **N T D A D R G K L P G K K L P P D V L R E L E A N A R R A G C T R G**
1701 **CTGCCTCATTTGCCTCTCCACATTAAGTGCACCCCTAAGATGAAGAAATTTATCCCTGGCAGGTGCCACACTTATGAAGGTGAAAAGGAGTCTGCTCAG**
95▶ **C L I C L S H I K C T P K M K K F I P G R C H T Y E G E K E S A Q**
1801 **GGAGGATTGGAGAGGCAATTTGATATCCCAGAGATTCCTGGCTCAAGGATAAGGAGCCACTGGACCAGTTTATTGCTCAAGTGGACCTCTGTGCTG**
129▶ **G G I G E A I V D I P E I P G F K D K E P L D Q F I A Q V D L C A**
1901 **ATTGCACCACTGGCTGTCTGAAGGGCCTTGCCAATGTCCAGTGTCTGACCTCTGAAGAAGTGGCTTCCCAGAGGTGTACCACTTTTGCAGCAAGAT**
162▶ **D C T T G C L K G L A N V Q C S D L L K K W L P Q R C T T F A S K I**
NheI (2049)
2001 **TCAGGGTAGGGTGGACAAAATCAAGGGTCTGGCTGGGGACAGATGATAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAAACCACAA**
195▶ **Q G R V D K I K G L A G D R •**
2101 **CTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAACAAGTTAACAAACAA**
2201 **TTGCATTCATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTAATTCTAAAATAC**
SspI (2528)
2301 **AGCATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATT**
2401 **AGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGGTTTGAAGTACTCTTCATTCTTTATGTTTTAAATGCACTG**
SgrAI (2956)
2501 **ACCTCCCACATTCCTTTTTAGTAAAATTTAGAAAATTTAAATACATCATTGCAATGAAAATAATGTTTTTATTAGCAGAATCCAGATGCTCA**
2601 **AGGCCCTCATAATATCCCCAGTTTGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTTGGACAGCAAGAAAGCGAGCTCTAGCTTATCC**
2701 **TCAGTCTGCTCCTCTGCCACAAAGTGCACGCAGTTGCCGGCCGGTTCGCGCAGGGCGAACTCCCCCCCCACGGCTGCTCGCCGATCTCGGTGATGGCC**
125▶ **• D Q E E A V F H V C N G A P D R L A F E R G W P Q E G I E T M A**
2801 **GGCCCGGAGGCGTCCCGAAGTTCGTGGACACGACCTCCGACCACTCGGCGTACAGCTCGTCCAGGCCGCGCACCCACACCAGGCCAGGGTGTGTCGG**
91▶ **P G S A D R F N T S V V E S W E A Y L E D L G R V W V W A L T N D P**
2901 **GCACCACCTGGTCTGGACCGCTGATGAACAGGGTCACGTCGTCGCCGACCACACCGGCGAAGTCTCCTCCACGAAGTCCCGGGAGAACCCGAGCCG**
58▶ **V V Q D Q V A S I F L T V D D R V V G A F D D E V F D R S F G L R**

3001 GTCGGTCCAGAACTCGACCGCTCCGGCGACGTCGCGCGGGTGAGCACGGAAACGGCACTGGTCAACTGGCCATGATGGCTCCTCCTGTCAGGAGAGGA
25 D T W F E V A G A V D R A T L V P V A S T L K A M
3101 AAGAGAAGAAGGTTAGTACAATTGCTATAGTGAGTTGTATTATACTATGCAGATATACTATGCCAATGATTAATTGTCAAAGTAGGGCTGCAGGTTAATT
3201 AAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACA
3301 AAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCGCTCCTCTGTTCCGAC
3401 CCTGCCGCTTACCGGATACCTGTCCGCCTTCTCCCTTCGGGAAGCGTGGCGCTTTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTT
3501 CGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACT
3601 TATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACAC
3701 TAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGC
3801 GGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAAACG
3901 AAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCA