



2601 **SspI (2623)** **Swal (2637)** **EcoO109I**
CTCCACATTCCCTTTTATAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAAATAATGTTTTTATTAGGCAGAATCCAGATGCTCAAG

2701 **SacI (2898)**
GCCCTTCATAATATCCCCAGTTTGTAGTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTC
141 • N

2801 **BstXI (2927)**
CTGGTGTACTTGAGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGA
138 R T Y 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

2901 **StuI (3062)**
GCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTG
105 E R 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

3001 **BbsI (3208)**
CCCCTTGCTCACAGCAGACCCAATGGCAATGGCTTACGACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAAGT
72 G N 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

3101 **XmnI (3204)** **AseI (3270)**
TTGGTCTGATGGCCGCCCGACATGGTGTGTTGTCCCTCATAGAGCATGGTATCTTCTCAGTGGCAGCTCCACCAGCTCCAGATCTGCTGAGAGA
38 K T R 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

3201 **SacI (3327)**
TGTTGAAGGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTC
5 N F T K M

3301 **SpeI (3425)**
TCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACG

3401 **SnaBI (3553)**
ACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAACAAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCGTGAGTCAAACCGCTATCCA

3501 **NdeI (3658)**
CGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACT

3601 **PacI (3844)**
GGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGCGTACTTGGCATATGATACACTTGTACTGCCAAGTGGGCAGTTTACCG

3701 **PstI (3837)** **SdaI (3836)** **BspLU11I (3854)**
TAAATACTCCACCATTGACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCAATTATTGACGTCAATGGGGGGGGTCTTGGGGCGT

3801 **EagI (4604)** **NotI (4603)**
CAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCAGGTTAAATAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGC

3901 **ApaLI (4168)**
CGCGTTGCTGGCGTTTTTCCATAGGCTCCGCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGA

4001 **EagI (4604)** **NotI (4603)**
TACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGGCTCTCTGTTCGACCCTGCCGTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGG

4101 **PacI (4584)** **Swal (4593)**
CGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCCGGTGTAGGTGCTTCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTACGCCGACCGCTG

4201 **EagI (4604)** **NotI (4603)**
CGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTAT

4301 **EagI (4604)** **NotI (4603)**
GTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCG

4401 **EagI (4604)** **NotI (4603)**
GAAAAAGAGTTGGTAGTCTTGATCCGGCAACAAACCACCGTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATC

4501 **EagI (4604)** **NotI (4603)**
TCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAA

4601 **EagI (4604)** **NotI (4603)**
TCAGCGCCGCAATAAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAA

4701 **EagI (4604)** **NotI (4603)**
ACAAAACAAACTAGCAAAATAGGCTGTCCCAAGTCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA