



2401 **SacI** (2406) **BstXI** (2435)
AGAGATGAGCTCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAG
108 S I L 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

2501 **StuI** (2570)
TCCTTCTGCCCGTTGCTCACAGCAGACCAATGGCAATGGCTTCCAGCAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCT
74 D K Q 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

2601 **BbsI** (2716)
CCCGAGTCTTGGTCTGATGGCCGCCCGACATGGTGTCTGTGTCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTG
41 G T 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

2701 **XmnI** (2712) **AseI** (2778)
CTGAGAGATGTTGAAGGCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGG
8 Q S I N F T K M

2801 **SacI** (2835)
ATGGCGTCTCCAGCTTATCTGACGGTCTACTAAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAG

2901 **SpeI** (2933)
TTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAACAACTCCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACC

3001 **SnaBI** (3061)
GCTATCCACGCCATTGATGTACTGCCAAAACGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGT

3101 **NdeI** (3166)
CATGTAAGTGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTACTGCAAGTGGGCA

3201 **PacI** (3352)
GTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGTCGT

3301 **PstI** (3345) **SdaI** (3344) **BspLU11I** (3362)
TGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAAITTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGT

3401 **EagI** (4112)
AAAAAGGCCGCGTTGCTGGCGTTTTTTCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCAAACCCGACAGGAC

3501 **SwaI** (4101) **NotI** (4111)
TATAAAGATACCAGGCGTTTCCCCTGGAAGTCCCTCGTGGCTCTCTGTTCGACCTGCCGTTACCGGATACCTGTCCGCTTCTCCCTTCGGG

3601 **ApaLI** (3676)
AAGCGTGGCGTTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTGTTGCTCCAAGCTGGGTGTGTGCACGAACCCCGTTCAGCCC

3701 **EagI** (4112)
GACCGTGGCGCTTATCCGTAAGTATCGTCTTGTAGTCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAG

3801 **EagI** (4112)
CGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGT

3901 **EagI** (4112)
TACCTTCGAAAAAGAGTTGGTAGTCTTGTATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAA

4001 **PacI** (4092)
AAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAA

4101 **EagI** (4112)
CATTTAAATCAGCGGCCGAATAAAATATCTTTATTTTATTACATCTGTGTGGTTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAC

4201 **EagI** (4112)
AAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCAGTCAAGTGCAGGTGCCAGAACATTTCTATCGAA