



100



2201 ACCTCCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAAACTC **SpeI (2281)**

2300 CCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTA CTG CCAAAACCGCATCATCATGGTAATAG

**SnaBI (2409)**  
**Eco105I (2409)**

2400 CGATGACTAATACGTAGATGTACTGCCAAGTAGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAAT

**NdeI (2514)**

2500 AGGGGGCGTACTTGGCATATGATACACTTGATGTA CTG CCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGC

**PstI (2693)**  
**SdaI (2692)**

2600 GTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCCTGCA

**PacI (2700)** **BspLU11I (2710)**

2699 **G G T T A A T T A A G A A C A T G T G A G C A A A A G G C C A G C A A A A G G C C A G G A A C C G T A A A A A G G C C G C G T T G C T G G C G T T T T C C A T A G G C T C C G C C C C C T G A C C**

2798 **A G C A T C A C A A A A A T C G A C G C T C A A G T C A G A G G T G G C G A A A C C C G A C A G G A C T A T A A A G A T A C C A G G C G T T T C C C C T G G A A G C T C C C T C G T G C G C T C T C C**

2898 **T G T T C C G A C C C T G C C G T T A C C G G A T A C C T G T C C G C T T T C T C C C T T C G G A A G C G T G G C G C T T T C T C A T A G C T C A C G C T G T A G G T A T C T C A G T T C G G T G**

**ApaLI (3024)**

2998 **T A G G T C G T T C G C T C C A A G C T G G G C T G T G T G C A C G A A C C C C C G T T C A G C C C G A C C G C T G C G C T T A T C C G G T A A C T A T C G T C T T G A G T C C A A C C C G G T A A**

3098 **G A C A C G A C T T A T C G C C A C T G G C A G C A G C C A C T G G T A A C A G G A T T A G C A G A G C G A G G T A T G T A G G C G G T G C T A C A G A G T T C T T G A A G T G G T G G C C T A A C T A**

3198 **C G G C T A C T A G A A G A A C A G T A T T T G G T A T C T G C G C T C T G C T G A A G C C A G T T A C C T T C G A A A A A G A G T T G G T A G C T C T T G A T C C G G C A A A C A A A C C A C C**

3298 **G C T G G T A G C G G T G G T T T T T T T T G T T T G C A A G C A G C A G A T T A C G C G C A G A A A A A A A G G A T C T C A A G A A G A T C C T T T G A T C T T T T C T A C G G G G T C T G A C G C T C**

**EagI (3460)**  
**PacI (3440)** **Swal (3449)** **NotI (3459)**

3398 **A G T G G A A C G A A A A C T C A C G T T A A G G G A T T T T G G T C A T G G C T A G T T A A T T A A C A T T T A A A T C A G C G G C C G C A A T A A A A T A T C T T T A T T T T C A T T A C A T C T G**

3498 **T G T G T T G G T T T T T T G T G T G A A T C G T A A C T A A C A T A C G C T C C C A T C A A A A C A A A A C G A A A C A A A A C A A A C T A G C A A A A T A G G C T G T C C C C A G T G C A A G T G**

3598 **C A G G T G C C A G A A C A T T T C T A T C G A A**