



2301 CTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCTTCTTTCATGGAG

2401 TTTAAGATATAGTGATTTTTCCAAGGTTTGAAGTCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCTTTTTAGTAAATATTC **SspI (2493)**

2501 AGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGT **Swal (2507)** **EcoO109I (2568)**

2601 TGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGAGGGGGATGAGTTCCTCA **141** • N R T Y K L P I L E E

2701 ATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGCTGACCA **SacI (2768)** **BstXI (2797)**

128 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 C M G C P S V V

2801 CCCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAAT **95** 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 S V A S G I A I

2901 GGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCCTCAATGTGGACAGCAGAGATGATCTCCCCAGTCTTGGTCTGTAGGCCGCCCGACATGGTGC **StuI (2932)**

62 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 I A G V H H

3001 TTGTTGCTCATAGAGCATGGTGTCTCTCAGTGGCGACCTCCACCAGTCCAGATCCTGTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTAT **BbsI (3078)** **XmnI (3074)**

28 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 T K M

3101 AGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCCTAAACGAG **AseI (3140)** **SacI (3197)**

3201 CTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAG **SpeI (3295)**

3301 TCAAAAACAACTCCCATTGACGTCATGGGGTGGAGACTTGAAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATC

3401 ATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAAGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGT **SnaBI (3423)**

3501 CATTGACGTCATAGGGGGCTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCGAGTTTACCGTAAATACTCCACCATTGACGTCATGGAA **NdeI (3528)**

3601 AGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCATGGGGGGGGTCTTGGGCGGTGAGCCAGGCGGGCCATTTACCGTAAAGTTAT

3701 GTAACGCTGCAGGTTAAATAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCG **PacI (3714)** **PstI (3707)** **SdaI (3706)** **BspLU11I (3724)**

3801 CCCCCGTGACGAGCATCAGAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTC

3901 GTGCGCTCTCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTCGGAAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATC

4001 TCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTGAGCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTC **ApaI (4038)**

4101 CAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGG

4201 TGGCTAACTACGGTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCA

4301 AACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGG **EagI (4474)** **NotI (4473)**

4401 GTCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTT **PacI (4454)** **Swal (4463)**

4501 CATTACATCTGTGTGGTTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAATAGGCTGTCCC

4601 CAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA