





3401 CAGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCCTGATGGCCGCCCGACATGGTGCTTGTTCCTCATA  
57↓I Thr Val Arg Gl y I e Tyr Al a Gl u l e Hi s Va l Al a Ser I l e l l e Gl u Gl y Thr Lys Thr Arg l e Al a Al a Gl y Va l Hi s Hi s Lys Asn Asp Gl u Tyr  
BspHI (3569)  
XmnI (3565)  
3501 GAGCATGGTGATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCTGCTGAGAGATGTTGAAGGCTTTCATGATGGCCCTCTATAGTGAGTCGTATTA  
24↓Leu Me t Thr l l e Lys Gl u Thr Al a Va l Gl u Va l Leu Gl u Leu Asp Gl n Gl n Ser l l e Asn Phe Thr Lys Me t  
AseI (3628)  
3601 TACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACATAACGAGCTCTGCTTATATAG  
SpeI (3782)  
3701 ACCTCCACCCTACACGCCCTACCGCCATTTGCGTCAATGGGGCGGAGTGTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCC  
3801 CATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGC  
3901 GATGACTAATACGTAGATGTACTGCCAAGTAGAAAGTCCATAAGGTCATGTAAGTGGCATAATGCCAGGCGGGCATTACCGTCAATTGACGTCAATA  
4001 GGGGGCGTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGCGAGTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCG  
SdaI (4198)  
4101 TTA CTATGGGAACATACGTCATTATTGACGTCAATGGCGGGGGTCTGGTGGCGGTCAGCCAGGCGGGCCATTACCGTAAAGTTATGTAACGCCTGCAGG  
PaeI (4205) BspLU11I (4211)  
4201 TTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCTTGCTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGC  
4301 ATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTGTCGCTCTCCTGT  
4401 TCCGACCCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAG  
4501 GTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTCAGCCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGTAAGAC  
4601 ACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGG  
4701 CTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTTTGATCCGGCAAAACAAACCACCGCT  
4801 GGTAGCGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGT  
PacI (4945) SmaI (4953) NotI (4961)  
4901 GGAACGAAAACCTACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGT  
5001 GTTGGTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCAGTGAAGTGCAG  
5101 GTGCCAGAACATTTCTCTATCGAA