



2701 CCATGAAGAAGATGATGAAGAAACTATGAGAATGTACCACGTGTATTACTGGCCTCAGACCTAGCCCTTACCCAGAGTGGCCACAGGACTAGTCC
 713▶ H E E D D E E N Y E N V P R V L L A S D H •
 2801 TAGCTCGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTTATTGTGAAATTTGTGATGCTATTGCTT

2901 TATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAAACAAGTTAACAAACAACAAATTGCATTCATTTTATGTTTCAGGTT
 HpaI (2961) MfeI (2972)
 3001 CAGGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACAAATGTGGTATGGAATTCATAAATACAGCATAGCAAACTTTAACCTCCAATCA
 3101 AGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATG

3201 GAGTTTAAGATATAGTGATTTTCCCAAGGTTTGAAGTACTCTTCATTTCTTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTATGAAAAATA
 SapI (3239) SspI (3296)
 3301 TTCAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCCAGTTTAGT
 SwaI (3310)
 3401 AGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCCTGGTACTTGAGGGGGATGAGTTCC
 141◀ • N R T Y K L P I L E
 3501 TCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTCTGCACATGCCACAGGGGCTGA
 129▶ 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 E R C M G C P S V
 BstXI (3600)
 3601 CCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCCTTCTGCCGTTGCTCACAGCAGACCAATGGC
 96▶ V 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
 3701 AATGGCTTCAGCACAGACAGTACCTGCCAATGTAGGCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTGGTCTGATGGCCGCCCGACATGG
 63▶ I 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 A G V H
 XmnI (3877)
 3801 TGCTTGTTCCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTATGGTGGCCCTCC
 29▶ 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 N F T K M ◀◀

3901 TATAGTGAGTCGTATTATACTATGCCGATATACTATGCGGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCACAAAC
 AseI (3943)
 4001 GAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCCTGTTGATTTAC
 SacI (4000) SpeI (4098)
 4101 TAGTCAAAACAACTCCATTGACGTCAATGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGTATCCACGCCATTGATGTAAGTCCAAAACCGC

4201 ATCATCATGGTAATAGCGATGACTAATACGTAGATGACTGCCAAGTAGGAAAGTCCATAAGGTCATGTAAGTGGCATAATGCCAGGGGGCCATTAC
 SnaBI (4226)
 4301 CGTCATTGACGTCAATAGGGGGGCTACTTGGCATATGATACACTTGTACTGCCAAGTGGGCGAGTTACCGTAAATACTCCACCCATTGACGTCAATG
 NdeI (4331)
 4401 GAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGGGCGGGGCTGTTGGGCGGTGAGCCAGGCGGGCCATTACCGTAAGT

4501 TATGTAACGCTGACAGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGTTGCTGGCGTTTTTCCATAGGCT
 SdaI (4509) PacI (4517) BspLU11I (4527)
 4601 CCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCTGGAAGCTCC
 4701 CTCGTGCGCTCTCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGT

4801 ATCTCAGTTCGGTGTAGGTCGTTGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCGCTTATCCGGTAACTATCGTCTTGA
 ApaLI (4841)
 4901 GTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGCGGTGTACAGAGTCTTGAAG
 5001 TGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCG
 5101 GCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTAC

5201 GGGGCTGACGCTCAGTGAACGAAAACCTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC AGCGGCCCAATAAAATATCTTTAT
 Pacl (5257) SwaI (5266) EagI (5277) NotI (5276)
 5301 TTTTATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGT
 5401 CCCAGTGAAGTGCAGGTGCCAGAACATTTCTTATCGAA