



2701 AGGCCAGGTGTAGCTCTGAGTGTGGGCACTTTGCCCTGGACAGTGGGGCAGGTTTCAAGGCAGTGGCACTGCCACTCCTTACGCCCTTATTACCACC
713▶ G P G V A L S V G T L P L D S G A G S E G S G T A T P S A L I T T
2801 AATATGGTAGCCATGGAGCCATCTGTCCAGAGGCATTGCCCGTCTTCCAACAGTGGCATCAACGTATGCAGGTGGCAGATCTGCAGTCCATTAATA
747▶ N M V A M E A I C P E G I A R L A N S G I N V M Q V A D L Q S I N
BglIII (2880) AseI (2893)
MscI (2938)
2901 TCAGTGGCAATGGCTTCTGAGATCAGGCACCCGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAA
780▶ I S G N G F •
HpaI (3070) MfeI (3081)
3001 AAAATGCTTTATTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAAGTTAAACAACAACATTGCATTCAATTTTATG
EcoRI (3166)
3101 TTTCAGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAGCATAGCAAACCTTTAAC
3201 TCCAAATCAAGCCTCTACTTGAATCCTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTGCAGCCTCACCT
SapI (3348)
3301 TCTTTCATGGAGTTAAGATATAGTGTATTTTCCCAAGTTTGAAGTACTGCTCTTCATTCTTTATGTTTTAAATGCAGTACCTCCACATTCCTTTTT
SspI (3405) SmaI (3419)
3401 AGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCC
3501 CAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTTAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTCTGGTGTACTTGAGGGGG
141▶ • N R T Y K L P
3601 ATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCAC
132▶ I L E 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
3701 AGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCCGTTGCTCACAGCAGA
99▶ P S V 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
3801 CCCAATGGCAATGGCTTACGACAGACAGTGCACCTGCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGGTCTGATGGCCGCC
66▶ 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 G T K T R I A A
BbsI (3990)
XmnI (3986)
3901 CCGACATGGTGTGTTGTCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGG
32▶ 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 Q S I N F T K M
AseI (4052)
4001 TGGCCCTCTATAGTGAGTCTGATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGT
4101 TCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCCCG
SpeI (4207)
4201 TTGATTTACTAGTCAAAACAAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGC
SnaBI (4335)
4301 CAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGG
NdeI (4440)
4401 GCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGATGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTG
4501 ACGTCAATGGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTACGCCAGGCGGGCCATT
PacI (4626)
SdaI (4618) BspLU11I (4636)
4601 ACCGTAAGTTATGTAACGCCTGCAGGTTAATTAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTT
4701 CCATAGGCTCCGCCCCCTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCT
4801 GGAAGCTCCCTCGTGGCTCTCTGTTCCGACCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCAC
ApaLI (4950)
4901 GCTGTAGGTATCTCAGTTCGGTGTAGTCTGTTCCGCTCAAGTGGGCTGTGTGACGAAACCCCGTTCAGCCCAGCCGTGCGCTTATCCGGTAACTA
5001 TCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAG
5101 TTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCT
5201 CTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGAT
EagI (5386)
PacI (5366) SmaI (5375) NotI (5385)
5301 CTTTTCTACGGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCAATGGCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAA
5401 TATCTTTATTTTCAATACATCTGTGTGGTTTTTTTGTGTAATCGTAACTAACATACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCAAA
5501 ATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA