





MscI (2978)

2901 ATGTTAAGCTGGTTAACCCAGGCAGAAAGAACGTGTTACACACAGCAATAAGAATATCCACCTCAATGATGTTGCTAGCTGGCCAGACATGATAAGATACA  
780 N V K L V N Q A E G T C Y T Q Q •

3001 TTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTG

---

HpaI (3110) MfeI (3121)

3101 CAATAAACCAAGTTAACCAACAACAAATTGCATTCATTTTTATGTTTCAGGTTCAAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGT

3201 GGTATGGAATTCTAAAATACAGCATAGCAAACCTTAACTCCAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGG

---

SapI (3388)

3301 GGCTGTTGCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTAAGATATAGTGTATTTTCCAAGGTTTGAACCTAGCTCTTCATTTTC

---

SwaI (3459)

3401 TTTATGTTTTAAATGCACCTGCCACATTCCCTTTTTAGTAAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATT

3501 AGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAA

3601 AGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCATTTCATCTCAATGAGCACAAAGCAG  
141 • N R T Y K L P I L E E I T T K V L K G N M E I L V F C

SacI (3720)

3701 TCAGGAGCATAGTCAGAGATGAGCTCTGACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCA  
112 D P A Y D S I L E R C M G C P S V V R I S R D V E D S Y P H R V A V

StuI (3884)

3801 CAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGAC  
79 I T D F D K Q G N S V A S G I A I A E A C V T V R G I Y A E I H V

3901 AGCAGAGATGATCTCCCAAGTCTTGGTCTGATGGCCGCCGACATGGTCTTGTTCCTCATAGAGCATGGTGTCTTCTCAGTGGCGACCTCCACC  
46 A S I I E G T K T R I A A G V H H K N D E Y L M T I K E T A V E V

4001 AGCTCCAGATCCTGCTGAGAGATGTTGAAGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCGGATATACTATGCCGATGATTAATTG  
12 L E L D Q Q S I N F T K M

---

SacI (4149)

4101 TCAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTCACTAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTCGG

---

SpeI (4247)

4201 TCAATGGGGCGGAGTTGTTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGGAAATCC

---

SnaBI (4375)

4301 CCGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGA

---

NdeI (4480)

4401 AAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTA

4501 CTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCATTATTGACGTCAA

---

PstI (4659)  
SdaI (4658) PacI (4666)

4601 TGGGCGGGGTCGTTGGGCGTCCAGCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAGAACATGTGAGCAAAAGGCCAGCAAA

4701 AGGCCAGGAACCGTAAAAAGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCTTGACGAGCATCACAAAAATCGACGCTCAAGTCAGAGGTGGCG

4801 AAACCCGACAGGACTATAAGATACCAGGCTTTCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCC

---

ApaLI (4990)

4901 TTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGATGATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAGCTGGGCTGTGTGCACGAAC

5001 CCCCCTTCAGCCCGACCCTGCGCTTATCCGTAACATATCGTCTTGTAGTCCAAACCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAA

5101 CAGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTCTTGAAGTGGTGGCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCT

5201 CTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGTATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGA

5301 TTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTCAT

---

EagI (5426)  
PacI (5406) SwaI (5415) NotI (5425)

5401 GGCTAGTTAATTAACATTTAAATC AGCGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTTGTGTAATCGTAACTAACATACG

5501 CTCTCCATCAAAACAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA