



SgfI (11) EcoRI (16)  
1 GGATCTGCGATCGCTGAATTC TGGGGACTTTCCACTGGGGACTTTCCACTGGGGACTTTCCACTGGGGACTTTCCACTCCTGCAGC  
101 AGTGGATATTTCCAGAAAAC TTTTTGGATGCAGTTGGGGATTCTCTTTACTGGATGTGGACAATATCTCTATTATTACAGGAAGCAATCCCTCCT  
HindIII (256) NcoI (283)  
201 A T A A A A G G G C C T C A G C A G A A G T A G T G T T C A G C T G T T C T T G G C T G A C T T C A C A T C A A A G C T T C T A T A C T G A C C T G A G A C A G A G C C A T G G A A A T C A A G G T G C  
1 M E I K V BglIII (398)  
301 T G T T T G C C C T C A T C T G T A T T G C T G T T G C T G A G G C A A A C C C A C T G A A A T C A A T G A A G A C C T C A A T A T A G T G C T G T G G C C C A A C T T T G C C A C C A C A G A  
6 L F A L I C I A V A E A K P T E I N E D L N I A A V A S N F A T T D  
401 T C T T G A G A C T G A C C T G T T C A C C A A C T G G G A G A C C A T G A A T G T G A T T A G C A C T G A C A C A G A G C A G G T G A A C A C A G A T G C T G A C A G G G C A A G C T G C C T G G C  
39 L E T D L F T N W E T M N V I S T D T E Q V N T D A D R G K L P G  
501 A A A A A C T C C C C C A G A T G T C C T G A G G G A G C T G G A G G C C A A T G C C A G A A G G G C T G G T T G C A C A A G A G G C T G C C T C A T T T G C C T C T C C C A C A T T A A G T G C A  
73 K K L P P D V L R E L E A N A R R A G C T R G C L I C L S H I K C  
EcoRV (695)  
601 C C C T A A G A T G A A G A A T T T A T C C C T G G C A G G T G C C A C A C T T A T G A A G G T G A A A A G G A G T C T G C T C A G G G A G G G A T T G G A G A G G C A A T T G T T G A T A T C C C  
106 T P K M K K F I P G R C H T Y E G E K E S A Q G G I G E A I V D I P  
701 A G A G A T T C C T G G C T T C A A G G A T A A G G A G C C A C T G G A C C A G T T T A T T G C T C A A G T G G A C C T C T G T G C T G A T T G C A C C A C T G G C T G T C T G A A G G G C C T T G C C  
139 E I P G F K D K E P L D Q F I A Q V D L C A D C T T G C L K G L A  
801 A A T G T C C A G T G C T C T G A C C T C C T G A A G A A G T G G C T T C C C C A G A G G T G T A C C A C T T T T G C C A G C A A G A T T C A G G G T A G G G T G G C A A A A T C A A G G G T C T G G  
173 N V Q C S D L L K K W L P Q R C T T F A S K I Q G R V D K I K G L  
NheI (917)  
901 C T G G G G A C A G A T G A T A G C T A G C T G G C C A G A C A T G A T A A G A T A C A T T G A T G A G T T T G G C A A A C C A C A A C T A G A A T G C A G T G A A A A A A T G C T T T A T T T G T  
206 A G D R •  
HpaI (1057)  
1001 G A A A T T T G T G A T G C T A T T G C T T A T T T G T A A C C A T T A A A G C T G C A A T A A A C A A G T T A A C A C A A C A A C A A T T G C A T T C A T T T T A T G T T T C A G G T T C A G G G G  
EcoRI (1151)  
1101 A G T G T G G G A G G T T T T T A A A G C A A G T A A A A C C T C T A C A A A T G T G G T A T G G A A T T C T A A A A T A C A G C A T A G C A A A A C T T T A A C T C C A A A T C A A G C C T C T  
1201 A C T T G A A T C C T T T T C T G A G G G A T G A A T A A G G C A T A G G C A T C A G G G G C T G T T G C C A A T G T G C A T T A G C T G T T T G C A G C C T C A C T T C T T T C A T G G A G T T A  
1301 A G A T A T A G T G T A T T T T C C A A G G T T T G A A C T A G C T C T C A T T T C T T A T G T T T A A A T G C A C T G A C C T C C C A C A T T C C C T T T T A G T A A A A T A T T C A G A A  
1401 A T A A T T T A A A T A C A T C A T T G C A A T G A A A A T A A A T G T T T T T A T T A G G C A G A A T C C A G A T G C T C A A G G C C C T T C A T A A T A T C C C C C A G T T T A G T A G T T G G A  
1501 C T T A G G G A C A A A G G A A C C T T T A A T A G A A A T T G G A C A G C A A G A A A G C G A G C T T C T A G C T T T A G T T C C T G G T G A C T T G A G G G G G A T G A G T T C C T C A A T G G  
141 • N R T Y K L P I L E E I T  
SacI (1669)  
1601 T G T T T T G A C C A G C T T G C C A T T C A T C T C A A T G A G C A C A A A G C A G T C A G G A G C A T A G T C A G A G A T G A G C T C T C G C A C A T G C C A C A G G G C T G A C C A C C T  
127 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 R  
1701 G A T G G A T C T G C C A C C T C A T C A G A G T A G G G G T G C C T G A C A G C C A C A A T G G T G T C A A A G T C C T T C T G C C G T T G C T C A C A G C A G A C C C A A T G G C A A T G G C T  
94 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 A  
StuI (1831)  
1801 T C A G C A C A G A C A G T G A C C C T G C C A A T G T A G G C C T C A A T G T G G A C A G C A G A G A T G A T C T C C C C A G T C T T G G T C C T G A T G G C C G C C C C A C A T G G T G C T T G T  
60 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 H K N  
1901 T G T C C T C A T A G A G C A T G G T G A T C T T C A G T G G C G A C C T C C A C C A G C T C C A G A T C C T G C T G A G A G A T G T T G A A G G T C T C A T G G T G G C C C T C C T A T A G T G  
27 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 (2038)  
2001 A G T C G T A T T A T A C T A T G C C G A T A T A C T A T G C C G A T G A T T A A T T G T C A A C T A C T G T T T G T A G G C C G G T C A C A G C T T G A T C T G T A A C G G C G C A G A A C A G  
2101 A A A A C G A A A C A A A G A C G T A G A G T T G A G C A A G C A G G G T C A G G C A A A G C G T G G A G A G C C G G C T G A G T C T A G G T A G G C T C C A A G G G A G C C G G A C A A A G G C C  
2201 C G G T C T C G A C C T G A G C T T T A A A C T T A C C T A G A C G G C G G A C G C A G T T C A G G A G G C A C C A C A G G C G G G A G G C G G C A G A A C G C G A C T C A A C C G G C T G G A T G G  
HindIII (2351) Psp1406I (2394)  
2301 C G G C C T C A G G T A G G G C G G C G G G C G T G A A G G A G A G A T G C G A G C C C C T C G A A G C T T C A G C T G T G T T C T G G C G G C A A C C C G T T G C G A A A A A G A A C G T T C A  
2401 C G G C G A C T A C T G C A C T T A T A T A C G G T T C T C C C C A C C C T C G G G A A A A G G C G G A G C C A G T A C A C G A C A T C A C T T T C C A G T T A C C C C G C C A C C T T C T  
AgeI (2507) SpeI (2576)  
2501 C T A G G C A C C G G T T C A A T T G C C G A C C C C T C C C C C A A C T T C T C G G G A C T G T G G G C A T G T G C G C T C T G C C C A C T G A C T A G T G G G C C C T G C A G G T T A A T T A

**BspLU11I (2604)**

2601 AGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGCGTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAA  
2701 AAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCTGTCCGACC  
2801 CTGCCGCTTACCGGATACCTGTCCGCCTTTCTCCCTTCGGGAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTT  
2901 GCTCCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCCCGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTT  
3001 ATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACT  
3101 AGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCG  
3201 GTGGTTTTTTGTTTGAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGAACGA

**NotI (3354)**

3301 AA~~ACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCG~~CAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTT  
3401 TTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACAAAACAAACTAGCAAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAG  
3501 AACATTTCTCTATCGAA