



**PvuI (7)**  
**SgfI (6)**
**MfeI (82)**

1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA  
 101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGATGTCGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

---

**Psp1406I (203)**
**HindIII (245)**

201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCCTTACGCGCCCGCCCTACCTGAGGCC  
 301 GCCATCCACGCGGGTTGAGTCGCGTTCTGCGCCCTCCCGCTGTGGTGCTCCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC  
 401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCACTCTACGTCTTTGTTTCGTTT

---

*NcoI (560)*

**BstEII (555)**

501 TCTGTTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCCCTA **KasI (535)** **AgeI (552)** *NcoI (579)* **BsrBI (594)**  
 601 CAAGCGGAGGAGCTGCTTCCAGGACAGGCAAGGACGGGAGCTTCTCGTGGTGGCCAGCGAGTCCATCTCCGGGCATACGCGCTCTGCGTGTGTAT  
 13▶r LysAl aGl uGl uLeuLeuSer ArgThr Gl yLysAspGl ySer PheLeuVal ArgAl aSer Gl uSer I l eSer ArgAl aTyrAl aLeuCysVal LeuTyr

**EcoRI (722)**

701 CGGAATTGCGTTTACACTTACAGAATTCTGCCAATGAAGATGATAAATCTACTGTTCAAGCATCCGAAGCGCTCTCCATGAGTTTCTCACCAAGCTGG  
 47▶ArgAsnCysVal TyrThr TyrArgI l eLeuProAsnGl uAspAspLysPheThr Val Gl nAl aSer Gl uGl yVal Ser Me tArgPhePheThr LysLeuA

**BstEII (840)**

801 ACCAGTCTCATCGAGTTTACAAGAAGGAAACATGGGGCTGGTACCCATCTGCAATACCCTGTGCCGCTGGAGGAAGAGGACACAGGCGACGCCCTGA  
 80▶spGl nLeuI l eGl uPheTyrLysLysGl uAsnMe tGl yLeuVal Thr Hi sLeuGl nTyrP roVal P roLeuGl uGl uGl uAspThr Gl yAspAspP roGl  
 901 GGAGGACACAGAAAGTGTGCTGTCTCCACCCGAGCTGCCCCAAAGAAACATCCCGCTGACTGCCAGCTCCTGTGAGGCCAAGGAGGTTCTTTTTCAAAC  
 113▶uGl uAspThr Gl uSer Val Val Ser P roP roGl uLeuP roP roArgAsnI l eP roLeuThr Al aSer Ser CysGl uAl aLysGl uVal P roPheSerAsn  
 1001 GAGAATCCCCGAGCGACCAGACAGCCGGCCGAGCCTCTCCGAGACATTGTTCCAGCGACTGCCAAAGCATGGACACCAGTGGCTTCCAGAAGAGCATC  
 147▶Gl uAsnP roArgAl aThr Gl uThr Ser ArgP roSer LeuSer Gl uThr LeuPheGl nArgLeuGl nSer Me tAspThr Ser Gl yLeuP roGl uGl uHi sL  
 1101 TTAAGGCCATCCAAGATTATTTAAGCACTCAGCTCGCCAGGACTCTGAATTTGTGAAGACAGGGTCCAGCAGTCTTCTCACCTGAAGAAACTGACCCAC  
 180▶euLysAl a l l eGl nAspTyrLeuSer Thr Gl nLeuAl aGl nAspSer Gl uPheVal LysThr Gl ySer Ser Ser LeuP roHi sLeuLysLysLeuThr Th

**BspEI (1232)**

1201 ACTGCTCTGCAAGGAGCTCTATGGAGAAGTCATCCGGACCCTCCCATCCCTGGAGTCTCTGCAGAGGTTATTTGACCAGCAGCTCTCCCCGGGCTCCGT  
 213▶r LeuLeuCysLysGl uLeuTyrGl yGl uVal I l eArgThr LeuP roSer LeuGl uSer LeuGl nArgLeuPheAspGl nGl nLeuSer P roGl yLeuArg

**SruI (1397)**

1301 CCACGTCTCAGGTTCTGGTGAGGCAATCCCATCAACATGGTGTCCAAGCTCAGCCAAGTACAAGCTGTGTGCTCCATTGAAGACAAGGTCAAGG  
 247▶P roArgP roGl nVal P roGl yGl uAl aAsnP ro l l eAsnMe tVal Ser LysLeuSer Gl nLeuThr Ser LeuLeuSer Ser I l eGl uAspLysVal LysA  
 1401 CCTTGCTGCACGAGGGTCTGAGTCTCCGCACCCGGCCCTCCCTATCCCTCCAGTACCTTTGAGGTGAAGGCAGAGTCTCTGGGATTCCTCAGAAAAT  
 280▶I l aLeuLeuHi sGl uGl yP roGl uSer P roHi sArgP roSer LeuI l eP roP roVal Thr PheGl uVal LysAl aGl uSer LeuGl yI l eP roGl nLysMe

**SaII (1510)**

1501 GCAGCTCAAAGTCGACGTTGAGTCTGGGAAACTGATCATTAAAGAGTCCAAGGATGGTCTGAGGACAAGTCTACAGCCACAAGAAAATCCTGCAGCTG  
 313▶tGl nLeuLysVal AspVal Gl uSer Gl yLysLeuI l eI l eLysLysSer LysAspGl ySer Gl uAspLysPheTyrSer Hi sLysLysI l eLeuGl nLeu  
 1601 ATTAAGTCACAGAAATTTCTGAATAAGTTGGTGATCTTGGTGGAAACAGAGAAGGAGAAGATCCTGCGGAAGGAATATGTTTTGCTGACTCCAAAAGA  
 347▶I l eLysSer Gl nLysPheLeuAsnLysLeuVal I l eLeuVal Gl uThr Gl uLysGl uLysI l eLeuArgLysGl uTyrVal PheAl aAspSer LysLysA

**BsaBI (1767)**

1701 GAGAAGGCTTCTGCCAGCTCCTGCAGCAGATGAAGAACAAGCACTCAGAGCAGCCGGAGCCGACATGATCACCATCTTCATCGGCACCTGGAACATGGG  
 380▶r gGl uGl yPheCysGl nLeuLeuGl nGl nMe tLysAsnLysHi sSer Gl uGl nP roGl uP roAspMe tI l eThr I l ePheI l eGl yThr TrpAsnMe tGl  
 1801 TAACGCCCCCTCCCAAGAAGTACAGTCTCGTCTTCTCCAAGGGCAGGAAAGACGCGGGACGACTCTGCGGACTACATCCCCATGACATTTAC  
 413▶yAsnAl aP roP roLysLysI l eThr Ser TrpPheLeuSer LysGl yGl nGl yLysThr ArgAspAspSer Al aAspTyrI l eP roHi sAspI l eTyr  
 1901 GTGATCGGCACCAAGAGGACCCCTGAGTGAGAAGGAGTGGCTGGAGATCCTCAAACACTCCCTGCAAGAAATCACCAGTGTGACTTTTAAACAGTCG  
 447▶Val I l eGl yThr Gl nGl uAspP roLeuSer Gl uLysGl uTrpLeuGl uI l eLeuLysHi sSer LeuGl nGl uI l eThr Ser Val Thr PheLysThr ValA

**MscI (2036)**

2001 CCATCCACACGCTCTGGAACATCCGCATCGTGGTGGCCAGCCTGAGCAGGAGAACCAGGATCAGCCACATCTGTACTGACAACGTGAAGACAGGCAT  
 480▶I al l eHi sThr LeuTrpAsnI l eArgI l eVal Val LeuAl aLysP roGl uHi sGl uAsnArgI l eSer Hi sI l eCysThr AspAsnVal LysThr Gl yI l  
 2101 TGCAAAACACTGGGGAACAAGGAGCCGTGGGGGTGTCGTTTCATGTTCAATGGAACCTCCTTAGGGTTCGTCACAGCCACTTGACTTCAGGAAGTGAA  
 513▶eAl aAsnThr LeuGl yAsnLysGl yAl aVal Gl yVal Ser PheMe tPheAsnGl yThr Ser LeuGl yPheVal AsnSer Hi sLeuThr Ser Gl ySer Gl u  
 2201 AAGAAACTCAGGCAAAACAAAATATATGAACATTCTCCGGTCTCGGCCCTGGGCGACAAGAAGCTGAGTCCCTTTAACATCACTCACCGCTTACGC  
 547▶LysLysLeuArgArgAsnGl nAsnTyrMe tAsnI l eLeuArgPheLeuAl aLeuGl yAspLysLysLeuSer P roPheAsnI l eThr Hi sArgPheThr H  
 2301 ACCTCTTCTGGTTTGGGATCTTAACACTACCGTGTGGATCTGCCTACCTGGGAGGCAGAAACCATCATCCAGAAAATCAAGCAGCAGCAGTACCGAGACCT  
 580▶i sLeuPheTrpPheGl yAspLeuAsnTyrArgVal AspLeuP roThr TrpGl uAl aGl uThr I l eI l eGl nLysI l eLysGl nGl nGl nTyrAl aAspLe

**XmnI (2439)**

2401 CCTGTCCCACGACAGCTGCTCACAGAGAGGAGGAGCAGAAGGTTCTTCTACACTTCGAGGAGGAAGAAATCACGTTTGCCCCAACCTACCGTTTTGAG  
 613▶uLeuSer Hi sAspGl nLeuLeuThr Gl uArgArgGl uGl nLysVal PheLeuHi sPheGl uGl uGl uI l eThr PheAl aP roThr TyrArgPheGl u  
 2501 AGACTGACTCGGGACAAAATACGCCTACACCAAGCAGAAAGCAGACGGATGAAGTACAACCTGCCTTCTGGTGTGACCGAGTCTCTGGAAGTCTTATC  
 647▶ArgLeuThr ArgAspLysTyrAl aTyrThr LysGl nLysAl aThr Gl yMe tLysTyrAsnLeuP roSer TrpCysAspArgVal LeuTrpLysSer TyrP

**DraIII (2608)**

**ApaLI (2605)**
**BspHI (2645)**

2601 CCCTGGTGACGTTGGTGTGTCAGTCTTATGGCAGTACCAGCGACATCATGACGAGTGACCACAGCCCTGTCTTTGCCACATTTGAGGCAGGAGTCACTTC  
 680▶r oLeuVal Hi sVal Val CysGl nSer TyrGl ySer Thr SerAspI l eMe tThr SerAspHi sSer P roVal PheAl aThr PheGl uAl aGl yVal Thr Se

**XcmI (2795)**

2701 CCAGTTTGTCTCCAAGAACGGTCCCGGACTGTTGACAGCCAAGGACAGATTGAGTTTCTCAGGTGCTATGCCACATTGAAGACCAAGTCCCAGACAAA  
 713▶r Gl nPheVal Ser LysAsnGl yP roGl yThr Val AspSer Gl nGl yGl nI l eGl uPheLeuArgCysTyrAl aThr LeuLysThr LysSer Gl nThr Lys

**XhoI (2817)**

2801 TTCTACCTGGAGTTCACCTCGAGCTGCTTGGAGAGTTTTGTCAAGAGTCAGGAAGGAGAAAATGAAGAAGGAAGTGAGGGGGAGCTGGTGGTGAAGTTTG  
747>PheTyrLeuGI uPheHi sSer Ser CysLeuGI uSer PheVal I LysSer GI nGI uGI yGI uAsnGI uGI uGI ySer GI uGI yGI uLeuVal I Val LysPheG  
2901 GTGAGACTCTTCCAAAGCTGAAGCCATTATCTCTGACCTGAGTACCTGCTAGACCAGCACATCCTCATCAGCATCAAGTCCCTGACAGCGACGAATC  
780>IyGI uThrLeuP roLysLeuLysP roI l eI l eSer AspP roGI uTyrLeuLeuAspGI nHi s l l eLeuI l eSer l l eLysSer Ser AspSer AspGI uSe

**NcoI (3075)**

3001 CTATGGCGAGGGCTGCATTGCCCTTCGGTTAGAGGCCACAGAAACGCAGCTGCCCATCTACACGCCTCTCACCCACCATGGGGAGTTGACAGGCCACTTC  
813>r TyrGI yGI uGI yCysI l eAl aLeuArgLeuGI uAl aThr GI uThr GI nLeuP roI l eTyrThr P roLeuThr Hi sHi sGI yGI uLeuThr GI yHi sPhe  
3101 CAGGGGGAGATCAAGCTGCAGACCTCTCAGGGCAAGACGAGGGAGAAGCTCTATGACTTTGTGAAGACGGAGCGTGATGAATCCAGTGGGCCAAAGACCC  
847>GI nGI yGI uI l eLysLeuGI nThr Ser GI nGI yLysThr ArgGI uLysLeuTyrAspPheVal I LysThr GI uArgAspGI uSer Ser GI yProLysThr L  
3201 TGAAGAGCCTCACAGCCACGACCCATGAAGCAGTGGGAAGTCACTAGCAGGGCCCCCTCCGTGACGTGGCTCCAGCATCACTGAAATCATCAACCCCAA  
880>euLysSer LeuThr Ser Hi sAspP roMe tLysGI nTrpGI uVal Thr Ser ArgAl aP roP roCysSer GI ySer Ser l l eThr GI uI l eI l eAsnP roAs

**SfiI (3314)**

3301 CTACATGGGAGTGGGGCCCTTTGGGCCACCAATGCCCTGCACGTGAAGCAGACCTTGTCCCCTGACCAGCAGCCACAGCCTGGAGCTACGACCAGCCG  
913>nTyrMet tGI yVal GI yProPheGI yProP roMe tP roLeuHi sVal I LysGI nThr LeuSer P roAspGI nGI nP roThr Al aTrpSer TyrAspGI nP ro

**SdaI (3422)**

3401 CCCAAGGACTCCCCGCTGGGGCCCTGCAGGGGAGAAAAGTCTCCGACACCTCCCGCCAGCCGCCATATCACCCAAGAAGTTTTTACCCTCAACAGCAA  
947>P roLysAspSer P roLeuGI yP roCysArgGI yGI uSer P roP roThr P roP roGI yGI nP roP roI l eSer P roLysLysPheLeuP roSer Thr Al aA  
3501 ACCGGGTCTCCCTCCAGGACACAGGAGTCAAGGCCAGTGACCTGGGGAAGAACGCAGGGGACACGCTGCCTCAGGAGGACCTGCCCTGACGAAGCC  
980>snArgGI yLeuP roP roArgThr GI nGI uSer ArgP roSer AspLeuGI yLysAsnAl aGI yAspThr LeuP roGI nGI uAspLeuP roLeuThr LysP r

**SandI (3625)**

**SacII (3685)**

3601 CGAGATGTTTGAGAACCCCTGTATGGTCCCTGAGTTCCTCCCTAAGCCTGCTCCCAGGAAGGACCAGGAATCCCCAAAATGCCCGGAAGAACCC  
1013>oGI uMe tPheGI uAsnP roLeuTyrGI ySer LeuSer Ser PheP roLysP roAl aP roArgLysAspGI nGI uSer P roLysMe tP roArgLysGI uP ro  
3701 CCGCCTGCCCGAACCCGGCATTTGTGCGCCAGCATCGTGTCAACAAAGCCAGGAGGCTGATCGCGCGAGGGGCCCGCAAGCAGGTGCCCGCGC  
1047>P roP roCysP roGI uP roGI yI l eLeuSer P roSer l l eVal LeuThr LysAl aGI nGI uAl aAspArgGI yGI uGI yP roGI yLysGI nVal I P roAl aP

**PshAI (3879)**

3801 CCCGGCTGCGCTCCTTACGTGCTCATCTCTGCGAGGGCAGGGCGCGGGGGACAAGAGCCAAGGGAAGCCCAAGACCCGGTACGCTCCAGGC  
1080>r oArgLeuArgSer PheThr CysSer Ser Ser Al aGI uGI yA rgAl aAl aGI yGI yAspLysSer GI nGI yLysP roLysThr P roVal Ser Ser GI nAl

**NotI (3975)**

**SacII (3973)**

3901 CCCGGTCCCGCCCAAGAGGCCCATCAAGCCTTCCAGATCGAAATCAACCAGCAGACCCCGCCACCCCGACGCCGGCGCCGCTGCCAGTCAAGAGC  
1113>aP roVal P roAl aLysArgP roI l eLysP roSer ArgSer GI uI l eAsnGI nGI nThr P roP roThr P roThr P roArgP roP roLeuP roVal I LysSer

**SfiI (4091)**

4001 CCGCGGTGCTGCACCTCCAGCACTCCAAGGGCCGCGACTACCGCGACAACACCGAGCTCCCGCATCACGGCAAGCACCGGCCGAGGAGGGGCCACCAAG  
1147>P roAl aVal LeuHi sLeuGI nHi sSer LysGI yArgAspTyrArgAspAsnThr GI uLeuP roHi sHi sGI yLysHi sArgP roGI uGI uGI yP roP roG  
4101 GGCCTCTAGGCAGGACTGCCATGCAGTGAAGCCTCAGTGAGCTGCCACTGAGTCGGGAGCCCAGAGGAACGGGTGAAGCCACTGGACCCTCTCCCGGG  
1180>IyP roLeuGI yA rgThr Al aMe tGI n●●●

**MscI (4250)**

**NheI (4244)**

4201 ACCTCCTGCTGGCTCCTCCTGCCAGCTTCTATGCAAGGCTTTGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAG

**HpaI (4382)**

**MfeI (4393)**

4301 AATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAACCAATTGC

**EcoRI (4478)**

4401 ATTCATTTTATGTTTCAGGTTTCAGGGGGAGGTGTGGGAGGTTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTTCTAAAATACAGCATAGC

4501 AAAACTTTAACCTCCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAGCTGTTT

4601 GCAGCCTCACCTTCTTTCATGGAGTTTAAGATATAGTGTATTTTCCAAAGTTTGAAGTACTGCTTTCATTTCTTTATGTTTTAAATGCACTGACCTCCCA

**SspI (4717)**

**SwaI (4731)**

4701 CATTCCCTTTTTAGTAAAATATTAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTT

4801 CATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTTGTTG

141<●●●AsnArgThr T

4901 TACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTC

136<y rLysLeuP roI l eLeuGI uGI uI l eThr Thr LysVal I LeuLysGI yAsnMetGI uI l eLeuVal I PheCysAspP roAl aTyrAspSer l l eLeuGI uAr

5001 TGCACATGCCACAGGGGTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTT

103<gCysMetGI yCysP roSer Val I Val I ArgI l eSer ArgAspVal I uAspSer TyrP roHi sArgVal I Al aVal I l eThr AspPheAspLysGI nGI yAsn

**StuI (5156)**

5101 GCTCACAGCAGACCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAAGTCTTGCTC

70<Ser Val I Al aSer GI yI l eAl aI l eAl aGI uAl aCysVal I Thr Val I ArgGI yI l eTyrAl aGI uI l eHi sVal I Al aSer l l eI l eGI uGI yThr LysThr A

**XmnI (5298)**

5201 CTGATGGCCGCCCGACATGGTCTTGTTCCTCATAGACATGGTATCTTCTCAGTGGCAGCTCCACCAGCTCCAGATCCTGCTGAGAGATGTTGA

36<r gl l eAl aAl aGI yVal Hi sHi sLysAsnAspGI uTyrLeuMe tThr l l eLysGI uThr Al aVal I GI uVal I LeuGI uLeuAspGI nGI nSer l l eAsnPh

**BspHI (5306)**

**AseI (5364)**

5301 AGGCTTTCATGATGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACAGCGTGGATGGCGTCTCCAGC

3<eThr LysMet

5401 T TATCTGACGGTTCACTAAACGAGCTCTGCTTATATAGACCTCCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGAGTTGTTACGACATTT

---

5501 TGGAAAGTCCCCTTGATTTACTAGTCAAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCCGTGAGTCAAACCGCTATCCACGCC

SpeI (5519) ←

---

5600 ATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCA

SnaBI (5647)

---

5700 TAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACACTTGTGTACTGCCAAGTGGGCAGTTTACCGTAAAT

NdeI (5752)

---

5800 ACTCCACCATTGACGTCAATGGAAAGTCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAATGGGCGGGGTCGTTGGGCGGTCAGCC

---

5900 AGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAA TTAAGAACATGTGAGCAAAAGGCCAGCAAAAGGCCAGGAACCGTAAAAAGGCCGC

SdaI (5930) PacI (5938) BspLU11I (5948) ←

---

5998 GTTGCTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATAC

---

6098 CAGGCGTTTCCCCTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGC

---

6198 TTTTCATAGCTCAGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCCAAGCTGGGCTGTGTGCACGAACCCCCGTTACGCCGACCGCTGCCG

ApaLI (6262)

---

6298 CTTATCCGGTAACTATCGTCTTGAGTCCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTA

---

6398 GCGGGTGTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGGAA

---

6498 AAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCCTGGTAGCGGTGGTTTTTTTGTGTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCA

---

6598 AGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAACGAAAACCTACGTTAAGGGATTTTGGTCATGGCTAGTTAATTAACATTTAAATC A

PacI (6678) SmaI (6687)

---

6698 GCGGCCGAATAAAATATCTTTATTTTATTACATCTGTGTGTTGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAACAAAACGAAACA

NotI (6697)

6798 AAACAACTAGCAAAATAGGCTGTCCCCAGTGAAGTGCAAGTGCCAGGTCAGAACATTTCTCTATCGAA