



Bsp120I (7) ScaI (18)
1 **SdaI (6)** **SpeI (13)** MfeI (98)
CCTGCAGGGCCCCTAGTACTATGCTGCACAAGCAATTTAAAAACCAACAGCAAAAAATACACTTCTCTGAAAAAGTCTTGGTCTAGGACCTAAACAA

101 **NaeI (187)**
TTGCTGAAACTGGGTAGACTTACACCAATGAGAGGCAGATAAAGAGATTAAGATTGAGGGAGTAGGGCAGGGCTTGCAATGGTGCCGGCCAGGATGTGG

201 CTGAGGGGGTGTGGGTGCCTGCCGTGGATGCTAGGGTAGAAGACGACTCTATTAAGTGGTGGCTGTAAGCAGTACCCAGGTCAATGCCTTTCATCTTCT

301 ACAACCTCGACGTTGCTGGAATCTAAATCTTTTTCTTCACTTAACAAACATCACCTCTGCTCAAATCTGCAACTGCTTTGATATCACACTGCCTTTTT

401 CACCCCTCTATTATAGATGGCATTATTTACTTACATGTTTTTCCCTACTAGACTATACTCCTTGAGAACAGCGATTGTGTCTTATTTATTTCTGAATC

501 **SphI (519)**
ACCAATTCAGACAGGCATGCAAACACTTGTGAACCAATGCACAAATATATTTTGTCTTCTTCATAGATTCTCCGGCCTCAGATGACCAGGCACCACT

601 **BspHI (659)**
AGATACAGAACACTGTGCTTTCCTTCTCCAAGTAAAGGAATAAATATCTGTTCCCTTCATGAAGTGTACTGTTGGGCCTTTATGCCATCCTGAAGCC

701 Bsp120I (793)
EcoNI (788)
ACCAGGATGTGGAACAGATCAGGGAGGTCACAGTTACAACCCCTTGTATCTGTAACACCAGCAGGACATTATCTACAGAGTCTGCTGCAGGGCCCCG

801 **NsiI (856)**
AATGAAGACAGCATTGCTGCTTTGTAGCGTGAGCAGTGCTGTAACAGTGATGCATGGATGTTCTCTGCTGCTGAAAGAATGTAGGTGCTTCTTGA

901 **AseI (918)**
AAGCTCTCTGCAACTTATTAATTGGGAGTGATTATGCGATGGAGAAAACAGAGTCCCATCACCCCTCAGTCTTCCCTGGGAAATCACAAGAGGGCTGAT

1001 **XmnI (1015)**
AGCTCTCTGTGAGGTGAACCGTTTCTAGAATCCCACCGTCTCGTCTGTTCTTCCGCCACCAGTTCTCAAGATAGCCCCTGTGGCTTCTGATGAA

1101 GTCACCACACCCTGGCTAATGAAGTAGATAAACAGAACAGTTTGGTTAACATTTAAGGTCAGAAACAGGAACCTTTCTAGAGGAGAAATCAAAAAAGC

1201 **PshAI (1277)**
AAAAGAAGTATAAGGGCAGCCCTCAACCAGTCAGAATACCGTGACCACCTGAGAGGCCGTGGCCAGCGGACACGGACGCATGTCAACTCTGGAGCAG

1301 ATATCTTCAGCGCAGCATCTGACCTGGGAGTACAGCCACATACCCTCATTCTAAACGGCAGATTGACTACTGCAGCCACACACAGTCTCCGGGCAATGT

1401 **SspI (1415)**
GGAGACATGTCTAATATTTAGTCAACATAAAGTCAAGGTGCCACAGTCTTCACTGTTGTGAGCACTTGGAGATGCTGCATTTGAAGATAGGAATTTGC

1501 CCTCAAGCATCTGGGGTTTGGGTACAGAACAGAGCTTCCCTGCCACCACCTGCTAATTTTATAAAATGTGCATTCAAAAAAAATCCTGCCTGTAAGAA

1601 **SwaI (1619)**
GGAATTAAGCTACCCATTTAAATATAACAGCTGCCTGTGCAATCTACTGCTGCTTTTATAGGAAACGCTTAAATAATTGAGATACTTAATTGGGTTAAA

1701 GAGATCCCTAGCACATAGATGTTCTATAAATAAAGAATGAGTAAATAATCTAGTAACCTTCTTTTCATGTCCTTCACTTAAAGAGATCGTTCTGTTTT

1801 GTTTGCACCAATAAGATCACTGTTAGAGGACTCCAGAGAGTTTATTTCAGGTGGGGTGGGGCTTCCCAAGGAAGTCCCTTTTCATTTGTTTCAGGTG

1901 TACTGCCACCTTTTTCCCTGGCTCTTCACTAAAAATGAAAAATTTGTTGATCTTTGCTGTAAGTAGGTAGGCATCTGGGCTTTGCTTTTCACTAGAG

2001 TCAAAGAAGTCAAGTTATCAGGCTGATCTTGCCTTGTATCTAGAATCAGAAAGTTAAGTAGCCAGGGACTACTCAAAGACAGCTGGAGGAGAAAGG

2101 **Tth111I (2145)**
GAGAGAGAAAAATGCTTATAAAGAGGTGGGCAAAAGAGCGGGACCTTGTCTCAAAAAAAAAAAAAAAAAAGAGGAAGTGGTAGGAGGTCTGAATTTCA

2201 CTGTGACCTGTTCTGTCAGGTGATTTTTGGTGGGGCGGGACATGAAAAAAAGTTAAATGTCCTTATAAGACAAAATCTTTTTCTTCTGGCTGAT

2301 XmnI (2375) ScaI (2401)
GATTTGTCATTCTAGTCACTTCTGCTTGTGACCACACACCCAGGCTTGACAAAGCTGTTCTGCAGATCAGAAAGAAGGGTCTCTGCTACACACCAG

2401 **NcoI (2406)** **NheI (2444)** **Acc65I**
TACTACCATGGGGGTTCTCATCATCATCATCATGGTATGGCTAGCATGACTGGTGGACAGCAAATGGGTCGGGATCTGTACGACGATGACGATAAG
32 ▶ Me tGI yGI ySer Hi sHi sHi sHi sHi sGI yMe tAl aSer Met Thr GI yGI yGI nGI nMe tGI yAr gAspLeuTyrAspAspAspAspLys

2501 **Bsu36I (2505)**
GTACCTAAGGATCAGCTTGGAGTTGATCCCGTCGTTTTACAACGTCGTGACTGGGAAAACCTGGCGTTACCCAACCTTAATCGCCTTGCAGCACATCCCC
32 ▶ Val ProLysAspGI nLeuGI yVal AspP roVal Val LeuGI nArgArgAspT rpGI uAsnProGI yVal Thr GI nLeuAsnArgLeuAl aAl aHi sP roP

2601 **FspI (2660)**
CTTTCGCCAGCTGGCGTAATAGCGAAGAGGCCCGACCGATCGCCCTTCCAACAGTTGCGCAGCCTGAATGGCGAATGGCGCTTTGCCTGGTTCCGGC
65 ▶ r oPheAl aSer T rpArgAsnSer GI uGI uAl aArgThrAspArgP roSer GI nGI nLeuArgSer LeuAsnGI yGI uT rpArgPheAl aT rpPheP roAl

2701 **Bsu36I (2742)**
ACCAGAAGCGGTGCCGAAAGCTGGCTGGAGTGCATCTTCCCTGAGCGGATACTGTCGTCGCTCCCTCAAACCTGGCAGATGCACGGTTACGATGCGCCC
98 ▶ aP roGI uAl aVal P roGI uSer T rpLeuGI uCvsAspLeuP roGI uAl aAspThr Val Val Val P roSer AsnT rpGI nMe tHi sGI vT vrAspAl aP ro

2801 ATCTACACCAACGTAACCTATCCATTACGGTCAATCCGCCGTTTGTCCACGGAGAATCCGACGGGTTGTTACTCGCTCACATTTAATGTTGATGAAA
132▶ I l e Tyr Thr Asn Val Thr Tyr Pro l e Thr Val Asn Pro P ro P he Val P ro Thr Gl y Asn P ro Thr Gl y Cys Tyr Ser Leu Thr P he Asn Val Asp Gl u S
2901 GCTGGCTACAGGAAGCCAGACGCGAATTATTTTTGATGGCGTAACTCGCGGTTTCATCTGTGGTGCACGGGCGCTGGGTCGGTACGCGCCAGGACAG
165▶ er T rp Leu Gl n Gl u Gl y Gl n Thr Arg l l e l e P he Arg Asp Val l e Asn Ser Al a P he Hi s Leu T rp Cys Asn Gl y A r g T rp Val Gl y T yr Gl y Gl n Asp Se
3001 TCGTTTGGCGTCTGAATTTGACCTGAGCGCATTTTTACGCGCCGAGAAAACCGCTCGCGGTGATGGTGCTGCGTTGGAGTGACGGCAGTTATCTGGAA
198▶ r Arg Leu P ro Ser Gl u P he Asp Leu Ser Al a P he Leu Arg Al a Gl y Gl u Asn Arg Leu Al a Val Met Val l e u Arg T rp Ser Asp Gl y Ser Tyr Leu Gl u

AatII (3141)

3101 GATCAGGATATGTGGCGGATGAGCGGCATTTCCGTGACGTCGTTGCTGCATAAACCGACTACACAAATCAGCGATTTCATGTTGCCACTCGCTTTA
232▶ Asp Gl n Asp Met T rp Arg Met Ser Gl y l e P he Arg Asp Val Ser Leu Leu Hi s Lys P ro Thr Thr Gl n l e Ser Asp P he Hi s Val Al a Thr Arg P he A
3201 ATGATGATTTACGCGCGCTGTACTGGAGGCTGAAGTTCAGATGTGCGCGAGTTGCGTGACTACCTACGGGTAACAGTTTCTTTATGGCAGGGTGAAC
265▶ sn Asp Asp P he Ser Arg Al a Val Leu Gl u Al a Gl u Val Gl n Met Cys Gl y Gl u Leu Arg Asp Tyr Leu Arg Val Thr Val Ser Leu T rp Gl n Gl y Gl u Th

ClaI (3342)

3301 GCAGGTGCGCAGCGCACCGCCCTTCGCGGTGAAATTCATGATGAGCGTGGTGGTTATGCCGATCGCGTCACACTACGTCTGAACGTCGAAAACCCG
298▶ r Gl n Val Al a Ser Gl y Thr Al a P ro P he Gl y Gl y Gl u l e l e l e Asp Gl u Arg Gl y Gl y Tyr Al a Asp Arg Val Thr Leu Arg Leu Asn Val Gl u Asn P ro
3401 AAAGTGTGGAGCGCCGAAATCCCGAATCTCTATCTGCGGTGGTTGAAGTGCACACCGCCGACGGCAGCGTATTGAAGCAGAAGCCTCGCATGTCGGTT
332▶ Lys Leu T rp Ser Al a Gl u l e P ro Asn Leu Tyr Arg Al a Val Val Gl u Leu Hi s Thr Al a Asp Gl y Thr Leu l e Gl u Al a Gl u Al a Cys Asp Val Gl y P
3501 TCCGCGAGGTGCGGATTGAAATGGTCTGCTGCTGCTGAACGGCAAGCCGTTGCTGATTCGAGGCGTTAACCGTCACGAGCATCATCTCTGCATGGTCA
365▶ he Arg Gl u Val l e Arg l l e Gl u Asn Gl y Leu Leu Leu Leu Asn Gl y Lys P ro Leu Leu l e Arg Gl y Val Asn Arg Hi s Gl u Hi s Hi s P ro Leu Hi s Gl y Gl
3601 GGTATGATGAGCAGACGATGGTGCAGGATATCTGCTGATGAAGCAGAACAACCTTTAACGCGTGCCTGTTGCGATTATCCGAACCATCCGCTGTGG
398▶ n Val Met Asp Gl u Gl n Thr Met Val Gl n Asp l e Leu Leu Met Lys Gl n Asn Asn P he Asn Al a Val l e Arg Cys Ser Hi s Tyr P ro Asn Hi s P ro Leu T rp

DraIII (3708)

SspI (3748)

3701 TACACGCTGTGCGACCGCTACGGCTGTATGTGGTGGATGAAGCCAATTTGAAACCCACGGCATGGTCCAATGAATCGTCTGACCGATGATCCGCGCT
432▶ Tyr Thr Leu Cys Asp Arg Tyr Gl y Leu Tyr Val Val Asp Gl u Al a Asn l e Gl u Thr Hi s Gl y Met Val P ro Met Asn Arg Leu Thr Asp Asp P ro Arg T

BsaBI (3844)

3801 GGCTACCGCGATGAGCGAACGCGTAACGCGAATGGTGACGCGCATGTAATCACCCGAGTGTGATCATCTGGTCGCTGGGGAATGAATCAGGCCACGG
465▶ r P ro Leu P ro Al a Me t Ser Gl u Arg Val Thr Arg Me t Val Gl n Arg Asp Arg Asn Hi s P ro Ser Val l e l e l e T rp Ser Leu Gl y Asn Gl u Ser Gl y Hi s Gl
3901 CGCTAATCAGCAGCGCTGTATCTGCTGGATCAAATCTGTGATCTTCCCGCGGTGAGTGAAGCGCGGAGCCGACCCAGCCACCGATTT
498▶ y Al a Asn Hi s Asp Al a Leu Tyr Arg T rp l e Lys Ser Val Asp P ro Ser Arg P ro Val Gl n Tyr Gl u Gl y Gl y Al a Asp Thr Thr Al a Thr Asp l e

BssHIII (4016)

4001 ATTTGCCCGATGTACGCGCGTGGATGAAGACAGCCCTTCCCGGCTGTGCCGAAATGGTCCATCAAAAAATGGCTTTCGCTACCTGGAGAGACGCGCC
532▶ I l e Cys P ro Me t Tyr Al a Arg Val Asp Gl u Asp Gl n P ro P he P ro Al a Val P ro Lys T rp Ser l l e Lys Lys T rp Leu Ser Leu P ro Gl y l e Th r Arg P
4101 CGCTGATCCTTTGCGAATACGCCACGCGATGGTAACAGCTTTGCGGCTTTCGCTAAATACTGGCAGGCGTTTCGTCAGTATCCCGGTTTACAGGGCGG
565▶ r o Leu l e Leu Cys Gl u Tyr Al a Hi s Al a Me t Gl y Asn Ser Leu Gl y Gl y P he Al a Lys Tyr T rp Gl n Al a P he Arg Gl n Tyr P ro Arg Leu Gl n Gl y Gl
4201 CTTCGCTGGGACTGGTGGATCAGTCGCTGATTAATATGATGAAAACGGCAACCCGTTGGTGGCTTACGGCGGTGATTTTGGCGATACGCCAACGAT
598▶ y P he Val T rp Asp T rp Val Asp Gl n Ser Leu l e Lys Tyr Asp Gl u Asn Gl y Asn P ro T rp Ser Al a Tyr Gl y Gl y Asp P he Gl y Asp Thr P ro Asn Asp

Eco47III (4353)

4301 CGCAGTTCTGTATGAACGGTCTGGTCTTTCGCGACCGCACGCGCATCCAGCGCTGACGGAAGCAAAAACACCAGCAGAGTTTTTCCAGTTCGCTTTAT
632▶ Arg Gl n P he Cys Me t Asn Gl y Leu Val P he Al a Asp Arg Thr P ro Hi s P ro Al a Leu Thr Gl u Al a Lys Hi s Gl n Gl n P he P he Gl n P he Arg Leu S

SacI (4458)

4401 CCGGCAAAACATCGAAGTGACCAGCGAATACCTGTTCCGTCATAGCGATAACGAGCTCCTGCCTGGATGGTGGCGCTGGATGGTAAGCCGCTGGCAAG
665▶ er Gl y Gl n Thr l l e Gl u Val Thr Ser Gl u Tyr Leu P he Arg Hi s Ser Asp Asn Gl u Leu Leu Hi s T rp Me t Val Al a Leu Asp Gl y Lys P ro Leu Al a Se
4501 CCGTGAAGTGCCTCTGGATGTCGCTCCACAAGGTAACAGTTGATTGAACTGCCTGAACTACCGCAGCCGAGAGCGCCGGCAACTCTGGCTCACAGTA
698▶ r Gl y Gl u Val P ro Leu Asp Val Al a P ro Gl n Gl y Lys Gl n Leu l e Gl u Leu P ro Gl u Leu P ro Gl n P ro Gl u Ser Al a Gl y Gl n Leu T rp Leu Thr Val
4601 CGGTAGTGCAACCGAACGCGACCGCATGGTCAGAAGCCGGGCACATACGCGCTGGCAGCAGTGGCGCTGGCGGAAAACCTCAGTGTGACGCTCCCGG
732▶ Arg Val Val Gl n P ro Asn Al a Thr Al a T rp Ser Gl u Al a Gl y Hi s l l e Ser Al a T rp Gl n Gl n T rp Arg Leu Al a Gl u Asn Leu Ser Val Thr Leu P ro A
4701 CCGCTCCACGCGCATCCCGCATCTGACCACCGCAATGGATTTTTGCATCGAGTGGGTAATAAGCGTTGGCAATTTAACCGCCAGTCAAGCTTTCT
765▶ l a Al a Ser Hi s Al a l e P ro Hi s Leu Thr Thr Ser Gl u Me t Asp P he Cys l l e Gl u Leu Gl y Asn Lys Arg T rp Gl n P he Asn Arg Gl n Ser Gl y P he Le
4801 TTCACAGATGTGGATTGGCGATAAAAAACAACCTGCTGACGCGCTGCGCATCAGTTACCCGTCACCGCTGGATAACGACATTGGCGTAAGTGAAGCG
798▶ u Ser Gl n Me t T rp l e Gl y Asp Lys Lys Gl n Leu Leu Thr P ro Leu Arg Asp Gl n P he Thr Arg Al a P ro Leu Asp Asn Asp l e Gl y Val l e Ser Gl u Al a
4901 ACCCGATTGACCCTAACGCTGGTGAACGCTGGAAGGCGGGCCATTACCAGGCCAAGCAGCGTTGTTGAGTGCACGGCAGATACACTTGCTG
832▶ Thr Arg l l e Asp P ro Asn Al a T rp Val Gl u Arg T rp Lys Al a Al a Gl y Hi s Tyr Gl n Al a Gl u Al a Leu Leu Gl n Cys Thr Al a Asp Thr Leu Al a A
5001 ATGCGGTGCTGATTACGACCGCTCAGCGTGGCAGCATCAGGGGAAAACCTTATTTATCAGCCGAAAACCTACCGGATTGATGGTAGTGGTCAAATGGC
865▶ sp Al a Val Leu l e Thr Thr Al a Hi s Al a T rp Gl n Hi s Gl n Gl y Lys Thr Leu P he l e Ser Arg Lys Thr Tyr Arg l l e Asp Gl y Ser Gl y Gl n Me t Al
5101 GATTACCGTTGATGTTGAAGTGGCGAGCGATACCCGATCCGCGCGGATTGGCCTGAACTGCCAGTGGCGCAGGTAGCAGAGCGGTAAACTGGCTC
898▶ al l e Thr Val Asp Val Gl u Val Al a Ser Asp Thr P ro Hi s P ro Al a Arg l l e Gl y Leu Asn Cys Gl n Leu Al a Gl n Val Al a Gl u Arg Val l e Asn T rp Leu

BsiWI (5288)

5201 GGATTAGGGCCGCAAGAAAATATCCCGACCGCTTACTGCCGCTGTTTACCAGCTGGGATCTGCCATTGTAGACATGTATACCCGTACGTCTTCC
932▶ Gl y Leu Gl y Pro Gl n Gl u Asn Tyr P ro Asp Arg Leu Thr Al a Al a Cys P he Asp Arg T rp Asp Leu P ro Leu Ser Asp Me t Tyr Thr P ro Tyr Val P he P
5301 CGAGCGAAAACGGTCTGCGTGGCGGACGCGCAATTTGAATTTAGCCACACAGTGGCGCGGCGACTTCCAGTTCAACATCAGCCGCTACAGTCAACA
965▶ r o Ser Gl u Asn Gl y Leu Arg Cys Gl y Thr Arg Gl u Leu Asn Tyr Gl y P ro Hi s Gl n T rp Arg Gl y Asp P he Gl n P he Asn l e Ser Arg Tyr Ser Gl n Gl

NdeI (5475)

5401 GCAACTGATGAAACAGCCATCGCATCTGCTGACGCGGAAGAACGACATGGCTGAATATCGACGTTTCCATATGGGGATTGGTGGCGACGACTCC
998▶ n Gl n Leu Me t Gl u Thr Ser Hi s Arg Hi s Leu Leu Hi s Al a Gl u Gl u Gl y Thr T rp Leu Asn l e Asp Gl y P he Hi s Me t Gl y l l e Gl y Gl y Asp Asp Ser

NheI (5597)

EcoRI (5591)

5501 TGGAGCCGTCAGTATCGCGGAATTACAGCTGAGCGCGGTCGCTACCATACCAGTTGGTCTGGTGTCAAAAAATAAATCTAGTCGAGAATTCGCTA
1032▶ T rp Ser P ro Ser Val Ser Al a Gl u Leu Gl n Leu Ser Al a Gl y Arg Tyr Hi s Tyr Gl n Leu Val T rp Cys Gl n Lys ●●●
5601 GCTCGACATGATAAGATACATTGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTA

MfeI (5771)

5701 TTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTAAGCTGCAATAAACAAGTTAACAACAACAATTGCATTCTTTATGTTTCAGGTTCA

SwaI (5862)

5801 GGGGGAGGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTAGATCCATTTAAATGTTAATTAAGTACGCGATGACCAAAATCCCTTAAC

5901 GTGAGTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTCTGCGCGTAATCTGCTGCTTGCAAACAAA
6001 AAAACCACCGCTACCAGCGGTGGTTTGTGGCCGATCAAGAGCTACCAACTCTTTTTCCGAAGGTAAGTGGCTTCAGCAGAGCGCAGATACCAATACT
6101 GTTCTTCTAGTGTAGCCGTAGTTAGGCCACCACCTCAAGAAGCTGTAGCACCCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCA
6201 GTGGCGATAAGTCGTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGGATAAGGCCAGCGGTGGGCTGAACGGGGGTTCTGTGCACACAGCCAG
6301 CTTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGAGCTATGAGAAAAGCCACGCTTCCCGAAGGGAGAAAAGCGGACAGGTATCCGGTA
6401 AGCGGCAGGGTCGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGAAACGCCTGGTATCTTTATAGTCCTGTCGGGTTTCGCCACCTCTGACTTGAGC
6501 GTCGATTTTTGTGATGCTCGTCAGGGGGCGGAGCCTATGAAAAACGCCAGCAACGCGGCCTTTTTACGGTTCCTGGCCTTTTGCTGGCCTTTTGCTCA

6601 CATGTTCTTAATTAATTTTCAAAGTAGTTGACAATTAATCATCGGCATAGTATATCGGCATAGTATAATACGACTCACTATAGGAGGGCCATCATG ^{AseI (6638)} ^{SfiI (6689)} **MseI (6700)**
6700 GCCAAGTTGACCAGTGCTGTCCAGTGCTCACAGCCAGGGATGTGGCTGGAGCTGTTGAGTTCTGGACTGACAGGTTGGGGTCTCCAGAGATTTTGTGG ¹Met
2 AlaLysLeuThr SerAl aVal P roVal LeuThr Al aArgAspValAl aGl yAl aVal Gl uPheTrpThrAspArgLeuGl yPheSer ArgAspPheVal G
6800 AGGATGACTTTGCAGGTGGTTCAGAGATGATGCACCCTGTTTCATCTCAGCAGTCCAGGACCAGGTGGTGCCTGACAACACCCTGGCTTGGGTGTGGGT
35 I uAspAspPheAl aGl yVal Val ArgAspAspVal Thr LeuPheI l eSer Al aVal Gl nAspGl nVal Val P roAspAsnThr LeuAl aTrpVal TrpVa
6900 GAGAGGACTGGATGAGCTGTATGCTGAGTGGAGTGAGGTGGTCTCCACCAACTTCAGGGATGCCAGTGGCCCTGCCATGACAGAGATTGGAGAGCAGCC
68 I ArgGl yLeuAspGl uLeuTyrAl aGl uTrpSer Gl uVal Val Ser ThrAsnPheArgAspAl aSer Gl yProAl aMe tThr Gl uI l eGl yGl uGl nP ro
7000 TGGGGAGAGAGTTTGCCTGAGAGACCCAGCAGGCAACTGTGTGCACCTTTGTGGCAGAGGAGCAGGACTGAGGATAAGAATTGAGTTTCAGAAAAGGGG ^{DraIII (7050)} ^{SfiI (7098)}
102 TrpGl yA rgGl uPheAl aLeuArgAspP roAl aGl yAsnCysVal Hi sPheValAl aGl uGl uGl nAsp•••
7100 GCCTGAGTGGCCCTTTTTCAACTTAATTAA