



2401 **SapI (2450)**
CTCTTTTCATGGAGTTTAAAGATATAGTGATTTTCCCAAGTTTGAAGCTAGCTCTTCATTTCTTTATGTTTTAAATGCACTGACCTCCACATTCCCTTT

2501 **SspI (2507)** **SwaI (2521)**
TTAGTAAAATATTCAGAAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTTATTAGGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCC

2601 **SacI (2782)**
CCCAGTTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTCTGGTGTACTTGAGGG
141 • N R T Y K L P

2701 **BstXI (2811)** **SacI (2782)**
GGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGTCAGAGATGAGCTCTGCACATGCC
133 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

2801 **StuI (2946)**
ACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTGCCGTTGCTCACAGCA
100 C 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

2901 **BbsI (3092)** **XmnI (3088)**
GACCCAATGGCAATGGCTTCAGCACAGACAGTGACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATCTCCCAGTCTTGGTCTGATGGCCG
66 S 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

3001 **AseI (3154)**
CCCCGACATGGTCTTGTTCCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGTCCAGATCTGCTGAGAGATGTTGAAGTCTTCAT
33 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

3101 **SacI (3211)**
GGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGTCAAACACAGCGTGGATGGCGTCTCCAGCTTATCTGACC
←

3201 **SpeI (3309)**
GTTCACTAAACGAGCTCTGCTTATATAGACCTCCACCCTACACGCCTACCGCCATTTCGCTCAATGGGGCGGAGTTGTTACGACATTTTGAAAGTCC

3301 **SnaBI (3437)**
CGTTGATTTACTAGTCAAACAAACTCCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCAAACCGCTATCCACGCCATTGATGTACT
←

3401 **NdeI (3542)**
GCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAAAGTCCATAAGGTCATGTACTGGGCATAATGCCAGGC

3501 **SdaI (3720)** **PacI (3728)** **BspLU11I (3738)**
GGGCCATTTACCGTCATTGACGTCAATAGGGGGCTACTTGGCATATGATACTTGTACTGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCAT
3601 TGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTATTATTGACGTCAATGGGCGGGGTCTTGGGCGGTCAGCCAGGCGGGCCAT

3701 **EagI (4488)**
TTACCGTAAGTTATGTAACGCCTGCAGGTTAAATAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCGTAAAAAGGCCGCTTCTGGCGTTT
←

3801 **PacI (4468)** **SwaI (4477)** **NotI (4487)**
TTCCATAGGCTCCGCCCCCTGACGAGCATCAAAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGACTATAAAGATACCAGGCGTTTCCCC

3901 **EagI (4488)**
CTGGAAGCTCCCTCGTGCCTCTCCTGTTCCGACCCTGCCGTTACCGGATACCTGTCCGCTTTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTC

4001 **ApaLI (4052)**
ACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTCCAGCCGACCGCTGCGCCTTATCCGGTAAC

4101 **EagI (4488)**
TATCGTCTTGTAGTCAACCCGGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGAGCGAGGTATGTAGGCGGTGTACAG

4201 **EagI (4488)**
AGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAG

4301 **EagI (4488)**
CTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGTTTTTTTTGTTTGCAAGCAGCAGATTACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTG

4401 **EagI (4488)**
ATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACCTCACGTTAAGGGATTTTGGTTCATGGCTAGTTAATTAACATTTAAATCAGCGGCCCAATAA
PacI (4468) SwaI (4477) NotI (4487)

4501 **EagI (4488)**
AATATCTTTATTTTACATCTGTGTGTTGTTTTTTGTGTGAATCGTAACATAACGCTCTCCATCAAACAAAACGAAACAAAACAACTAGCA

4601 **EagI (4488)**
AAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTATCGAA