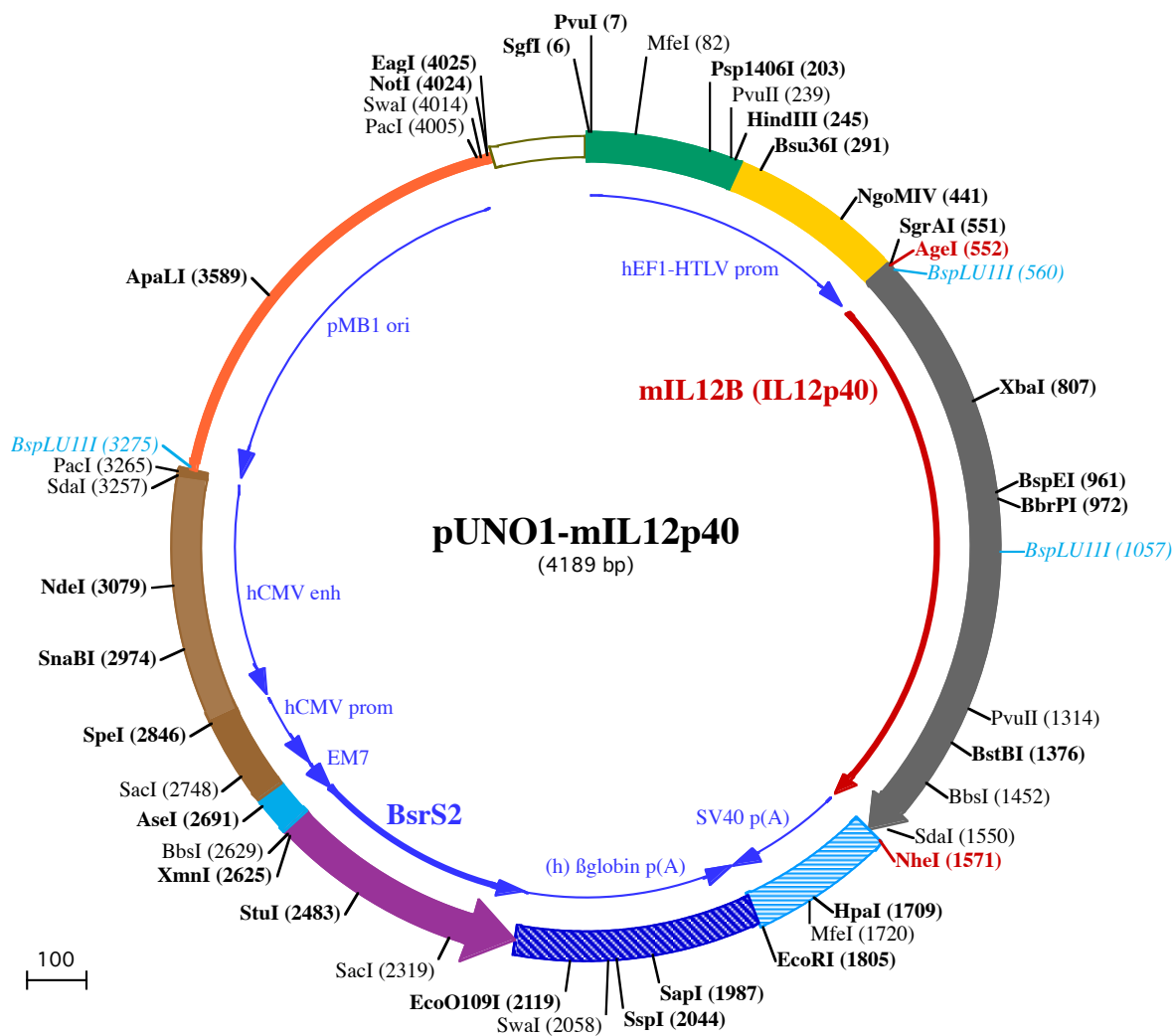




PvuI (7)
SgfI (6) MfeI (82)

1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGCAGAGCGCACATCGCCACAGTCCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA

101 GAGAAGGTGGCGCGGGGTAAACTGGGAAAGTATGTCGTGTACTGGCTCCGCCTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC

HindIII (245)

Psp1406I (203) PvuII (239) Bsu36I (291)

201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCGCGCCCTACCTGAGGCC

301 GCCATCCACGCCGGTTGAGTCGCGTTTCTGCCGCTCCCGCCTGTGGTGCTCCTGAACTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)

401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTGCCTGACCCTGCTTGCTCAACTCTACGCTTTGTTTCGTTT

AgeI (552)
SgrAI (551) BspLU11I (560)

501 TCTGTTCTGCGCGGTTACAGATCCAAGCTGTGACCGCGCCTACCTGAGATCACCGGTGAACATGTGTCCTCAGAAGCTAACCATCTCCTGGTTTGCCAT

601 CGTTTTGCTGGTGTCTCCACTCATGGCCATGTGGGAGCTGGAGAAAGACGTTTATGTTGTAGAGGTGGACTGGACTCCCGATGCCCTGGAGAAACAGTG

701 AACCTCACCTGTGACACGCTGAAGAAGATGACATCACCTGGACCTCAGACCAGAGACATGGAGTCATAGGCTCTGAAAAGACCTGACCATCACTGTCA

801 AAGAGTTTCTAGATGCTGGCCAGTACACCTGCCACAAAGGAGCGAGACTCTGAGCCACTCACATCTGCTGCTCCACAAGAAGGAAAAATGGAATTTGGTC

901 CACTGAAATTTTAAAAATTTCAAAAACAAGACTTTCTGAAGTGTGAAGCACAAATTACTCCGGACGGTTCACGTGCTCATGGCTGGTGCAGAAAGAAC

1001 ATGGACTTGAAGTTCAACATCAAGAGCAGTAGCAGTCCCCCGACTCTCGGGCAGTGACATGTGGAATGGCGTCTCTGTCTGCAGAGAAGGTCACTGG

1101 ACCAAAGGGACTATGAGAAGTATTCAGTGTCTGCCAGGAGGTGTACCTGCCAACTGCCGAGGAGACCCTGCCATTGAACTGGCGTTGGAAGCAGC

1201 GCAGCAGAATAAATATGAGAATACAGCACCAGCTTCTTCATCAGGGACATCATCAAAACCAGACCCGCCAAGAACTGCAGATGAAGCCTTTGAAGAAC

1301 TCACAGGTGGAGGTGAGCTGGGAGTACCTGACTCCTGGAGCACTCCCATTCCTACTTCTCCTCAAGTTCTTTGTTCCGAATCCAGCGCAAGAAAGAAA

1401 AGATGAAGGAGACAGAGGAGGGGTGTAACCAGAAAGGTGCGTTCCTCGTAGAGAAGACATCTACCGAAGTCCAATGCAAAGGCGGGAATGTCTGCGTGCA

1501 AGCTCAGGATCGCTATTACAATTCCTCATGCAGCAAGTGGGATGTGTTCCCTCGAGGTCGATCCATAAGCTAGCTGGCCAGACATGATAAGATACAT

1601 TGATGAGTTTGGACAAACCACAACCTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGC

HpaI (1709) MfeI (1720)

1701 AATAAACAAAGTTAAACAACAATTGCATTATTTTATGTTTCAGGTTTCAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAACCTCTACAAATGTG

EcoRI (1805)

1801 GTATGGAATTCATAAATACAGCATAGCAAACTTTAACCTCCAAATCAAGCCTCTACTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGG

SapI (1987)

1901 GCTGTTGCCAATGTGCATTAGCTGTTTGCAGCCTCACCTTCTTTCATGGAGTTTAAGATATAGTGTATTTTCCAAGGTTTGAAGTAGCTCTTCATTCT

SspI (2044) SmaI (2058)

2001 TTATGTTTTAAATGCACTGACCTCCACATTCCCTTTTGTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAATAAATGTTTTTTATTA

EcoO109I (2119)

2101 GGCAGAATCCAGATGCTCAAGGCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAAGAACCTTTAATAGAAATTGGACAGCAAGAAA

2201 GCGAGCTTCTAGCTTTAGTTCCTGGTGTACTTGAGGGGGATGAGTTTCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCAATGAGCACAAGCAGT

2301 CAGGAGCATAGTCAGAGATGAGCTCTGCACATGCCACAGGGGTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCTGACAGCCAC

SacI (2319)

112 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 (2483)

2401 AATGGTGTCAAAGTCCTTCTGCCGTTGCTCACAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTGACCTGCCAATGTAGGCCTCAATGTGGACA
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
2501 GCAGAGATGATCTCCCCAGTCTTGGTCCTGATGGCGCCCCGACATGGTGCTTGTGCTCATAGAGCATGGTGATCTTCTCAGTGGCGACCTCCACCA
45 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 L

BbsI (2629)

XmnI (2625)

AseI (2691)

2601 GCTCCAGATCCTGCTGAGAGATGTTGAAGGTCTTCATGGTGGCCCTCTATAGTGAGTCGTATTATACTATGCCGATATACTATGCCGATGATTAATTGT
12 E L D Q Q S I N F T K M

SacI (2748)

2701 CAAACAGCGTGGATGGCGTCTCCAGCTTATCTGACGGTTCATAAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCCATTTGCGT

SpeI (2846)

2801 CAATGGGCGGAGTTGTTACGACATTTTGAAAGTCCCGTTGATTTACTAGTCAAAACAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCC

SnaBI (2974)

2901 CGTGAGTCAAACCGCTATCCACGCCATTGATGTACTGCCAAAACCGCATCATCATGGTAATAGCGATGACTAATACGTAGATGTACTGCCAAGTAGGAA

NdeI (3079)

3001 AGTCCATAAGGTCATGTACTGGGCATAATGCCAGGCGGGCCATTTACCGTCATTGACGTCAATAGGGGGCGTACTTGGCATATGATACACTTGATGTAC

3101 TGCCAAGTGGGCAGTTTACCGTAAATACTCCACCCATTGACGTCAATGAAAGTCCCTATTGGCGTTACTATGGGAACATACGTCATTATTGACGTCAAT

SdaI (3257) PacI (3265) BspLU11I (3275)

3201 GGGCGGGGTCGTTGGGCGTCCAGGCGGGCCATTTACCGTAAGTTATGTAACGCTGCAGGTTAATTAAAGAACATGTGAGCAAAAGGCCAGCAAAA

3301 GGCCAGGAACCGTAAAAAGGCCGCTTGTGGCGTTTTTCCATAGGCTCCGCCCCCTGACGAGCATCACAAAATCGACGCTCAAGTCAGAGGTGGCGA

3401 AACCCGACAGGACTATAAAGATACCAGCGCTTCCCCCTGGAAGCTCCCTCGTGCGCTCTCCTGTTCCGACCTGCCGCTTACCGGATACCTGTCCGCT

ApaLI (3589)

3501 TTCTCCCTTCGGGAAGCGTGGCGCTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTGCTTCCGCTCAAGCTGGGCTGTGTGCACGAACC

3601 CCCGTTTCAGCCGACCGCTGCGCTTATCCGGTAACATCGTCTTGAGTCCAACCCGTAAGACACGACTTATCGCCACTGGCAGCAGCCACTGGTAAC

3701 AGGATTAGCAGAGCGAGGTATGTAGGCGGTGCTACAGAGTTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTC

3801 TGCTGAAGCCAGTTACCTTCGAAAAAGAGTTGGTAGCTCTTGATCCGGCAAACAAACCACCGCTGGTAGCGGTGGTTTTTTGTTTGCAAGCAGCAGAT

3901 TACGCGCAGAAAAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGTCTGACGCTCAGTGGAAACGAAAACTCACGTTAAGGGATTTTGGTCATG

EagI (4025)

PacI (4005) SmaI (4014) NotI (4024)

4001 GCTAGTTAATTAACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGC

4101 TCTCCATCAAAACAAACGAAACAAACAACTAGCAAAATAGGCTGTCCCAGTGCAAGTGCAGGTGCCAGAACATTTCTCTATCGAA