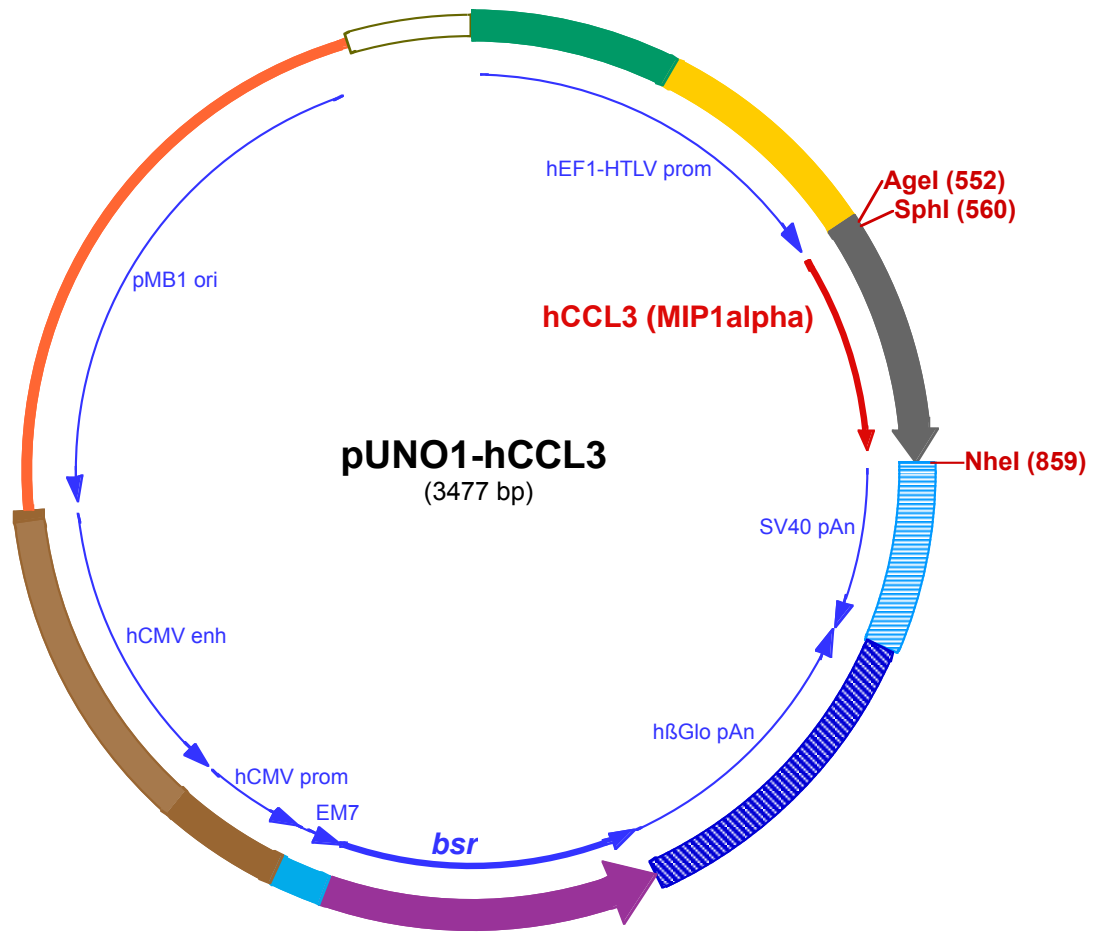


75



1 GGATCTGCGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATGCCACAGTCCCGAGAAGTTGGGGGAGGGGTCGGCAATTGAACGGGTGCCTA  
101 GAGAAGGTGGCGCGGGGTAAACTGGGAAAGTGATGTCGTGACTGGCTCCGCCTTTTTCCCGAGGGTGGGGGAGAACCGTATATAAGTGCAGTAGTCGCC  
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTCACGCGCCCGCCCTACCTGAGGCC  
301 GCCATCCACGCCGTTGAGTCGCGTTCTGCCGCTCCCGCTGTGGTGCCTCCTGAAGCTGCGTCCGCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC  
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCCGGCTCTCCACGCTTTGCCTGACCTGCTTGTCTCAACTCTACGCTTTGTTTCGTTT

**SphI (560)**  
**AgeI (552)**

501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCCTACCTGAGATCACCGGTCAGCATGCAGGTCTCCACTGCTGCCCTGTGTCCTCCTCTG  
1 M Q V S T A A L A V L L C  
601 CACCATGGCTCTCTGCAACCAAGTCTCTGCATCACTTGTGCTGACACGCCGACCGCTGCTGCTTACAGTACACCTCCCGCAGATTCCACAGAATTC  
13 T M A L C N Q F S A S L A A D T P T A C C F S Y T S R Q I P Q N F  
701 ATAGCTGACTACTTTGAGACGAGCAGCCAGTGTCCAAGCCTGGTGCATCTTCTAACCAGCGAAGCCGGCAGGTCTGTGCTGACCCAGTGAGGAGT  
47 I A D Y F E T S S Q C S K P G V I F L T K R S R Q V C A D P S E E

**NheI (859)**

801 GGGTCCAGAAATATGTCAGCGACCTGGAGCTGAGTGCCTGAGGGTCCAGAAGCTTCGAGCTAGTGGCCAGACATGATAAGATACATTGATGAGTTGG  
80 W V Q K Y V S D L E L S A •  
901 ACAAACCAACTAGAATGCAGTGAAAAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGTGAATAAACAAGTT

1001 AACAAACA AATTGCATTCA TTTTATGTTTCAGGTT CAGGGGAGGTGTGGGAGGTTTTTAAAGCAAGTAAAACCTCTACA AATGTGGTATGGAATTCT  
1101 AAAATACAGCATAGCAAACTTAACTCCAATCAAGCCTCTACTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTGCCAAT  
1201 GTGCATTAGCTGTTTGCAGCCTCACCTTCTTCATGGAGTTAAGATATAGTGTATTTCCCAAGGTTTGAAGTACTGCTTTCATTTCTTTATGTTTAAA  
1301 TGCACTGACCTCCACATTCCCTTTTAGTAAATATTCAGAAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAG  
1401 ATGCTCAAGGCCCTTCATAATATCCCCAGTTAGTAGTTGGACTTAGGGAACAAAGGAACCTTAAATAGAAATTGGACAGCAAGAAAGCGAGCTTCTAG  
1501 CTTAGTTCCTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCATCTCAATGAGCACAAAGCAGTCAGGAGCATAGT  
141 • N R T Y K L P 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  
1601 CAGAGTAGCTCTCTGCACATGCCACAGGGGCTGACCACCCTGATGGATCTGTCCACCTCATCAGAGTAGGGTGCCTGACAGCCACAATGGTGTCAA  
108 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 I T D F  
1701 GTCCTTCTGCCGTTGCTCAGCAGACCCAATGGCAATGGCTTCAGCACAGACAGTACCCTGCCAATGTAGGCCTCAATGTGGACAGCAGAGATGATC  
75 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 A S I I  
1801 TCCCCAGTCTGGTCTGATGGCCGCCCGACATGGTCTTGTGTCCTCATAGAGCATGGTATCTTCTCAGTGGCGACCTCCACCAGTCCAGATCCT  
41 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 E L D Q  
1901 GCTGAGAGATGTTGAAGTCTTCATGCTGGCCCTCTATAGTGAAGTCTGATTACTATGCGGATATACTATGCCGATGATTAATTGTCAAACAGCGTG  
8 Q S I N F T K M  
2001 GATGGCGTCTCCAGCTTATCTGACGGTTCACATAACGAGCTCTGCTTATATAGACCTCCACCGTACACGCCTACCGCCATTTGCGTCAATGGGGCGGA  
2101 GTTGTACGACATTTTGGAAAGTCCCGTTGATTTACTAGTCAAACAAACTCCATTGACGTCAATGGGGTGGAGACTTGAAATCCCGTGAGTCAAAC  
2201 CGCTATCCAGCCATTGATGACTGCCAAAACCGCATCATGGAATAGCGATGACTAATACGTAGATGACTGCAAGTAGGAAAGTCCATAAGG  
2301 TCATGACTGGGCATAATGCCAGGCGGGCATTACCCTGATTGACGTCAATAGGGGGGCTACTTGGCATATGATACACTTGTGACTGCAAGTGGGC  
2401 AGTTTACCGTAAATACTCCACCCATTGACGTCAATGGAAAGTCCCTATTGGCGTACTATGGGAACATACGTCAATTATTGACGTCAATGGCGGGGGTCCG  
2501 TTGGGCGTCCAGCAGCGGGCCATTACCCTAAGTTATGTAACCGCTGCAGGTTAATAAGAACATGTGAGCAAAGGCCAGCAAAGGCCAGGAACCG  
2601 TAAAAGCGCCGTTGCTGGCGTTTTCCATAGGCTCCGCCCCCTGACGAGCATCAGAAAATCGACGCTCAAGTCAGAGGTGGCGAAACCCGACAGGA  
2701 CTATAAGATACCAGGCGTTTTCCCTGGAAGCTCCCTCGTGCCTCTCTGTTCCGACCCTGCCGCTTACCGGATACCTGTCCGCTTTCTCCCTCCG  
2801 GAAGCGTGGCGCTTTCTCATAGCTCACGCTGTAGGTATCTCAGTTCGGTGTAGGTCGTTCCGCTCAAGCTGGGCTGTGTGCACGAACCCCGTTAGCC  
2901 CGACCGCTGCGCCTTATCCGGTAACTATCGTCTTGTGAGTCAACCCGGTAAAGACAGACTTATCGCACTGGCAGCAGCCACTGGTAACAGGATTAGCAGA  
3001 GCGAGGTATGTAGCGGTGCTACAGAGTCTTGAAGTGGTGGCCTAACTACGGCTACACTAGAAGAACAGTATTTGGTATCTGCGCTCTGCTGAAGCCAG  
3101 TTACCTTCGAAAAAGAGTTGGTAGCTCTTGTATCCGGCAAACAAACCCTGGTGGTGGTGGTTTTTTTGTGCAAGCAGCAGATTACGCGCAGAAA

3201 AAAAGGATCTCAAGAAGATCCTTTGATCTTTTCTACGGGGTCTGACGCTCAGTGGAACGAAACTCACGTTAAGGGATTTTGGTCATGGCTAGTTAATTA

---

3301 ACATTTAAATCAGCGGCCGCAATAAAATATCTTTATTTTCATTACATCTGTGTGTTGGTTTTTGTGTGAATCGTAACTAACATACGCTCTCCATCAAAA

3401 CAAAACGAAACAAAACAACTAGCAAAATAGGCTGTCCCAGTGCAAGTGCAAGTGCCAGAACATTTCTCTATCGAA