

pE-SUM0pro3 Kan

>pE-SUM0pro3 Kan (5610 bp)

AGATCTCGATCCCAGCAAATTAATACGACTACTATAGGGGAATTGTGAGCGGATAACAATTCCCCTCTAGAAATAATTT
TGTTTAACTTTAAGAAGGAGATATACCATGGGTATCACCATCATCATCACGGGTCCCTGCAGGAGGAGAAGCCCAAGGA
GGGTGTGAAGACAGAGAATGACCACATCAACCTGAAGGTGGCCGGCAGGACGGCTCCGTGGTGCAGTTCAAGATCAAGA
GGCACACGCCGCTGAGCAAGCTGATGAAGGCCTACTGCGAGAGGCAGGGCTTGTCAATGAGGCAGATCAGATTCAGGTTT
GACGGGCAGCCAATCAATGAACTGACACTCCAGCACAGCTGGAGATGGAGGACGAGGACACCATCGACGTGTTCCAGCA
GCAGACGGGAGGTTGAGACCACTAGTGGTACCGGTCTCACTAGAGGATCCGAATTCGAGCTCCGTGACAAGCTTGCGGC
CGCACTCGAGCACCACCACCACCACCCTGAGATCCGGCTGCTAACAAAGCCCGAAAGGAAGCTGAGTTGGCTGCTGCCA
CCGCTGAGCAATAACTAGCATAACCCCTTGGGGCCTCTAAACGGGTCTTGAGGGGTTTTTGTGAAAGGAGGAACTATA
TCCGGATTGGCGAATGGGACGCGCCCTGTAGCGGCGCATTAAAGCGCGCGGGTGTGGTGGTTACGCGCAGCGTGACCGCT
ACACTTGCCAGCGCCCTAGCGCCCGCTCCTTTGCTTTCTTCCCTTCTTTCTCGCCACGTTTCGCCGGCTTTCCCCGTCA
AGCTCTAAATCGGGGGCTCCCTTTAGGGTTCGATTTAGTGCTTTACGGCACCTCGACCCCAAAAACTTGATTAGGGTG
ATGGTTCACGTAGTGGGCCATCGCCCTGATAGACGGTTTTTCGCCCTTGACGTTGGAGTCCACGTTCTTTAATAGTGGA
CTCTTGTTCCAACTGGAACAACACTCAACCCTATCTCGGTCTATTCTTTTATTTATAAGGGATTTTGCCGATTTCCGC
CTATTGGTTAAAAATGAGCTGATTTAACAAAAATTTAACGCGAATTTTAAACAAATATTAACGTTTACAATTTAGGTG
GCACTTTTGGGGAAATGTGCGCGGAACCCCTATTTGTTATTTTTCTAAATACATTCAAATATGTATCCGCTCATGAAT
TAATCTTAGAAAACTCATCGAGCATCAAATGAACTGCAATTTATTCATATCAGGATTATCAATACCATATTTTTGAA
AAAGCCGTTTCTGTAATGAAGGAGAAAACCTACCGAGGCAGTTCATAGGATGGCAAGATCCTGGTATCGGTCTGCGATT
CCGACTCGTCCAACATCAATAACAACCTATTAATTTCCCTCGTCAAAAATAAGTTATCAAGTGAAGAAATCACCATGAGT
GACGACTGAATCCGGTGAGAATGGCAAAAGTTTATGCATTTCTTTCCAGACTTGTTCAACAGGCCAGCCATTACGCTCGT
CATCAAAATCACTCGCATCAACCAAACCGTTATTCATTCTGATGTCGCTGAGCGGAGACGAAATACGCGATCGCTGTTA
AAAGGACAATTACAAACAGGAATCGAATGCAACCGGCGCAGGAACACTGCCAGCGCATCAACAATATTTTACCTGAATC
AGGATATTCTTCTAATACCTGGAATGCTGTTTTCCCGGGGATCGCAGTGGTGAATACCATGCATCATCAGGAGTACGGA
TAAATGCTTGATGGTGGGAAGAGGCATAAATCCGTGAGCCAGTTTAGTCTGACCATCTCATCTGTAACATCATTGGCA
ACGCTACCTTTGCCATGTTTCAAGAAACAACCTCTGGCGCATCGGGCTTCCCATACAATCGATAGATTGTGCGACCTGATTG
CCCAGATTATCGCGAGCCATTTATACCCATATAAATCAGCATCCATGTTGGAATTTAATCGCGGCCTAGAGCAAGACG
TTTCCCGTTGAATATGGCTCATAACACCCCTTGTATTACTGTTTATGTAAGCAGACAGTTTTATTGTTTATGACCAAAAT
CCCTTAACGTGAGTTTTCGTTCCACTGAGCGTCAAGCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTTC
TGCGCGTAATCTGCTGCTTGCAAACAAAAAACCACCGCTACCAGCGGTGGTTTGTGTTGCCGGATCAAGAGCTACCACT
CTTTTTCCGAAGGTAACCTGGCTTCAAGCAGAGCGCAGATACCAAAACTGTCTTCTAGTGTAGCCGTAGTTAGGCCACCA
CTTCAAGAACTCTGTAGCACCGCCTACATACCTCGCTCTGCTAATCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGT
CGTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTGGGGCTGAACGGGGGGTTCTGTGCACA
CAGCCAGCTTGGAGCGAACGACCTACACCGAAGTGAATACCTACAGCGTGAAGTATGAGAAAGCGCCACGCTTCCCGA
AGGGAGAAAGGCGGACAGGTATCCGGTAAGCGGCAGGGTCGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGGAAACG
CCTGGTATCTTTATAGTCTGTGCGGTTTTCGCCACCTCTGACTTGAGCGTGCATTTTTGTGATGCTCGTCAAGGGGGCGG
AGCCTATGGAAAAACGCCAGCAACGCGGCCTTTTTACGGTTCCTGGCCTTTTGTGCGCCTTTTGTGCTCACATGTTCTTCC
TGCGTTATCCCCTGATTCTGTGGATAACCGTATTACCGCCTTTGAGTGAAGTGTATACCGCTCGCCGCAGCCGAACGACCG
AGCGCAGCGAGTCAAGTGAAGCGAGGAAGCGGAAGAGCGCCTGATGCGGTATTTTTCTCTTACGCATCTGTGCGGTATTTCA
CACCGCATATATGGTGCCTCTCAGTACAATCTGCTCTGATGCCGCATAGTTAAGCCAGTATACACTCCGCTATCGCTAC
GTGACTGGGTATGGCTGCGCCCCGACACCCGCCAACACCCGCTGACGCGCCCTGACGGGCTTGTCTGCTCCCGGCATCC
GTTACAGACAAGCTGTGACCGTCTCCGGGAGCTGCATGTGTGAGAGTTTTACCGTCAACCGAAACGCGCGAGGCA
GCTGCGGTAAAGCTCATCAGCGTGGTCTGAAGCGATTACAGATGTCTGCCTGTTTATCCGCGTCCAGCTCGTTGAGTT
TCTCCAGAAGCGTTAATGTCTGGCTTCTGATAAAGCGGGCCATGTTAAGGGCGTTTTTTTCTGTTTGGTCACTGATGCC
TCCGTGTAAGGGGGATTTCTGTTTATGGGGGTAATGATACCGATGAAACGAGAGAGGATGCTCACGATACGGGTTACTGA
TGATGAACATGCCCGTTACTGGAACGTTGTGAGGGTAAACAACCTGGCGGTATGGATGCGGGCGGGACCAGAGAAAAATCA
CTCAGGGTCAATGCCAGCGCTTCTGTTAATACAGATGTAGGTGTTCCACAGGGTAGCCAGCAGCATCTGCGATGCAGATC
CGGAACATAATGGTGCAGGGCGCTGACTTCCGCGTTTTCCAGACTTTACGAAACACGGAAACCGAAGACCATTATGTTGT
TGCTCAGGTCGAGACGTTTTGCAGCAGCAGTCTGTTTACGTTTCTGCTCGCGTATCGGTGATTCACTTCTGCTAACCAAGTAA

pE-SUMOpro3 Kan

GGCAACCCCGCCAGCCTAGCCGGGTCCTCAACGACAGGAGCACGATCATGCGCACCCGTGGGGCCGCCATGCCGGCGATA
ATGGCCTGCTTCTCGCCGAAACGTTTGGTGGCGGGACCAGTGACGAAGGCTTGAGCGAGGGCGTGCAAGATTCCGAATAC
CGCAAGCGACAGGCCGATCATCGTCGCGCTCCAGCGAAAGCGGTCTCGCCGAAAATGACCCAGAGCGCTGCCGGCACCT
GTCCTACGAGTTGCATGATAAAGAAGACAGTCATAAGTGCGGCGACGATAGTCATGCCCCGCGCCACCGGAAGGAGCTG
ACTGGGTTGAAGGCTCTCAAGGGCATCGGTGAGATCCCGGTGCCTAATGAGTGAGCTAACTTACATTAATTGCGTTGCG
CTCACTGCCCGCTTTCCAGTCGGGAAACCTGTCGTGCCAGCTGCATTAATGAATCGGCCAACGCGCGGGGAGAGGCGGTT
TGCGTATTGGGCGCCAGGGTGGTTTTTCTTTTACCAGTGAGACGGGCAACAGCTGATTGCCCTTACCAGCCTGGCCCTG
AGAGAGTTGCAGCAAGCGGTCCACGCTGGTTTGGCCAGCAGGCGAAAATCCTGTTTGATGGTGGTTAACGGCGGGATAT
AACATGAGCTGTCTTCGGTATCGTCGTATCCCACTACCGAGATATCCGCACCAACGCGCAGCCCGGACTCGGTAATGGCG
CGCATTGCGCCAGCGCCATCTGATCGTTGGCAACCAGCATCGCAGTGGGAACGATGCCCTCATTACGATTTGCATGGT
TTGTTGAAAACCGGACATGGCACTCCAGTCGCCTTCCCGTTCGGCTATCGGCTGAATTTGATTGCGAGTGAGATATTTAT
GCCAGCCAGCCAGACGCAGACGCGCCGAGACAGAACTTAATGGGCCCGCTAACAGCGCGATTTGCTGGTGACCCAATGCG
ACCAGATGCTCCACGCCCAGTCGCGTACCGTCTTCATGGGAGAAAATAATACTGTTGATGGGTGCTGGTCAGAGACATC
AAGAAATAACGCCGGAACATTAGTGACGGCAGCTTCCACAGCAATGGCATCCTGGTCATCCAGCGGATAGTTAATGATCA
GCCCACTGACGCGTTGCGCGAGAAGATTGTGCACCGCCGCTTTACAGGCTTCGACGCGCTTCGTTCTACCATCGACACC
ACCACGCTGGCACCCAGTTGATCGGCGCGAGATTTAATCGCCGCGACAATTTGCGACGGCGCGTGACGGCCAGACTGGA
GGTGGCAACGCCAATCAGCAACGACTGTTTGGCCGCCAGTTGTTGTGCCACGCGGTTGGGAATGTAATTCAGCTCCGCCA
TCGCCGCTTCCACTTTTTCCCGCTTTTTCGAGAAACGTGGCTGGCCTGGTTACCACGCGGGAAACGGTCTGATAAGAG
ACACCGGCATACTCTGCGACATCGTATAACGTTACTGGTTTTACATTACCACCCTGAATTGACTCTCTTCCGGGCGCTA
TCATGCCATAACCGCGAAAGTTTTGCGCCATTCGATGGTGTCCGGGATCTCGACGCTCTCCCTTATGCGACTCCTGCATT
AGGAAGCAGCCAGTAGTAGGTTGAGGCCGTTGAGCACCGCCGCCGCAAGGAATGGTGCATGCAAGGAGATGGCGCCCAA
CAGTCCCCCGGCCACGGGGCTGCCACCATACCCACGCCGAAACAAGCGCTCATGAGCCCGAAGTGGCGAGCCCGATCTT
CCCCATCGGTGATGTCGGCGATATAGGCGCCAGCAACCGCACCTGTGGCGCCGGTATGCCGGCCACGATGCGTCCGGCG
TAGAGGATCG