

pE-SUMOstar Amp

>pE-SUMOstar Amp (5761 bp)

AGATCTCGATCCCGCAAATTAATACGACTACTATAGGGGAATTGTGAGCGGATAACAATTCCCCTCTAGAAATAATTT  
TGTTTAACTTTAAGAAGGAGATATACCATGGGTATCACCATCATCATCACGGGTCCTGCAGGACTCAGAAGTCAATCA  
AGAAGCTAAGCCAGAGGTCAAGCCAGAAGTCAAGCCTGAGACTCACATCAATTTAAAGGTGTCCGATGGATCTTCAGAGA  
TCTTCTTCAAGATCAAAAAGACCACTCCTTTAAGAAGGCTGATGGAAGCGTTTCGCTAAAAGACAGGGTAAGGAAATGGAC  
TCCTTAAACGTTCTTGTACGACGGTATTGAAATTCAGCTGATCAGACCCCTGAAGATTTGGACATGGAGGATAACGATAT  
TATTGAGGCTCACCGCGAACAGATTGGAGGTTGAGACCACTAGTGGTACCGGTCTCACTAGAGGATCCGAATTCGAGCTC  
CGTCGACAAGCTTGCGGCCGCACTCGAGCACCACCACCACCACCCTGAGATCCGGCTGCTAACAAAGCCCCGAAAGGAAG  
CTGAGTTGGCTGCTGCCACCGCTGAGCAATAACTAGCATAACCCCTTGGGGCCTCTAAACGGGTCTTGAGGGGTTTTTGTG  
CTGAAAGGAGGAATAATCCGGATTGGCGAATGGGACGCGCCCTGTAGCGGCGCATTAAAGCGCGGCGGGTGTGGTGGTT  
ACGCGCAGCGTGACCGCTACACTTGCCAGCGCCCTAGCGCCCGCTCCTTTCGCTTCTTCCCTTCTTCTCGCCACGTT  
CGCCGGCTTTCCTCGTCAAGCTCTAAATCGGGGGCTCCTTTAGGGTTCGATTTAGTGCTTACGGCACCTCGACCCCA  
AAAACTTGATTAGGGTGTGGTTCACGTAGTGGGCCATCGCCCTGATAGACGGTTTTTCGCCCTTGACGTTGGAGTCC  
ACGTTCTTTAATAGTGGACTCTTGTCCAACTGGAACAACACTCAACCCTATCTCGGTCTATTCTTTTGATTTATAAGG  
GATTTTGCCGATTTGCGCCTATTGGTTAAAAATGAGCTGATTTAACAAAAATTTAACGCGAATTTTAACAAAATATTA  
CGTTTACAATTTAGGTGGCACTTTTCGGGAAATGTGCGCGGAACCCCTATTTGTTATTTTTCTAAATACATTCAAT  
ATGTATCCGCTCATGAGACAATAACCCTGATAAATGCTTCAATAATATTGAAAAAGGAAGAGTATGAGTATTCAACATTT  
CCGTGTCGCCCTTATTCCCTTTTTTTCGGCATTTTTGCCTTCTGTTTTTGTCTACCCAGAAACGCTGGTGAAGTAAAAG  
ATGCTGAAGATCAGTTGGGTGCACGAGTGGGTACATCGAAGTGGATCTCAACAGCGGTAAGATCCTTGAGAGTTTTCGC  
CCCGAAGAACGTTTTCCAATGATGAGCACTTTTAAAGTTCGCTATGTGGCGCGGTATTATCCCGTATTGACGCCGGCA  
AGAGCAACTCGGTGCGCCGATACACTATTCTCAGAATGACTTGGTTGAGTACTCACCAGTACAGAAAAGCATCTTACGG  
ATGGCATGACAGTAAGAGAATTATGCAGTGTGCCATAACCATGAGTGATAAAGTACTGCGGCAACTTACTTCTGACAACG  
ATCGGAGGACCGAAGGAGCTAACCCTTTTTTGCACAACATGGGGGATCATGTAAGTACTCGCTTGATCGTTGGGAACCGGA  
GCTGAATGAAGCCATACCAACGACGAGCGTGACACCACGATGCCTGCAGCAATGGCAACAACGTTGCGCAAATATTA  
CTGGCGAACTACTTACTCTAGCTTCCCGCAACAATTAATAGACTGGATGGAGGCGGATAAAGTTGCAGGACCACTTCTG  
CGCTCGGCCCTTCCGGCTGGCTGGTTTTATTGCTGATAAATCTGGAGCCGGTGGAGCGTGGTTCTCGCGGTATCATTGCAGC  
ACTGGGGCCAGATGGTAAGCCCTCCCGTATCGTAGTTATCTACACGACGGGGAGTCAAGCAACTATGGATGAACGAAATA  
GACAGATCGCTGAGATAGGTGCCTCACTGATTAAGCATTGGTAAGTGTGACACCAAGTTTACTCATATATACTTTAGATT  
GATTTAAAACCTTCAATTTTTAATTTAAAAGGATCTAGGTGAAGATCCTTTTTGATAATCTCATGACCAAAATCCCTAACG  
TGAGTTTTCGTTCCTGAGCGTCAAGCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTTCTGCGCGTAA  
TCTGCTGCTTGCAACAAAAAAACCACCGCTACCAGCGGTGGTTTTGTTTGCAGGATCAAGAGCTACCAACTTTTTTCCG  
AAGGTAAGTGGCTTCAAGCAGAGCGCAGATACCAATACTGTCCTTCTAGTGTAGCCGTAGTTAGGCCACCACTTCAAGAA  
CTCTGTAGCACCGCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTGCTGCCAGTGGCGATAAGTCTGTCTTA  
CCGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTGGGCTGAACGGGGGGTTCGTGCACACAGCCCAGC  
TTGGAGCGAACGACCTACACCGAACTGAGATACCTACAGCGTGAAGTATGAGAAAGCGCCACGCTTCCGAAGGGAGAAA  
GGCGGACAGGTATCCGGTAAGCGGCAGGGTGGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGGAAACGCCTGGTATC  
TTTATAGTCTGTGCGGTTTTCGCCACCTCTGACTTGAGCGTCAATTTTTGTGATGCTCGTCAAGGGGGGCGGAGCCTATGG  
AAAAACGCCAGCAACGCGGCTTTTTACGGTTCTGGCCTTTTGTGTCCTTTGCTCACATGTTCTTCTGCGTTATC  
CCCTGATTCTGTGGATAACCGTATTACCGCTTTGAGTGAAGTGTATACCGCTCGCCGAGCCGAACGACCGAGCGCAGCG  
AGTCAGTGAAGCAGGAAGCGGAAGAGCGCCTGATGCGGTATTTTCTCCTTACGCATCTGTGCGGTATTTACACCCGATA  
TATGGTGCCTCTCAGTACAATCTGCTCTGATGCCGCATAGTTAAGCCAGTATACACTCCGCTATCGCTACGTGACTGGG  
TCATGGCTGCGCCCCGACACCCGCCAACACCCGCTGACGCGCCCTGACGGGCTTGTCTGCTCCCGCATCCGCTTACAGA  
CAAGCTGTGACCGTCTCCGGGAGCTGCATGTGTGAGAGGTTTTACCAGTATCACCGAAACGCGCAGGAGGAGCTGCGGTA  
AAGCTCATCAGCGTGGTCTGAAGCGATTACAGATGTCTGCCTGTTTATCCGCTCCAGCTCGTTGAGTTTCTCCAGAA  
GCGTTAATGTCTGGCTTCTGATAAAGCGGGCCATGTTAAGGGCGGTTTTTCTGTTTGGTCACTGATGCCTCCGTGTAA  
GGGGGATTTCTGTTTATGGGGTAATGATACCGATGAAACGAGAGAGGATGCTCACGATACGGGTTACTGATGATGAACA  
TGCCCGGTTACTGGAACGTTGTGAGGGTAACAACCTGGCGGTATGGATGCGGCGGGACCAGAGAAAAATCACTCAGGGTC  
AATGCCAGCGCTTCGTTAATACAGATGTAGGTGTTCCACAGGGTAGCCAGCAGCATCCTGCGATGCAGATCCGGAACATA

pE-SUMOstar Amp

ATGGTGCAGGGCGCTGACTTCCGCGTTTTCCAGACTTTACGAAACACGGAAACCGAAGACCATTCATGTTGTTGCTCAGGT  
CGCAGACGTTTTGCAGCAGCAGTCGCTTACGTTTCGCTCGCGTATCGGTGATTATTCTGCTAACCAGTAAGGCAACCCC  
GCCAGCCTAGCCGGGTCCTCAACGACAGGAGCACGATCATGCGCACCCGTGGGGCCGCCATGCCGGCGATAATGGCCTGC  
TTCTCGCCGAAACGTTTTGGTGGCGGGACCAGTGACGAAGGCTTGAGCGAGGGCGTGCAAGATTCCGAATACCGCAAGCGA  
CAGGCCGATCATCGTCGCGCTCCAGCGAAAGCGGTCTCGCCGAAAATGACCCAGAGCGCTGCCGGCACCTGTCCTACGA  
GTTGCATGATAAAGAAGACAGTCATAAGTGCGGCGACGATAGTCATGCCCCGCGCCACCGGAAGGAGCTGACTGGGTTG  
AAGGCTCTCAAGGGCATCGGTGAGATCCCGGTGCCTAATGAGTGAGCTAACTTACATTAATTGCGTTGCGCTCACTGCC  
CGTTTTCCAGTCGGGAAACCTGTCGTGCCAGCTGCATTAATGAATCGGCCAACGCGCGGGGAGAGGGCGGTTTTGCGTATTG  
GGCGCCAGGGTGGTTTTTTCTTTTACCAGTGAGACGGGCAACAGCTGATTGCCCTTACCAGCCTGGCCCTGAGAGAGTTG  
CAGCAAGCGGTCCACGCTGGTTTTGCCCCAGCAGGCGAAAATCCTGTTTGATGGTGGTTAACGGCGGGATATAACATGAGC  
TGTCTTCGGTATCGTCGTATCCCCTACCAGATATCCGCACCAACGCGCAGCCCGGACTCGGTAATGGCGCGCATTGCG  
CCCAGCGCCATCTGATCGTTGGCAACCAGCATCGCAGTGGGAACGATGCCCTCATTACGATTTGCATGGTTTTGTTGAAA  
ACCGGACATGGCACTCCAGTCGCCTTCCCGTTCCGCTATCGGCTGAATTTGATTGCGAGTGAGATATTTATGCCAGCCAG  
CCAGACGCAGACGCGCCGAGACAGAACTTAATGGGCCCGCTAACAGCGCGATTTGCTGGTGACCCAATGCGACCAGATGC  
TCCACGCCCAGTCGCGTACCGTCTTCATGGGAGAAAATAATACTGTTGATGGGTGTCTGGTCAGAGACATCAAGAAATAA  
CGCCGGAACATTAGTGCAGGCAGCTTCCACAGCAATGGCATCCTGGTCATCCAGCGGATAGTTAATGATCAGCCCACTGA  
CGCGTTGCGCGAGAAGATTGTGCACCGCCGCTTTACAGGCTTCGACGCGCTTCGTTCTACCATCGACACCACCACGCTG  
GCACCCAGTTGATCGGCGCGAGATTTAATCGCCGCGACAATTTGCGACGGCGCGTGCAGGGCCAGACTGGAGGTGGCAAC  
GCCAATCAGCAACGACTGTTTTGCCCGCCAGTTGTTGTGCCACGCGGTTGGGAATGTAATTCAGCTCCGCCATCGCCGCTT  
CCACTTTTTCCCGCTTTTTCGCAGAAACGTGGCTGGCCTGGTTACCACGCGGGAAACGGTCTGATAAGAGACACCGGCA  
TACTCTGCGACATCGTATAACGTTACTGGTTTTCACATTCACCACCCTGAATTGACTCTCTTCCGGGCGCTATCATGCCAT  
ACCGCGAAAGGTTTTGCGCCATTCGATGGTGTCCGGGATCTCGACGCTCTCCCTTATGCGACTCCTGCATTAGGAAGCAG  
CCCAGTAGTAGGTTGAGGCCGTTGAGCACCGCCGCGCAAGGAATGGTGCATGCAAGGAGATGGCGCCCAACAGTCCCCC  
GGCCACGGGGCCTGCCACCATACCCACGCCGAAACAAGCGCTCATGAGCCCCGAAGTGGCGAGCCCCGATCTTCCCCATCGG  
TGATGTGGCGATATAGGCGCCAGCAACCGCACCTGTGGCGCCGGTGATGCCGGCCACGATGCGTCCGGCGTAGAGGATC  
G