

>pE-SUMOpro Amp

AGATCTCGATCCCGCAAATTAATACGACTCACTATAGGGGAATTGTGAGCGGATAACAATCCCCTCTA  
GAAATAATTTTGTTTAACTTTAAGAAGGAGATATACCATGGGTCATCACCATCATCATCAGGGTCCCTG  
CAGGACTCAGAAGTCAATCAAGAAGCTAAGCCAGAGGTCAAGCCAGAAGTCAAGCCTGAGACTCACATCA  
ATTTAAAGGTGTCCGATGGATCTTCAGAGATCTTCTTCAAGATCAAAAAGACCACTCCTTTAAGAAGGCT  
GATGGAAGCGTTCGCTAAAAGACAGGGTAAGGAAATGGACTCCTTAAGATTCTTGTACGACGGTATTAGA  
ATTCAAGCTGATCAGGCCCTGAAGATTTGGACATGGAGGATAACGATATTATTGAGGCTCACCGCGAAC  
AGATTGGAGGTTGAGACCACTAGTGGTACCGGTCTCACTAGAGGATCCGAATTCGAGCTCCGTCGACAAG  
CTTGCGGCCGCACTCGAGCACCACCACCACCACCCTGAGATCCGGCTGCTAACAAAGCCCGAAAGGAAG  
CTGAGTTGGCTGCTGCCACCGCTGAGCAATAACTAGCATAACCCCTTGGGGCCTCTAAACGGGTCTTGAG  
GGTTTTTTGCTGAAAGGAGGAATAATCCGGATTGGCGAATGGGACGCGCCCTGTAGCGGCGCATTAA  
GCGCGGCGGGTGTGGTGGTTACGCGCAGCGTGACCGCTACACTTGCCAGCGCCCTAGCGCCCGCTCCTTT  
CGCTTTCTCCCTCCTTTCTCGCCACGTTCCGCGGCTTTCCCCGTCAAGCTCTAAATCGGGGGCTCCCT  
TTAGGGTCCGATTTAGTGCTTTACGGCACCTCGACCCAAAAAATTGATTAGGGTGATGGTTCACGTA  
GTGGGCCATCGCCCTGATAGACGGTTTTTCGCCCTTTGACGTTGGAGTCCACGTTCTTAATAGTGGACT  
CTTGTTCCAACTGGAACAACACTCAACCCTATCTCGGTCTATTCTTTTGATTTATAAGGGATTTTGCCG  
ATTTCCGCCTATTGGTTAAAAAATGAGCTGATTTAACAAAAATTAACGCGAATTTAACAAAAATATAA  
CGTTTACAATTCAGGTGGCACTTTTCGGGGAAATGTGCGCGGAACCCCTATTTGTTATTTTTCTAAAT  
ACATTCAAATATGTATCCGCTCATGAGACAATAACCCTGATAAATGCTTCAATAATATTGAAAAAGGAAG  
AGTATGAGTATTCAACATTTCCGTGTCGCCCTTATCCCTTTTTGCGGCATTTGCTTCTGTTTTG  
CTCACCAGAAACGCTGGTGAAAGTAAAAGATGCTGAAGATCAGTTGGGTGCACGAGTGGGTTACATCGA  
ACTGGATCTCAACAGCGGTAAGATCCTTGAGAGTTTTCGCCCCGAAGAACGTTTTCCAATGATGAGCACT  
TTAAAGTTCTGCTATGTGGCGCGGTATTATCCCGTATTGACGCCGGGCAAGAGCAACTCGGTCGCCGCA  
TACACTATTCTCAGAATGACTTGGTTGAGTACTCACCAGTCACAGAAAAGCATCTTACGGATGGCATGAC  
AGTAAGAGAATTATGCAGTGCTGCCATAACCATGAGTGATAAACTGCGGCCAATTACTTCTGACAACG  
ATCGGAGGACCGAAGGAGCTAACCCTTTTTGCAACAATGCGGGGATCATGTAACCTGCCTTGATCGTT  
GGGAACCGGAGCTGAATGAAGCCATACCAAACGACGAGCGTGACACCACGATGCCTGCAGCAATGGCAAC  
AACGTTGCGCAAATTAATACTGGCGAACTACTTACTCTAGCTTCCCGCAACAATTAATAGACTGGATG  
GAGGCGGATAAAGTTGCAGGACCACTTCTGCGCTCGGCCCTCCGGCTGGCTGGTTATTGCTGATAAAT  
CTGGAGCCGTGAGCGTGGTTCTCGCGGTATCATTGCAGCACTGGGGCCAGATGGTAAGCCCTCCCGTAT  
CGTAGTTATCTACACGACGGGGAGTCAGGCAACTATGGATGAACGAAATAGACAGATCGCTGAGATAGGT  
GCCTCACTGATTAAGCATTGGTAACCTGTCAGACCAAGTTTACTCATATATACTTTAGATTGATTTAAAAC  
TTCATTTTTAATTTAAAAGGATCTAGGTGAAGATCCTTTTTGATAATCTCATGACCAAATCCCTAACG  
TGAGTTTTCGTTCCACTGAGCGTCAGACCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTT  
CTGCGCGTAATCTGCTGCTTCAAACAAAAAACCCCGCTACCAGCGGTGGTTTGTGGCCGGATCAAG  
AGCTACCAACTCTTTTTCCGAAGGTAACCTGCTTCAAGCAAGCGCAGATACCAATACTGCTTCTAGT  
GTAGCCGTAGTTAGGCCACCACTTCAAGAACTCTGTAGCACCGCTACATACCTCGCTCTGCTAATCCTG  
TTACCAGTGGCTGCTGCCAGTGCCGATAAGTCGTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGG  
ATAAGGCGCAGCGTCCGGCTGAACGGGGGGTTCGTGCACACAGCCAGCTTGGAGCGAACGACCTACAC  
CGAACTGAGATACTACAGCGTGAGCTATGAGAAAGCGCCACGCTTCCCGAAGGGAGAAAGGCGGACAGG  
TATCCGGTAAGCGGCAGGGTCCGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGGAAACGCCTGGTATC  
TTTATAGTCTGTGGGTTTTGCCACCTCTGACTTGAGCGTCGATTTTTGTGATGCTCGTCAGGGGGGCG  
GAGCCTATGGAAAACGCCAGCAACGCGGCTTTTTACGGTTCCTGGCCTTTTGTGCTGCTTTTGTCTCAC  
ATGTTCTTCTGCTTATCCCTGATTCTGTGGATAACCGTATTACCGCTTTGAGTGAGCTGATACCG  
CTCGCCGACGCCAACGACCGAGCGCAGCGAGTCACTGAGCGAGGAAGCGGAAGAGCGCCTGATGCGGTA  
TTTTCTCCTTACGCATCTGTGCGGATTTTACACCGCATATATGGTGCACTCTCAGTACAATCTGCTCTG  
ATGCCGCATAGTTAAGCCAGTATACTCCGCTATCGCTACGTGACTGGGTCATGGCTGCGCCCCGACAC  
CCGCCAACCCCGCTGACGCGCCCTGACGGGCTTGTCTGCTCCCGCATCCGCTTACAGACAAGCTGTGA

CCGTCTCCGGGAGCTGCATGTGTCAGAGGTTTTACCCTCATCACCGAAACGCGCGAGGCAGCTGCGGTA  
AAGCTCATCAGCGTGGTCGTGAAGCGATTACAGATGTCTGCTGTTCCATCCGCGTCCAGCTCGTTGAGT  
TTCTCCAGAAGCGTTAATGTCTGGCTTCTGATAAAGCGGGCCATGTTAAGGGCGGTTTTTCTGTTTGG  
TCACTGATGCCTCCGTGTAAGGGGGATTCTGTTTCATGGGGTAATGATACCGATGAAACGAGAGAGGAT  
GCTCACGATACGGGTTACTGATGATGAACATGCCCCGTTACTGGAACGTTGTGAGGGTAACAACACTGGCG  
GTATGGATGCGGCGGGACCAGAGAAAAATCACTCAGGGTCAATGCCAGCGCTTCGTTAATACAGATGTAG  
GTGTTCCACAGGGTAGCCAGCAGCATCCTGCGATGCAGATCCGGAACATAATGGTGCAGGGCGCTGACTT  
CCGCGTTTTCCAGACTTTACGAAACACGGAACCGAAGACCATTTCATGTTGTTGCTCAGGTGCGAGACGTT  
TTGCAGCAGCAGTCGTTACGTTGCTCGGTATCGGTGATTTCATTCTGCTAACAGTAAGGCAACCCC  
GCCAGCCTAGCCGGGTCTCAACGACAGGAGCAGATCATGCGCACCCGTGGGGCCGCCATGCCGGCGAT  
AATGGCCTGCTTCTCGCCGAAACGTTTTGGTGGCGGGACCAGTGACGAAGGCTTGAGCGAGGGCGTGCAAG  
ATCCGAATACCGCAAGCGACAGGCCGATCATCGTCGCGCTCCAGCGAAAGCGGTCCTCGCCGAAAATGA  
CCCAGAGCGCTGCCGGCACCTGTCCTACGAGTTGCATGATAAAGAAGACAGTCATAAGTGCGGCGACGAT  
AGTCATGCCCCGCGCCACCGGAAGGAGCTGACTGGGTTGAAGGCTCTCAAGGGCATCGGTGCGAGATCCC  
GGTGCTAATGAGTGAGCTAATTACATTAATTGCGTTGCGCTCACTGCCCGCTTCCAGTCGGGAAACC  
TGTCGTGCCAGCTGCATTAATGAATCGCCAACGCGCGGGGAGAGGCGGTTTTGCGTATTGGGCGCCAGGG  
TGTTTTTCTTTTACCAGTGAGACGGGCAACAGCTGATTGCCCTTACCAGCCTGGCCCTGAGAGAGTTG  
CAGCAAGCGGTCCACGCTGGTTTGCCCCAGCAGGCGAAAATCTGTTTGATGGTGGTTAACGGCGGGATA  
TAACATGAGCTGTCTTCGGTATCGTCGTATCCACTACCGAGATATCCGCACCAACGCGCAGCCCCGACT  
CGGTAATGGCGCGCATTGCGCCAGCGCCATCTGATCGTTGGCAACCAGCATCGCAGTGGGAACGATGCC  
CTCATTACGATTTGCATGGTTTGTGAAAACCGGACATGGCACTCCAGTCGCCTTCCCGTTCCGCTATC  
GGCTGAATTTGATTGCGAGTGAGATATTTATGCCAGCCAGCCAGACGCAGACGCGCCGAGACAGAACTTA  
ATGGGCCCCGTAACAGCGCGATTTGCTGGTGACCAATGCGACCAGATGCTCCACGCCCAGTCGCGTACC  
GTCTTCATGGGAGAAAATAATACTGTTGATGGGTGTCTGGTCAGAGACATCAAGAAATAACGCCGGAACA  
TTAGTGAGGCAGCTTCCACAGCAATGGCATCCTGGTCATCCAGCGGATAGTTAATGATCAGCCCACTGA  
CGCGTTGCGCGAGAAGATTGTGCACCGCCGCTTACAGGCTTCGACGCCGCTTCGTTCTACCATCGACAC  
CACCACGCTGGCACCCAGTTGATCGGCGCGAGATTTAATCGCCGCGACAATTTGCGACGGCGCGTGCAGG  
GCCAGACTGGAGGTGGCAACGCCAATCAGCAACGACTGTTGCCCGCCAGTTGTTGTGCCACGCGGTTGG  
GAATGTAATTCAGCTCCGCCATCGCCGCTTCCACTTTTTCCCGCTTTTCGAGAAACGTGGCTGGCCTG  
GTTACCACGCGGGAAACGGTCTGATAAGAGACACCGGCATACTCTGCGACATCGTATAACGTTACTGGT  
TTCACATTCACCACCCTGAATTGACTCTTCCGGGCGCTATCATGCCATACCGCGAAAGTTTTGCGCC  
ATTCGATGGTGTCCGGATCTCGACGCTCTCCCTTATGCGACTCCTGCATTAGGAAGCAGCCCAGTAGTA  
GGTTGAGGCCGTTGAGCACCGCCGCGCAAGGAATGGTGCATGCAAGGAGATGGCGCCCAACAGTCCCCC  
GGCCACGGGGCCTGCCACCATACCCACGCCGAAACAAGCGCTCATGAGCCCGAAGTGCGGAGCCCCGATCT  
TCCCCATCGGTGATGTCGGCGATATAGGCGCCAGCAACCGCACCTGTGGCGCCGGTGTATGCCGGCCACGA  
TTCGTCCGGCGTAGAGGATCG