

>pE-SUMO Amp (direct)

AGATCTCGATCCC GCGAAATTAATACGACTCACTATAGGGGAATTGTGAGCGGATAACAATTCCCCTCTA  
GAAATAATTTTTGTTTAACTTTAAGAAGGAGATATACCATGGGTCATCACCATCATCATCACGGGTCCCTG  
CAGGACTCAGAAGTCAATCAAGAAGCTAAGCCAGAGGTCAAGCCAGAAGTCAAGCCTGAGACTCACATCA  
ATTTAAAGGTGTCCGATGGATCTTCAGAGATCTTCTTCAAGATCAAAAAGACCACTCCTTTAAGAAGGCT  
GATGGAAGCGTTTCGCTAAAAGACAGGGTAAGGAAATGGACTCCTTAAGATTCTTGTACGACGGTATTAGA  
ATTTCAAGCTGATCAGGCCCTGAAGATTTGGACATGGAGGATAACGATATTATTGAGGCTCACC GCGAAC  
AGATTGGAGGTTGAGACCACTAGTGGTACCGGTCTCACTAGAGGATCCGAATTCGAGCTCCGTGACAAG  
CTTGCGGCCGCACTCGAGCACCACCACCACCACCCTGAGATCCGGCTGCTAACAAAGCCCCGAAAGGAAG  
CTGAGTTGGCTGCTGCCACCGCTGAGCAATAACTAGCATAACCCCTTGGGGCCTCTAAACGGGTCTTGAG  
GGGTTTTTTGCTGAAAGGAGGAATATATCCGGATTTGGCGAATGGGACGCGCCCTGTAGCGGCGCATTA  
GCGCGGCGGGTGTGGTGGTTACGCGCAGCGTGACCCTACTTACCAGCGCCCTAGCGCCCGCTCCTTT  
CGCTTTCTTCCCTTCTTCTCGCCACGTTTCGCCGGCTTTCCCGTCAAGCTCTAAATCGGGGGCTCCCT  
TTAGGGTTCCGATTTAGTGCTTTACGGCACCTCGACCCCAAAAACTTGATTAGGGTGATGGTTCACGTA  
GTGGGCCATCGCCCTGATAGACGGTTTTTCGCCCTTTGACGTTGGAGTCCACGTTCTTTAATAGTGGACT  
CTTGTTCCAACTGGAACAACACTCAACCCTATCTCGGTCTATTCTTTTGATTTATAAGGGATTTTGCCG  
ATTTCCGCCCTATTGGTTAAAAAATGAGCTGATTTAACAAAAATTTAACGCGAATTTTAACAAAATATTAA  
CGTTTTACAATTTCCAGGTGGCACTTTTCGGGAAATGTGCGCGGAACCCCTATTTGTTTTATTTTCTAAAT  
ACATTTCAAATATGTATCCGCTCATGAGACAATAACCCTGATAAATGCTTCAATAATATTGAAAAAGGAAG  
AGTATGAGTATTCAACATTTCCGTGTCGCCCTTATTCCCTTTTTTTCGGGCATTTTGCCTTCCTGTTTTG  
CTCACCCAGAAACGCTGGTGAAGTAAAAGATGCTGAAGATCAGTTGGGTGCACGAGTGGGTTACATCGA  
ACTGGATCTCAACAGCGGTAAGATCCTTGAGAGTTTTCGCCCCGAAGAACGTTTTCCAATGATGAGCACT  
TTTAAAGTTCTGCTATGTGGCGCGGTATTATCCCGTATTGACGCCGGGCAAGAGCAACTCGGTGCGCCGA  
TACACTATTCTCAGAATGACTTGGTTGAGTACTACCAGTCACAGAAAAGCATCTTACGGATGGCATGAC  
AGTAAGAGAATTATGCAGTGTGCCATAACCATGAGTGATAACACTGCGGCCAACTTACTTCTGACAACG  
ATCGGAGGACCGAAGGAGCTAACCCTTTTTTGACAACATGGGGGATCATGTAACCTCGCCTTGATCGTT  
GGGAACCGGAGCTGAATGAAGCCATACCAAACGACGAGCGTGACACCACGATGCCTGCAGCAATGGCAAC  
AACGTTGCGCAAACCTAATACTGGCGAACTACTTACTCTAGCTTCCCGGCAACAATTAATAGACTGGATG  
GAGGCGGATAAAGTTGCAGGACCACTTCTGCGCTCGGCCCTTCCGGCTGGCTGGTTTTATTGCTGATAAAT  
CTGGAGCCGGTGAGCGTGGTTCTCGCGGTATCATTGCAGCACTGGGGCCAGATGGTAAGCCCTCCCGTAT  
CGTAGTTATCTACACGACGGGGAGTCAGGCAACTATGGATGAACGAAATAGACAGATCGCTGAGATAGGT  
GCCTCACTGATTAAGCATTGGTAACTGTCAGACCAAGTTTACTCATATATACTTTAGATTGATTTAAAC  
TTCATTTTTAATTTAAAAGGATCTAGGTGAAGATCCTTTTTGATAATCTCATGACCAAAATCCCTTAACG  
TGAGTTTTCGTTCCTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTTCTTGAGATCCTTTTTTT  
CTGCGCGTAATCTGCTGCTTGCAAACAAAAAACCACCGCTACCAGCGGTGGTTTTGTTTGCCGGATCAAG  
AGCTACCAACTCTTTTTCCGAAGGTAACCTGGCTTCAGCAGAGCGCAGATACCAAATACTGTCCTTCTAGT  
GTAGCCGTAGTTAGGCCACCCTTCAAGAACTCTGTAGCACC GCCTACATACTCGCTCTGCTAATCCTG  
TTACCAGTGGCTGCTGCCAGTGGCGATAAGTCGTGTCTTACC GGGTTGGACTCAAGACGATAGTTACCGG  
ATAAGGCGCAGCGGTGCGGCTGAACGGGGGTTCTGTGCACACAGCCCAGCTTGGAGCGAACGACCTACAC  
CGAACTGAGATACCTACAGCGTGAGCTATGAGAAAGCGCCACGCTTCCCGAAGGGAGAAAGGCGGACAGG  
TATCCGGTAAGCGGCAGGGTCCGGAACAGGAGAGCGCACGAGGGAGCTTCCAGGGGGAAACGCCTGGTATC  
TTTATAGTCCTGTCCGGTTTTCGCCACCTCTGACTTGAGCGTCGATTTTTTGTGATGCTCGTCAGGGGGCG  
GAGCCTATGAAAAACGCCAGCAACGCGGCCTTTTTACGGTTTCTGGCCTTTTGTGCTGGCCTTTTGTCTAC  
ATGTTCTTTCTGCGTTATCCCTGATTCTGTGGATAACCGTATTACC GCTTTGAGTGAGCTGATACCG  
CTCGCCGACCGAACGACCGAGCGCAGCGAGTCAGTGAGCGAGGAAGCGGAAGAGCGCCTGATGCGGTA  
TTTTCTCCTTACGCATCTGTGCGGTATTTACACCCGATATATGGTGC ACTCTCAGTACAATCTGCTCTG  
ATGCCCATAGTTAAGCCAGTATACACTCCGCTATCGCTACGTGACTGGGT CATGGCTGCGCCCCGACAC  
CCGCCAACACCCGCTGACGCGCCCTGACGGGCTTGTCTGCTCCCGGCATCCGCTTACAGACAAGCTGTGA  
CCGTCTCCGGGAGCTGCATGTGTGAGAGTTTTTACC GTCATACCGAAACGCGCGAGGCAGCTGCGGTA  
AAGCTCATCAGCGTGGTGTGAGCGATTACAGATGTCTGCCTGTT CATCCGCGTCCAGCTCGTTGAGT  
TTCTCCAGAAGCGTTAATGTCTGGCTTCTGATAAAGCGGGCCATGTTAAGGGCGGTTTTTTCTGTTTGG  
TCACTGATGCCTCCGTGTAAGGGGGATTTCTGTT CATGGGGTAATGATAACCGATGAAACGAGAGAGGAT  
GCTCACGATACGGTTACTGATGATGAACATGCCCGGTTACTGGAACGTTGTGAGGGTAAACA ACTGGCG  
GTATGGATGCGGCGGGACCAGAGAAAAATCACTCAGGGTCAATGCCAGCGCTTCGTTAATACAGATGTAG  
GTGTTCCACAGGGTAGCCAGCAGCATCCTGCGATGCAGATCCGGAACATAATGGTGCAGGGCGCTGACTT  
CCGCGTTTTCCAGACTTTACGAAACACGGAAACCGAAGACCATT CATGTTGTTGCTCAGGTGCGAGACGTT  
TTGAGCAGCAGTCGCTTACGTTTCGCTCGCGTATCCGGT GATTATTCTGCTAACCAGTAAGGCAACCC

GCCAGCCTAGCCGGGTCTCAACGACAGGAGCACGATCATGCGCACCCGTGGGGCCGCCATGCCGGCGAT  
AATGGCCTGCTTCTCGCCGAAACGTTTTGGTGGCGGGACCAGTGACGAAGGCTTGAGCGAGGGCGTGCAAG  
ATTCCGAATACCGCAAGCGACAGGCCGATCATCGTCGCGCTCCAGCGAAAGCGGTCTCGCCGAAAATGA  
CCCAGAGCGCTGCCGGCACCTGTCCTACGAGTTGCATGATAAAGAAGACAGTCATAAGTGCGGGCGACGAT  
AGTCATGCCCCGCGCCACCAGGAGGAGCTGACTGGGTTGAAGGCTCTCAAGGGCATCGGTGAGATCCC  
GGTGCCTAATGAGTGAGCTAACTTACATTAATTGCGTTGCGCTCACTGCCCGCTTTCCAGTCGGGAAACC  
TGTCGTGCCAGCTGCATTAATGAATCGGCCAACGCGCGGGGAGAGGCGGTTTGCGTATTGGGCGCCAGGG  
TGGTTTTTCTTTTACCAGTGAGACGGGCAACAGCTGATTGCCCTTACCAGCTGGCCCTGAGAGAGTTG  
CAGCAAGCGGTCCACGCTGGTTTTGCCCCAGCAGGCGAAAATCCTGTTTGATGGTGGTTAACGGCGGGATA  
TAACATGAGCTGTCTTCGGTATCGTCGTATCCCCTACCAGATATCCGCACCAACGCGCAGCCCCGGACT  
CGGTAATGGCGCGCATTGCGCCAGCGCCATCTGATCGTTGGCAACCAGCATCGCAGTGGGAACGATGCC  
CTCATTACGATTTGCATGGTTTTGTTGAAAACCGGACATGGCACTCCAGTCGCCTTCCCGTTCCGCTATC  
GGCTGAATTTGATTGCGAGTGAGATATTTATGCCAGCCAGCCAGACGCAGACGCGCCGAGACAGAACTTA  
ATGGGCCCCGCTAACAGCGCGATTTGCTGGTGACCCAATGCGACCAGATGCTCCACGCCCAGTCGCGTACC  
GTCTTCATGGGAGAAAATAATACTGTTGATGGGTGTCTGGTCAGAGACATCAAGAAAATAACGCCGGAACA  
TTAGTGCAGGCAGCTTCCACAGCAATGGCATCCTGGTCATCCAGCGGATAGTTAATGATCAGCCCCACTGA  
CGCGTTGCGCGAGAAGATTGTGCACCGCCGCTTTACAGGCTTCGACGCCGCTTCGTTCTACCATCGACAC  
CACCACGCTGGCACCCAGTTGATCGGCGCGAGATTTAATCGCCGCGACAATTTGCGACGGCGCGTGCAGG  
GCCAGACTGGAGGTGGCAACGCCAATCAGCAACGACTGTTTGGCCGCCAGTTGTTGTGCCACGCGGTTGG  
GAATGTAATTCAGCTCCGCCATCGCCGCTTCCACTTTTTCCCGGTTTTTCGAGAAACGTGGCTGGCCTG  
GTTACCACGCGGGAAACGGTCTGATAAGAGACACCGGCATACTCTGCGACATCGTATAACGTTACTGGT  
TTCACATTCACCACCCTGAATTGACTCTCTTCCGGGCGCTATCATGCCATACCGCGAAAGGTTTTGCGCC  
ATTCGATGGTGTCCGGGATCTCGACGCTCTCCCTTATGCGACTCCTGCATTAGGAAGCAGCCCAGTAGTA  
GGTTGAGGCCGTTGAGCACCGCCGCCGCAAGGAATGGTGCATGCAAGGAGATGGCGCCCAACAGTCCCCC  
GGCCACGGGGCCTGCCACCATACCCACGCCGAAACAAGCGCTCATGAGCCCCAAGTGGCGAGCCCCGATCT  
TCCCCATCGGTGATGTCGGCGATATAGGCGCCAGCAACCGCACCTGTGGCGCCGGTATGCCGGCCACGA  
TGCGTCCGGCGTAGAGGATCG