

pE-SUMOstar Kan

>pE-SUMOstar Kan (5628 bp)

AGATCTCGATCCCGCAAATTAATACGACTACTATAGGGGAATTGTGAGCGGATAACAATTCCCCTCTAGAAATAATTT
TGTTTAACTTTAAGAAGGAGATATACCATGGGTCATCACCATCATCATCACGGGTCCTGCAGGACTCAGAAGTCAATCA
AGAAGCTAAGCCAGAGGTCAAGCCAGAAGTCAAGCCTGAGACTCACATCAATTTAAAGGTGTCCGATGGATCTTCAGAGA
TCTTCTTCAAGATCAAAAAGACCACTCCTTTAAGAAGGCTGATGGAAGCGTTTCGCTAAAAGACAGGGTAAGGAAATGGAC
TCCTTAAACGTTCTTGTACGACGGTATTGAAATTCAAGCTGATCAGACCCCTGAAGATTTGGACATGGAGGATAACGATAT
TATTGAGGCTCACCGCGAACAGATTGGAGGTTGAGACCACTAGTGGTACCGGTCTCACTAGAGGATCCGAATTCGAGCTC
CGTCGACAAGCTTGCGGCCGCACTCGAGCACCACCACCACCACCCTGAGATCCGGCTGCTAACAAAGCCCCGAAAGGAAG
CTGAGTTGGCTGCTGCCACCGCTGAGCAATAACTAGCATAACCCCTTGGGGCCTCTAAACGGGTCTTGAGGGGTTTTTGTG
CTGAAAGGAGGAATAATCCGGATTGGCGAATGGGACGCGCCCTGTAGCGGCGCATTAAAGCGCGCGGGTGTGGTGGTT
ACGCGCAGCGTGACCGCTACACTTGCCAGCGCCCTAGCGCCCGCTCCTTTCGCTTCTTCCCTTCTTCTCGCCACGTT
CGCCGGCTTTCCTCGTCAAGCTCTAAATCGGGGGCTCCTTTAGGGTTCGATTTAGTGCTTTACGGCACCTCGACCCCA
AAAACTTGATTAGGGTGTGGTTCACGTAGTGGGCCATCGCCCTGATAGACGGTTTTTCGCCCTTTCGAGTTGGAGTCC
ACGTTCTTTAATAGTGGACTCTTGTCCAACTGGAACAACACTCAACCCTATCTCGGTCTATTCTTTTGATTTATAAGG
GATTTTGCCGATTTGCGCCTATTGGTTAAAAAATGAGCTGATTTAACAAAAATTTAACGCGAATTTTAAACAAAATATTA
CGTTTACAATTTAGGTGGCACTTTTCGGGAAATGTGCGCGGAACCCCTATTTGTTATTTTTCTAAATACATTCAAT
ATGTATCCGCTCATGAATTAATTCTTAGAAAACTCATCGAGCATCAAATGAAACTGCAATTTATTATATCAGGATTAT
CAATACCATATTTTTGAAAAAGCCGTTTCTGTAATGAAGGAGAAAACCTCACCGAGGCAGTTCATAGGATGGCAAGATCC
TGGTATCGGTCTGCGATTCCGACTCGTCCAACATCAATAACCTATTAATTTCCCTCGTCAAAAATAAGGTTATCAAG
TGAGAAATCACCATGAGTGACGACTGAATCCGGTGAGAATGGCAAAAGTTTATGCATTTCTTCCAGACTTGTCAACAG
GCCAGCCATTACGCTCGTCATCAAAATCACTCGCATCAACCAAACCGTTATTCATTCTGATTGCGCCTGAGCGAGACGA
AATACGCGATCGCTGTTAAAAGGACAATTACAAACAGGAATCGAATGCAACCGGCGCAGGAACACTGCCAGCGCATCAAC
AATATTTTACCTGAATCAGGATATTCTTCTAATACCTGGAATGCTGTTTTCCCGGGGATCGCAGTGGTGAATACCATG
CATCATCAGGAGTACGGATAAAATGCTTATGGTTCGGAAGAGGCATAAATCCGTCAGCCAGTTTAGTCTGACCATCTCA
TCTGTAACATCATTGGCAACGCTACCTTGGCATGTTTTCAGAAACAACCTCTGGCGCATCGGGCTTCCCATACAATCGATA
GATTGTGCGACCTGATTGCCCGACATTATCGCGAGCCATTTATACCCATATAAATCAGCATCCATGTTGGAATTTAATC
GCGGCCTAGAGCAAGACGTTTTCCCGTTGAATATGGCTCATAACACCCCTTGTATTACTGTTTATGTAAGCAGACAGTTTT
ATTGTTTCATGACCAAAATCCCTTAACTGAGTTTTCGTTCCACTGAGCGTCAGACCCCGTAGAAAAGATCAAAGGATCTT
CTTGAGATCCTTTTTTCTGCGCGTAATCTGCTGCTTGCAAACAAAAAACCACCGCTACCAGCGGTGGTTGTTTGCCG
GATCAAGAGCTACCAACTCTTTTTCCGAAGGTAACCTGGCTTTCAGCAGAGCGCAGATACCAAATACTGTCCTTCTAGTGTA
GCCGTAGTTAGGCCACCACTTCAAGAACTCTGTAGCACCGCCTACATACCTCGCTCTGCTAATCCTGTTACCAGTGGCTG
CTGCCAGTGGCGATAAGTCGTGTCTTACCGGGTTGGACTCAAGACGATAGTTACCGGATAAGGCGCAGCGGTGGGCTGA
ACGGGGGGTTCGTGCACACAGCCAGCTTGGAGCGAACGACCTACACCGAAGTACCTACAGCGTGAGCTATGAGA
AAGCGCCACGCTTCCCGAAGGGAGAAAGGCGGACAGGTATCCGGTAAGCGGCAGGGTCGGAACAGGAGAGCGCACGAGGG
AGCTTCCAGGGGGAAACGCCTGGTATCTTTATAGTCTGTGGGTTTTCGCCACCTCTGACTTGAGCGTCGATTTTTGTGA
TGCTCGTCAGGGGGGCGGAGCCTATGGAAAAACGCCAGCAACGCGGCCTTTTTACGGTTCCTGGCCTTTTGTGGCCTTT
TGCTCACATGTTCTTCTGCGTTATCCCCTGATTCTGTGGATAACCGTATTACCGCCTTTGAGTGAGCTGATACCGCTC
GCCGACCCGAACGACCGAGCGCAGCGAGTCACTGAGCGAGGAAGCGGAAGAGCGCCTGATGCGGTATTTTCTCCTTACG
CATCTGTGCGGTATTTACACCCGCATATATGGTGCCTCTCAGTACAATCTGCTCTGATGCCGCATAGTTAAGCCAGTAT
ACACTCCGCTATCGCTACGTGACTGGGTATGGCTGCGCCCCGACACCCGCCAACACCCGCTGACGCGCCCTGACGGGCT
TGTCTGCTCCCGGCATCCGCTTACAGACAAGCTGTGACCGTCTCCGGGAGCTGCATGTGTGAGAGTTTTACCCTCATC
ACCGAAACGCGCGAGGACGCTGCGGTAAGCTCATCAGCGTGGTGTGAAGCGATTACAGATGCTGCCTGTTTATCCG
CGTCCAGCTCGTTGAGTTTTCTCCAGAAGCGTTAATGTCTGGCTTCTGATAAAGCGGGCCATGTTAAGGGCGGTTTTTCC
TGTTTTGGTCACTGATGCCTCCGTGTAAGGGGGATTTCTGTTTATGGGGGTAATGATACCGATGAAACGAGAGAGGATGCT
CACGATACGGGTTACTGATGATGAACATGCCCGGTTACTGGAACGTTGTGAGGGTAAACAACCTGGCGGTATGGATGCGGC
GGGACCAGAGAAAAATCACTCAGGGTCAATGCCAGCGCTTCTGTTAATACAGATGTAGGTGTTCCACAGGGTAGCCAGCAG
CATCTGCGATGCAGATCCGGAACATAATGGTGCAGGGCGCTGACTTCCGCGTTTTCCAGACTTTACGAAACACGGAACCC
GAAGACCATTATGTTGTTGCTCAGGTCGACAGCGTTTTGCAGCAGCAGTCGTTTACGTTTCGCTCGCGTATCGGTGATT

pE-SUMOstar Kan

CATTCTGCTAACCAGTAAGGCAACCCCGCCAGCCTAGCCGGTCTCAACGACAGGAGCACGATCATGCGCACCCGTGGG
GCCGCCATGCCGGCGATAATGGCCTGCTTCTCGCCGAAACGTTTGGTGGCGGGACCAGTGACGAAGGCTTGAGCGAGGGC
GTGCAAGATTCCGAATACCGCAAGCGACAGGCCGATCATCGTCGCGCTCCAGCGAAAGCGGTCTCGCCGAAAATGACCC
AGAGCGCTGCCGGCACCTGTCCTACGAGTTGCATGATAAAGAAGACAGTCATAAGTGCGGGCAGCATAGTCATGCCCCGC
GCCACCGGAAGGAGCTGACTGGGTTGAAGGCTCTCAAGGGCATCGGTGCGAGATCCCGGTGCCTAATGAGTGAGCTAACT
TACATTAATTGCGTTGCGCTCACTGCCCCGTTTTCCAGTCGGGAAACCTGTCGTGCCAGCTGCATTAATGAATCGGCCAAC
GCGCGGGGAGAGGCGGTTTTGCGTATTGGGCGCCAGGGTGGTTTTTCTTTTACCAGTGAGACGGGCAACAGCTGATTGCC
CTTACCAGCCTGGCCCTGAGAGAGTTGCAGCAAGCGGTCCACGCTGGTTTGGCCCAGCAGGCGAAAATCCTGTTTGATGG
TGGTTAACGGCGGGATATAACATGAGCTGTCTTCGGTATCGTCGTATCCCACTACCGAGATATCCGCACCAACGCGCAGC
CCGACTCGGTAATGGCGCGCATTGCGCCAGCGCCATCTGATCGTTGGCAACCAGCATCGCAGTGGGAACGATGCCCTC
ATTCAGCATTTGCATGGTTTTGTTGAAAACCGGACATGGCACTCCAGTCGCCTTCCCGTTCGGCTATCGGCTGAATTTGAT
TGCGAGTGAGATATTTATGCCAGCCAGCCAGACGCAGACGCGCCGAGACAGAATTAATGGGCCCGCTAACAGCGCGATT
TGCTGGTGACCCAATGCGACCAGATGCTCCACGCCAGTCGCGTACCGTCTTCATGGGAGAAAATAATACTGTTGATGGG
TGTCTGGTCAGAGACATCAAGAAATAACGCCGGAACATTAGTGACGGCAGCTTCCACAGCAATGGCATCCTGGTCATCCA
GCGGATAGTTAATGATCAGCCCACTGACGCGTTGCGCGAGAAGATTGTGCACCGCCGCTTTACAGGCTTCGACGCGCTT
CGTTCTACCATCGACACCACCGCTGGCACCCAGTTGATCGGCGCGAGATTTAATCGCCGCGACAATTTGCGACGGCGC
GTGCAGGGCCAGACTGGAGGTGGCAACGCCAATCAGCAACGACTGTTTGGCCGCGAGTTGTTGTGCCACGCGGTTGGGAA
TGTAATTCAGCTCCGCCATCGCCGCTTCCACTTTTTCCCGGTTTTTCGAGAAACGTGGCTGGCCTGGTTACCACGCGG
GAAACGGTCTGATAAGAGACACCGGCATACTCTGCGACATCGTATAACGTTACTGGTTTTACATTCACCACCCTGAATTG
ACTCTTTCCGGGCGCTATCATGCCATACCGCGAAAGTTTTGCGCCATTCGATGGTGTCCGGGATCTCGACGCTCTCCC
TTATGCGACTCCTGCATTAGGAAGCAGCCAGTAGTAGGTTGAGGCCGTTGAGCACCGCCGCGCAAGGAATGGTGCATG
CAAGGAGATGGCGCCCAACAGTCCCCCGGCCACGGGGCCTGCCACCATACCCACGCCGAAACAAGCGCTCATGAGCCCGA
AGTGGCGAGCCGATCTTCCCCATCGGTGATGTCGGCGATATAGGCGCCAGCAACCGCACCTGTGGCGCCGGTGATGCCG
GCCACGATGCGTCCGGCGTAGAGGATCG