

>EGFR wt sequence

atgcgaccctccgggacggccggggcagcgcctcctggcgctgctggctgctgctctgcccggcgagtcgggctctggaggaaaagaaagtgtccaag
gcacgagtaacaagctcacgcagttgggcacttttgaagatcattttctcagcctccagaggatgttcaataactgtgaggtggccttgggaatttggga
aattacatgtgagaggaattatgatctttccttctaaagaccatccaggaggtggctggttatgtcctcattgcctcaacacagtgaggcgaatt
cctttggaaaacctgcagatcatcagaggaaatgtactacgaaaattcctatgccttagcagcttatctaactatgatgcaataaaaacgggactg
aaggagctgcccagagaaatttacaggaaatcctgcatggcgccgtgcggttcagcaacaacctgcctgtgcaacgtggagagcatccagtggc
gggacatagtcagcagtgactttctcagcaaatgctgatggacttccagaacctgggagcagctgcaaaaagtgtgatccaagctgtccaatggg
agctgctggggtgcaggagaggagaactgccagaaactgacaaaatcatctgtgccagcagtgctccggcgctgcccgtggcaagtccccagtg
actgctgccacaaccagtgtgctgcaggctgcacagggccccgggagagcgactgcctggctgcccgaattccgagacgaagccacgtgcaagga
cacctgccccactcatgcttcaaccccaccacgtaccagatggatgtgaaccccaggggcaatacagctttggtgccacctgctgaagaagt
gtccccgaattatggtgacagatcacggctcgtgctccgagcctgtggggccgacagctatgagatggaggaagacggcgtccgcaagtgtaa
gaagtgcgaaggccttgcgcaaaagtgtgaacggaataggtattggtgaatttaaagactcactctcataaatgctacgaatattaacacttca
aaaactgcacctcatcagtgccgatctccacatcctgcccgtggcatttaggggtgactccttcacacatactcctcctctggatccacaggaactgg
atattctgaaaaccgtaaaggaaatcacagggttttctgattcaggcttggcctgaaaacaggacggacctccatgccttgagaacctgaaatc
atacggcgaggaccaagcaacatggctcagtttctctgagctcgtcagcctgaacataacatccttgggattacgctccctcaaggagataagt
ggagatgtgataattcaggaacaaaaatttgtgctatgcaatacaataaactgaaaaaactgtttgggacctccggtcagaaaacaaaaattat
aagcaacagaggtgaaaacagctgcaaggccacaggccaggtctgcatgcttgtgctccccgagggtgctgctggggccggagcccagggactg
cgtctcttgcggaatgtcagccgaggcagggaatgctggacaagtcaaccttctggagggtgagccaaggaggttgggagaactctgagtg
atacagtgccaccagagtgctgcctcaggccatgaacatcacctgcacaggacggggaccagacaactgtatccagtggtcccactacattgacg
gccccactgctcaagacctgcccggcaggagtcagggagaaaacaacacctggctggaagtacgcagacgcccggccatgtgtgccacctgtg
ccatcaaaactgcacctacggatgactgggacaggtctgaaggcttcaacgaatgggctaaaatcccgtccatgcctcaggatggtggggg
ccctccttctgctgctggtggtggccctggggatcgccctctcatgcaaggcgcacatcgttcggaagcgcacgctgaggaggtgctgcaggaga
gggagcttggagcctctacaccagtgagaaactccaaccaagctctctgaggatctgaaggaaactgaattcaaaaagatcaaatgctg
ggctccggtgcttggcagcgtgtataaggactctggatcccagaaggtgagaaagttaaattcccgtcgtatcaaggaattaagagaagcaa
catctcgaagcaacaaggaaatcctcgatgaagcctacgtgatggcagcgtggacaacccccacgtgtgcccctgctgggcatctgctcacc
tccacctgcaactcatcacgcagctcatgccctcggctgctcctgactatgtccgggaacacaagacaatattggctcccagtacctgctcaac
tgggtgtgagatgcaagggcatgaactacttgaggagacctgcttgggtgaccgcgacctggggccaggaaactgactggtgaaaacaccg
agcatgtcaagatcacagatttgggctggccaaactgctgggtgcggaagagaaagaataccatgcagaaggaggcaaagtcctatcaagtga
tggcattggaatcaattctacacagaatctatacccaccagagtgatgtctggagctacggggtgacctttgggagttgatgaccttggatccaagc
catatgacggaatccctgccagcagatctcctcatcctggagaaaggagaacgctccctcagccaccatattgacctgatgtctacatgatca
tggtaagtgtggatgatagcagatagtcgcccgaagtccgtgagttgatcatgaattctcaaaaatggcccagacccccagcgtaccttg
tattcaggggatgaaagaatgattgccaagtctacagactccaacttctaccgtgcctgatggatgaagaagacatggacgacgtggtggat
gccgacgagtacctatcccacagcagggtcttcagcagcccctcacgtcacggactcccctcctgagctctctgagtgcaaccagcaacaattcc
accgtggcttgattgatagaatgggctgcaagctgtccatcaaggaagacagcttctgagcgatacagctcagacccccagggccttgac
tggagacagcatagacacacctcctcccagtgctgaatacataaaccagtccgttccaaaaggcccgtggctctgtgcagaatcctgtctatca
caatcagccttgaaccccgcgccagcagagaccacactaccaggacccccagcactgcagtggggcaaccccagatctcaacactgtccag
cccactgtgcaacagcattcgacagccctgccactgggcccagaaaggcagccaccaaattagcctggacaacctgactaccagcaggact
tctttccaaggaagccaagccaatggcatcttaagggtccacagctgaaaatgcagaatacctaagggtcgcccaaaagcagtgaaattatt
ggagca