

>EGFR del 19

atcgaccctccgggacggccggggcagcgcctcctggcgtgctggctgcgctctgccggcgagtgggctctggaggaaaagaaagttgccaag
gcacgagtaacaagctcacgcagttgggcacttttgaagatcattttctcagcctccagaggatgttcaataactgtgaggtggtccttgggaatttggga
aattacatgtgagaggaattatgatctttccttctaaagaccatccaggaggtggctggttatgtcctcattgcctcaacacagtgagcgaatt
cctttggaaaacctgcagatcatcagaggaaatgtactacgaaaattcctatgccttagcagcttatctaactatgatgcaataaaaacgggactg
aaggagctgcccatagagaaatttacaggaaatcctgcatggcgcgctgcggttcagcaacaacctgcccctgtgcaacgtggagagcatccagtggc
gggacatagtcagcagtgactttctcagcaacatgctgatggacttcagaaccacctgggagcagctgcaaaaagtgtgatccaagctgtccaatggg
agctgctggggtgcaggagaggagaactgccagaaactgacaaaatcatctgtgccagcagtgctccgggctgctccgtggcaagtccccagtg
actgctgccacaaccagtgtgctgcaggctgcacagggccccgggagagcgcactgctggctgcccgaattccgagacgaagccacgtgcaagga
cacctgccccactcatgcttcaaccccaccacgtaccagatggatgtgaaccccaggggcaatacagctttggtgccacctgctgaagaagt
gtccccgtaattatgtggtgacagatcacggctcgtgctccgagcctgtggggccgacagctatgagatggaggaagacggcgtccgcaagtgtaa
gaagtgcgaaggccttgcgcaaaagtgtgaacggaataggatgtggaatttaagactcactctcataaatgctacgaatattaacacttca
aaaactgcacctcatcagtggtgctcctcactcctgcccgtggcatttaggggtgactccttcaacatactcctcctctggtaccacaggaactgg
atattctgaaaaccgtaaaggaaatcacagggttttctgctgattcaggcttggcctgaaaacaggacggacctccatgccttgagaacctgaaatc
atacggcgaggaccaagcaacatggctcagtttctctgagctcgtcagcctgaacataacatccttgggattacgctccctcaaggagataagt
ggagatgtgataattcaggaacaaaaatttctgctatgcaatacaataaactgaaaaaactgtttgggacctccggtcagaaaacaaaaattat
aagcaacagaggtgaaaacagctgcaaggccacaggccaggtctgcatgcttctgctccccgagggtgctgctggggccggagcccaggggactg
cgtctcttgcggaatgtcagccgaggcagggaatgctggacaagtcaaccttctggagggtgagccaaggaggttgggagaactctgagtg
atacagtgccaccagagtgctgctcaggccatgaacatcacctgcacaggacggggaccagacaactgtatccagtggtgcccactacattgacg
gccccactgctcaagacctgcccggcaggagtcagggagaaaacaacacctggctggaagtacgcagacgcccggccatgtgtgccacctgtg
ccatcaaaactgcacctacggatgactgggacaggtctgaaggctgccaacgaatgggctcaaaatcccgtccatgcccactgggatggtggggg
cctcctcttctgctgctggtggtggccctggggatcgccctcttcatgcaaggcgcacatcgttccggaagcgcaagctgctggaggtgctg
gggagcttggagcctcttacaccagtgagaaactccaaccaagctctctgaggatctgaaggaaactgaattcaaaaagatcaaatgctg
ggctccggtgcttggcagcgggtgataaggactctggatcccagaaggtgagaaagttaaattcccgtcgtatcaagacatctccgaaagcaa
caaggaaatcctgatgaagcctacgtgatggccagcgtggacaacccccacgtgctgcccctgctgggcatctgctcactccacctgcaactca
tcacgcagctcatgccctcggctgctcctggactatgtcgggaaacaaaagacaatattggctcccagctgctcaactggtgtgtgagatcg
caaaggcatgaactactggaggacctgcttgggtgaccgcgacctggcggccaggaacgtactggtgaaaaaccgcagcatgtcaagatca
cagatgggctggccaaactgctgggtgcggaagagaaagaataccatgcagaaggaggcaaagtcctatcaagtggatggcattggaatcaa
ttctacacagaatctatacccaccagagtgatgtctggagctacggggtgacctttgggagttgatgacctttggatccaagccatgatgaggaatcc
ctgcccagcagatctcctcatcctggagaaaggagaaacgctccctcagccaccatgtaccatcagatgtctacatgatcatggtcaagtgtgga
tgatagacgcagatagtcgcccaggtccgtgagttgatcatgaattctcaaaatggcccagaccccagcgtacctgtcattcagggggatg
aaagaatgcatgttccaagtctacagactccaacttctaccgtgcccctgatggatgaagaagacatggacgacgtggtggatgccgacgagtacct
atcccacagcagggtcttccagcagcccctccacgtcacggactcccctcctgagctctctgagtgcaaccagcaacaattccacctggtctgattg
atagaaatgggctgcaaagctgtcccataaggaagacagcttctgagcgatagctcagaccccacaggcgcctgactgaggacagcataga
cgacacctctcccagtgctgaatacataaacagctcgttccaaaaggcccgtgctgctgagaaatcctgtctatcaaatcagcctctgaa
ccccgcccagcagagaccacactaccagaccccacagcactgagtgggcaaccccagatctcaacactgtccagcccactgtgtcaac
agcattcagacgcctgcccactgggcccagaaaggcagccaccaaatagcctggacaacctgactaccagcaggactcttcccaaggaag
ccaagccaaatggcatcttaagggtccacagctgaaaatgcagaatacctaagggtcgcgcacaaaagcagtgaaattattggagca