

>EGFR del19 T790M sequence

atgcgacctccgggacggccggggcagcgcctcctggcgtgctggctgcgctctgccggcgagtcgggctctggaggaaaagaaagttagccaag
gcacgagtaacaagctcacgcagttgggcacttttgaagatcattttctcagcctccagaggatgttaataactgtgaggtggccttgggaatttggga
aattacctatgtgcagaggaattatgatctttccttctaaagaccatccaggaggtggctggttatgtcctcattgcctcaacacagtgagcgaatt
cctttggaaaacctgcagatcatcagaggaaatgtactacgaaaattcctatgccttagcagcttatctaactatgatgcaataaaaacgggactg
aaggagctgcccagagaaatttacaggaaatcctgcatggcgcgctgcggttcagcaacaacctgcctgtgcaacgtggagagcatccagtggc
gggacatagtcagcagtgactttctcagcaacatgctgatggacttccagaaccctgggagcagctgcaaaaagtgtgatccaagctgtccaatggg
agctgctggggtgcaggagaggagaactgccagaaactgacaaaatcatctgtgccagcagtgctccgggctgctccgtggcaagtccccagtg
actgctgccacaaccagtgtgctgcaggctgcacagggccccgggagagcagctgctggctgcccgaattccgagacgaagccacgtgcaagga
cacctgccccactcatgcttcaaccccaccacgtaccagatggatgtgaaccccaggggcaatacagcttgggtgccacctgctgaagaagt
gtccccgtaattatgtggtgacagatcacggctcgtgctccgagcctgtggggccgacagctatgagatggaggaagacggcgtccgcaagttaa
gaagtgcgaaggccttccgcaaaagtgtgaacggaataggtattgtgaatttaaagactcactctcataaatgctacgaatattaacacttca
aaaactgcacctccatcagtggtgctcctccatcctgcccgtggcatttaggggtgactccttcaacatactcctcctctggatccacaggaactgg
atattctgaaaaccgtaaaggaaatcacagggttttctgattcaggcttggcctgaaaacaggacggacctccatgccttgagaacctagaaatc
atcgcggcaggaccaagcaacatggctcagtttctctgagctcgtcagcctgaacataacatccttgggattacgctccctcaaggagataagt
ggagatgtgataattcaggaacaaaaatttgtgctatgcaatacaataaactggaaaaaactgtttgggacctccggtcagaaaacaaaaattat
aagcaacagaggtgaaaacagctgcaaggccacaggccaggtctgcatgcttctgctccccgagggtgctgctggggccggagcccaggggactg
cgtctcttccggaatgtcagccgaggcagggaatgctggacaagtcaaccttctggagggtgagccaaggaggttgggagaactctgagtg
atacagtgccaccagagtgctgctcaggccatgaacatcacctgcacaggacggggaccagacaactgtatccagtggtgcccactacattgacg
gccccactgctcaagacctgcccggcaggagtcagggagaaaacaacacctggtctggaagtacgcagacgcccggccatgtgtgccacctgtg
ccatcaaaactgcacctacggatgactgggcccaggtctgaaggctgccaacgaatgggcccataatcccgtccatgcacctgggatggtggggg
cctcctcttctgctgctgggtggccctggggatcgcccttctatgcgaaggcgcacatcgttccggaagcgcaagctgctggaggtgctgcaggaga
gggagcttggagcctcttacaccagtgagaaactccaaccaagctctctgaggatctgaaggaaactgaattcaaaaagatcaaagtgtg
ggctccggtgcttccgacgggtgtataaggactctggatcccagaaggtgagaaagttaaattcccgtcgtatcaagacatctccgaaagccaa
caaggaaatcctcagatgaagcctacgtgatggccagcgtggacaacccccacgtgctgcccgtgctgggcatctgctcactccacctgcaactca
tcagtcagctcatgccctcggctgctcctggactatgtccgggaacacaaagacaatattggctcccagtacctgctcaactgggtgtgctcagatgc
aaaggcatgaactacttgaggacctgctgctggtgaccgcgacctggcggccaggaaactgactggtgaaaacaccgcagatgtcaagatcac
agatttgggctggccaaactgctgggtgcggaagagaaagaataccatgcagaaggaggcaagtgctatcaagtggatggcattggaatcaatt
ctacacagaatctatacccaccagagtgatgtctggagctacggggtgaccttgggagttgatgaccttggatccaagccatgatgacggaatcct
gccagcagatctcctcatcctggagaaaggagaacgcctcctcagccaccatattgaccatcagatgtctacatgatcatggtcaagtgtggatg
atagacgcagatagtcgcccaggtccgtgagttgatcatcgaattctccaaaatggcccagagacccccagcgtaccttgcattcagggggatga
aagaatgcatttgcgaagtctacagactccaacttctaccgtgccctgatggatgaagaagacatggacgacgtggtggatgccgacgagtagctca
tcccacagcagggtcttctcagcagcccctccacgtcacggactcccctcctgagctctctgagtgcaaccagcaacaattccaccgtggcttgcattga
tagaaatgggctgcaaagctgtcccataaggaagacagcttcttgcagcgtacagctcagacccccagggccttgcagtaggacagcatagac
gacaccttctcccagtgctgaatacataaaccagtccgttccaaaaggcccgtggctctgtgcaaatcctgtctatcacaatcagcctctgaac
cccgcgccagcagagaccacactaccaggacccccacagcactgcagtgggcaacccccgagtatctcaactgtccagcccactgtgtcaaca
gcacattcgacagccctgcccactgggcccagaaaggcagccaccaaattagcctggacaacctgactaccagcaggacttcttcccaaggaagc
caagccaaatggcatcttaagggtccacagctgaaaatgcagaatacctaagggtcgcgcccaaaagcagtgaaattattggagca