

>EGFR T790M

atcgaccctccgggacggccgggagcgcctcctggcgtgctggctgctgctctgcccggcgagtcgggctctggaggaaaagaaagtgtccaag
gcacgagtaacaagctcacgcagttgggcacttttgaagatcattttctcagcctccagaggatgttcaataactgtgaggtggtccttgggaatttgga
aattacctatgtgcagaggaattatgatctttccttctaaagaccatccaggaggtggctggttatgtcctcattgcctcaacacagtgagcgaatt
cctttggaaaacctgcagatcatcagaggaaatagtactacgaaaattcctatgccttagcagcttatctaactatgatgcaataaaaacgggactg
aaggagctgcccagagaaatctacaggaaatcctgcatggcgcgctgcggttcagcaacaacctgcctgtgcaacgtggagagcatccagtggc
gggacatagtcagcagtgactttctcagcaaatgctgatggacttccagaacctgggagcagctgcaaaaagtgtgatccaagctgtccaatggg
agctgctggggtgcaggagaggagaactgccagaaactgacaaaatcatctgtgccagcagtgctccgggctgctgctggcaagtccccagtg
actgctccacaaccagtgtgctgcaggctgcacagggccccgggagagcagctgctggctgcccgaattccgagacgaagccacgtgcaagga
cacctgccccactcatgcttcaacccaccacgtaccagatggatgtgaacccgaggggcaatacagctttgggtgccacctgctgaagaagt
gtccccgtaattatgtggtgacagatcacggctcgtgctccagcctgtggggccgacagctatgagatggaggaagacggcgtccgcaagtgtaa
gaagtgcgaaggccttccgcaaaagtgtgaacggaataggtattggtgaatttaaagactcactctcataaatgctacgaatattaacacttca
aaaactgcacctccatcagtggtgcatctccacatcctgcccgtggcatttaggggtgactccttcacacatactcctcctctggatccacaggaactgg
atattctgaaaaccgtaaaggaaatcacagggttttctgattcaggcttggcctgaaaacaggacggacctccatgcctttgagaacctgaaatc
atcgcggcaggaccaagcaacatggtcagtttctctgagctcgtcagcctgaacataacatccttgggattacgctccctcaaggagataagtat
ggagatgtgataattcaggaacaaaaattgtgctatgcaatacaataaactgaaaaaactgtttgggacctccggtcagaaaacaaaaattat
aagcaacagaggtgaaaacagctgcaaggccacaggccaggtctgcatgcttctgctccccgagggtgctgctggggccggagcccaggggactg
cgtctcttccggaatgtcagccgaggcagggaatgctggacaagtcaaccttctggagggtgagccaaggaggttgggagaactctgagtg
atcagtgccaccagagtgctgctcaggccatgaacatcacctgcacaggacggggaccagacaactgtatccagtggtcccactacattgacg
gccccactgctcaagacctgcccggcaggagtcagggagaaaacaacacctggtctggaagtacgcagacgcccagccatgtgtgccacctgtg
ccatcaaaactgcacctacggatgactgggacaggtctgaaggctgccaacgaatgggctcaaaatcccgtccatgcacctgggatggtggggg
cctcctcttctgctgctggtggtggccctggggatcgccctctcatgcaaggcgcacatcgttcggaagcgcacgctgaggaggtgctgcaggaga
gggagcttggagcctcttacaccagtgagaaactccaaccaagctctcttggagatctgaaggaaactgaattcaaaaagatcaaatgtctg
ggctccggtgcttggcagcgggtgataagggactctggatcccagaaggtgagaaagttaaattcccgtcgtatcaaggaattaagagaagcaa
catctcgaagcaacaaggaaatcctcgatgaagcctacgtgatggcagcgtggacaacccccacgtgtgcccctgctgggcatctgctcacc
tccacctgcaactcatcagcagctcatgcccttggctgctcctggactatgcccgggaacaaaagacaatattggctcccagctacgtctcaact
ggtgtgtcagatcgaaggatgaactactggaggaccgtcgttgggtgaccgagacctggcggccaggaactactggtgaaaaaccgca
gcatgtcaagatcacagatttgggctggcaaaactgctgggtgcggaagagaaagaataccatgcagaaggaggcaaatgctctatcaagtggat
ggcattggaatcaattctacagaaatctatacccaccagagtgatgtctggagctacggggtgaccttgggagttgatgacctttggatccaagcc
atatgacggaatccctgccagcgagatctcctcatcctggagaaaggagaacgcctccctcagccacctatgtaccatcgatgtctacatgatcat
ggtcaagtgtggatgatagacgcagatagtcgcccgaagttccgtgagttgatcatcgaattctcaaaatggcccagacccccagcgtacctgt
cattcagggggatgaaagaatgcaatttccaagtctacagactccaacttctaccgtgccctgatggatgaagaagacatggacgacgtggtggatg
ccgacgagtacctcatcccacagcagggtcttctcagcagcccctccacgtcacggactcccctcctgagctctctgagtgcaaccagcaacaattcca
ccgtggcttgattgatagaatgggctgcaaaactgtccatcaaggaagacagcttctgacgatacagctcagacccccacaggcgccttgact
gaggacagcatagacgacaccttctccagtcctgaatacataaacagtccttccaaaaggccgctggctctgtgcaaatcctgtctatcac
aatcagcctctgaaccccgcccagcagacccacactaccaggacccccacagcactgcagtgggcaacccccagatctcaacactgtccagc
ccacctgttcaacagcattcgcagccctgcccactgggcccagaaggcagccaccaaattagcctggacaacctgactaccagcaggactt
ctttccaaggaaagcaagcaaatggcatctttaaagggtccacagctgaaaatgcagaatacctaagggtcgcccaaaagcagtgatatttattg
gagca