

> EGFR (Y764\_V765insHH)

ATGCGACCCTCCGGGACGGCCGGGGCAGCGCTCCTGGCGCTGCTGGCTGCGCTCTGCCCGGCGAGTCGGGCTCT  
GGAGGAAAAGAAAGTTTTGCCAAGGCACGAGTAACAAGCTCACGCAGTTGGGCACTTTTGAAGATCATTTTCTCAG  
CCTCCAGAGGATGTTCAATAACTGTGAGGTGGTCTTGGGAATTTGGAAATTACCTATGTGCAGAGGAATTATGAT  
CTTTCCTTCTTAAAGACCATCCAGGAGGTGGCTGGTTATGTCCTCATTGCCCTCAACACAGTGGAGCGAATTCCTT  
GGAAAACCTGCAGATCATCAGAGGAAATATGTACTACGAAAATTCCTATGCCTTAGCAGTCTTATCTAACTATGAT  
GCAAATAAAACCGGACTGAAGGAGCTGCCATGAGAAATTTACAGGAAATCCTGCATGGCGCCGTGCGGTTCCAGC  
AACAACCTGCCCTGTGCAACGTGGAGAGCATCCAGTGGCGGGACATAGTCAGCAGTGACTTTCTCAGCAACATG  
TCGATGGACTTCCAGAACCACCTGGGCAGCTGCCAAAAGTGTGATCCAAGCTGTCCAATGGGAGCTGCTGGGGT  
GCAGGAGAGGAGAACTGCCAGAACTGACCAAATCATCTGTGCCAGCAGTGTCCGGGCGCTGCCGTGGCAA  
GTCCCCAGTGACTGCTGCCACAACCAGTGTGCTGCAGGCTGCACAGGCCCGGGAGAGCGACTGCCTGGTCTG  
CCGCAAATTCGAGACGAAGCCACGTGCAAGGACACCTGCCCCCACTCATGCTCTACAACCCACCACGTACCAG  
ATGGATGTGAACCCCGAGGGCAAATACAGCTTTGGTGCCACCTGCGTGAAGAAGTGTCCCCGTAATTATGTGGTG  
ACAGATCACGGCTCGTGCGTCCGAGCCTGTGGGGCCGACAGCTATGAGATGGAGGAAGACGGCGTCCGCAAGTG  
TAAGAAGTGCGAAGGGCCTTGCCGAAAAGTGTGTAACGGAATAGGTATTGGTGAATTTAAAGACTCACTCTCCAT  
AAATGCTACGAATATTAACACTTCAAAAACCTGCACCTCCATCAGTGGCGATCTCCACATCCTGCCGGTGGCATTTA  
GGGGTGACTCCTTCACACATACTCCTCCTGTGATCCACAGGAACTGGATATTCTGAAAACCGTAAAGGAAATCAC  
AGGGTTTTTGTGATTGAGGCTTGGCCTGAAAACAGGACGGACCTCCATGCCTTTGAGAACCTAGAAATCATACGC  
GGCAGGACCAAGCAACATGGTCAGTTTTCTTGCAGTCGTCAGCCTGAACATAACATCCTTGGGATTACGCTCCC  
TCAAGGAGATAAGTGATGGAGATGTGATAATTTAGGAAACAAAATTTGTGCTATGCAAATACAATAAACTGGA  
AAAAACTGTTTGGGACCTCCGGTCAGAAAACCAAATATAAGCAACAGAGGTGAAAACAGCTGCAAGGCCACA  
GGCCAGGTCTGCCATGCCTTGTGCTCCCCGAGGGCTGCTGGGGCCCGGAGCCCAGGGACTGCGTCTCTTGGCCG  
AATGTCAGCCGAGGCAGGGAATGCGTGGACAAGTGAACCTTCTGGAGGGTGAAGCAAGGGAGTTTGTGGAGA  
ACTCTGAGTGCATACAGTGCCACCCAGAGTGCCTGCCTCAGGCCATGAACATCACCTGCACAGGACGGGGACCAG  
ACAAGTGTATCCAGTGTGCCCACTACATTGACGGCCCCCACTGCGTCAAGACCTGCCCGGACAGGAGTCATGGGAG  
AAAACAACACCCCTGGTCTGGAAGTACGCAGACGCCGCCATGTGTGCCACCTGTGCCATCCAACTGCACCTACG  
GATGCACTGGGCCAGGTCTTGAAGGCTGTCCAACGAATGGGCCTAAGATCCCGTCCATCGCCACTGGGATGGTGG  
GGGCCCTCCTCTTGTGCTGGTGGTGGCCCTGGGGATCGGCCTCTTATGCGAAGGCGCCACATCGTTCGGAAGC  
GCACGCTGCGGAGGCTGCTGCAGGAGAGGGAGCTTGTGGAGCCTTTACACCCAGTGGAGAAGCTCCCAACCAA  
GCTCTCTTGAAGGATCTTGAAGGAACTGAATCAAAAAGATCAAAGTGTGGGCTCCGGTGCCTTCGGCACGGTG  
TATAAGGGACTCTGGATCCCAGAAGGTGAGAAAAGTTAAATTTCCCGTCGCTATCAAGGAATTAAGAGAAGCAACA  
TCTCCGAAAGCCAACAAGGAAATCCTCGATGAAGCCTACACCACTGATGGCCAGCGTGGACAACCCCCACGTG  
TGCCGCCTGCTGGGCATCTGCCTCACCTCCACCGTGCAACTCATCACGCAGCTCATGCCCTTCGGCTGCCTCCTGGA  
CTATGTCCGGGAACACAAAGACAATATTGGCTCCAGTACCTGCTCAACTGGTGTGTGCAGATCGCAAAGGGCAT  
GAACTACTTGGAGACCGTCTGCTTGGTGCACCGCGACCTGGCAGCCAGGAACGTACTGGTAAAACACCCGACG  
ATGTCAAGATCACAGATTTTGGGCTGGCCAACTGCTGGGTGCGGAAGAGAAAGAATACCATGCAGAAGGAGGC  
AAAGTGCCTATCAAGTGGATGGCATTGGAATCAATTTTACACAGAATCTATAACCCACCAGAGTGATGCTGGAGCT  
ACGGGGTACTGTTTGGGAGTTGATGACCTTTGGATCCAAGCCATATGACGGAATCCCTGCCAGCGAGATCTCCTC  
CATCCTGGAGAAAGGAGAACGCCTCCCTCAGCCACCCATATGTACCATCGATGTCTACATGATCATGGTCAAGTGC  
TGGATGATAGACGCAGATAGTCGCCAAAGTTCCGTGAGTTGATCATCGAATTCTCCAAAATGGCCCGAGACCCCC  
AGCGCTACCTTGTATTAGGGGGATGAAAGAATGCATTTGCCAAGTCTACAGACTCCAACTTCTACCGTGCCCT  
GATGGATGAAGAAGACATGGACGACGTGGTGGATGCCGACGAGTACCTCATCCACAGCAGGGCTTCTTACGCA  
GCCCCCTCACGTACGGACTCCCCCTGAGCTCTCTGAGTGCAACCAGCAACAATTCACCGTGGCTTGCATTGAT  
AGAAATGGGCTGCAAAGCTGTCCATCAAGGAAGACAGCTTCTTGCAGCGATACAGCTCAGACCCCCACAGGCGCC

TTGACTGAGGACAGCATAGACGACACCTTCCTCCCAGTGCCTGAATACATAAACCAGTCCGTTCCAAAAGGCCCG  
CTGGCTCTGTGCAGAATCCTGTCTATCACAAATCAGCCTCTGAACCCCGCACCCAGCAGAGACCCACACTACCAGGA  
CCCCACAGCACTGCAGTGGGCAACCCCGAGTATCTCAACTGTCCAGCCCACCTGTGTCAACAGCACATTGAC  
AGCCCTGCCCACTGGGCCAGAAAGGCAGCCACCAAATTAGCCTGGACAACCCTGACTACCAGCAGGACTTCTTTC  
CCAAGGAAGCCAAGCCAAATGGCATCTTTAAGGGCTCCACAGCTGAAAATGCAGAATACCTAAGGGTCGCGCCAC  
AAAGCAGTGAATTTATTGGAGCA