

>EGFR M766\_A767insA1

ATGCGACCCTCCGGGACGGCCGGGCAGCGCTCCTGGCGCTGCTGGCTGCGCTCTGCCCGGCGAGTCGGGCTCT  
GGAGGAAAAGAAAGTTTTGCCAAGGCACGAGTAACAAGCTCACGCAGTTGGGCACTTTTGAAGATCATTCTCAG  
CCTCCAGAGGATGTTCAATAACTGTGAGGTGGTCTTGGGAATTTGGAAATTACCTATGTGCAGAGGAATTATGAT  
CTTTCCTTCTTAAAGACCATCCAGGAGGTGGCTGGTTATGTCCTCATTGCCCTCAACACAGTGGAGCGAATTCCTT  
GGAAAACCTGCAGATCATCAGAGGAAATATGTACTACGAAAATTCCTATGCCTTAGCAGTCTTATCTAACTATGAT  
GCAAATAAAACCGGACTGAAGGAGCTGCCATGAGAAATTTACAGGAAATCCTGCATGGCGCCGTGCGGTTCCAGC  
AACAACCTGCCCTGTGCAACGTGGAGAGCATCCAGTGGCGGGACATAGTCAGCAGTGACTTTCTCAGCAACATG  
TCGATGGACTTCCAGAACCACCTGGGCAGCTGCCAAAAGTGTGATCCAAGCTGTCCAATGGGAGCTGCTGGGGT  
GCAGGAGAGGAGAACTGCCAGAACTGACCAAATCATCTGTGCCAGCAGTGTCCGGGCGCTGCCGTGGCAA  
GTCCCCAGTGACTGCTGCCACAACCAGTGTGCTGCAGGCTGCACAGGCCCCCGGGAGAGCGACTGCCTGGTCTG  
CCGCAAATTCGAGACGAAGCCACGTGCAAGGACACCTGCCCCCACTCATGCTCTACAACCCACCACGTACCAG  
ATGGATGTGAACCCCGAGGGCAAATACAGCTTTGGTGCCACCTGCGTGAAGAAGTGTCCCCGTAATTATGTGGTG  
ACAGATCACGGCTCGTGCGTCCGAGCCTGTGGGGCCGACAGCTATGAGATGGAGGAAGACGGCGTCCGCAAGTG  
TAAGAAGTGCGAAGGGCCTTGCCGAAAAGTGTGAACGGAATAGGTATTGGTGAATTTAAAGACTCACTCTCCAT  
AAATGCTACGAATATTAACACTTCAAAAACCTGCACCTCCATCAGTGGCGATCTCCACATCCTGCCGGTGGCATTTA  
GGGGTGACTCCTTCACACATACTCCTCCTGTGATCCACAGGAACTGGATATTCTGAAAACCGTAAAGGAAATCAC  
AGGGTTTTTGTGATTGAGGCTTGGCCTGAAAACAGGACGGACCTCCATGCCTTTGAGAACCTAGAAATCATACGC  
GGCAGGACCAAGCAACATGGTCAGTTTTCTTGCAGTCGTCAGCCTGAACATAACATCCTTGGGATTACGCTCCC  
TCAAGGAGATAAGTGATGGAGATGTGATAATTTAGGAAACAAAATTTGTGCTATGCAAATACAATAAACTGGA  
AAAAACTGTTTGGGACCTCCGGTCAGAAAACCAAATATAAGCAACAGAGGTGAAAACAGCTGCAAGGCCACA  
GGCCAGGTCTGCCATGCCTGTGCTCCCCGAGGGCTGCTGGGGCCCGGAGCCCAGGGACTGCGTCTCTTGGCCG  
AATGTCAGCCGAGGCAGGGAATGCGTGGACAAGTGAACCTTCTGGAGGGTGAGCCAAGGGAGTTTGTGGAGA  
ACTCTGAGTGCATACAGTGCCACCCAGAGTGCCTGCCTCAGGCCATGAACATCACCTGCACAGGACGGGGACCAG  
ACAACGTATCCAGTGTGCCCACTACATTGACGGCCCCCACTGCGTCAAGACCTGCCCGGACAGGAGTCATGGGAG  
AAAACAACACCCCTGGTCTGGAAGTACGCAGACGCCGCCATGTGTGCCACCTGTGCCATCCAACTGCACCTACG  
GATGCACTGGGCCAGGTCTTGAAGGCTGTCCAACGAATGGGCCTAAGATCCCGTCCATCGCCACTGGGATGGTGG  
GGGCCCTCCTCTTGTGCTGGTGGTGGCCCTGGGGATCGGCCTCTTATGCGAAGGCGCCACATCGTTCGGAAGC  
GCACGCTGCGGAGGCTGCTGCAGGAGAGGGAGCTTGTGGAGCCTTTACACCCAGTGGAGAAGCTCCCAACCAA  
GCTCTCTTGAAGGATCTTGAAGGAACTGAATCAAAAAGATCAAAGTGTGGGCTCCGGTGCCTTCGGCACGGTG  
TATAAGGGACTCTGGATCCCAGAAGGTGAGAAAAGTTAAATTCCTGTCGCTATCAAGGAATTAAGAGAAGCAACA  
TCTCCGAAAGCCAACAAGGAAATCCTCGATGAAGCCTACGTGATG**GCCATA**GCCAGCGTGGACAACCCCCACGTG  
TGCCGCCTGCTGGGCATCTGCCTCACCTCCACCGTGCAACTCATCACGCAGCTCATGCCCTTCGGTGCCTCCTGGA  
CTATGTCCGGGAACACAAAGACAATATTGGCTCCAGTACCTGCTCAACTGGTGTGTGCAGATCGCAAAGGGCAT  
GAACTACTTGGAGACCGTCTGCTTGGTGCACCGCGACCTGGCAGCCAGGAACGTAAGTGGTAAAACACCCGACG  
ATGTCAAGATCACAGATTTTGGGCTGGCCAACTGCTGGGTGCGGAAGAGAAAGAATACCATGCAGAAGGAGGC  
AAAGTGCCTATCAAGTGGATGGCATTGGAATCAATTTACACAGAATCTATAACCCACCAGAGTGATGCTGGAGCT  
ACGGGGTACTGTTTGGGAGTTGATGACCTTTGGATCCAAGCCATATGACGGAATCCCTGCCAGCGAGATCTCCTC  
CATCCTGGAGAAAGGAGAACGCCTCCCTCAGCCACCCATATGTACCATCGATGTCTACATGATCATGGTCAAGTGC  
TGGATGATAGACGAGATAGTCGCCAAAGTTCCGTGAGTTGATCATCGAATTCTCCAAAATGGCCCGAGACCCCC  
AGCGCTACCTGTCAATCAGGGGGATGAAAGAATGCATTTGCCAAGTCTACAGACTCCAACTTCTACCGTGCCCT  
GATGGATGAAGAAGACATGGACGACGTGGTGGATGCCGACGAGTACCTCATCCACAGCAGGGCTTCTTACGCA  
GCCCCCTCACGTACGGACTCCCCCTGAGCTCTCTGAGTGCAACCAGCAACAATCCACCGTGGCTTGCATTGAT  
AGAAATGGGCTGCAAAGCTGTCCATCAAGGAAGACAGCTTCTTGCAGCGATACAGCTCAGACCCCCACAGGCGCC

TTGACTGAGGACAGCATAGACGACACCTTCCTCCCAGTGCCTGAATACATAAACCCAGTCCGTTCCAAAAGGCCCG  
CTGGCTCTGTGCAGAATCCTGTCTATCACAAATCAGCCTCTGAACCCCGCACCCAGCAGAGACCCACACTACCAGGA  
CCCCACAGCACTGCAGTGGGCAACCCCGAGTATCTCAACTGTCCAGCCCACCTGTGTCAACAGCACATTGAC  
AGCCCTGCCCACTGGGCCAGAAAGGCAGCCACCAAATTAGCCTGGACAACCCTGACTACCAGCAGGACTTCTTTC  
CCAAGGAAGCCAAGCCAAATGGCATCTTTAAGGGCTCCACAGCTGAAAATGCAGAATACCTAAGGGTCGCGCCAC  
AAAGCAGTGAATTTATTGGAGCA