

>EGFR (A767\_dupASV)

ATGCGACCCTCCGGGACGGCCGGGCAGCGCTCCTGGCGCTGCTGGCTGCGCTCTGCCCGGCGAGTCGGGCTCT  
GGAGGAAAAGAAAAGTTTGGCAAGGCACGAGTAACAAGCTCACGCAGTTGGGCACTTTTGAAGATCATTTTCTCAG  
CCTCCAGAGGATGTTCAATAACTGTGAGGTGGTCTTGGGAATTTGGAAATTACCTATGTGCAGAGGAATTATGAT  
CTTTCCTTCTTAAAGACCATCCAGGAGGTGGCTGGTTATGTCCTCATTGCCCTCAACACAGTGGAGCGAATTCCTT  
GGAAAACCTGCAGATCATCAGAGGAAATATGTACTACGAAAATTCCTATGCCTTAGCAGTCTTATCTAACTATGAT  
GCAAATAAAACCGGACTGAAGGAGCTGCCATGAGAAATTTACAGGAAATCCTGCATGGCGCCGTGCGGTTCCAGC  
AACAACCTGCCCTGTGCAACGTGGAGAGCATCCAGTGGCGGGACATAGTCAGCAGTGACTTTCTCAGCAACATG  
TCGATGGACTTCCAGAACCACCTGGGCAGCTGCCAAAAGTGTGATCCAAGCTGTCCAATGGGAGCTGCTGGGGT  
GCAGGAGAGGAGAACTGCCAGAACTGACCAAATCATCTGTGCCAGCAGTGTCCGGGCGCTGCCGTGGCAA  
GTCCCCAGTGACTGCTGCCACAACCAGTGTGCTGCAGGCTGCACAGGCCCCCGGGAGAGCGACTGCCTGGTCTG  
CCGCAAATTCGAGACGAAGCCACGTGCAAGGACACCTGCCCCCACTCATGCTCTACAACCCACCACGTACCAG  
ATGGATGTGAACCCCGAGGGCAAATACAGCTTTGGTGCCACCTGCGTGAAGAAGTGTCCCCGTAATTATGTGGTG  
ACAGATCACGGCTCGTGCGTCCGAGCCTGTGGGGCCGACAGCTATGAGATGGAGGAAGACGGCGTCCGCAAGTG  
TAAGAAGTGCGAAGGGCCTTGCCGAAAAGTGTGAACGGAATAGGTATTGGTGAATTTAAAGACTCACTCTCCAT  
AAATGCTACGAATATTAACACTTCAAAAATGCACCTCCATCAGTGGCGATCTCCACATCCTGCCGGTGGCATTTA  
GGGGTGACTCCTTCACACATACTCCTCCTTGATCCACAGGAACTGGATATTCTGAAAACCGTAAAGGAAATCAC  
AGGGTTTTTGTGATTGAGGCTTGGCCTGAAAACAGGACGGACCTCCATGCCTTTGAGAACCTAGAAATCATACGC  
GGCAGGACCAAGCAACATGGTCAGTTTTCTTGCAGTCGTCAGCCTGAACATAACATCCTTGGGATTACGCTCCC  
TCAAGGAGATAAGTGATGGAGATGTGATAATTTAGGAAACAAAATTTGTGCTATGCAAATACAATAAACTGGA  
AAAAACTGTTTGGGACCTCCGTCAGAAAACCAAATATAAGCAACAGAGGTGAAAACAGCTGCAAGGCCACA  
GGCCAGGTCTGCCATGCCTTGTGCTCCCCGAGGGCTGCTGGGGCCCGGAGCCCAGGGACTGCGTCTCTTGGCCG  
AATGTCAGCCGAGGCAGGGAATGCGTGGACAAGTGAACCTTCTGGAGGGTGAGCCAAGGGAGTTTGTGGAGA  
ACTCTGAGTGCATACAGTGCCACCCAGAGTGCCTGCCTCAGGCCATGAACATCACCTGCACAGGACGGGGACCAG  
ACAAGTGTATCCAGTGTGCCACTACATTGACGGCCCCACTGCGTCAAGACCTGCCCGGAGGAGTATGGGAG  
AAAACAACACCCTGGTCTGGAAGTACGCAGACGCCGCCATGTGTGCCACCTGTGCCATCCAACTGCACCTACG  
GATGCACTGGGCCAGGTCTTGAAGGCTGTCCAACGAATGGGCCTAAGATCCCGTCCATCGCCACTGGGATGGTG  
GGGCCCTCCTCTTGTGCTGGTGGTGGCCCTGGGGATCGGCCTCTTATGCGAAGGCGCCACATCGTTCGGAAGC  
GCACGCTGCGGAGGCTGCTGCAGGAGAGGGAGCTTGTGGAGCCTTTACACCCAGTGGAGAAGCTCCCAACCA  
GCTCTCTTGAAGGATCTTGAAGGAACTGAATCAAAAAGATCAAAGTGTGGGCTCCGGTGCCTTGGGCACGGTG  
TATAAGGGACTCTGGATCCCAGAAGGTGAGAAAAGTTAAATTTCCCGTCGCTATCAAGGAATTAAGAGAAGCAACA  
TCTCCGAAAGCCAACAAGGAAATCCTCGATGAAGCCTACGTGATGGCCAGCGTGGCCAGCGTGGACAACCCCCAC  
GTGTGCCGCTGCTGGGCATCTGCCTCACCTCCACCGTGCAGCTCATCACGCAGCTCATGCCCTTCGGTGCCTCCT  
GGACTATGTCCGGGAACACAAAGACAATATTGGCTCCAGTACCTGCTCAACTGGTGTGTGCAGATCGCAAAGGG  
CATGAACTACTTGGAGACCGTGCCTTGGTGCACCCGACCTGGCAGCCAGGAACGTAAGTGGTGAACACCCGCA  
GCATGTCAAGATCACAGATTTTGGGCTGGCCAACTGCTGGGTGCGGAAGAGAAAGAATACCATGCAGAAGGAG  
GCAAAGTGCCTATCAAGTGGATGGCATTGGAATCAATTTACACAGAATCTATAACCCACCAGAGTGTGTCTGGAG  
CTACGGGGTACTGTTTGGGAGTTGATGACCTTTGGATCCAAGCCATATGACGGAATCCCTGCCAGCGAGATCTCC  
TCCATCCTGGAGAAAGGAGAACGCCTCCCTCAGCCACCCATATGTACCATCGATGTCTACATGATCATGGTCAAGT  
GCTGGATGATAGACGCAGATAGTCGCCAAAAGTTCCGTGAGTTGATCATCGAATTCCTCAAAAATGGCCGAGACC  
CCCAGCGCTACCTTGTCAATCAGGGGGATGAAAGAATGCATTTGCCAAGTCTACAGACTCCAATTCTACCGTGC  
CCTGATGGATGAAGAAGACATGGACGACGTGGTGGATGCCGACGAGTACCTCATCCCACAGCAGGGCTTCTCAG  
CAGCCCCTCCACGTACGGACTCCCCTCCTGAGCTCTCTGAGTGCAACCAGCAACAATTCACCGTGGCTTGCATTG  
ATAGAAATGGGCTGCAAAGCTGTCCATCAAGGAAGACAGCTTCTTGCAGCGATACAGCTCAGACCCACAGGCC

CCTTGACTGAGGACAGCATAGACGACACCTTCCTCCCAGTGCCTGAATACATAAACCAGTCCGTTCCCAAAGGCC  
CGCTGGCTCTGTGCAGAATCCTGTCTATCACAAATCAGCCTCTGAACCCCGCACCCAGCAGAGACCCACACTACCAG  
GACCCCAACAGCACTGCAGTGGGCAACCCCGAGTATCTCAAACTGTCCAGCCACCTGTGTCAACAGCACATTCG  
ACAGCCCTGCCCACTGGGCCCAGAAAGGCAGCCACCAAATTAGCCTGGACAACCCTGACTACCAGCAGGACTTCTT  
TCCAAGGAAGCCAAGCCAAATGGCATCTTTAAGGGCTCCACAGCTGAAAATGCAGAATACCTAAGGGTCGCGCC  
ACAAAGCAGTGAATTTATTGGAGCA