

>EGFR A763_Y764insFQEA

ATGCGACCCTCCGGGACGGCCGGGGCAGCGCTCCTGGCGCTGCTGGCTGCGCTCTGCCCGGCGAGTCGGGCTCT
GGAGGAAAAGAAAGTTTGGCAAGGCACGAGTAACAAGCTCACGCAGTTGGGCACTTTTGAAGATCATTTTCTCAG
CCTCCAGAGGATGTTCAATAACTGTGAGGTGGTCTTGGGAATTTGGAAATTACCTATGTGCAGAGGAATTATGAT
CTTTCCTTCTTAAAGACCATCCAGGAGGTGGCTGGTTATGTCCTCATTGCCCTCAACACAGTGGAGCGAATTCCTT
GGAAAACCTGCAGATCATCAGAGGAAATATGTACTACGAAAATTCCTATGCCTTAGCAGTCTTATCTAACTATGAT
GCAAATAAAACCGGACTGAAGGAGCTGCCCATGAGAAATTTACAGGAAATCCTGCATGGCGCCGTGCGGTTCCAGC
AACAACCTGCCCTGTGCAACGTGGAGAGCATCCAGTGGCGGGACATAGTCAGCAGTGACTTTCTCAGCAACATG
TCGATGGACTTCCAGAACCACCTGGGCAGCTGCCAAAAGTGTGATCCAAGCTGTCCAATGGGAGCTGCTGGGGT
GCAGGAGAGGAGAACTGCCAGAACTGACCAAATCATCTGTGCCAGCAGTGTCCGGGCGCTGCCGTGGCAA
GTCCCCAGTGACTGCTGCCACAACCAGTGTGCTGCAGGCTGCACAGGCCCCCGGGAGAGCGACTGCCTGGTCTG
CCGCAAATTCGAGACGAAGCCACGTGCAAGGACACCTGCCCCCACTCATGCTCTACAACCCACCACGTACCAG
ATGGATGTGAACCCCGAGGGCAAATACAGCTTTGGTGCCACCTGCGTGAAGAAGTGTCCCCGTAATTATGTGGTG
ACAGATCACGGCTCGTGCGTCCGAGCCTGTGGGGCCGACAGCTATGAGATGGAGGAAGACGGCGTCCGCAAGTG
TAAGAAGTGCGAAGGGCCTTGCCGAAAAGTGTGTAACGGAATAGGTATTGGTGAATTTAAAGACTCACTCTCCAT
AAATGCTACGAATATTAACACTTCAAAAACCTGCACCTCCATCAGTGGCGATCTCCACATCCTGCCGGTGGCATTTA
GGGGTGACTCCTTCACACATACTCCTCCTCTGGATCCACAGGAACTGGATATTCTGAAAACCGTAAAGGAAATCAC
AGGGTTTTTGTGATTGAGGCTTGGCCTGAAAACAGGACGGACCTCCATGCCTTTGAGAACCTAGAAATCATACGC
GGCAGGACCAAGCAACATGGTCAGTTTTCTTGCAGTCGTCAGCCTGAACATAACATCCTTGGGATTACGCTCCC
TCAAGGAGATAAGTGATGGAGATGTGATAATTTAGGAAACAAAATTTGTGCTATGCAAATACAATAAACTGGA
AAAAACTGTTTGGGACCTCCGGTCAGAAAACCAAATATAAGCAACAGAGGTGAAAACAGCTGCAAGGCCACA
GGCCAGGTCTGCCATGCCTTGTGCTCCCCGAGGGCTGCTGGGGCCCGGAGCCCAGGGACTGCGTCTCTTGGCCG
AATGTCAGCCGAGGCAGGGAATGCGTGGACAAGTGAACCTTCTGGAGGGTGAAGCAAGGGAGTTTGTGGAGA
ACTCTGAGTGCATACAGTGCCACCCAGAGTGCCTGCCTCAGGCCATGAACATCACCTGCACAGGACGGGGACCAG
ACAAGTGTATCCAGTGTGCCCACTACATTGACGGCCCCCACTGCGTCAAGACCTGCCCGGACAGGAGTCATGGGAG
AAAACAACACCCCTGGTCTGGAAGTACGCAGACGCCGCCATGTGTGCCACCTGTGCCATCCAACTGCACCTACG
GATGCACTGGGCCAGGTCTTGAAGGCTGTCCAACGAATGGGCCTAAGATCCCCTCCATCGCCACTGGGATGGTG
GGGCCCTCCTCTTGTGCTGGTGGTGGCCCTGGGGATCGGCCTCTTATGCGAAGGCGCCACATCGTTCGGAAGC
GCACGCTGCGGAGGCTGCTGCAGGAGAGGGAGCTTGTGGAGCCTTTACACCCAGTGGAGAAGCTCCCAACCAA
GCTCTCTTGAAGGATCTTGAAGGAACTGAATCAAAAAGATCAAAGTGTGGGCTCCGGTGCCTTGGGCACGGTG
TATAAGGGACTCTGGATCCCAGAAGGTGAGAAAAGTAAAATCCCCTCGCTATCAAGGAATTAAGAGAAGCAACA
TCTCCGAAAGCCAACAAGGAATCCTCGATGAAGCCTTCCAGGAAGCCTACGTGATGGCCAGCGGCATCTGCCTC
ACCTCCACCGTGCAGCTCATCACGCAGCTCATGCCCTTCGGCTGCCTCCTGGACTATGTCCGGGAACACAAAGACA
ATATTGGCTCCAGTACCTGCTCAACTGGTGTGTGCAGATCGCAAAGGGCATGAACTACTTGGAGGACCGTGCCTT
GGTGACCCGCGACCTGGCAGCCAGGAACGTACTGGTAAAACACCCGACGATGTCAAGATCACAGATTTTGGGCT
GGCAAACCTGCTGGGTGCGGAAGAGAAAGAATACCATGCAGAAGGAGCAAAGTGCCTATCAAGTGGATGGCAT
TGGAATCAATTTTACACAGAATCTATAACCACAGAGTGTGCTGGAGCTACGGGGTACTGTTTGGGAGTTGAT
GACCTTTGGATCCAAGCCATATGACGGAATCCCTGCCAGCGAGATCTCCTCCATCCTGGAGAAAGGAGAACGCCT
CCCTCAGCCACCCATATGTACCATCGATGTCTACATGATCATGGTCAAGTGTGGATGATAGACGCAGATAGTCGC
CCAAAGTTCGGTGAAGTTCATCGAATTTCCAAAATGGCCCAGACCCCCAGCGCTACCTTGTCAATTCAGGGGG
ATGAAAGAATGCATTTGCCAAGTCTACAGACTCCAATTCTACCGTGCCTGATGGATGAAGAAGACATGGACG
ACGTGGTGGATGCCGACGAGTACCTCATCCCACAGCAGGGCTTCTTACAGAGCCCCTCCACGTACGGACTCCCCT
CCTGAGCTCTCTGAGTGAACCCAGCAACAATCCACCGTGGCTTGCATTGATAGAAATGGGCTGCAAAGCTGTCCC
ATCAAGGAAGACAGCTTCTTGCAGCGATACAGCTCAGACCCACAGGGCGCCTTACTGAGGACAGCATAGACGAC

ACCTTCCTCCCAGTGCCTGAATACATAAACCAGTCCGTTCCCAAAGGCCCGCTGGCTCTGTGCAGAATCCTGTCTA
TCACAATCAGCCTCTGAACCCCGCACCCAGCAGAGACCCACACTACCAGGACCCCCACAGCACTGCAGTGGGCAA
CCCCGAGTATCTCAACACTGTCCAGCCCACCTGTGTCAACAGCACATTCGACAGCCCTGCCCACTGGGCCAGAAA
GGCAGCCACCAAATTAGCCTGGACAACCCTGACTACCAGCAGGACTTCTTTCCAAGGAAGCCAAGCCAAATGGC
ATCTTTAAGGGCTCCACAGCTGAAAATGCAGAATACCTAAGGGTCGCGCCACAAAGCAGTGAATTTATTGGAGCA