

>EGFR H773_V774insH

ATGCGACCCTCCGGGACGGCCGGGCAGCGCTCCTGGCGCTGCTGGCTGCGCTCTGCCCGGCGAGTCGGGCTCT
GGAGGAAAAGAAAGTTTTGCCAAGGCACGAGTAACAAGCTCACGCAGTTGGGCACTTTTGAAGATCATTTTCTCAG
CCTCCAGAGGATGTTCAATAACTGTGAGGTGGTCTTGGGAATTTGGAAATTACCTATGTGCAGAGGAATTATGAT
CTTTCCTTCTTAAAGACCATCCAGGAGGTGGCTGGTTATGTCCTCATTGCCCTCAACACAGTGGAGCGAATTCCTT
GGAAAACCTGCAGATCATCAGAGGAAATATGTACTACGAAAATTCCTATGCCTTAGCAGTCTTATCTAACTATGAT
GCAAATAAAACCGGACTGAAGGAGCTGCCCATGAGAAATTTACAGGAAATCCTGCATGGCGCCGTGCGGTTCCAGC
AACAACCTGCCCTGTGCAACGTGGAGAGCATCCAGTGGCGGGACATAGTCAGCAGTGACTTTCTCAGCAACATG
TCGATGGACTTCCAGAACCACCTGGGCAGCTGCCAAAAGTGTGATCCAAGCTGTCCAATGGGAGCTGCTGGGGT
GCAGGAGAGGAGAACTGCCAGAACTGACCAAATCATCTGTGCCAGCAGTGTCCGGGCGCTGCCGTGGCAA
GTCCCCAGTGACTGCTGCCACAACCAGTGTGCTGCAGGCTGCACAGGCCCGGGAGAGCGACTGCCTGGTCTG
CCGCAAATTCGAGACGAAGCCACGTGCAAGGACACCTGCCCCCACTCATGCTCTACAACCCACCACGTACCAG
ATGGATGTGAACCCCGAGGGCAAATACAGCTTTGGTGCCACCTGCGTGAAGAAGTGTCCCCGTAATTATGTGGTG
ACAGATCACGGCTCGTGCGTCCGAGCCTGTGGGGCCGACAGCTATGAGATGGAGGAAGACGGCGTCCGCAAGTG
TAAGAAGTGCGAAGGGCCTTGCCGAAAAGTGTGTAACGGAATAGGTATTGGTGAATTTAAAGACTCACTCTCCAT
AAATGCTACGAATATTAACACTTCAAAAACCTGCACCTCCATCAGTGGCGATCTCCACATCCTGCCGGTGGCATTTA
GGGGTGACTCCTTCACACATACTCCTCCTCTGGATCCACAGGAACTGGATATTCTGAAAACCGTAAAGGAAATCAC
AGGGTTTTTGTGATTGAGGCTTGGCCTGAAAACAGGACGGACCTCCATGCCTTTGAGAACCTAGAAATCATACGC
GGCAGGACCAAGCAACATGGTCAGTTTTCTTGCAGTCGTCAGCCTGAACATAACATCCTTGGGATTACGCTCCC
TCAAGGAGATAAGTGATGGAGATGTGATAATTTAGGAAACAAAATTTGTGCTATGCAAATACAATAAACTGGA
AAAAACTGTTTGGGACCTCCGGTCAGAAAACCAAATATAAGCAACAGAGGTGAAAACAGCTGCAAGGCCACA
GGCCAGGTCTGCCATGCCTTGTGCTCCCCGAGGGCTGCTGGGGCCCGGAGCCCAGGGACTGCGTCTCTTGGCCG
AATGTCAGCCGAGGCAGGGAATGCGTGGACAAGTGAACCTTCTGGAGGGTGAGCCAAGGGAGTTTGTGGAGA
ACTCTGAGTGATACAGTGCCACCCAGAGTGCCTGCCTCAGGCCATGAACATCACCTGCACAGGACGGGGACCAG
ACAACGTATCCAGTGTGCCCACTACATTGACGGCCCCCACTGCGTCAAGACCTGCCCGGACAGGAGTATGGGAG
AAAACAACACCCTGGTCTGGAAGTACGCAGACGCCGCCATGTGTGCCACCTGTGCCATCCAACTGCACCTACG
GATGCACTGGGCCAGGTCTTGAAGGCTGTCCAACGAATGGGCCTAAGATCCCGTCCATCGCCACTGGGATGGTGG
GGGCCCTCCTCTTGTGCTGGTGGTGGCCCTGGGGATCGGCCTCTTATGCGAAGGCGCCACATCGTTCGGAAGC
GCACGCTGCGGAGGCTGCTGCAGGAGAGGGAGCTTGTGGAGCCTTTACACCCAGTGGAGAAGCTCCCAACCAA
GCTCTCTTGAAGGATCTTGAAGGAACTGAATCAAAAAGATCAAAGTGTGGGCTCCGGTGCCTTGGGCACGGTG
TATAAGGGACTCTGGATCCAGAAGGTGAGAAAAGTTAAATTCCTGTCGCTATCAAGGAATTAAGAGAAGCAACA
TCTCCGAAAGCCAACAAGGAAATCCTCGATGAAGCCTACGTGATGGCCAGCGTGGACAACCCCCACACCGTGTGC
CGCCTGCTGGGCATCTGCCTCACCTCCACCGTGAACCTCATCACGCAGCTCATGCCCTTCGGCTGCCTCCTGGACTA
TGTCCGGGAACACAAAGACAATATTGGCTCCAGTACCTGCTCAACTGGTGTGTGCAGATCGCAAAGGGCATGAA
CTACTTGGAGGACCGTGCCTTGGTGCACCCGACCTGGCAGCCAGGAACGTAAGTGGTAAAACACCCGACGATGT
CAAGATCACAGATTTTGGGCTGGCCAACTGCTGGGTGCGGAAGAGAAAGAATACCATGCAGAAGGAGGCAAAG
TGCCTATCAAGTGGATGGCATTGGAATCAATTTTACACAGAATCTATACCCACCAGAGTGTGCTGGAGTACGG
GGTGACTGTTTGGGAGTTGATGACCTTTGGATCCAAGCCATATGACGGAATCCCTGCCAGCGAGATCTCCTCCATC
CTGGAGAAAGGAGAACGCCTCCCTCAGCCACCCATATGTACCATCGATGTCTACATGATCATGGTCAAGTGTGGA
TGATAGACGCAGATAGTCGCCAAAGTTCCGTGAGTTGATCATCGAATTTCTCAAATGGCCCCGAGACCCCAAGC
GCTACCTTGTCAATCAGGGGGATGAAAGAATGCATTTGCCAAGTCTACAGACTCCAACCTTCTACCGTGCCTGAT
GGATGAAGAAGACATGGACGACGTGGTGGATGCCGACGAGTACCTCATCCACAGCAGGGCTTCTTACGACGCC
CCTCCACGTACGGACTCCCCTCCTGAGCTCTCTGAGTGAACCCAGCAACAATCCACCGTGGCTTGCATTGATAGA
AATGGGCTGCAAAGCTGTCCATCAAGGAAGACAGCTTCTTGCAGCGATACAGCTCAGACCCACAGGGCGCCTTG

ACTGAGGACAGCATAGACGACACCTTCCTCCCAGTGCCTGAATACATAAACCAGTCCGTTCCCAAAGGCCCGCTG
GCTCTGTGCAGAATCCTGTCTATCACAATCAGCCTCTGAACCCCGCACCCAGCAGAGACCCACACTACCAGGACCC
CCACAGCACTGCAGTGGGCAACCCCGAGTATCTCAACACTGTCCAGCCCACCTGTGTCAACAGCACATTCGACAGC
CCTGCCCACTGGGCCCAGAAAGGCAGCCACCAAATTAGCCTGGACAACCCTGACTACCAGCAGGACTTCTTTCCCA
AGGAAGCCAAGCAAATGGCATCTTTAAGGGCTCCACAGCTGAAAATGCAGAATACCTAAGGGTCGCGCCACAAA
GCAGTGAATTTATTGGAGCA