

>EGFR L858R sequence

ATGCGACCCTCCGGGACGGCCGGGCAGCGCTCCTGGCGCTGCTGGCTGCGCTCTGCCCGCGAGTCGGGCTCT
GGAGGAAAAGAAAGTTTGC AAGGCACGAGTAACAAGCTCACGCAGTTGGGCACTTTTGAAGATCATTTTCTCAG
CCTCCAGAGGATGTTCAATAACTGTGAGGTGGTCTTGGGAATTTGGAAATTACCTATGTGCAGAGGAATTATGAT
CTTTCCTTCTTAAAGACCATCCAGGAGGTGGCTGGTTATGTCCTCATTGCCCTCAACACAGTGGAGCGAATTCCTT
GGAAAACCTGCAGATCATCAGAGGAAATATGTACTACGAAAATTCCTATGCCTTAGCAGTCTTATCTAACTATGAT
GCAAATAAAAACCGGACTGAAGGAGCTGCCATGAGAAATTTACAGGAAATCCTGCATGGCGCCGTGCGGTTCCAGC
AACAACCTGCCCTGTGCAACGTGGAGAGCATCCAGTGGCGGGACATAGTCAGCAGTGACTTTCTCAGCAACATG
TCGATGGACTTCCAGAACCACCTGGGCAGCTGCCAAAAGTGTGATCCAAGCTGTCCAATGGGAGCTGCTGGGGT
GCAGGAGAGGAGAACTGCCAGAACTGACCAAATCATCTGTGCCAGCAGTGTCCGGGCGCTGCCGTGGCAA
GTCCCCAGTGACTGCTGCCACAACCAGTGTGCTGCAGGCTGCACAGGCCCGGGAGAGCGACTGCCTGGTCTG
CCGCAAATTCGAGACGAAGCCACGTGCAAGGACACCTGCCCCCACTCATGCTCTACAACCCACCACGTACCAG
ATGGATGTGAACCCCGAGGGCAAATACAGCTTTGGTGCCACCTGCGTGAAGAAGTGTCCCCGTAATTATGTGGTG
ACAGATCACGGCTCGTGCGTCCGAGCCTGTGGGGCCGACAGCTATGAGATGGAGGAAGACGGCGTCCGCAAGTG
TAAGAAGTGCGAAGGGCCTTGCCGAAAAGTGTGTAACGGAATAGGTATTGGTGAATTTAAAGACTCACTCTCCAT
AAATGCTACGAATATTAACACTTCAAAAACCTGCACCTCCATCAGTGGCGATCTCCACATCCTGCCGGTGGCATTTA
GGGGTGACTCCTTCACACATACTCCTCCTCTGGATCCACAGGAACTGGATATTCTGAAAACCGTAAAGGAAATCAC
AGGGTTTTTGTGATTGAGGCTTGGCCTGAAAACAGGACGGACCTCCATGCCTTTGAGAACCTAGAAATCATACGC
GGCAGGACCAAGCAACATGGTCAGTTTTCTTGCAGTCGTCAGCCTGAACATAACATCCTTGGGATTACGCTCCC
TCAAGGAGATAAGTGATGGAGATGTGATAATTTAGGAAACAAAATTTGTGCTATGCAAATACAATAAACTGGA
AAAAACTGTTTGGGACCTCCGGTCAGAAAACCAAATATAAGCAACAGAGGTGAAAACAGCTGCAAGGCCACA
GGCCAGGTCTGCCATGCCTTGTGCTCCCCGAGGGCTGCTGGGGCCCGAGGCCAGGGACTGCGTCTCTTGGCCG
AATGTCAGCCGAGGCAGGGAATGCGTGGACAAGTGAACCTTCTGGAGGGTGAAGCAAGGGAGTTTGTGGAGA
ACTCTGAGTGATACAGTGCACCCAGAGTGCCTGCCTCAGGCCATGAACATCACCTGCACAGGACGGGGACCAG
ACA ACTGTATCCAGTGTGCCACTACATTGACGGCCCCACTGCGTCAAGACCTGCCCGCAGGAGTATGGGAG
AAAACAACACCCTGGTCTGGAAGTACGCAGACGCCGCCATGTGTGCCACCTGTGCCATCCAACTGCACCTACG
GATGCACTGGGCCAGGTCTTGAAGGCTGTCCAACGAATGGGCCTAAGATCCCGTCCATCGCCACTGGGATGGTG
GGGCCCTCCTCTTGTGCTGGTGGTGGCCCTGGGGATCGGCCTCTTATGCGAAGGCGCCACATCGTTCGGAAGC
GCACGCTGCGGAGGCTGCTGCAGGAGAGGGAGCTTGTGGAGCCTTTACACCCAGTGGAGAAGCTCCCAACCAA
GCTCTCTTGAAGGATCTTGAAGGAACTGAATCAAAAAGATCAAAGTGTGGGCTCCGGTGCCTTCGGCACGGTG
TATAAGGGACTCTGGATCCCAGAAGGTGAGAAAAGTTAAATTCCTGTCGCTATCAAGGAATTAAGAGAAGCAACA
TCTCCGAAAGCCAACAAGGAAATCCTCGATGAAGCCTACGTGATGGCCAGCGTGGACAACCCCCACGTGTGCCGC
CTGCTGGGCATCTGCCTCACCTCCACCGTGCAGCTCATCACGCAGCTCATGCCCTTCGGTGCCTCCTGGACTATGT
CCGGGAACACAAAGACAATATTGGCTCCAGTACCTGCTCAACTGGTGTGTGCAGATCGCAAAGGGCATGAACTA
CTTGGAGGACCGTGCCTTGGTGCACCGCGACCTGGCAGCCAGGAACGTACTGGTAAAACACCGCAGCATGTCAA
GATCACAGATTTTGGGCGGCCAACTGCTGGGTGCGGAAGAGAAAAGAATACCATGCAGAAGGAGGCAAAGTG
CCTATCAAGTGGATGGCATTGGAATCAATTTTACACAGAATCTATACCCACCAGAGTGATGTCTGGAGCTACGGGG
TGACTGTTTGGGAGTTGATGACCTTTGGATCCAAGCCATATGACGGAATCCCTGCCAGCGAGATCTCCTCCATCCT
GGAGAAAGGAGAACGCCTCCCTCAGCCACCCATATGTACCATCGATGTCTACATGATCATGGTCAAGTGTGGAT
GATAGACGCAGATAGTCGCCAAAGTTCCTGAGTTGATCATCGAATTTCCAAAATGGCCGAGACCCCCAGCG
CTACCTTGTATTAGGGGGATGAAAGAATGCATTTGCCAAGTCTACAGACTCCAATTCTACCGTGCCTGATG
GATGAAGAAGACATGGACGACGTGGTGGATGCCGACGAGTACCTCATCCACAGCAGGGCTTCTCAGCAGCCCC
TCCACGTACGGACTCCCCTCCTGAGCTCTCTGAGTGCAACCAGCAACAATCCACCGTGGCTTGCATTGATAGAA
ATGGGCTGCAAAGCTGTCCCATCAAGGAAGACAGTCTTTCGAGCGATACAGCTCAGACCCACAGGCGCCTTGA

CTGAGGACAGCATAGACGACACCTTCCTCCCAGTGCCTGAATACATAAACCAGTCCGTTCCAAAAGGCCCGCTGG
CTCTGTGCAGAATCCTGTCTATCACAAATCAGCCTCTGAACCCCGCACCCAGCAGAGACCCACACTACCAGGACCCC
CACAGCACTGCAGTGGGCAACCCCGAGTATCTCAACACTGTCCAGCCCACCTGTGTCAACAGCACATTGACAGCC
CTGCCCACTGGGCCCAGAAAGGCAGCCACCAAATTAGCCTGGACAACCCTGACTACCAGCAGGACTTCTTTCCCAA
GGAAGCCAAGCCAAATGGCATCTTTAAGGGCTCCACAGCTGAAAATGCAGAATACCTAAGGGTCGCGCCACAAAG
CAGTGAATTTATTGGAGCA