

# Active Kinase Mutants



## Technical Bulletin

### Essential Tools for Developing Targeted Drugs

Mutations in protein kinases are often implicated in the primary causes of various diseases. These include autoimmune and metabolic disorders, as well as cancer. Mutations in kinases can also confer drug resistance, which poses an increasing challenge to development of effective therapeutics.

SignalChem offers the most extensive range of medically relevant human recombinant Active Kinase Mutants, specifically to support research and development for the next generation of kinase drugs. Our research and development team focus on expanding this particular line of products, as to provide the best tools for understanding and curing human diseases.

#### The Largest Number of Mutant Kinases

- Disease Relevant Mutants
- 200 Active Kinase Mutants and Growing
- Representing Over 50 Different Proteins

#### Dependable Performance

- >80% Average Purity
- Ready for Kinase Activity Assays
- Lot-to-lot Consistency

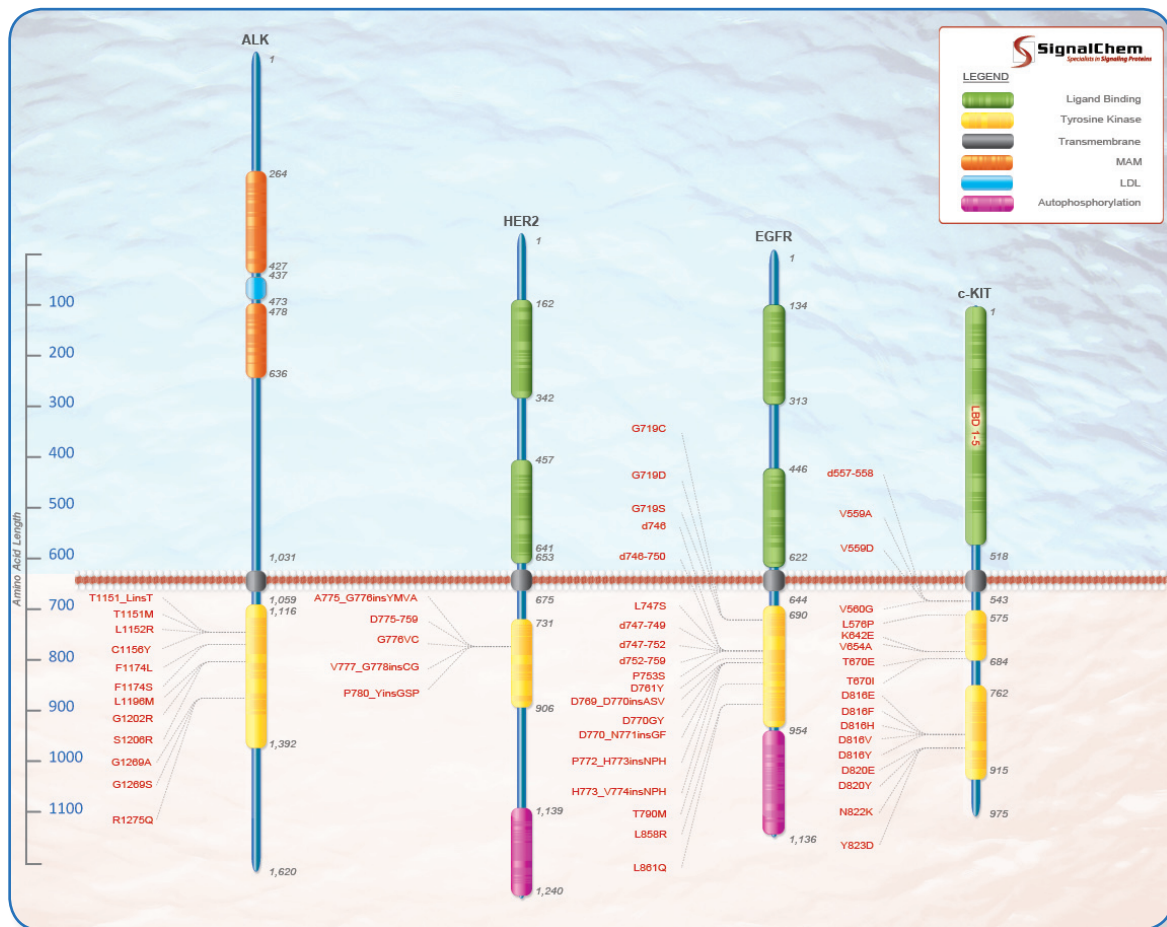


Figure: Sample of available Active Kinase Mutants from SignalChem (Specific Mutations in Red)