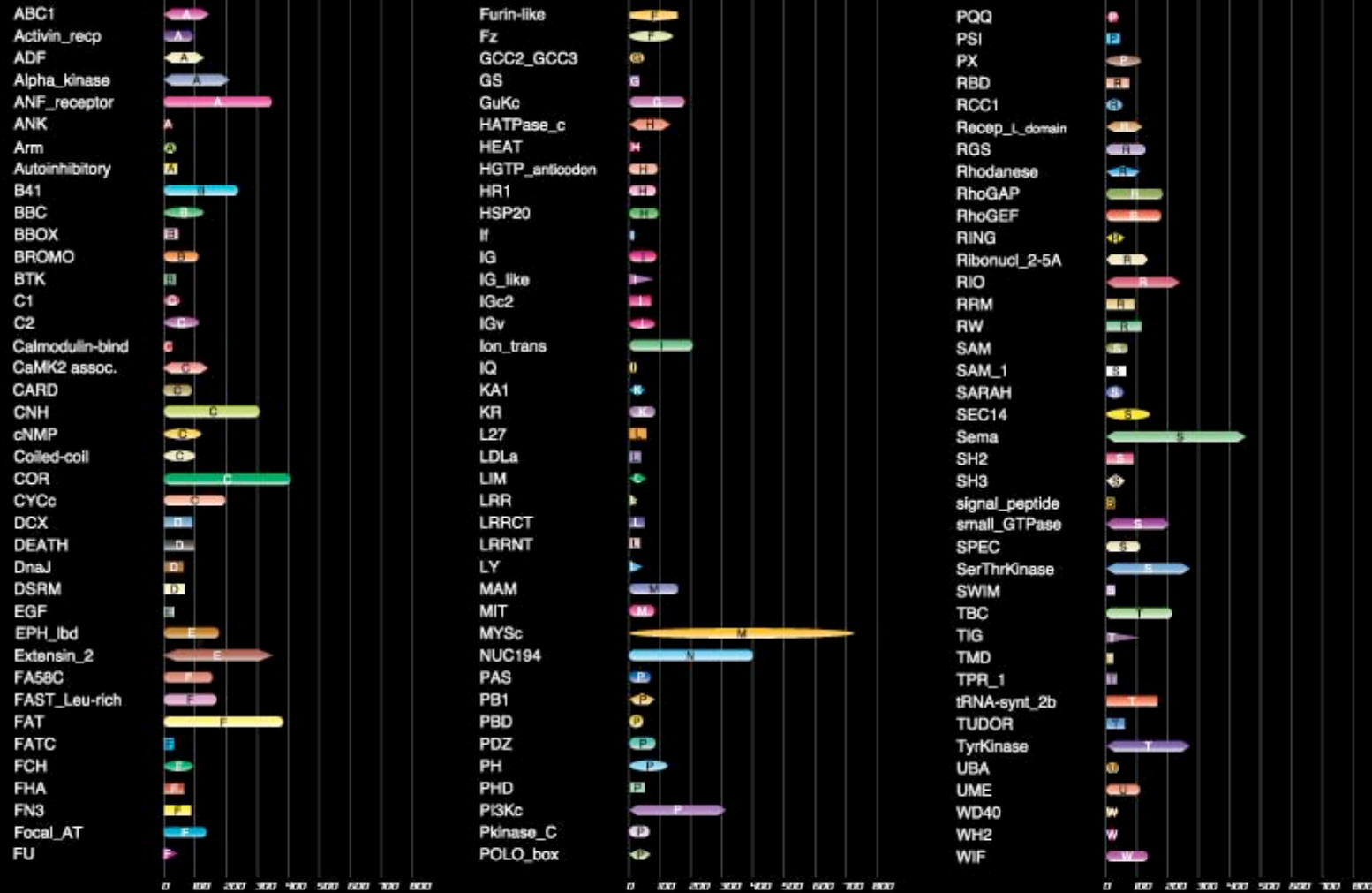


Legend for Domains in Human Protein Kinases



Phosphorylation site type: ● Phospho-Serine ■ Phospho-Threonine ▼ Phospho-Tyrosine

Protein kinases are identified with their Kinbase_Name (www.kinase.com) and HGNC_Name (www.gene.ucl.ac.uk/nomenclature/index.html).
 Further domain information can be obtained from <http://pfam.wustl.edu/browse.shtml> and <http://smart.embl-heidelberg.de/browse.shtml>.
 Further phosphorylation site information can be retrieved from <http://www.phosphosite.org>.

