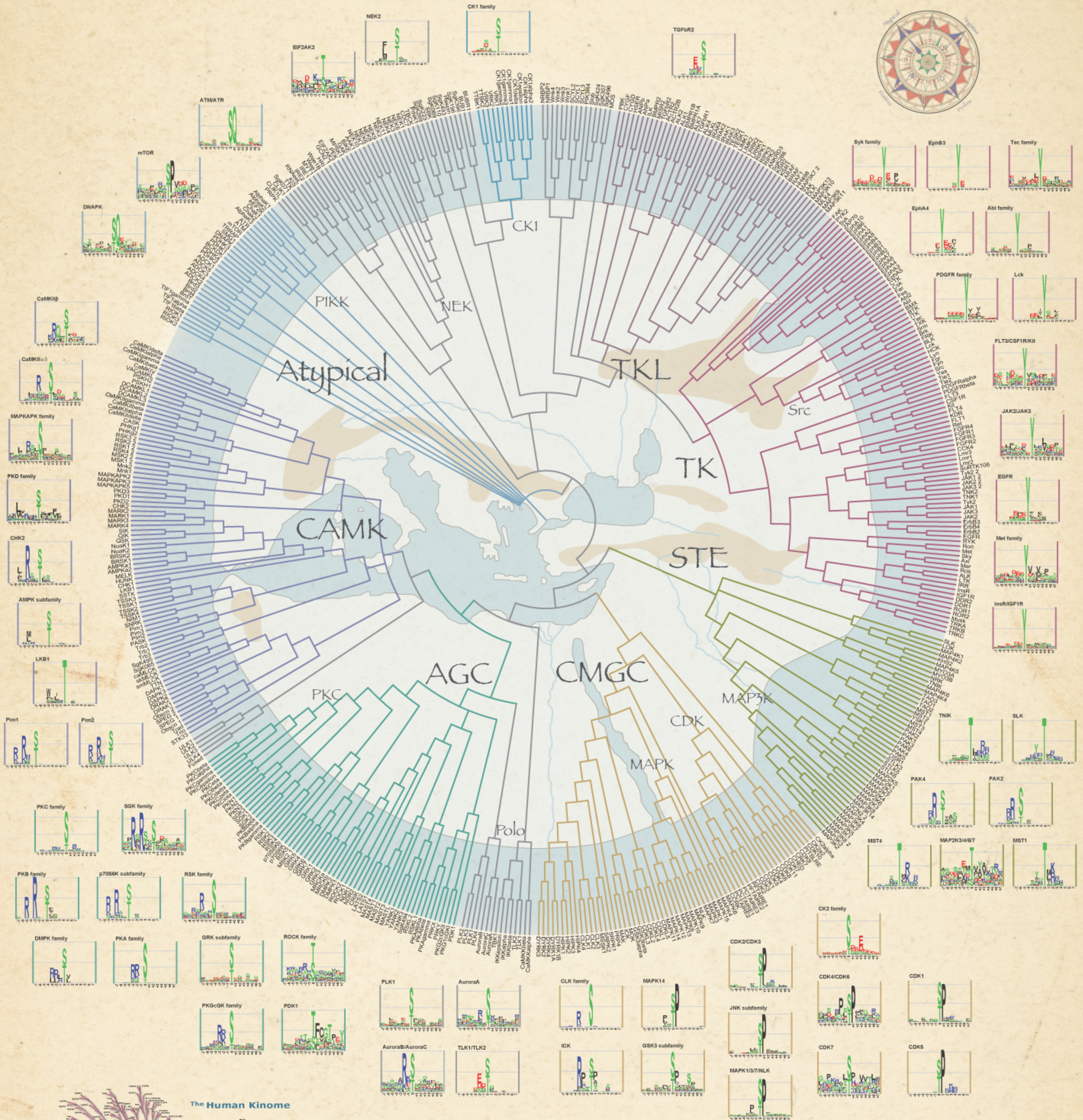


A Sequence-Specificity Atlas of the Kinase World



Finding the Targets of the Human Kinome

Protein kinases are enzymes that mediate cellular decision processes by catalyzing the addition of a negatively charged phosphate group to protein substrates, which can subsequently be recognized by phosphorylation-dependent binding domains in other proteins. The human genome encodes 518 protein kinases, which have been organized into a tree that represents evolutionary relationships. The substrate specificities of kinases are in part determined by the amino acid sequence of the phosphorylation sites, and comprehensively mapping the consensus sequence motif recognized by each kinase catalytic domain is thus crucial for understanding phosphorylation-mediated signaling networks.

NetPhorest (<http://netphorest.info>) is a community resource that uses phylogenetic trees to organize data from both in vivo and in vitro experiments to derive sequence specificities for individual kinase or phosphorylation-

dependent binding domains or families of closely related domains. This poster presents the specificities of kinases as sequence logo plots around a phylogenetic map of the human protein kinases. The height of a stack of letters represents the information content of a particular position relative to a phosphorylation site, and the height of the individual letters is proportional to the frequency of the amino acid residue in question. The serine, threonine, or tyrosine phosphoacceptor residues are placed at the center of each logo plot. Closely related kinases typically have similar substrate specificities; however, distantly related kinases may not have divergent substrate specificities. The compass summarizes the major trends of the specificity atlas in the center; for example, most kinases that prefer positively charged residues are located in the lower left-hand part of the phylogenetic map.

Poster authors: Martin Lee Miller, Soren Brunak, Lars Juhl Jensen, Michael B. Yaffe, Rune Linding
Science Signaling coordinator: Nancy R. Gosugh
Design and production: Lars Juhl Jensen, Rune Linding, Dave Comb, Aileen Taylor, Marcus Spiegler
Copyeditor: Harry Jach
Sponsored by: Cell Signaling Technology, Inc.
 This poster accompanies the Science Signaling paper "Linear Motif Atlas for Phosphorylation-Dependent Signaling" by Martin Lee Miller, Lars Juhl Jensen, Francesca Diella, Claus Jorgensen, Michele Timi, Lei Li, Marilyn Huang, Sirkka A. Parker, Jennifer Bordeaux, Thomas Schaeuble, Martina Orlowski, Adam Pasculic, Jes Alexander, Stefan Knapp, Nikolaj Blom, Peer Bork, Sharon Li, Gianni Cesareni, Tony Pawson, Benjamin E. Turk, Michael B. Yaffe, Soren Brunak, Rune Linding.

1 Center for Biological Sequence Analysis, Technical University of Denmark, 2800 Lyngby, Denmark.
 2 The Novo Nordisk Foundation Center for Protein Research, University of Copenhagen, 2200 Copenhagen, Denmark.
 3 European Molecular Biology Laboratory, 65117 Heidelberg, Germany.
 4 Samuel Lunenfeld Research Institute, Mount Sinai Hospital, Toronto, Ontario, Canada M5G 1X5.
 5 University of Rome, Tor Vergata, 00133 Rome, Italy.
 6 University of Western Ontario, London, Ontario, Canada N6A 5S1.
 7 Department of Pharmacology, Yale University School of Medicine, New Haven, CT 06520 USA.
 8 Center for Cancer Research, Massachusetts Institute of Technology, Cambridge, MA 02139, USA.
 9 Structural Genomics Consortium, University of Oxford, Oxford OX3 7DG, UK.
 10 Max-Delbrück Center for Molecular Medicine, 13082 Berlin, Germany.
 11 Cellular & Molecular Logic Team, The Institute of Cancer Research, London SW3 6JB, UK.
 *These authors contributed equally to this work.
 †To whom correspondence should be addressed. E-mail: brunak@dbu.dk (S.B.), myaffe@mit.edu (M.B.Y.), and rune.linding@gmail.com (R.L.)

