

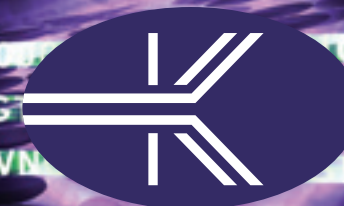
Kinexus is revolutionizing the discovery-based research paradigm so all researchers can access + benefit from systems biology approaches to investigate their experimental model systems.

We have created a powerful discovery engine that is unparalleled for the generation + analysis of quantitative proteomics data.

Kinexus Bioinformatics Corporation
Suite 402, 6190 Agronomy Road
Vancouver, British Columbia, Canada V6T 1Z3

Mapping cell signalling systems + identifying their linkages to normal + disease-related processes is the most challenging + promising biomedical research endeavor of this decade.

Kinexus supports this critical research effort through our unrivalled suite of antibody-base proteomics + bioinformatics services.



KINEXUS™
LISTENING TO CELLS TO SILENCE DISEASE

website	toll free phone	facsimile	e-mail
www.kinexus.ca	866-KINEXUS	604-822-9964	info@kinexus.ca

Researchers usually guess which signalling proteins may be relevant in their work, gamble on which antibodies to use + trust that their experiments are performed without other technical issues.

Kinexus minimizes the risks + maximizes the prospects for research success by providing services that are unmatched for breadth, reliability, speed, convenience + affordability.

KiNET™

The KiNET™ bioinformatics services provide website access to our unique functional proteomics database. For a low subscription fee, clients have unlimited Internet use of KiNET™ to query about the expression levels and phosphorylation states of hundreds of target proteins in hundreds of well defined experimental model systems. Using KiNET™, clients can compare their own Kinetworks™ results with the data from thousands of other experiments analyzed with the very same reagents + methodology by our scientists and technicians.

Our standard Kinetworks™ multi-immunoblotting services track panels of 25 to 78 thematically-related signal transduction protein targets for their expression + phosphorylation with high accuracy. With our custom Kinetworks™ services, clients can choose any 18 from over 500 different pan- + phospho-site-specific antibodies for analysis of 1 tissue/cell sample or up to 3 targets of distinct molecular mass in 8 different samples. The custom Kinetworks™ services are ideal for validation of results obtained from our standard Kinetworks™ and Kinex™ services.

KiNEX™

Our Kinex™ antibody microarray services monitor the binding of over 500 target proteins to the same antibodies used in our Kinetworks™ services. There are many reasons why antibody microarrays can generate false positive and negative data, including cross-reactivity, protein complexes and masked epitopes. However, the cost, speed + low specimen material requirement for these services makes this a high value proposition for our clients. The Kinex™ approach is an ideal first step for novel biomarker discovery.

KiNETXWORKS™

KINETTM

Text about the service

KINEXTM

Text about the service

KINET
KIRKSTM

Text about the service



KINEXUSTM
LISTENING TO CELLS TO SILENCE DISEASE