



**KINEXUS**

## **KINEXUS BIOINFORMATICS CORPORATION HONORED AS A RECIPIENT OF RED HERRING'S TOP 100 PRIVATE COMPANIES**

*Award recognizes 100 private companies in North America that drive the technology industry's future*

**PRESS RELEASE - 2005 MAY 13**

**FOR IMMEDIATE RELEASE MAY 13, 2005**

**VANCOUVER, British Columbia** – Kinexus Bioinformatics Corporation, a Canadian proteomics company, was recently honored by Red Herring as one of the Top 100 Private Companies in North America. The award winning Silicon Valley media company selected Kinexus from over 900 companies based on its demonstrated critical contributions in novel technologies and services and its innovative business strategy.

“Five years after the dot-com bubble burst, innovation and entrepreneurship are thriving. The companies on this list are proof,” said Red Herring editor-in-chief Joel Dreyfuss. Red Herring’s lists of private companies are an important part of the magazine’s tradition of identifying new and innovative technology companies and entrepreneurs. Companies like Google and eBay were spotted in their early days by Red Herring editors as those that would change the way we live and think. Since 1996, the *Red Herring 100* has served as the place to look for emerging leaders; young companies that retain the capacity to disrupt existing markets and create entirely new ones.

“We are thrilled to be honored by Red Herring” said Dr. Steven Pelech, President and Chief Scientific Officer for Kinexus, “especially being one of only two Canadian companies chosen. Kinexus has a unique business model and proprietary technologies that generate revenues and expedites the drug discovery process for academia and biopharmaceutical companies. The accumulation of 6 years of data generated from our proteomics services into a query able, Internet subscription database, called KiNET, is the realization of the Kinexus vision”.

Kinexus is poised to become a major proteomics and bioinformatics player with its focus on protein kinases and other cell signalling proteins for diagnostic and therapeutic applications. The world market for kinase profiling is estimated at around \$200 million and climbing. Kinexus currently offers 13 unique signal transduction protein profiling services for tracking more than 200 protein kinases, 200 known phosphorylation site targets of kinases and 110 other signalling proteins that regulate cell proliferation, stress and death. The application of these methods in characterizing cell signalling networks provides Kinexus and its clients with powerful information for disease diagnosis and drug development for personalized therapies to improve human health.

The winners of the Top 100 Private Companies in North America will be recognized at the Red Herring conference from May 17-19, 2005 in Monterey, California and will be published in the May 23, 2005 issue of Red Herring Magazine. A list of the winning companies can be viewed on the Red Herring website at: [www.redherring.com](http://www.redherring.com).

**For further information, please contact Kinexus Bioinformatics Corporation toll free at  
1-866-KINEXUS or visit our website at [www.kinexus.ca](http://www.kinexus.ca).**