



KINEXUS

RED HERRING SELECTS KINEXUS BIOINFORMATICS CORPORATION FOR PRESTIGIOUS RED HERRING TOP 100 INNOVATORS AWARD

The Award Recognizes 100 Companies That Embody Innovation

PRESS RELEASE - 2004 DECEMBER 7

FOR IMMEDIATE RELEASE DECEMBER 7, 2004

VANCOUVER, British Columbia – Kinexus Bioinformatics Corporation, a Canadian proteomics company is pleased to announce that it has recently been selected as a Top 100 Innovator by Red Herring, the award winning Silicon Valley media company that covers technology, innovation, entrepreneurship and business. The award honors 100 international public and private companies that have made their mark with critical contributions in technologies, products and services and have demonstrated creativity and innovation in marketing, sales, financing, human relations and business strategies.

The editors worked hard to identify the top companies who will benefit from the important technology trends in the coming year,” said Joel Dreyfuss, Editor-in-Chief of Red Herring. Red Herring’s editorial staff evaluated over 1200 submissions from 900 companies. The editors used a careful analysis of financial data as well as subjective criteria, including quality of management, execution of strategy, and dedication to research and development. Kinexus was 1 of just 12 life sciences companies chosen for this award, while the other 88 companies largely dominated the information technology, communications, and energy industries. The Red Herring Award places Kinexus among the top 100 most innovative technology companies in the world, which includes such technology giants as IBM, Sun Microsystems, Hewlett Packard, Amazon.Com and Google.

“We are honored to be chosen as 1 of only 3 Canadian companies by the editors of Red Herring magazine and recognized as an industry leader amongst a prominent and elite group of 100 international companies,” said Dr. Steven Pelech, President and Chief Scientific Officer of Kinexus and a Professor in the Department of Medicine at the University of British Columbia, Canada. “This distinction acknowledges our continuous innovation in the health care industry and our dedication to putting that innovation to work for our customers, our shareholders and ultimately for the benefit of the public. We look forward to expanding our strategic vision into 2005 and beyond.”

The Top 100 Innovators were announced during the Red Herring Fall 2004 conference, December 6-8, 2004 at the Monterey Plaza Hotel in Monterey, California and will be published in the December 13, 2004 issue of Red Herring Magazine. A list of the winning companies can be viewed on the Red Herring website at: 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.

About Kinexus

Kinexus Bioinformatics Corporation is a private, biopharmaceutical company engaged in the research and development of innovative proteomics methods to discover and exploit the relationships of signalling proteins in cellular communication networks. The application of this knowledge positions Kinexus and its clients in disease diagnosis, drug development and personalized therapies to improve human health. The Kinetworks™ screening services are a cost-effective solution for the identification of disease diagnostic markers and drug targets, delineation of mechanisms of drug action and toxicity, and validation of animal models for testing drug leads. Kinexus currently tracks over 180 protein kinases, 250 phospho-sites in kinase targets, and 120 other signalling proteins that regulate cell reproduction, stress responses and cell death. For further information, please contact Kinexus Bioinformatics Corporation toll free at 1-866-KINEXUS or visit our website at www.kinexus.ca.



About Red Herring

Red Herring covers technology innovation, venture financing, and the deals that make a difference. Its award-winning journalists go deeper, providing a comprehensive, critical analysis of what's new and why it matters. Red Herring covers the business trends with insight that counts, serving today's entrepreneurs and the technology of tomorrow. Visit www.redherring.com for penetrating articles, newsletters, and research reports.