



**KINEXUS**

## **Kinexus Bioinformatics Corporation is Pleased to Announce the Appointment of Ms. Catherine Sutter as Director of Human Resources**

*HR Director to Implement Employment Contracts and New Policies for Employees*

**FOR IMMEDIATE RELEASE SEPTEMBER 1, 1999**

**VANCOUVER, British Columbia** - Dr. Steven Pelech, President and C.E.O. of Kinexus Bioinformatics Corporation, is pleased to announce the appointment of Ms. Catherine Sutter as the Company's Director of Human Resources effective September 1, 1999. Her new position involves the development and implementation of policies and procedures affecting all employees including organization design and development, performance management, training and education, recruitment and retention, compensation, benefits and incentives, equity programs, and health and safety.

Ms. Sutter has over ten years experience in human resources and administration including three years as the Director of Human Resources for Kinetek Pharmaceuticals, Inc. and four years as the Director of Biotechnology for Kinetek Biotechnology Corporation. Prior to that, she worked as a Laboratory Manager at the University of British Columbia and Office Administrator for Gulf Shark Research in Bayou la Batre, Alabama. She is also a founding principal and life-time member of the Board of Directors of the Children's Research Foundation of British Columbia, a charitable Foundation dedicated to the advancement of research and knowledge of life-threatening and disabling diseases affecting children.

Kinexus Bioinformatics Corporation capitalizes on protein kinases to revolutionize medicine by improving the efficiency of drug discovery in support of personalized healthcare. Kinexus is harnessing the powerful synergies of genomics, proteomics, bioinformatics, and the Internet to advance drug discovery, disease diagnosis and global biomedical research. Protein kinases, key proteins responsible for communication and control inside cells, operate within complex networks of functionally interconnected signalling pathways. Their malfunction has been linked to hundred of diseases. They regulate all cellular processes by catalyzing the reversible phosphorylation of thousands of different target proteins. The importance of protein kinases as critical targets for the diagnosis and treatment of diseases is well recognized in the pharmaceutical industry.

**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).**

**PRESS RELEASE - 1999 SEPTEMBER 1**