



**KINEXUS**

## **Kinexus Bioinformatics Corporation Appoints Mr. Kevin McDuffie as Director of Corporate Development and Mr. Jimmy Wong as Controller**

*Two Key Players Recently Appointed to the Management Team of Kinexus*

**PRESS RELEASE - 2000 JULY 10**

### **FOR IMMEDIATE RELEASE JULY 10, 2000**

**VANCOUVER, British Columbia** - Dr. Steven Pelech, President and CEO of Kinexus Bioinformatics Corporation is pleased to announce the appointment of Kevin McDuffie to the position of Director of Corporate Development, and Jimmy Wong, as Controller.

Kevin McDuffie has an extensive background in licensing, contract negotiations and sales and marketing. His experience includes five years with University Technologies International Inc. as the Manager of Technology Transfer at the University of Calgary, and ten years with Kodak, Bio-Rad and Pharmacia. He completed his Bachelor of Science in Biochemistry (honors) at the University of Saskatchewan.

Jim Wong contributes fifteen years of international financial management expertise to Kinexus, most recently as President and Controller of Leader Direct Response Services Ltd. Prior to that, he held positions as Controller and Financial Accountant with several U.S. and Canadian firms. He obtained his Bachelor of Science from the University of British Columbia, a Diploma of Technology (honors) in Financial Management and is currently completing his MBA.

Kinexus is a private proteomics/bioinformatics company specializing in protein kinases for the molecular diagnosis of disease and drug discovery. Kinexus has developed unique and robust screening methods of high stringency to facilitate the validation of drug targets and the rapid exclusion of inappropriate drug leads in the early stages of the drug discovery process. The application of these methods in characterizing cell signalling networks provides Kinexus, its clients and associates with powerful knowledge for drug development, rational drug design, disease diagnosis, and personalized therapies.

**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).**