



KINEXUS

## **Kinexus Bioinformatics Corporation Announces the Launch of its Fourth Proteomics Analysis Service**

*Stress and Heat Shock Protein Screen Available to the Public*

**FOR IMMEDIATE RELEASE JANUARY 30, 2002**

**VANCOUVER, British Columbia** - Kinexus Bioinformatics Corporation is pleased to announce the release of a new proteomics service with the Kinetworks™ Stress and Heat Shock Protein Screen, or KHSP-1.0. The KHSP screen is designed to track up to 25 stress-regulated and heat shock proteins on a single gel and is the fourth signal transduction protein screen offered by Kinexus. Heat shock proteins (HSP) are currently the focus of intense work by scientists worldwide as markers of stress and therapeutic adjuvants to treat cancer and other diseases with immune system activation. The KHSP Screen is complementary to the three existing services available from Kinexus including the Kinetworks™ Protein Phosphatase Screen (KPPS), Phospho-Site Screen (KPSS) and Protein Kinase Screen (KPKS). Sample preparation instructions, service agreements and order forms are available online at [www.kinexus.ca](http://www.kinexus.ca).

“These screens offer biomedical researchers in academia and the biopharmaceutical industry the ability to simultaneously track the presence and activation states of hundreds of different regulatory and cell signalling proteins to assist in the discovery of new drug targets, new drugs, and the molecular diagnosis of disease” said Dr. Steven Pelech, President and Chief Executive Officer of Kinexus, and a Professor of Medicine at the University of British Columbia. “Kinexus continues to exceed its milestones in product development, with apoptosis and cell cycle protein screens expected to be released in the second quarter of 2002”.

Kinexus Bioinformatics Corporation is a private, biopharmaceutical company engaged in the research and development of innovative methods to understand the relationships of signalling proteins in cellular communication networks. The application of this knowledge positions Kinexus and its clients in drug development, rational drug design, disease diagnosis and personalized therapies to improve human health.

**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 - 2002 JANUARY 30**