



KINEXUS

New Kinexus Website with Technical & Educational Resources for Cell Signalling Research

New Knowledge Based Website for Biomedical Research

FOR IMMEDIATE RELEASE OCTOBER 26, 2007

VANCOUVER, British Columbia – Kinexus Bioinformatics Corporation, a world leader in functional proteomics research, is pleased to announce the launch of its newly redesigned corporate website located at www.kinexus.ca. In keeping with Kinexus' commitment to customer service excellence and technological innovation, the new website offers scientists and other interested individuals additional ease and flexibility in accessing information, new features and greatly expanded content. The navigation on the site has also been enhanced and the entire site was built using internal resources, tools, and expertise.

“Kinexus is always looking for ways to improve and expand the range of services we provide to our clients. As we revamped the website, we took every opportunity to add features that would make it easier to use and to incorporate much more detailed information for the expert and the novice on cell signalling systems” said Dr. Steven Pelech, President and Chief Scientific Officer of Kinexus. “We hope to make our website one of the world’s leading cell signalling resources for information and education.”

In addition to its new look and feel, the newly redesigned website contains many valuable scientific resources and free educational downloads to facilitate signal transduction research and teaching. Key features of the new site include:

- Detailed information about the suite of unique proteomics and bioinformatics services offered by Kinexus and links to nearly 150 scientific publications using these services;
- Over 100 charts and figures related to cell signalling proteins;
- Nearly 100 movies and animations, including animations of 3D crystallographic structures of signal transduction proteins in complexes;
- Links to about 200 other useful genomics, proteomics and bioinformatics websites as well as other fascinating and entertaining websites;
- Links to over 200 published scientific reviews about cell signalling;
- Galleries with science-related movies, animations, still art, and photographs as well as links to other website galleries that feature compelling imagery;
- Glossary of scientific terms; and
- Visual tours of the Kinexus facilities and photographs of our staff.

Kinexus currently has service agreements with over 1000 laboratories in life science companies, universities, government institutions, and hospitals in over 33 different countries. Kinexus is a private, biotechnology company engaged in the research and development of innovative methods to track hundreds of signalling proteins in cellular communication networks. The application of this knowledge positions Kinexus and its clients in disease biomarker discovery, rational drug target identification, and drug development, to permit delivery of personalized therapies to improve human welfare.

**For further information, please contact Kinexus Bioinformatics Corporation
toll free at 1-866-KINEXUS or visit our website at www.kinexus.ca**

PRESS RELEASE - 2007 OCTOBER 26