



KINEXUS

Kinexus Announces the Launch of New Peptide Production Services

New Peptide Array Facility Designed to Expedite the Drug Discovery Process

PRESS RELEASE - 2010 MARCH 16

FOR IMMEDIATE RELEASE MARCH 16, 2010

VANCOUVER, British Columbia – Kinexus Bioinformatics Corporation, a world leader in functional proteomics research, announced the launch of a new peptide synthesis facility to provide researchers with the latest in peptide array technologies. These new services will facilitate and expedite access to peptide platforms for innovative approaches to proteomic research and discovery. A variety of peptide services are being made available by Kinexus to the research community including:

- Protein labeling,
- Screening of peptide arrays like Epitope Mapping, Substitution Analysis, and
- Peptide Library Screening,
- Custom Peptide Synthesis, including modified peptides like phosphopeptides, histones, cyclized and labeled peptides
- CelluSpots™ Custom Peptide Arrays on Glass Slides, and
- CelluSpots™ Kinase Substrate Peptide Arrays on Glass Slides.

“Kinexus’ peptide-based products strongly complement our existing portfolio of cell signaling, proteomics and bioinformatics services,” stated Dr. Steven Pelech, president and Chief Scientific Officer of Kinexus Bioinformatics Corporation. “These new products will give Kinexus customers access to high-throughput peptide synthesis and screening platforms, which will accelerate research in the field of signal transduction. The Peptide Array Facility is the latest addition to a unique integrated suite of proteomics services offered by Kinexus which delivers solutions and services for pharmaceutical, biotechnology, academic, life science, clinical and government research laboratories from around the world.

To learn more about the new Peptide Production Services or any of the other proteomics services available, please visit www.kinexus.ca. Kinexus Bioinformatics Corporation is a privately held, biotechnology company engaged in the research and development of innovative methods to understand the relationships of signaling proteins in molecular 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