



KINEXUS

Kinexus Bioinformatics Corporation Appoints Mr. Lyle Arnason as Director of Sales & Marketing

Former Bayer, Inc. Executive to Oversee Kinetworks™ Screening Services

PRESS RELEASE - 2001 JUNE 1

FOR IMMEDIATE RELEASE JUNE 1, 2001

VANCOUVER, British Columbia - Dr. Steven Pelech, President and C.E.O. of Kinexus Bioinformatics Corporation, is pleased to announce the appointment of Mr. Lyle Arnason as the Company's new Director of Sales and Marketing, effective June 1, 2001.

Mr. Arnason has over ten years of experience in the life science field most recently with Bayer Inc. as the Western Regional Business Manager. He has extensive experience in international sales, marketing and contract negotiations. Prior to Bayer Inc., Mr. Arnason held various positions with Becton Dickinson, Roche (formerly Boehringer Mannheim) and Fisher Scientific. He completed his B.Sc. at the University of Alberta and has an M.B.A degree with a Marketing specialization.

Kinexus Bioinformatics Corporation is a Vancouver-based biotechnology/information technology company striving to emerge as a world-leader in providing knowledge about intracellular communication systems that involve protein kinases. Defective cell signalling through these enzymes has been linked to hundreds of diseases in humans. Kinexus plans to deduce the composition and architecture of kinase networks and apply this knowledge for human health and welfare with diversified partners and clients by:

- providing cost effective screening services to track the presence and regulation of hundreds of different kinases and their targets;
- developing and marketing proprietary kinase-focused databases;
- modeling disease states in signal transduction pathways; and
- identifying drug-pathway interactions by 3D molecular modeling of kinases for rational design of drug candidates.

Kinexus is strategically positioned to take full advantage of the success of the human genome project 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.