



**KINEXUS**

## **Kinexus Announced as Third Winner of the ADP Small Business Grant Contest**

*Win will help Vancouver-based bioinformatics firm hire university students*

**Toronto, March 11, 2013** – ADP Canada announced today that Vancouver-based Kinexus, a biotechnology company dedicated to improving diagnostics and therapeutics for treating diseases including cancer and diabetes among others, is the third monthly winner of ADP Canada’s Small Business Grant Contest. As part of the contest, ADP Canada is giving six entrepreneurs \$10,000 cash plus free payroll services for a year. One winner is chosen and announced each month during the contest period.

In its winning entry, Kinexus described how the company’s work focuses on understanding a group of important enzymes that regulate all cellular processes. The malfunction of these proteins has been linked to more than 400 human diseases. Having a better understanding of how these enzymes work is key to finding the right diagnostics and medications to recognize and treat disease.

“One of the biggest challenges facing Kinexus is research funding,” said Catherine Sutter, Director of Sales and Marketing for Kinexus. “There have been a lot of cuts to government research grants to our clients and it is increasingly difficult to get private investors to invest in biotech companies.”

As a consistent source of funding, Kinexus has developed a business model that enables it to provide service contracts and analyses to more than 750 academic labs and 200 industrial labs for more than 100 companies around the world, including most leading pharmaceutical and biotech companies. With the \$10,000, Kinexus plans to hire local university students to assist with the design, development and programming of a database that will play an important part in mapping all of the 22,000 proteins encoded by the human genome and their interconnections with each other. Upon completion the database would be made available to the scientific community worldwide with open-access to accelerate the development of better medications to treat disease and improve patient care.

Through its partnerships with more than 400,000 small businesses across North America, ADP has unique insights into the challenges that confront owners every day. ADP’s innovative small business solutions help reduce risks associated with running an establishment—from human resources to cash flow to compliance with complex rules and regulations—so owners can focus on growing their business.

“Kinexus is a compelling demonstration of the critical role entrepreneurs play in medical R&D that can help improve the health and lives of Canadians and people around the world,” said Greg Secord, President, ADP Canada, Employer Services Division. “We hope that innovative businesses like Kinexus will continue to come forward and share their inspiring stories in the contest entries.

**PRESS RELEASE - 2013 MARCH 11**

Kinexus currently has agreements with over 1700 research laboratories in companies, universities, government institutions and hospitals in over 35 different countries. To learn more about the diverse proteomics and bioinformatics services offered by Kinexus, please visit <http://www.kinexus.ca>. Kinexus is a private, biotechnology company engaged in the research and development of innovative methods to map, track and manipulate 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 email [csutter@kinexus.ca](mailto:csutter@kinexus.ca) or visit our  
website at [www.kinexus.ca](http://www.kinexus.ca)**