



# KINEXUS

## **Kinexus a Finalists for the "Red Herring 100 Global" Awards 2007**

*The Most Promising Private Technology Ventures in the World to be Celebrated at Event in Seattle, Washington*

**FOR IMMEDIATE RELEASE NOVEMBER 13, 2007**

**San Francisco, CA, November 13, 2007** – Red Herring Magazine announced today that Kinexus Bioinformatics Corporation has been short listed for the Red Herring 100 Global 2007 Awards. Red Herring 100 winners and finalists from North America, Europe and Asia of the last three years were eligible for this award. The Red Herring editorial team used a very competitive process to whittle down this pool of 1800 eligible promising companies to the 200 finalists of this first-time award. Evaluations were made on both quantitative and qualitative criteria such as financial performance, innovation, management, global strategy, and ecosystem integration. The names of the 200 companies short-listed as finalists can be found online at <http://www.herringevents.com/global07/index.html>.

"The 200 finalists we selected from around the world are all excellent contenders," said Joel Dreyfuss, Editor-in-Chief of Red Herring. "They are exceptional companies with smart and innovative platforms that attest to the importance of technology in today's global economy." Kinexus made the final group because of its outstanding achievements, and Red Herring Magazine is honored to announce it is one of the most promising technology firms in the world. In 2004, Red Herring chose Kinexus as one of the top 100 most innovative technology companies in the world, along with such technology giants as IBM, Sun Microsystems, Hewlett Packard, Amazon.Com and Google, and in 2005 as one of the top 100 Private Companies in North America.

The finalists companies will be celebrated at the Red Herring Global event in Seattle, Washington on December 3-5. The 100 winners will be announced at the Award Ceremony on December 4. Red Herring Global is an exclusive event honoring cutting-edge private technology companies from across the world. The event brings together an elite roster of entrepreneurial and global venture investment firms to showcase excellence in innovation. It features keynote discussions, roundtable panels and networking events designed for participants to make connections and learn what it takes to achieve the best worldwide in business and technology. Distinguished speakers at Red Herring Global include Bruce Jaffe, Corporate VP at Microsoft; Promod Haque at Norwest Venture Partners and David Chao, Cofounder and General Partner, DCM; among others.

### **About Red Herring ([www.redherring.com](http://www.redherring.com))**

Red Herring is a global media company uniting the world's best high technology innovators, venture investors and business decision makers in a variety of forums: a leading innovation magazine; an online daily technology news service; technology research, and major events for technology leaders around the globe. Red Herring provides an insider's access to the global innovation economy, featuring unparalleled insights on the emerging technologies driving the economy.

### **About Kinexus Bioinformatics Corporation ([www.kinexus.ca](http://www.kinexus.ca))**

Kinexus Bioinformatics Corporation is a private, biotechnology 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 - 2007 NOVEMBER 13**