



Kinexus Bioinformatics Corporation
www.kinexus.ca call toll free: 1-866-546-3987

KINEXUS



Kinexus Bioinformatics Corporation
maps cell communications systems to
enable diagnostics and therapeutics
discovery and development.
Our antibody-based platform technologies
permit us to track hundreds of key cell
signalling proteins for our industrial and
academic clients.

KINEXUS BIOINFORMATICS

www.kinexus.ca

cell signalling

post genomics

bioinformatics

listening to cells
to silence disease

KINEXUS

SC-300TOP

