

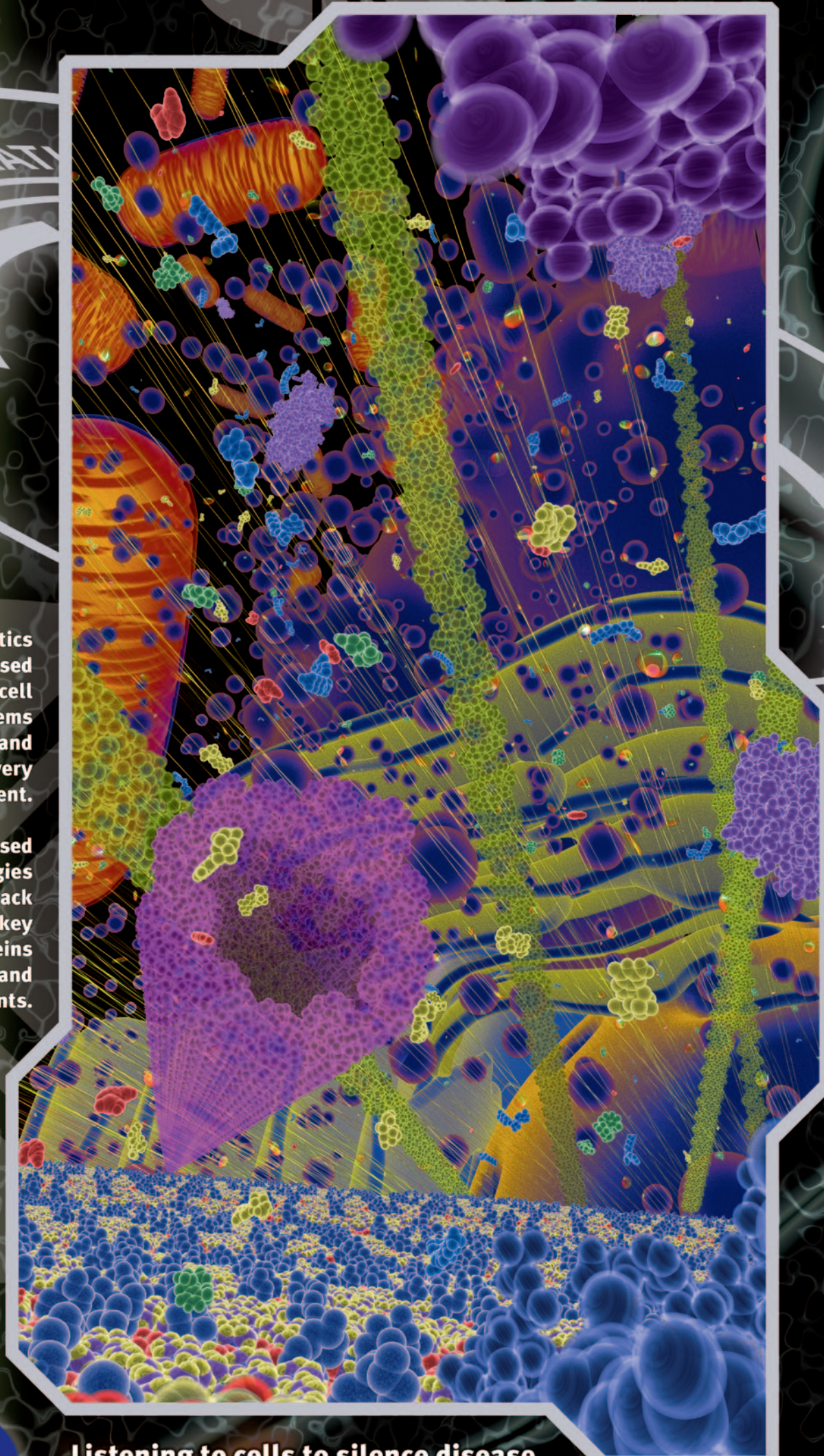
KINEXUS

KINEXUS BIOINFORMATICS

Kinexus Bioinformatics Corporation is focused on mapping 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.

Call toll free:
866-546-3987
info@kinexus.ca



Listening to cells to silence disease.



www.kinexus.ca

Poster design by Cameron Bowyer www.carbon-digital.com. Printed in Canada.