

Crystallography Services Transitioning to New Partner

Carna's collaborative agreement with the German company, Crelux GmbH, will expire on June 30, 2016 after a seven year business relationship.

We are pleased to inform you we are finalizing an Agreement with a new crystallography service provider, enabling us to provide new and expanded services to you our valued customers. With this new collaboration, we look forward to meeting and exceeding your expectations. In selecting our new partner, we have looked for a company who could provide their services at the same high quality on which our reputation has been built.

Should you have any questions or inquiries during this interim period, please feel welcome to contact :

Carna Biosciences Inc.

Marketing and Sales Department

at info@carnabio.com or orders@carnabio.com (US and Canada).