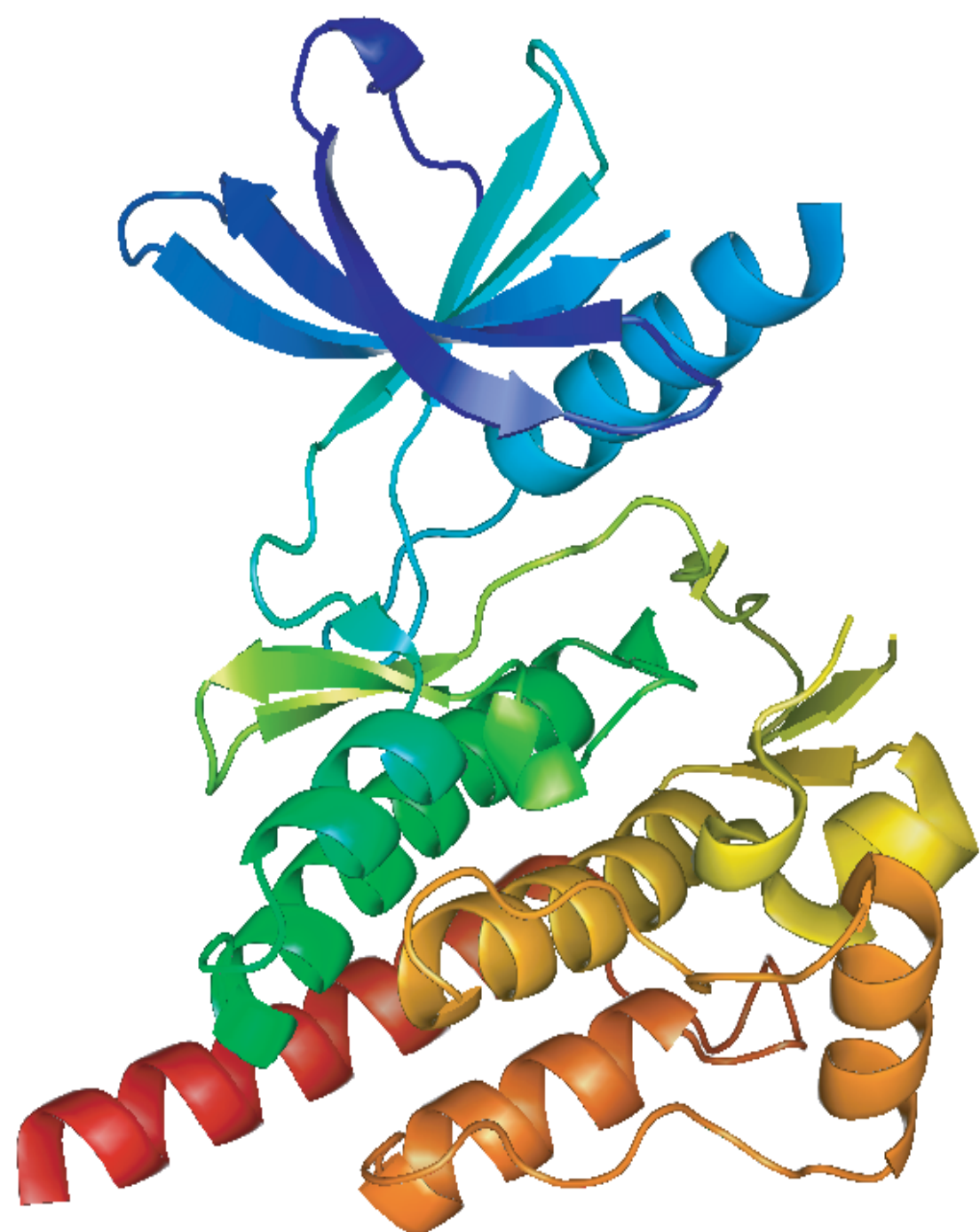


Crystallography

products and services

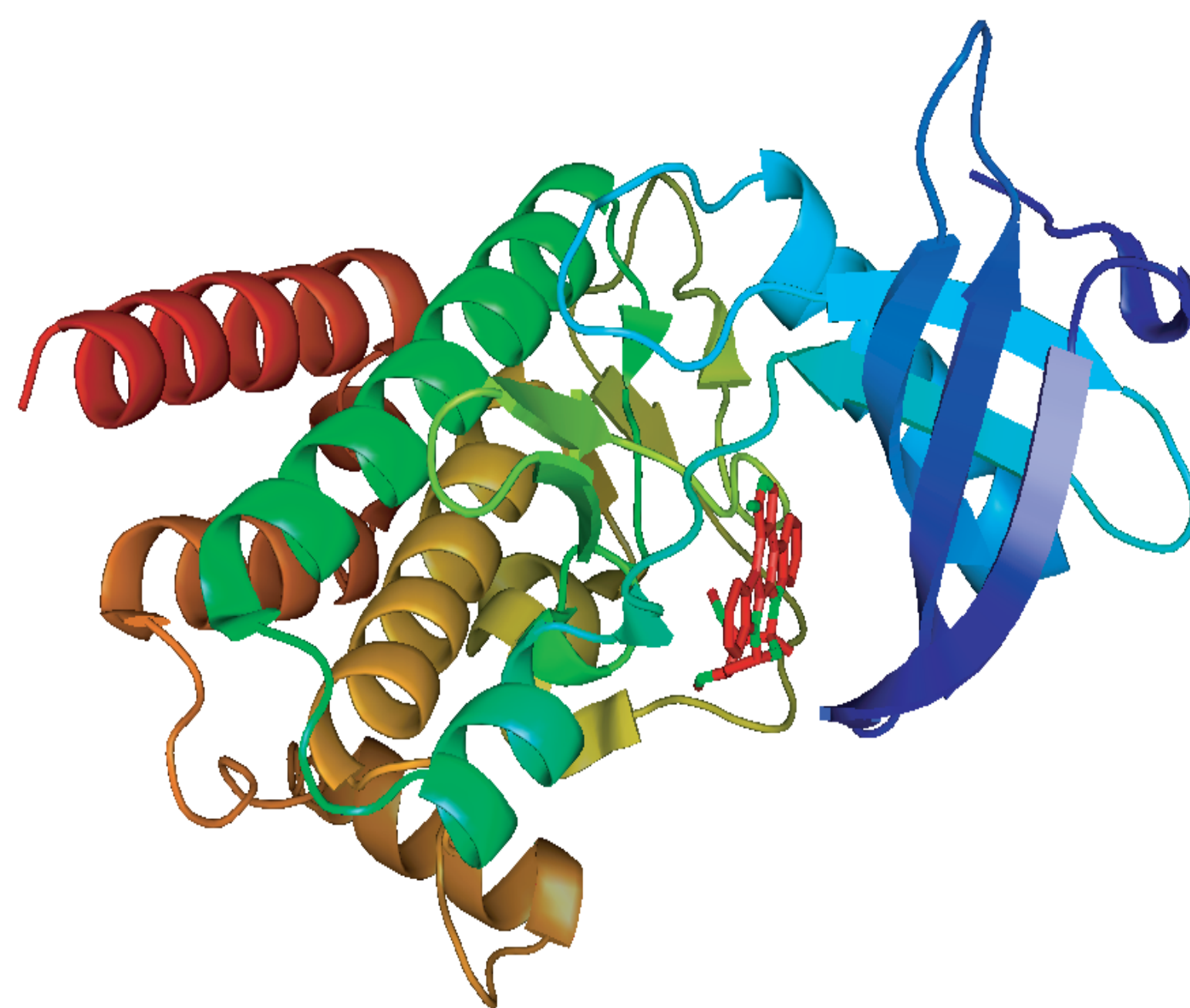
Carna Biosciences is please to announce the availability of crystallization-related products and service through our alliance with **CrystalGenomics, Inc.**



Detailed structural analysis of disease target kinases, in the presence or absence of their putative therapeutic compounds, has become essential for the SBDD (structure-based drug design) processes now favored by many biopharmaceutical companies, particularly those investigating the modulation of known protein kinases. The 3-D structural information of kinases, crystallized alone or co-crystallized with such compounds, provides these biopharmaceutical companies the ability to optimize their compounds rapidly, rationally, and efficiently using our crystallization related products and services.

"Go Co-crystal" service

Detailed 3-D structural information is currently available for 8 protein kinases. With this information, knowledge of the conditions necessary for appropriate crystal growth, and the unique co-crystallization services provided by CrystalGenomics, we can rapidly provide similarly detailed 3-D structural information on any of these protein kinases co-crystallized with your compounds. These procedures can significantly reduce lead optimization-related time and cost requirements, for our biopharmaceutical customers involved in SBDD activities.



Available Kinase List

AurA
Erk1
Erk2

GSK3 β
JNK1
JNK2

JNK3 -available soon-
KDR
CDK4 surrogate

