



Carna Biosciences focuses on kinases. We have substantial know-how covering the areas of gene cloning, expression of active protein kinases, identification of substrates for kinase assays, and developing assay systems for protein kinases suitable for high-throughput screening ("HTS") as well as Kinase Profiling. Through our alliance with CrystalGenomics, we can now offer additional crystallization related products and services.

Our highly experienced group of researchers are taking place in Kobe, Japan, a city known for its state-of-the-art medical technology and industry. We use newest biotechnologies, in order to contribute to the realization of human desire.



Our highly advanced technology can provide a wide variety of drug discovery tools and services to reduce the time and cost significantly, for biopharmaceutical company's drug discovery processes.

Profiling Service

Carna Biosciences offers a broad range of profiling services using our extensive panel of high quality kinase proteins to rapidly and efficiently determine the efficacy, specificity and selectivity of your kinase activity modulator. Also, we are providing custom profiling service, such as IC₅₀ determinations.

Assay Development for HTS

Carna Biosciences will work with biopharmaceutical customers to develop high-throughput screening (HTS) assays to meet their specific requirements. Carna Biosciences' services include cloning the specified protein kinase, expressing and purifying the resulting protein in large quantity, and optimizing assay conditions in a timely fashion (typically 2-to-6 months).

Bulk Kinase Protein

Carna Biosciences has developed expression vectors for more than 500 protein kinases. These vectors are being used to clone and produce bulk quantities for immediate purchase (for list of currently available protein kinases, please contact us).

Crystallography products and services

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

