

## News Release

February 4, 2019  
Carna Biosciences, Inc.

### **Carna Announces Opening of New Clinical Development Office in South San Francisco**

Carna Biosciences, a biopharmaceutical company focusing on the discovery and development of innovative drugs targeting kinases to address serious unmet medical needs in oncology and autoimmune diseases, announces today that CarnaBio USA, Inc, a wholly owned subsidiary of the company, has opened a new office in South San Francisco for its newly established Clinical Development Division to implement overseas clinical trials of drug candidates discovered from our drug discovery programs. Carna Biosciences will commence clinical studies of our drug candidates in the U.S. in cooperation with CarnaBio USA to accelerate the clinical development of new therapies worldwide.

“We are excited to announce the opening of the new office for clinical development in South San Francisco. This marks the important step to achieve our aim since the foundation of the company to conduct clinical studies by ourselves. We believe that the U.S. is the best country to establish a key clinical development function to accelerate the development of our breakthrough drug candidates to treat cancers,” said Dr. Kohichiro Yoshino, President and CEO of Carna Biosciences.

Dr. Masaaki Sawa, Chief Scientific Officer of Carna Biosciences noted, “The San Francisco Bay Area is an ideal location bringing advantage of pharmaceutical industry to conduct clinical trials of our drug candidates. Our new clinical development team will gear up to move AS-1763, a next generation non-covalent BTK inhibitor targeting blood cancer, into first-in-human clinical trial in the U.S.”

#### **About CarnaBio USA Clinical Development Division**

Address: 400 Oyster Point Boulevard, South San Francisco, CA 94080, USA

Date of establishment: February 1, 2019

Purpose: Clinical Development of Carna’s proprietary drug candidates

#### Contact:

Corporate Planning

Carna Biosciences, Inc.

TEL: +81-78-302-7075

<https://www.carnabio.com/english/>