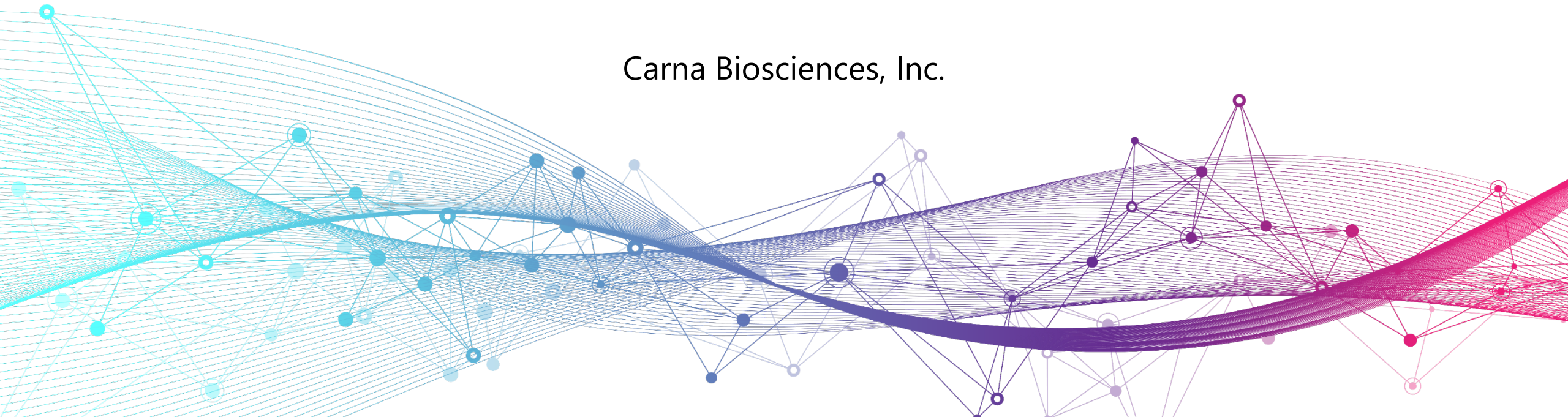


# Financial Results

## Q2 FY2024

(January to June 2024)

Carna Biosciences, Inc.



# AGENDA

1

FY2024 Q2 Key Highlights

2

Updates on Pipelines in Clinical Development

3

Updates on Licensed Pipelines

4

Drug Discovery Support Business

5

FY2024 Q2 Results

6

Appendix



## BTK inhibitor AS-1763

- Phase 1b study is ongoing in the U.S.

Multi-center clinical study

Study Lead : Prof. Nitin Jain, MD, Department of Leukemia,  
University of Texas MD Anderson Cancer Center.

- Safe and well tolerated at the first four dose levels and moved to the dose level 5 in June 2024.
- Encouraging preliminary data from ongoing Phase 1b study was presented at EHA 2024 in June. Ref. P.12-P.13

A favorable safety and PK profile as well as promising efficacy in patients with CLL who have been heavily pretreated with systemic therapies, including covalent BTK inhibitors and BCL2 inhibitor.

- Obtained a patent in China on combination therapy with AS-1763 and BCL-2 inhibitor to strengthen intellectual property protection.



## CDC7 inhibitor monzosertib (AS-0141)

- Phase 1 study in cancer patients is in progress in Japan.

Clinical trial site : National Cancer Center Hospital and National Cancer Center Hospital East

- Presented new preclinical data on monzosertib at AACR in April 2024.

The preclinical data presented in the poster demonstrate the antitumor efficacy of monzosertib alone and in combination with current therapies against human AML cell lines in vitro and in vivo. The new preclinical data support advancement into blood cancer indications.

## DGK $\alpha$ inhibitor : GS-9911 (Out-licensed to Gilead Sciences, Inc.)

- Phase 1 study in patients with solid tumor is ongoing.

## Financing

- Fundraising through the 3rd party allotment of common shares (May)





# Updates on Pipelines in Clinical Development

- 1 AS-1763**
- 2 sofno Brutinib (AS-0871)**
- 3 monzosertib (AS-0141)**

International Nonproprietary Name (INN) : sofno Brutinib, Code name : AS-0871

International Nonproprietary Name (INN) : monzosertib, Code name : AS-0141



Compound	Target	Indication	Status
AS-1763	BTK	Blood Cancer	<ul style="list-style-type: none"><li>Phase 1b dose escalation part is ongoing in the U.S.</li><li>Safe and well tolerated at the first four dose levels and moved to the dose level 5 in June 2024.</li><li>Encouraging preliminary data was presented at EHA 2024 in June.</li><li>Plan to initiate the dose expansion part by the end of 2024, to accelerate development timelines.</li></ul> <div>Multi-center clinical study Study Lead : Prof. Nitin Jain, MD, Department of Leukemia, University of Texas MD Anderson Cancer Center.</div>
sofnobrutinib (AS-0871)	BTK	Immune-inflammatory diseases	<ul style="list-style-type: none"><li>Completed Phase 1 clinical trial in healthy volunteers in the Netherlands, in which a favorable safety and tolerability profile as well as a promising PK/PD profile were confirmed.</li><li>Conducting preclinical studies necessary for Phase 2 trial initiation.</li><li>Seeking a strategic partner for further development.</li></ul>
monzosertib (AS-0141)	CDC7/ ASK	Cancer	<ul style="list-style-type: none"><li>Dose escalation part of Phase 1 study in cancer patients is ongoing in Japan.</li></ul> <div>Clinical trial site : National Cancer Center Hospital and National Cancer Center Hospital East</div>



## Mechanism/ Indication

**Orally available** small molecule inhibitor of Bruton's Tyrosine Kinase (BTK) targeting B-cell malignancies

## To overcome drug resistance

AS-1763 is a **non-covalent inhibitor** that reversibly inhibits BTK, **having a potential to be effective for patients who have developed resistance to the existing BTK inhibitors.**

## To minimize a risk of side effects

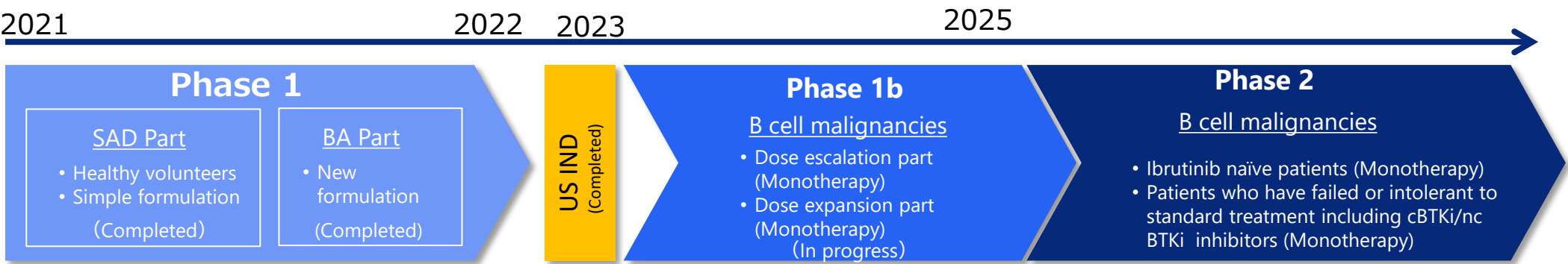
AS-1763 is designed to selectively inhibit BTK **to reduce a risk of potential side effects.**



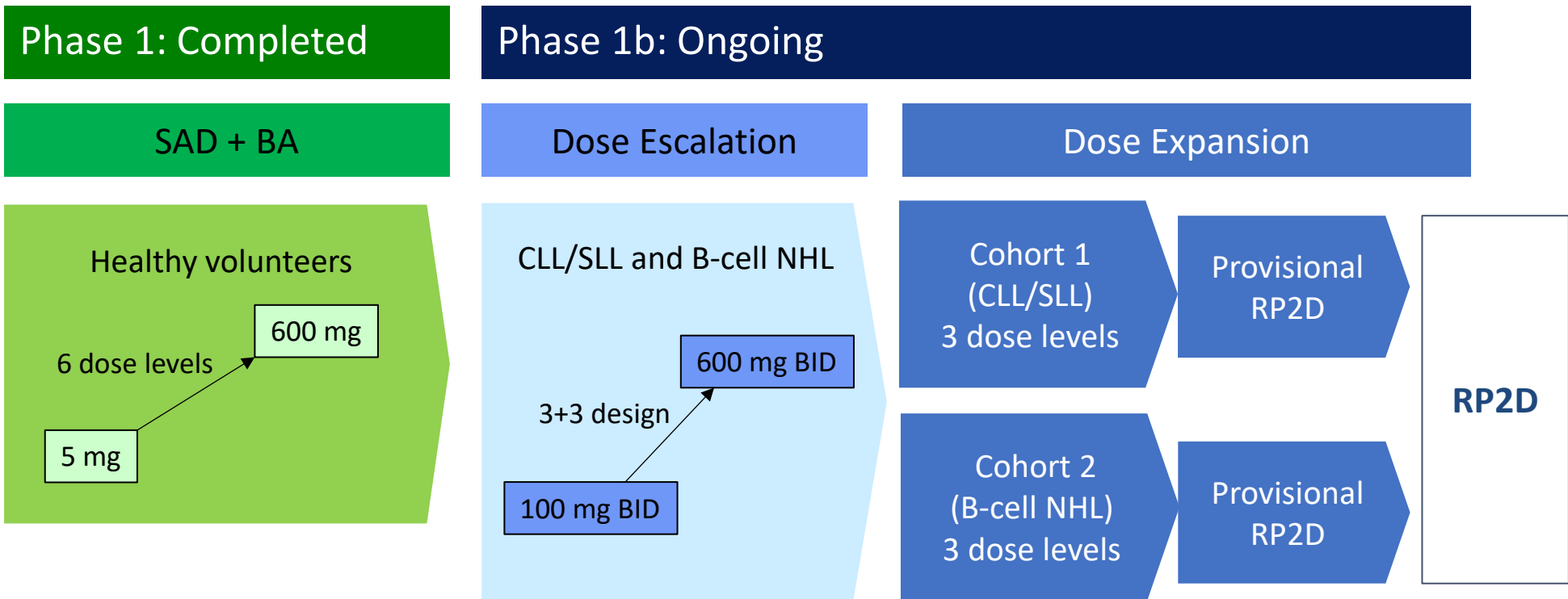
### AS-1763 : Targeting Blood Cancer

- Small molecule BTK inhibitor
- Non-covalent/reversible
- High kinase selectivity
- Pan-mutant BTK inhibitor
- Orally available

- Active against covalent/non-covalent BTK inhibitor-resistant mutations found in patients (C481x, T474x, T316A, L528x)
- The first patient was dosed in August 2023 in Ph 1b study in the U.S
- Dose escalation part is ongoing at the dose level 5 (500 mg BID)



IND application: Investigational New Drug application  
BID: Twice a day  
SAD: Single Ascending Dose  
BA: Bioavailability  
B-cell malignancies: Chronic Lymphocytic Leukemia (CLL), Small Lymphocytic Lymphoma (SLL), and B-cell non-Hodgkin Lymphoma (B-cell NHL), etc.  
cBTKi: covalent BTK inhibitor  
ncBTKi: noncovalent BTK inhibitor



- ◆ Double-blinded, randomized, and placebo-controlled Phase 1 study in the Netherlands
- ◆ Healthy volunteers

- ◆ Open-label, multi-center, Phase 1b study in the U.S.
- ◆ Patients with CLL/SLL or B-cell NHL who have failed or intolerant to at least two lines of systemic therapy
- ◆ Prior therapy with a covalent BTKi is permitted

CLL: Chronic lymphocytic leukemia  
SLL: Small lymphocytic lymphoma  
B-cell NHL: B-cell non-Hodgkin lymphoma  
RP2D: recommended phase 2





## Objectives of the study

**A single dose of AS-1763 was administered orally to healthy volunteers to evaluate:**

- plasma concentration of AS-1763
- safety assessments (clinical laboratory, ECGs, or vital signs, etc.)
- inhibitory potency on B cell activation

## Result of the study

**After a single-dose administration of 5 mg, 25 mg, 100 mg, 300 mg, 500 mg, and 600 mg of AS-1763:**

- plasma concentrations of AS-1763 were increased in a dose dependent manner
- no clinically meaningful issues were reported in all safety assessments
- dose-dependent inhibition in B-cell activation was observed

## New tablet formulation for Phase 1b study

**Relative oral bioavailability was evaluated after administering newly developed tablet formulation containing 100 mg of AS-1763 to healthy volunteers.**

- the new tablet formulation demonstrated good oral bioavailability, supporting that the tablet formulation can be used for Phase 1b study.



**Multi-center clinical study**  
**Study Lead : Prof. Nitin Jain, MD, Department of Leukemia, University of Texas MD Anderson Cancer Center.**

**Clinical trails in progress**

Phase 1b dose escalation part was initiated in the U.S.

**Indication**

Patients with CLL(Chronic Lymphocytic Leukemia), SLL(Small Lymphocytic Leukemia), and B-cell NHL(B-cell non-Hodgkin Lymphoma).

**Status**

- Opened 9 clinical sites.
- Safe and well tolerated at the first four dose levels.
- Moved to the dose level 5 (500 mg BID).

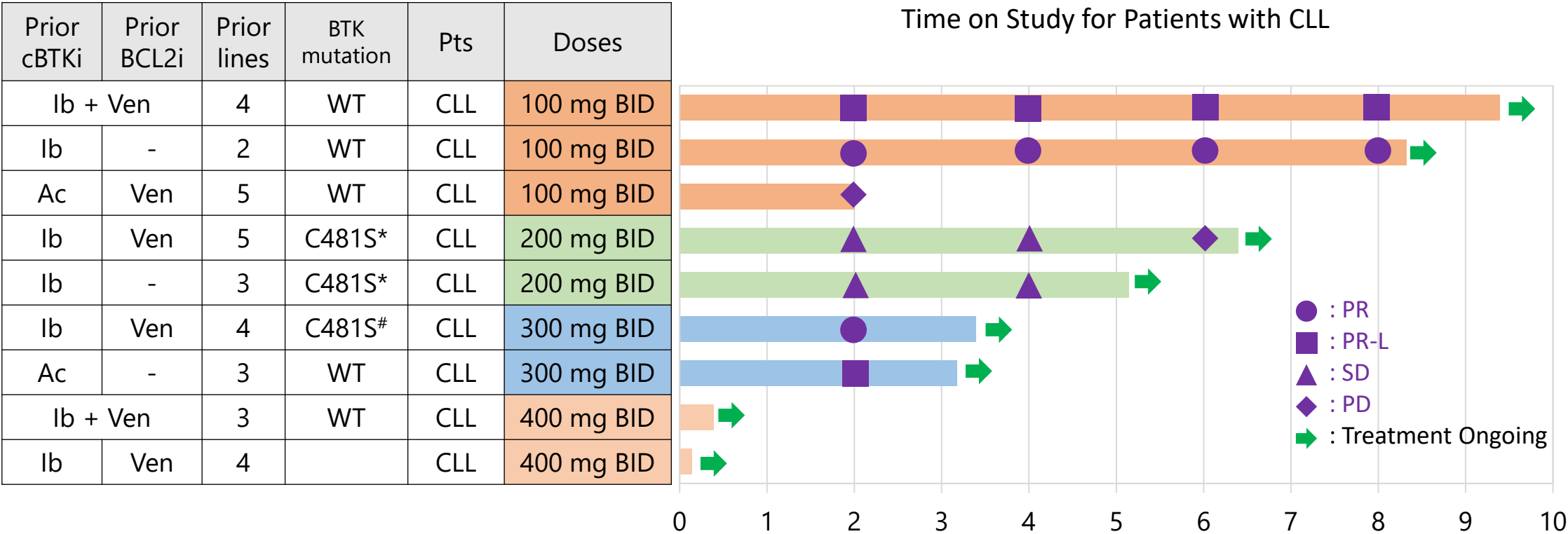


# Preliminary Phase 1b Data in Patients with CLL



AS-1763 demonstrates impressive clinical responses in heavily pretreated patients with CLL, including covalent BTK inhibitors and BCL2 inhibitors.

Data cutoff: 19 April 2024



\*BTK C481S detected centrally by next generation sequencing (CLL mutation panel)

#BTK C481S detected locally by Sanger sequencing

Ib: ibrutinib, Ac: acalabrutinib, Ven: venetoclax, Ib + Ven: combo



- The first patient was dosed in Phase 1b study in August 2023
- As of April 19, 2024, 12 patients (9 CLL, 2 FL, 1 MCL) were enrolled to 4 dose levels (100 mg BID - 400 mg BID)
- No DLTs were observed, and the maximum tolerated dose was not reached yet
- No drug-related atrial fibrillation or bleeding-related events were reported
- At the time of the data cut-off (April 19, 2024), AS-1763 demonstrated an overall response rate of 57% among 7 patients with CLL who have been heavily pretreated with systemic therapies, including cBTKi and BCL2i
  - ✓ 2 of 3 patients (67%) for 100 mg BID and 2 of 2 patients (100%) for 300 mg BID achieved PR or PR-L
  - ✓ 2 patients with CLL at 200 mg BID experienced SD with 16-45% reduction in tumor size
- The dose escalation with higher doses is ongoing

*Based on the promising results, we plan to initiate the dose expansion part later this year*



## Competitors: other non-covalent BTK inhibitors in clinical development

Compound	Company	Development Phase
pirtobrutinib (LOXO-305)	Lilly (Loxo)	Approved/P3
nemtabrutinib (ARQ 531)	Merck (ArQule)	P3
TT-01488	TransThera	P1
HMPL-760	HutchMed	P1

- FDA granted an accelerated approval to pirtobrutinib, the most advanced competitor to AS-1763, in the U.S early 2023.
- Focused differentiation strategy from pirtobrutinib is the key to the success of AS-1763.
- In preclinical studies, AS-1763 demonstrated to be effective for mutant BTKs that confer resistance to pirtobrutinib\* in addition to ibrutinib \*\*.

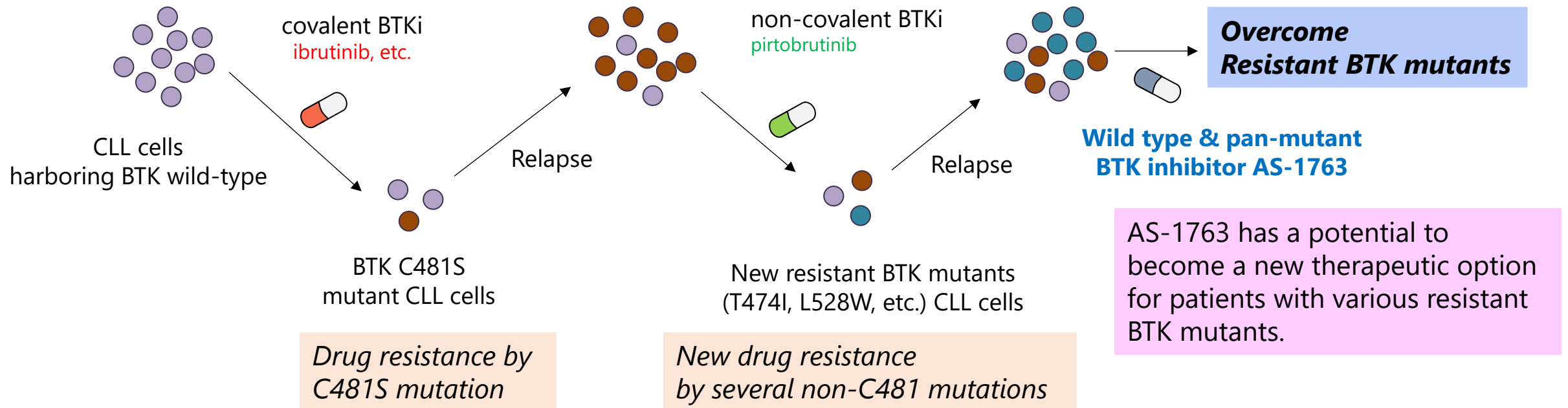
\*N. Engl. J. Med. 2022;386(8):735–743.

\*\*Poster presentation at ASH 2023



# AS-1763: Wild-type & Pan-mutant BTK Inhibitor

- Covalent BTK inhibitors have been widely appreciated as a promising targeted therapy for patients with B-cell malignancies.
- However, patients are reported to develop resistance during the treatment due to substitution of cysteine residue at 481 position with serine (C481S mutation) in BTK, which reduces the efficacy of the covalent BTK inhibitors.
- In addition, the emergence of other types of resistance mutations to non-covalent BTK inhibitor, pirtobrutinib approved in 2023, has been reported.
- AS-1763 potently inhibited both wild type and those mutant BTKs, strongly suggesting that AS-1763 will be a new therapeutic option for treating patients with B-cell malignancies both having wild type and resistance mutations in BTK.





Present

Covalent BTK inhibitors including ibrutinib, acalabrutinib, and zanubrutinib are key therapeutic options.

Opportunity

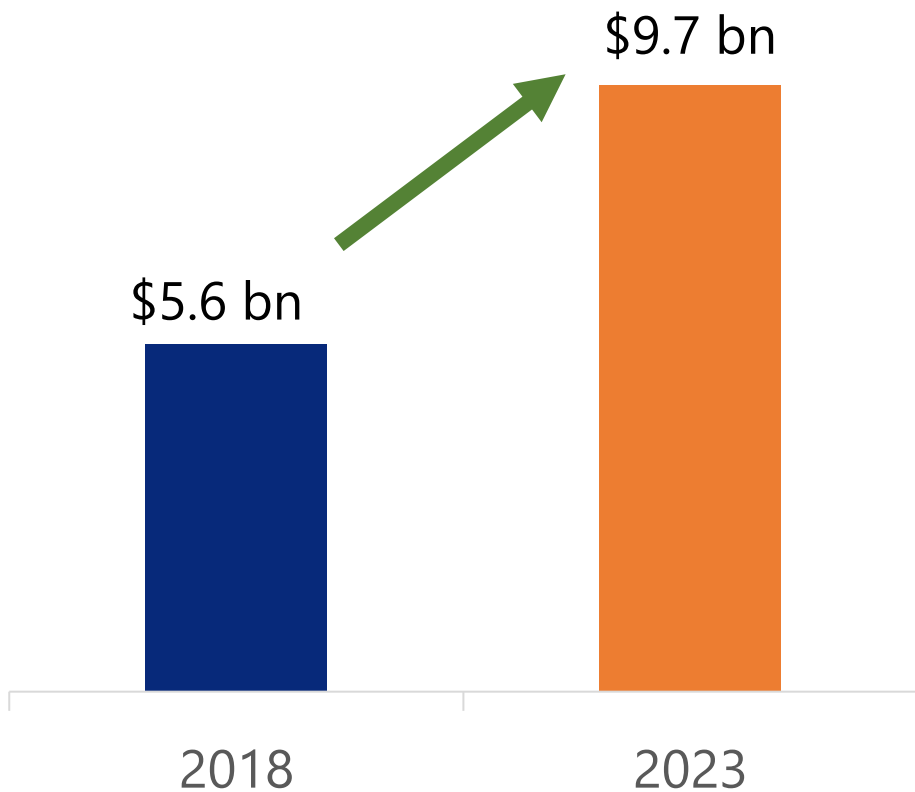
Patients treated with covalent BTK inhibitors are reported to develop resistance to the drugs



High unmet needs for new therapeutic options to treat patients with B-cell malignancies who have developed resistance to covalent BTK inhibitors



- Next generation non-covalent BTK inhibitors including AS-1763 are expected to be promising therapeutic options for the patients who have developed resistance to covalent BTK inhibitors. Our goal is to take a certain share of this potential large market, making AS-1763 a blockbuster drug.



BTK Inhibitors WW sales  
(1st generation/covalent type)

The sales growth of first-generation BTK inhibitors has slowed due to potential side effects and drug-resistance by C481S mutation.



Next generation BTK inhibitors including AS-1763 are under development to overcome side effects and drug-resistance.

*Global BTK inhibitors market size is still expanding significantly and expected to reach > \$22Bn by 2030!*

<https://www.insightaceanalytic.com>



## Mechanism/ Indication

Sofnobrutinib is an **orally available** Bruton's Tyrosine Kinase (BTK) inhibitor to treat autoimmune diseases by inhibiting activation of immune cells such as B cells, macrophages, and mast cells.

## To minimize a risk of side effects

Sofnobrutinib is designed to selectively inhibit BTK **to reduce a risk of potential side effects.**

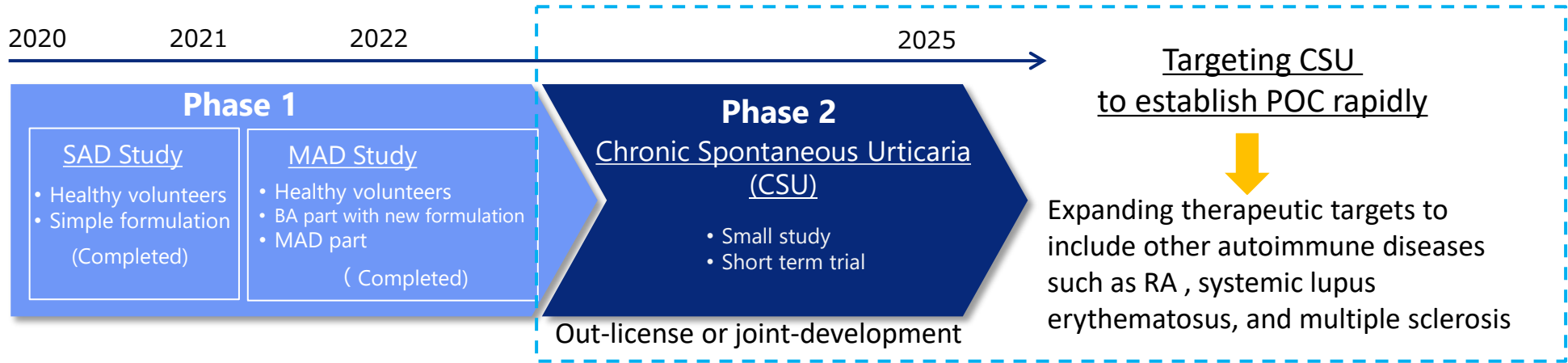
## Characteristics

Sofnobrutinib is a **non-covalent BTK inhibitor** that reversibly inhibits BTKs to reduce safety concerns associated with covalent inhibitors.



## Sofnobrutinib (AS-0871) : Targeting Immune-inflammatory diseases

- |   |  |
|---|--|
| <ul style="list-style-type: none"><li>• Small molecule BTK inhibitor</li><li>• Non-covalent/reversible</li><li>• High kinase selectivity</li><li>• Orally available</li></ul> | <ul style="list-style-type: none"><li>• Demonstrated significant efficacies in arthritis models</li><li>• Showed efficacy in systemic lupus erythematosus model</li><li>• Phase 1 Clinical Trial was completed</li><li>• Find a partner to conduct further development</li></ul> |
|---|--|



SAD: Single Ascending Dose  
MAD: Multiple Ascending Dose  
BA: Bioavailability  
POC: Proof of Concept





Completed

## Phase 1 in the Netherlands SAD study (Healthy volunteers)

- Safe and well-tolerated at all dose levels
- Favorable Pharmacokinetic (PK) Profile
- Promising Pharmacodynamic(PD) effects
- Conducted using simple formulation



Developed multiple new formulations



## Phase 1 in the Netherlands MAD study (Healthy volunteers)

**BA Part**

Evaluate the relative bioavailability of multiple new formulations to select the best formulation

Completed



**MAD Part**

Evaluate the safety, tolerability, PK and PD in the 2-week multiple ascending dose of sofno Brutinib .

Completed



## Objectives of Single Ascending Dose (SAD) study

Ref. P.53-P.55

**A single dose of Sofnobrutinib was administered orally to healthy volunteers to evaluate:**

- plasma concentration of sofno Brutinib
- safety assessments (clinical laboratory, ECGs, or vital signs, etc.)
- inhibitory potency on B cell and basophil activation

## Result of SAD study

**After a single-dose administration of 5 mg, 25 mg, 100 mg, 300 mg, 600 mg, and 900 mg of sofno Brutinib:**

- plasma concentration of AS-0871 were increased in a dose dependent manner.
- no clinically meaningful issues were reported in all safety assessments.
- sufficient inhibition of B cell and basophil activations was observed at 100 mg or above.

## New formulations for Multiple Ascending Dose (MAD) study

**Plasma concentration was evaluated after administering newly developed capsule formulation and tablet formulation containing 50 mg of sofno Brutinib to healthy volunteers:**

- the new tablet formulation was shown to be safe and demonstrated good oral bioavailability, and the tablet formulation was selected for the MAD part.



### Multiple Ascending Dose (MAD) study MAD part design

- Double blinded, placebo-controlled, randomized multiple ascending dose study in healthy volunteers
- 14-day multiple oral doses of sofno Brutinib tablets in 3 cohorts (50, 150 or 300 mg twice daily)
- The safety, tolerability, PK and PD were evaluated.

### Result of MAD study MAD part

- Well tolerated with no dose-limiting adverse events (AEs): AEs were mostly mild.
- Favorable safety profile up to 300 mg twice daily
- Favorable PK profile: approximately dose proportional with increasing doses
- Robust PD effect was observed: over 90% inhibition for basophils activation on Day 14 at 150 and 300 mg BID doses.

**These results from the Phase 1 studies of sofno Brutinib support to advance sofno Brutinib into Phase 2 clinical development for further investigations.**



# Sofnobrutinib: CSU is a skin disease with unmet medical needs

**Chronic Spontaneous Urticaria (CSU)** is a distressing skin disorder that is characterized by itching and hives lasting for more than 4 weeks with unknown causes. The symptoms can last months or years, affecting QoL of patients.

## Challenges of CSU

- A significant number of patients having uncontrolled CSU by existing drugs.
- High socio-economic costs for patients with high disease activity\*
- Large number of patients; approximately 1% of the global population is affected\*

**High unmet medical needs with potential large market**

\* Br J Dermatol 2021;184:226-36.

## Competitors

Compound	Company	Development Phase
Remibrutinib (LOU064)	Novartis	P3

The Phase III trials met the primary endpoints and showed rapid symptom control in CSU, supporting the potential of BTK inhibitors as a new treatment option for those uncontrolled by first-line H1-antihistamines.\*

\* <https://www.novartis.com/news>

## Opportunity

- Approval of new treatment options may trigger the expansion of CSU market.
- We plan to pursue the clinical implications of sofno Brutinib (non-covalent BTK inhibitor) to provide clinical benefits by minimizing potential adverse events associated with covalent BTK inhibitors including remibrutinib.

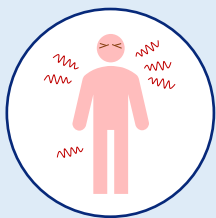


# Chronic Spontaneous Urticaria (CSU)

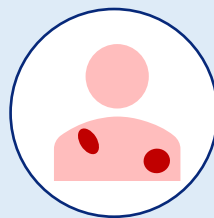
CSU is a debilitating disease of chronic itch, hives and angioedema, lasting six weeks or more.

## Symptoms

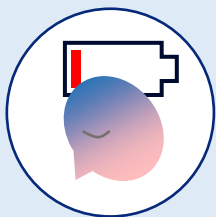
There is no specific external trigger for CSU, but the autoimmune system may play a role.



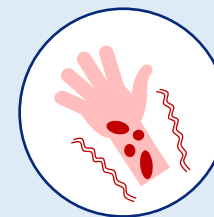
Spontaneously present & re-occur



Red swollen hives



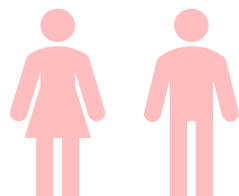
Lack of Energy  
Depression/Anxiety  
Chronic (Lasting for at least six weeks)



Itch

- ✓ Approximately 50% of CSU patients don't respond to H1-antihistamine.
- ✓ Curative treatment is not available.
- ✓ High socio-economic costs for patients with high disease activity.

## Number of Patients

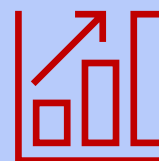


**2.8 million**

diagnosed prevalent cases in major seven markets

- ✓ Approximately 1% of the population worldwide is affected.

## Market Size



**\$1,315 million**

in 2021 in seven major markets

- ✓ The market size of CSU in major seven countries is expected to reach \$7,664 mn by 2032 growing at a CAGR of 14.96% from 2019.

<https://www.delveinsight.com/>

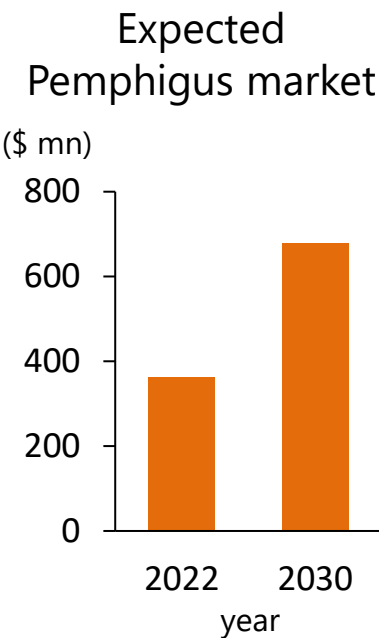
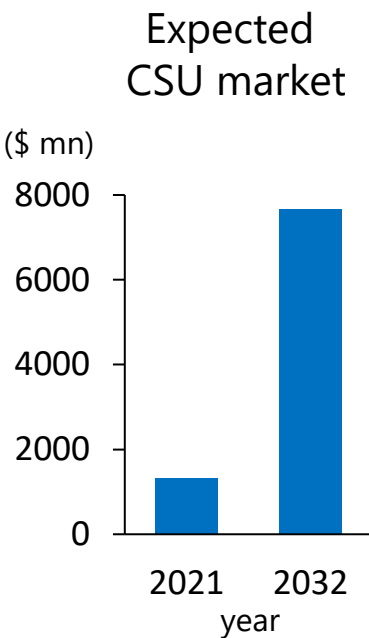




Initial focus

Diseases	Number of patients
CSU	<ul style="list-style-type: none"><li>Diagnosed prevalent cases : 2.8 mn*</li><li>WW population affected: 76 mn</li></ul>
Pemphigus	<ul style="list-style-type: none"><li>Diagnosed prevalent cases : 40,000*</li></ul>

\*in major 7 markets



Other potential therapeutic area

Diseases	Number of patients	Market size in value
Systemic lupus erythematosus (SLE)	Global SLE prevalence is estimated to be 15.87 to 108.92 per 100,000 people	expected to reach \$3,517 mn by 2030
Multiple sclerosis (MS)	In 2016, an estimated 2.2 million people worldwide had MS, corresponding to a prevalence of 30.1 cases per 100,000 population	expected to reach \$34 bn by 2031
Rheumatoid arthritis (RA)	18 million people worldwide were living with RA	expected to reach \$70 bn by 2030

<https://www.delveinsight.com/>  
<https://www.databridgemarketresearch.com/>  
<https://ard.bmj.com/>  
<https://straitresearch.com/>  
<https://www.skyquestt.com/>  
<https://www.who.int/>  
Ann Rheum Dis 2023;82:351–356  
Lancet Neurol 2019 ; 18: 269–85



## Mechanism/ Indication

Monzosertib is an **orally available** CDC7 kinase inhibitor targeting cancers.

## To minimize a risk of side effects

Monzosertib is designed to selectively inhibit CDC7 kinase **to reduce a risk of potential side effects.**

## Potentially effective for various cancers

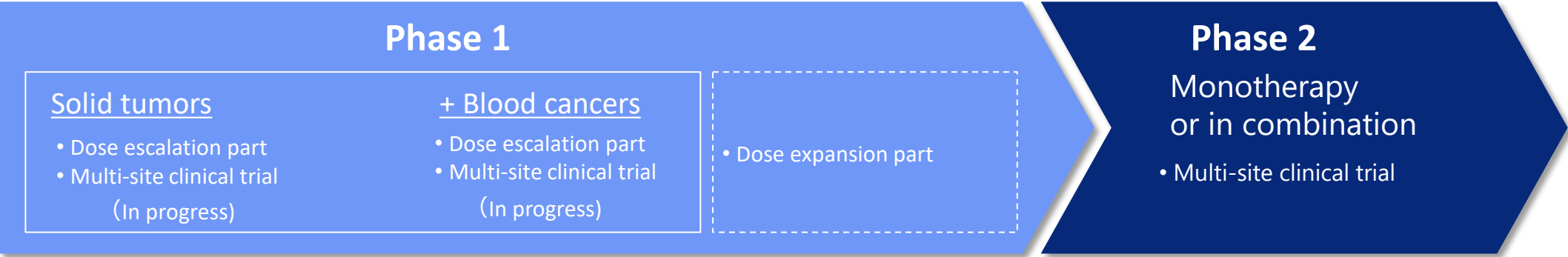
Monzosertib exhibited **a potent anti-proliferative activity against a wide range of cancer cell lines in preclinical studies.**

## Potential first-in-class molecule

Monzosertib has a potential to become a first-in-class drug as no CDC7 inhibitors have been approved.



Monzosertib (AS-0141) : Targeting Cancer	
<ul style="list-style-type: none"><li>● Small molecule CDC7 inhibitor</li><li>● High kinase selectivity</li><li>● Potential First-in-class drug</li><li>● Orally available</li></ul>	<ul style="list-style-type: none"><li>● Potent anti-proliferative activity against various cancer cell lines</li><li>● Demonstrated strong anti-tumor activity in several human tumor xenograft models</li><li>● Conducting Phase 1 study in Japan targeting solid tumors and blood cancers</li></ul>





# Monzosertib (AS-0141): Phase 1 Clinical Trials



Clinical trial sites : National Cancer Center Hospital and National Cancer Center Hospital East

**Clinical trails in progress**

Phase 1 dose escalation study targeting cancer patients is ongoing.

**Objectives of the study**

The primary objectives of the dose escalation study is to assess safety, tolerability, maximum tolerated dose (MTD), preliminary anti-tumor activity, and pharmacokinetics (plasma concentration, duration) of monzosertib.

**Dosage**

Oral administration, twice a day

**Status**

- 80 mg BID (5d on/2d off) was well-tolerated and safe.
- Switched to a continuous dosing schedule (without drug holiday) to maximize efficacy. Safety and tolerability were confirmed at the first dose level (50 mg BID) and dosing at the second dose level (80 mg BID) is ongoing.
- Expanded the target indications to blood cancers to exploratorily investigate the efficacy.



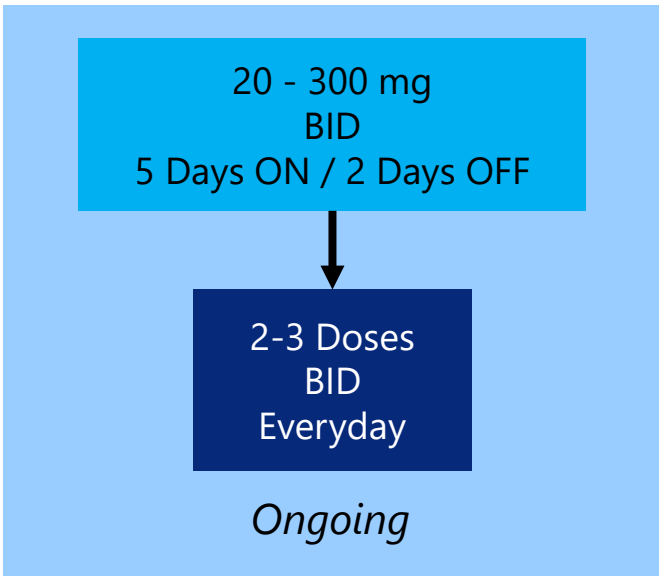
# Monzosertib (AS-0141): Phase 1 Clinical Trial



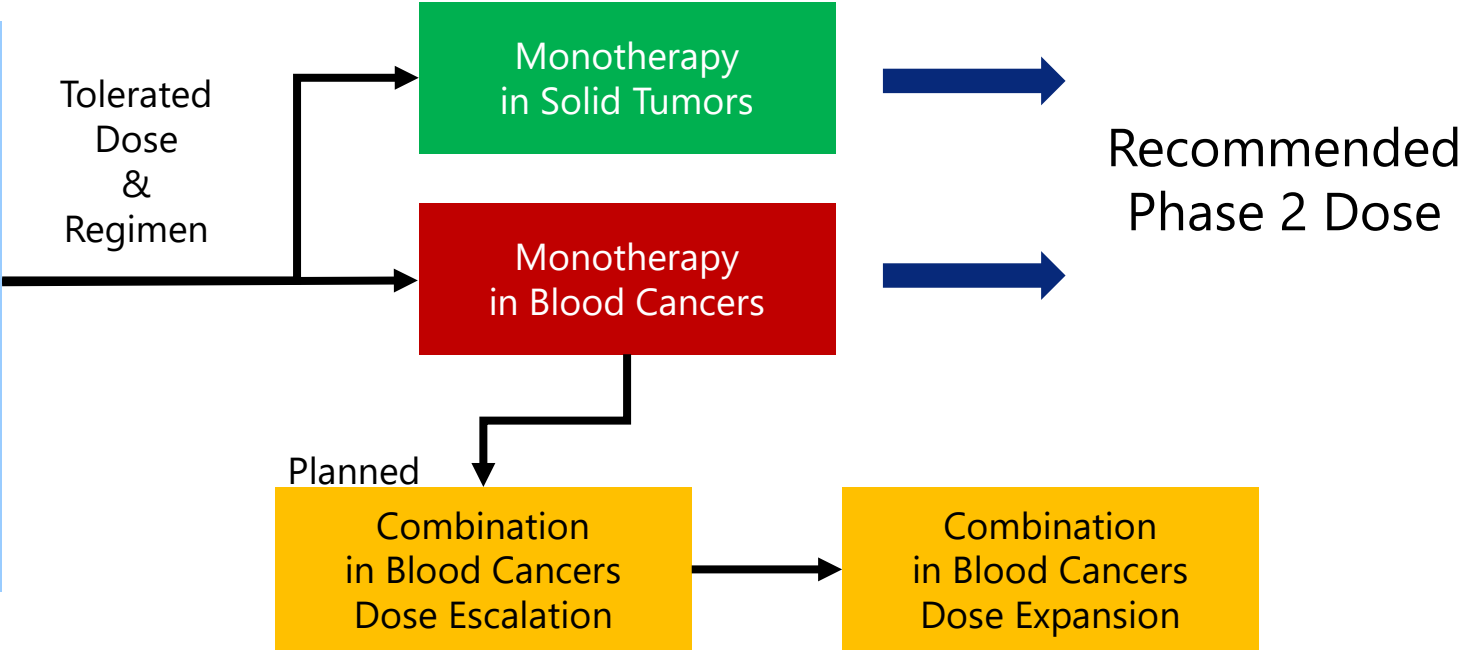
## Phase 1 study targeting cancer patients

- The Phase 1 study in patients with advanced, metastatic, relapsed or refractory malignancies. The protocol has been changed to include patients with blood cancers.
- The study consists of two parts, a dose escalation and an expansion.
- The primary objective is to assess safety, tolerability, maximum tolerated dose (MTD), preliminary anti-tumor activity, and PK / PD as well as to determine RP2D.
- Pharmacodynamic effect of monzosertib was confirmed at 40 mg BID and above.

### Dose Escalation Part



### Dose Expansion Part





*Switched to a continuous dosing schedule (without drug holiday) allowing persistent inhibition of CDC7 to maximize efficacy*

- 50 mg BID (every day) was well-tolerated and safe
- 3 of 4 efficacy-evaluable patients (50 mg BID and 80 mg BID) achieved SD
- One patient receiving 50 mg BID (every day) is still on treatment (>200 days)

Doses	Regimen	n	DLT
50 mg BID	every day	3	0
80 mg BID	every day	3	DLT assessment period
100 mg BID	every day	planned	

Dose Escalation with “every day” regimen is on-going



# Updates on Licensed Pipelines

- 1 DGK $\alpha$  Inhibitor** (Gilead Sciences, Inc.)
- 2 Joint Research with Sumitomo Pharma**



# Out-licensed Programs

Program/ Partner	Compound (Target)	Status	Upfront payment	Total milestone payments expected	Royalty	Region	Contract date	Milestones received
<b>DGK<math>\alpha</math> inhibitor</b> Gilead Sciences (Out-license)	GS-9911 (Immuno- oncology)	Phase 1	\$20M	\$450M	Undisclosed	Worldwide	Jun. 2019	Received milestones twice, totaling \$15M
Joint Research with Sumitomo Pharma	Kinase inhibitor (Psychiatric and neurological disorders)	Late discovery	JPY80M (including research milestone)	JPY10.6B	Undisclosed	Worldwide	Mar. 2018	





Partner



GILEAD

**Gilead Sciences, Inc.**

**Out-licensed in June 2019** (worldwide rights)

Deal size

- Upfront payment \$20 million
- Maximum of \$450 million potential milestone payments upon achievement of certain development and commercial milestones

Royalties

- Royalties on future net sales

## 1. Investigational DGK $\alpha$ inhibitor: GS-9911

## 2. Indication: Cancer (immunotherapy)

- At JP Morgan Annual Healthcare Conference held in January 2024, Gilead introduced DGK $\alpha$  as next generation target in oncology and presented GS-9911 as the DGK $\alpha$  inhibitor in Phase 1 trials.

## 3. The Phase 1 study in patients with solid tumors is ongoing.

Refer Gilead's website for details of the study.

<https://www.gileadclinicaltrials.com/study?nctid=NCT06082960>



Partner



**Sumitomo Pharma Co., Ltd.**  
**Joint Research Agreement in March 2018**  
(worldwide rights)

Deal size	<ul style="list-style-type: none"><li>• Upfront payment + Research milestone JPY80 million</li><li>• Maximum of JPY10.6 billion potential milestone payments upon achievement of certain development and commercial milestones</li></ul>
Royalties	<ul style="list-style-type: none"><li>• Royalties on future net sales</li></ul>

- 1. Joint research to discover novel kinase inhibitors to treat psychiatric and neurological disorders.**
- 2. The term of the joint research was extended in December 2021.**
- 3. Joint research is ongoing to identify preclinical candidates.**



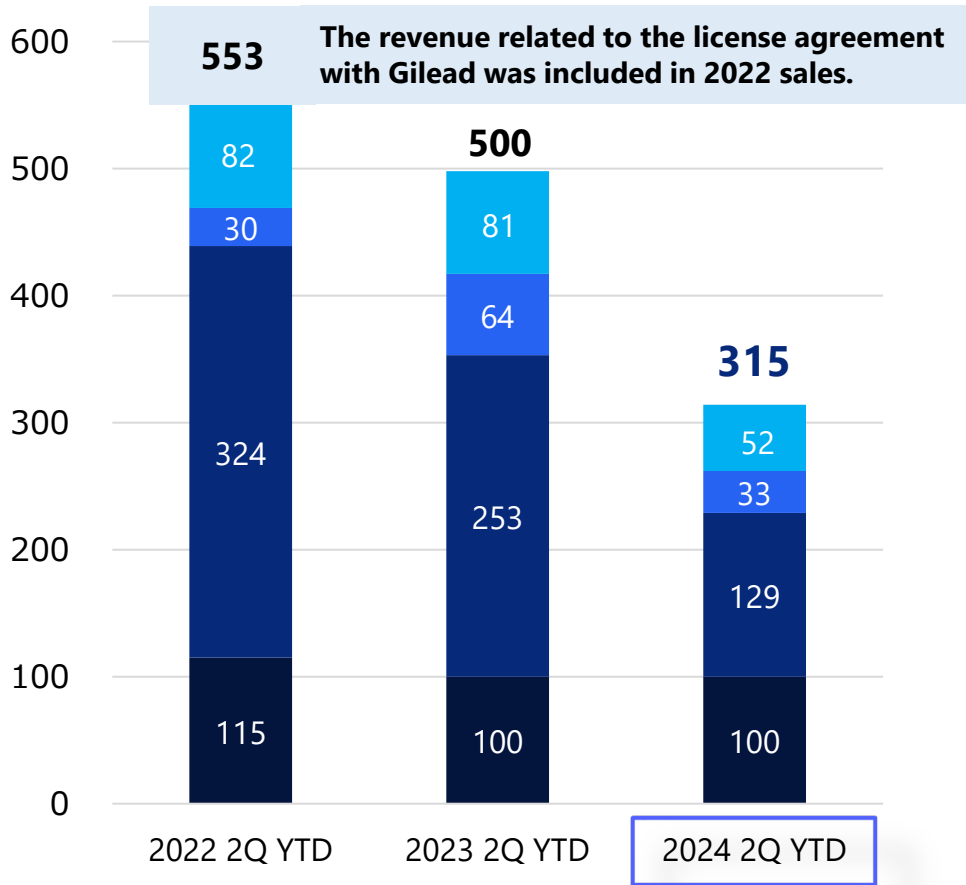
# **Drug Discovery Support (ddSP) Business**



# FY2024 Q2 YTD Drug Discovery Support Business Sales Trend

Drug Discovery Support Business Sales Trend by Region  
(Consolidated)

(JPY million)    Other   Europe   North America   Japan



Japan

## Decreased 0.2% YoY

- Sales of proteins to pharmaceutical companies and distributors were robust, overall sales remained at the same level YoY.

North America

## Decreased 49.1% YoY

- Sales decreased YoY. Accompanied with our major customers' projects progress, the needs of kinase protein declined.
- Sales of profiling services to AI-driven drug discovery companies remained strong.

Europe

## Decreased 48.6% YoY

- Overall sales declined YoY due to the progress of research of our major customers.

Other

## Decreased 35.3% YoY

- The economic recession and the possibility of restraint of trade with Chinese biotech companies by the U.S. government had negative impacts on Chinese market and decreased demand from CROs.



## Profiling Services

- We have successfully developed unique profiling system using the combination of Sciex BioPhase8800 of 8 capillary electrophoresis device, and the robot arm with stacker. We launched this profiling service in May 2024.
  - As the sole provider capable of offering kinase assay services using the reliable Mobility Shift Assay system, we are committed to providing continuous high-quality data.
  - With this cutting-edge system, we aim to attract more customers.
- 
- Currently, we are focusing on promoting awareness of our new system.
  - We have released a promotional video at conferences, on our website, and through social media.
  - It is now available at : [https://www.carnabio.com/common/images/product/video\\_top\\_msa.mp4?240725](https://www.carnabio.com/common/images/product/video_top_msa.mp4?240725)
  - Our customers highly appreciate our continuous service provision with mobility shift assay system and data continuity.

\*At the end of 2024, the support for current mobility shift assay system (EZReader) will be discontinued.



Figure1 BioPhase8800 and the robot arm with stacker

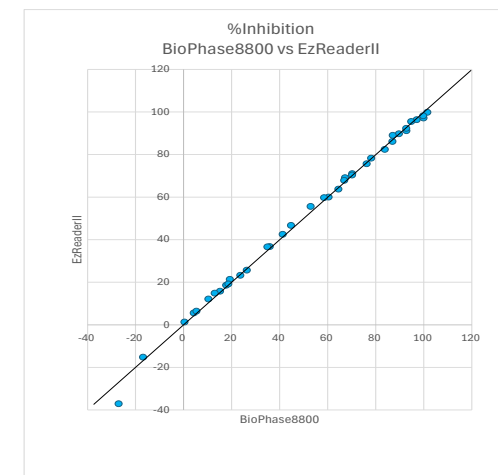


Figure2 This plot shows the correlation between BioPhase8800 and EzReaderII





### Biotinylated Protein Kinases

- The combination of LSA<sup>XT</sup>, a novel system developed by Carterra, for efficient screening of antibodies or small molecules binding to proteins, great compatibility with Carna's biotinylated kinase proteins, has been demonstrated to allow simultaneous measurement of binding interactions between approximately 100 kinases and small molecules in a short time. Furthermore, this system enables quantitative analysis of binding reaction rates and selectivity between small molecules and proteins. In February, we presented these findings at SLAS 2024, leading to the acquisition of new customers.
- Currently, we are conducting experiments using novel combinations of kinases and small molecules. In addition to traditional kinase inhibitors, we are evaluating PROTAC, a protein degradation inducer that has recently gained attention, as well as assessing our biotinylated kinase proteins for even smaller fragment compounds.
- We plan to present these results at a conference scheduled for this fall.



As the use of Carterra LSA<sup>XT</sup> for screening becomes more widespread in the market, we anticipate an expansion in sales of our biotinylated kinase proteins.



# FY2024 Q2 Results





# FY2024 Q2 Results by Business Segment



(JPY million)	Q2FY2023 Actual	Q2FY2024 Actual	YoY Change	FY2024 Plan	
Total Sales	500	315	-184 -36.9%	925	
ddSP business	500	315	-184 -36.9%	925	• Overall sales declined due to weaker than expected overseas sales while sales in Japan remained solid.
ddRD business	—	—	—	—	
Total Operating Profit/Loss	-863	-1,095	-231	-2,201	
ddSP business	169	-24	-193 -114.4%	229	
ddRD business	-1,033	-1,070	-37	-2,431	• Continued investment in the clinical-stage programs.
Ordinary Loss	-868	-1,087	-218	-2,208	
Net Loss	-885	-1,094	-208	-2,225	
R&D cost	959	992	+32	2,309	• Continued investment in the clinical-stage programs including costs related to clinical studies and manufacturing of investigational new drugs for AS-1763 and monzosertib (AS-0141)

Business plan for FY2024 dose not include potential milestone payments or upfront payments as the timing or the amounts are difficult to predict.

Note : Rounded down to the nearest million yen



# Consolidated Balance Sheet

(JPN million)	As of Dec. 31,2023	As of Jun. 30,2024	Change	Reason for changes
Current assets	4,191	3,544	-646	<ul style="list-style-type: none"><li>Cash and deposits +136</li><li>Accounts receivable-trade -689</li></ul>
Cash and deposits	2,889	<b>3,026</b>	+136	
Non-current Assets	158	141	-17	
Total assets	4,349	3,685	-664	
Current liabilities	375	370	-5	<ul style="list-style-type: none"><li>Long-term loans payable within 1 year -49</li><li>Accounts payable +67</li></ul>
Non-current liabilities	96	99	+3	
Total liabilities	472	470	-2	
Total net assets	3,877	3,215	-662	<ul style="list-style-type: none"><li>Capital stock and capital surplus +364</li><li>Retained earnings -1,094</li></ul>
Total liabilities and net assets	4,349	3,685	-664	
Shareholders' equity ratio	89.1%	87.2%		
BPS	226.16yen	178.16yen		
PBR	2.3x	2.2x		
Share price of Carna	522yen	390yen		

Note: Share price is the closing price of the term end.



## Conducted a fundraising through the 3rd party allotment of common shares (May)

- Settlement date      May 31<sup>st</sup>, 2024
- Allotee                Athos Asia Event Driven Master Fund
- Net proceeds        JPY338,570,000
- Use of proceeds     Investment in development of AS-1763 and monzosertib (AS-0141)

### Financing

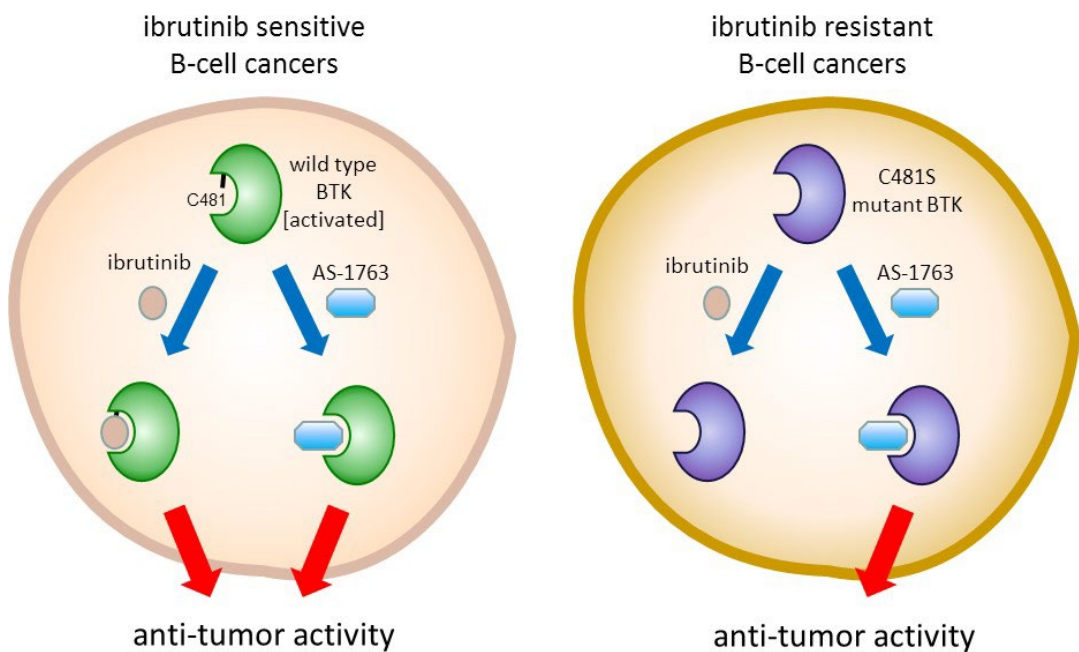
Additional financing may be considered as necessary in order to accelerate the clinical trials of AS-1763, our most important asset.



# Appendix



# AS-1763: Potent Inhibitor of C481S mutant BTK



## Discovery of AS-1763: A Potent, Selective, Noncovalent, and Orally Available Inhibitor of Bruton's Tyrosine Kinase

Wataru Kawahata,\* Tokiko Asami, Takao Kiyoi, Takayuki Irie, Shigeki Kashimoto, Hatsuo Furuichi, and Masaaki Sawa

Cite This: *J. Med. Chem.* 2021, 64, 14129–14141

Read Online

### ◆ IC<sub>50</sub> values of AS-1763 against wild-type and C481S-mutant BTK

	IC <sub>50</sub> (nM)	
	BTK[A]	BTK <sup>C481S</sup>
AS-1763	0.85	0.99

*J Med Chem.* 2021 Oct 14;64(19):14129-14141.





# AS-1763: Strong Cellular Activity and High Kinase Selectivity

## ◆ In vitro pharmacological activities of AS-1763

	IC <sub>50</sub> (nM)	
	AS-1763	Ibrutinib
Autophosphorylation BTK (Ramos)	1.4	1.1
CD69 activation (Human whole blood)	11	8.1
Cancer cell growth OCI-Ly10 cells	1.8	0.75
Cancer cell growth OCI-Ly10 [BTK C481S] cells	20	1030
Normal cell growth HEL299 cells	6370	6870

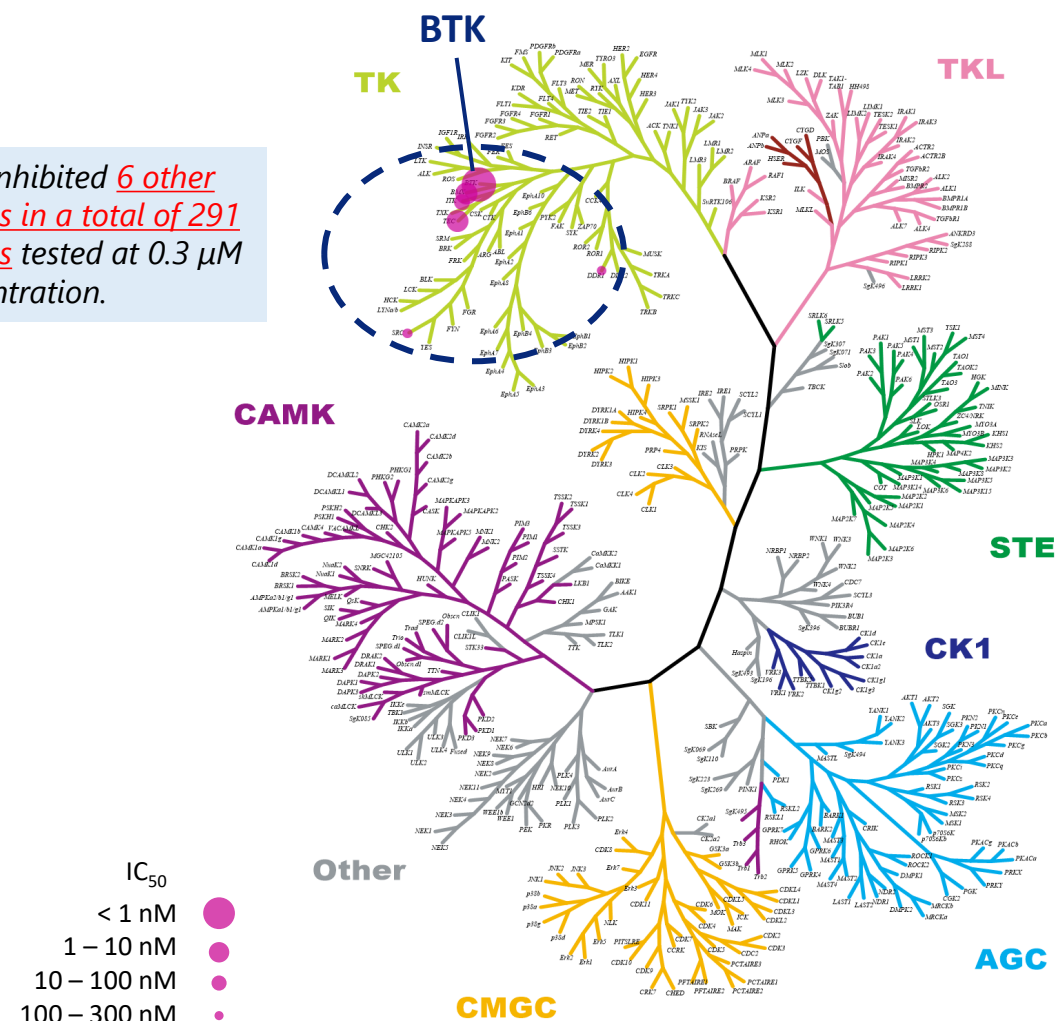
50-fold Stronger activity

Ramos: human Burkitt lymphoma cell line  
OCI-Ly10: human B-cell non-Hodgkin lymphoma cell line  
OCI-Ly10 [BTK C481S]: BTK[C481S] knock-in OCI-Ly10 cells  
HEL299: human embryo lung cell line

J Med Chem. 2021 Oct 14;64(19):14129-14141.

## ◆ Kinase selectivity profiling

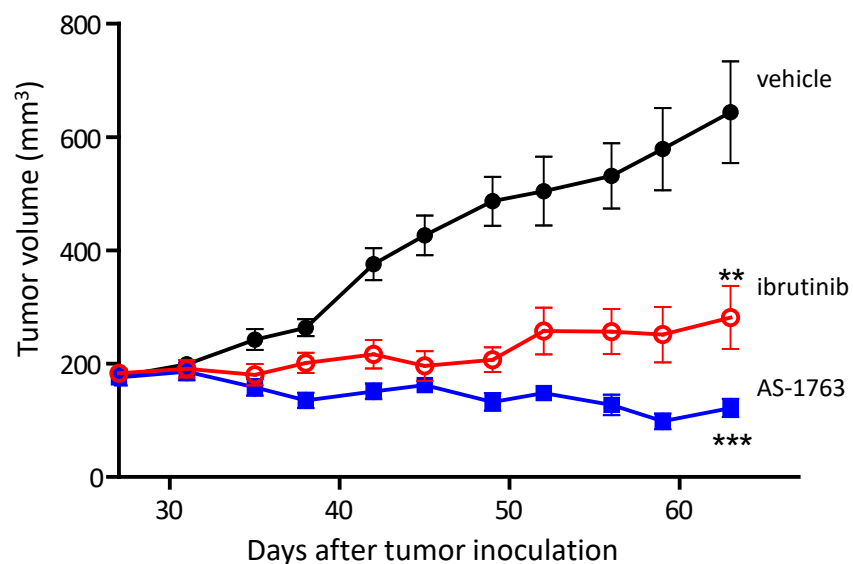
Only inhibited 6 other kinases in a total of 291 kinases tested at 0.3  $\mu$ M concentration.





# AS-1763: In Vivo Antitumor Effect against BTK<sup>C481S</sup> Mutant

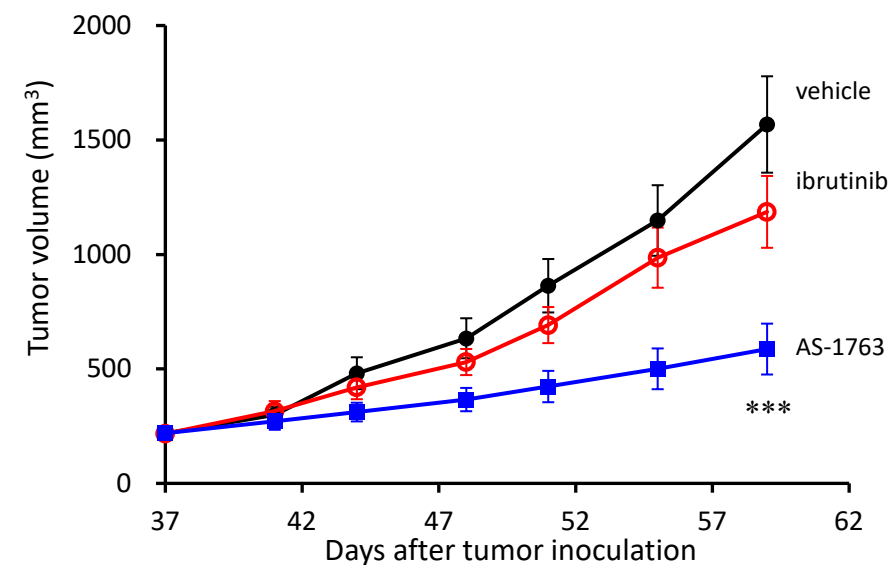
- ◆ In vivo antitumor effects of AS-1763 on human B-cell non-Hodgkin lymphoma cell line, OCI-LY10 tumor xenograft mouse model (n=8-10)



Ibrutinib: 25 mg/kg QD  
AS-1763: 60 mg/kg BID

\*\*: p < 0.01  
\*\*\*: p < 0.001

- ◆ In vivo antitumor effects of AS-1763 on ibrutinib-resistant BTK<sup>C481S</sup> knock-in OCI-LY10 tumor xenograft mouse model (n=11)



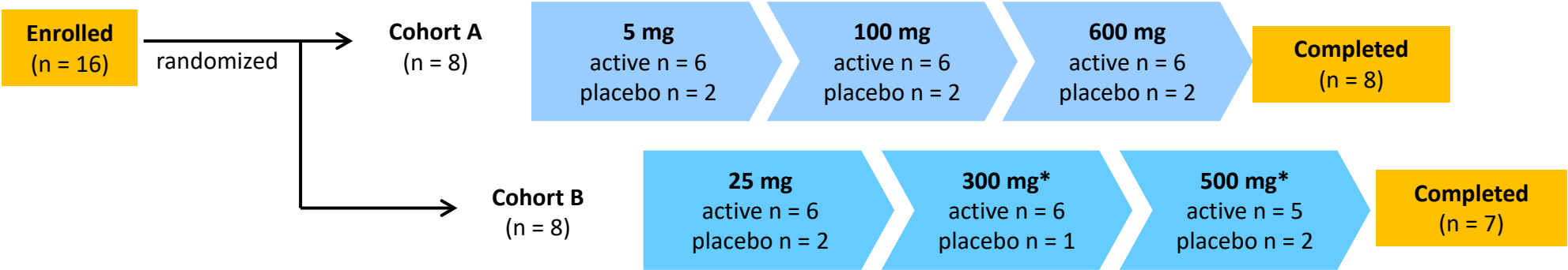
Ibrutinib: 25 mg/kg QD  
AS-1763: 60 mg/kg BID

\*\*\*: p < 0.001



## Study Design

Step 1 Single Ascending Dose (SAD) Part	Step 2 Relative Bioavailability (BA) Part
<ul style="list-style-type: none"><li>• Double-blind, placebo-controlled, randomized FIH study</li><li>• Simple formulation (solution)</li><li>• 6 dose levels (8 subjects/cohort A, 8 subjects/cohort B)</li><li>• 6 active / 2 placebo for each dose level</li><li>• Safety and tolerability</li><li>• Pharmacokinetics and pharmacodynamics (PD; CD69 upregulation on naïve B cells )</li></ul>	<ul style="list-style-type: none"><li>• Open label study</li><li>• Another cohort of 8 subjects</li><li>• The subjects were dosed with a single dose of AS-1763 100-mg tablet, and relative bioavailability with simple formulation was evaluated</li></ul>



\*One subject was withdrawn from the study on Day 1 of 300-mg period before the intake of treatment medication (placebo) by physician's decision. This subject showed AEs (Grade 2 lymphocytosis and Grade 2 neutropenia) which were considered treatment-emergent but not trial medication-related. No replacement was done at 300-mg and the following 500-mg periods.





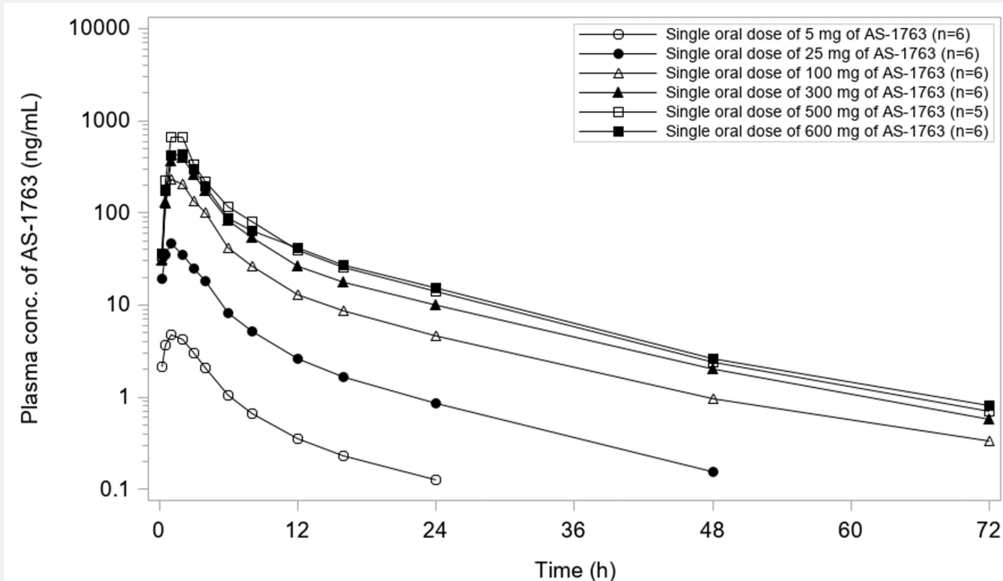
## Safety and tolerability

- AS-1763 was well-tolerated after single dose administration up to the maximum dose level (600 mg).
- No serious adverse events (AEs) were reported during the trial.
- Two Grade 2 AEs were reported in one subject, which were considered not related to trial medication.
- Other AEs reported were of mild intensity and showed no apparent dose-relationship in frequency.
- No clinically relevant changes from baseline were observed in all other safety parameters assessed (clinical laboratory, 12-lead safety ECGs, vital signs, or physical examinations).

## Pharmacokinetics

- After a single-dose oral administration, plasma concentration of AS-1763 rapidly reached the maximum and then declined in a biphasic manner across the dose range (median  $t_{max}$  between 0.5 and 1.5 hours; mean  $t_{1/2}$  between 8.4 and 12.1 hours).
- Mean AS-1763 exposures generally increased with dose up to 500 mg.

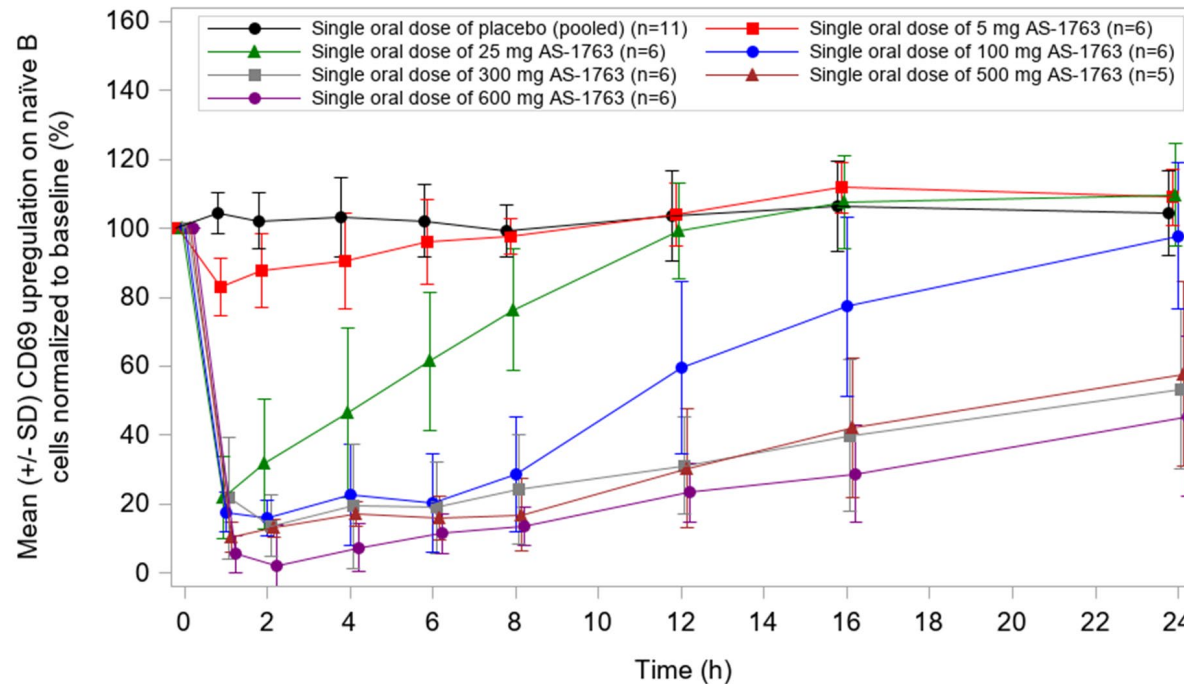
### < Plasma concentration of a single oral dose of AS-1763 >



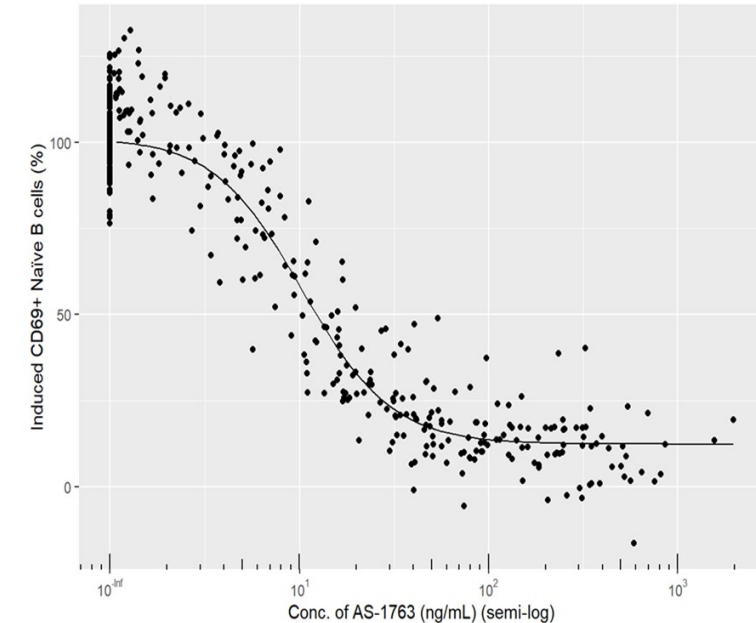


- Inhibition of B cell CD69 upregulation was observed for 5 mg onwards.
- Maximum inhibition (arbitrarily defined as  $\geq 80\%$ ) was observed at 1-2 hours post-dose from 100 mg to 600 mg, and the duration of inhibition was dose-dependent with values of 2, 6, 8 and 8 hours for 100, 300, 500, and 600 mg, respectively.
- Based on a PK/PD correlation analysis, the IC<sub>50</sub> value of AS-1763 on CD69 upregulation was calculated to be 10.5 ng/mL.

## < B cell CD69 upregulation after a single oral dose of AS-1763 >



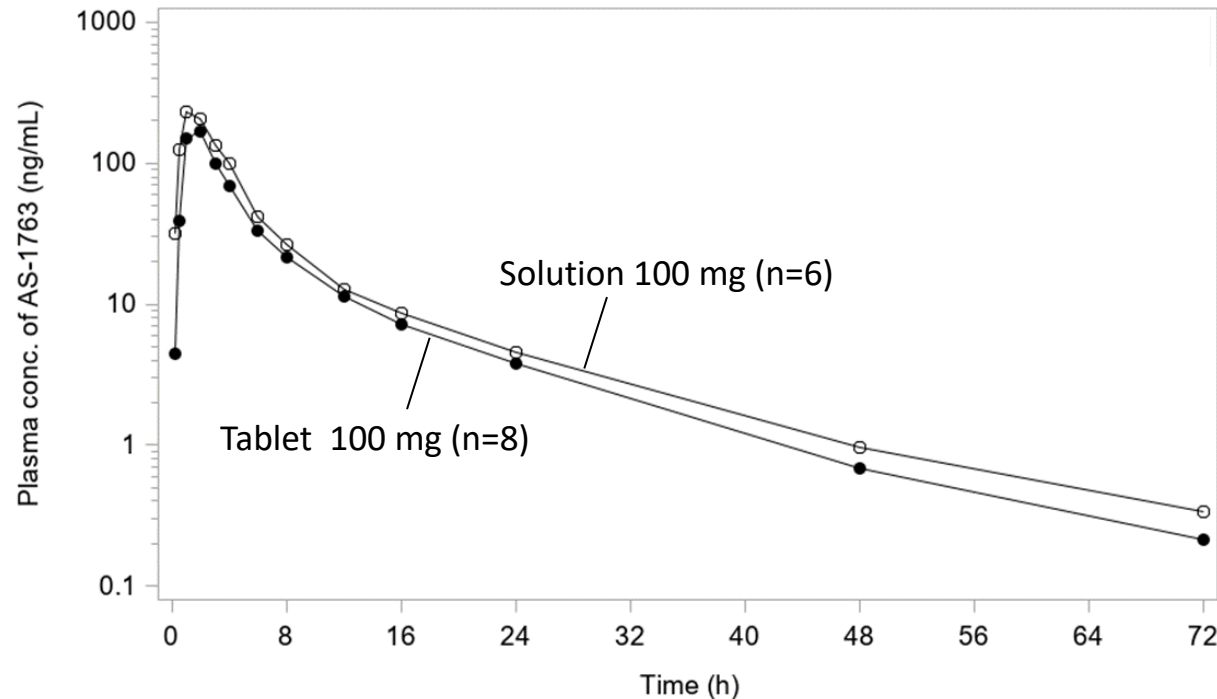
## < PK/PD correlation analysis >





- In the BA part, 100 mg tablet and the solution showed almost similar PK profile while the exposure of 100 mg tablet was slightly lower than the that of the solution.
- The PK/PD data and favorable safety profile in healthy volunteers support a planned Phase 1b clinical study with AS-1763 tablet twice daily dosing in relapsed/refractory CLL and B-cell NHL.

## <PK of Tablet vs Solution after a Single oral dose AS-1763>





# Sofnobrutinib (AS-0871): Excellent Kinase Selectivity

## ◆ Targeting Inactive Conformation of BTK



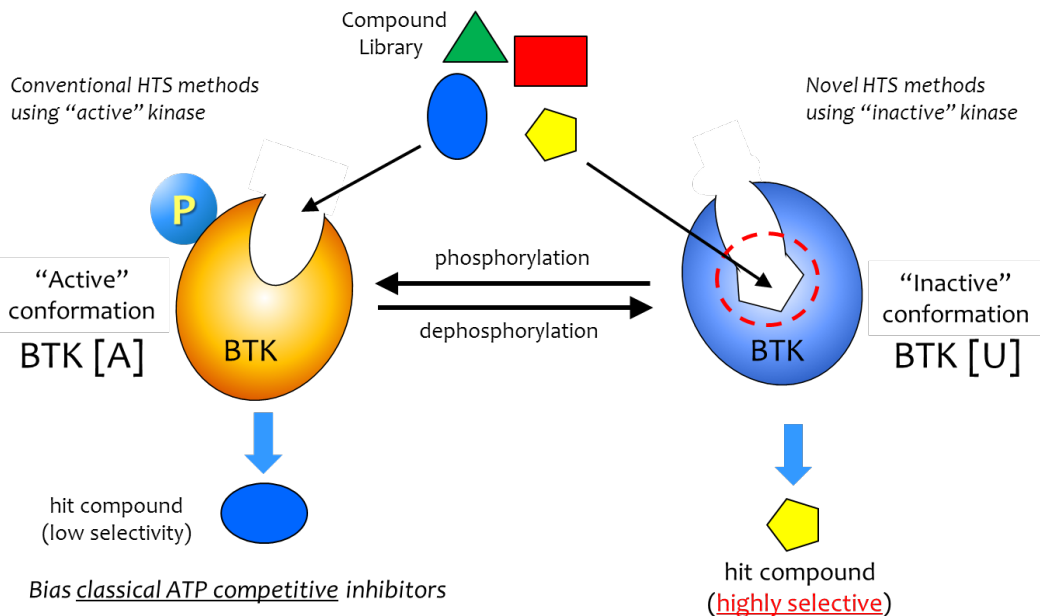
### TR-FRET binding assay targeting unactivated form of Bruton's tyrosine kinase

Tokiko Asami\*, Wataru Kawahata, Masaaki Sawa  
Carna Biosciences, Inc., BMA 3F, 1-5-5 Minatogino-Minamimachi, Chuo-ku, Kobe 650-0047, Japan



### Design and Synthesis of Novel Amino-triazine Analogues as Selective Bruton's Tyrosine Kinase Inhibitors for Treatment of Rheumatoid Arthritis

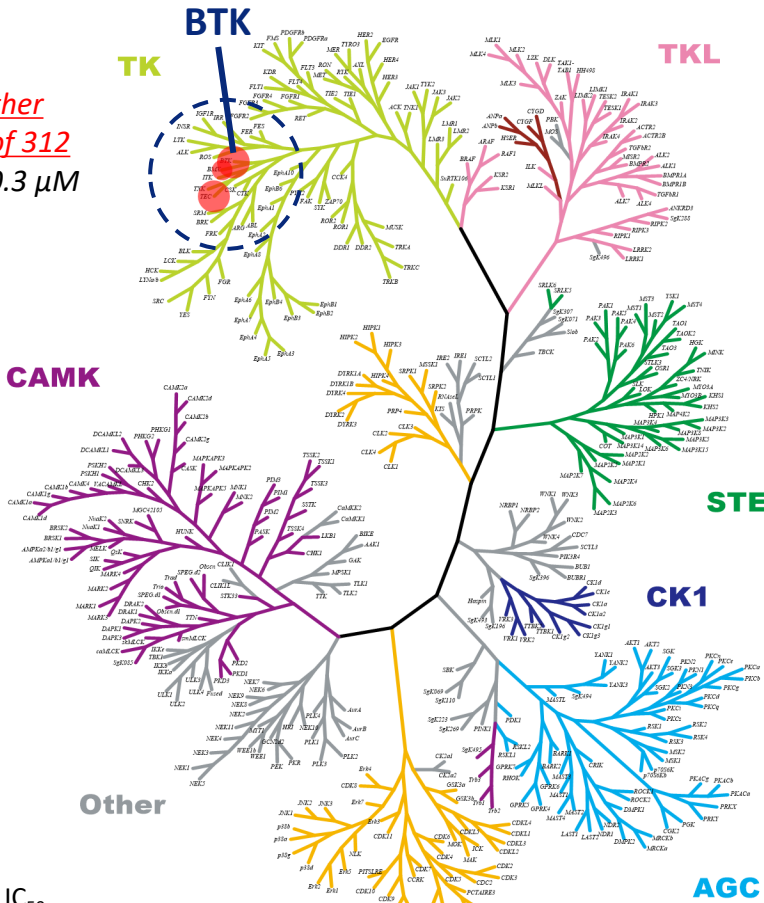
Wataru Kawahata\*, Tokiko Asami, Takao Kiyoi, Takayuki Irie, Haruka Taniguchi, Yuko Asamitsu, Tomoko Inoue, Takahiro Miyake, and Masaaki Sawa  
Research and Development, Carna Biosciences, Inc., 3rd Floor, BMA, 1-5-5 Minatogino-Minamimachi, Chuo-ku, Kobe 650-0047, Japan



	BTK IC <sub>50</sub> (nM)	
	BTK[A]	BTK[U]
sofnobrutinib	3.4	0.3

## ◆ Kinase Selectivity Profiling

Only inhibited 2 other kinases in a total of 312 kinases tested at 0.3  $\mu$ M concentration.



IC<sub>50</sub>  
< 10 nM  
10 – 100 nM  
100 – 1000 nM

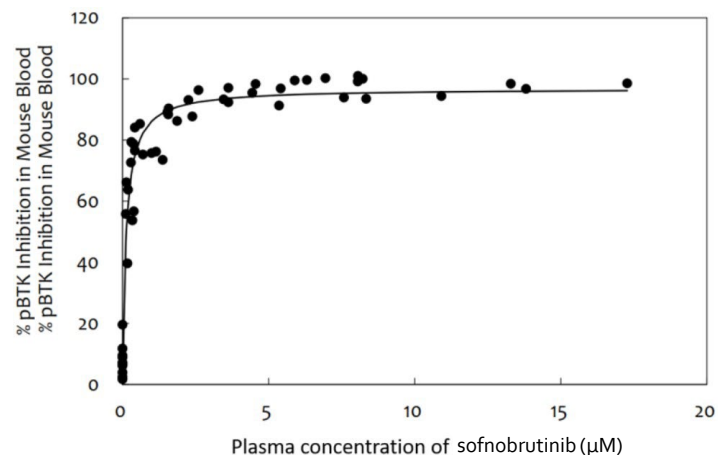
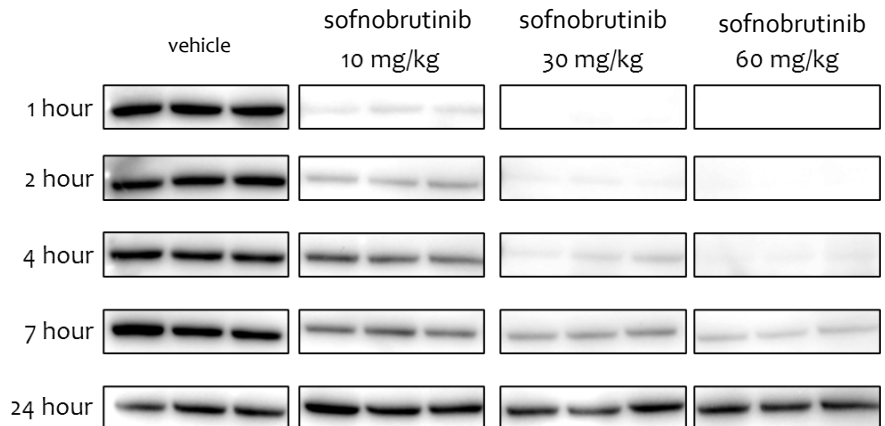
CMGC



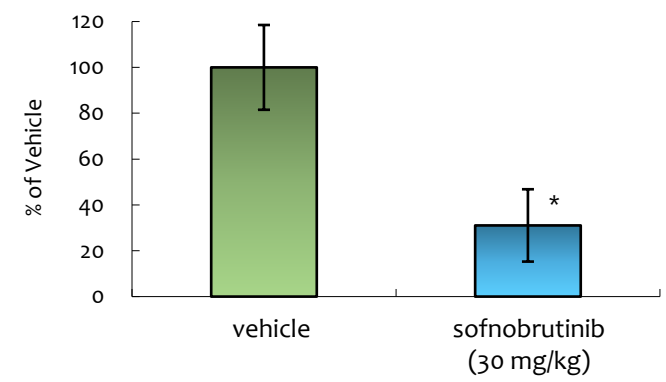
# Sofnobrutinib (AS-0871): In Vivo Therapeutic Efficacy

## ◆ PK/PD Study

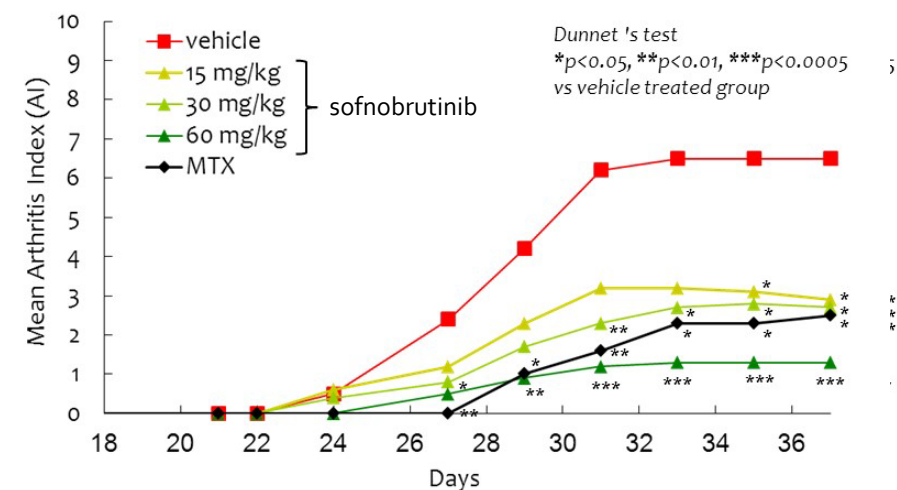
Auto-phosphorylation status of BTK was measured following oral single administration of sofno Brutinib .



## ◆ Passive cutaneous anaphylaxis (PCA) mouse model (n=5)



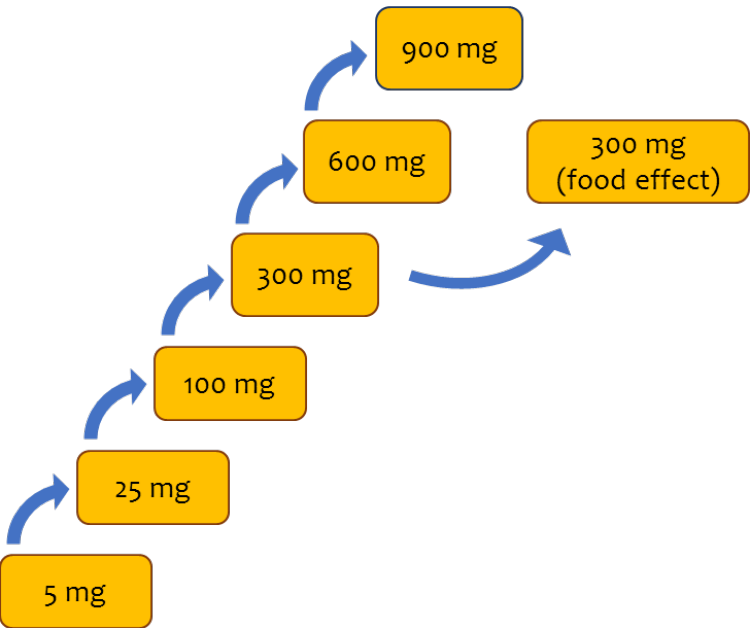
## ◆ Collagen-induced arthritis (CIA) mouse model (n=10)





## SAD Part (Completed)

Step 1 Single Ascending Dose (SAD)	Step 2
<ul style="list-style-type: none"><li>• 6 dose levels (8 subjects/cohort)</li><li>• Placebo controlled (6 active / 2 placebo)</li><li>• Safety and tolerability</li><li>• Pharmacokinetics and pharmacodynamics</li></ul>	<ul style="list-style-type: none"><li>• Food effect</li></ul>



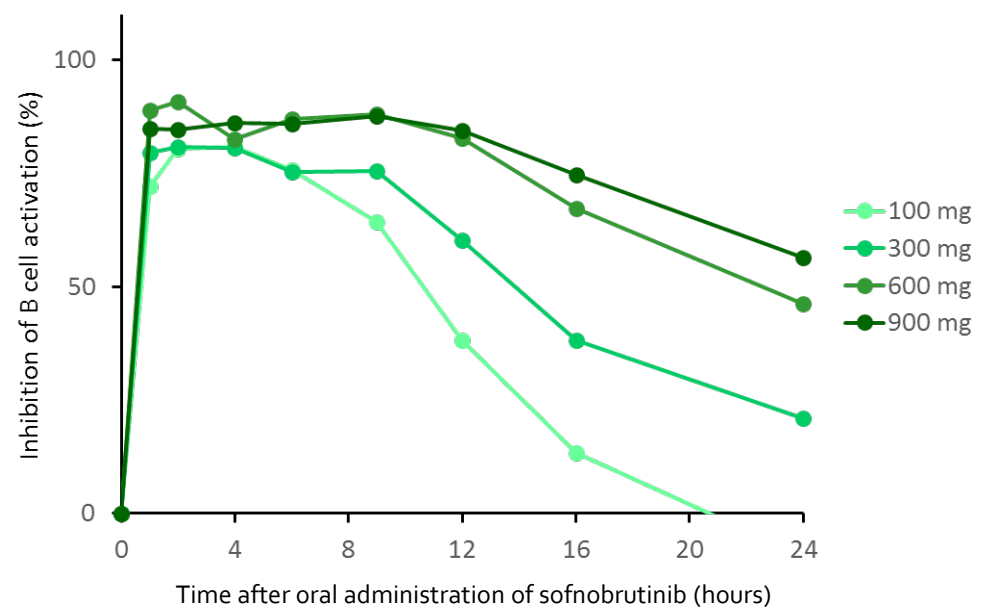
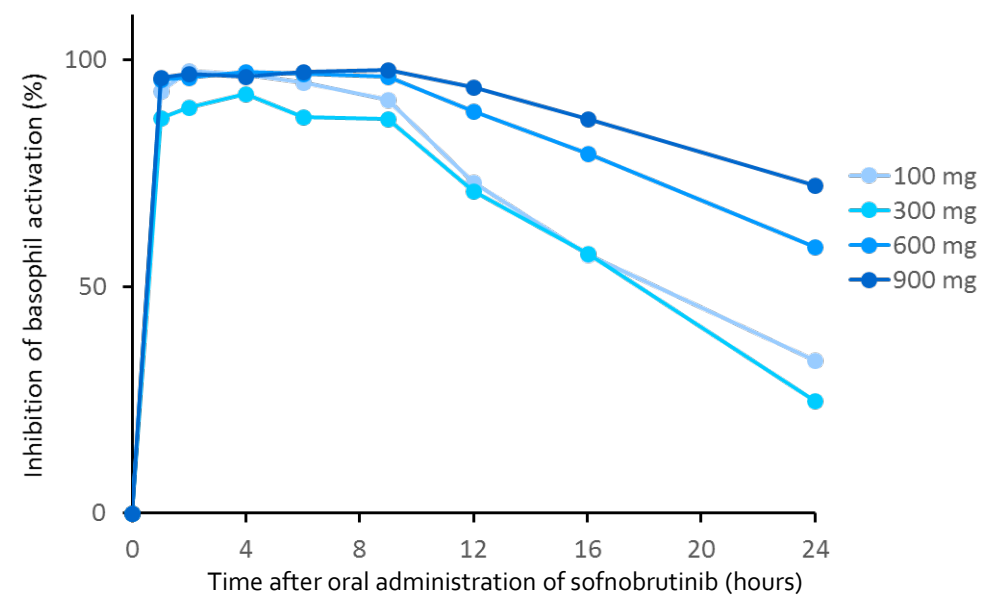
- Sofnobrutinib is well-tolerated without any safety concerns.
- Favorable pharmacokinetic profile.
- Blood samples to assess PD effects were analyzed for evaluation of the B-cell and basophil responses. Administration of sofno Brutinib at 100mg or above resulted in strong inhibition of B-cell and basophil activation.
- Switching to a new formulation in the MAD study.



# Pharmacodynamics of sofno Brutinib (AS-0871)



- Pharmacodynamic study demonstrated that subjects who received sofno Brutinib showed dose proportional inhibitions in basophil and B-cell activations, and significant and sustained inhibitory effects were observed at 100 mg and above.
- Oral administration of sofno Brutinib achieved therapeutic plasma levels needed to inhibit B cells and basophils activation, suggesting that sofno Brutinib has a potential to become a new treatment option for inflammatory diseases.



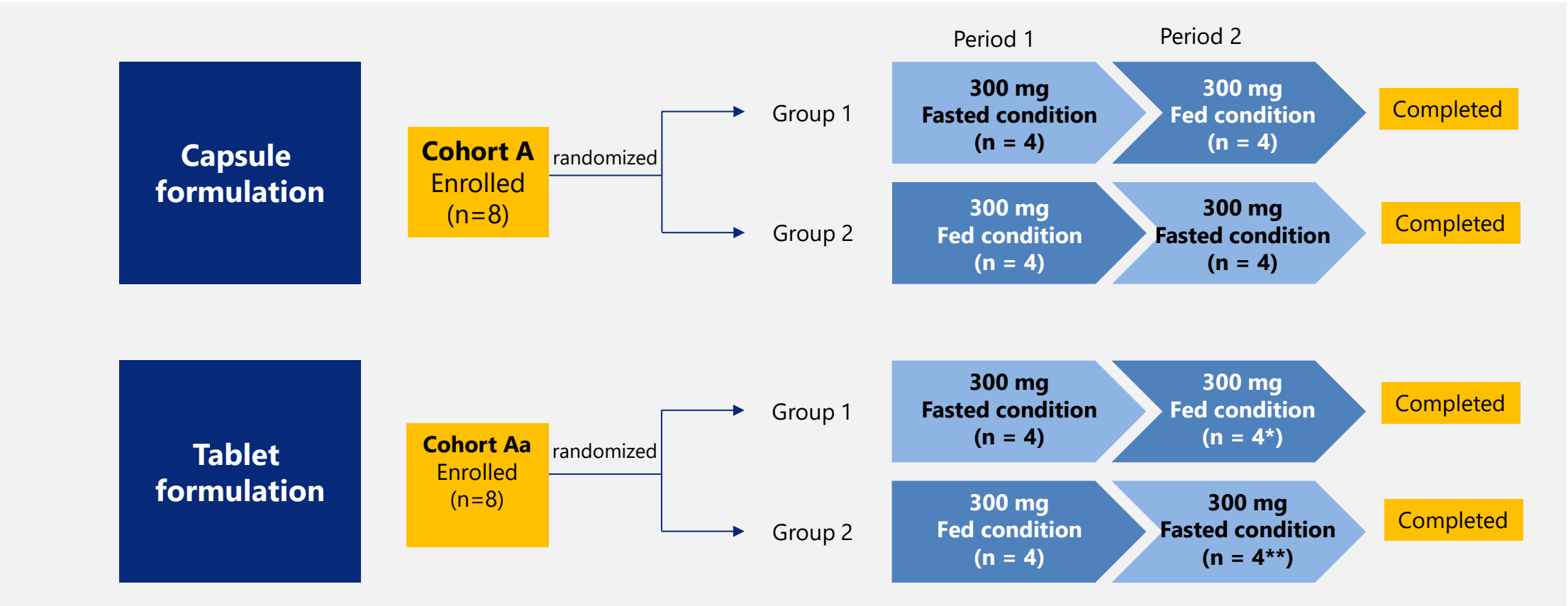




# Sofnobrutinib (AS-0871): Phase 1 MAD Study BA part

## Study Design of rBA/FE part

PK, safety, and tolerability after single-dose oral administration of sofnobrutinib, formulated as capsules or tablets, were evaluated under fasted and fed conditions in an open-label, randomized, 2-period crossover design. Eight healthy subjects (Cohort A or Cohort Aa) were randomized to either Group 1 or Group 2 (4 subjects per group).



\*One subject vomited after dosing (considered not related to study drug), excluded from the PK analysis.

\*\*One subject withdrew from the study due to personal reasons before dosing.

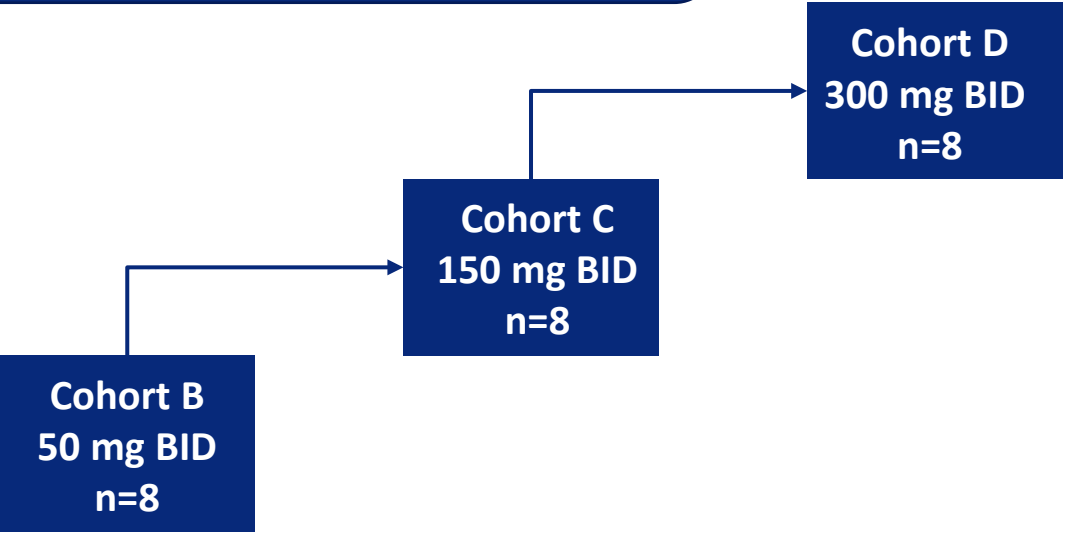




## Study Design of MAD part

- In the MAD part, safety, tolerability, PK, and PD of 3 multiple ascending doses of sofno Brutinib, following 14-day multiple dose oral administration of sofno Brutinib, will be investigated using a double blind, placebo-controlled, randomized design in 3 cohorts of 8 healthy subjects each.
- ◆ Dosing was completed as planned.
- ◆ The final Clinical Study Report (CSR) for the Phase 1 MAD study has been finalized in November 2023.

14-days dosing for each cohort



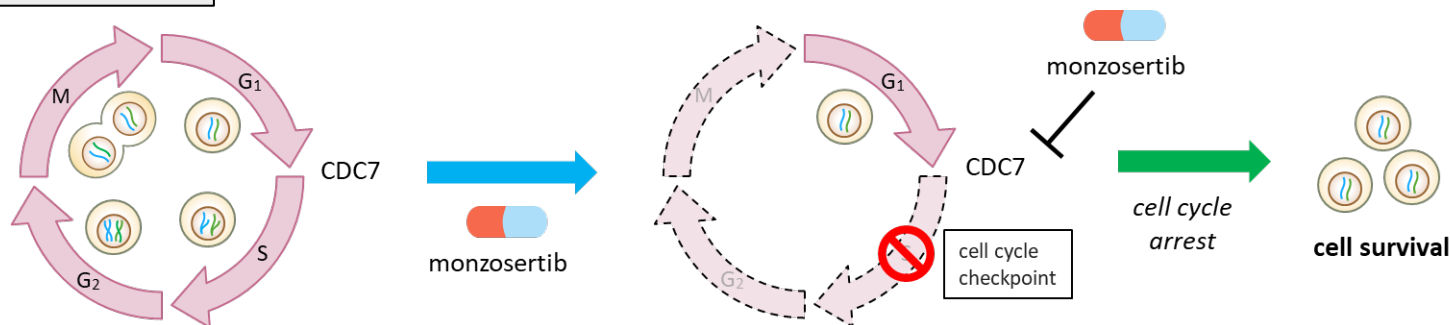
- Well tolerated with no dose-limiting AEs.
- Favorable safety profile up to 300 mg BID.
- The exposure levels increased dose-dependently, and favorable PK profile with tablet formulation was confirmed.
- Achieved almost complete inhibition of basophil activation (PD marker) at 150 mg and 300 mg BID at a steady state



## CDC7 Kinase Inhibitor: MoA of monzosertib

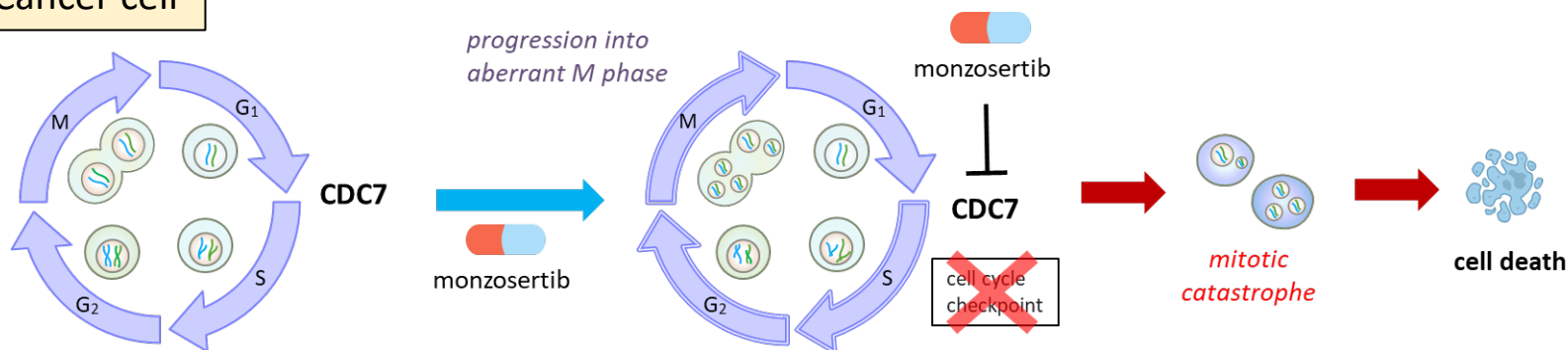
- ◆ CDC7 (cell division cycle 7) is a serine/threonine kinase that facilitates DNA replication during DNA synthesis
- ◆ Over expression of CDC7 has been reported to cause uncontrolled proliferation of many cancer types

### Normal cell



*Inhibition of CDC7 in normal cells does not cause cell death, only leads to cell cycle arrest at the DNA replication checkpoint*

### Cancer cell



*Inhibition of CDC7 in cancer cells causes lethal S phase or M phase progression when the cancer cells have a defective cell cycle checkpoint*

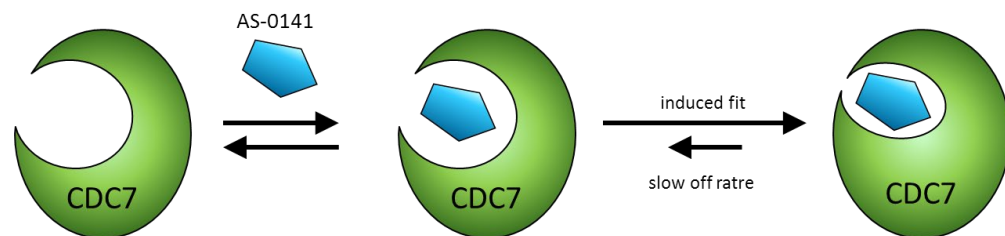


## ◆ Monzosertib has a unique inhibitory mechanism for CDC7 kinase (time-dependent inhibition)



Research paper  
Discovery of novel furanone derivatives as potent Cdc7 kinase inhibitors  
Takayuki Irie<sup>a,\*</sup>, Tokiko Asami<sup>a</sup>, Ayako Sawa<sup>a</sup>, Yuko Uno<sup>a</sup>, Mitsuharu Hanada<sup>a</sup>, Chika Taniyama<sup>b</sup>, Yoko Funakoshi<sup>b</sup>, Hisao Masai<sup>c</sup>, Masaaki Sawa<sup>a</sup>  
<sup>a</sup> Research and Development, Carina Biosciences, Inc., 3F BMA, 1-5-5 Minatogawa-Minamimachi, Chuo-ku, Kobe, 650-0047, Japan  
<sup>b</sup> Research and Development Department, SRI Biotech Co., Ltd., Ezumi Garden Tower 18F, 1-6-1 Rappongi, Minato-ku, Tokyo 106-6018, Japan  
<sup>c</sup> Department of Genome Medicine, Tokyo Metropolitan Institute of Medical Science, 2-1-6 Konishikazawa, Setagaya-ku, Tokyo 158-8506, Japan

pubs.acs.org/jmc Drug Annotation  
**Discovery of AS-0141, a Potent and Selective Inhibitor of CDC7 Kinase for the Treatment of Solid Cancers**  
Takayuki Irie<sup>a,\*</sup>, Tokiko Asami<sup>a</sup>, Ayako Sawa<sup>a</sup>, Yuko Uno<sup>a</sup>, Chika Taniyama<sup>b</sup>, Yoko Funakoshi<sup>b</sup>, Hisao Masai<sup>c</sup>, and Masaaki Sawa<sup>a</sup>  
Cite This: *J. Med. Chem.* 2021, 64, 14153–14164 Read Online



Inhibitory potency (IC<sub>50</sub>) for CDC7 in the presence of 1 mM ATP

Without  
Preincubation

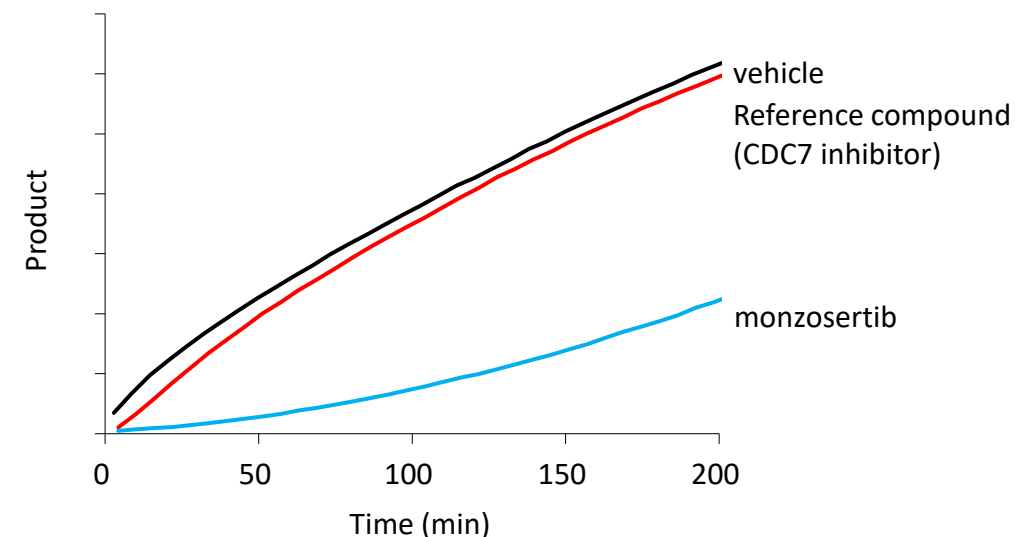
503 nM

With  
Preincubation

2.4 nM

## ◆ Monzosertib inhibits CDC7 in a reversible fashion but has a very slow off-rate

Rapid dilution assay for Cdc7 inhibitors. Recovery of enzymatic activity was monitored by formation of the phosphorylated product.



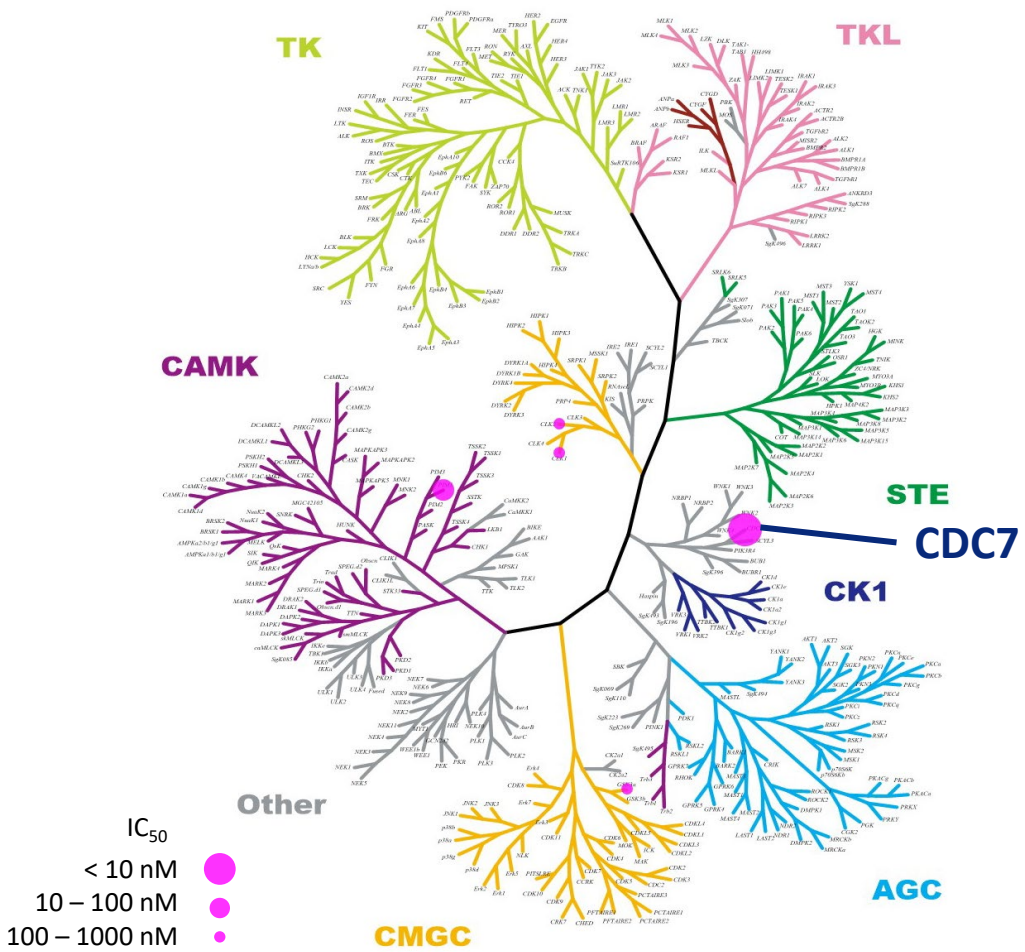


# Monzosertib (AS-0141): High Kinase Selectivity



## Kinase Selectivity Profiling

In the presence of 1 mM ATP with preincubation



## IC<sub>50</sub> values of hit kinases (at 1 mM ATP)

	IC <sub>50</sub> (nM)	
	Preincubation	
	-	+
CDC7	503	2.4
PIM1	30	34
CLK1	212	206
CLK2	270	227
GSK3α	189	251

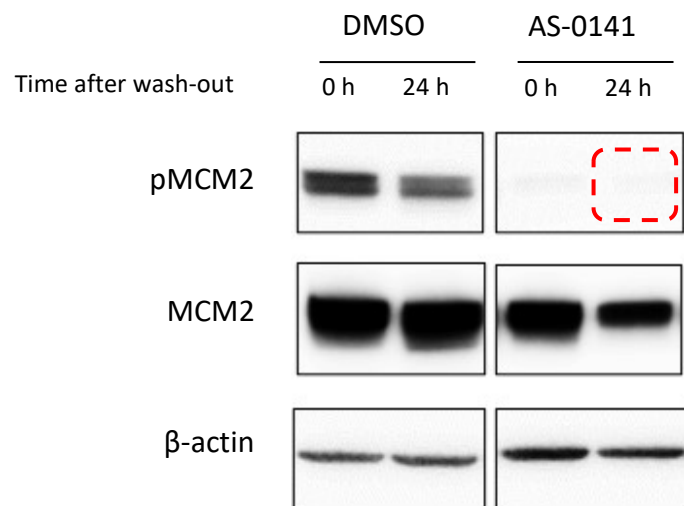
CDC7 is the only kinase that shows preincubation effect



# Monzosertib (AS-0141): Strong Cellular Activity

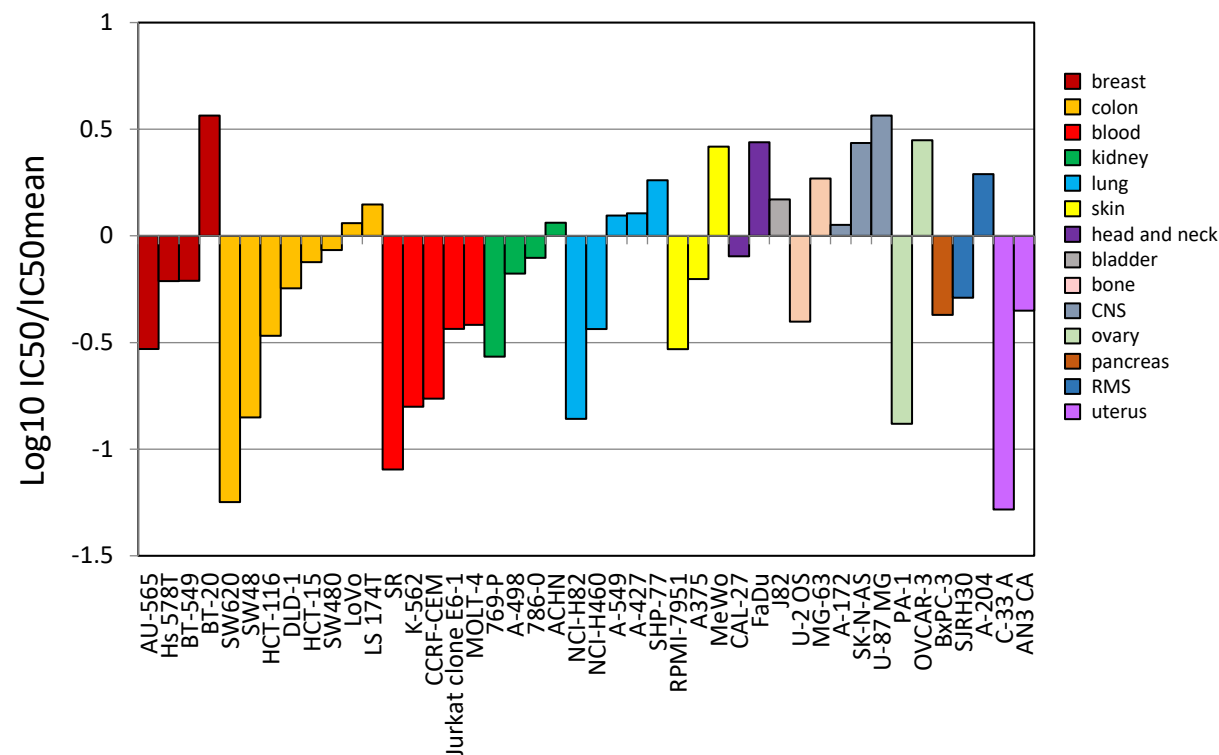
## ◆ Prolonged inhibition in cells

Human colon cancer cell line, Colo-205 cells were treated with DMSO control or monzosertib. After washout of the inhibitor, the cells were further incubated in the same media for 0 or 24 h and subjected to western blot analysis.



the inhibitory effect of monzosertib on the phosphorylation of MCM2 in cells continued up to 24 h after washing out

## ◆ Monzosertib potently inhibited growth in a wide range of tumor cell lines, including solid and hematological tumors

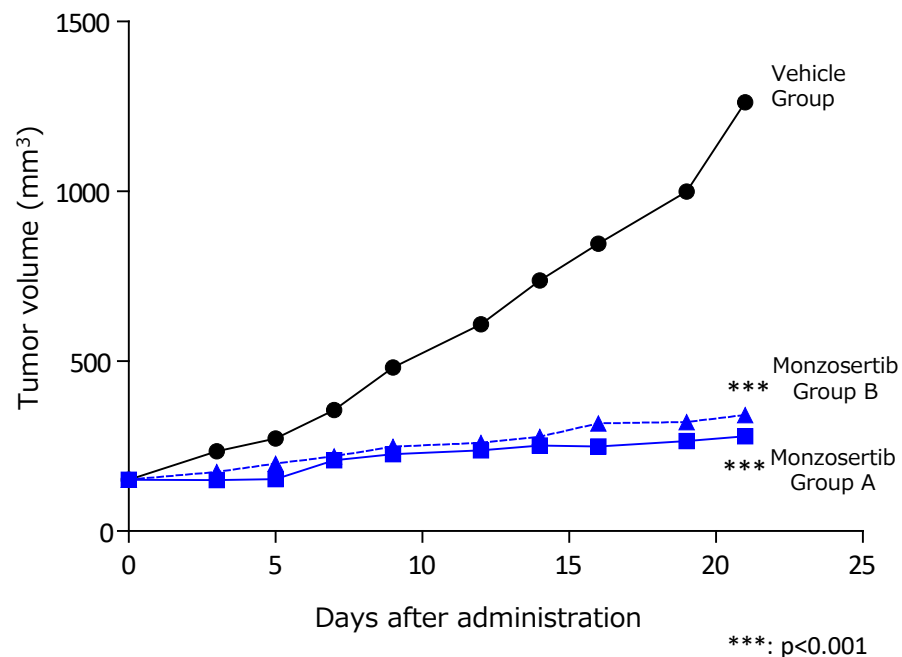


44 Cancer cell lines (Oncolines at NTRC)

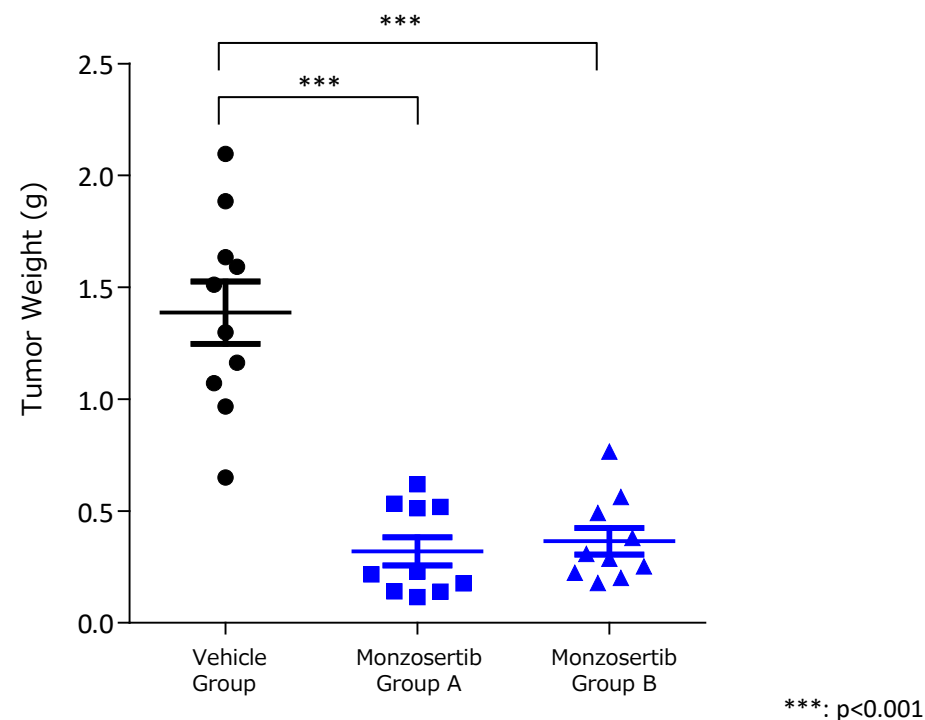


## In vivo antitumor efficacy of monzosertib in a SW620 (human colon cancer) xenograft mouse model

Tumor Growth Curve (Mean, n = 10)



Final Tumor Weight of Each Mouse



Monzosertib group A: 60 mg/kg TID, 4d ON/2d OFF  
Monzosertib group B: 120 mg/kg QD



“Carna” is a goddess of Roman mythology who takes care of human health, protecting the human heart and other organs as well as everyday life, and is said to be the root for the word “cardiac.”

The word “biosciences” is derived from the words 'biology' and 'life sciences.'

Carna Biosciences has created contemporary Carna goddess with protein kinase.

## **Carna Biosciences, Inc.**

### **Corporate Planning**

BMA3F 1-5-5 Minatojia-Minaimachi,

Chuo-ku, Kobe 650-0047

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

ir-team@carnabio.com

---

This document was prepared for the sole purpose of providing information to investors and is not intended as a solicitation for investment.

The forward-looking statements contained in this document are based on our plans and estimation and do not imply a commitment or guarantee of actual outcomes.

Investors should aware that the actual performance of the company could be materially different from our current forecasts.

The statements on the industry and other information were prepared based on the data assumed to be reliable. However, no guarantee is given regarding the accuracy or completeness of the information.

This document is presented on the assumption that all investors will make use of this document on their own judgment and responsibilities regardless of their purposes. Therefore, we do not assume no responsibility for any consequence caused by using this document.