

Financial Results

Q3 FY2023

(January to September 2023)

Carna Biosciences, Inc.



AGENDA

- 1** FY2023 Q3 Results
- 2** Updates on Pipelines in Clinical Development
- 3** Updates on Licensed Pipelines
- 4** Appendix



Q3 FY2023 Key Highlights

BTK inhibitor: AS-1763

- The first patient was dosed in Phase 1b study in the U.S in August 2023.
Multi-center clinical study
Principal Investigator : Dr. Nitin Jain, Department of Leukemia,
University of Texas MD Anderson Cancer Center.

BTK inhibitor: sofnobrutinib (AS-0871)

- The final Clinical Study Report (CSR) for the Phase 1 MAD study has been finalized in November 2023.
- Demonstrated a favorable safety and tolerability profile as well as a promising pharmacokinetic (PK) and pharmacodynamic (PD) profile in the MAD study.

International Nonproprietary Name (INN) : sofnobrutinib, Code name : AS-0871
MAD : Multiple Ascending Dose

DGK α inhibitor: GS-9911 (Out-licensed to Gilead Sciences, Inc.)

- The Phase 1 study information has been published on ClinicalTrials.gov in October 2023.
- A patent application filed jointly with Gilead: Received a notice of allowance in the U.S. in July 2023.



Q3 FY2023 Key Highlights

STING inhibitor

- Fresh Tracks Therapeutics (FRTX), the licensee of STING inhibitor, announced that its Board of Directors has approved a plan of liquidation and dissolution of the company in September 2023.

ddSP

- Sales were JPY711 mn, 78% progress to FY plan of JPY902 mn.
- Sales of kinase protein were JPY448 mn, reached a record-high for Q3 YTD.

Others

- All the unexercised series 19th subscription rights to shares expired at the end of the exercise period in August 2023.

ddSP : Drug Discovery Support business



FY2023 Q3 YTD Results by Business Segment

(JPY million)	FY2022 Q3YTD Actual	FY2023 Q3YTD Actual	YoY Change	FY2023 Plan as of Feb 10	vs. FY Plan	
Total Sales	1,095	711	-384 -35.1%	902	78.8%	
ddSP business	809	711	-98 -12.1%	902	78.8%	<ul style="list-style-type: none"> Sales were on track to achieve FY plan. While sales of kinase proteins were solid, overall sales declined YoY as revenue related to the license agreement with Gilead contributed to FY2022 sales.
ddRD business	286	—	-286	—	—	<ul style="list-style-type: none"> Received an upfront payment from FRTX and a milestone payment from BioNova in FY2022.
Total Operating Profit/Loss	-753	-1,201	-448	-1,890	—	
ddSP business	335	218	-117 -34.9%	221	98.8%	<ul style="list-style-type: none"> Strong sales of highly profitable protein accelerated the operating profit growth toward FY plan achievement. Operating profit declined YoY due to a decrease in sales.
ddRD business	-1,089	-1,420	-331	-2,111	—	<ul style="list-style-type: none"> Received an upfront payment and a milestone payment in 2022. Operating loss increased YoY due to the increase in R&D expenses.
Ordinary Profit/Loss	-735	-1,203	-467	-1,911	—	
Net Profit/Loss	-795	-1,230	-435	-1,936	—	
R&D Cost	1,267	1,323	+55	1,968		<ul style="list-style-type: none"> Continued investment in the programs in clinical trials.

Business plan for FY2023 in ddRD does not include potential milestone payments or upfront payments as the timing or the amounts are difficult to predict.

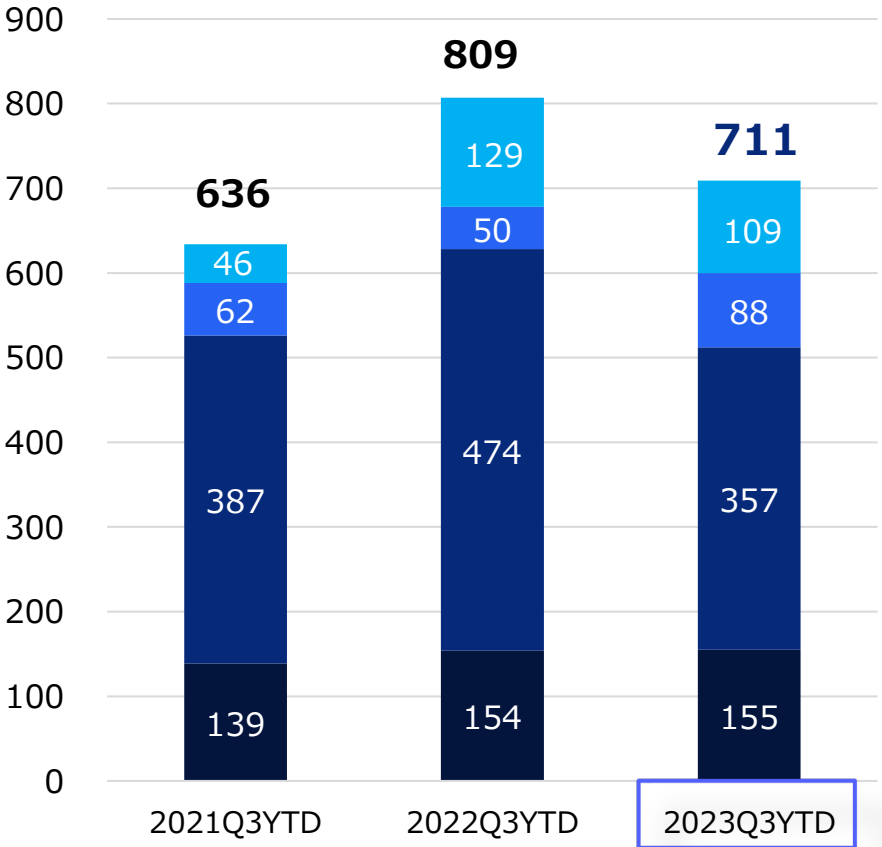
FRTX : Fresh Tracks Therapeutics, Inc.

Note : Rounded down to the nearest million yen



Drug Discovery Support Business Sales Trend by Region
(Consolidated)

(JPY million) Other Europe North America Japan



Japan

Increased 0.1% YoY

- While sales in 1H decreased due to a lack of large orders, the sales of profiling service were solid in Q3 and contributed to a marginal increase in sales YoY.

North America

Decreased 24.6% YoY

- Sales of kinase proteins increased significantly.
- Demand from biotech companies including AI-driven drug discovery companies remained strong.
- Overall sales declined YoY as sales from Gilead contributed to FY2022 sales.

Europe

Increased 74.4% YoY

- Thanks to the continuous order from AI-driven drug discovery companies, sales of kinase proteins were very strong.

Other

Decreased 15.2% YoY

- In China, the overall demand is still robust but sales declined YoY compared to the same period last year when sales were very strong.



Consolidated Balance Sheet



(JPY million)	As of Dec.31, 2022	As of Sep.30, 2023	Change	Reason for changes
Current assets	4,104	4,050	-53	Cash and deposits +11
Cash and deposits	3,379	3,390	+11	Amount raised from exercised share options +1,349
Non-current assets	162	141	-20	
Total assets	4,266	4,192	-74	
Current liabilities	436	277	-159	Accounts payable -127 Current portion of long-term debt -32
Non-current liabilities	188	98	-89	Long term loans payable-89
Total liabilities	624	375	-248	
Total net assets	3,641	3,816	+174	Capital stock and capital surplus +1,388 Retained earnings -1,230
Total liabilities and net assets	4,266	4,192	-74	

Shareholders' equity ratio	85.0%	91.0%
BPS	255.0yen	222.61yen
PBR	2.0x	3.3 x
Share price of Carna	520yen	727yen

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



<Exercise of Subscription Rights to Shares for Q3 YTD FY2023>

	19 th Subscription Rights to Shares	20 th Subscription Rights to Shares	Total
Amount raised (JPY)	47mil.	1,302mil.	1,349mil.
No. of shares exercised (Shares)	50,000	2,836,500	2,886,500

<20th Subscription Rights to Shares >

	DEC. 2022	JAN.-APR.2023	Total
Amount raised (JPY)	300 mil.	1,302mil.	1,602mil.
No. of shares exercised (Shares)	550,000	2,836,500	3,386,500
No. of Exercised rights / No. of total rights issued	16.2 %	83.8%	100%

Cash and deposits as of the end of September 2023 : 3,390million JPY.

Financing

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



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">Phase 1 SAD and BA part in healthy volunteers were completed in the Netherlands.In August 2023, the first patient was dosed in Phase 1b study in the U.S. <div>Multi-center clinical study Principal Investigator : Dr. Nitin Jain, Department of Leukemia, University of Texas MD Anderson Cancer Center.</div>
sofno Brutinib (AS-0871)	BTK	Immune-inflammatory diseases	<ul style="list-style-type: none">Phase 1 clinical trials (SAD study and MAD study) in healthy volunteers has been completed.The Clinical Study Report (CSR) for the Phase 1 MAD study has been finalized in November 2023 .Demonstrated a favorable safety and tolerability profile as well as a promising PK/PD profile in the MAD study.
monzosertib (AS-0141)	CDC7/ASK	Cancer	<ul style="list-style-type: none">Phase 1 study in cancer patients is in progress in Japan.Dose escalation part is on going. <div>Clinical trial site : National Cancer Center Hospital and National Cancer Center Hospital East</div>

SAD : Single Ascending Dose
MAD : Multiple Ascending Dose
BA : Bioavailability



Mechanism/ Indication

AS-1763 is an **orally available** Bruton's Tyrosine Kinase (BTK) inhibitor 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 ibrutinib**, the first BTK inhibitor in the market.

To minimize a risk of side effects

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



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 assays 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
Principal Investigator : Dr. Nitin Jain, Department of Leukemia, University of Texas MD Anderson Cancer Center.

Clinical trails in progress

Phase 1b study in the U.S.: Open and enrolling.

Indication

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

Status

- Opened 6 clinical sites.
- First patient dosed in August.
- Observed safety and tolerability at the first dose level.
- Proceed to the next higher dose level in October.



AS-1763: Potential Market Size (B-cell Malignancies)



Present

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

Est. market size

Estimated sales of covalent BTK inhibitors (ibrutinib, acalabrutinib, zanubrutinib)

Est. sales for 2021 \$9 billion

Est. sales for 2029 \$17 billion ※Expected CAGR is 8%

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.



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.



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.



Objectives of Single Ascending Dose (SAD) study

Ref. P.38, P.41-P.43

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 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

Novartis announced a positive Ph 2b result that remibrutinib rapidly and effectively improved the symptoms of patients whose CSU was inadequately controlled. *

*J Allergy Clin Immunol.2022;150:1498-506.

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.



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.



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

- Open-label Phase I study in patients with unresectable advanced, recurrent, or metastatic solid tumors in Japan
- Dose Escalation on-going



Updates on Licensed Pipelines

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



Out-licensed Programs

Program/ Partner	Compound (Target)	Upfront payment	Total milestone payments expected	Royalty	Region	Contract date	Milestones received
DGKα inhibitor Gilead Sciences (Out-license)	GS-9911 (Immuno- oncology)	\$20M	\$450M	Undisclosed	Worldwide	Jun. 2019	\$10M (Dec. 2021)
Joint Research with Sumitomo Pharma	Kinase inhibitor (Psychiatric and neurological disorders)	JPY80M (including research milestone)	JPY10.6B	Undisclosed	Worldwide	Mar. 2018	

※Deleted the description on STING inhibitor from the list above due to the announcement of board approval of a plan of liquidation and dissolution by Fresh Tracks Therapeutics in September 2023.



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 α inhibitor: GS-9911

2. Indication: Cancer (immunotherapy)

- GS-9911 was introduced by Gilead in its presentation “Oncology Deep Dive” held in April 2022. The presentation suggested that GS-9911 potentiated anti-tumor immunity by activating T cell function and demonstrated anti-tumor immune responses in mouse models in combination with anti-PD-1 antibody.

3. The Phase 1 study information has been published on [ClinicalTrials.gov](https://clinicaltrials.gov/ct2/show/study/NCT06082960) in October 2023.

Official Title	A Phase 1 Study to Evaluate the Safety and Tolerability of GS-9911 as Monotherapy and in Combination With an Anti-PD-1 Monoclonal Antibody in Adults With Advanced Solid Tumors
ClinicalTrials.gov ID	NCT06082960



Partner



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

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

- 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.**

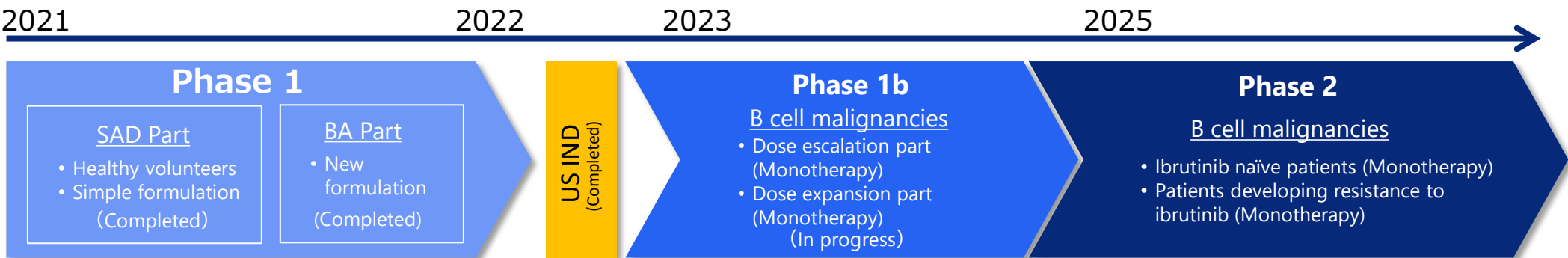


Appendix



AS-1763 : Targeting Blood Cancer

- Small molecule BTK inhibitor
- Non-covalent/reversible
- High kinase selectivity
- Inhibits both BTK wild type and ibrutinib resistant BTK C481S mutants
- Orally available
- Displayed strong anti-tumor effects in lymphoma model with both wild type and C481S mutant BTK
- Displayed efficacy in immuno-oncology model
- The first patient was dosed in August 2023 in Ph 1b study in the U.S.



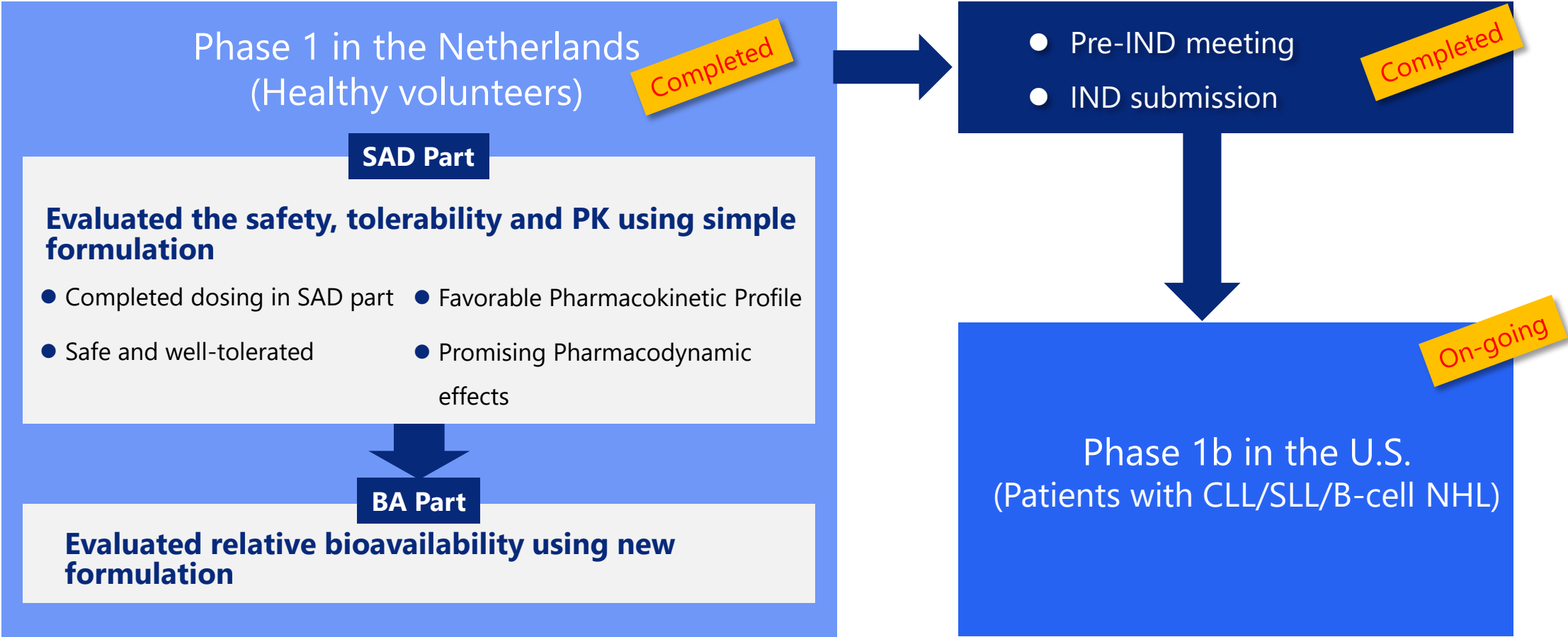
IND application: Investigational New Drug application

FPI: First Patient In

SAD: Single Ascending Dose

BA: Bioavailability

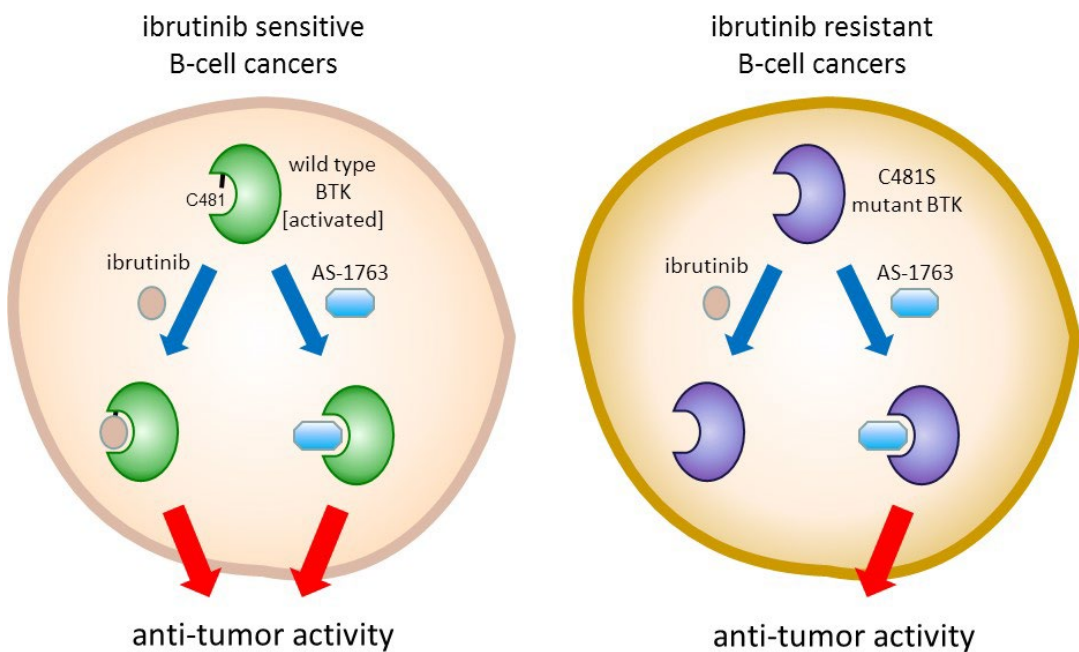
B-cell malignancies: Chronic Lymphocytic Leukemia (CLL), Small Lymphocytic Lymphoma (SLL), and B-cell non-Hodgkin Lymphoma (B-cell NHL), etc.



◆ The first patient was dosed in Phase 1b study in August 2023.



AS-1763: Potent Inhibitor of C481S mutant BTK



Journal of
**Medicinal
Chemistry**

pubs.acs.org/jmc

Drug Annotation

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

◆ IC₅₀ values of AS-1763 against wild-type and C481S-mutant BTK

	IC ₅₀ (nM)	
	BTK[A]	BTK ^{C481S}
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 ₅₀ (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

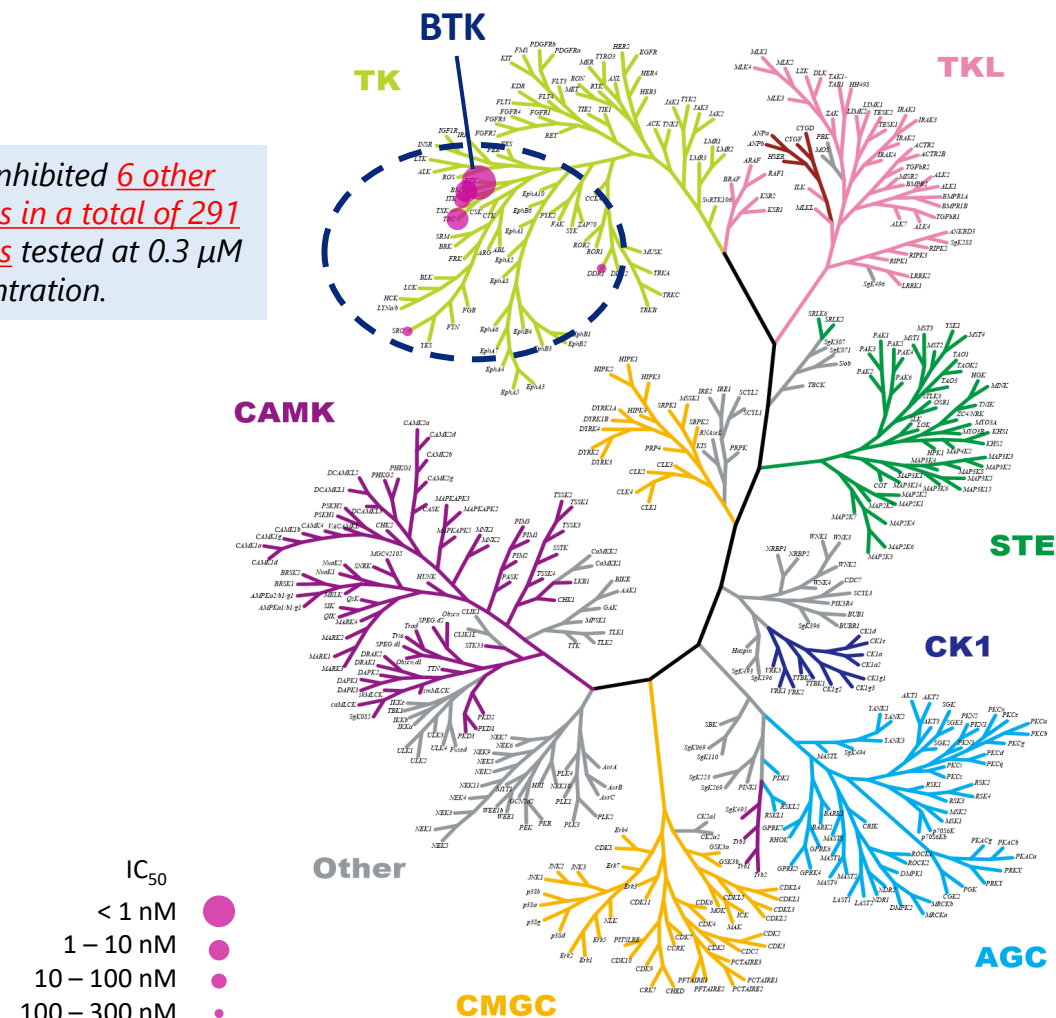
50-fold Stronger activity

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

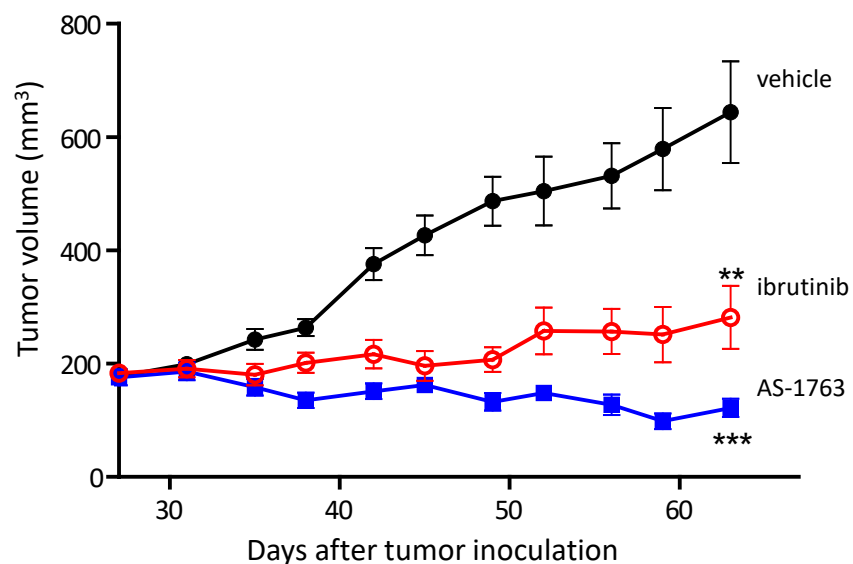
Only inhibited 6 other kinases in a total of 291 kinases tested at 0.3 μ M concentration.





AS-1763: In Vivo Antitumor Effect against BTK^{C481S} Mutant

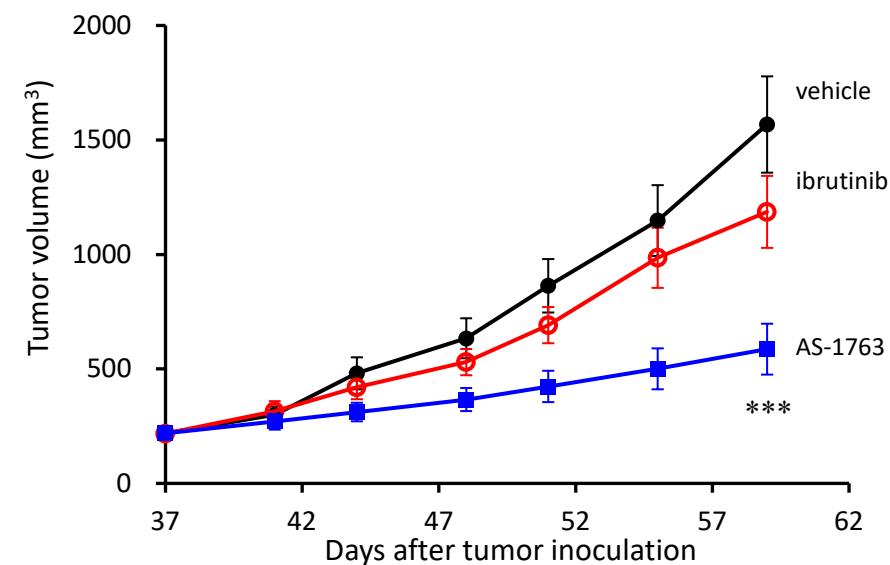
- ◆ In vivo antitumor effects of AS-1763 on human B-cell non-Hodgkin lymphoma cell line, OCI-LY10 tumor xenograft mouse model (n=8-10)



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^{C481S} knock-in OCI-LY10 tumor xenograft mouse model (n=11)



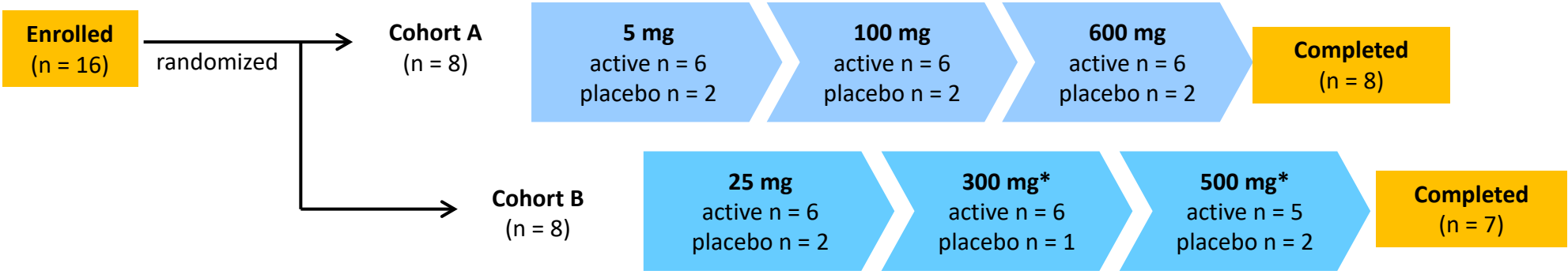
Ibrutinib: 25 mg/kg QD
AS-1763: 60 mg/kg BID

***: p<0.001



Study Design

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



*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.



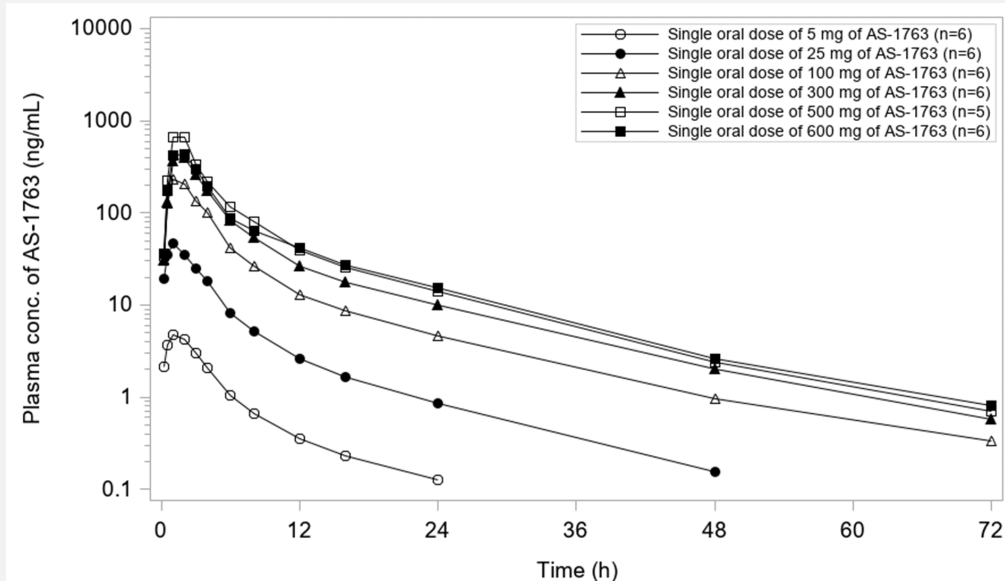
Safety and tolerability

- 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 >

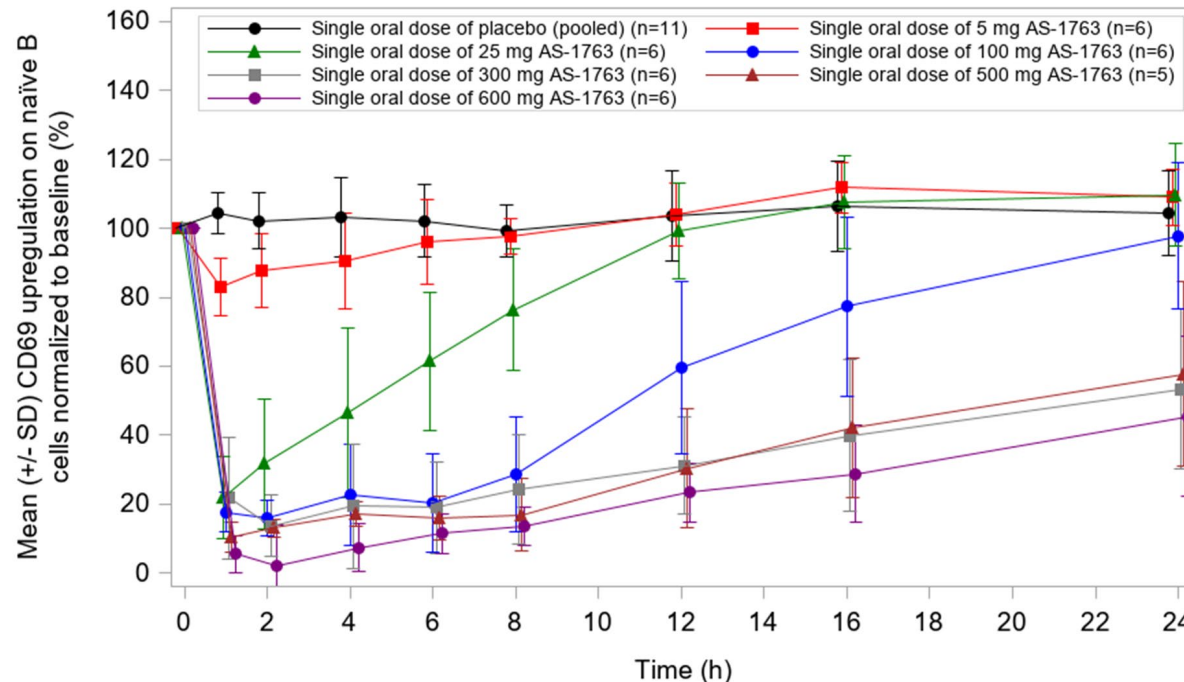




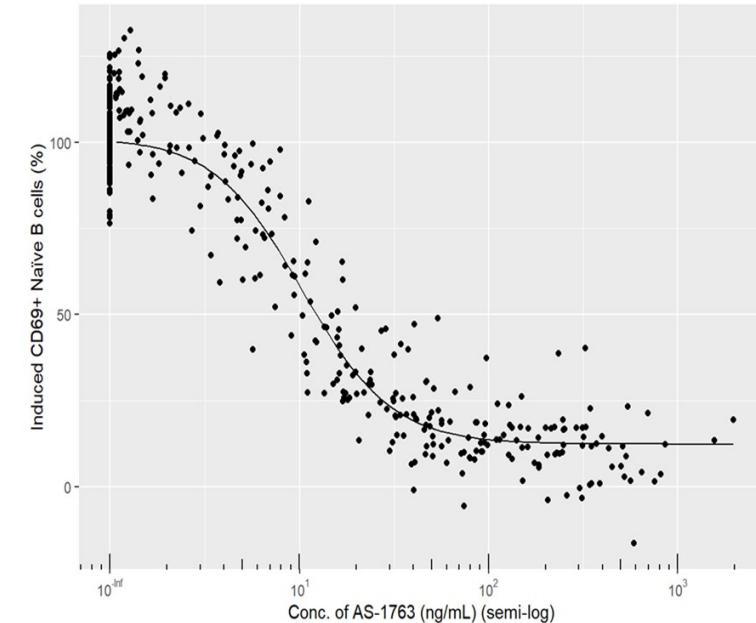
Pharmacodynamics of AS-1763

- 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 IC₅₀ 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 >

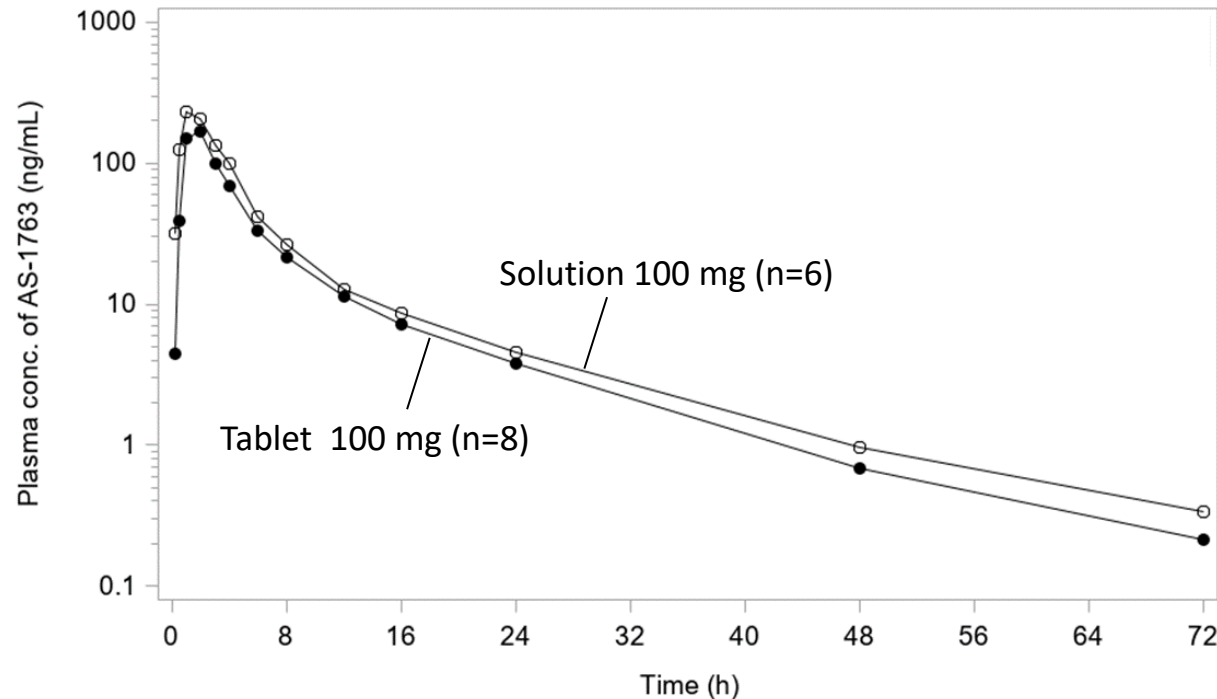




AS-1763: BA Part

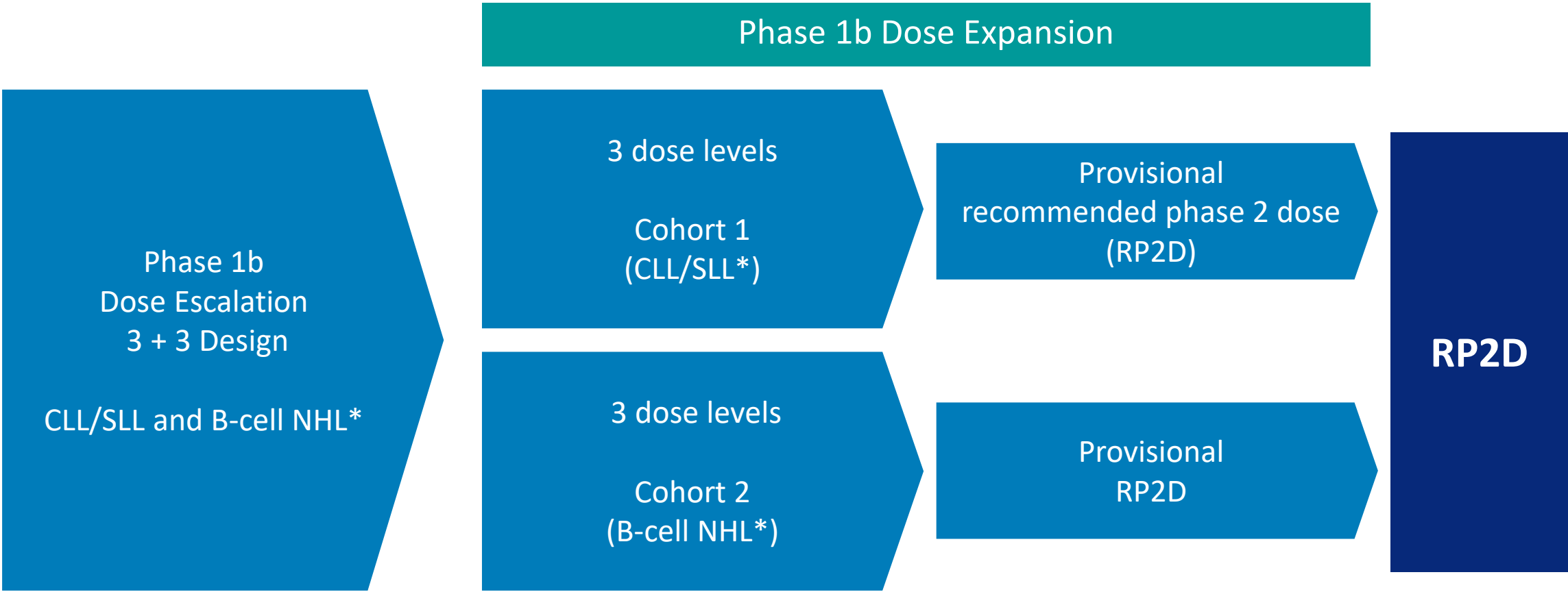
- In the BA part, 100 mg tablet and the solution showed almost similar PK profile while the exposure of 100 mg tablet was slightly lower than the that of the solution.
- The PK/PD data and favorable safety profile in healthy volunteers support a planned Phase 1b clinical study with AS-1763 tablet twice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

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





AS-1763: Phase 1b Schema (US)



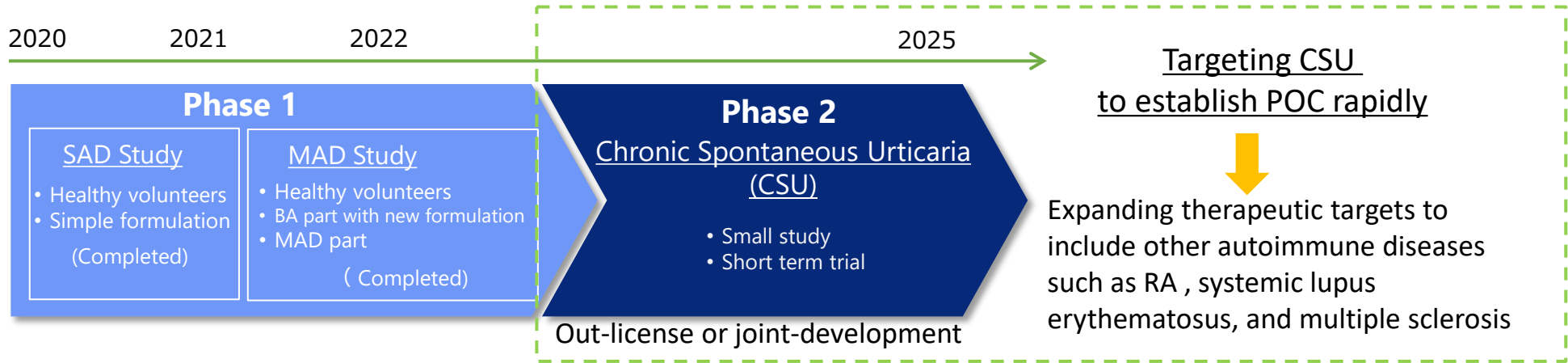
*Previously treated patients

◆ The first patient was dosed in August 2023.



AS-0871 : Targeting Immune-inflammatory diseases

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



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



Sofnobrutinib (AS-0871): Phase 1 Clinical Trial in Progress

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

MAD Part

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

Completed

Completed



Sofnobrutinib (AS-0871): Excellent Kinase Selectivity

◆ Targeting Inactive Conformation of BTK



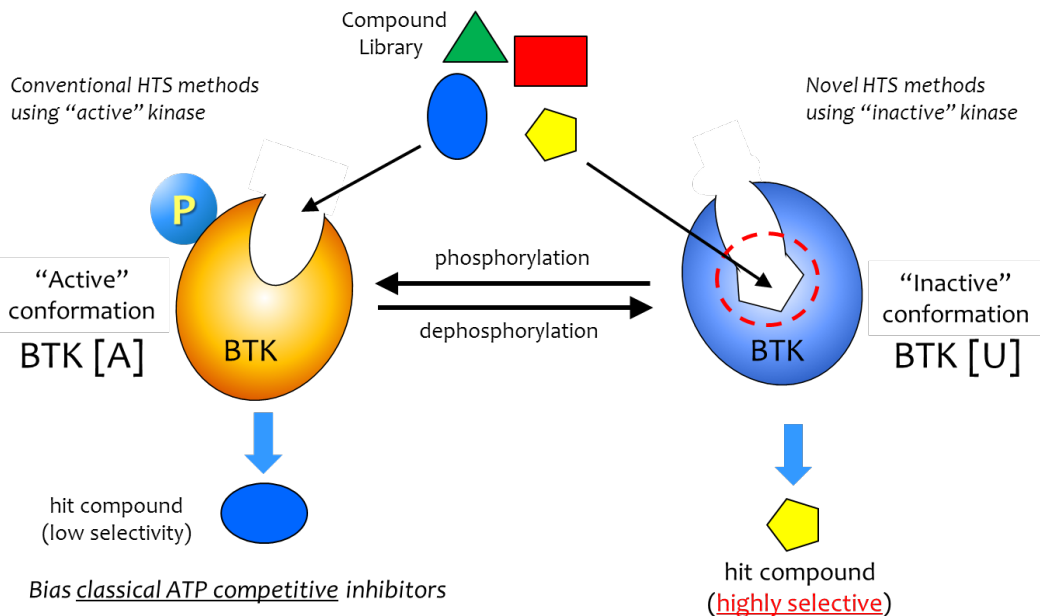
TR-FRET binding assay targeting unactivated form of Bruton's tyrosine kinase

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

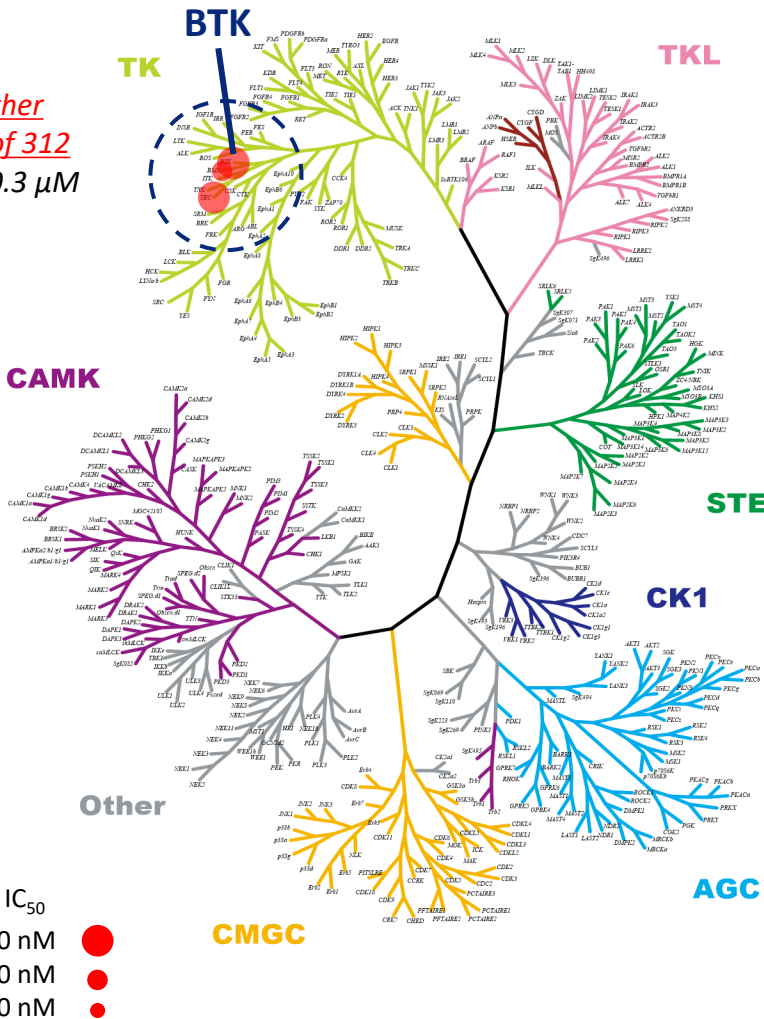
Wataru Kawahata*, Tokiko Asami, Takao Kiyoi, Takayuki Irie, Haruka Taniguchi, Yuko Asamitsu, Tomoko Inoue, Takahiro Miyake, and Masaaki Sawa
Research and Development, Carna Biosciences, Inc., 3rd Floor, BMA, 1-5-5 Minatogino-Minamimachi, Chuo-ku, Kobe 650-0047, Japan



	BTK IC ₅₀ (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 μ M concentration.

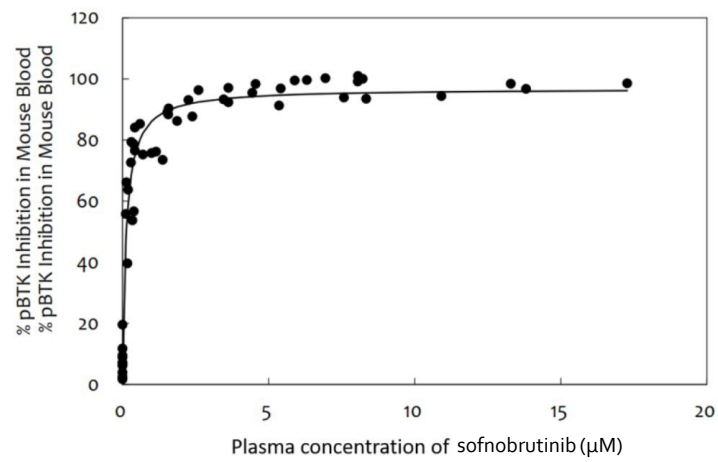
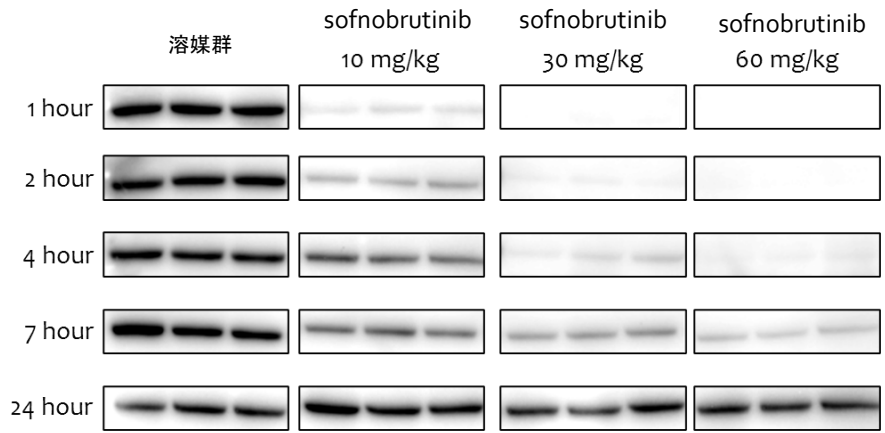




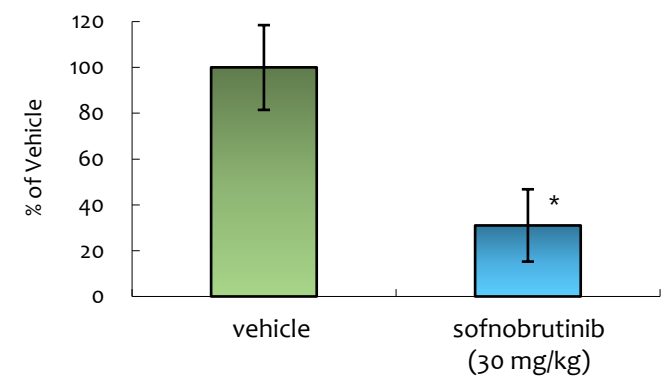
Sofnobrutinib (AS-0871): In Vivo Therapeutic Efficacy

◆ PK/PD Study

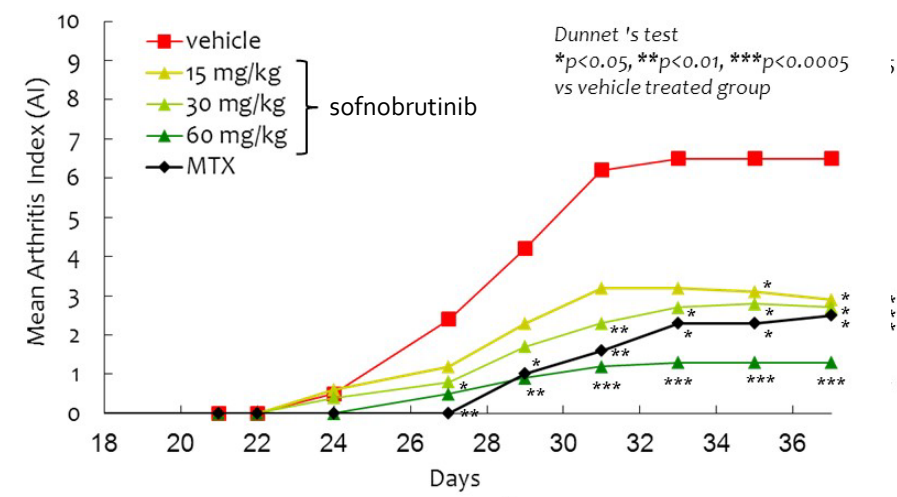
Auto-phosphorylation status of BTK was measured following oral single administration of sofno Brutinib .



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



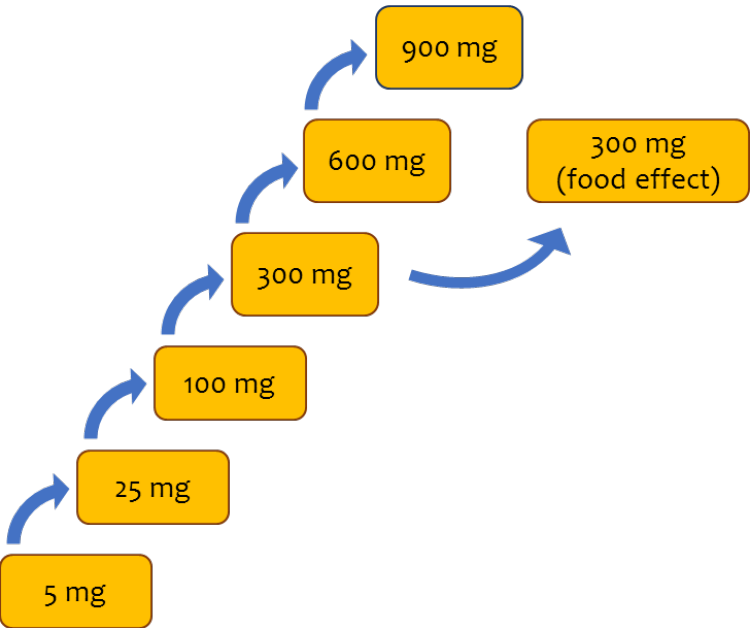
◆ Collagen-induced arthritis (CIA) mouse model (n=10)





SAD Part (Completed)

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



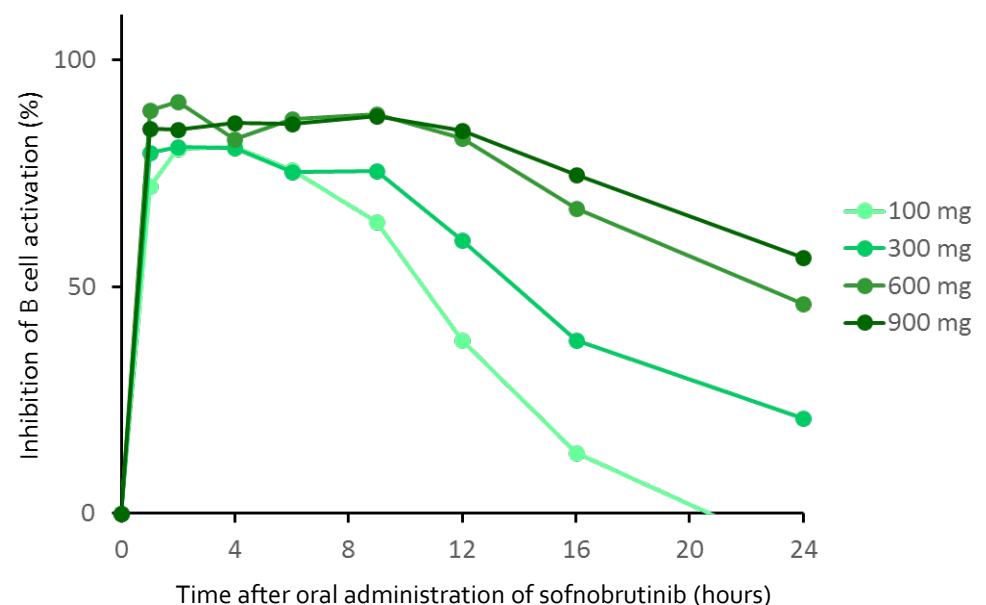
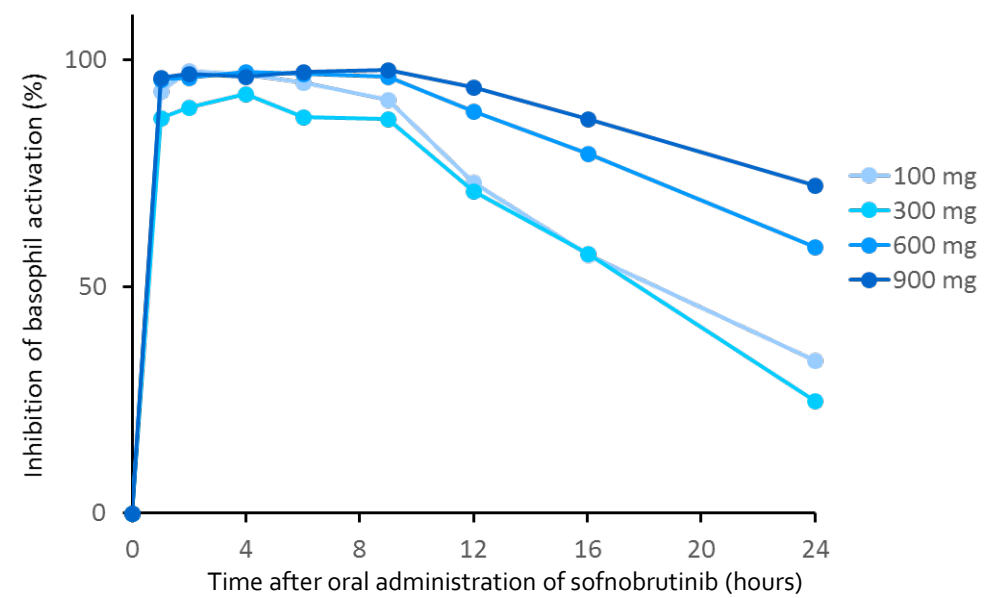
- Sofnobrutinib is well-tolerated without any safety concerns.
- Favorable pharmacokinetic profile.
- Blood samples to assess PD effects were analyzed for evaluation of the B-cell and basophil responses. Administration of sofno Brutinib at 100mg or above resulted in strong inhibition of B-cell and basophil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

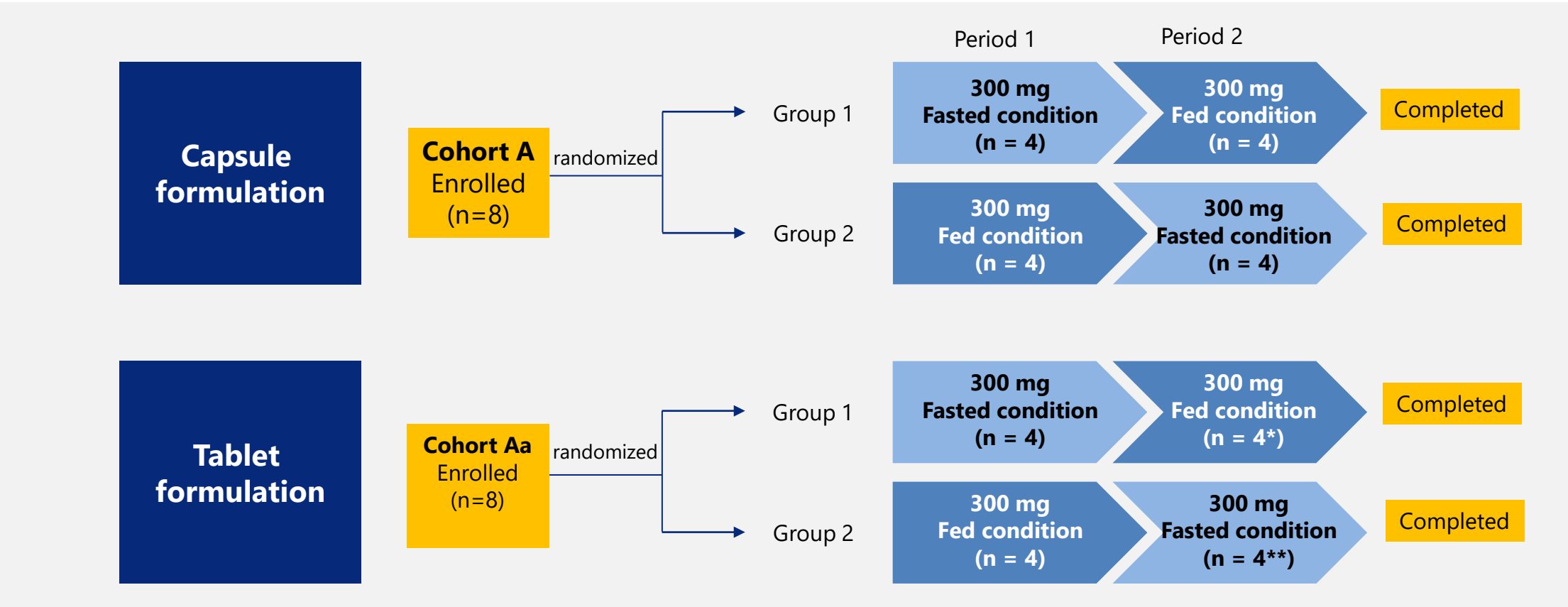




Sofnobrutinib (AS-0871): Phase 1 MAD Study BA part

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.

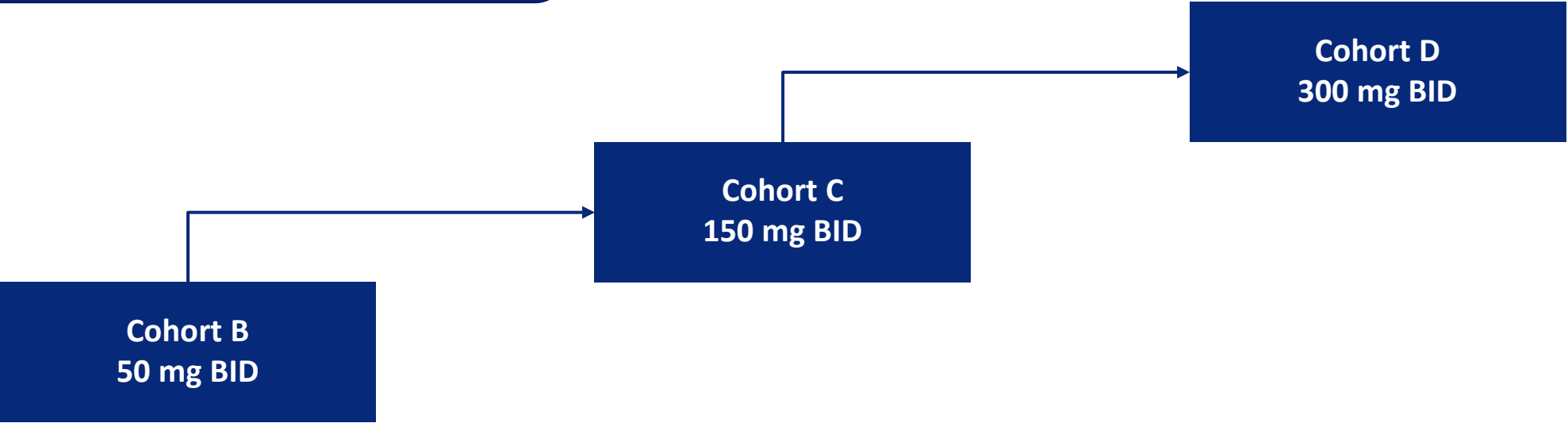
**One subject withdrew from the study due to personal reasons before dosing.



Study Design of MAD part

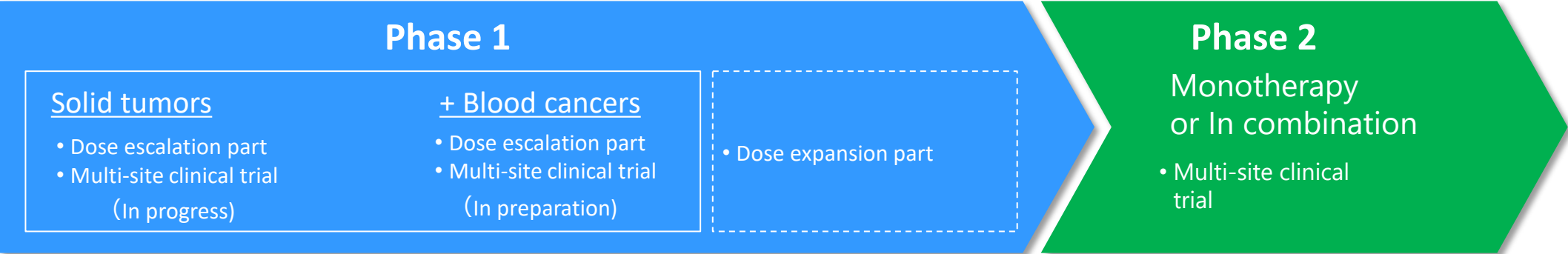
- In the MAD part, safety, tolerability, PK, and PD of 3 multiple ascending doses of sofno Brutinib, following 14-day multiple dose oral administration of sofno Brutinib, will be investigated using a double blind, placebo-controlled, randomized design in 3 cohorts of 8 healthy subjects each.
- ◆ Dosing 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





monzosertib : Targeting Cancer	
<ul style="list-style-type: none">● Small molecule CDC7 inhibitor● High kinase selectivity● Potential First-in-class drug● Orally available	<ul style="list-style-type: none">● Potent anti-proliferative activity against various cancer cell lines● Demonstrated strong anti-tumor activity in several human tumor xenograft models● Conducting Phase 1 study in Japan targeting solid tumors● Planning to expand to blood cancers as monotherapy or in combination with other drugs





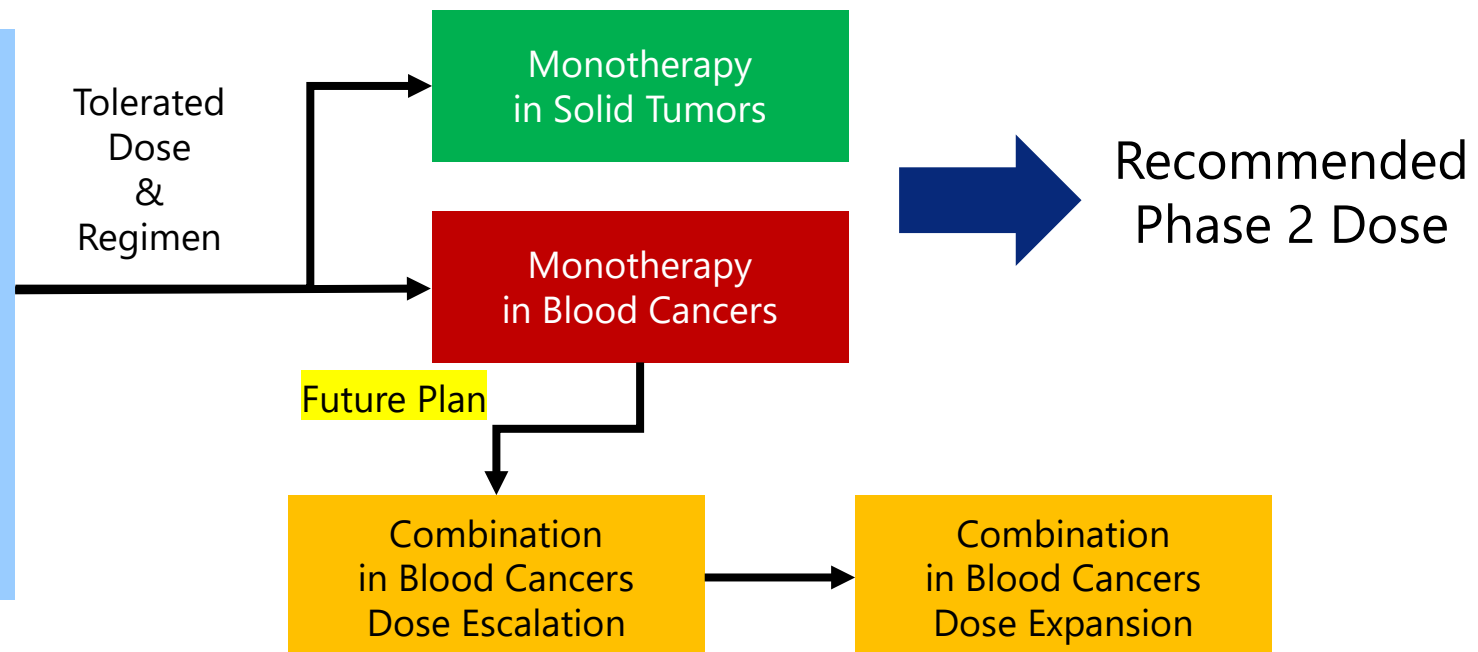
Phase 1 study targeting cancer patients

- Phase 1 study in patients with unresectable, advanced, recurrent, or metastatic solid tumors was initiated in Japan in 2021.
- 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-80 mg BID and above.

Dose Escalation Part in Advanced Solid Tumors



Dose Expansion Part



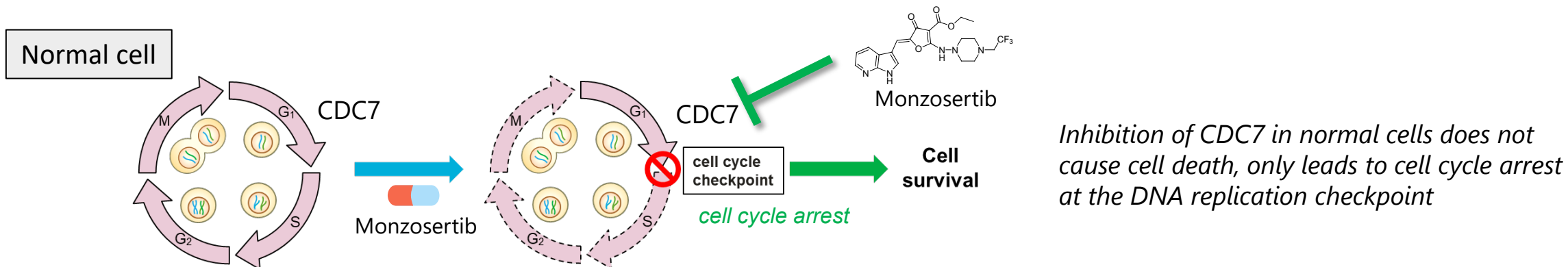


Monzosertib (AS-0141)

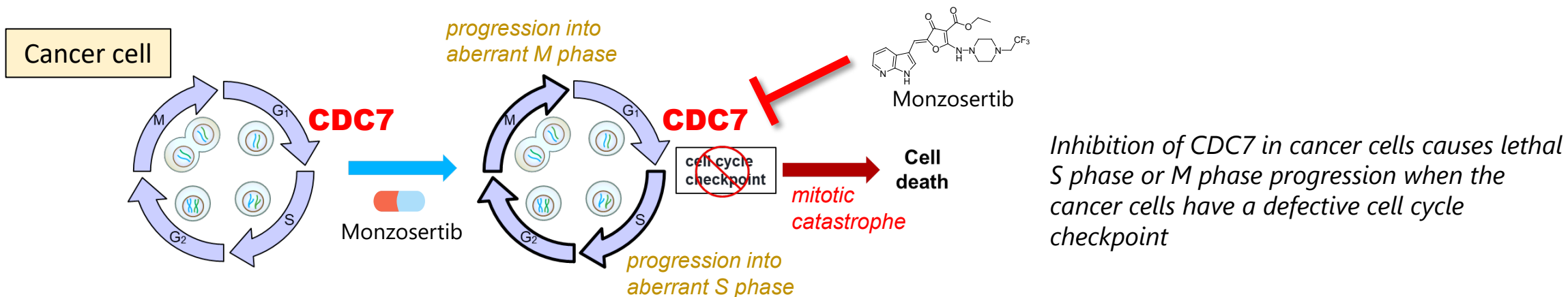
Potential First-in-class CDC7 Inhibitor Targeting Cancer

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





◆ 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^{a,*}, Tokiko Asami^a, Ayako Sawa^a, Yuko Uno^a, Mitsuharu Hanada^a, Chika Taniyama^b, Yoko Funakoshi^b, Hisao Masai^c, Masaaki Sawa^a

^a Research and Development, Carina Biosciences, Inc., 3F BMS, 1-5-5 Minatogawa-Minamimachi, Chuo-ku, Kobe, 650-0047, Japan
^b Research and Development Department, SRI Biotech Co., Ltd., Izumi Garden Tower B5F, 1-6-1 Rappongi, Minato-ku, Tokyo 106-8018, Japan
^c Department of Genome Medicine, Tokyo Metropolitan Institute of Medical Science, 2-1-6 Minamishinboku, Setagaya-ku, Tokyo 158-8506, Japan

Journal of
Medicinal
Chemistry

pubs.acs.org/jmc

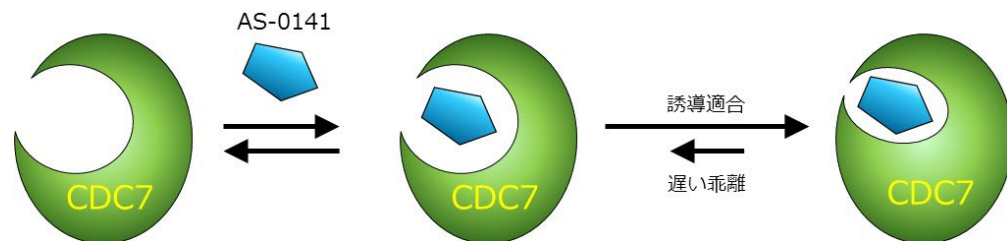
Drug Annotation

Discovery of AS-0141, a Potent and Selective Inhibitor of CDC7 Kinase for the Treatment of Solid Cancers

Takayuki Irie^{a,*}, Tokiko Asami, Ayako Sawa, Yuko Uno, Chika Taniyama, Yoko Funakoshi, Hisao Masai, and Masaaki Sawa

Cite This: *J. Med. Chem.* 2021, 64, 14153–14164

Read Online



Inhibitory potency (IC₅₀) for CDC7 in the presence of 1 mM ATP

Without
Preincubation

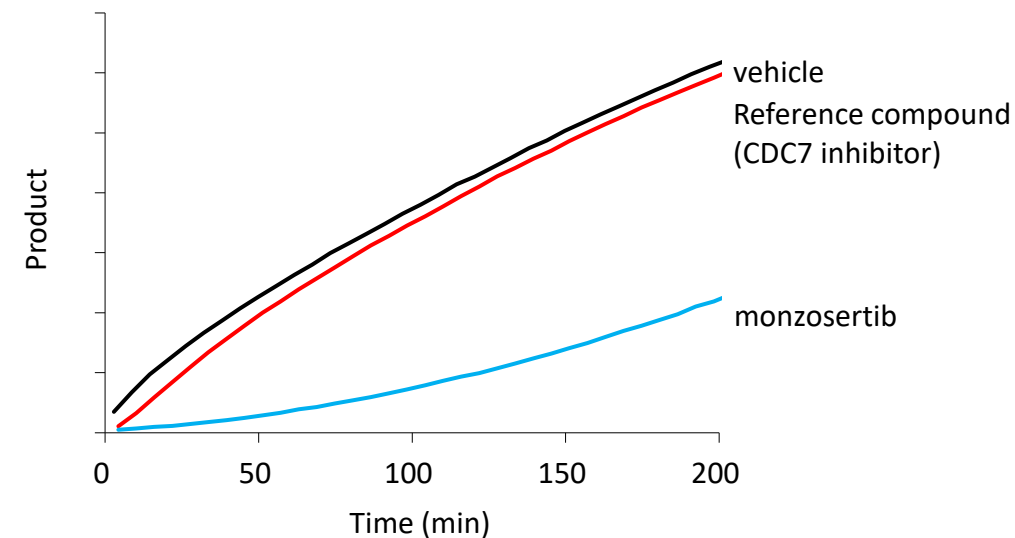
503 nM

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.



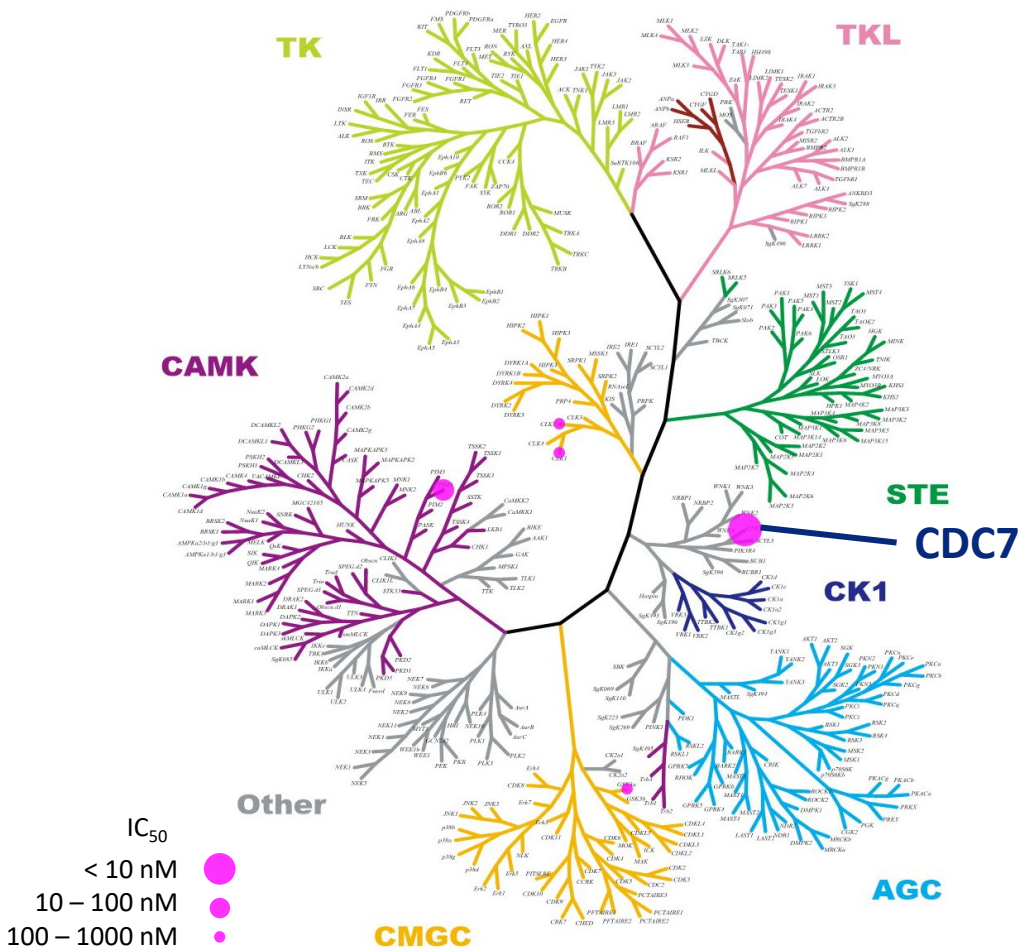


Monzosertib (AS-0141): High Kinase Selectivity



Kinase Selectivity Profiling

In the presence of 1 mM ATP with preincubation



IC50 values of hit kinases (at 1 mM ATP)

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

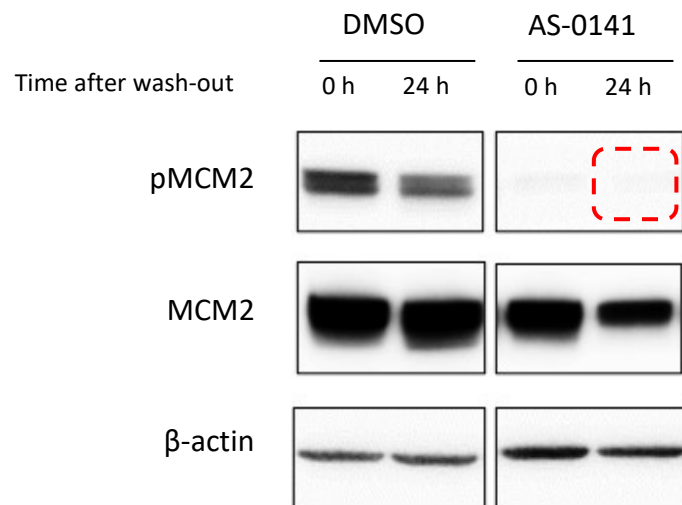
CDC7 is the only kinase that shows preincubation effect



Monzosertib (AS-0141): Strong Cellular Activity

◆ Prolonged inhibition in cells

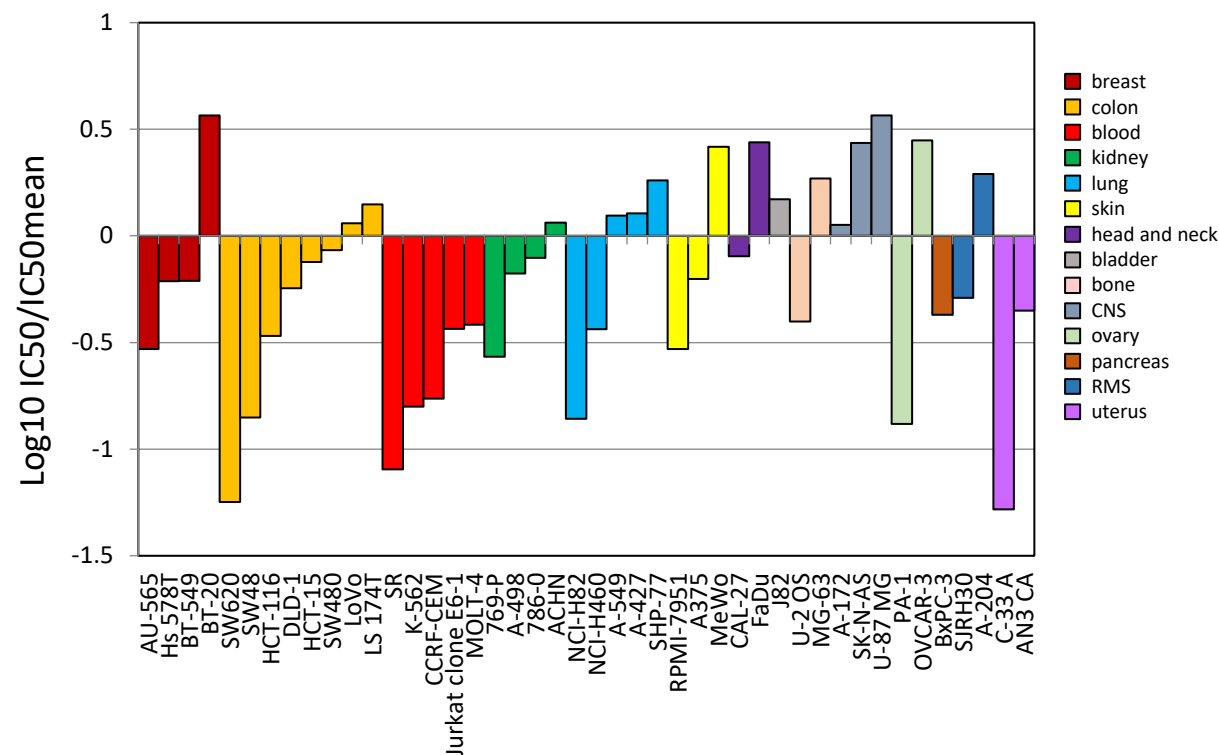
Human colon cancer cell line, Colo-205 cells were treated with DMSO control or monzosertib. After washout of the inhibitor, the cells were further incubated in the same media for 0 or 24 h and subjected to western blot analysis.



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

J Med Chem. 2021 Oct 14;64(19):14153-14164.

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

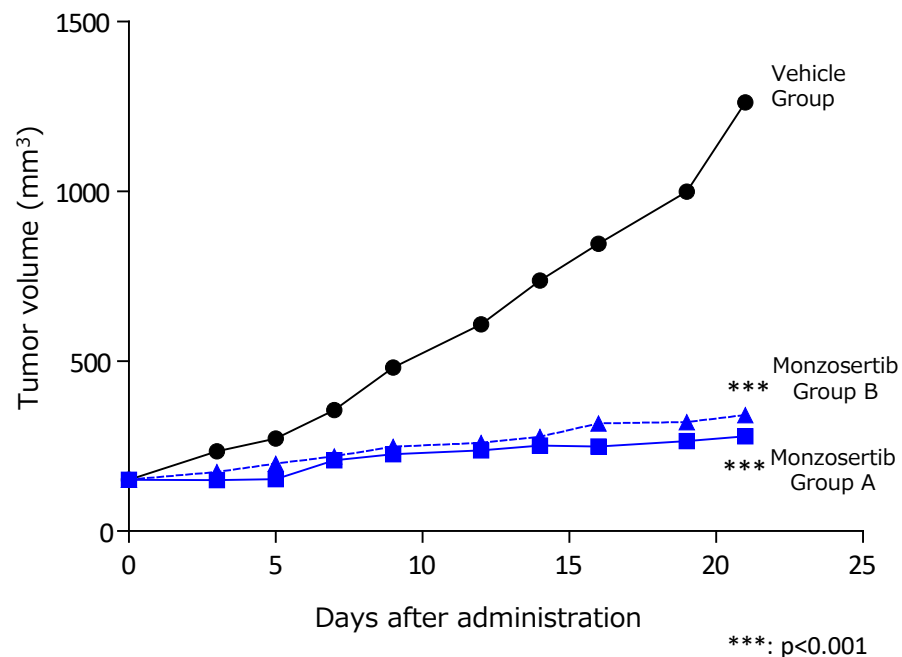


44 Cancer cell lines (Oncolines at NTRC)

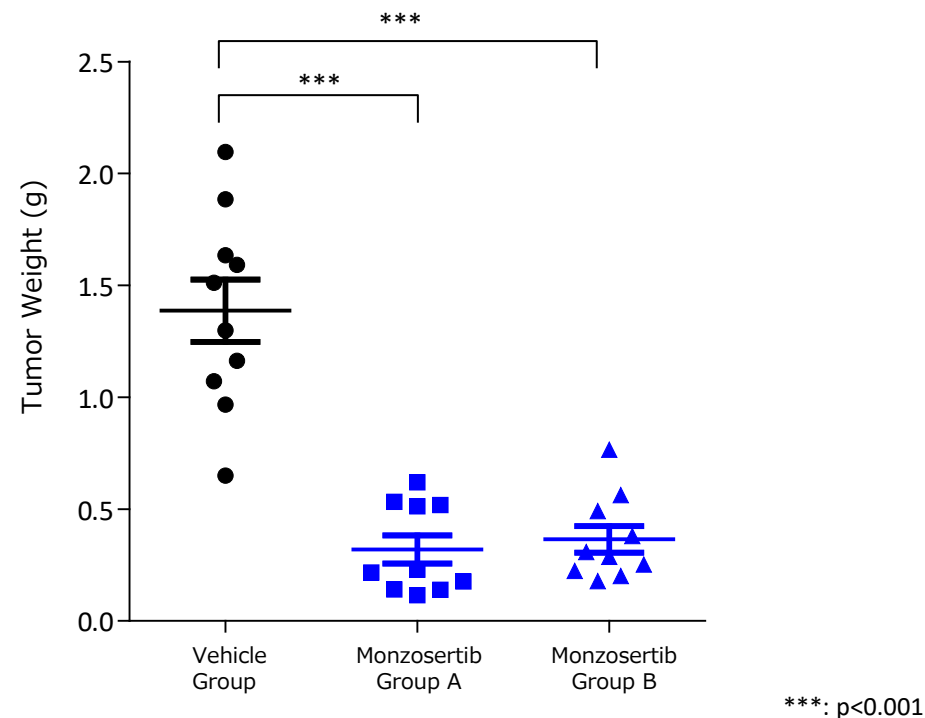


In vivo antitumor efficacy of monzosertib in a SW620 (human colon cancer) xenograft mouse model

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.