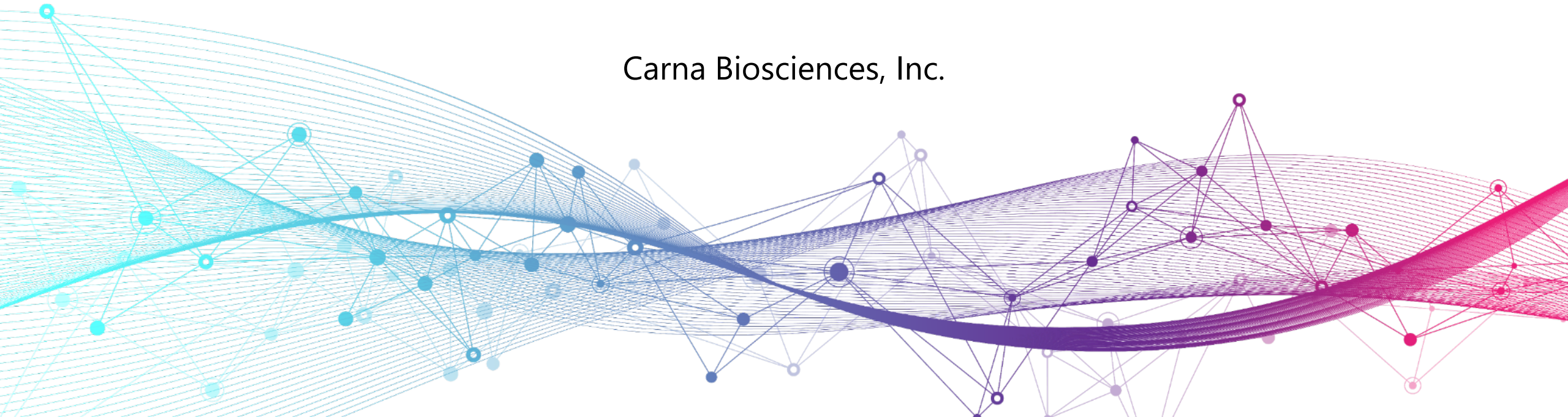


Financial Results FY2024

(January to December 2024)

Carna Biosciences, Inc.



AGENDA

1

Updates on Pipelines in Clinical Development

2

Updates on Licensed Pipelines

3

FY2024 Results

4

Business Plan for FY2025

5

Appendix



Updates on Pipelines in Clinical Development

- 1 docirbrutinib (AS-1763)**
- 2 sofnobrutinib (AS-0871)**
- 3 monzosertib (AS-0141)**

International Nonproprietary Name (INN) : docirbrutinib, Code name : AS-1763
International Nonproprietary Name (INN) : sofnobrutinib, Code name : AS-0871
International Nonproprietary Name (INN) : monzosertib, Code name : AS-0141



Compound	Target	Indication	Status
docirbrutinib (AS-1763)	BTK	Blood Cancer	<ul style="list-style-type: none">Phase 1b clinical trial is ongoing in the U.S.Dose expansion part was initiated in October to accelerate the development timeline.Dose escalation part was completed in December.Encouraging preliminary data was presented at ASH 2024 in December. <div><p>Multi-center clinical study Study Lead : Prof. Nitin Jain, MD, Department of Leukemia, University of Texas MD Anderson Cancer Center.</p></div>
sofnobrutinib (AS-0871)	BTK	Immune-inflammatory diseases	<ul style="list-style-type: none">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.Performed a preclinical study (EFD study) to establish a best-in-class status; potential advantages of sofnobrutinib over other BTK inhibitors.Seeking a strategic partner for further development.
monzosertib (AS-0141)	CDC7/ ASK	Cancer	<ul style="list-style-type: none">Phase 1 clinical trial in cancer patients is ongoing in Japan.Solid tumor : Completed dose escalation part Initiated patient recruitment in dose expansion part in January 2025.Blood cancer : Dose escalation part is ongoing. <div><p>Clinical trial sites</p><ul style="list-style-type: none">- National Cancer Center Hospital and National Cancer Center Hospital East- The Cancer Institute Hospital of JFCR</div>



Docirbrutinib (AS-1763): Highlights



Key Highlights

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

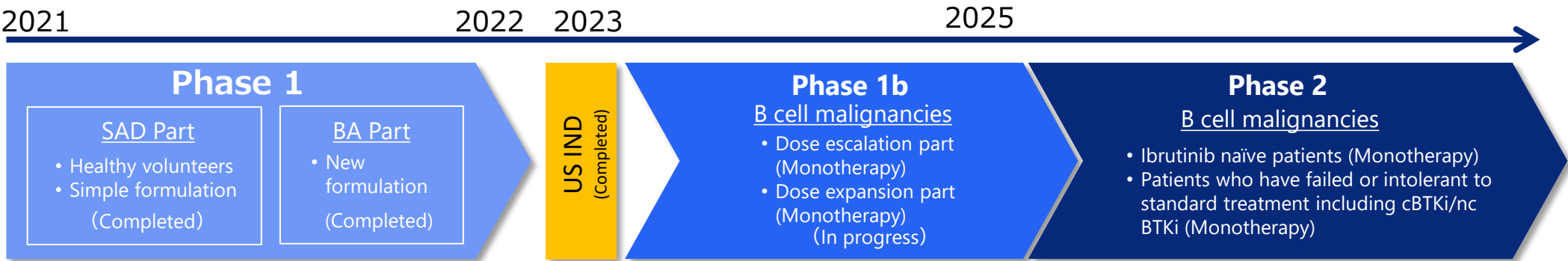
- ✓ Indication : CLL/SLL and B-cell NHL
- ✓ Non-covalent BTK inhibitor
- ✓ Docirbrutinib has a potential to be effective for patients who have developed resistance to the existing BTK inhibitors.

Status

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

- ✓ Dose escalation part was completed in December 2024.
- ✓ Dose expansion part was initiated to accelerate the development timeline in October 2024.



IND application : Investigational New Drug application 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: non-covalent BTK inhibitor



Docirbrutinib (AS-1763): Phase 1b Trial Design



Dose Escalation

Completed

CLL/SLL and B-cell NHL

3+3 design

100 mg BID

500 mg BID

RP2D: Recommended phase 2
BID: Twice a day

Dose Expansion

Cohort 1 (CLL/SLL)

High dose 500 mg BID

Medium dose 400 mg BID

Low dose 300 mg BID

In progress

Cohort 2 (B-cell NHL)

High dose 500 mg BID

Medium dose 400 mg BID

Low dose 300 mg BID

In progress

Cohort 3 (Prior therapy with pirtobrutinib for CLL/SLL or B-cell NHL)

High dose 500 mg BID

Medium dose 400 mg BID

In progress

RP2D

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.



Clinical sites (As of January 2025)

- UC Irvine Health
- Mount Sinai Comprehensive Cancer Center
- Moffitt Cancer Center
- Northwestern Memorial Hospital
- University of Maryland Medical Center-Greenebaum Comprehensive Cancer Center
- University of Massachusetts Memorial Medical Center
- Clinical Research Alliance, Inc.
- University of Texas MD Anderson Cancer Center
- The Medical College of Wisconsin

- ✓ **Phase 1b study is ongoing at nine top clinical sites in the US.**
- ✓ **Planning to activate additional clinical sites to accelerate the development timeline.**



CLL Society

CLL Society is the world's leading authority for chronic lymphocytic leukemia and small lymphocytic lymphoma cancer patients.

Mission

CLL Society is an inclusive, patient-centric, physician-curated nonprofit organization that addresses the unmet needs of the chronic lymphocytic leukemia and small lymphocytic lymphoma (CLL / SLL) community through patient education, advocacy, support, and research.

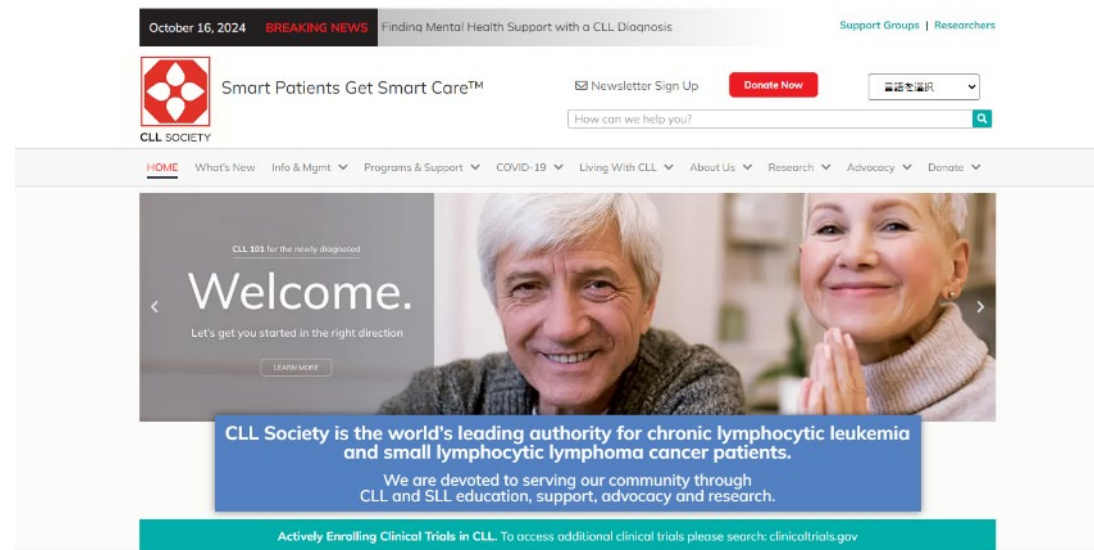
Vision

We envision a world in which the entire CLL / SLL community can equitably access quality education, support, and care, to lead healthier and richer lives.

– Cited from the website of CLL Society –

Carna offered sponsorship at the request of CLL Society.

<https://cllsociety.org/>



A clinical study for BTK inhibitors resistance
Docirbrutinib: an Investigational Oral BTK Inhibitor

Currently enrolling patients with CLL/SLL and B-cell NHL

LEARN MORE

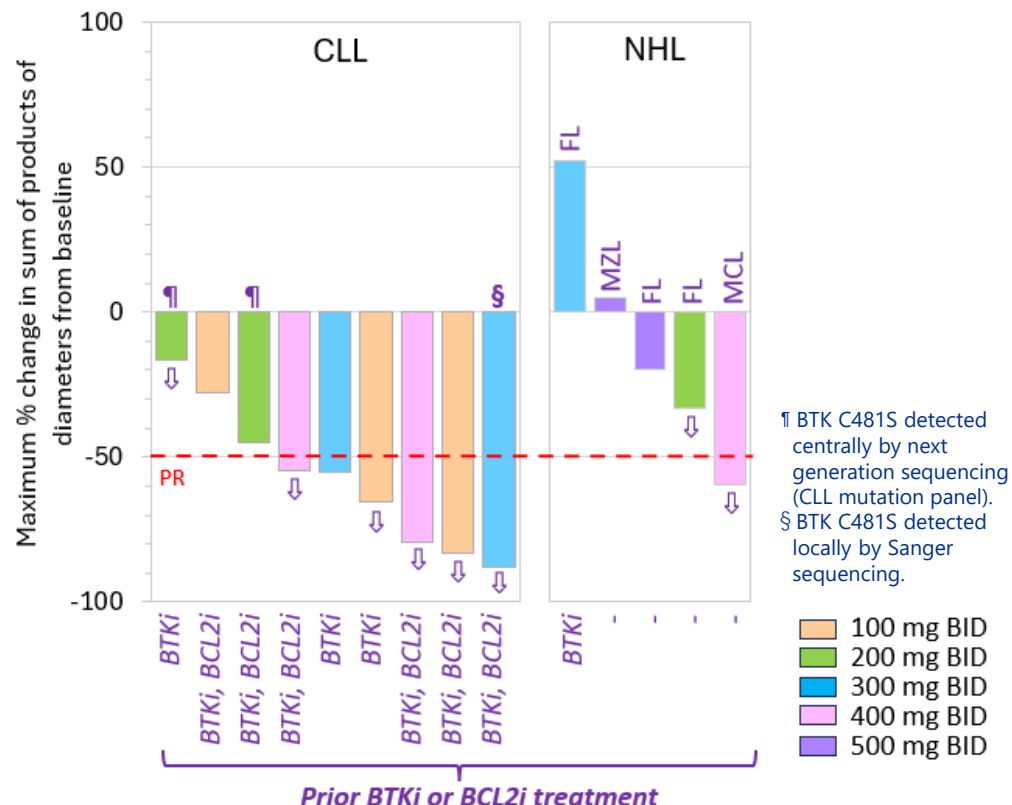


A clinical trial ad for docirbrutinib is on CLL Society's website to promote patient enrollment.

Preliminary Phase 1b data (1)

Data from ASH2024 poster presentation
The 66th American Society of Hematology Annual Meeting & Exposition, December 7-10, 2024

Tumor response (lymph node size)



CLL: Chronic Lymphocytic Leukemia, NHL: non-Hodgkin Lymphoma, FL: Follicular Lymphoma, MCL: Mantle cell lymphoma, MZL: Marginal zone lymphoma, BTKi: BTK inhibitor, BCL2i: BCL2 inhibitor, BID: Twice a day

Efficacy of docirbrutinib

CLL: All patients experienced lymph node size reduction. 6 out of 9 evaluable patients (67%) with CLL achieved PR or PR-L with 50% reduction in lymph node size. The exposures at ≥ 300 mg BID exceeded the IC_{90} throughout the dosing interval, and all 4 CLL patients (100%) receiving ≥ 300 mg BID achieved PR or PR-L.

NHL: 3 out of 5 patients with NHL experienced lymph node size reduction. One MCL patient experienced PR with over 50% reduction in lymph node size.

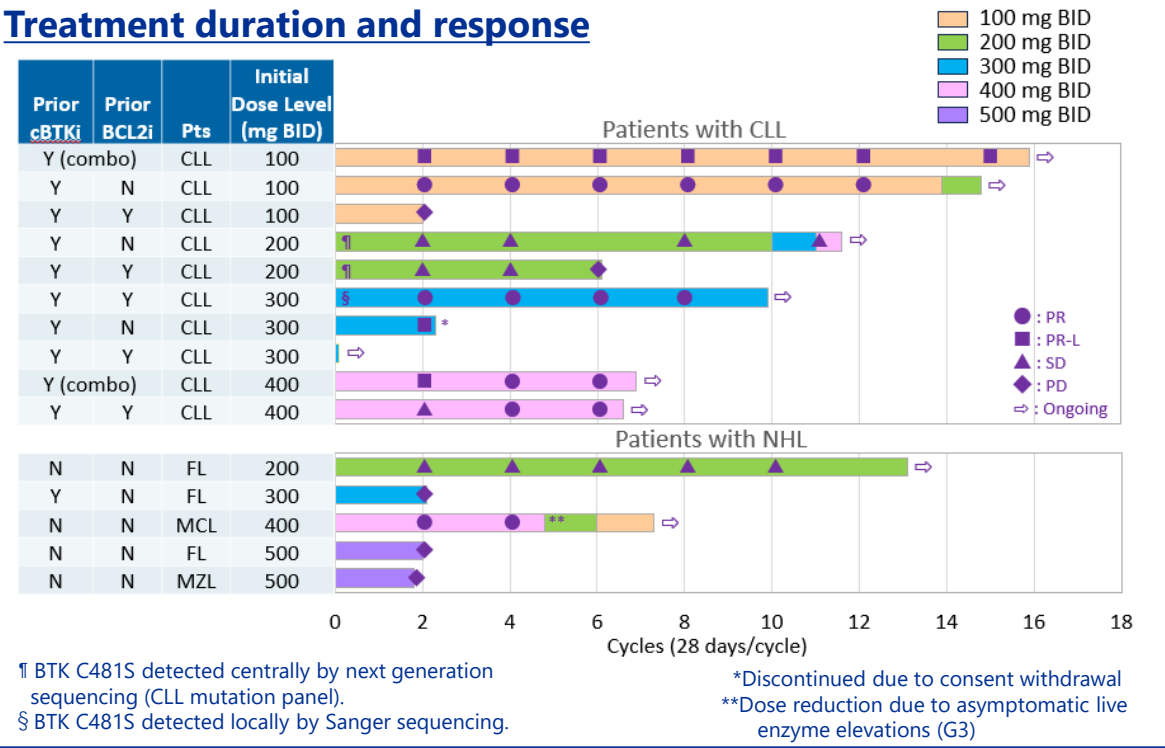
These preliminary data indicate clinical responses in patients with B-cell malignancies.



Preliminary Phase 1b data (2)

Data from ASH2024 poster presentation

Treatment duration and response



- The bar charts indicate the duration of treatment.
- The color of the bar charts indicates dose levels. The change in the color indicates that the dose level was changed during the treatment.
- Arrows indicate that the treatment is ongoing.
- PR (Partial Response) indicates that the size of lymph nodes has decreased $\geq 50\%$ and other parameters for PR, e.g. reduction in the number of lymphocytes, have been met.
- PR-L (PR-Lymphocytosis) indicates that the size of lymph nodes has decreased $\geq 50\%$ but the reduction in the number of lymphocytes has not met the criteria for PR.
- SD (Stable Disease) indicates that the disease remains stable.
- PD (Progression Disease) indicates that the disease has progressed.

CLL: Chronic Lymphocytic Leukemia, NHL: non-Hodgkin Lymphoma, FL: Follicular Lymphoma
MCL: Mantle cell lymphoma, MZL: Marginal zone lymphoma, BTKi: BTK inhibitor, BCL2i: BCL2 inhibitor, BID: Twice a day

Safety profile of docirbrutinib

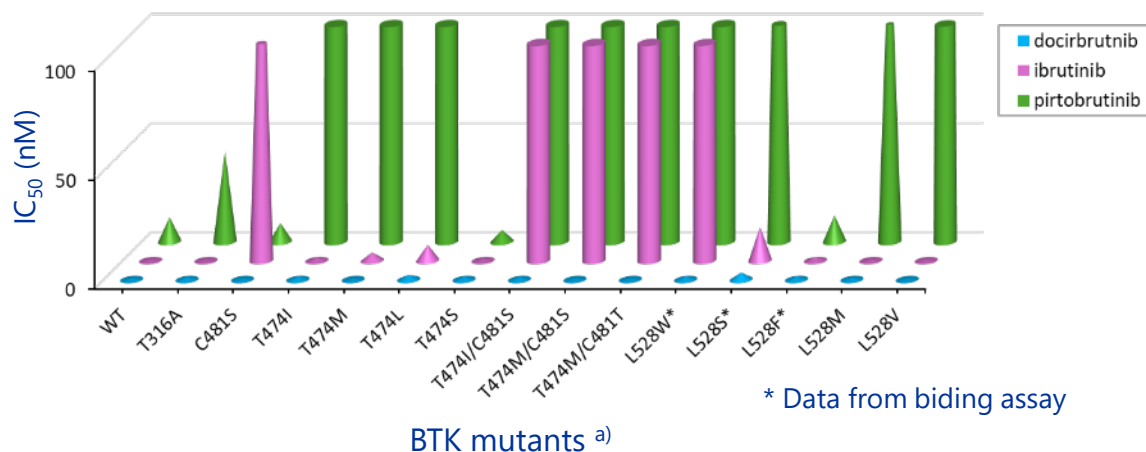
- No dose-limiting toxicities were observed. No treatment discontinuation due to AEs and no drug-related atrial fibrillation or hypertension were reported at doses of 100-500 mg BID.
- Asymptomatic ALT/AST elevations (G3) were reported in one patient (7%) as the only drug-related $\geq G3$ AEs other than neutrophil count decrease, indicating a favorable safety profile of docirbrutinib.

These preliminary data indicate a favorable safety profile and clinical responses in heavily pretreated patients with B-cell malignancies.



Preclinical study

Inhibitory potency of BTK inhibitors against BTK mutants



^{a)} Patients are reported to develop resistance during the treatment with covalent BTK inhibitors including ibrutinib due to substitution of cysteine residue at 481 position with serine (C481S mutation) in BTK, which reduces the efficacy of covalent BTK inhibitors. In addition, the emergence of other types of mutations, such as T474x and L528x, has been reported during the treatment with pirtobrutinib.

In vitro study using recombinant BTK mutant proteins

The bar charts show the comparison of the inhibitory potency of BTK inhibitors against various BTK mutants. A shorter bar indicates stronger potency.

- Ibrutinib and pirtobrutinib showed weak inhibitory potency against many resistant BTK mutants.
- Docirbrutinib showed strong inhibitory potency against all resistant BTK mutants.

Docirbrutinib is expected to be effective against patients who have developed resistance to the existing BTK inhibitors.

Drug resistance: the reduction in effectiveness of a drug during targeted therapies due to alterations of drug targets including the mutation of the target proteins.



Safety profile of docirbrutinib

Preliminary data from Phase 1b study

Data from ASH2024 poster presentation

Treatment-Emergent Adverse Event (TEAE)	All Doses and Pts (n=15)			
	Any		Treatment-related	
	Any Grades n (%)	Grade ≥3 n (%)	Any Grades n (%)	Grade ≥3 n (%)
≥15%				
Dizziness	9 (60%)	0	2 (13%)	0
Headache	6 (40%)	0	1 (7%)	0
Nausea	5 (33%)	0	2 (13%)	0
Neutrophil count decreased	5 (33%)	2 (13%)	4 (27%)	2 (13%)
Blood creatinine increased	4 (27%)	0	0	0
Fatigue	4 (27%)	0	1 (7%)	0
Abdominal pain	3 (20%)	0	0	0
Anemia	3 (20%)	0	1 (7%)	0
Constipation	3 (20%)	0	1 (7%)	0
Cough	3 (20%)	0	0	0
Fever	3 (20%)	0	0	0
Myalgia	3 (20%)	0	0	0
TEAEs of Special Interest				
Bruising ^a	2 (13%)	0	1 (7%)	0
Hemorrhage ^b	1 (7%)	0	1 (7%)	0

The preliminary data from the Phase 1b study indicates impressive safety profile of docirbrutinib.

^a Contusion (G1), ^b Epistaxis (G1).



Selected BTK inhibitors in clinical development

Compound	Type	Company (Originator)	Development Phase
pirtobrutinib (LOXO-305)	Non-covalent BTK inhibitor	Lilly (Loxo)	Approved/P3
nemtabrutinib (ARQ 531)	Non-covalent BTK inhibitor	Merck (ArQule)	P3
NX-5948	BTK degrader	Nurix	P1
BGB-16673	BTK degrader	BeiGene	P1
docirbrutinib (AS-1763)	Non-covalent BTK inhibitor	Carna	P1

Carna plans to accelerate the enrollment in the dose expansion part of the Phase 1b study and accumulate the clinical data to demonstrate:

Best-in-Class BTKi

- ✓ Safer profile compared with other BTKi
- ✓ Effective against resistant mutants



***Discontinuation of ibrutinib treatment
is commonly due to intolerance***

The discontinuation rate of ibrutinib treatment was
41% in the US, and the majority reasons of
discontinuation was AEs.

Reasons for ibrutinib discontinuation	Ibrutinib in front-line	Ibrutinib in relapse
Toxicity	63.1%	50.2%
CLL Progression	15.8%	20.9%
Others	21.1%	28.9%

Data cited from Mato AR, et al., Haematologica. 2018;103(5):874-879

Docirbrutinib (AS-1763) in Phase 1b :

No discontinuation due to adverse events so far



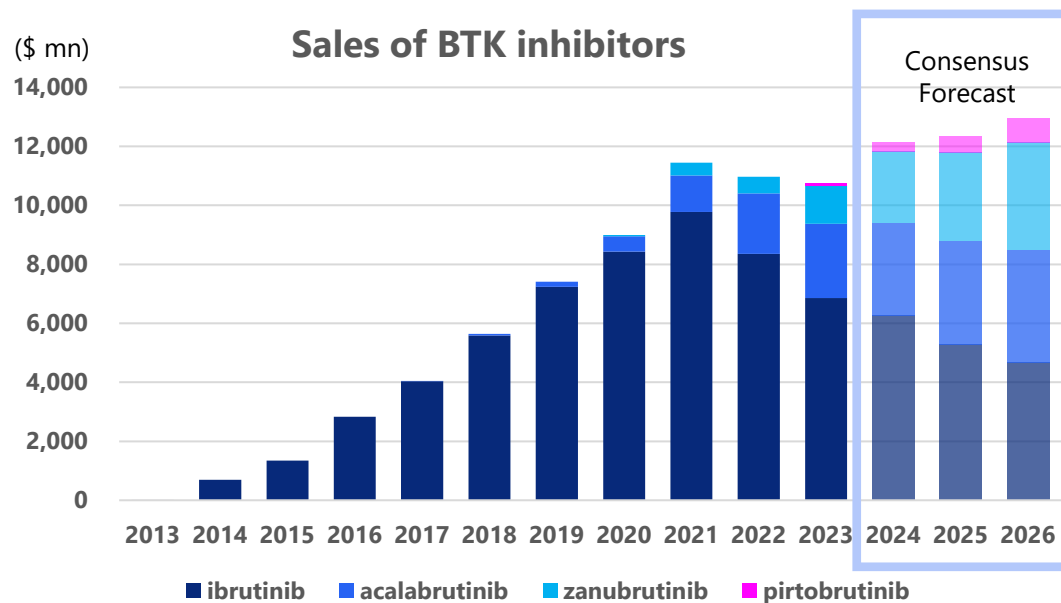
**Docirbrutinib demonstrates safer profiles,
suggesting better option for effective therapy**



BTK inhibitors form a large market

Sales of the currently approved BTK inhibitors, ibrutinib, acalabrutinib, zanubrutinib, and pirtobrutinib, exceed \$10 bn.

- Acalabrutinib, zanubrutinib, and pirtobrutinib are taking a market share from ibrutinib as safer BTK inhibitors.
- Acquired resistance against ibrutinib, acalabrutinib, zanubrutinib, and pirtobrutinib present a need for new therapeutic options



Source: Clarivate

Product Positioning of docirbrutinib

Offer a new therapeutic option to:

- *patients who discontinued the existing BTK inhibitors due to adverse events as a safer BTK inhibitor*
- *patients who have developed resistance to the existing BTK inhibitors as a pan-mutant BTK inhibitor*

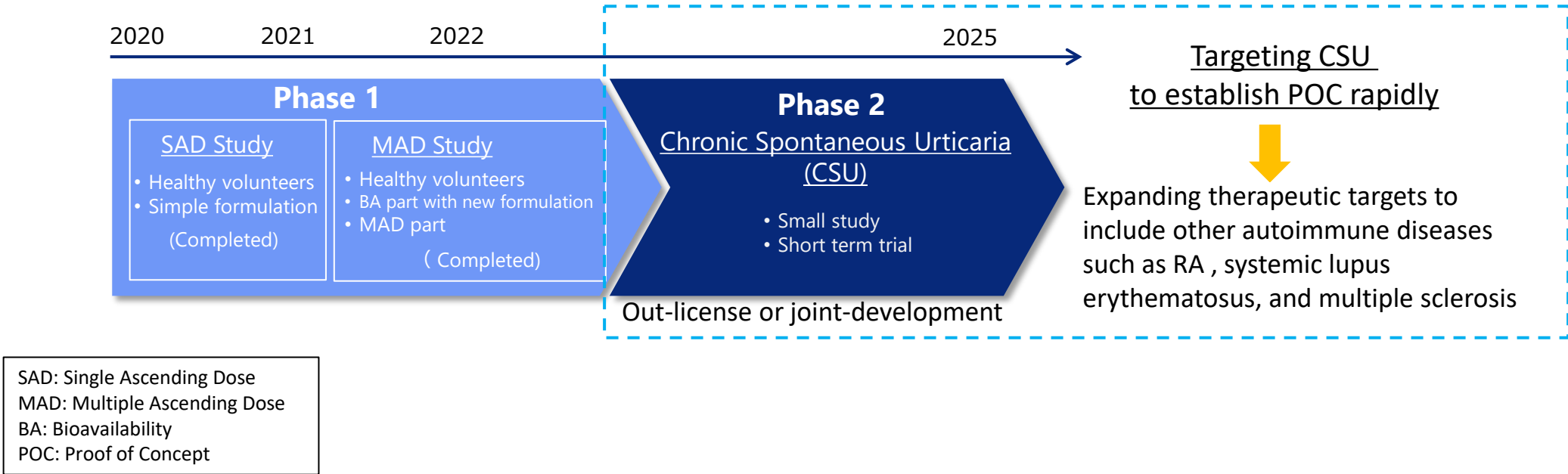


Mechanism/ Indication	Sofnobrutinib is an orally available Bruton's Tyrosine Kinase (BTK) inhibitor to treat autoimmune diseases by inhibiting activation of immune cells such as B cells, macrophages, and mast cells.
Clinical trials	<u>Completed Phase 1 clinical trial in healthy volunteers in the Netherlands.</u> A favorable safety and tolerability profile as well as a promising PK/PD profile were confirmed and these results support to advance sofno Brutinib into Phase 2 clinical development.
Status	<ul style="list-style-type: none">✓ Performed a preclinical study to establish a best-in-class status; potential advantages of sofno Brutinib over other BTK inhibitors.✓ Seeking a strategic partner for further development.



Sofnobrutinib : Targeting Immune-inflammatory diseases

- Small molecule BTK inhibitor
- Non-covalent/reversible
- High kinase selectivity
- Orally available
- Demonstrated significant efficacies in arthritis models
- Showed efficacy in systemic lupus erythematosus model
- Phase 1 Clinical Trial was completed
- Find a partner to conduct further development
- Performed a preclinical study to establish a best-in-class status





Embryo-Fetal Development (EFD) toxicity study was performed to prove potential advantages of sofnobrutinib over other BTK inhibitors.

Sofnobrutinib showed "No Teratogenic Effect" in the EFD study, suggesting it is suitable for the treatment of dermatologic diseases including CSU.

As most BTK inhibitors approved are teratogenic, their use should be limited especially for women.

Sofnobrutinib is confirmed to be non-teratogenic in the EFD toxicity study, providing a treatment option for a wider range of patients.

Sofnobrutinib is the only BTK inhibitor having a non-covalent inhibitory mechanism of action with no teratogenic effect.



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.

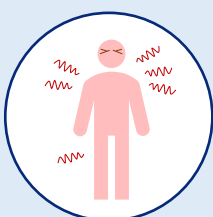


Chronic Spontaneous Urticaria (CSU)

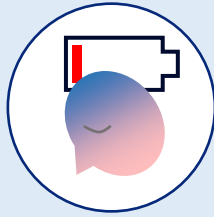
CSU is a debilitating disease of chronic itch, hives and angioedema, lasting six weeks or more.

Symptoms

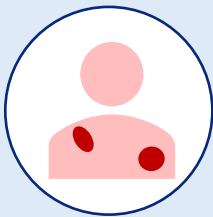
There is no specific external trigger for CSU, but the autoimmune system may play a role



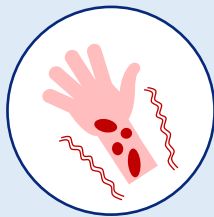
Spontaneously present & re-occur



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

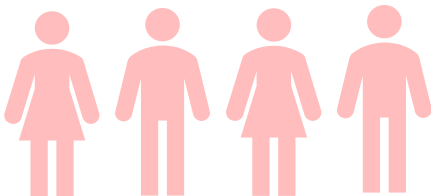


Red swollen hives



Itch

Number of Patients



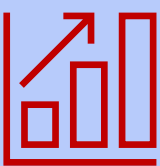
2.8 million

diagnosed prevalent cases in major seven markets

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

- ✓ 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.

Market Size



\$2,844 million

in 2020 in seven major markets

- ✓ The market size of CSU in major seven countries is expected to reach \$8,043 mn by 2030

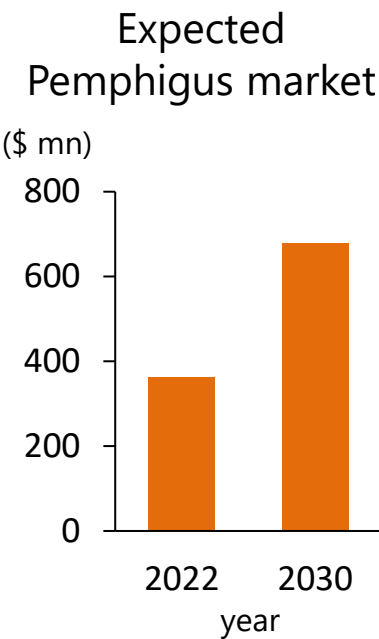
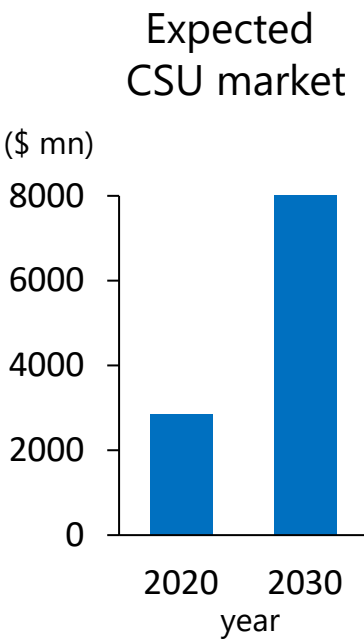
<https://www.delveinsight.com/>
Source : Clarivate



Initial focus

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

*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
Source : Clarivate



Highlights

Monzosertib is an orally available CDC7 kinase inhibitor targeting cancer

Clinical trials in progress

Conducting Phase 1 study in Japan targeting solid tumors and blood cancers

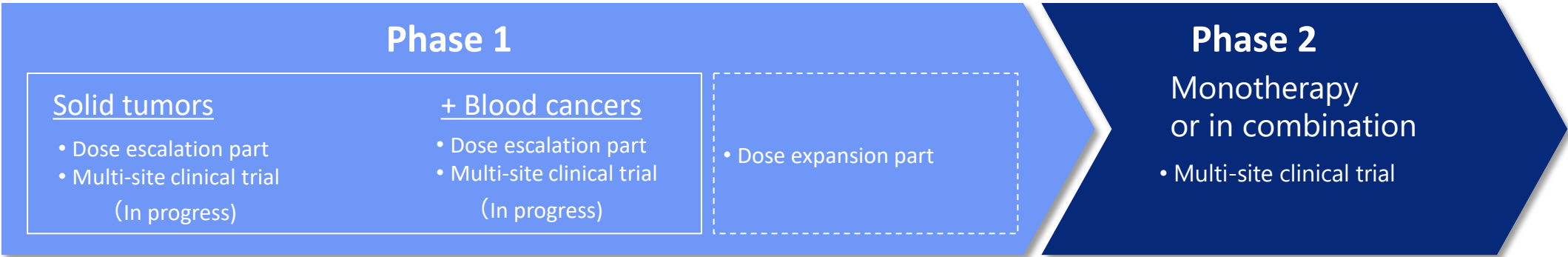
- ✓ Solid tumor : Completed dose escalation part.
Initiated patient recruitment in dose expansion part in January 2025.
- ✓ Blood cancer : Dose escalation part is ongoing.

Clinical trial sites

- National Cancer Center Hospital and National Cancer Center Hospital East
- The Cancer Institute Hospital of JFCR will be activated in dose expansion part.



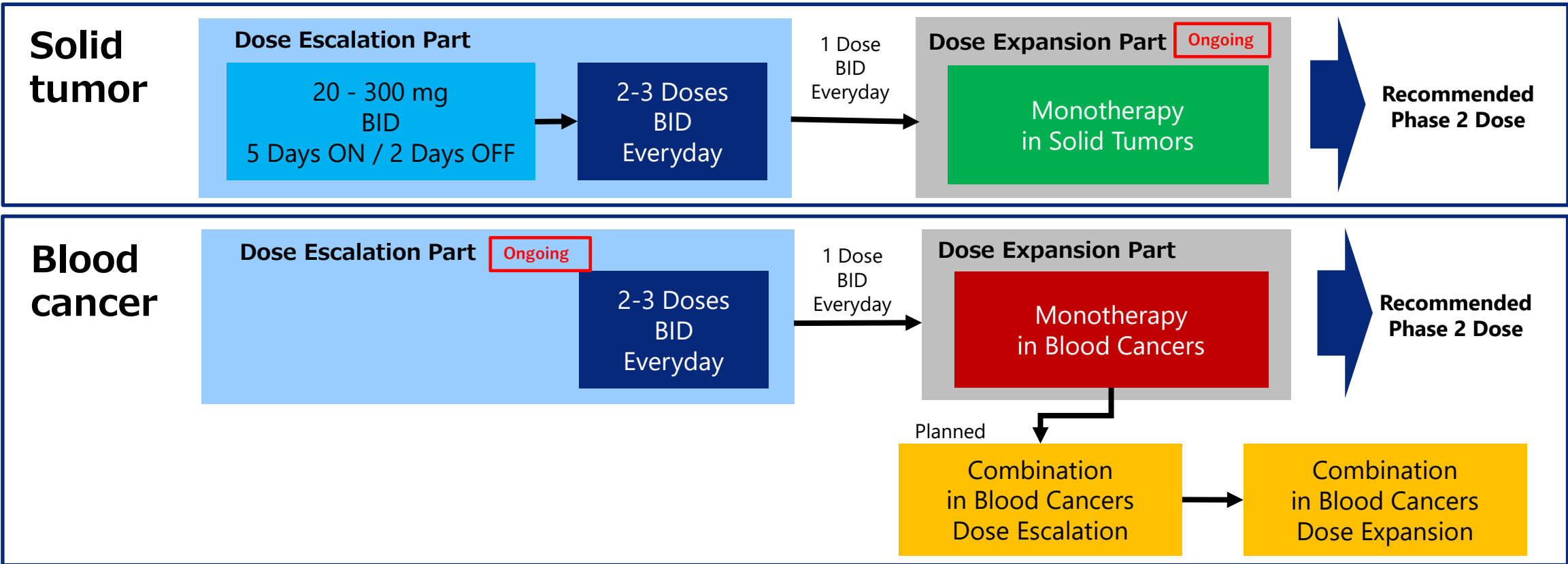
Monzosertib : Targeting Cancer	
<ul style="list-style-type: none">● Small molecule CDC7 inhibitor● High kinase selectivity● Potential First-in-class drug● Orally available	<ul style="list-style-type: none">● Potent anti-proliferative activity against various cancer cell lines● Demonstrated strong anti-tumor activity in several human tumor xenograft models● Conducting Phase 1 study in Japan targeting solid tumors and blood cancers<ul style="list-style-type: none">Solid tumor : Initiated patient recruitment in dose expansion part in January 2025.Blood cancer : Dose escalation part is ongoing.





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.





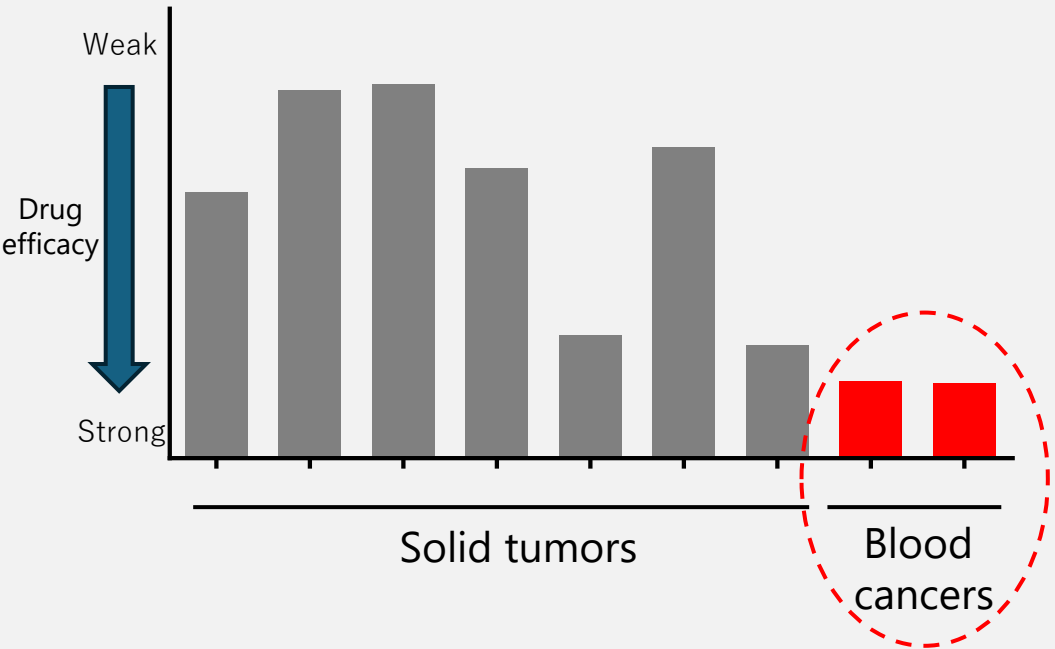
Monzosertib (AS-0141): Dose escalation part targeting blood cancer is ongoing



Non-clinical study



Antiproliferative effects of monzosertib on 35 human cancer cell lines
(Each bar are presented as mean of different cell lines (N = 1 – 7))



✓ Monzosertib demonstrated robust tumor growth inhibition in a human AML xenograft mouse model.

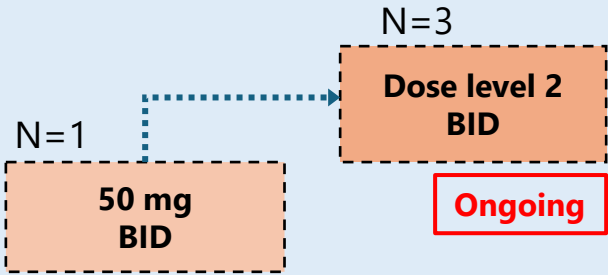
Phase 1 study



Patient Population: advanced, relapsed, refractory or distant metastasis malignant tumors

Initiated dose escalation part targeting blood cancer patients in 2024.

✓ Safety and tolerability were confirmed in a patient dosed 50 mg BID, and dosing at dose level 2 has started.

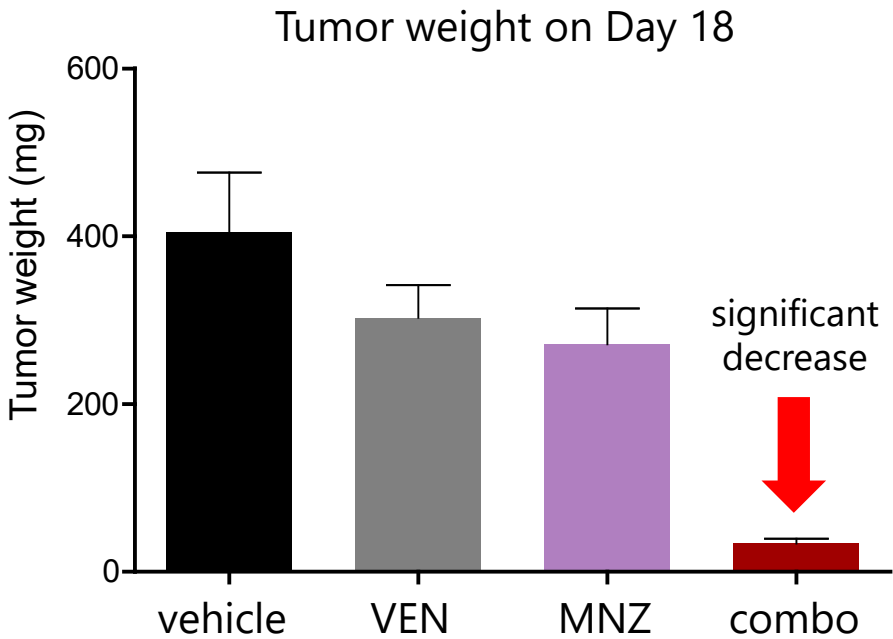




Synergistic effect of monzosertib with BCL2 inhibitor venetoclax was assessed using a MV-4-11 human AML mouse model.

AACR2024 Poster Presentation
AACR2024 : American Association for Cancer Research
Annual Meeting 2024

In vivo antitumor efficacy in MV-4-11 xenograft model mice



VEN: venetoclax (10 mg/kg, QD)
MNZ: monzosertib (60mg/kg, BID)
5 days/week, dosed orally

- In combination with venetoclax, monzosertib demonstrated robust tumor growth inhibition in a MV-4-11 human AML xenograft model.



Phase 1 dose escalation part targeting solid tumor

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

- 100 mg BID (twice a day) was well-tolerated and safe and MTD was determined to be this dose level.
- Dose expansion part was initiated at 80 mg BID , one dose level lower than MTD
- 6 of 10 efficacy-evaluable patients achieved SD
- One patient receiving 50 mg BID (every day) achieved long SD (>6 months).

Doses	Regimen	n	DLT
50 mg BID	every day	3	0
80 mg BID	every day	4	0
100 mg BID	every day	3	0

Completed dose escalation part, and dose expansion part is ongoing.

MTD: Maximum Tolerated Dose

DLT: Dose Limiting Toxicity



Updates on Licensed Pipelines

- 1 DGK α 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
DGKα inhibitor 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 α inhibitor: GS-9911

2. Indication: Cancer (immunotherapy)

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

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

In December 2024, Carna received a progress report on the development of GS-9911 from Gilead and confirmed the progress is on track.

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">• 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.**



FY2024 Results



FY2024 Consolidated Financial Results

(JPY million)	FY2023 Actual	FY2024 Actual	YoY Change	FY2024 Plan	FY2024 Dec.19 Revised Plan	
Total Sales	1,625	636	-989 -60.8%	925	630	
ddSP business	918	636	-224 -31.5%	925	630	<ul style="list-style-type: none"> Overall sales declined due to drastic decrease of overseas sales than expected including in the U.S. and China while sales of proteins and profiling services in Japan remained solid.
ddRD business	707	—	-707	—	—	
Total Operating Profit/Loss	-1,116	-2,076	-959	-2,201	-2,153	
ddSP business	225	-34	-259	229	-38	
ddRD business	-1,342	-2,041	-699	-2,431	-2,115	<ul style="list-style-type: none"> Continued investment in the clinical-stage programs.
Ordinary Loss	-1,126	-2,080	-954	-2,208	-2,168	
Net Loss	-1,152	-2,178	-1,025	-2,225	-2,180	
R&D cost	1,903	1,886	-16	2,309	1,965	<ul style="list-style-type: none"> Phase 1b study of docirbrutinib (AS-1763) is on track. Continued investment in the clinical-stage programs including costs related to clinical studies and manufacturing of investigational new drugs for docirbrutinib (AS-1763) and monzosertib (AS-0141).

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

ddRD: Drug Discovery R&D business ddSP: Drug Discovery Support business

Note : Rounded down to the nearest million yen



Consolidated Balance Sheet

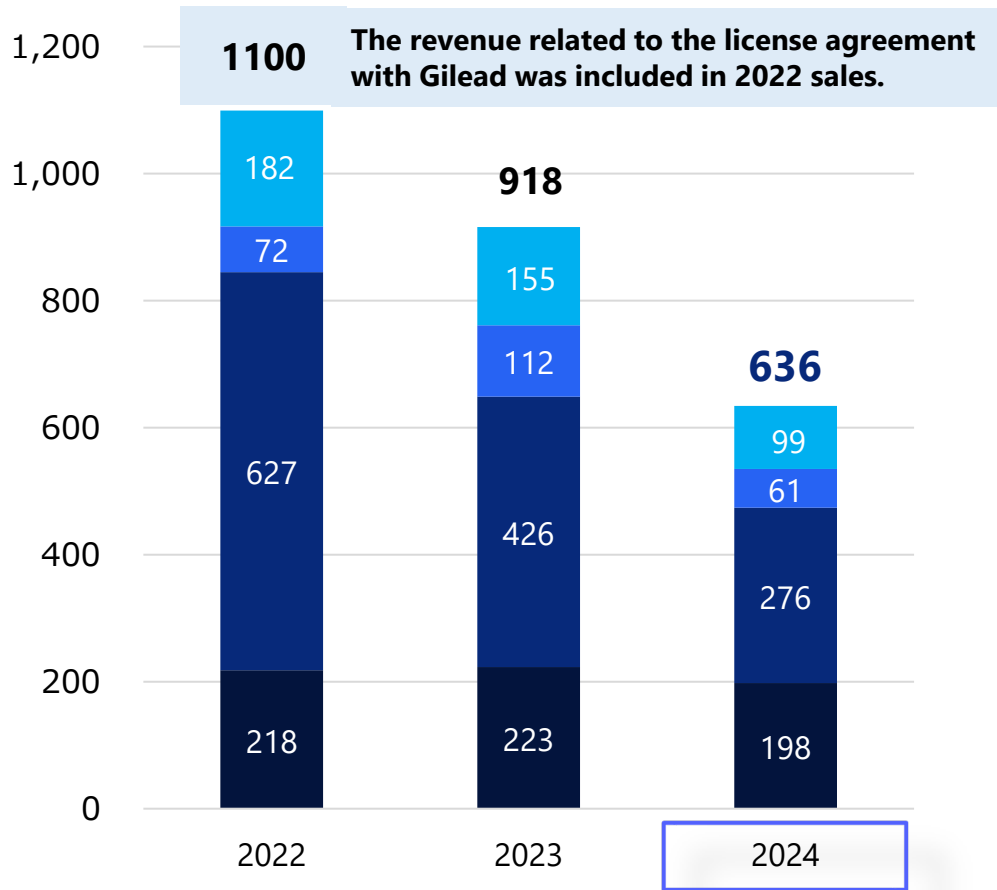
(JPN million)	As of Dec. 31,2023	As of Dec. 31,2024	Change	Reason for changes
Current assets	4,191	2,737	-1,453	<ul style="list-style-type: none">Cash and deposits -780Accounts receivable-trade -701
Cash and deposits	2,889	2,108	-780	
Non-current Assets	158	34	-124	
Total assets	4,349	2,772	-1,577	
Current liabilities	375	222	-152	<ul style="list-style-type: none">Long-term loans payable within 1 year -100Accounts payable -37
Non-current liabilities	96	73	-22	
Total liabilities	472	296	-175	
Total net assets	3,877	2,475	-1,402	<ul style="list-style-type: none">Capital stock and capital surplus +371Retained earnings -2,178
Total liabilities and net assets	4,349	2,772	-1,577	
Shareholders' equity ratio	89.1%	89.2%		
BPS	226.16yen	129.62yen		
PBR	2.3x	2.3x		
Share price of Carna	522yen	300yen		

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



Drug Discovery Support Business Sales Trend by Region (Consolidated)

(JPY million) Other Europe North America Japan



Japan

Decreased 10.9% YoY

- Sales of proteins to pharmaceutical companies and distributors were robust.
- Sales of profiling services remained solid while sales of cell-based assay services declined.

North America

Decreased 35.1% 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 45.7% YoY

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

Other

Decreased 35.9% 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 the start of production of kinase protein by some Chinese CROs also caused the continuous reduction in demand in China.



Raised additional capital through the 3rd party allotment of common shares in May and in October.

● Allotee Athos Asia Event Driven Master Fund

Settlement date	May 31 st , 2024	October 11 th , 2024
Net proceeds (JPN million)	362	378
Use of proceeds	Investment in development of docirbrutinib (AS-1763) and monzosertib (AS-0141)	

Financing

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



Business Plan



Key Milestones for 2025

Business		Key Milestones		
		Milestones for 2024	Achievement in 2024	Milestones for 2025
ddRD	Docirbrutinib (AS-1763)	<input type="checkbox"/> Present interim clinical data from ongoing Ph1b study	<input checked="" type="checkbox"/> Present interim clinical data from ongoing Ph1b study • Presented the preliminary data of Ph1b study at EHA2024 in June 2024 and ASH in December 2024	<input type="checkbox"/> Promote Ph1b dose expansion part and present interim clinical data from the study
	Sofnibrutinib (AS-0871)	<input type="checkbox"/> Promote partnering activity to find a strategic partner	<input checked="" type="checkbox"/> Promote partnering activity to find a strategic partner	<input type="checkbox"/> Out-license or initiate Ph2 by joint-development
	Monzosertib (AS-0141)	<input type="checkbox"/> Enroll patients with blood cancers <input type="checkbox"/> Initiate Ph1 dose expansion part	<input checked="" type="checkbox"/> Enroll patients with blood cancers <input type="checkbox"/> Initiate Ph1 dose expansion part Solid tumor : Initiated patient recruitment in January 2025	<input type="checkbox"/> Promote Ph1 study and choose cancer types for the clinical development
ddSP		<input type="checkbox"/> Expand sales of in-house developed products and services in North America and Asia <input type="checkbox"/> Increase line-up of protein kinase products <input type="checkbox"/> Expand sales of cell-based assay	<input type="checkbox"/> Expand sales of in-house developed products and services in North America and Asia <input checked="" type="checkbox"/> Increase line-up of protein kinase products <input type="checkbox"/> Expand sales of cell-based assay	<input type="checkbox"/> Expand sales of in-house developed products and services in North America, Europe and Asia <input type="checkbox"/> Increase line-up of protein kinase products <input type="checkbox"/> Expand sales of cell-based assay

EHA : European Hematology Association
 ASH : American Society of Hematology
 ddRD: Drug Discovery R&D business
 ddSP: Drug Discovery Support Business

☒ Achieved
☐ Plan or to be achieved



(JPY million)	FY2024 Actual	FY2025 Plan	Outlook for 2026 – 2029
Total Sales	636	722	
ddSP business	636	722	Maintain stable sales
ddRD business	—	—	Revenue from milestone payments and upfront payments
Total Operating Loss	(2,076)	(2,133)	
ddSP business	(34)	83	Maintain stable profit while investing in product developments
ddRD business	(2,041)	(2,216)	Continue to invest in R&D and deliver profits depending on the size of milestone payments and upfront payments
Ordinary Loss	(2,080)	(2,137)	
Net Loss	(2,178)	(2,147)	

(JPY million)	FY2024 Actual	FY2025 Plan	Outlook for 2026 – 2029
R&D Cost	1,886	2,059	Continue to invest in R&D for the future growth. The R&D cost may vary from JPY 1 bn to 2.5 bn, depending on the size of clinical studies.
Capex	13	7	Invest in equipment for R&D and IT system (JPY20 mn to 100 mn)

- Business plan for FY2025 dose not include potential milestone payments or upfront payments as the timing or the amounts are difficult to predict.
- Numerical targets for 2026-2029 are not disclosed for the same reason.

ddRD: Drug Discovery R&D business
ddSP: Drug Discovery Support Business



➤ Advance clinical trials of our innovative pipelines to maximize corporate value

	Started internal drug discovery activity	Demonstrated strong capabilities in drug discovery	Maximize the value of pipelines	Continue delivering profits
	2010-2015	2016-2020	2021-2025 (Plan)	2026-2030 (Plan)
PLAN	<ul style="list-style-type: none">Established in-house research capabilityEstablished pipeline	<ul style="list-style-type: none">Out-licensed multiple programsInitiated clinical trials	<ul style="list-style-type: none">Advance clinical trials of docirbrutinib (AS-1763), sofnobrutinib (AS-0871) and monzosertib (AS-0141)Earn revenue from new license dealsReceive milestone payments from the out-licensed programs and deliver profitsInitiate pre-clinical and clinical studies of new pipelines	<ul style="list-style-type: none">Receive milestone payments and royalty income from the out-licensed programs and expand profitsEarn revenue from new license dealsInitiate pre-clinical and clinical studies of new pipelines
ACTUAL	<ul style="list-style-type: none">✓ Out-licensed : J&J (2015)	<ul style="list-style-type: none">✓ Out-licensed : Sierra Oncology (2016)✓ Joint research : Sumitomo Pharma (2018)✓ Out-licensed : Gilead (2019)✓ Out-licensed : BioNova (2020)✓ Initiated clinical trials of sofnobrutinib (2020)	<ul style="list-style-type: none">✓ Out-licensed : Fresh Tracks Therapeutics (2022)✓ 1st Milestone : BioNova (2022)✓ 1st Milestone : Gilead (2021)✓ 2nd Milestone : Gilead (2023)✓ Initiated clinical trials of docirbrutinib (2021)✓ Initiated clinical trials of monzosertib (2021)	

<ddRD>

- ✓ Advance clinical trials of docirbrutinib (AS-1763) and monzosertib(AS-0141)
- ✓ Create next wave of pipeline
- ✓ Receive milestone payments and royalty income from out-licensed programs

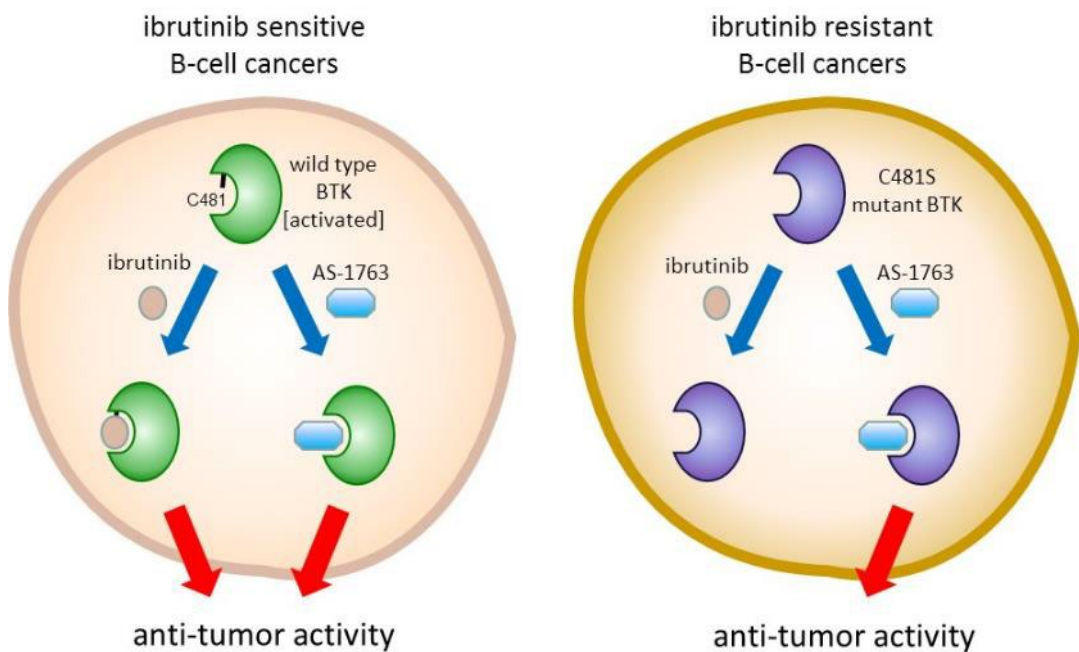
<ddSP>

- Expand sales of in-house developed products and services in North America and Asia
- Secure sustainable sales growth by launching new products and services and reaching out to new customers
- Generate cash to invest in ddRD

ddRD: Drug Discovery R&D business
ddSP: Drug Discovery Support Business



Appendix 1



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 docirbrutinib against wild-type and C481S-mutant BTK

	IC ₅₀ (nM)	
	BTK[A]	BTK ^{C481S}
docirbrutinib	0.85	0.99

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

◆ In vitro pharmacological activities of docirbrutinib

	IC ₅₀ (nM)	
	docirbrutinib	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

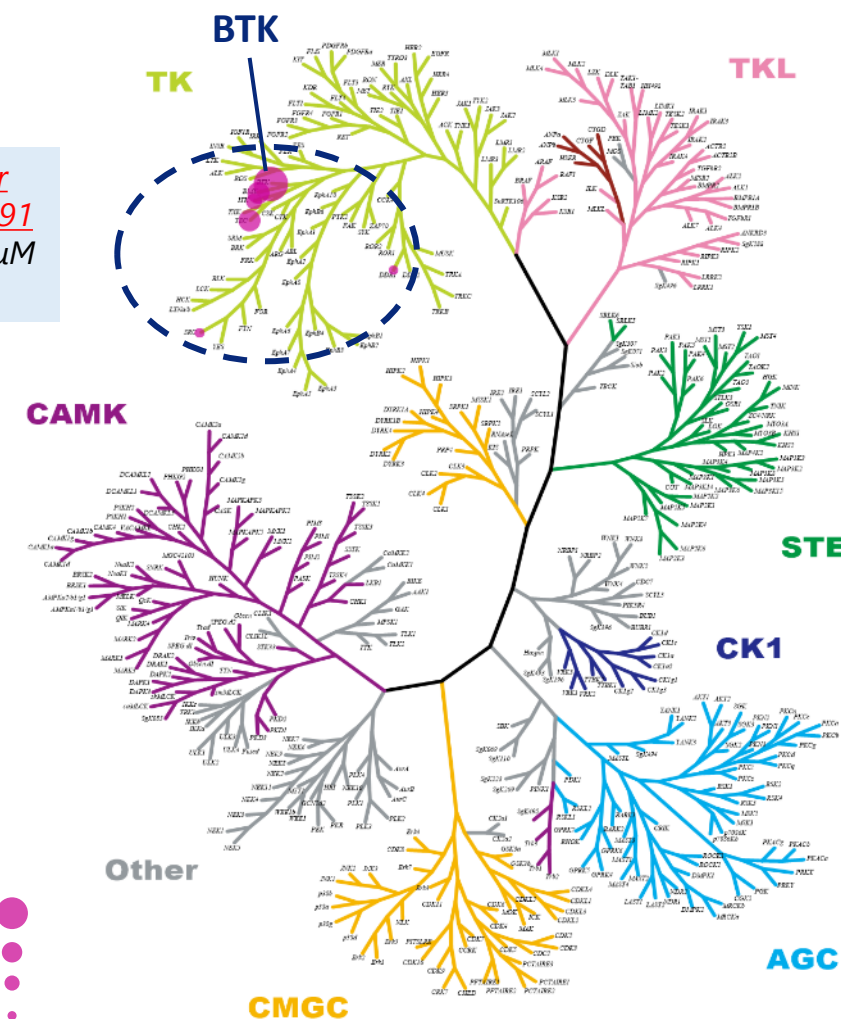
50-fold Stronger activity

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.

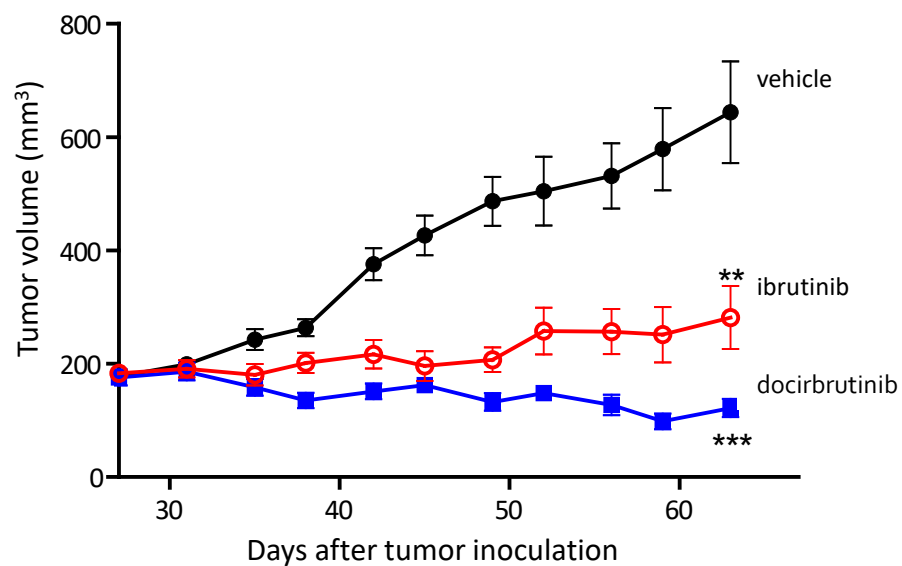
◆ Kinase selectivity profiling

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





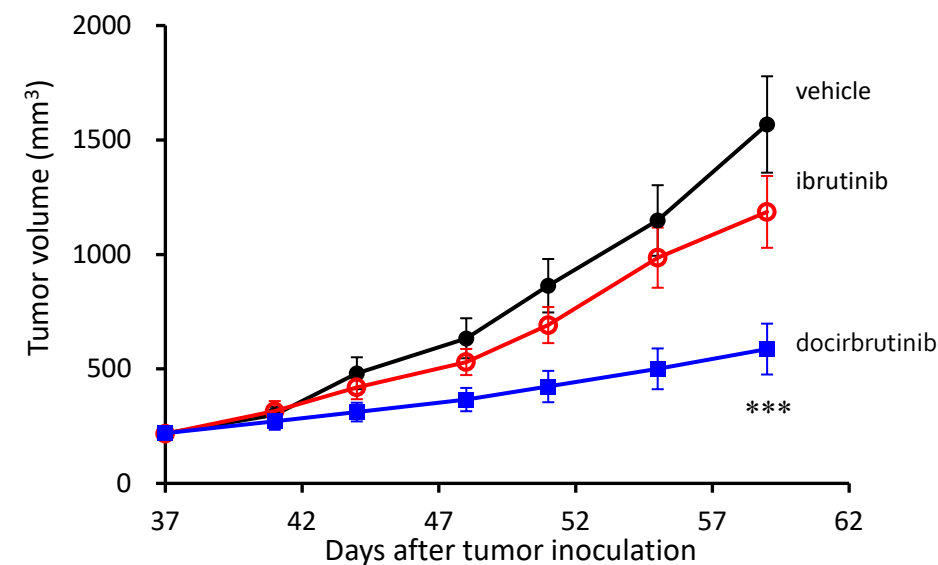
- ◆ In vivo antitumor effects of docirbrutinib on human B-cell non-Hodgkin lymphoma cell line, OCI-LY10 tumor xenograft mouse model (n=8-10)



Ibrutinib: 25 mg/kg QD
Docirbrutinib : 60 mg/kg BID

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

- ◆ In vivo antitumor effects of docirbrutinib on ibrutinib-resistant BTK^{C481S} knock-in OCI-LY10 tumor xenograft mouse model (n=11)



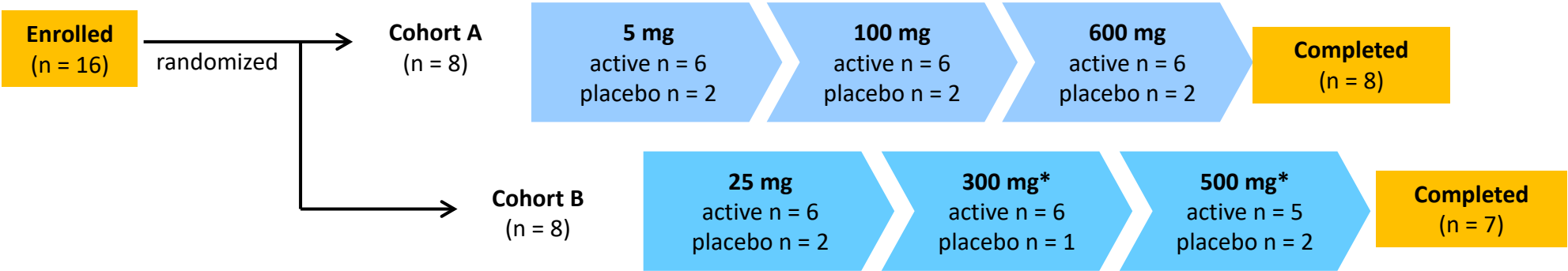
Ibrutinib: 25 mg/kg QD
Docirbrutinib : 60 mg/kg BID

***: 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 docirbrutinib 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.



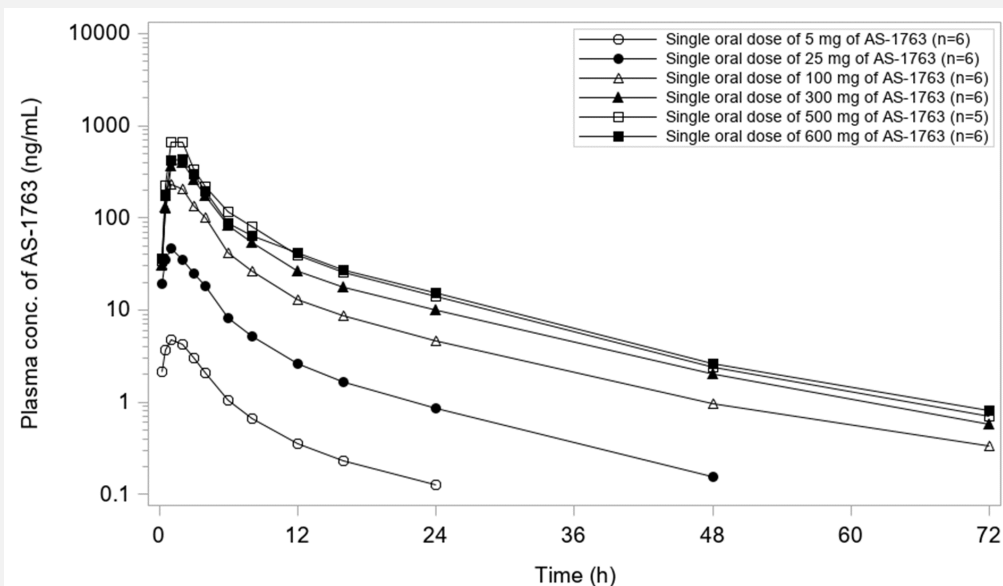
Safety and tolerability

- Docirbrutinib 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 docirbrutinib 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 docirbrutinib exposures generally increased with dose up to 500 mg.

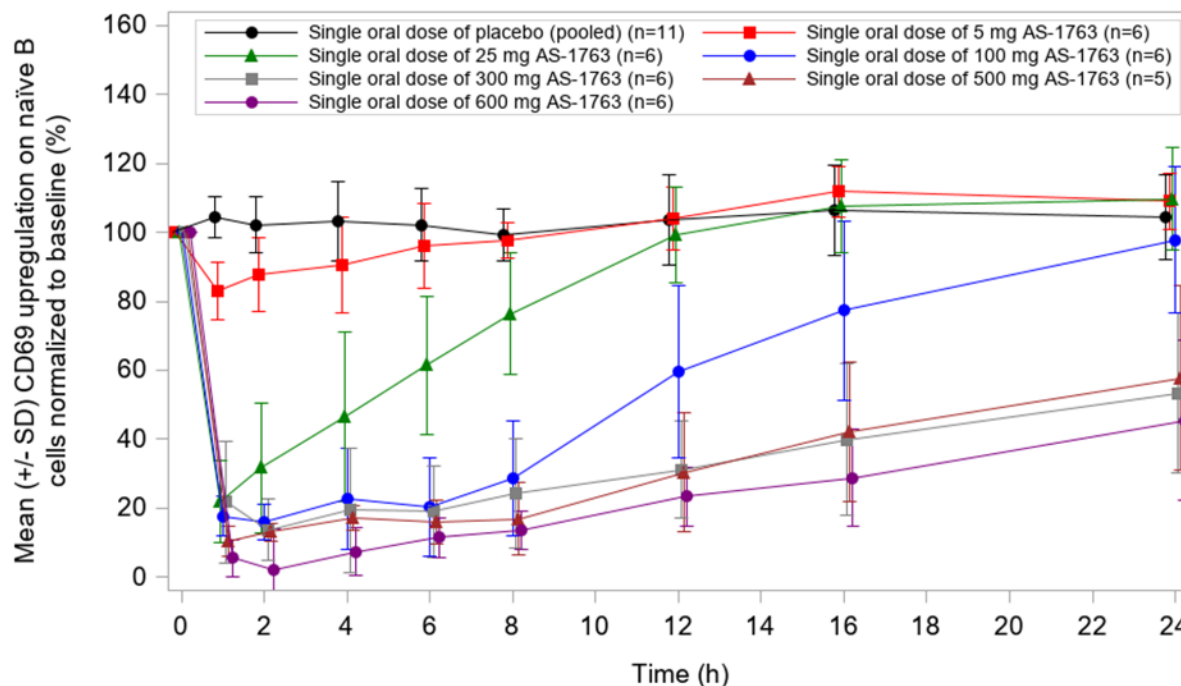
< Plasma concentration of a single oral dose of docirbrutinib >



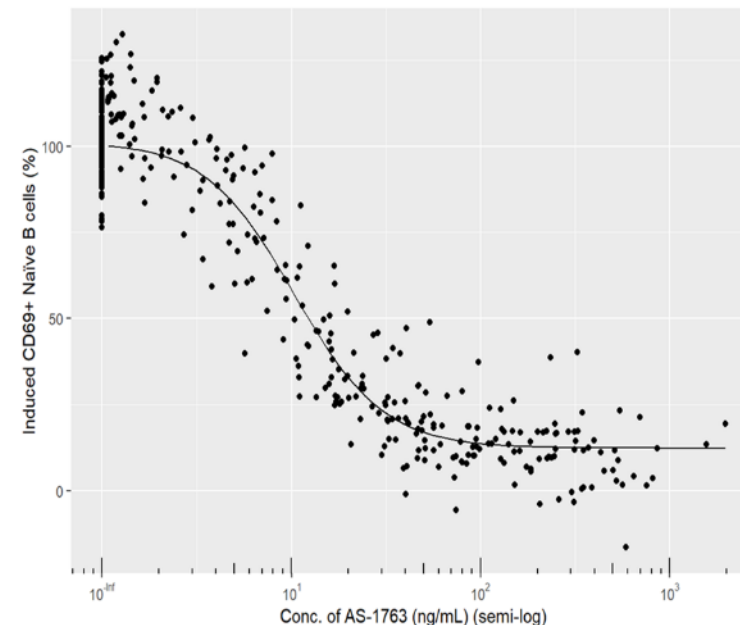


- Inhibition of B cell CD69 upregulation was observed for 5 mg onwards.
- Maximum inhibition (arbitrarily defined as $\geq 80\%$) was observed at 1-2 hours post-dose from 100 mg to 600 mg, and the duration of inhibition was dose-dependent with values of 2, 6, 8 and 8 hours for 100, 300, 500, and 600 mg, respectively.
- Based on a PK/PD correlation analysis, the IC₅₀ value of docirbrutinib on CD69 upregulation was calculated to be 10.5 ng/mL.

< B cell CD69 upregulation after a single oral dose of docirbrutinib >



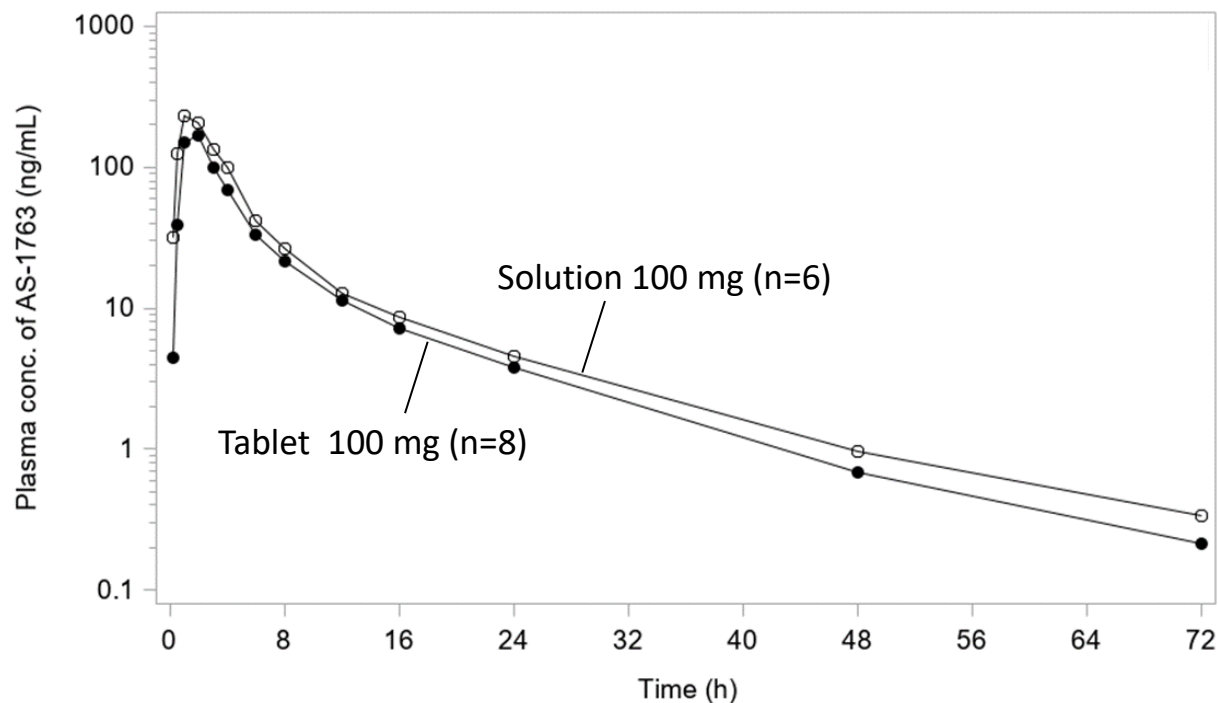
< PK/PD correlation analysis >





- 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 docirbrutinib tablet twice daily dosing in relapsed/refractory CLL and B-cell NHL.

<PK of Tablet vs Solution after a Single oral dose docirbrutinib>





Objectives of the study

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

- plasma concentration of docirbrutinib
- 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 docirbrutinib :

- plasma concentrations of docirbrutinib 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 docirbrutinib to healthy volunteers.

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

◆ Targeting Inactive Conformation of BTK



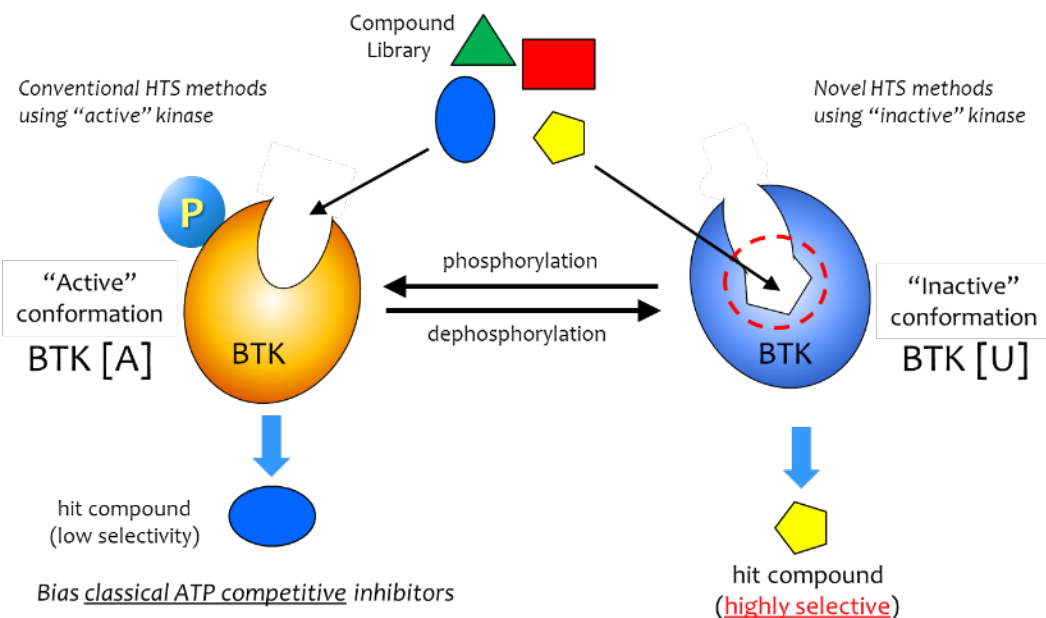
TR-FRET binding assay targeting unactivated form of Bruton's tyrosine kinase

Tokiko Asami¹, Wataru Kawahata, Masaaki Sawa
Carna Biosciences, Inc., 4-1-15 Minamigoshi, Minamikuwa, 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

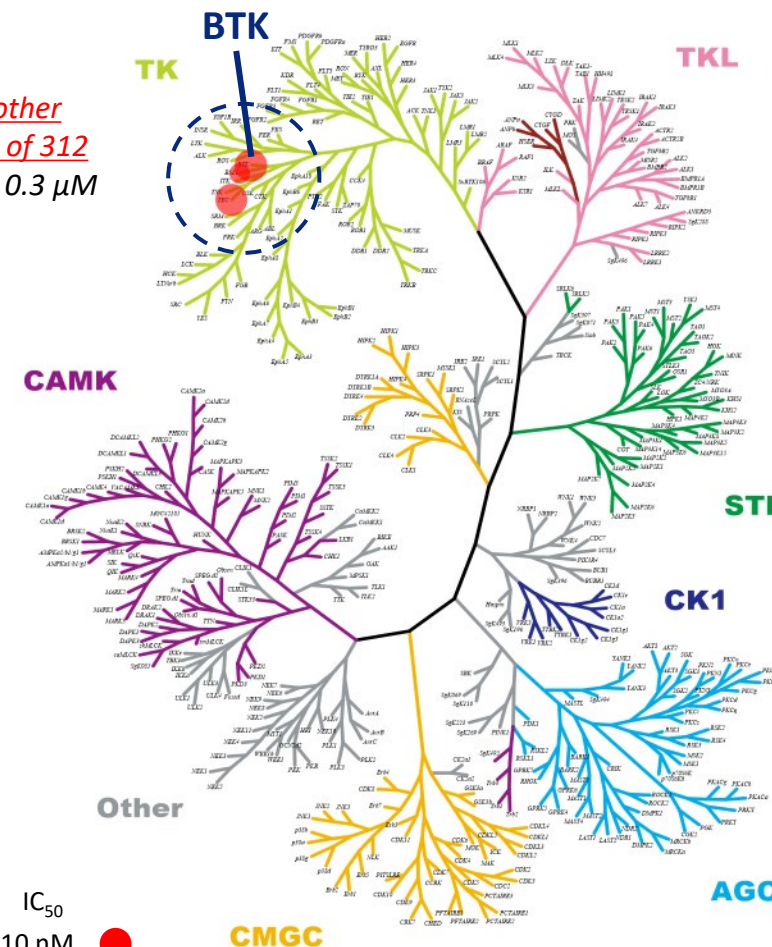
Wataru Kawahata¹, Tokiko Asami, Takao Kiyoi, Takayuki Irie, Haruka Taniguchi, Yuko Asanishi, Tomoe Inoue, Takahiro Miyake, and Masaaki Sawa¹
Research and Development, Carna Biosciences, Inc., 4-1-15 Minamigoshi, Minamikuwa, Chuo-ku, Kobe 650-0047, Japan



	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.

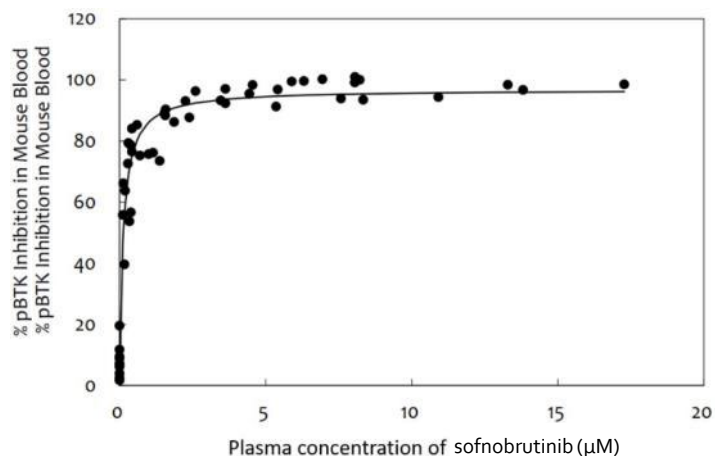
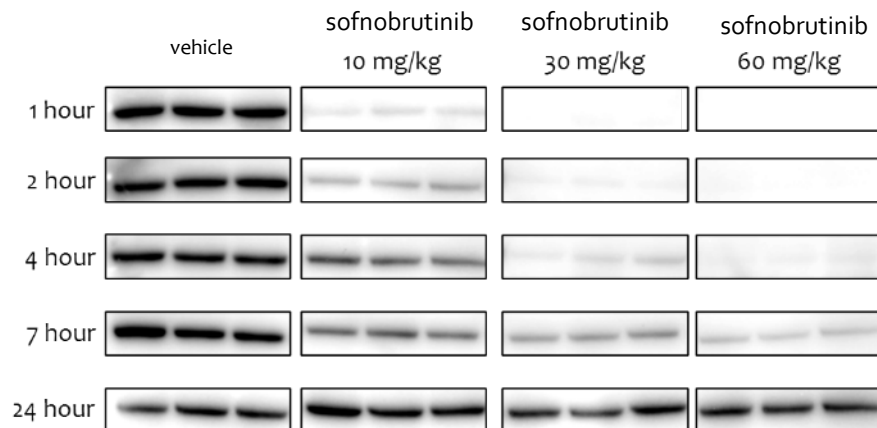


IC₅₀
< 10 nM
10 - 100 nM
100 - 1000 nM

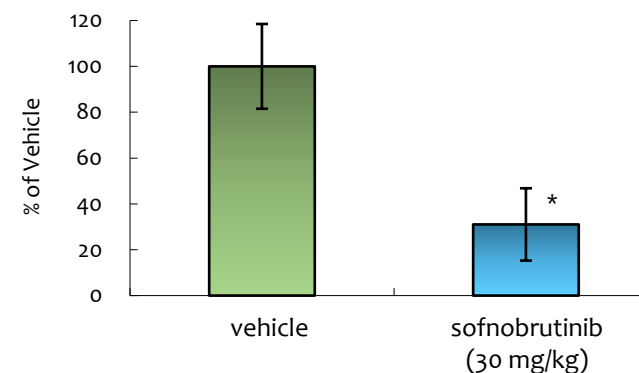


◆ PK/PD Study

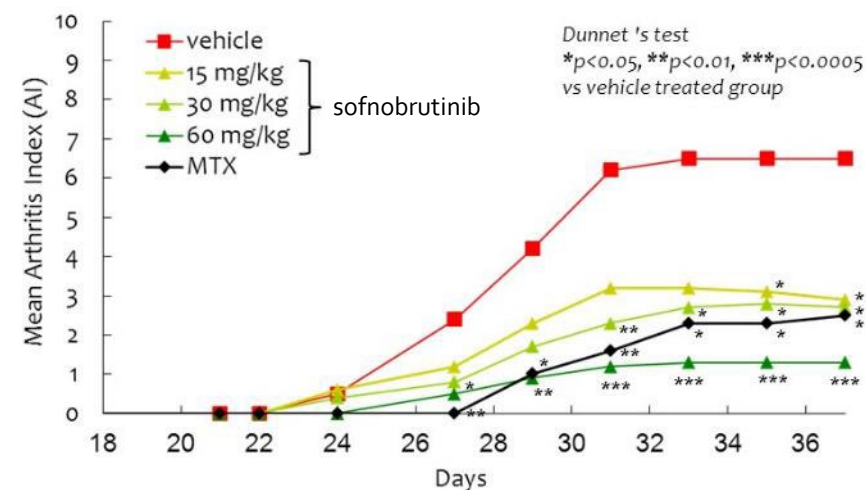
Auto-phosphorylation status of BTK was measured following oral single administration of sofno Brutinib.



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



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





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

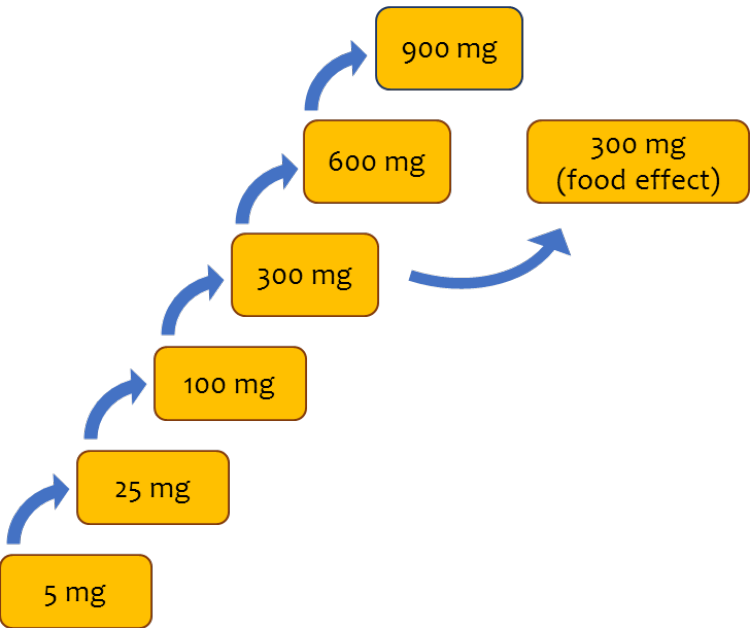
Evaluate the safety, tolerability, PK and PD in the 2-week multiple ascending dose of sofnobrutinib.

Completed



SAD Part (Completed)

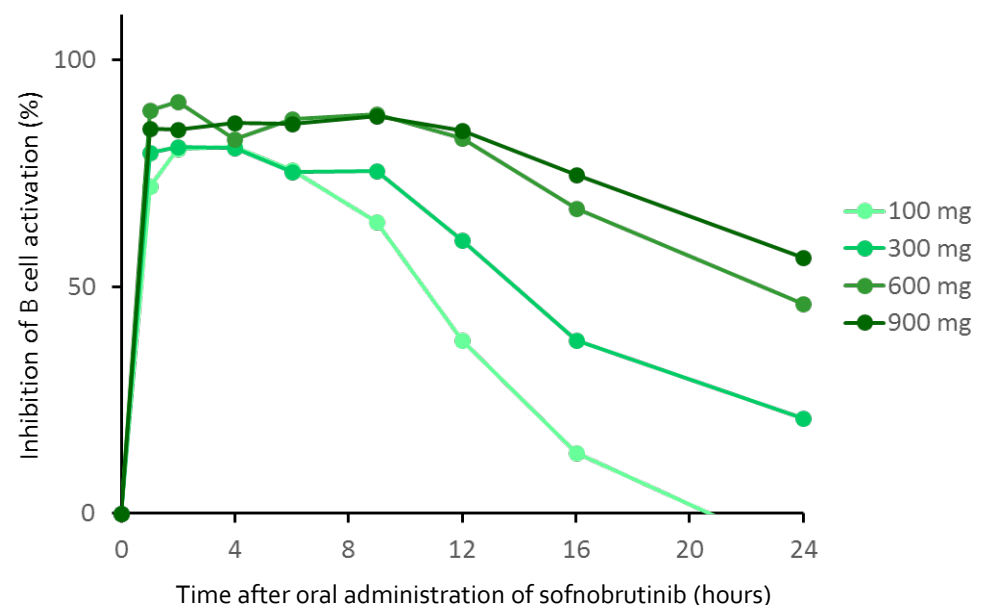
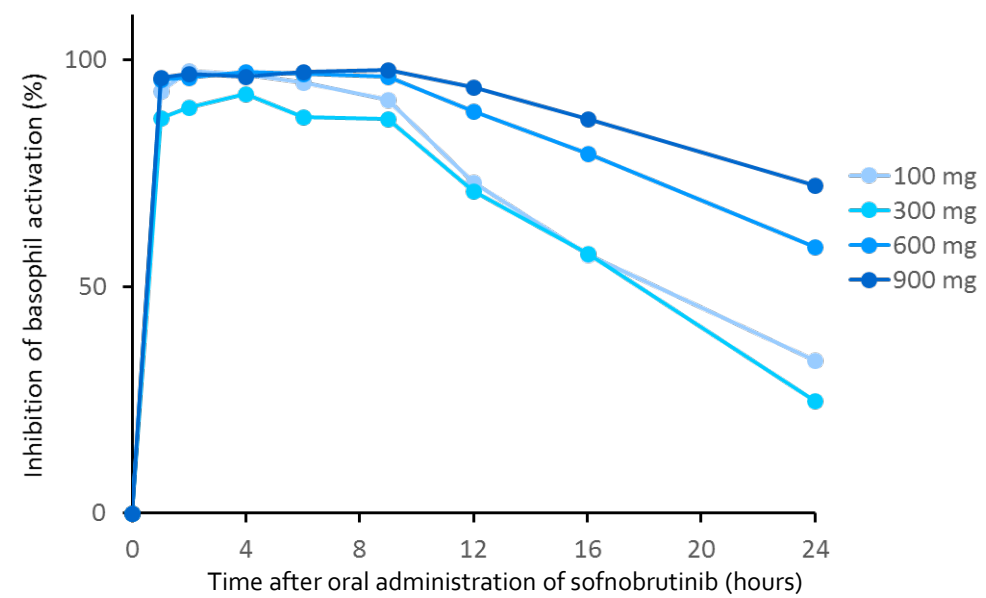
Step 1 Single Ascending Dose (SAD)	Step 2
<ul style="list-style-type: none">• 6 dose levels (8 subjects/cohort)• Placebo controlled (6 active / 2 placebo)• Safety and tolerability• Pharmacokinetics and pharmacodynamics	<ul style="list-style-type: none">• Food effect



- 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.



- 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.

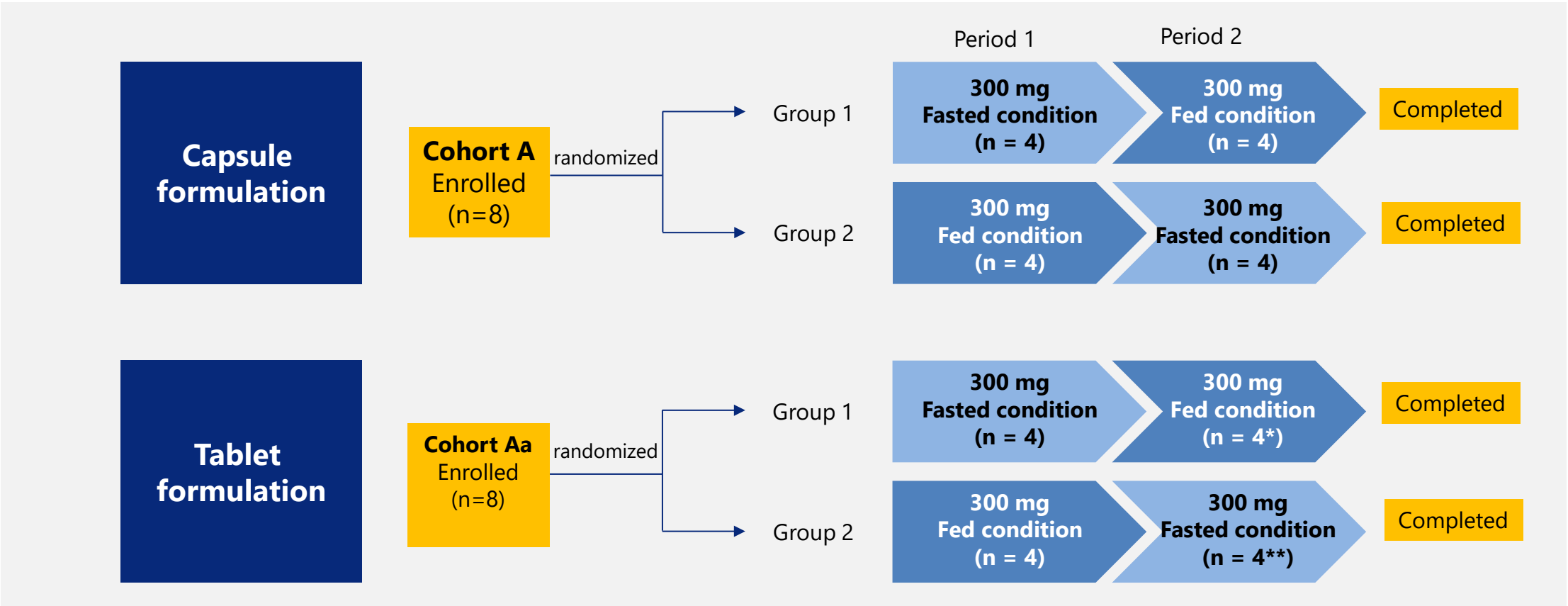




Sofnobrutinib (AS-0871) : Phase 1 MAD Study BA part

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).



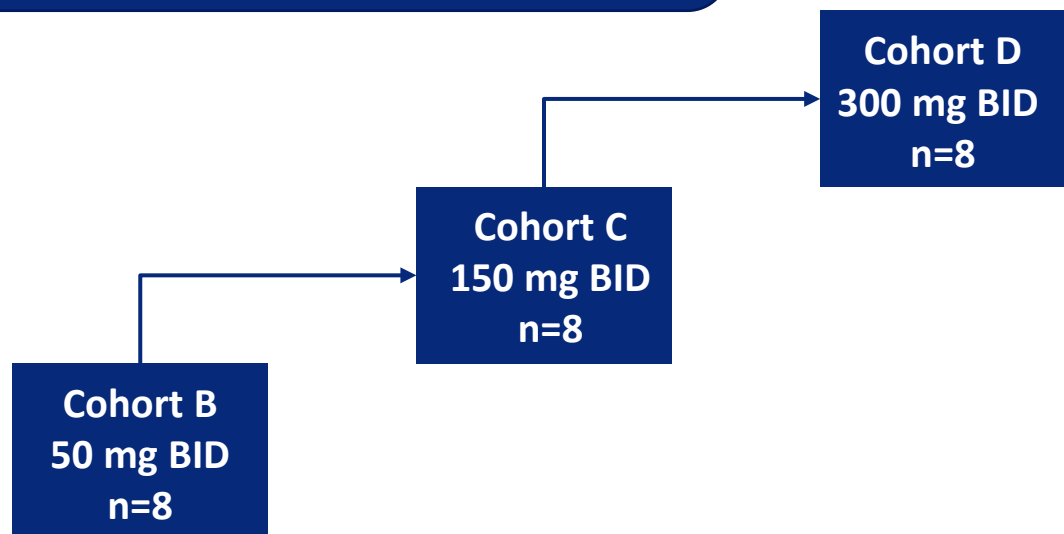
*One subject vomited after dosing (considered not related to study drug), excluded from the PK analysis.
**One subject withdrew from the study due to personal reasons before dosing.



Study Design of MAD part

- In the MAD part, safety, tolerability, PK, and PD of 3 multiple ascending doses of sofnobrutinib, following 14-day multiple dose oral administration of sofnobrutinib, 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



Objectives of Single Ascending Dose (SAD) study

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 sofno Brutinib 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.

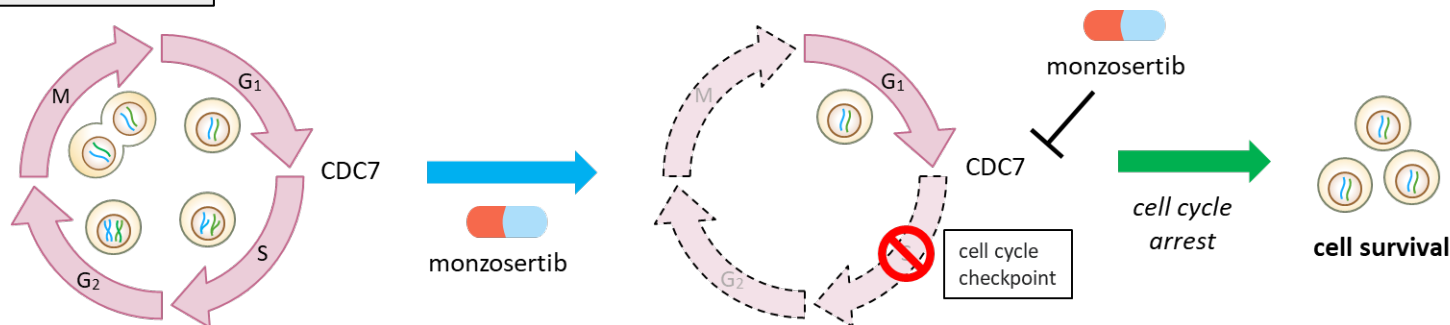
These results from the Phase 1 studies of sofno Brutinib support to advance sofno Brutinib into Phase 2 clinical development for further investigations.



CDC7 Kinase Inhibitor: MoA of monzosertib

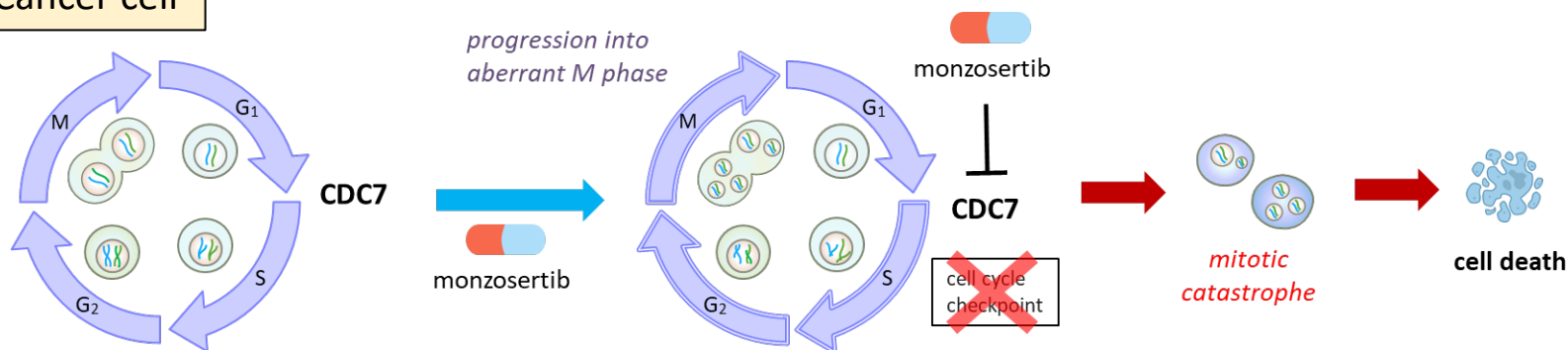
- ◆ CDC7 (cell division cycle 7) is a serine/threonine kinase that facilitates DNA replication during DNA synthesis
- ◆ Over expression of CDC7 has been reported to cause uncontrolled proliferation of many cancer types

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

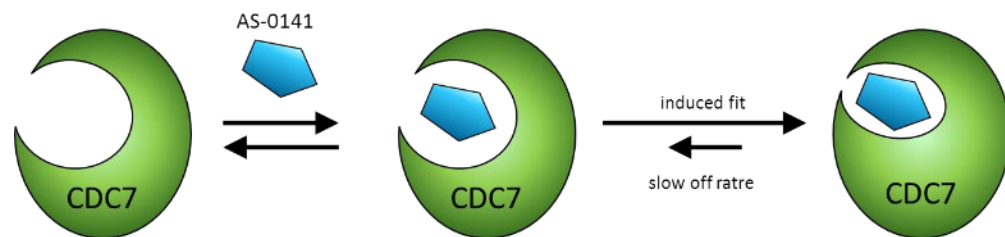


Inhibition of CDC7 in cancer cells causes lethal S phase or M phase progression when the cancer cells have a defective cell cycle checkpoint

◆ 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^{a,*}, Tokiko Asami^a, Ayako Sawa^a, Yuko Uno^a, Mitsuhiro Hanada^a, Chika Taniyama^a, Yoko Funakoshi^a, Hisao Masai^b, Masaki Sawa^c
^a Research and Development, Carina Biosciences, Inc., 2F Bldg., 1-5-2 Minamigino Minamigino-ku, Kyoto 604-0047, Japan
^b Research and Development Department, Otsuka Pharmaceutical Co., Ltd., 18-18-1 Hongo, Bunkyo-ku, Tokyo 112-8604, Japan
^c Department of Cellular Medicine, Tokyo Metropolitan Institute of Medical Sciences, 2-1-6, Minamishinjyuku, Shinjyuku-ku, Tokyo 162-8602, Japan



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

Without
Preincubation

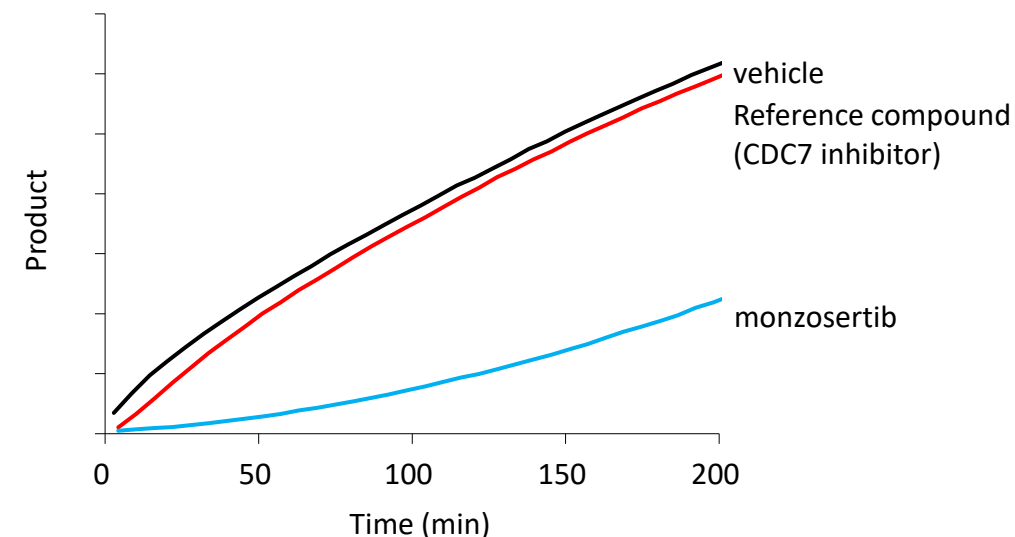
503 nM

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.

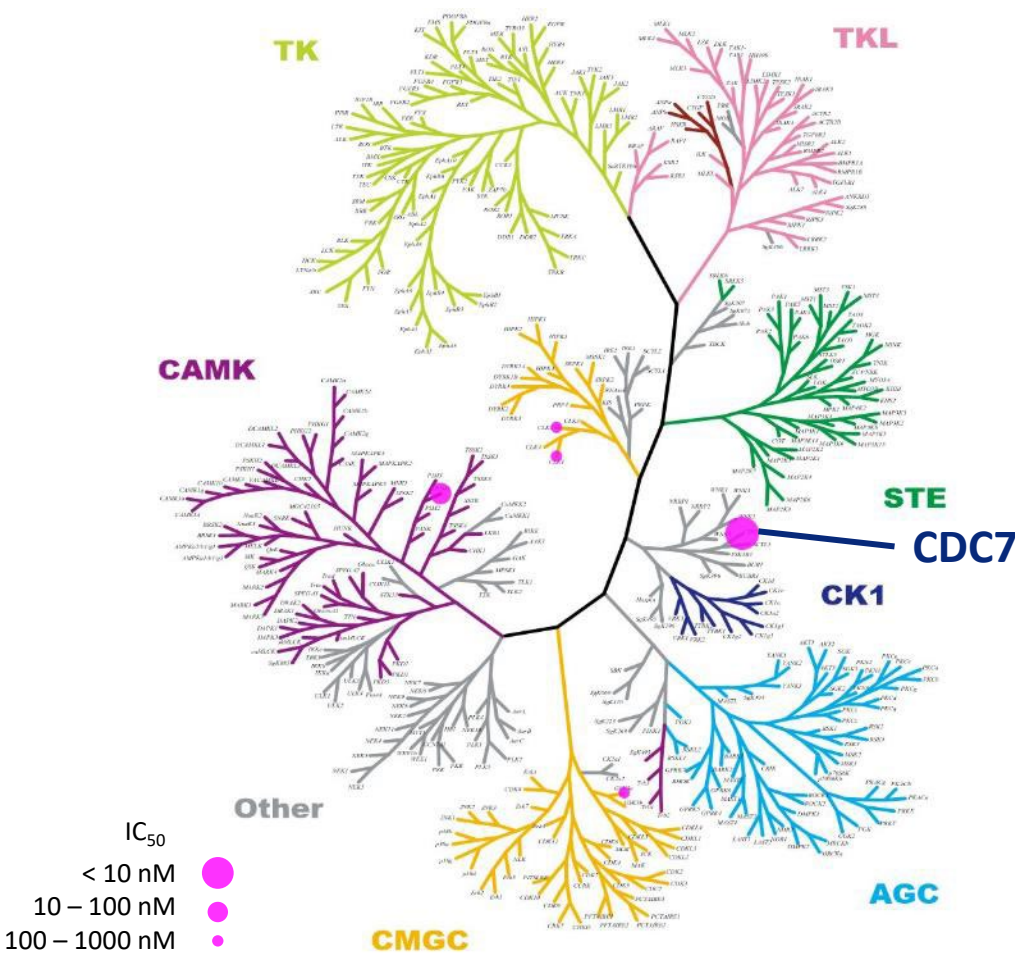




Monzosertib (AS-0141) : High Kinase Selectivity

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

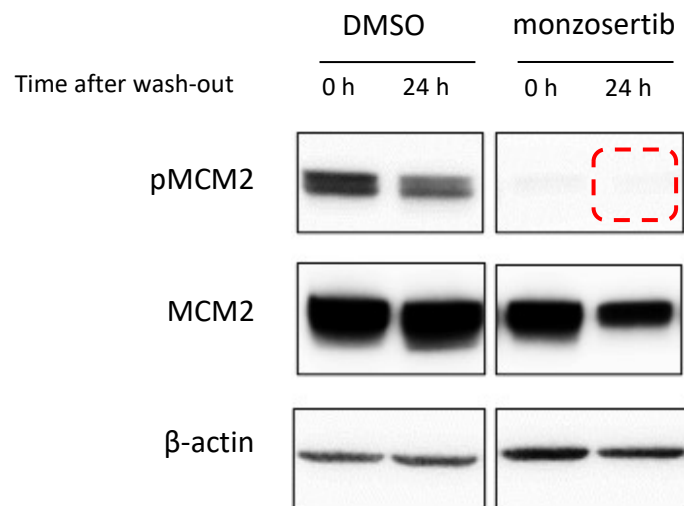


Monzosertib (AS-0141) : Strong Cellular Activity



◆ Prolonged inhibition in cells

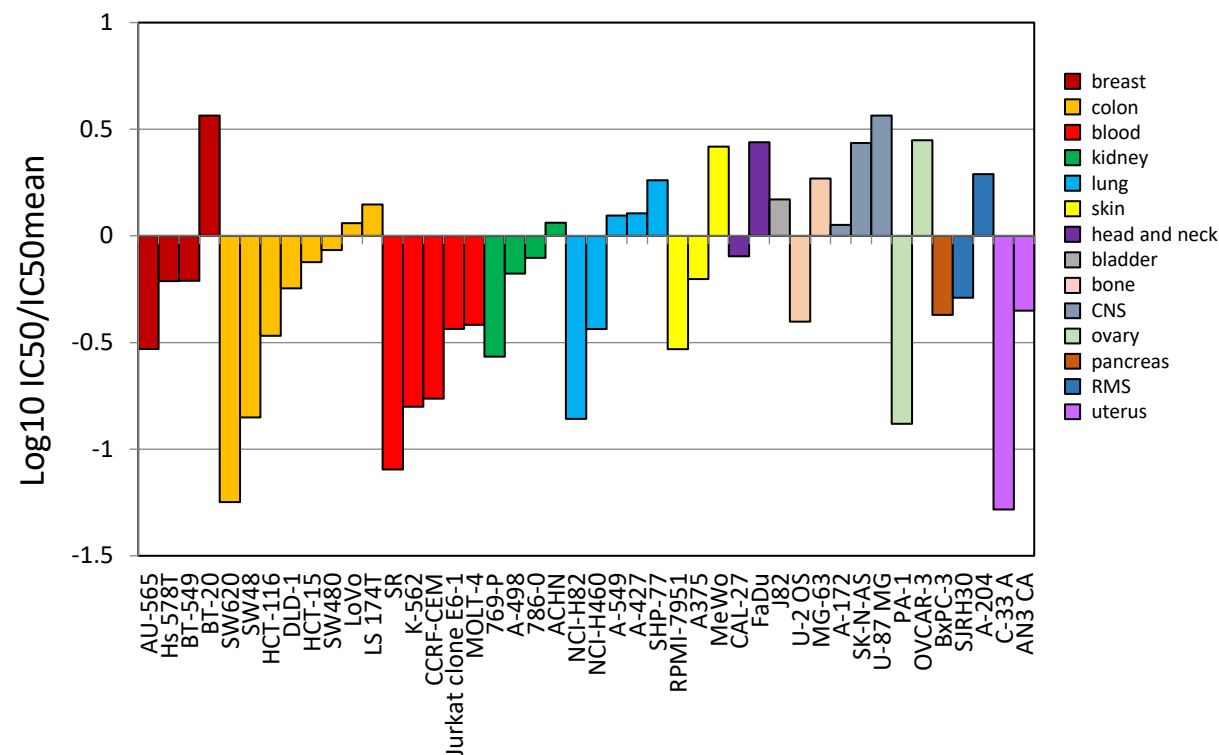
Human colon cancer cell line, Colo-205 cells were treated with DMSO control or monzosertib. After washout of the inhibitor, the cells were further incubated in the same media for 0 or 24 h and subjected to western blot analysis.



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

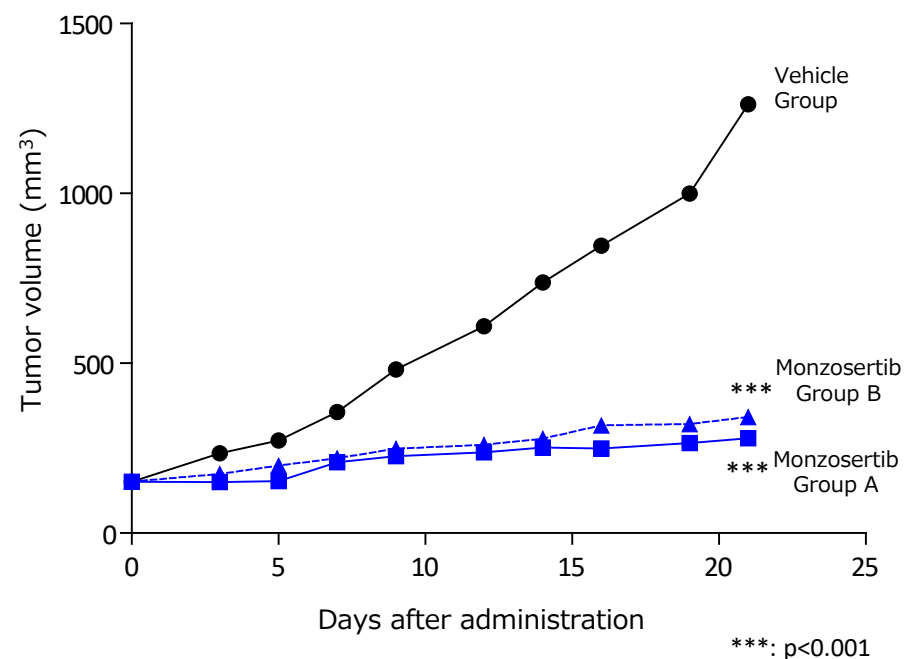


44 Cancer cell lines (Oncolines at NTRC)

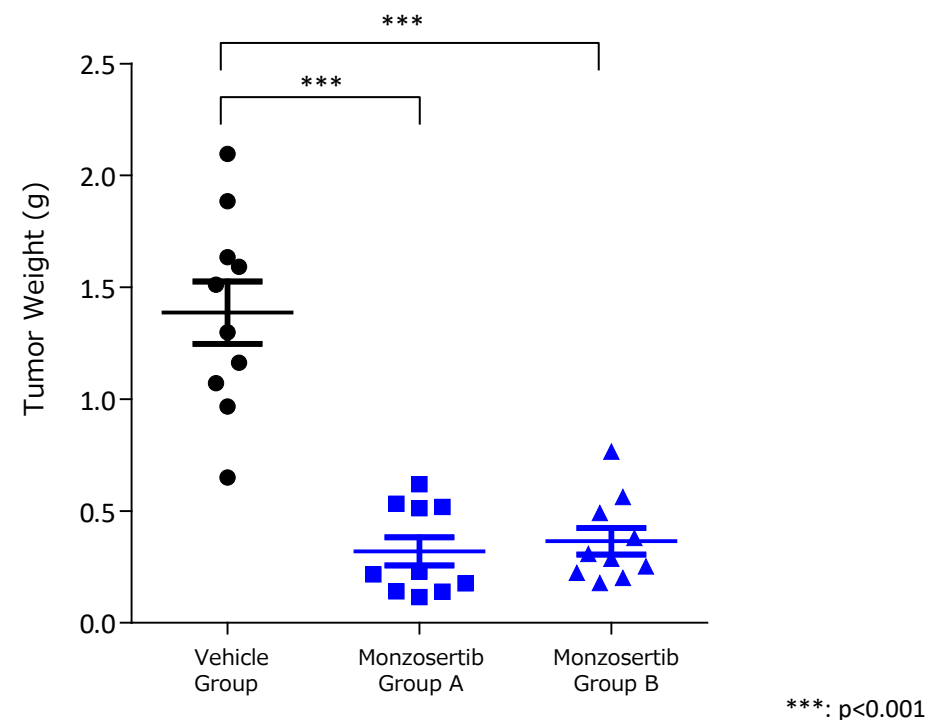


In vivo antitumor efficacy of monzosertib in a SW620 (human colon cancer) xenograft mouse model

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



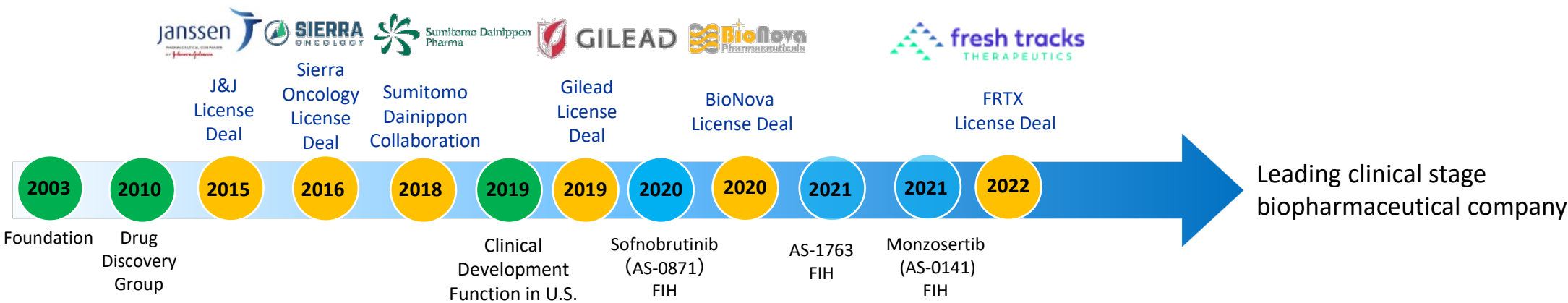
Appendix 2



Building Long-Term Value



Our goal is to deliver innovative therapies for patients suffering from serious diseases



2003 – 2024	2025 Plan	Long term plan
<ul style="list-style-type: none">● Founding members who had expertise in kinase drug discovery technology spun out from Nippon Organon and established Carna.● Started offering kinase proteins and screening services to pharmaceutical companies for kinase inhibitor drug discovery.● In 2010, Drug Discovery Group was established to conduct internal drug discovery.● Entered into five license agreements and one joint-development agreement with pharmaceutical companies.● Initiated FIH study of BTK inhibitor docirbrutinib (AS-1763), sofnobrutinib (AS-0871) and CDC7 inhibitor monzosertib(AS-0141).● Completed Phase 1 study of BTK inhibitor sofnobrutinib(AS-0871) and started partnering activities.	<ul style="list-style-type: none">● Actively seek a strategic partner to bring sofnobrutinib into late clinical development stages.● Advance Phase 1 studies of BTK inhibitor docirbrutinib (AS-1763) and CDC7 inhibitor monzosertib (AS-0141).● Strengthen global clinical development capability.● Create next wave of pipeline.	<ul style="list-style-type: none">● Advance clinical studies of docirbrutinib (AS-1763) and monzosertib (AS-0141).● Find strategic partners for late stage development and commercialization.● Receive milestone payments and royalties from licensees and strengthen financial position.● Create next wave of pipeline.

FIH: First in Human



“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.