

Financial Results

FY2021 Q3

(January to September 2021)












Carna Biosciences, Inc.



Stock Code: 4572








- Announced positive results for AS-0871 Phase 1 Single Ascending Dose study. (July)
- Completed dosing in Phase 1 Single Ascending Dose study of BTK inhibitor AS-1763. (July)
- Bought back and canceled Series 18th Subscription Rights to Shares. (July)
- Issued Series 19th Subscription Rights Shares. (July)
- The board resolved to apply for the listing on the Growth Market under the new market segment of Tokyo Stock Exchange.(September)
- The research paper on the discovery of AS-1763 was published in *Journal of Medicinal Chemistry*. (September)
- The research paper on the discovery of AS-0141 was published in *Journal of Medicinal Chemistry*. (October)

<Oncology>

| Compound | Target | Indication | Discovery/Preclinical | Clinical | Partner |
|----------------|----------|-----------------|---|---|---|
| AS-0141 | CDC7/ASK | Cancer |   |  | |
| Small Molecule | Kinase | Immuno-Oncology |  | |  |
| AS-1763 | BTK | Blood Cancer |   |  |  |
| Small Molecule | ALK5 | Immuno-Oncology |  | | |
| Small Molecule | CDK1 | Cancer |  | | |

*Greater China only

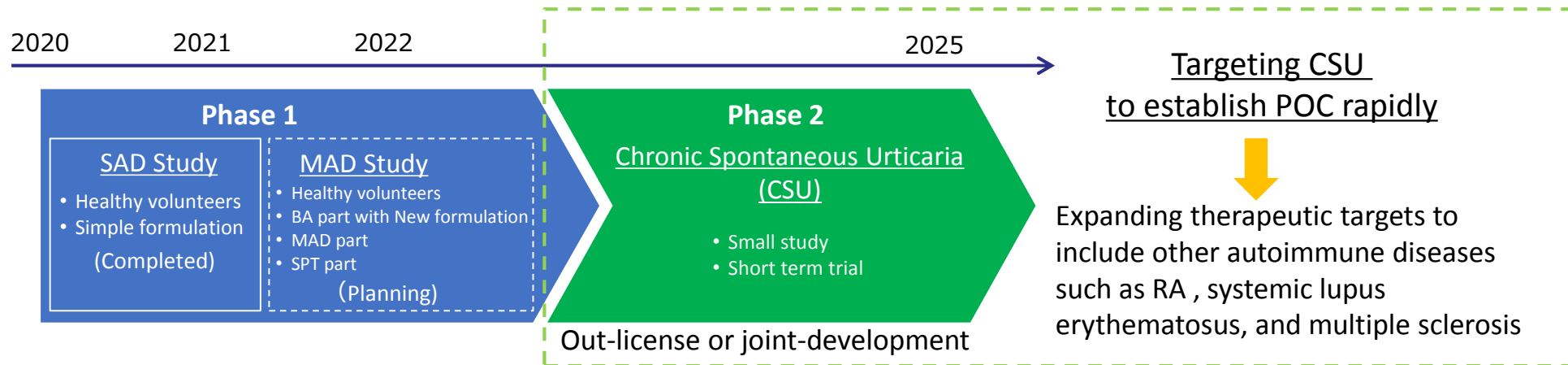
<Other Therapeutic Areas>

| Compound | Target | Indication | Discovery/Preclinical | Clinical | Partner |
|----------------|--------|------------------------------|---|---|---|
| Small Molecule | Kinase | Psychiatry & neurology |  | |  |
| AS-0871 | BTK | Immune-inflammatory diseases |   |  | |
| Small Molecule | N/A | Malaria |  | | |
| Small Molecule | STING | Immune-inflammatory diseases |  | | |

✓ We are actively pursuing early discovery programs to create next wave of pipeline.

AS-0871 : 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
- Find a partner to conduct further development after completing Phase 1 study



■ Phase 1 Single Ascending Dose (SAD) study: *Completed*

- ✓ The SAD study in healthy volunteers was initiated in H2 2020 in the Netherlands
- ✓ Safe and well-tolerated at all dose levels tested from 5 mg to 900 mg

■ Favorable Pharmacokinetic Profile

- ✓ Demonstrated a favorable pharmacokinetic profile with dose-dependent exposures (C_{max} and AUC)
- ✓ Elimination half-life ranging 7-9 hours after administration at 100 mg or above

■ Promising Pharmacodynamic effects

- ✓ Demonstrated a dose-dependent and marked inhibition of basophil and B-cell activation

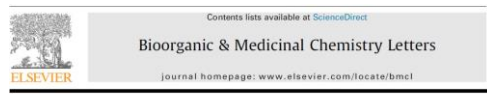
■ Clinical Development Plan

- ✓ Plan to initiate a Multiple Ascending Dose (MAD) study consisted of three parts with a new formulation
 - BA part: To evaluate the bioavailability (BA) of capsule and tablet formulations
 - MAD part: To evaluate the safety, tolerability, PK and PD in the 2-week multiple ascending dose (MAD) part
 - SPT part: To evaluate the effect on allergen-induced skin reaction in the skin prick test (SPT) to assess the potential of AS-0871 for the treatment of Chronic Spontaneous Urticaria (CSU), a disease with high unmet needs
- ✓ The bioavailability (BA) part in human with the capsule formulation will be conducted in H2 2021 and the GMP production of the drug product is underway
- ✓ New tablet formulation is under development, and the BA part is also planned
- ✓ Plan to initiate the MAD part with the new formulation based on the BA studies

AS-0871: Excellent Kinase Selectivity

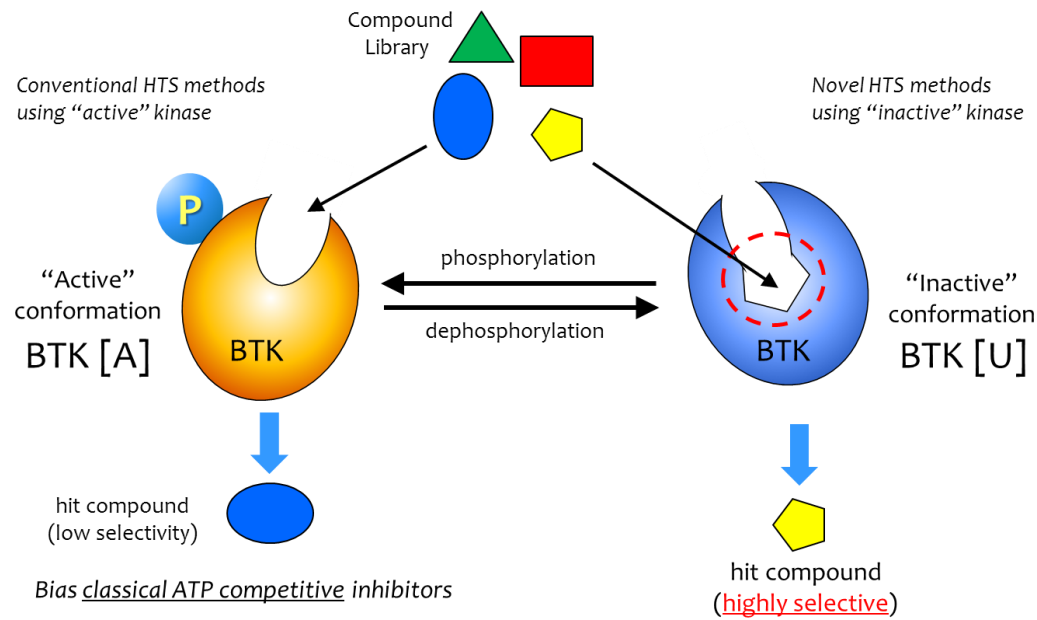


◆ Targeting Inactive Conformation of BTK



Journal of Medicinal Chemistry
Design and Synthesis of Novel Amino-triazine Analogues as Selective Bruton's Tyrosine Kinase Inhibitors for Treatment of Rheumatoid Arthritis
Wataru Kawahata,*¹ Tokiko Asami, Takao Kiyoi, Takayuki Irie, Haruka Taniguchi, Yuko Asamitsu, Tomoko Inoue, Takahiro Miyake, and Masaaki Sawa²
Research and Development, Carna Biosciences, Inc., 3rd Floor, RMA, 1-5-5 Minatojima Minamimachi, Chuo-ku, Kobe 650-0047, Japan

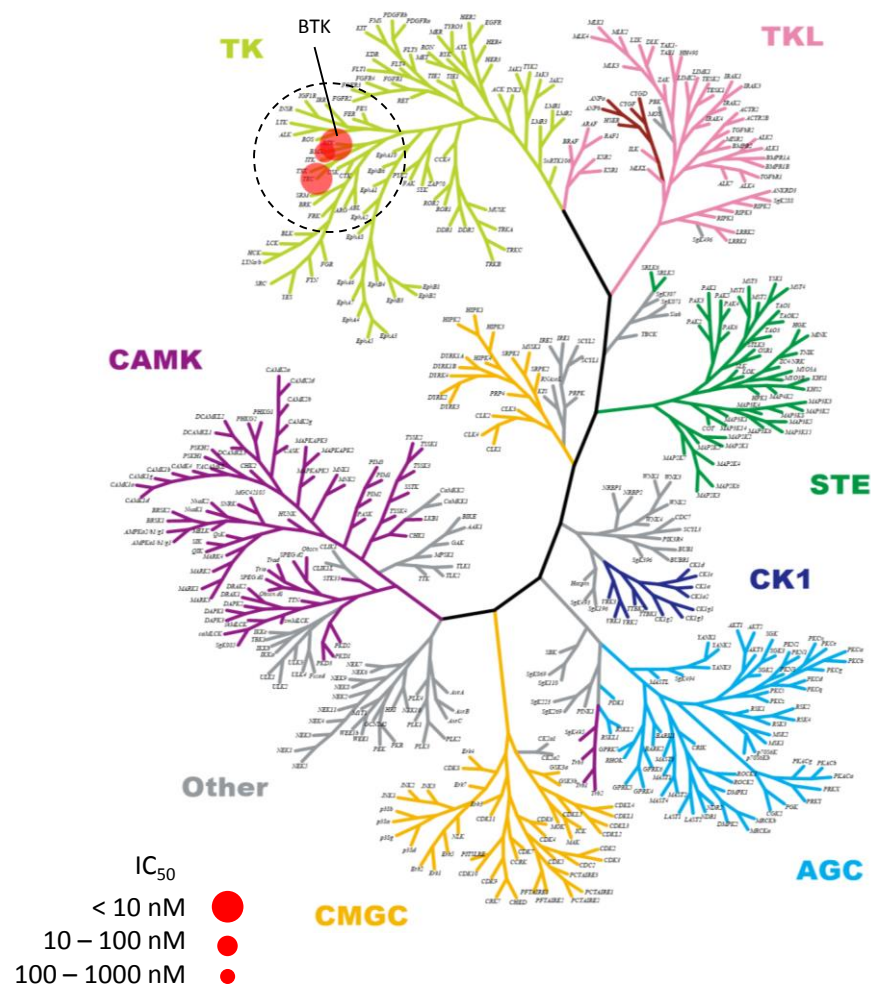
TR-FRET binding assay targeting unactivated form of Bruton's tyrosine kinase
Tokiko Asami¹, Wataru Kawahata, Masaaki Sawa
Carina Biosciences, Inc., RMA 3F, 1-5-5 Minatojima Minamimachi, Chuo-ku, Kobe 650-0047, Japan



| | BTK IC ₅₀ (nM) | |
|---------|---------------------------|---------|
| | BTK [A] | BTK [U] |
| AS-0871 | 3.4 | 0.3 |

◆ Kinase Selectivity Profiling

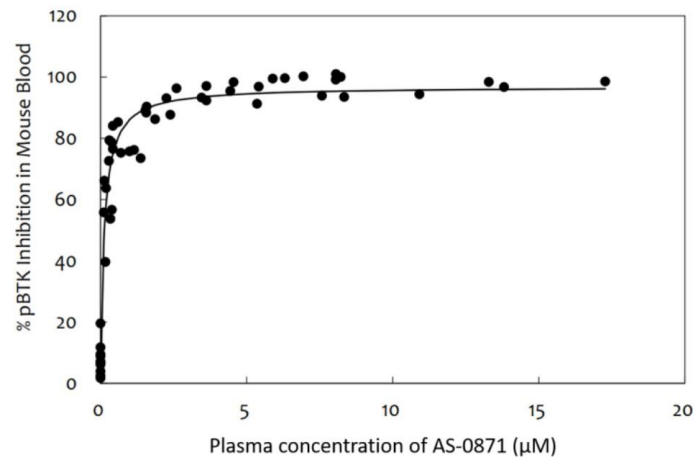
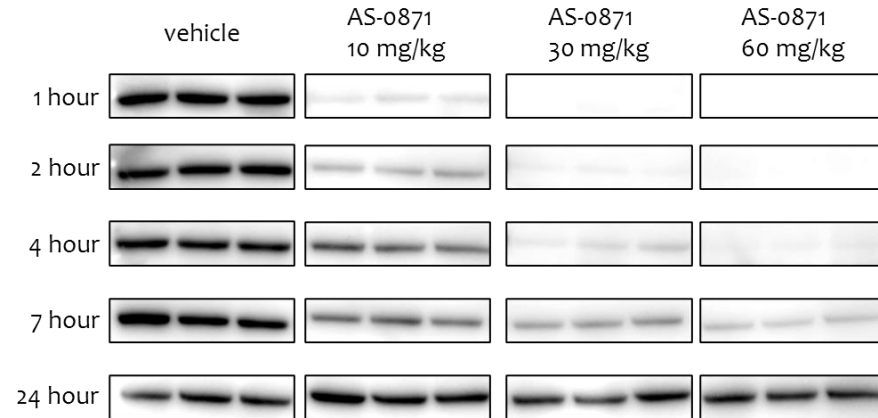
Only inhibited 2 other kinases in a total of 312 kinases tested at 0.3 μ M concentration.



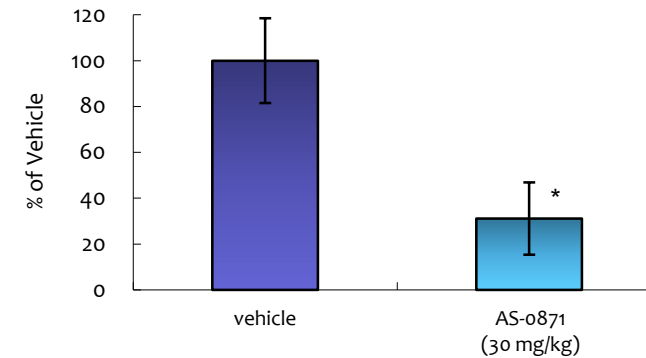
AS-0871: In Vivo Therapeutic Efficacy

◆ PK/PD Study

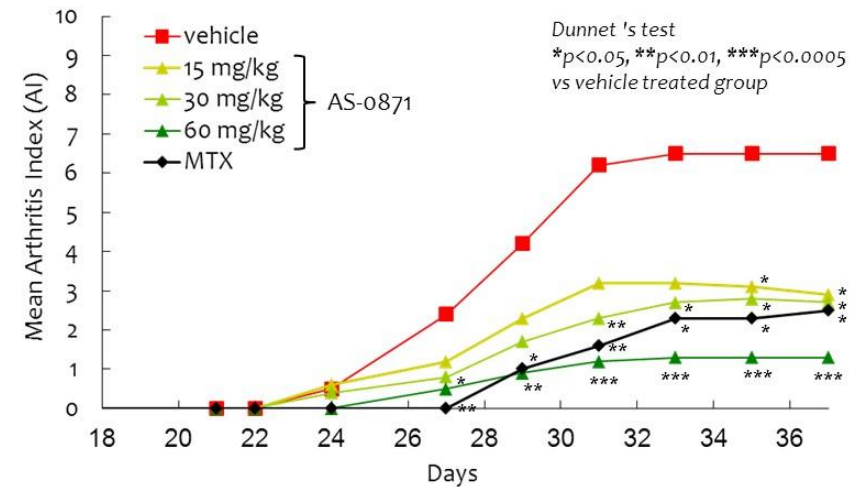
Auto-phosphorylation status of BTK was measured following oral single administration of AS-0871



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

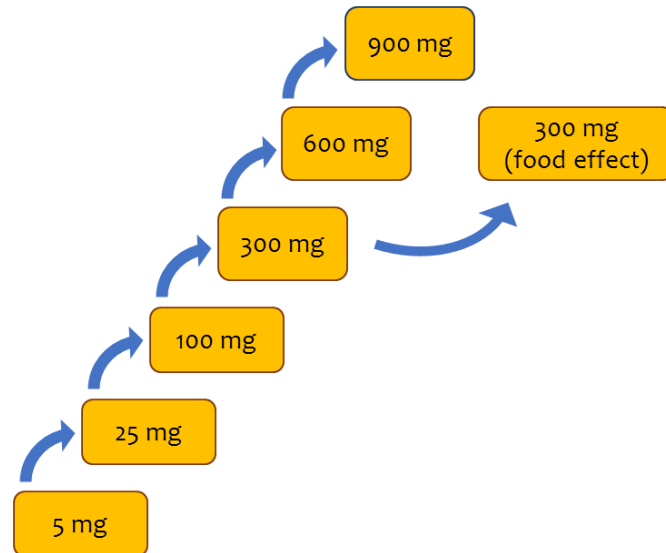


◆ Collagen-induced arthritis (CIA) mouse model (n=10)



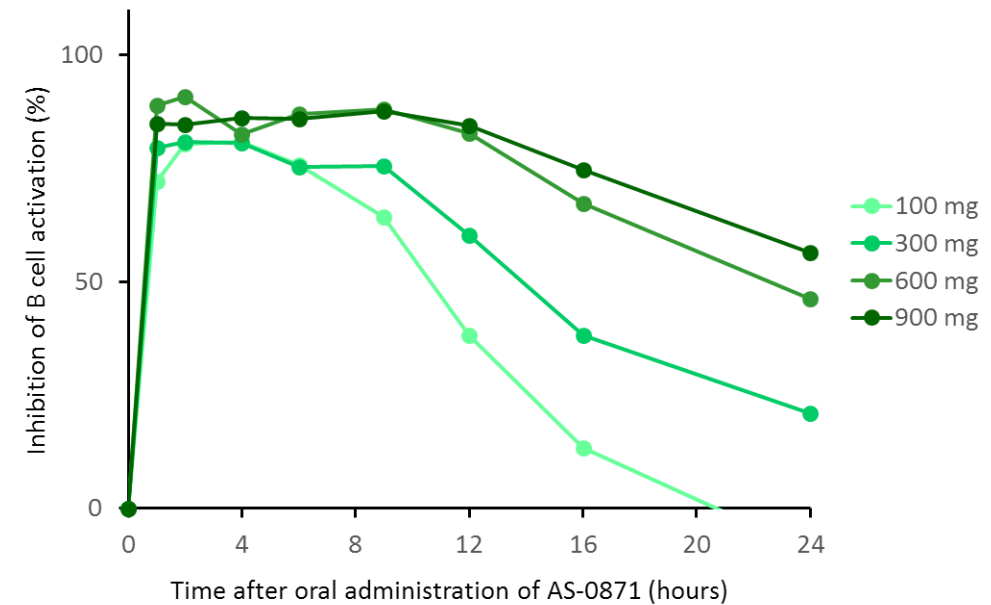
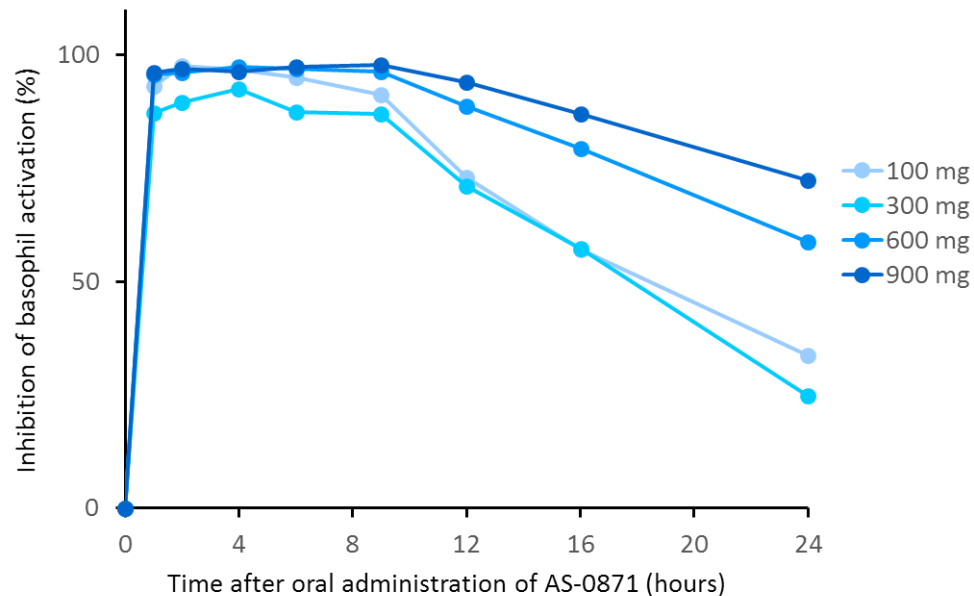
SAD Part (Completed)

| Step 1 Single Ascending Dose Study (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 |



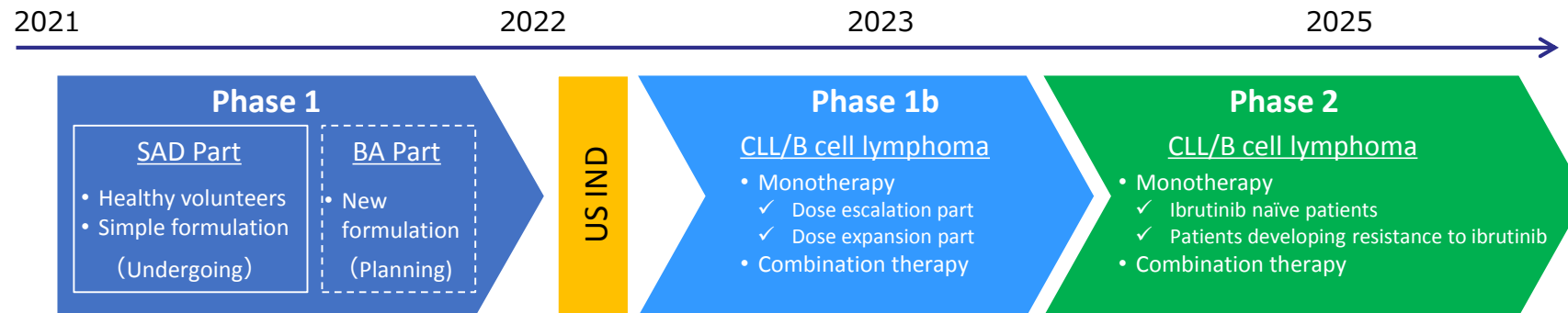
- ✓ AS-0871 is well-tolerated without any safety concerns.
- ✓ Favorable pharmacokinetic profile.
- ✓ Blood samples to assess PD effects were analyzed for evaluation of the B-cell and basophil responses. Administration of AS-0871 at 100mg or above resulted in strong inhibition of B-cell and basophil activation.
- ✓ Switching to a new formulation in the MAD study.

- ✓ Pharmacodynamic study demonstrated that subjects who received AS-0871 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 AS-0871 achieved therapeutic plasma levels needed to inhibit B cells and basophils activation, suggesting that AS-0871 has a potential to become a new treatment option for inflammatory diseases.



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
- Potential applications for autoimmune diseases
- Plan to accelerate the clinical studies utilizing the clinical data of BioNova, the licensee in Greater China



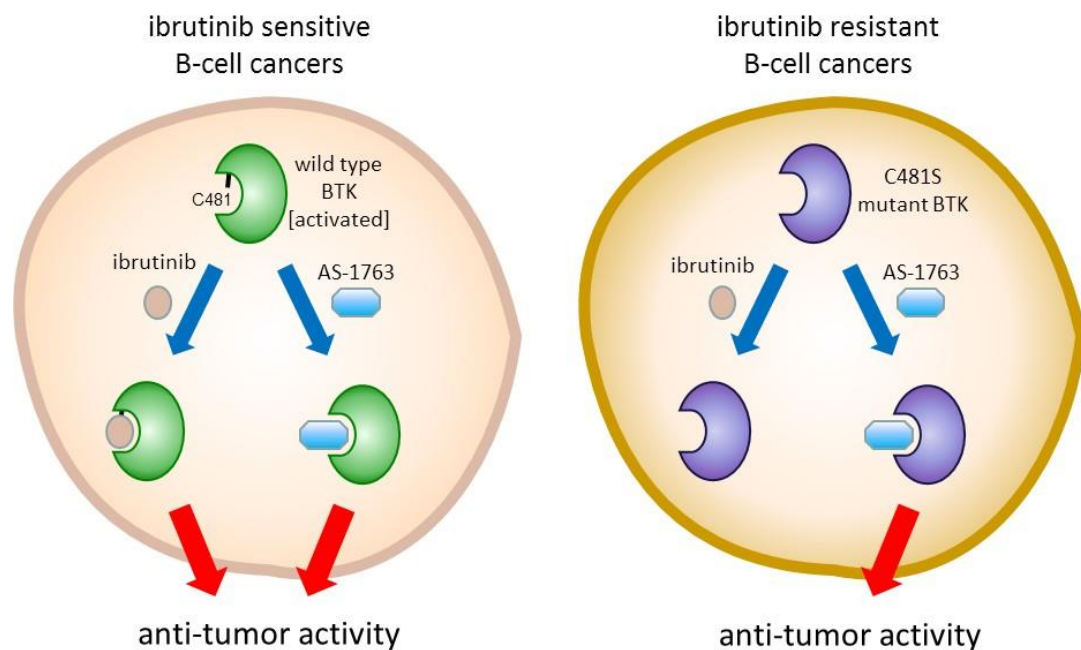
BA: Bioavailability
IND: Investigational New Drug application

■ Phase 1 Study in healthy volunteers

- ✓ The Phase 1 study in healthy volunteers was initiated in the Netherlands in H1 2021
- ✓ Dosing in the SAD part was completed
- ✓ Well-tolerated without any safety concerns
- ✓ Pharmacokinetic profile was also found favorable

■ Clinical Development Plan

- ✓ New formulation has been developed and the GMP production is underway
- ✓ The BA part in the Phase 1 using new formulation will be conducted in H2 2021
- ✓ Plan to initiate Phase 1b study in patients with chronic lymphocytic leukemia (CLL)/B cell lymphoma in the U.S. in 2022
- ✓ Request a pre-IND (Investigational Drug Application) meeting with FDA
- ✓ Aim to conduct clinical studies efficiently, collaborating with BioNova



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

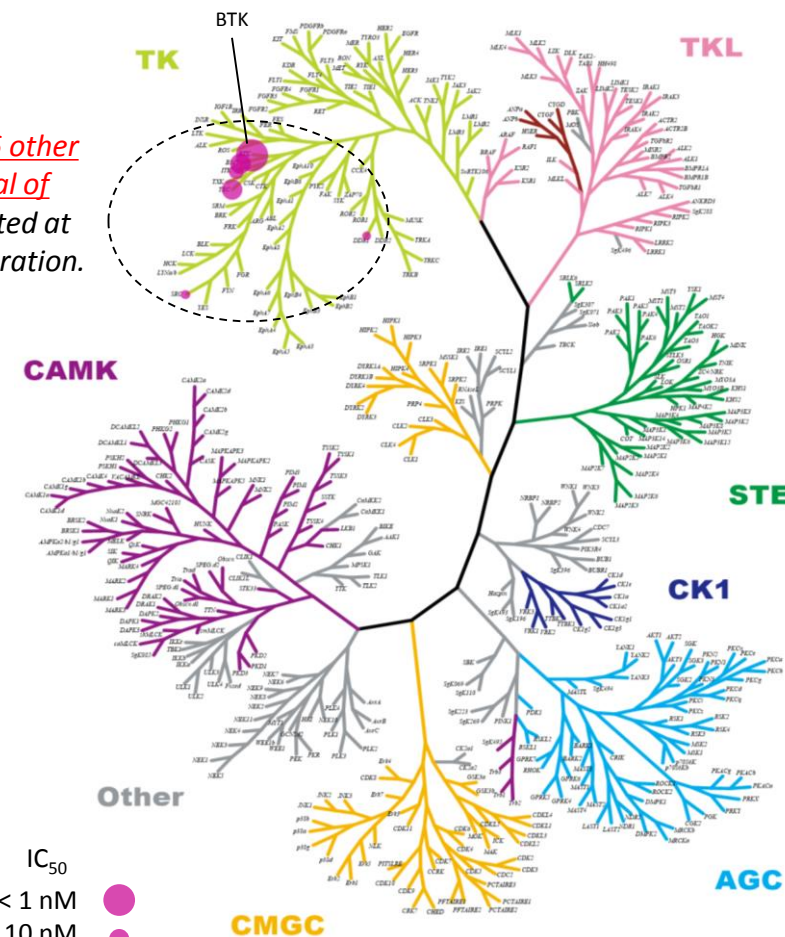
| | IC ₅₀ (nM) | |
|---|-----------------------|-----------|
| | AS-1763 | ibrutinib |
| Autophosphorylation BTK (Ramos) | 1.4 | 1.1 |
| CD69 activation (Human whole blood) | 11 | 8.1 |
| Cancer cell growth OCI-Ly10 cells | 1.8 | 0.75 |
| Cancer cell growth OCI-Ly10 [BTK C481S] cells | 20 | 1030 |
| 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

◆ Kinase selectivity profiling

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

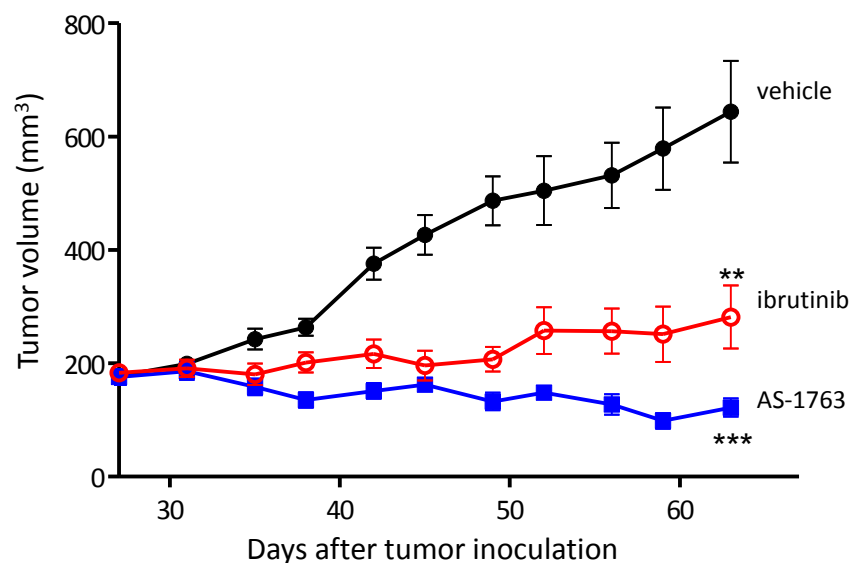


IC₅₀
 < 1 nM
 1 – 10 nM
 10 – 100 nM
 100 – 300 nM

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



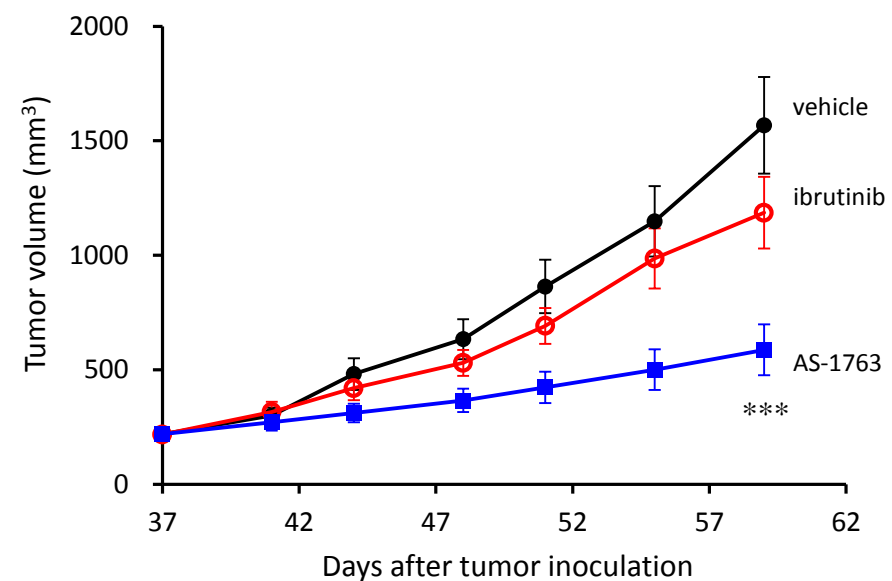
- ◆ In vivo antitumor effects of AS-1763 on human B-cell non-Hodgkin lymphoma cell line, OCI-LY10 tumor xenograft mouse model (n=8-10)



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)

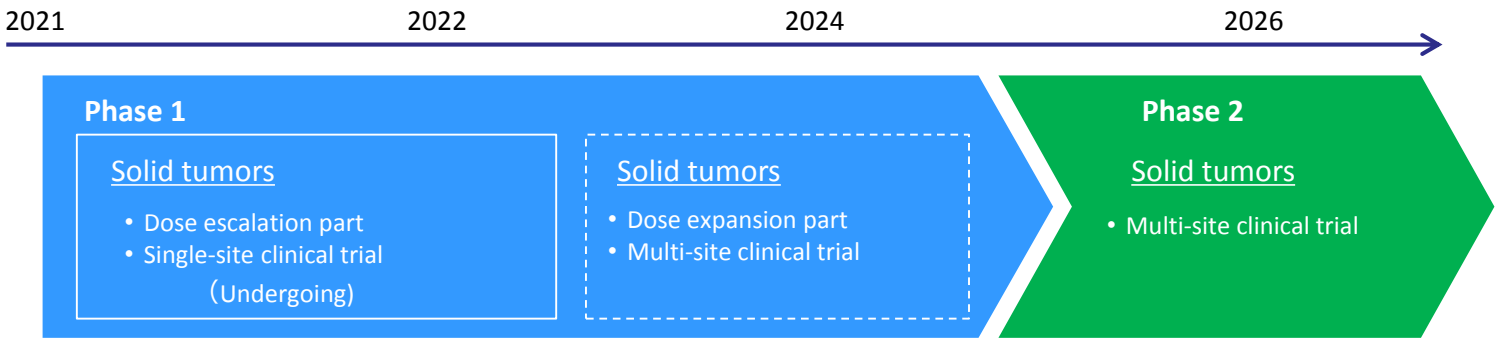


Ibrutinib: 25 mg/kg QD
AS-1763: 60 mg/kg BID

***: p<0.001

AS-0141 : Targeting Cancer

- Small molecule CDC7 inhibitor
- High kinase selectivity
- Potential First-in-class drug
- Orally available
- 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



■ Phase 1 Study in patients

- ✓ The Phase 1 study in patients with unresectable, advanced, recurrent, or metastatic solid tumors was initiated in Japan in H1 2021
- ✓ The study consists of two parts, a dose escalation and an expansion
- ✓ The dose escalation part is ongoing
- ✓ No dose-limiting toxicity (DLT) has been observed so far
- ✓ Approved dose escalation to Cohort 3 (dose level 3)

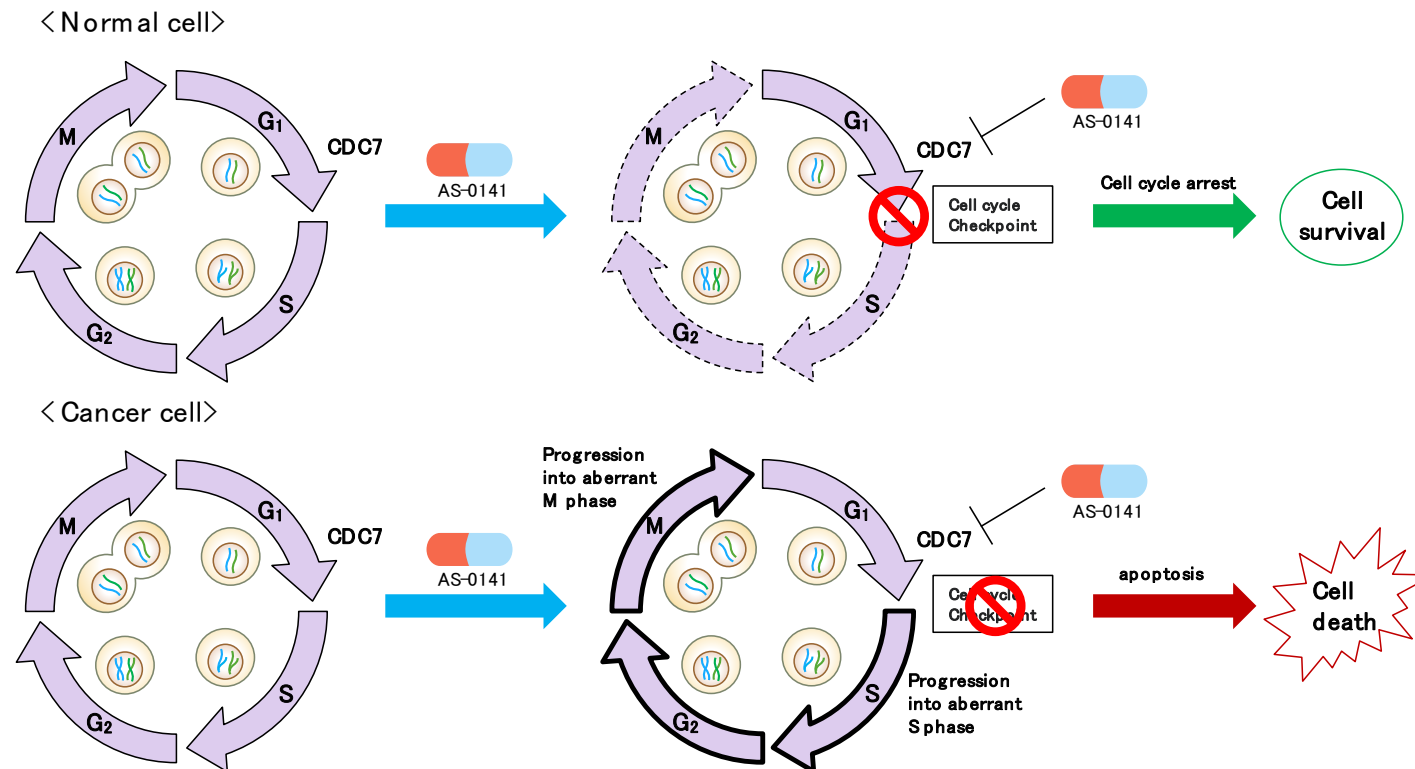
■ Clinical Development Plan

- ✓ Continuing dose escalation
- ✓ Select dose level for expansion phase
- ✓ Exploring multiple types of solid tumors

AS-0141: Highly Selective CDC7 Inhibitor

■ CDC7 kinase inhibitor

CDC7 (cell division cycle 7) is a serine-threonine kinase that plays a critical role in DNA synthesis and is required for the activation of DNA replication origins throughout the S phase of the cell cycle. Inhibition of CDC7 in cancer cells causes lethal S phase or M phase progression, whereas normal cells survive, most likely through induction of cell cycle arrest at the DNA replication checkpoint. It has been reported in the literature that CDC7 is overexpressed in many cancers. Therefore, CDC7 is an attractive target for cancer drug development.



AS-0141: Time-Dependent Inhibitor of CDC7

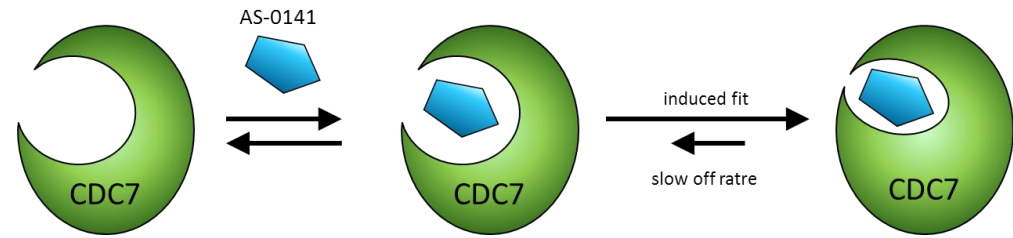
◆ AS-0141 has a unique inhibitory mechanism for CDC7 kinase (time-dependent inhibition)

◆ AS-0141 inhibits CDC7 in a reversible fashion but has a very slow off-rate



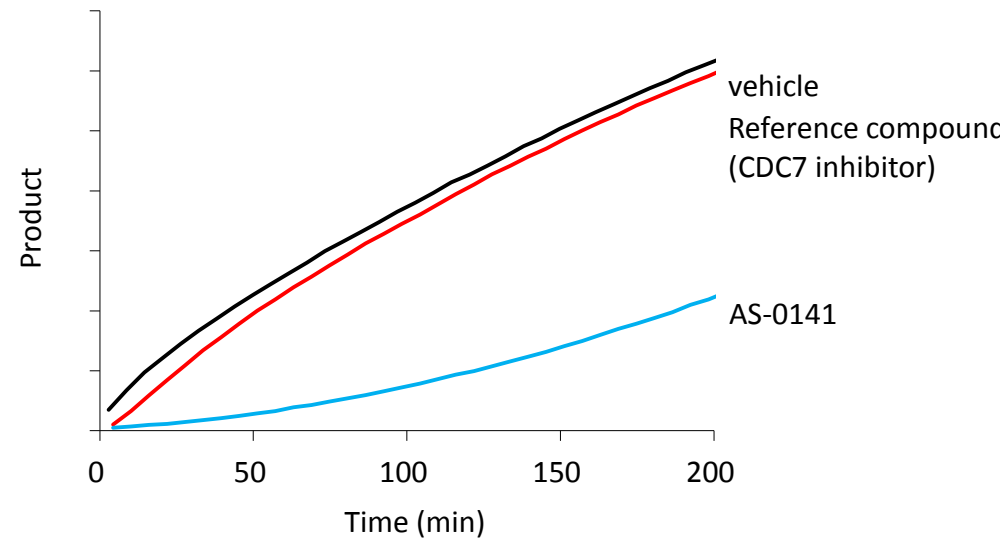
Research paper
Discovery of novel furanone derivatives as potent Cdc7 kinase inhibitors
Takayuki Irie^{a,*}, Tokiko Asami^a, Ayako Sawa^a, Yuko Uno^a, Mitsuharu Hanada^a, Chika Taniyama^b, Yoko Funakoshi^c, Hisao Masai^c, Masaaki Sawa^b
^a Research and Development, Carina Biosciences, Inc., 3F BMA, 1-5-5 Minamijima Minamimachi, Chuo-ku, Kobe, 650-0047, Japan
^b Research and Development Department, SRI Biotech Co., Ltd., Izumi Garden Tower 3F, 1-6-1 Akappongi, Minato-ku, Tokyo 106-6018, Japan
^c Department of Genome Medicine, Tokyo Metropolitan Institute of Medical Science, 2-1-6 Kamikiazawa, Setagaya-ku, Tokyo 158-8501, 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



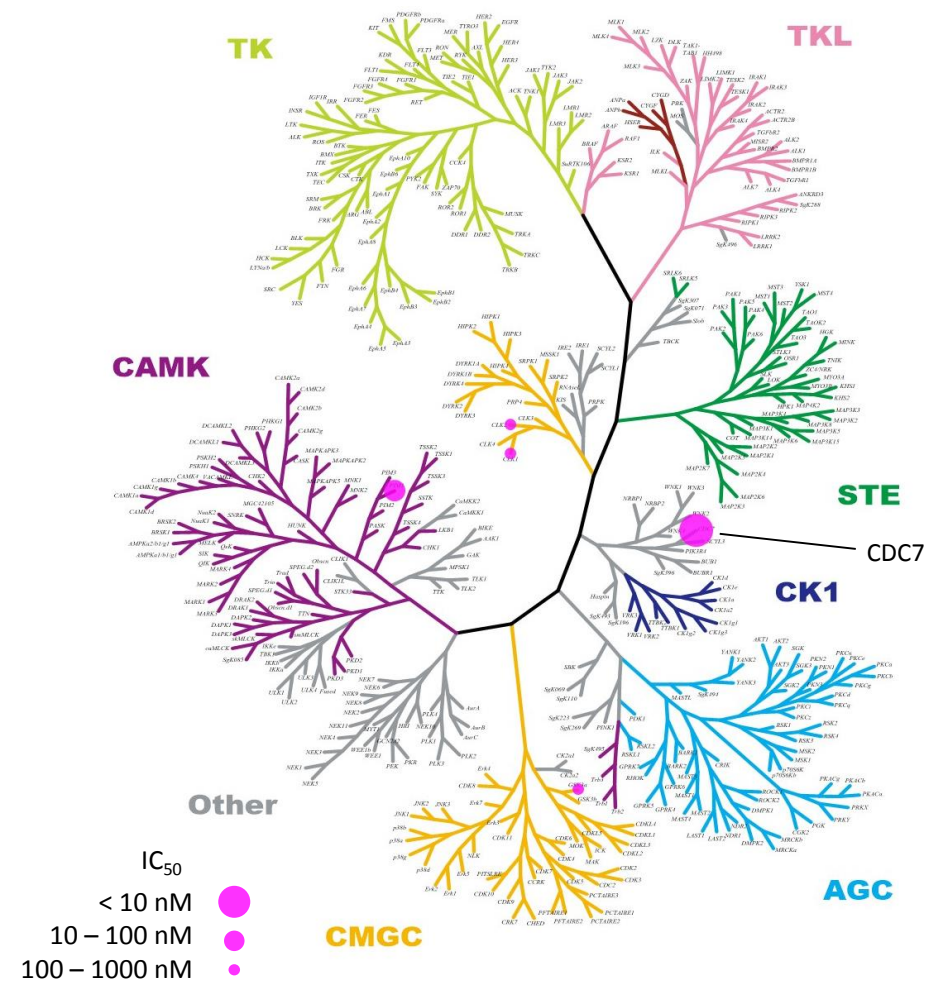
| Inhibitory potency (IC ₅₀) for CDC7 in the presence of 1 mM ATP | |
|---|--------------------|
| Without Preincubation | With Preincubation |
| 503 nM | 2.4 nM |

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



◆ Kinase Selectivity Profiling

In the presence of 1 mM ATP with preincubation



◆ IC₅₀ values of hit kinases (at 1 mM ATP)

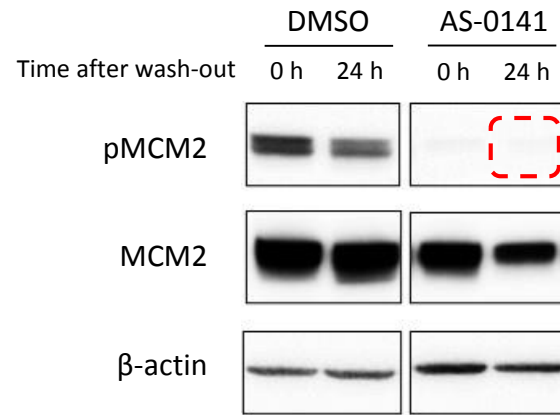
| | IC ₅₀ (nM) | |
|-------|-----------------------|-----|
| | Preincubation | |
| | - | + |
| CDC7 | 503 | 2.4 |
| PIM1 | 30 | 34 |
| CLK1 | 212 | 206 |
| CLK2 | 270 | 227 |
| GSK3a | 189 | 251 |

CDC7 is the only kinase that shows preincubation effect

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

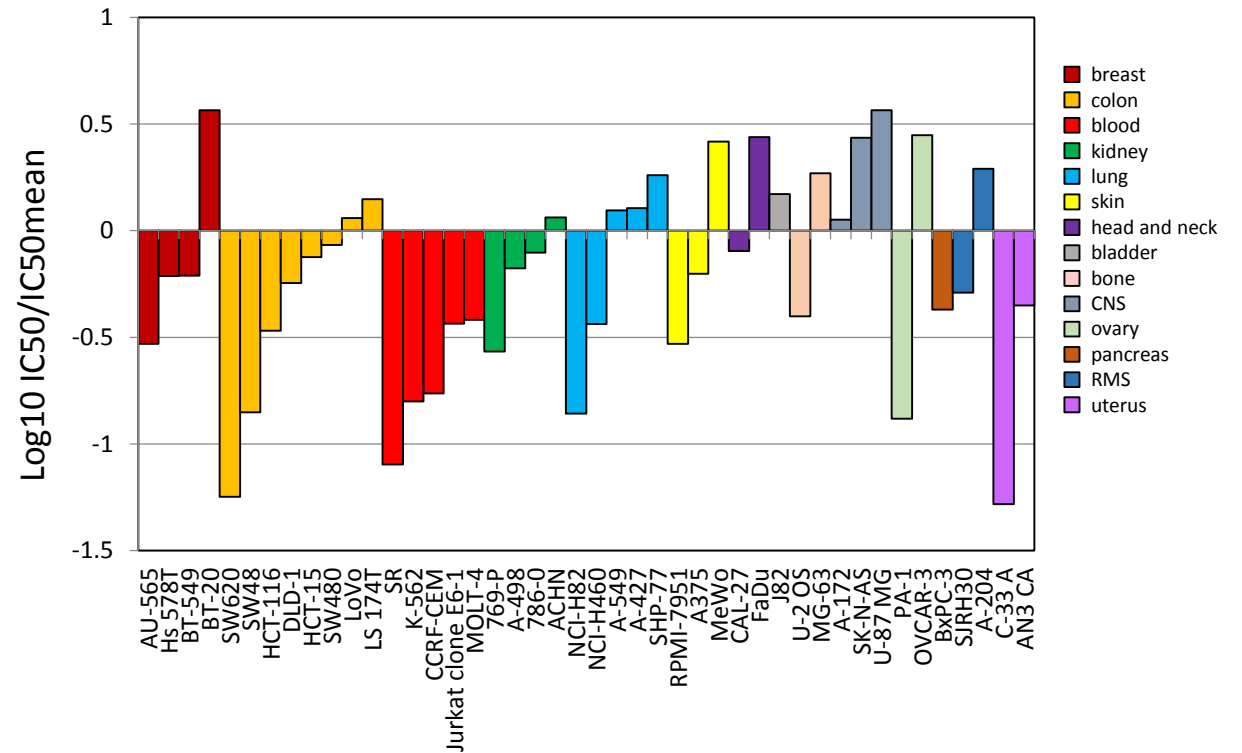
◆ Prolonged inhibition in cells

Human colon cancer cell line, Colo-205 cells were treated with DMSO control or AS-0141. 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 AS-014 on the phosphorylation of MCM2 in cells continued up to 24 h after washing out

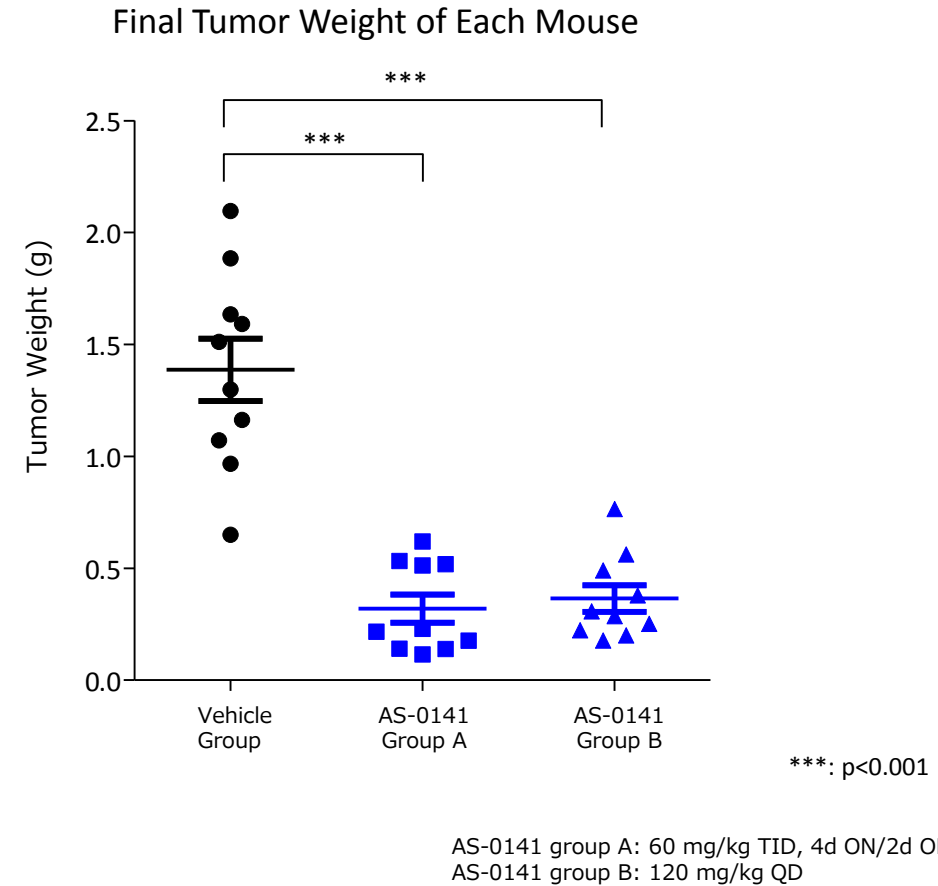
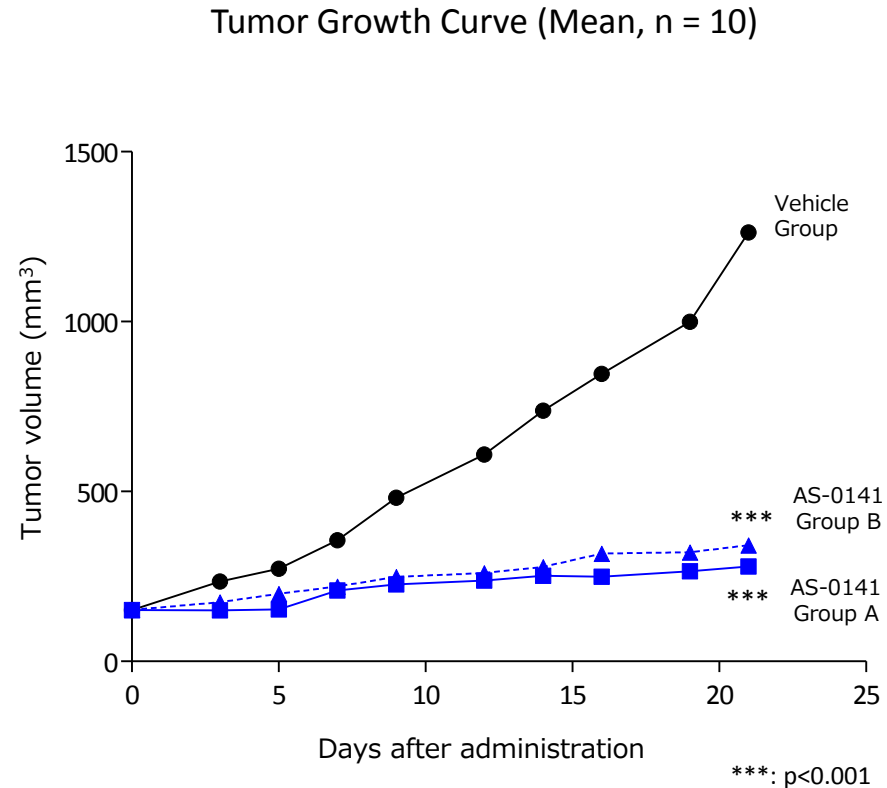
◆ AS-0141 potently inhibited growth in a wide range of tumor cell lines, including solid and hematological tumors



44 Cancer cell lines (Oncolines at NTRC)

AS-0141: Robust In Vivo Antitumor Efficacy

- ◆ In vivo antitumor efficacy of AS-0141 in a SW620 (human colon cancer) xenograft mouse model



- At Drug Discovery Support business, sales in Q1-3 were slightly behind the full year plan while operating profit was stronger than expected.
 - ✓ In North America, sales declined yoy but they were stronger than expected thanks to sales to Gilead.
 - ✓ In Japan, sales of kinase proteins were strong while profiling service and agent business (cell-based assay and X-ray crystallography) were weaker than expected. Weak agent business sales had limited impact on operating profit as the profitability of the agent business is relatively low.
 - ✓ Sales of Cell-based assay service using NanoBRET™ technology were strong. Launched a full-panel assay service (192 kinds of kinases) aiming to boost sales further.
 - ✓ In China, orders from the agent were strong. However, sales in September and October were temporarily affected by the delay in the import custom clearance due to COVID-19. Working closely with the agent to deliver products as soon as possible.
- Focusing on the growth areas to achieve full year sales target.
 - ✓ Increase sales from newly established biotech companies in the U.S.
 - ✓ Increase sales of custom kinase proteins only Carna can offer.
 - ✓ Conduct wet lab work for AI drug discovery companies to increase profiling sales.

FY2021 Q3 Consolidated Financial Results



| (JPY million) | FY2020 Q3 Actual | FY2021 Q3 Actual | YoY Change | FY2021 Plan | |
|-----------------------|---------------------|---------------------|----------------|----------------|---|
| Sales | 847 | 636 | -211 -24.9% | 923 | -Slightly behind the FY sales plan due to weaker than expected sales in Japan. -Received an upfront payment from licensing in Q1 FY2020. |
| Operating Profit/Loss | (615) | (1,169) | -553 | (1,811) | |
| Ordinary Profit/Loss | (625) | (1,171) | -546 | (1,816) | |
| Net Profit/Loss | (649) | (1,178) | -528 | (1,825) | |
| R&D Cost | 941 | 1,310 | +368 +39.1% | 1,981 | -Investment in clinical studies. |

Note 1: Rounded down to the nearest million yen.

Note 2: YoY change % for Operating Profit/Loss, Ordinary Profit/Loss, and Net Profit/Loss are not presented since losses were recorded.

Note 3: FY2021 plan was disclosed on February 12, 2021.

FY2021 Q3 Results by Business Segment



| (JPY million) | FY2020 Q3 Actual | FY2021 Q3 Actual | YoY Change | FY2021 Plan | vs. FY Plan | |
|-----------------------------|---------------------|---------------------|----------------|----------------|----------------|---|
| Total Sales | 847 | 636 | -211 -24.9% | 923 | 69.0% | |
| ddSP business | 794 | 636 | -158 -19.9% | 923 | 69.0% | Slightly behind the FY sales plan due to weaker than expected sales in Japan. |
| ddRD business | 53 | — | -53 | — | — | Received an upfront payment from licensing in Q1 2020. |
| Total Operating Profit/Loss | (615) | (1,169) | -553 | (1,811) | — | |
| ddSP business | 347 | 199 | -148 -42.7% | 207 | 95.8% | Sales of high-margin kinase proteins were robust. |
| ddRD business | (963) | (1,368) | -404 | (2,019) | — | Investment in clinical studies. |

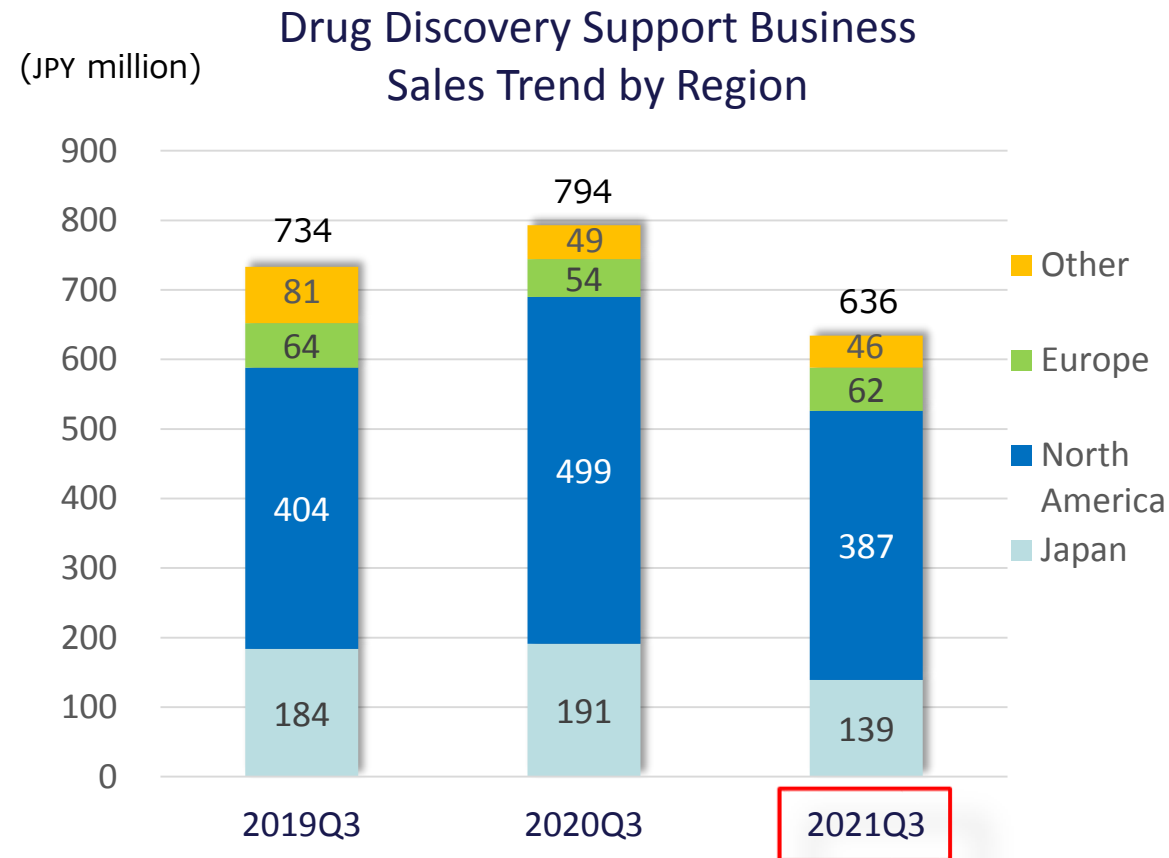
Note 1: Rounded down to the nearest million yen.

Note 2: YoY change % for consolidated operating profit/loss and ddRD operating profit/loss are not presented since losses were recorded.

Note 3: FY2021 plan was disclosed on February 12, 2021.

Note 4: ddRD: Drug Discovery R&D business, ddSP: Drug Discovery Support Business

FY2021 Q3 Discovery Support (ddSP) Business Sales by Region



- ▣ Japan: Decreased 26.8% YoY
 - Sales of profitable kinase proteins were robust, offsetting weak profiling sales.
 - Agent business (cell-based assay service and X-ray crystallography) were weak but the impact on the operating profit was limited as its profitability is relatively low.
- ▣ North America: Decreased 22.4% YoY
 - Contribution from sales to Gilead continued.
 - NanoBRET assay service showed strong growth.
 - Agent business (cell-based assay service) were weak.
- ▣ Europe: Increased 13.6% YoY
 - Kinase proteins, profiling service and NanoBRET assay service were robust.
- ▣ Other: Decreased 5.2% YoY
 - Orders from Chinese agent were robust.
 - Sales in September and October were temporarily affected by the delay in the import custom clearance in China.

Balance Sheet

| (JPY million) | As of Dec. 31, 2020 | As of Sep. 30, 2021 | Change | Reason for changes |
|----------------------------------|------------------------|------------------------|--------|---|
| Current assets | 4,708 | 4,084 | -624 | |
| Cash and deposits | 4,299 | 3,790 | -508 | |
| Non-current Assets | 127 | 127 | +0 | |
| Total assets | 4,835 | 4,211 | -624 | |
| Current liabilities | 727 | 584 | -142 | |
| Non-current liabilities | 284 | 147 | -136 | Long term loans payable -105 Bonds payable -28 |
| Total liabilities | 1,011 | 731 | -279 | |
| Total net assets | 3,824 | 3,479 | -344 | Capital stock and capital surplus +810, Retained earnings -1,178 |
| Total liabilities and net assets | 4,835 | 4,211 | -624 | |

| | | |
|----------------------------|-----------|-----------|
| Shareholders' equity ratio | 79.0% | 82.4% |
| BPS | 308.0 yen | 264.4 yen |
| PBR | 3.9x | 4.2x |
| Share price of Carna | 1,212 yen | 1,099 yen |

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



“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.