

Financial Results FY2021

(January to December 2021)

Carna Biosciences, Inc.



February 10, 2022

Stock Code: 4572



- AS-0871 (BTK inhibitor) : Initiated Multiple Ascending Dose (MAD) study of Phase 1 study in Europe.
- AS-1763 (BTK inhibitor) : Completed dosing in Single Ascending Dose (SAD) part of Phase 1 study in Europe and initiated bioavailability (BA) part.
- AS-0141 (CDC7 inhibitor) : Initiated Phase 1 study in Japan.
- Received the first milestone payment from Gilead regarding novel immuno-oncology program licensed in 2019.
- STING antagonist: Advanced to preclinical stage.
- Issued Series 19th Subscription Rights Shares.

Key Milestones for 2021



Business		Milestones for 2021	Achievement in 2021
	AS-0871	Initiate Ph1 MAD part (H2 2021)	Initiated Ph1 MAD part in Dec. 2021
ddRD	AS-1763	Initiate Ph1 (H1 2021)	Initiated Ph1 in Apr. 2021
	AS-0141	□Initiate Ph1 (H1 2021)	☑Initiated Ph1 in June 2021
	Research program	Bring one or more programs in preclinical stage	STING antagonist was advanced to preclinical stage
ddSP		 Achieve sales target of JPY920 mn Launch new products Expand NanoBRET service Propose project-based service to collaborate with clients, leveraging Carna's drug discovery technology 	 Actual sales were JPY889 mn, 96% of initial target Launched 31 new kinase protein products and 4 new EGFR mutant targets to profiling service Expand NanoBRET service Propose project-based service to collaborate with clients, leveraging Carna's drug discovery technology



Drug Discovery R&D (ddRD) Business



- Clinical development
 - ✓ AS-0871: Completed Phase 1 SAD study and initiated MAD study in the Netherlands.
 - ✓ AS-1763: Completed dosing in Phase 1 SAD part and initiated BA part in the Netherlands.
 - ✓ AS-0141: Initiated Phase 1 study in Japan.
- Research
 - ✓ STING antagonist was advanced to preclinical stage.
- \checkmark Extended the term of joint research with Sumitomo Dainippon Pharma.
- Other highlights
 - Received the first milestone payment from Gilead under the agreement to develop and commercialize small molecule compounds in immuno-oncology.
 - BioNova Pharmaceuticals, the licensee of AS-1763 in Greater China, submitted an Initial New Drug(IND) application in China for AS-1763 (BN102) in January 2022.
 - ✓ The patents for AS-1763 were registered in Japan and China.
 - ✓ Published research papers in Journal of Medicinal Chemistry on the discovery of AS-1763 and AS-0141.







♥ BrickellBio

In Feb. 2022, Carna and Brickell Biotech entered into a license agreement to grant Brickell the exclusive, worldwide rights to develop and commercialize Carna's portfolio of novel, potent, and orally available STING antagonists.

Deal size	 ✓ Upfront payment of \$2 million ✓ Success-based development, regulatory, and sales milestone payment of up to \$258
Royalty	 Tiered royalty payments ranging up to 10% of net sales

- Carna initiated the STING antagonist program in 2019 to create next wave of pipeline.
- The program was advanced to preclinical stage in Q4 2021.
- Carna can continue its own research on STING modulator (agonist/antagonist).
 (The antagonist research is limited to a new chemical scaffold in the filed of oncology.)

Successfully licensed non-kinase target project in a short period of time.





Robust Drug Pipeline (as of Q1 2022)





<Oncology>

Compound	Target	Indication	Discovery/Preclinical	Clinical	Partner
AS-0141	CDC7/ASK	Cancer			
Small Molecule	Kinase	Immuno-Oncology	Out-license	ed to Gilead	🧭 GILEAD
AS-1763	ВТК	Blood Cancer			Bionova *
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	Joint rese Sumitomo Dair	arch with nippon Pharma	Sumitomo Dainippon Pharma
AS-0871	ВТК	Immune-inflammatory diseases			
Small Molecule	N/A	Malaria			
Small Molecule	STING	Immune-inflammatory diseases	Out-license	d to Gilead	⊖ BrickellBio

✓ 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 		



AS-0871 : Phase 1 Clinical Trial in Progress



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

- ✓ Safe and well-tolerated at all dose levels
- ✓ Favorable Pharmacokinetic Profile
- Promising Pharmacodynamic effects
- ✓ Conducted using simple formulation

Developing several new formulations



AS-0871: Excellent Kinase Selectivity



Targeting Inactive Conformation of BTK



Journal of

_	BIR 1650 (1111)	
	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)



ARNA BIOSCIENCES



SAD Part (Completed)

Step 1 Single Ascending Dose Study (SAD)	Step 2
 6 dose levels (8 subjects/cohort) Placebo controlled (6 active / 2 placebo) Safety and tolerability Pharmacokinetics and pharmacodynamics 	• 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.

Pharmacodynamics of AS-0871

- 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.



ARNA BIOSCIENCES



AS-1763: Targeting <u>Blood Cancer</u>			
 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

AS-1763 : Phase 1 Clinical Trial in Progress





- Plan to present the Phase 1 data at AACR2022
- ◆ Aim to conduct clinical studies efficiently, collaborating with BioNova







Journal of

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	50-fold Stronger activity
Normal cell growth HEL299 cells	6370	6870	

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



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

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) 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



ARNA BIOSCIENCES



AS-0141: Targeting <u>Cancer</u>			
 Small molecule CDC7 inhibitor High kinase selectivity Potential First-in-class drug 	 Potent anti-proliferative activity against various cancer cell lines Demonstrated strong anti-tumor activity in several human tumor xenograft models 		
Orally available	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 primary objective is to assess safety, tolerability, maximum tolerated dose, preliminary anti-tumor activity, and pharmacokinetics(PK)/pharmacodynamics(PD) as well as to determine recommended Phase 2 dose.
- ✓ The dose escalation part is ongoing.
- ✓ No dose-limiting toxicity (DLT) has been observed so far.
- ✓ Advanced to dose level 3 (Cohort 3) underway.

Clinical Development Plan

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

CARNA BIOSCIENCES

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



Preincubation	Preincubation	
503 nM	2.4 nM	

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

AS-0141 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.



AS-0141: High Kinase Selectivity



Kinase Selectivity Profiling

In the presence of 1 mM ATP with preincubation



• IC_{50} values of hit kinases (at 1 mM ATP)

	IC ₅₀ (nM)		
	Preincucabation		
	-	+	
CDC7	503 — 210	→ 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.

CARNA BIOSCIENCES

• 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

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

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





AS-0141: Robust In Vivo Antitumor Efficacy

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



AS-0141 group A: 60 mg/kg TID, 4d ON/2d OFF AS-0141 group B: 120 mg/kg QD

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

ARNA BIOSCIENCES

CARNA BIOSCIENCES

- ✓ The cyclic GMP-AMP synthase (cGAS) stimulator of interferon genes (STING) signaling pathway plays a central role in innate immunity.
- ✓ The cGAS-STING signaling is activated in response to the presence of cytosolic DNAs produced by microbial infection or cellular stress. The activated cGAS-STING signaling induces type I interferons (IFNs) production to trigger immune responses for host defense.
- ✓ It has been reported that aberrant cGAS-STING signaling is implicated in the pathogenesis of several diseases. Mutations in several genes have been identified that cause the constitutive activation of the cGAS-STING pathway, resulting in severe autoinflammation in lung, kidney, joint, etc.
- ✓ Moreover, numbers of non-genetic diseases are also suggested to be associated with the aberrant cGAS-STING signaling. These include a subset of systemic lupus erythematosus (SLE), rheumatoid arthritis (RA), and several cancers.
- There is a high unmet medical need to develop novel STING antagonists to treat these diseases. (ref.1)
- ✓ Recently, a research paper was published suggesting that STING antagonist has a potential to treat severe lung inflammation induced by SARS-Cov-2. (ref.2)

Decout A., et al. Nat Rev Immunol. 2021 Sep;21(9):548-569.
 Di Domizio J., et al. Nature. 2022 Jan 19. doi: 10.1038/s41586-022-04421-w.

Cytosolic DNAs produced by microbial infection or cellular stress





Discovery Support (ddSP) Business



- At ddSP, full year sales reached 96% of initial plan.
 - ✓ In North America, contribution from sales to Gilead continued.
 - 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.
 - ✓ In China, full year sales beat the initial plan although sales were temporarily affected in September and October by the delay in the import custom clearance due to COVID-19. Continue working closely with the agent to expand sales further.
- Robust start for the new service
 - ✓ Launched a full-panel assay service (192 kinds of kinases) for cell-based assay service using NanoBRET[™] technology developed by Promega.
- Expanding lineup of kinase proteins and profiling service
- ✓ 31 kinase protein products, including high-demand mutant kinase biotinylated kinases, have been newly added to the line-up.
- ✓ 4 EGFR mutant targets were added to the profiling service.



FY2021 Financial Results

FY2021 Consolidated Financial Results



(JPY million)	FY2020 Actual	FY2021 Actual	YoY Change	FY2021 Initial Plan	FY2021 Dec. 23 Revised Plan	
Sales	1,133	2,017	+884 +78%	923	2,006	Sales were strong compared to the initial plan and the last year thanks to the first milestone from Gilead.
Operating Profit/Loss	(1,057)	(531)	+525	(1,811)	(548)	Operating loss was smaller compared to the initial plan and the last year thanks to milestone payment.
Ordinary Profit/Loss	(1,077)	(522)	+554	(1,816)	(552)	
Net Profit/ Loss	(1,111)	(534)	+576	(1,825)	(560)	
R&D Cost	1,474	1,841	+367 +24.9%	1,981	1,860	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 plan was revised on Dec. 23, 2021.

FY2021 Results by Business Segment



(JPY million)	FY2020 Actual	FY2021 Actual	YoY Change	FY2021 Plan	FY2021 Dec. 23 Revised Plan		
Total Sales	1,133	2,017	+884 +78.0%	923	2,006		
ddSP business	1,080	889	(190) -17.7%	923	878	Slightly behind the FY sales plan due to weaker than expected sales in Japan.	
ddRD business	53	1,128	+1,074 +2027%	_	1,128	Received milestone payment from Gilead.	
Total Operating Profit/Loss	(1,057)	(531)	+525	(1,811)	(548)		
ddSP business	458	289	(169) -37.0%	207	288	Beat the initial plan thanks to robust sales of high-margin kinase proteins.	
ddRD business	(1,515)	(820)	+695	(2,019)	(836)	Investment in clinical studies continued. Loss were smaller than the initial plan thanks to the milestone payment.	

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 since losses were recorded.

Note 3: FY2021 plan was disclosed on February 12, 2021. FY2021 plan was revised on Dec. 23, 2021.

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





- □ Japan: Decreased 26.2% 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 21.9% YoY
- Contribution from sales to Gilead continued.
- NanoBRET assay service showed strong growth.
- Agent business (cell-based assay service) were weak.
- Europe: Increased 14.3% YoY
- Kinase proteins, profiling service and NanoBRET assay service were robust.
- Other: Increased 22.4% YoY
- Full year sales were robust although sales in September and October were temporarily affected by the delay in the import custom clearance in China.



(JPY million)	As of Dec. 31, 2020	As of Dec. 31, 2021	Change	Reason for changes
Current assets	4,708	5,318	+610	Accounts receivable-trade +1,114 Cash and deposits -481
Cash and deposits	4,299	3,817	-481	
Non-current Assets	127	114	-13	
Total assets	4,835	5,432	+597	
Current liabilities	727	774	+47	
Non-current liabilities	284	342	+58	Long term loans payable +88 Bonds payable -28
Total liabilities	1,011	1,116	+105	
Total net assets	3,824	4,315	+491	Capital stock and capital surplus +996, Retained earnings -534
Total liabilities and net assets	4,835	5,432	+597	

Shareholders' equity ratio	79.0%	79.3%
BPS	308.0 yen	323.5 yen
PBR	3.9x	3.4x
Share price of Carna	1,212 yen	1,102 yen

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



Business Plan for FY2022

Key Milestones for 2022



Achieved

Plan or to be achieved

Business		Key Milestones				
		Milestones for 2021	Achievement in 2021	Next Milestones for 2022		
	AS-0871	Initiate Ph1 MAD part (H2 2021)	Initiated Ph1 MAD part in Dec. 2021	Start partnering activity		
ddRD	AS-1763	□ Initiate Ph1 (H1 2021)	☑Initiated Ph1 in Apr. 2021	Initiate Ph1b (US)		
	AS-0141	Initiate Ph1 (H1 2021)	Initiated Ph1 in June 2021	Initiate Ph1 expansion part		
	Research program	Bring one or more programs in preclinical stage	STING antagonist was advanced to preclinical stage	Bring one or more programs in preclinical stage or license a program.		
ddSP		 Achieve sales target of JPY920 mn Launch new products Expand NanoBRET service Propose project-based service to collaborate with clients, leveraging Carna's drug discovery technology 	 Actual sales were JPY889 mn, 96% of initial target Launched 31 new kinase protein products and 4 new EGFR mutant targets to profiling service Expand NanoBRET service Propose project-based service to collaborate with clients, leveraging Carna's drug discovery technology 	 Expand sales of in-house developed products and services Expand line-up of protein kinase products Increase target kinases to expand profiling service Seek collaboration opportunities to boost Carna's 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)
 Established in-house research capability Established pipeline 	 Out-licensed multiple programs Initiated clinical trials 	 Advance clinical trials of AS-0871, AS-1763, and AS-0141 Receive milestone payments from the out- licensed programs and deliver profits Initiate pre-clinical and clinical studies of new pipelines 	 Receive milestone payments and royalty income from the out-licensed programs and expand profits Earn revenue from new license deals Initiate pre-clinical and clinical studies of new pipelines



<ddRD>

✓ Advance clinical trials of AS-0871, AS-1763, and AS-0141

- ✓ Create next wave of pipeline
- Receive milestone payments and royalty income from out-licensed programs

- 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

Business Plan



(JPY million)	FY2021 Actual	FY2022 Plan	Outlook for 2023 - 2026	
Total Sales	2,017	1,127		
ddSP business	889	900	Maintain stable sales	
ddRD business	1,128	227	Revenue from milestone payments and upfront payments	
Total Operating Loss	(531)	(1,730)		
ddSP business	289	300	Maintain stable profit while investing in product developments	
ddRD business	(820)	(2,031)	Continue to invest in R&D and deliver profits depending on the size of mileston payments and upfront payments	
Ordinary Loss	(522)	(1,744)		
Net Loss	(534)	(1,799)		
(JPY million)	FY2021 Actual	FY2022 Plan	Outlook for 2023 – 2026	
R&D Cost	1,841	2,166	Invest in R&D (JPY1 bn to 2.5 bn) for the future growth.	
Сарех	41	124	Invest in equipment for R&D and IT system (JPY20 mn to 100 mn)	

* Business plan for FY2022 includes an upfront payment from Brickell Biotech but does not include milestone payments and upfront payments related to other license agreements as the timing or the amounts are difficult to predict. Numerical targets for 2023-2026 are not disclosed for the same reason.

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



(JPY mn)	FY2020 Actual	FY20201 Actual	FY2022 Plan
ddSP	1,080	889	900
Kinase Proteins	276	323	372
Assay Development	433	274	184
Profiling & Screening	230	203	222
Cell-based Assay	46	57	96
Cell-based Assay (agent business)	67	12	10
Others	23	17	13

Exchange rate(US\$):	106.77 yen	109.90 yen	110 yen
% of Overseas sales:	74.3%	76.9%	75.5%



Note: Actual foreign exchange rate is average rate of the term





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