

# 2020年3月期 第3四半期決算補足資料

Supplement for the Consolidated Financial Statements  
for the Third Quarter Ended December 31, 2019

タカラバイオ株式会社

Takara Bio Inc.

2020年2月13日

February 13, 2020

## 第3四半期 業績（連結）

Consolidated Quarterly Financial Results for the Third Quarter ended December 31, 2019

(百万円： millions of yen)

	19/03期 第3四半期 FY2019 (3Q)	20/03期 第3四半期 FY2020 (3Q)	前 期 比 Year on year	
			増 減 Change	増 減 率 Ratio
売 上 高 Net sales	25,888	24,558	▲ 1,329	▲ 5.1%
売 上 原 価 Cost of sales	10,947	9,767	▲ 1,180	▲ 10.8%
売 上 総 利 益 Gross profit	14,940	14,791	▲ 149	▲ 1.0%
販 売 費 及 び 一 般 管 理 費 SG&A expenses	11,507	10,911	▲ 595	▲ 5.2%
営 業 利 益 Operating income	3,432	3,879	+446	+13.0%
経 常 利 益 Ordinary income	3,549	3,896	+346	+9.8%
親会社株主に帰属する 四半期純利益 Net income attributable to owners of the parent	2,088	2,708	+619	+29.7%
研 究 開 発 費 R&D expenses	3,164	2,812	▲ 352	▲ 11.1%

## 事業セグメント別売上高（連結）

Consolidated Results by Business Segment

(百万円：millions of yen)

	19/03期 第3四半期 FY2019 (3Q)	20/03期 第3四半期 FY2020 (3Q)	前 期 比 Year on year	
			増 減 Change	増 減 率 Ratio
バイオ産業支援 Bioindustry	21,984	22,388	+403	+1.8%
研究用試薬 Research reagents	17,093	17,750	+657	+3.8%
理化学機器 Scientific instruments	1,907	952	▲ 955	▲ 50.1%
受 託 Contract service	2,647	3,392	+745	+28.2%
そ の 他 Others	335	291	▲ 44	▲ 13.2%
遺伝子医療 Gene therapy	2,304	2,170	▲ 134	▲ 5.8%
そ の 他* Others	1,599	-	▲ 1,599	-
売上高合計 Net sales total	25,888	24,558	▲ 1,329	▲ 5.1%

\* 旧医食品バイオ Former AgriBio

3 2020年3月期第3四半期 決算補足資料：Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## 事業セグメント別業績

Results by Business Segment

(百万円：millions of yen)

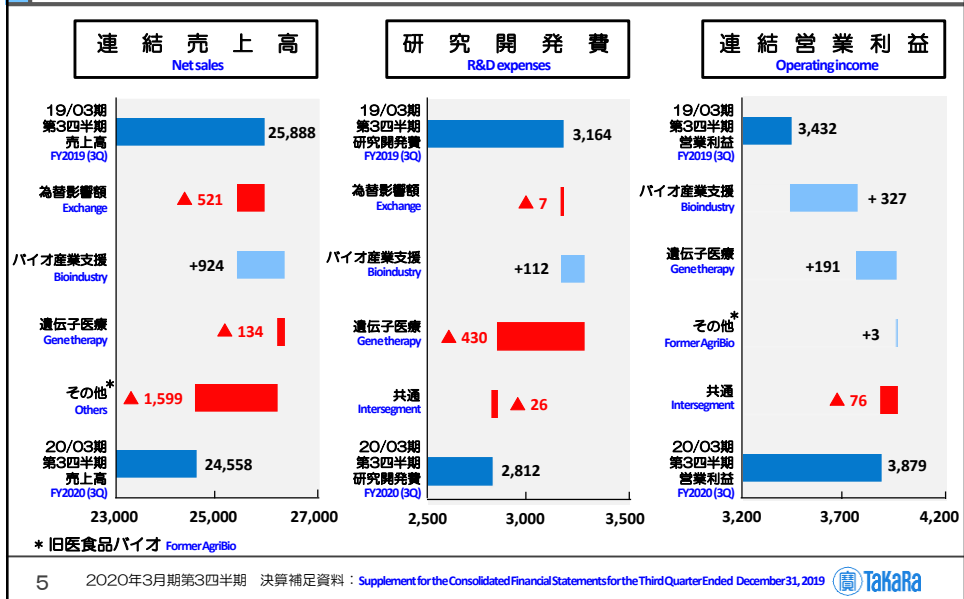
19/03期第3四半期 FY2019 (3Q)	バイオ産業支援 Bioindustry	遺伝子医療 Gene therapy	旧医食品バイオ Former AgriBio	共 通 Intersegment	計 Total
売上高 Net sales	21,984	2,304	1,599	0	25,888
売上総利益 Gross profit	12,691	1,886	361	0	14,940
販売費及び一般管理費 SG&A expenses	8,553	1,000	365	1,587	11,507
内、研究開発費 R&D expenses	2,021	988	0	155	3,164
営業利益 Operating income	4,138	885	▲ 3	▲ 1,587	3,432
20/03期第3四半期 FY2020 (3Q)	バイオ産業支援 Bioindustry	遺伝子医療 Gene therapy	そ の 他 Others	共 通 Intersegment	計 Total
売上高 Net sales	22,388	2,170	0	0	24,558
売上総利益 Gross profit	13,106	1,684	0	0	14,791
販売費及び一般管理費 SG&A expenses	8,640	607	0	1,663	10,911
内、研究開発費 R&D expenses	2,126	557	0	128	2,812
営業利益 Operating income	4,465	1,077	0	▲ 1,663	3,879

4 2020年3月期第3四半期 決算補足資料：Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## 事業セグメント情報<前期比増減>

Analysis of Results by Business Segment

(百万円 : millions of yen)



5

2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019



## 会社別業績

Results by Subsidiary

(百万円 : millions of yen)

	従業員数**	19/03期 第3四半期 FY2019 (3Q)		20/03期 第3四半期 FY2020 (3Q)	
		売上高	営業利益	売上高	営業利益
	Employee	Net sales	Operating income	Net sales	Operating income
タカラバイオ(単体) Takara Bio	523	15,597	1,645	15,289	1,685
Takara Bio Europe SAS(連結)*	73	2,319	126	2,304	73
宝生物工程(大連)* Takara Bio Biotechnology (Dalian)	520	2,881	662	2,455	722
宝日医生物技術(北京)* Takara Biomedical Technology (Beijing)	70	4,444	793	4,638	911
Takara Korea Biomedical *	25	759	123	761	117
DSS Takara India Pvt. Ltd. *	67	239	4	270	22
Takara Bio USA Inc. *	204	8,106	948	8,020	899

\* 決算期が1-12月であり、数値は1-9月実績 : Financial report from Jan. to Sep. \*\* 従業員数は、2019年12月31日現在 : Number of Employees as of Dec. 31, 2019

6

2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019



## 研究用試薬の外部売上高

Sales by Region for Research Reagents

(百万円：millions of yen)

	19/03期 第3四半期 FY2019 (3Q)	20/03期 第3四半期 FY2020 (3Q)	前 期 比 Year on year change	前 期 比 (為替の影響除く) Year on year change (Foreign exchange excluded)
日 本 Japan	3,880	4,171	+290	+290
米 国 USA	6,276	6,389	+113	+141
欧 州 Europe	2,107	2,010	▲ 97	+37
中 国 China	3,929	4,274	+344	+599
韓 国 Korea	669	646	▲ 22	+21
印 度 India	228	258	+29	+43

7 2020年3月期第3四半期 決算補足資料：Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## (通期予想) 通期業績 (連結)

Consolidated Forecast for the Year ending March 31, 2020

(百万円：millions of yen)

	20/03期 通期予想 FY2020 Forecast	前回予想比(11/12) To previous forecast as of Nov. 12		前 期 比 Year on year	
		増 減 Change	増減率 Ratio	増 減 Change	増減率 Ratio
売 上 高 Net sales	34,300	+400	+1.2%	▲ 1,541	▲ 4.3%
売 上 原 価 Cost of sales	13,169	+118	+0.9%	▲ 1,986	▲ 13.1%
売 上 総 利 益 Gross profit	21,130	+281	+1.4%	+445	+2.2%
販 売 費 及 び 一 般 管 理 費 SG&A expenses	14,930	+281	+1.9%	▲ 290	▲ 1.9%
営 業 利 益 Operating income	6,200	-	-	+736	+13.5%
経 常 利 益 Ordinary income	6,300	▲ 50	▲ 0.8%	+634	+11.2%
親会社株主に帰属する 当 期 純 利 益 Net income attributable to owners of the parent	3,700	▲ 550	▲ 12.9%	+42	+1.2%
研 究 開 発 費 R&D expenses	3,934	+20	+0.5%	▲ 402	▲ 9.3%

8 2020年3月期第3四半期 決算補足資料：Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## (通期予想) 事業セグメント別売上高 (連結)

Consolidated FY2020 Forecast by Business Segment

(百万円 : millions of yen)

	20/03期 通期予想 FY2020 Forecast	前回予想比(11/12) To previous forecast as of Nov. 12		前 期 比 Year on year	
		増 減 Change	増減率 Ratio	増 減 Change	増減率 Ratio
バイオ産業支援 Bioindustry	32,018	+487	+1.5%	+442	+1.4%
研究用試薬 Research reagents	24,358	+210	+0.9%	+757	+3.2%
理化学機器 Scientific Instruments	1,165	▲ 89	▲ 7.1%	▲ 1,405	▲ 54.7%
受 託 Contract service	6,082	+334	+5.8%	+1,128	+22.8%
そ の 他 Others	411	+31	+8.2%	▲ 38	▲ 8.5%
遺伝子医療 Gene therapy	2,281	▲ 87	▲ 3.7%	▲ 161	▲ 6.6%
そ の 他* Others	-	-	-	▲ 1,822	-
売上高合計 Net sales total	34,300	+400	+1.2%	▲ 1,541	▲ 4.3%

\* 旧医食品バイオ Former AgriBio

9 2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## (通期予想) 事業セグメント別業績

FY2020 Forecast by Business Segment

(百万円 : millions of yen)

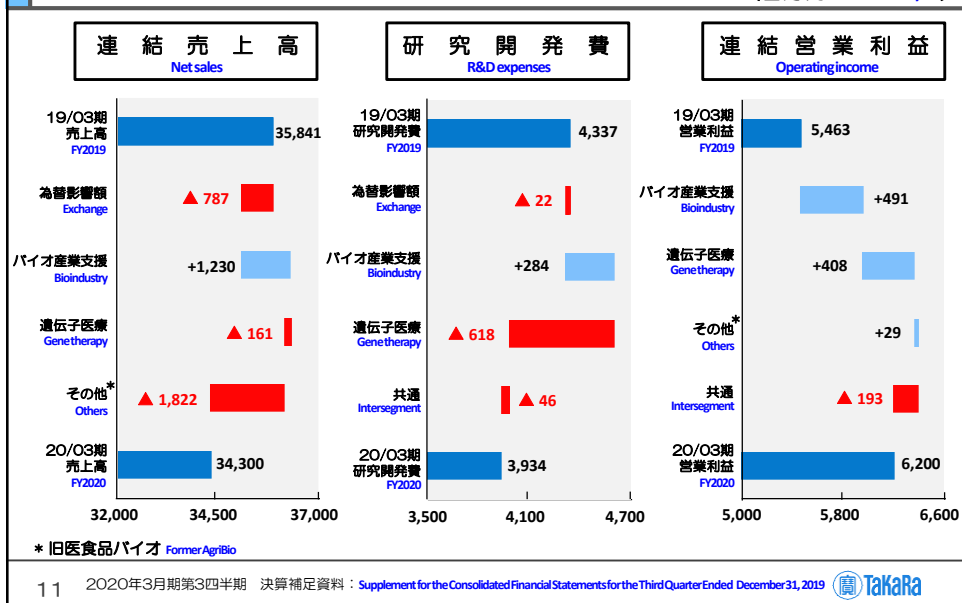
19/03期実績 FY2019 Actual	バイオ産業支援 Bioindustry	遺伝子医療 Gene therapy	旧医食品バイオ Former AgriBio	共 通 Intersegment	計 Total
売 上 高 Net sales	31,575	2,443	1,827	▲ 5	35,841
売 上 総 利 益 Gross profit	18,369	1,915	399	0	20,685
販売費及び一般管理費 SG&A expenses	11,269	1,409	429	2,114	15,221
内、研究開発費 R&D expenses	2,723	1,391	0	222	4,337
営 業 利 益 Operating income	7,100	506	▲ 29	▲ 2,114	5,463
20/03期通期予想 FY2020 Forecast	バイオ産業支援 Bioindustry	遺伝子医療 Gene therapy	そ の 他 Others	共 通 Intersegment	計 Total
売 上 高 Net sales	32,018	2,281	-	0	34,300
売 上 総 利 益 Gross profit	19,372	1,757	-	0	21,130
販売費及び一般管理費 SG&A expenses	11,780	843	-	2,307	14,930
内、研究開発費 R&D expenses	2,985	773	-	175	3,934
営 業 利 益 Operating income	7,592	914	-	▲ 2,307	6,200

10 2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## (通期予想) 事業セグメント情報<前期比増減>

Analysis of FY2020 Forecast by Business Segment

(百万円：millions of yen)



11 2020年3月期第3四半期 決算補足資料：Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019

## (通期予想) 会社別業績

FY2020 Forecast by Subsidiary

(百万円：millions of yen)

	19/03期 通期 FY2019 Actual		20/03期 通期予想 FY2020 Forecast	
	売上高 Net sales	営業利益 Operating income	売上高 Net sales	営業利益 Operating income
タカラバイオ(単体) Takara Bio	21,740	2,312	21,618	2,472
Takara Bio Europe SAS(連結)*	3,362	244	3,090	175
宝生物工程(大連)* Takara Bio Biotechnology (Dalian)	3,805	875	3,477	1,062
宝日医生物技术(北京)* Takara Biomedical Technology (Beijing)	6,121	1,057	6,378	1,224
Takara Korea Biomedical *	1,034	173	1,014	159
DSS Takara India Pvt. Ltd. *	335	5	365	32
Takara Bio USA Inc. *	11,224	1,533	11,217	1,625

\* 決算期が1-12月であり、数値は1-12月実績：Financial report from Jan. to Dec.

12 2020年3月期第3四半期 決算補足資料：Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019

## (通期予想) 研究用試薬の外部売上高

FY2020 Forecast by Region for Research Reagents

(百万円 : millions of yen)

	19/03期 通期 FY2019 Actual	20/03期 通期予想 FY2020 Forecast	前期比 Year on year change	前期比 (為替の影響除く) Year on year change (Foreign exchange excluded)
日本 Japan	5,715	6,000	+285	+285
米国 USA	8,230	8,532	+302	+413
欧州 Europe	2,948	2,703	▲ 245	▲ 60
中国 China	5,470	5,907	+437	+789
韓国 Korea	917	865	▲ 52	+10
印度 India	319	349	+29	+45

13 2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## 為替前提および実績 (海外子会社P/L換算用)

Exchange Rates for the Income Statement of Overseas Subsidiaries

(単位 : 円) (Unit : Yen)	19/03期 第3四半期 FY2019 (3Q)	19/03期 通期 FY2019 (Full year)	20/03期 第3四半期 FY2020 (3Q)	20/03期 通期 FY2020 (Full year)	20/03期 通期 FY2020 (Full year)
	実績 Actual	実績 Actual	実績 Actual	11/12予想 Forecast as of Nov. 12	2/13予想 Forecast as of Feb. 13
米ドル US dollar	109.61	110.45	109.13	108.30	109.03
ユーロ Euro	130.87	130.35	122.62	121.30	122.03
人民元 Yuan	16.84	16.71	15.89	15.60	15.77
100ウォン 100 Won	10.05	10.04	9.40	9.20	9.36
インドルピー Rupee	1.64	1.62	1.56	1.55	1.55
スウェーデン クローナ Sweden Krona	12.79	12.71	11.61	11.50	11.53

14 2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019 

## 遺伝子医療事業開発パイプラインの状況

Pipeline Status for developing in Gene Therapy Business

プロジェクト Projects		対象疾患 Target disease	現状 Status	提携先 Partnerships		
腫瘍溶解性ウイルス Oncolytic Virus	C-REV (TBI-1401)	日本 Japan	膵臓がん Pancreatic cancer	Phase I 進行中 Phase I in progress	大塚製薬(株) Otsuka Pharma	
	一般名: canerpaturev Nonproprietary name	韓国 Korea	全疾患 All indications	検討中 In preparation	東亜ST(株) Dong-AST Co., Ltd	
	商品名: EPLICANA® Trade name	米国 USA	メラノーマ Melanoma	医師主導治験 Phase II 進行中 Phase II Investigator initiated trial in progress	-	
遺伝子改変 T細胞療法 Engineered T cell Therapy	siTCR®	NY-ESO-1 (TBI-1301)	日本 Japan	滑膜肉腫 Synovial sarcoma	Phase I / II 進行中 2020年度申請予定 Phase I / II in progress Planned for application in FY2021	大塚製薬(株) Otsuka Pharma
		カナダ Canada	固形がん Solid cancer	医師主導治験 Phase I b 進行中 Phase I b Investigator initiated trial in progress	-	
	CAR	CD19 (TBI-1501)	日本 Japan	成人ALL* Adult ALL, etc	Phase I / II 進行中 Phase I / II in progress	大塚製薬(株) Otsuka Pharma

\* ALL : 急性リンパ芽球性白血病 Acute lymphoblastic leukemia

15

2020年3月期第3四半期 決算補足資料 : Supplement for the Consolidated Financial Statements for the Third Quarter Ended December 31, 2019



## 将来の見通しに関する注意事項

Forward-Looking Statements

資料中の当社による現在の計画、見通し、戦略、確信などのうち、歴史的事実でないものは、将来の業績に関する見通しであり、これらは現時点において入手可能な情報から得られた当社経営陣の判断に基づくものですが、重大なリスクや不確実性を含んでいる情報から得られた多くの仮定および考えに基づきなされたものであります。実際の業績は、さまざまな要素によりこれら予測とは大きく異なる結果となり得ることをご承知おきください。実際の業績に影響を与える要素には、経済情勢、特に消費動向、為替レートの変動、法律・行政制度の変化、競合会社の価格・製品戦略による圧力、当社の既存製品および新製品の販売力の低下、生産中断、当社の知的所有権に対する侵害、急速な技術革新、重大な訴訟における不利な判決等がありますが、業績に影響を与える要素はこれらに限定されるものではありません。

お問合せ先：広報・IR部

E-mail: bio-ir@takara-bio.co.jp

Statements in this news release, other than those based on historical fact, concerning the current plans, prospects, strategies and expectations of the Company and its Group represent forecasts of future results. While such statements are based on the conclusions of management according to information available at the time of writing, they reflect many assumptions and opinions derived from information that includes major risks and uncertainties. Actual results may vary significantly from these forecasts due to various factors. Factors that could influence actual results include, but are not limited to, economic conditions, especially trends in consumer spending, as well as exchange rate fluctuations, changes in laws and government systems, pressure from competitors' prices and product strategies, decline in selling power of the Company's existing and new products, disruptions to production, violations of our intellectual property rights, rapid advances in technology and unfavorable verdicts in major litigation.

Contact information : PR/IR Department

E-mail: bio-ir@takara-bio.co.jp

