

Summary of Financial Report for the FY ending December 2009

February 12, 2010

Company Name: Carna Biosciences Inc.

Listed Securities Exchange: JQ・NEO

Stock Code: 4572 JQ・NEO

URL <http://www.carnabio.com>

Representative: Kohichiro Yoshino, President

Contact: Norio, Aikawa, Director

TEL (078)302-7039

General meeting of shareholders scheduled: March 24, 2010

Payment of dividends scheduled: -

Release of the financial report scheduled: March 25, 2010

(Rounding down to the nearest million yen)

1. Consolidated results for the FY2009 (January 1, 2009 to December 31, 2009)

(1)Business Results(Consolidated)

(% indicates the increase / decrease from previous fiscal year)

| | Total Revenue | | Operating income | | Ordinary income | | Net income | |
|--------|---------------|------|------------------|---|-----------------|---|-------------|---|
| | million yen | % | million yen | % | million yen | % | million yen | % |
| FY2009 | 687 | 37.5 | -344 | - | -349 | - | -315 | - |
| FY2008 | 499 | - | -296 | - | -346 | - | -503 | - |

| | Net income per share | Net income per share (diluted) | Return on equity | Ordinary income on assets | Operating income on revenue |
|--------|----------------------|--------------------------------|------------------|---------------------------|-----------------------------|
| | yen | yen | % | % | % |
| FY2009 | -5,873.72 | - | -17.6 | -17.0 | -50.1 |
| FY2008 | -9,814.39 | - | -31.2 | -18.8 | -59.3 |

Note: Net income per share of FY2008 and 2009(Diluted): No statement - We had net loss per share even though we had diluted shares.

(2)Financial Condition(Consolidated)

| | Total assets | Net assets | Equity ratio | Net assets per share |
|--------|--------------|-------------|--------------|----------------------|
| | million yen | million yen | % | Yen |
| FY2009 | 2,043 | 1,801 | 88.1 | 30,680.71 |
| FY2008 | 2,070 | 1,789 | 86.4 | 33,587.30 |

(3)Cash Flow(Consolidated)

| | Cash flow from operating activities | Cash flow from investing activities | Cash flow from financing activities | Cash and cash equivalents at the end-of-term |
|--------|--|--|--|--|
| | million yen | million yen | million yen | million yen |
| FY2009 | -419 | 83 | 392 | 1,490 |
| FY2008 | -267 | -313 | 813 | 1,432 |

2. Cash Dividends

| (criterion) | Dividends per share | | | | |
|------------------|--|--|--|-----------|------|
| | End of 1 st quarter period | End of 2 nd quarter period | End of 3 rd quarter period | End of FY | FY |
| | yen | yen | yen | yen | yen |
| FY2008 | - | 0.00 | - | 0.00 | 0.00 |
| FY2009 | - | 0.00 | - | 0.00 | 0.00 |
| FY2010(forecast) | - | 0.00 | - | 0.00 | 0.00 |

3. Consolidated forecast for FY2010 (January 1, 2010 to December 31, 2010)

(%indicates the increase/decrease from previous fiscal year)

| | Total revenue | Operating income | Ordinary income | Net income | Net income per share |
|----------------|---------------|---------------------|-----------------|---------------|-------------------------|
| | million yen % | million yen % | million yen % | million yen % | Yen |
| Interim FY2010 | 385 21.1 | -194 - | -165 - | -173 - | -2,952.92 |
| FY2010 | 840 22.3 | -321 - | -294 - | -304 - | -5,182.64 |

*Disclaimer regarding appropriate use of forecasts and related points of note

1. The revised forecasts are based on the information available on the date of release of these materials and on various assumptions made on the date of the release of these material including uncertain factors and may affect future earnings. Actual operating results may differ significantly from those forecasts as a consequence of various factors.

2. The forecasted net income per share above is calculated based on average numbers of shares outstanding