

Combined Shareholders' Meeting
 7 April 2015, 1:30 pm, headquarters of the Company (Aubagne)

SARTORIUS STEDIM BIOTECH SA.

FIVE-YEAR FINANCIAL RESULTS

**Five-Year Financial Results of the Parent
 Company Sartorius Stedim Biotech S.A.**

| € in K | 2010 | 2011 | 2012 | 2013 | 2014 |
|--|------------|------------|------------|------------|------------|
| Share capital at end of period | | | | | |
| Share capital (capital stock) | 10,378 | 10,386 | 10,395 | 10,396 | 15,359 |
| Number of shares outstanding | 17,013,448 | 17,025,948 | 17,041,306 | 17,042,306 | 15,359,238 |
| Transactions and financial performance | | | | | |
| Sales revenue (excl. VAT) | 65,026 | 71,855 | 81,942 | 1,501 | 1,465 |
| Profit before tax, employee profit sharing plan, amortization, depreciation and provision expenses (and reversals) | 25,884 | 23,617 | 26,218 | 21,180 | 25,967 |
| Income tax | 1,185 | 1,069 | 678 | 292 | 468 |
| Contribution to employee profit-sharing plan | 0 | 0 | 0 | 0 | 0 |
| Net profit | 21,066 | 23,860 | 26,198 | 20,875 | 24,845 |
| Dividends paid or proposal of dividend | 13,783 | 15,327 | 16,877 | 16,878 | 18,412 |
| Earnings per share | | | | | |
| EPS after tax and employee profit-sharing, but before amortization, depreciation and provision expenses | 1.59 | 1.45 | 1.58 | 1.26 | 1.66 |
| EPS after tax and employee profit-sharing, amortization, depreciation and provision expenses | 1.24 | 1.40 | 1.54 | 1.22 | 1.62 |
| Dividend per share | 0.90 | 1.00 | 1.10 | 1.10 | 1.20 |
| Personnel | | | | | |
| Workforce size | 293 | 336 | 388 | 0 | 0 |
| Personnel costs | 11,177 | 11,843 | 14,171 | 0 | 0 |
| Social security costs | 6,007 | 6,574 | 7,969 | 0 | 0 |