

Corporate Calendar 2009 Sartorius AG

March 11, 2009	Annual press conference, Goettingen, Germany
April 23, 2009	Annual Shareholder's Meeting, Goettingen, Germany
April 29, 2009	Three-month report 2009
August 6, 2009	Six-month report 2009
October 31, 2009	Nine-month report 2009
November 9 – 11, 2009	Analysts' conference, German Equity Forum Fall 2009, Frankfurt am Main, Germany