

**Total number of shares and voting rights at the end of each calendar month**

Statement in compliance with Article 222-12-5 of the General Regulation of the *Autorité des Marchés Financiers* (AMF)

	<b>Share capital</b>	<b>Number of shares</b>	<b>Number of voting rights</b>	<b>Treasury shares</b>
<b>January 31, 2007</b>	EUR 3 783 562, 50	12,611,875	12,611,875	8,861
<b>January 31, 2008</b>	EUR 3 783 562, 50	12,611,875	12,611,875	8,861
<b>January 31, 2009</b>	EUR 3 783 562, 50	12,611,875	12,611,875	8,861