

**DECLARATION RELATIVE TO  
THE NUMBER OF SHARES AND VOTING RIGHTS  
MAKING UP THE ISSUED CAPITAL**

Article L 233-8-II of the Code of commerce  
and article 222-12-5 of the general regulations of the AMF

**Designation of issuer**

STEDIM SA  
BP 1051  
Z I des Paluds  
Avenue de Jouques  
13781 Aubagne cedex

<b>Deadline date for information</b>	<b>Total number of shares making up the issued capital</b>	<b>Total number of voting rights</b>
June 28 <sup>th</sup> , 2007	7 136 825	Gross total of voting rights : 10 618 050
		Net total of voting rights : 10 618 050

\* Net total = total number of voting rights attached to shares – private share voting right