

**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

Deadline date for information	Total number of shares making up the issued capital	Total number of voting rights
December 31 st , 2006	7 057 555	Gross total of voting rights : 10 539 130
		Net total of voting rights : 10 539 130

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