

Independent Auditors' Fees

Principal Independent Auditors

Ernst and Young Audit

408, avenue du Prado – BP 116 – 13267 Marseille Cedex 08 – France

Represented by Anis Nassif.

First commissioned by the Combined General Shareholders' Meeting on June 28, 1985. Date commission expires: 2015 Annual General Shareholders' Meeting to approve the 2014 financial statements.

Member of Compagnie régionale de Versailles.

Deloitte et Associés

10, Place de la Joliette – Les Docks – Atrium 10.4 – BP 64529 - 13567 Marseille Cedex 02 – France

Represented by Christophe Perrau.

First commissioned by the Annual General Shareholders' Meeting on May 19, 2006. Date commission expires: 2018 Annual General Shareholders' Meeting to approve the 2017 financial statements.

Member of Compagnie régionale de Versailles.

Independent Auditors' Fees

€ in K	Ernst & Young				Deloitte			
	2012		2011		2012		2011	
Audit								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	129	94.9%	115	92.0%	129	20.9%	115	17.3%
Subsidiaries	5	3.7%	5	4.0%	391	63.5%	332	49.8%
Services directly related to audit services								
Parent company								
Subsidiaries								
Subtotal	134	98.5%	120	96.0%	520	84.4%	447	67.1%
Other services								
Legal, tax, corporate					18	2.9%	50	7.5%
Information technology, other	2	1.5%	5	4.0%	78	12.7%	169	25.4%
Subtotal	2	1.5%	5	4.0%	96	15.6%	219	32.9%
Total	136	100%	125	100%	616	100%	666	100%

Substitute Independent Auditors

Auditex

Tour Ernst & Young – Faubourg de l'Arche – 92037 Paris-La Défense Cedex

First commissioned by the Annual General Shareholders' Meeting on April 21, 2009. Date commission expires: 2015 Annual General Shareholders' Meeting to approve the 2014 financial statements.

Member of Compagnie régionale de Versailles.

BEAS

7/9, Villa Houssay – 92200 Neuilly sur Seine – France
Represented by Alain Pons.

Commissioned by the Annual General Shareholders' Meeting on May 19, 2006. Date commission expires: 2018 Annual General Shareholders' Meeting to approve the 2017 financial statements.

Member of Compagnie régionale de Versailles.

2012		Other 2011		2012		Total 2011	
				258	22.9%	230	19.8%
273	72.4%	296	80.4%	669	59.3%	633	54.6%
273	72.4%	296	80.4%	927	82.1%	863	74.5%
58	15.4%	45	12.2%	76	6.7%	95	8.2%
46	12.2%	27		126	11.2%	201	17.3%
104	27.6%	72	19.6%	202	17.9%	296	25.5%
377	100%	368	100%	1,129	100%	1,159	100%