

Independent Auditors' Fees

Principal Independent Auditors

Ernst and Young Audit

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

Represented by Xavier Senent.

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			
	2013		2012		2013		2012	
Audit								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	80	34.9%	129	94.9%	110	23.2%	129	20.9%
Subsidiaries	50	21.8%	5	3.7%	340	71.6%	391	63.5%
Services directly related to audit services								
Parent company								
Subsidiaries								
Subtotal	130	56.8%	134	98.5%	450	94.7%	520	84.4%
Other services								
Legal, tax, corporate	0				17	3.6%	18	2.9%
Information technology, other	99	43.2%	2	1.5%	8	1.7%	78	12.7%
Subtotal	99	43.2%	2	1.5%	25	5.3%	96	15.6%
Total	229	100%	136	100%	475	100%	616	100%

Substitute Independent Auditors

Auditex

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

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.

2013		Other 2012		2013		Total 2012	
				190	18.8%	258	22.9%
277	89.9%	273	72.4%	667	65.9%	669	59.3%
277	89.9%	273	72.4%	857	84.7%	927	82.1%
25	8.1%	58	15.4%	42	4.2%	76	6.7%
6	1.9%	46	12.2%	113	11.2%	126	11.2%
31	10.1%	104	27.6%	155	15.3%	202	17.9%
308	100%	377	100%	1,012	100%	1,129	100%