

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			
	2014		2013		2014		2013	
Audit								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	55	45.8%	80	34.9%	87	17.1%	110	23.2%
Subsidiaries	61	50.8%	50	21.8%	398	78.3%	340	71.6%
Services directly related to audit services								
Parent company								
Subsidiaries								
Subtotal	116	96.7%	130	56.8%	485	95.5%	450	94.7%
Other services								
Legal, tax, corporate	4	3.3%			23	4.5%	17	3.6%
Information technology, other	0	0.0%	99	43.2%	0	0.0%	8	1.7%
Subtotal	4	3.3%	99	43.2%	23	4.5%	25	5.3%
Total	120	100.0%	229	100%	508	100.0%	475	100%

Substitute Independent Auditors

Auditex

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

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.

2014		Other 2013		2014		Total 2013	
				142	15.1%	190	18.8%
255	81.7%	277	89.9%	714	76.0%	667	65.9%
255	81.7%	277	89.9%	856	91.1%	857	84.7%
57	18.3%	25	8.1%	84	8.9%	42	4.2%
0	0.0%	6	1.9%	0	0.0%	113	11.2%
57	18.3%	31	10.1%	84	8.9%	155	15.3%
312	100.0%	308	100%	940	100.0%	1,012	100%