

Independent Auditors' Fees

Principal Independent Auditors

KPMG S.A.

480, avenue du Prado – CS 90021 – 13272 Marseille Cedex 08 – France

Represented by John Evans.

First commissioned by the Combined General Shareholders' Meeting on 7 April 2015.

Date commission expires: 2021 Annual General Shareholders' Meeting to approve the 2020 financial statements.

Member of the 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 19 May 2006.

Date commission expires: 2018 Annual General Shareholders' Meeting to approve the 2017 financial statements.

Member of the Compagnie régionale de Versailles.

Independent Auditors' Fees

€ in K	KPMG				Deloitte			
	2016		2015		2016		2015	
Audit								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	83	11.6%	56	7.8%	73	30.9%	75	27.0%
Subsidiaries	627	87.3%	510.8	71.1%	137	57.8%	161.4	58.1%
Services directly related to audit services								
Parent company								
Subsidiaries								
Subtotal	710	98.9%	566.8	78.9%	210	88.8%	236.4	85.0%
Other services								
Legal, tax, corporate	0	0.0%	1.8	0.3%	2	0.6%	10.5	3.8%
Information technology, other	8	1.1%	0	0.0%	25	10.6%	31.1	11.2%
Subtotal	8	1.1%	1.8	0.3%	27	11.2%	41.6	15.0%
Total	718	100.0%	568.6	79%	236	100.0%	278	100%

Substitute Independent Auditors

SALUSTRO REYDEL

3, cours du Triangle – Immeuble Le Palatin -92939
Paris-La Défense Cedex – France

Represented by Jean-Claude Reydel.

First commissioned by the Annual General Shareholders' Meeting on 7 April 2015.

Date commission expires: 2021 Annual General Shareholders' Meeting to approve the 2020 financial statements.

Member of the 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 19 May 2006.

Date commission expires: 2018 Annual General Shareholders' Meeting to approve the 2017 financial statements.

Member of the Compagnie régionale de Versailles.

		Ernst & Young		Other				Total			
		2016	2015	2016		2015		2016		2015	
		13	52.0%					156	14.1%	144	13.7%
		12	48.0%	118	76.3%	131.8	74.9%	881	79.5%	816	77.9%
		25	100.0%	118	76.3%	131.8	74.9%	1,037	93.6%	960	91.6%
		0		18	11.4%	17.7	10.1%	19	1.7%	30	2.9%
		0	0.0%	19	12.2%	26.5	15.1%	52	4.7%	57.6	5.5%
		0	0.0%	37	23.7%	44.2	25.1%	71	6.4%	87.6	8.4%
		25	100%	155	100.0%	176	100%	1,109	100.0%	1,048	100%