

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			
	2017		2016		2017		2016	
Audit								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	76	8.7%	83	9.5%	58	35.2%	73	30.9%
Subsidiaries	753	85.9%	627	71.6%	97	58.8%	137	57.9%
Services directly related to audit services								
Parent company								
Subsidiaries								
Subtotal	829	94.6%	710	81.0%	155	93.9%	210	88.8%
Other services								
Legal, tax, corporate	47	5.4%	0	0.0%	0	0.0%	2	0.6%
Information technology, other	0	0.0%	8	0.9%	10	6.1%	25	10.6%
Subtotal	47	5.4%	8	0.9%	10	6.1%	27	11.2%
Total	876	100.0%	718	82%	165	100.0%	236	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					
2017		2016		2017		2016		2017		2016	
								134	10.4%	156	14.1%
				88	35.6%	118	76.3%	938	72.7%	881	79.5%
		0		88	35.6%	118	76.3%	1,072	83.1%	1,037	93.6%
1	100.0%	0		140	56.4%	18	11.4%	188	14.6%	19	1.7%
		0		20	8.0%	19	12.2%	30	2.3%	52	4.7%
1		0		160	64.4%	36.6	23.7%	218	16.9%	71	6.4%
1	100.0%	0		248	100.0%	154.6	100%	1,291	100.0%	1,109	100%