

# 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		2015		2014	
		2015	2014					
<b>Audit</b>								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	56	9.8%		75	27.0%	87	17.1%	
Subsidiaries	511	89.8%		161	58.1%	398	78.3%	
Services directly related to audit services								
Parent company								
Subsidiaries								
<b>Subtotal</b>	<b>567</b>	<b>99.7%</b>		<b>236</b>	<b>85.0%</b>	<b>485</b>	<b>95.5%</b>	
<b>Other services</b>								
Legal, tax, corporate	2	0.3%		11	3.8%	23	4.5%	
Information technology, other	0	0.0%		31	11.2%	0	0.0%	
<b>Subtotal</b>	<b>2</b>	<b>0.3%</b>		<b>42</b>	<b>15.0%</b>	<b>23</b>	<b>4.5%</b>	
<b>Total</b>	<b>569</b>	<b>100.0%</b>		<b>278</b>	<b>100.0%</b>	<b>508</b>	<b>100%</b>	

