

# 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 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.

### Deloitte et Associés

10, Place de la Joliette – Les Docks – Atrium 10.4 – BP 64529 - 13567 Marseille Cedex 02 – France

Represented by Vincent Gros.

First commissioned by the Annual General Shareholders' Meeting on 19 May 2006.

Date commission expires: 2024 Annual General Shareholders' Meeting to approve the 2023 financial statements.

Member of the Compagnie régionale de Versailles.

## Independent Auditors' Fees

€ in K	KPMG				Deloitte			
	2018		2017		2018		2017	
<b>Audit</b>								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	61	7.4%	76	9.3%	54	37.2%	58	35.2%
Subsidiaries	754	92.0%	753	91.9%	91	62.8%	97	58.8%
Services directly related to audit services								
Parent company								
Subsidiaries								
<b>Subtotal</b>	<b>815</b>	<b>99.5%</b>	<b>829</b>	<b>101.1%</b>	<b>145</b>	<b>100.0%</b>	<b>155</b>	<b>93.9%</b>
<b>Other services</b>								
Legal, tax, corporate	5	0.5%	47	5.8%	0	0.0%	0	0.0%
Information technology, other	0	0.0%	0	0.0%	0	0.0%	10	6.1%
<b>Subtotal</b>	<b>5</b>	<b>0.5%</b>	<b>47</b>	<b>5.8%</b>	<b>0</b>	<b>0.0%</b>	<b>10</b>	<b>6.1%</b>
<b>Total</b>	<b>820</b>	<b>100.0%</b>	<b>876</b>	<b>107%</b>	<b>145</b>	<b>100.0%</b>	<b>165</b>	<b>100%</b>

## 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.

Ernst & Young		Other				Total			
2018	2017	2018		2017		2018		2017	
						115	10.2%	134	10.4%
		60	37.8%	88	35.7%	905	80.6%	938	72.7%
		<b>0</b>	<b>37.8%</b>	<b>88</b>	<b>35.5%</b>	<b>1,020</b>	<b>90.8%</b>	<b>1,072</b>	<b>83.1%</b>
		1	32.4%	140	56.5%	56	5.0%	188	14.6%
		0	29.7%	20	8.0%	47	4.2%	30	2.3%
		<b>1</b>	<b>62.2%</b>	<b>160</b>	<b>64.5%</b>	<b>103</b>	<b>9.2%</b>	<b>218</b>	<b>16.9%</b>
		<b>1</b>	<b>100.0%</b>	<b>248</b>	<b>100%</b>	<b>1,123</b>	<b>100.0%</b>	<b>1,291</b>	<b>100%</b>