

# Independent Auditors' Fees

## Principal Independent Auditors

### Ernst and Young Audit

400, Promenade des Anglais – BP 33124 – 06203 Nice Cedex 3 – France

Represented by Anis Nassif.

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 Vincent Gros.

First commissioned by the Annual General Shareholders' Meeting on May 19, 2006. Date commission expires: 2012 Annual General Shareholders' Meeting to approve the 2011 financial statements.

## Independent Auditors' Fees

€ in K	Ernst & Young				Deloitte			
	2009		2008		2009		2008	
<b>Audit</b>								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	105	75.0%	183	96.8%	110	21.6%	163	29.5%
Subsidiaries	31	22.1%	6	3.2%	370	72.5%	308.1	55.7%
Services directly related to audit services								
Parent company								
Subsidiaries				0.0%		0.0%		0.0%
<b>Subtotal</b>	<b>136</b>	<b>97.1%</b>	<b>189</b>	<b>100.0%</b>	<b>480</b>	<b>94.1%</b>	<b>471.1</b>	<b>85.2%</b>
<b>Other services</b>								
Legal, tax, corporate	2				30		82	
Information technology, other	2							
<b>Subtotal</b>	<b>4</b>	<b>3%</b>	<b>0</b>	<b>0%</b>	<b>30</b>	<b>6%</b>	<b>82</b>	<b>15%</b>
<b>Total</b>	<b>140</b>	<b>100%</b>	<b>189</b>	<b>100%</b>	<b>510</b>	<b>100%</b>	<b>553.1</b>	<b>100%</b>

## Substitute Independent Auditors

### Auditex

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

Member of Compagnie régionale de Versailles.

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.

### 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: 2012 Annual General Shareholders' Meeting to approve the 2011 financial statements.

	2009		Other 2008		2009		Total 2008
	0.0%		0.0%	215	19.1%	346	35.5%
249	52.1%	232.8	100.0%	650	57.6%	546.9	56.1%
				0		0	
	0.0%		0.0%	0	0.0%	0	0.0%
249	52.1%	232.8	100.0%	865	76.7%	892.9	91.6%
229				261		82	
				2		0	
229	48%	0	0%	263	23%	82	8%
478	100%	232.8	100%	1,128	100%	974.9	100%