

## Independent Auditors' Fees

### Principal Independent Auditors

#### Ernst and Young Audit

408, avenue du Prado – BP 116 – 13267 Marseille Cedex 08 – 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	2010		Ernst & Young 2009		2010		Deloitte 2009	
<b>Audit</b>								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	108	78.8%	105	75.0%	113	19.8%	110	21.6%
Subsidiaries	9	6.6%	31	22.1%	315	55.1%	370	72.5%
Services directly related to audit services								
Parent company								
Subsidiaries								
<b>Subtotal</b>	<b>117</b>	<b>85.4%</b>	<b>136</b>	<b>97.1%</b>	<b>428</b>	<b>74.8%</b>	<b>480</b>	<b>94.1%</b>
<b>Other services</b>								
Legal, tax, corporate	0		2	1.4%	71	12.4%	30	5.9%
Information technology, other	20	14.6%	2	1.4%	73	12.8%		
<b>Subtotal</b>	<b>20</b>	<b>14.6%</b>	<b>4</b>	<b>2.9%</b>	<b>144</b>	<b>25.2%</b>	<b>30</b>	<b>5.9%</b>
<b>Total</b>	<b>137</b>	<b>100%</b>	<b>140</b>	<b>100%</b>	<b>572</b>	<b>100%</b>	<b>510</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.

	2010		Other 2009		2010		Total 2009
	0.0%		0.0%	221	20.1%	215	19.1%
260	66.5%	249	52.1%	584	53.1%	650	57.6%
<b>260</b>	<b>66.5%</b>	<b>249</b>	<b>52.1%</b>	<b>805</b>	<b>73.2%</b>	<b>865</b>	<b>76.7%</b>
131	33.5%	229	47.9%	202	18.4%	261	23.1%
				93	8.5%	2	0.2%
<b>131</b>	<b>33.5%</b>	<b>229</b>	<b>47.9%</b>	<b>295</b>	<b>26.8%</b>	<b>263</b>	<b>23.3%</b>
<b>391</b>	<b>100%</b>	<b>478</b>	<b>100%</b>	<b>1,100</b>	<b>100%</b>	<b>1,128</b>	<b>100%</b>