

## 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.

Proposition to renew Deloitte et Associés at the 17 April 2012 Annual General Shareholders' meeting.

### Independent Auditors' Fees

€ in K	Ernst Et Young				Deloitte			
	2011		2010		2011		2010	
<b>Audit</b>								
Independent audit, certification, parent company & consolidated financial statements								
Parent company	115	92.0%	108	78.8%	115	17.3%	113	19.8%
Subsidiaries	5	4.0%	9	6.6%	332	49.8%	315	55.1%
Services directly related to audit services								
Parent company								
Subsidiaries								
<b>Subtotal</b>	<b>120</b>	<b>96.0%</b>	<b>117</b>	<b>85.4%</b>	<b>447</b>	<b>67.1%</b>	<b>428</b>	<b>74.8%</b>
<b>Other services</b>								
Legal, tax, corporate					50	7.5%	71	12.4%
Information technology, other	5	4.0%	20	14.6%	169	25.4%	73	12.8%
<b>Subtotal</b>	<b>5</b>	<b>4.0%</b>	<b>20</b>	<b>14.6%</b>	<b>219</b>	<b>32.9%</b>	<b>144</b>	<b>25.2%</b>
<b>Total</b>	<b>125</b>	<b>100%</b>	<b>137</b>	<b>100%</b>	<b>666</b>	<b>100%</b>	<b>572</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.

Proposition to renew BEAS at the 17 April 2012 Annual General Shareholders' meeting.

2011		Other 2010		2011		Total 2010	
				230	19.8%	221	20.1%
296	80.4%	260	66.5%	633	54.6%	584	53.1%
<b>296</b>	<b>80.4%</b>	<b>260</b>	<b>66.5%</b>	<b>863</b>	<b>74.5%</b>	<b>805</b>	<b>73.2%</b>
45	12.2%	131	33.5%	95	8.2%	202	18.4%
27	7.3%			201	17.3%	93	8.5%
<b>72</b>	<b>19.6%</b>	<b>131</b>	<b>33.5%</b>	<b>296</b>	<b>25.5%</b>	<b>295</b>	<b>26.8%</b>
<b>368</b>	<b>100%</b>	<b>391</b>	<b>100%</b>	<b>1,159</b>	<b>100%</b>	<b>1,100</b>	<b>100%</b>