

# Independent Auditors' Fees

## Principal Independent Auditors

### Ernst and Young Audit

408, Avenue du Prado – BP 116 –  
13267 Marseille Cedex 08 – France  
Represented by Jérôme Magnan.  
First commissioned by the Combined General  
Shareholders' Meeting on June 28, 1985.  
Date commission expires: 2009 Annual General  
Shareholders' Meeting to approve the 2008  
financial statements.  
Member of Compagnie régionale de Versailles.

### Deloitte et Associés

10, place de la Joliette – Les Docks – Atrium 10.4 –  
13002 Marseille – 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					
	2008		2007		2006		2008		2007		2006	
<b>Audit</b>												
Independent audit, certification, parent company & consolidated financial statements												
Parent company	<b>183</b>	<b>96.8%</b>	70	17.0%	104	87.4%	<b>163</b>	<b>29.5%</b>	64	10.7%	51	94.4%
Subsidiaries	<b>6</b>	<b>3.2%</b>	5	1.2%		0.0%	<b>308.1</b>	<b>55.7%</b>	320	53.4%		0.0%
Services directly related to audit services												
Parent company												
Subsidiaries			337	81.8%	15	12.6%	<b>0.0%</b>		94	15.7%	3	5.6%
Subtotal	<b>189</b>	<b>100.0%</b>	412	100.0%	119	100.0%	<b>471.1</b>	<b>85.2%</b>	478	79.8%	54	100.0%
<b>Other services</b>												
Legal, tax, corporate												
							<b>82</b>		121			
Information technology, other												
Subtotal	<b>0</b>	<b>0%</b>	0	0%	0	0%	<b>82</b>	<b>15%</b>	121	20%	0	0%
<b>Total</b>	<b>189</b>	<b>100%</b>	<b>412</b>	<b>100%</b>	<b>119</b>	<b>100%</b>	<b>553.1</b>	<b>100%</b>	<b>599</b>	<b>100%</b>	<b>54</b>	<b>100%</b>

## Substitute Independent Auditors

### Patrick Gounelle

Tour Ernst & Young – Faubourg de l'Arche –  
92037 Paris La Défense Cedex – France  
Member of Compagnie régionale de Versailles.  
First commissioned by the Annual General  
Shareholders' Meeting on March 22, 1991.  
Date commission expires: 2009 Annual General  
Shareholders' Meeting to approve the 2008  
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.

				Other						Total	
2008		2007		2006		2008		2007		2006	
	0.0%		0.0%		0.0%	<b>346</b>	<b>35.5%</b>	134	10.4%	155	63.5%
<b>232.8</b>	<b>100.0%</b>	208	73.5%	71	100.0%	<b>546.9</b>	<b>56.1%</b>	533	41.2%	71	29.1%
						<b>0</b>		<b>0</b>		<b>0</b>	
	0.0%		0.0%		0.0%	<b>0</b>	<b>0.0%</b>	431	33.3%	18	7.4%
<b>232.8</b>	<b>100.0%</b>	208	73.5%	71	100.0%	<b>892.9</b>	<b>91.6%</b>	1098	84.9%	244	100.0%
		75				<b>82</b>		196		<b>0</b>	
						<b>0</b>		<b>0</b>		<b>0</b>	
<b>0</b>	<b>0%</b>	75	27%	0	0%	<b>82</b>	<b>8%</b>	196	15%	<b>0</b>	<b>0%</b>
<b>232.8</b>	<b>100%</b>	<b>283</b>	<b>100%</b>	<b>71</b>	<b>100%</b>	<b>974.9</b>	<b>100%</b>	<b>1294</b>	<b>100%</b>	<b>244</b>	<b>100%</b>