

Fees from the Group's Statutory auditors and members of their networks (2007-2008)

	Ernst & Young				Groupe Y Audit			
	Amount (Eur excl. tax)		%		Amount (Eur excl. tax)		%	
	2008	2007	2008	2007	2008	2007	2008	2007
Audit								
Audit, duties and certification of annual and consolidated financial statements:								
. Cerep SA	83,753	60,298	100%	63%	58,038	47,438	100%	100%
. Fully consolidated subsidiaries	-	25,286	-	26%	-	-	-	-
Other assignments and services directly related to the Statutory auditors:								
. Cerep SA	-	10,750	-	11%	-	-	-	-
. Fully consolidated subsidiaries	-	-	-	-	-	-	-	-
Sub-total	83,753	96,334	100%	100%	58,038	47,438	100%	100%
Other services provided by the networks to fully consolidated subsidiaries								
Legal, tax, labor	-	-	-	-	-	-	-	-
Other	-	-	-	-	-	-	-	-
Sub-total	-	-	-	-	-	-	-	-
Total	83,753	96,334	100%	100%	58,038	47,438	100%	100%