

Report Parameters

Our sustainability progress is calculated on a calendar-year basis, with information in this report covering January 1, 2010, to December 31, 2011.

Reporting on Promega Corporate Responsibility progress is completed on a calendar year basis with information in this report covering January 1, 2010 to December 31, 2010. This is the third Promega report in this area following the initial report released in July of 2009. This process of reporting will continue on an annual basis. Corporate Responsibility reporting attempts to focus on the environmental and social impacts of Promega operations worldwide using the framework established by the Global Reporting Initiative's G3 Guidelines and the principles of the United Nations Global Compact.

By establishing a more timely process for gathering information worldwide, Promega has seen improved accuracy and transparency in its key indicators for environmental and social impacts. Increases in the scope, materiality, and comprehensiveness have been experienced but Promega recognizes that there is still significant room for growth. Information has been gathered from all Promega branches and subsidiaries worldwide, including Promega Korea and Promega Biosystems in Sunnyvale, California, both added in 2010. In some instances, additional or more accurate information has been gathered resulting in variations from reported indicators in previous reports. Estimations for

previous years' indicators has been made where information was unavailable using revenue as a factor.

Areas that have not been measured in this report due to lack of current information are:

- Packaging: By implementing a new system for gathering and tracking data, Promega will have a better understanding of gross packaging material usage by type and the impacts from these activities.
- Staff Commute
- Effluents to Water
- Supply Chain Analysis

Carbon footprint calculations have been made using the emission factors provided by the [World Resources Institute Greenhouse Gas Protocol](#) on energy and business travel. The reported emissions from distribution were calculated with the conversion factors provided by [Defra's 2010 Greenhouse Gas Conversion Factors](#). Lastly, the [Environmental Defense Fund's Paper Calculator](#) has been used for calculating the life cycle impacts due to paper usage. Current and previous years' carbon footprints have been calculated using the most updated information and current emission factors listed above.