



Promega Corporation
2800 Woods Hollow Road
Madison, WI 53711-5399
United States

tel: 608.274.4330 • fax: 608.277.2516 • www.promega.com

Promega REACH Policy

The Registration, Evaluation, Authorization and Restriction of Chemicals (REACH) is a European Community Regulation on chemicals and their safe use (Regulation (EC) No 1907/2006).

Promega is fully aware of the requirements of REACH and the dates by which various activities need to be implemented. It is our intention to assist our customers and suppliers to be in full compliance with the legal requirements of REACH.

Promega supports REACH in improving the protection of human health and the environment through the better and earlier identification of intrinsic properties of chemical substances. Promega meets all applicable REACH requirements and is committed to provide our customers with information about substances in its products as required.

A handwritten signature in black ink, appearing to read 'Ron Wheeler', is written over a horizontal line.

Ron Wheeler
Senior Director, Quality Assurance and Regulatory Affairs