



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

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

Promega Conflict Minerals Policy

Promega is committed to the responsible sourcing of material, products and components necessary to the functioning of the products that Promega manufactures or contracts to manufacture. Accordingly, Promega will collaborate with its suppliers to work towards a conflict-free mineral supply chain.

Promega expects its suppliers to adopt policies and management systems consistent with the Dodd-Frank Act (or similar regulations) to supply materials, products, and components that are conflict-free. Promega also requires their suppliers to adopt similar policies, systems and socially responsible sourcing practices. Therefore, we expect our suppliers to respond to our inquiries and provide declarations supporting their commitment to becoming conflict-free and documenting countries of origin for conflict mineral that they purchase.

Customers

Promega customers who are required to comply with the Dodd-Frank Act or similar regulations may request information regarding conflict mineral content in the products sold by Promega. Promega understands the critical need to respond to these requests in a timely manner.

Thank you for your cooperation and assistance in this matter.

A handwritten signature in black ink, appearing to read "Ron Wheeler".

Ron Wheeler
Senior Director, Quality Assurance and Regulatory Affairs