



Luminex Corporation Conflict Minerals Policy Statement

Luminex Corporation (Luminex) is committed to sound ethical practices and full compliance with all applicable laws and regulations. We strive to act in an environmentally and socially responsible manner, in compliance with applicable laws, while meeting the needs of our customers and shareholders.

In July 2010, the U.S. Congress passed Section 1502 of the Dodd-Frank Financial Reform and Consumer Protection Act (Dodd-Frank Act) and in 2012 the U.S. Securities and Exchange Commission (SEC) issued final rules implementing the "conflict mineral" disclosure provisions of Section 1502 of the Dodd-Frank Act. Section 1502 applies to manufactured goods containing the following minerals: tantalum, tin, gold, or tungsten, hereafter collectively referred to as "conflict minerals". Under these rules, publicly-traded companies, like Luminex, are required to make reasonable country of origin inquiries and, depending on the circumstances, undertake further due diligence to determine the source and chain of custody of "conflict minerals" that are contained in their products. Further, these companies must annually disclose and submit Conflict Mineral Reports to the SEC if their products contain "conflict minerals" originating from the Democratic Republic of Congo (DRC) or adjoining countries, where armed groups controlling local mines are using mining proceeds to fund regional conflicts.

Luminex is collaborating with its suppliers, other companies, service providers and consultants to track the origin of metals used in our products to ensure compliance with SEC rules. The international supply chain for these minerals is complex and tracking them is challenging. We are using our reasonable best efforts to ensure that we do not directly or indirectly support violence and human rights abuses in the DRC region.

Although Luminex does not directly purchase any conflict minerals from suppliers, Luminex purchases component parts from its suppliers that Luminex endeavors to ensure do not contain conflict minerals. Suppliers are expected to ensure that parts and products supplied to Luminex Corporation are DRC conflict-free and do not contain metals derived from "conflict minerals" or their derivatives such that they directly or indirectly finance or benefit armed groups through mining or mineral trading in the DRC region. Suppliers are expected to establish policies, due diligence frameworks, and management systems, consistent with the Organisation for Economic Co-operation and Development's Due Diligence Guidance for Responsible Supply Chains of Minerals from Conflict-Affected and High-Risk Areas, that are designed to accomplish this goal. We also request them to submit their sourcing information using the Electronics Industry Citizenship Coalition-Global e-Sustainability Initiative's Conflict Minerals Reporting Template as a standard.

Luminex will evaluate its relationships with its suppliers on an ongoing basis to ensure continued compliance with this policy. Luminex fully expects our suppliers to cooperate with us and to provide information to support these efforts. Suppliers that do not reasonably comply with Luminex's Conflict Mineral Policy will be reviewed by Luminex's supply chain management to assess whether Luminex will conduct business with those suppliers in the future.