



January 30, 2004

The Thermo Electron product Orion number listed below does not require a Material Safety Data Sheet per OSHA 29 CFR 1910.1200. It either does not contain any hazardous ingredients or it is composed of a mixture of chemicals and water with less than 1% hazardous chemical and less than 0.1% of a chemical that would be considered to be a carcinogen.

Orion Number-----700012

Sincerely,

A handwritten signature in black ink, appearing to read "R. Geis", written in a cursive style.

Ron Geis
Director of Continuous Improvement