

**G-Biosciences/Genotech:** St. Louis, MO, USA; [www.GBiosciences.com](http://www.GBiosciences.com) email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

**Product:** Assay Cuvettes (1ml)      **Catalog #:** 786-009 & 786-009A      [LU 01.15.10/JK-IA]

<u>Components</u>	<u>Catalog #</u>	<u>Quantity</u>
Assay Cuvettes (1ml)	786-009	5 x100
Assay Cuvettes (1ml)	786-009A	100

### **Material Safety Data Sheet**

Assay Cuvettes (1ml)

**Catalog #:** a component of product Cat # 786-009 &786-009A indicated above

**Composition:** Plastic

The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. Genotech shall not be held liable for any damage resulting from handling or from contact with the above product. See reverse side of invoice or packing slip for additional terms and conditions of sale.

***G-Biosciences/Genotech St. Louis, MO, USA***

**Web:** [www.GBiosciences.com](http://www.GBiosciences.com), email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)