

G-Biosciences/Genotech: St. Louis, MO, USA; www.GBiosciences.com email: technical@GBiosciences.com

Product: Assay Tubes (2ml) **Catalog #:** 786-008

[LU 01.15.10/JK-IA]

Components

Quantity

Assay Tubes (2ml)

500

Material Safety Data Sheet

Assay Tubes (2ml)

Catalog #: a component of product Cat # 786-008 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, email: technical@GBiosciences.com