

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

Product: Centrifuge Tube Adaptor

Catalog #: 786-145

[LU 5.19.09/JK]

Components

Quantity

Centrifuge Tube Adaptor

2

Material Safety Data Sheet

Composition/Information on Ingredients

Plastic material, believed to be non-hazardous and non-toxic.

Other Information

The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. G-Biosciences 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