

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

Product: Human Tumor Esophagus Single Tissue Blot **Catalog #:** TB91 [LU 4.27.09/MM]

Components

Quantity

Human Tumor Esophagus Single Tissue Blot

1 Blot

Material Safety Data Sheet

Section 1: Chemical Identification

for Research Use Only

Name: Human Tumor Esophagus Single Tissue Blot

Catalog #: TB91

Section 2: Composition/Information on Ingredients

Protein Transfer membrane; PVDF containing bands of denatured proteins from Human Tumor Esophagus

Section 3: Hazards Identification

Believed to be NON-Hazardous

Section 4: First-Aid Measures

None needed

Section 5: Fire Fighting Measures

None needed

Section 6: Accidental Release Measures

None needed

Section 7: Handling and Storage

Ambient to 4⁰C.

Section 8: Exposure Controls/Personal Protection

General Protective and Hygiene Measures: Avoid contact with eyes, skin and clothing. Wash hands thoroughly after handling.

Section 9: Physical and Chemical Properties

Thin membrane sandwiched in blotting papers.

Section 10: Stability and Reactivity

Stable at 4⁰C to Room Temp, Non-oxidizing agents

Section 11: Toxicological Information

Believed to be non-toxic

Section 12: Ecological Information

Data not yet available

Section 13: Disposal Considerations

Observe all federal, state and local environmental regulations.

Section 14: Transport Information

Contact *Genotech* for transportation information.

Section 15: Regulatory Information

Irritating to eyes, respiratory system and skin. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing.

Section 16: 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. 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