

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

**Product:** Normal Human Brain Single Tissue Blot      **Catalog #:** TB07      [LU 5.20.09/MM]

**Components**

**Quantity**

Normal Human Brain Single Tissue Blot

1 Blot

**Material Safety Data Sheet**

**Section 1: Chemical Identification**

*for Research Use Only*

**Name:** Normal Human Brain Single Tissue Blot

**Catalog #:** TB07

**Section 2: Composition/Information on Ingredients**

Protein Transfer membrane; PVDF containing bands of denatured proteins from Normal Human Brain tissue

**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 4<sup>0</sup>C.

**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 4<sup>0</sup>C 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](http://www.GBiosciences.com), email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)*