

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

**Product:** Tween<sup>®</sup> 80 [Polyethylene glycol sorbitan monooleate] [LU 5.11.09/MK]

<u>Components</u>	<u>Catalog No.</u>	<u>Quantity</u>
Tween <sup>®</sup> 80 [Polyethylene glycol sorbitan monooleate]	DG013	5 x 5ml
	DG014	10 x 10ml
	DG513	1 x 50ml
	DG520	1 x 100ml

## Material Safety Data Sheet

### Section 1: Chemical Identification

**Name:** Tween<sup>®</sup> 80 [Polyethylene glycol sorbitan monooleate]

**Catalog #:** a component of product Cat # DG013, DG014, DG513, and DG520 indicated above

### Section 2: Composition/Information on Ingredients

Tween<sup>®</sup> 80 [Polyethylene glycol sorbitan monooleate] 99%, [CAS# 9005-65-6]

### Section 3: Hazards Identification

Label precautionary statements. 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 4: First-Aid Measures

**After eye contact:** Immediately flush eyes with copious amounts of water for at least 15 minutes.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** Remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, give oxygen.

**After swallowing:** Wash out mouth with water provided person is conscious. Call a physician. Wash contaminated clothing before reuse.

### Section 5: Fire Fighting Measures

**Extinguishing Agents:** Water spray, Carbon dioxide, dry chemical powder or appropriate foam.

**Special fire fighting procedures:** None Needed

**Unusual fire and explosions hazards:** None known

### Section 6: Accidental Release Measures

**Human Safety Precautions:** Wear chemical safety goggles, rubber boots and heavy rubber gloves.

**Cleaning/Collecting Measures:** Sweep up, place in a bag and hold for waste disposal. Ventilate area and wash spill site after material pickup is complete.

### **Section 7: Handling and Storage**

Refer to section 8.

### **Section 8: Exposure Controls/Personal Protection**

**Personal protective equipment:** Chemical safety goggles, compatible chemical-respirator, NIOSH/MSHA-approved respirator, and safety shower and eye bath.

**General Protective and Hygiene Measures:** Avoid contact with eyes, skin and clothing. Wash hands thoroughly after handling. Irritant. Keep tightly closed. Store in a cool dry place.

### **Section 9: Physical and Chemical Properties**

**Form:** Viscous Liquid

**Color:** Yellow

**Odor:** None

### **Section 10: Stability and Reactivity**

**Thermal Decomposition/Conditions to be avoided:** No decomposition if used according to specifications. Avoid strong oxidizing, heavy metal salts, bases.

**Dangerous Reactions:** No dangerous reactions known.

**Dangerous Products of Decomposition:** Carbon oxides.

### **Section 11: Toxicological Information**

**Primary Irritant Effects:** Harmful if inhaled or ingested.

**On the skin:** Material is irritating to skin and mucous membranes.

**On the eye:** Causes eye irritation.

**Sensitization:** No sensitizing effects known.

### **Section 12: Ecological Information**

Data not yet available.

### **Section 13: Disposal Considerations**

#### **Recommendation:**

Must not be disposed of together with household garbage. Do not allow product to reach sewage system. Hand over to hazardous waste disposers. Observe all federal, state and local environmental regulations.

### **Section 14: Transport Information**

Not dangerous goods. Safe for all transport.

**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. G-Biosciences/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)*