

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

**Product:** HOOK™-Biotin Kit **Catalog #:** BK-01 to BK-20 [LU 5.1.09/MK]

**Components Quantity**

Hook Biotin Agent 25mg x 1

(Selected from the list BK-01 to BK-20)

Optimizer Buffer I to V [5X] 25ml x 2

(As per Biotin agent selected)

Tube-O-Reactor Tube-O-Dialyzer (4kDa, Micro) x 5

With 5 Floats, 5 Microdialyzer tanks and 30 Glass balls

Avidin 5mg x 2

HABA Dye Reagent 1ml x 1

BiotinQuant Assay Buffer 25ml x 1

## **Material Safety Data Sheet**

### **Section 1: Chemical Identification**

**Name:** Hook-Biotin Kit

**Catalog #:** BK-01 to BK20 as indicated above.

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

Contains one or more of the following items - Derivatives of Biotin, Dilute solution of Acid/Base/Salts, chelating agent and Sodium Azide (0.05%).

### **Section 3: Hazards Identification**

Irritating to Eyes and Skin. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing and protective Eye glasses.

### **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:**

Wear self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.

**Unusual fire and explosions hazards:** Emits toxic fumes under fire conditions.

**Section 6: Accidental Release Measures**

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

**Cleaning/Collecting Measures:** Sweep up, place in a bag and hold for waste disposal. Avoid raising dust. 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 Hygienic Measures:**

Mechanical exhaust required. Do not breathe in dust. Avoid contact with eyes, skin and clothing. Wash thoroughly after handling. Irritant. Keep tightly closed. Store in a cool dry place.

**Section 9: Physical and Chemical Properties**

Reagents are clear liquid and in solid powder form with no odor

**Section 10: Stability and Reactivity****Thermal Decomposition/Conditions to be Avoided:**

No decomposition if used according to specifications.

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

**Dangerous Products of Decomposition:** No known dangerous decomposition products.

**Section 11: Toxicological Information****Acute Effects/Toxicity:**

**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.

**Additional toxicological information:** None

**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**

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)