

Product: Hook PEO Iodoacetyl Biotin

Catalog #: BG-11

[LU 5.1.09/MK]

Components

Quantity

PEO Iodoacetyl Biotin

2 x 25mg

Material Safety Data Sheet

Section 1: Chemical Identification

Name: Biotinyl-Iodoacetamidyl-3,6-dioxaoctanediamine

Catalog #: BG-11

Section 2: Composition/Information on Ingredients

Formula: C₁₈H₃₁N₄O₅SI CAS #: NA

MW: 542.43

Section 3: Hazards Identification

Label precautionary statements. Harmful Toxic. May cause liver damage. Irritating to respiratory system. Risk of serious damage to eyes. Possible sensitizer. 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

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

In case of skin contact: Immediately wash skin with soap and copious amounts of water.

In case of Inhalation: Remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, give oxygen.

In case of 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 Hygenic 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

Physical Properties

White to off-white solid

Section 10: Stability and Reactivity

Stability

Stable

Incompatibilities

Strong oxidizing agents, strong acids, strong bases. Reactive with strong nucleophiles, such as amines and mercaptans. Decomposes with heat.

Hazardous Combustion or Decomposition Products

Carbon monoxide, Carbon dioxide, Sulfur oxide, Nitrogen oxides, Possibly NH₃.

Section 11: Toxicological Information

Acute Effects/Toxicity: May cause skin irritation.

Primary Irritant Effects: Harmful if absorbed through the skin, inhaled or ingested.

On the skin: Material is irritating to mucous membranes and upper respiratory tract.

On the eye: Causes eye irritation.

Toxicological properties have not been thoroughly investigated. Exercise caution when handling.

Section 12: Ecological Information

Data not yet available

Section 13: Disposal Considerations

Recommendation: Appropriate Method of Disposal of Substance or Preparation: Must not be disposed of together with household garbage. Do not allow product to reach sewage system. Hand over to hazardous waste disposers.

Contact a licensed professional waste disposal service to dispose of this material.. Dissolve or mix the material with combustible solvent and burn in a chemical incinerator equipped with an afterburner and scrubber.

Observe all federal, state and local environmental regulations.

Section 14: Transport Information

Contact *Genotech* for transportation information.

Section 15: Regulatory information

Substance not yet fully tested 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. Observe all federal, state and local environmental regulations

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