

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

Product: Iodoacetamide **Catalog #:** 786-228 [LU 7.16.09/JK]

Components **Quantity**

Iodoacetamide 5g

Material Safety Data Sheet

Section 1: Chemical Identification

Name: Iodoacetamide

Catalog #: a component of product Cat # 786-228 as indicated above

Section 2: Composition/Information on Ingredients

Iodoacetamide 98-100%, Formula: C₂H₄INO, CAS # 144-48-9

Section 3: Hazards Identification

Toxic if swallowed. Irritating to eyes, respiratory system, and, skin. Label precautionary statements.

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. 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 Hygiene Measures:

Mechanical exhaust required.

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: Solid

Color: Variable, slightly variable

Odor: Odorless

Section 10: Stability and Reactivity

Conditions / Materials to be avoided: Light, Strong acids, Strong bases, Strong oxidizing agents, and Strong reducing agents.

Hazardous Decomposition Products: Carbon monoxide, Carbon dioxide, Nitrogen oxide, Hydrogen iodide.

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: May cause allergic reaction.

Signs and Symptoms of Exposure: Symptoms of exposure may include burning sensation, coughing, wheezing, laryngitis, shortness of breath, headache, nausea, and vomiting.

Chronic Exposure – Mutagen

Species: Human, Dose: 10UMOL/L, Cell Type: Lymphocyte, Mutation Test: DNA Inhibition

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

Shipping Name: Toxic Solids, organic, n.o.s.

UN#: 2811, Class: 6.1, Packing Group: III, Hazard Label: Toxic substances

Section 15: Regulatory information

US Classification and Label Text: Indication of Danger: Toxic, Risk Statements: Toxic if swallowed. May cause sensitization by inhalation and skin contact. Safety Statements: Do not breathe dust. Wear suitable protective clothing and gloves. In case of accident or if you feel unwell, seek medical advice immediately.

European Union (EU) Additional Classification: Symbol of Danger: T, Indication of Danger: Toxic, R: 25 42/43, Risk Statements: Toxic if swallowed. May cause sensitization by inhalation and skin contact. S: 22 36/37 45. Safety Statements: Do not breathe dust. Wear suitable protective clothing and gloves. In case of accident or if you feel unwell, seek medical advice immediately.

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 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