

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

Product: Agarose 1 **Catalog #:** RC-005, RC-006 & RC-149 [LU 04.03.09/JK/IA]

<u>Components</u>	<u>Catalog #</u>	<u>Quantity</u>
Agarose 1	RC-005	25g
Agarose 1	RC-006	100g
Agarose 1	RC-149	500g

Material Safety Data Sheet

Section 1: Chemical Identification

Name: Agarose 1

Catalog #: a component of product RC-005, RC-006 & RC-149 indicated above

Section 2: Composition/Information on Ingredients

Agarose 1; CAS# 9012-36-6

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. Call a physician.

After skin contact: Immediately wash skin with soap and copious amounts of water. Remove contaminated clothing and shoes. Call a physician.

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 immediately.

Section 5: Fire Fighting Measures

Extinguishing Agents: 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 self-contained breathing apparatus, 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.

Do not breathe dust. Do not get in eyes, on skin or clothing. Avoid prolonged or repeated exposure. . Keep tightly closed. Store in a cool dry place. Light sensitive.

Section 9: Physical and Chemical Properties

Form: Solid

Color: White

Odor: None

Section 10: Stability and Reactivity

Thermal Decomposition/Conditions to be avoided: Air, light.

Dangerous Reactions: Avoid contact with acid, oxidizing agents, iron and iron salts, copper, brass, free radical initiators.

Dangerous Products of Decomposition: Carbon monoxide, carbon dioxide, ammonia.

Section 11: Toxicological Information

Primary Irritant Effects: Toxic if inhaled or ingested.

On the skin: Material is toxic to skin and mucous membranes.

On the eye: Causes eye irritation.

Sensitization: Sensitizer.

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. Geno Technology 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