

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

Product: MES SDS Running Buffer [20X] **Catalog #:** 786-531 [LU 5.04.09/JK]

<u>Components</u>	<u>Catalog No.</u>	<u>Quantity</u>
MES SDS Running Buffer [20X]	786-531	500ml

Material Safety Data Sheet

Section 1: Chemical Identification

Name: MES SDS Running Buffer [20X]

Catalog #: a component of product Cat # 786-531 as indicated above

Section 2: Composition/Information on Ingredients

Contains: Tris hydroxymethyl aminomethane [CAS# 77-86-1] <30%; MES [CAS#] <30%; EDTA [CAS# 60-00-4] <5%; SDS [CAS# 151-21-3] <5%

Section 3: Hazards Identification

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

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.

After Inhalation: Remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, call a physician.

After swallowing: Wash out mouth with water provided person is conscious. Call a physician.

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 needed.

Section 6: Accidental Release Measures

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

Cleaning/Collecting Measures: Mop up, place in a closed container 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: Mechanical exhaust required.

Avoid contact with eyes, skin and clothing. Wash thoroughly after handling. Keep tightly closed. Store in cool dry place.

Section 9: Physical and Chemical Properties

Form: Liquid

Color: None

Odor: None

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

On the skin: May be harmful if absorbed through the skin. May cause skin irritation.

On the eye: May cause eye irritation.

Inhalation: May be harmful if inhaled. Material may be irritating to mucous membranes and upper respiratory tract.

Ingestion: May be harmful if swallowed.

Section 12: Ecological Information

Data not yet available

Section 13: Disposal Considerations

Recommendation:

Observe all federal, state and local environmental regulations.

Section 14: Transport Information

UN# 2926 Proper shipping name: Toxic, organic, N.O.S.

Class: 4.1 (6.1) Packing group: II

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