

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

**Product:** SAM*fluoro*: SAM Methyltransferase Assay Kit      **Catalog #:** 786-431      [LU 09.04.09/CCH-IA]

**Components**

**Quantity**

|   |           |
|---|-----------|
| SAM Methyltransferase Assay Buffer (0.1M Tris·HCl, pH8.0) | 20ml      |
| Positive Control : Adenosylhomocysteine (1 mM)            | 0.2ml     |
| SAM <i>fluoro</i> Enzyme Mix                              | 5 x 300µl |
| S-Adenosylmethionine                                      | 5 x 200µl |
| SAM Methyl transferase Assay Buffer Additive              | 0.2ml     |
| SAM <i>fluoro</i> Fluorometric Mix                        | 3 vials   |
| Resorufin Standard  | 100µl     |
| DMSO  | 1 ml      |
| Resorufin Buffer  | 50 ml     |

**Material Safety Data Sheet**

**Section 1: Chemical Identification**

**Name:** SAM Methyltransferase Assay Buffer

**Catalog #:** a component of product Cat# 786-431 indicated above

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

Contains Tris hydroxymethyl aminomethane [CAS# 77-86-1];  $\text{NH}_2\text{C}(\text{CH}_2\text{OH})_3$ , <0.2M

**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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

**Section 6: Accidental Release Measures**

**Human Safety Precautions:** None needed

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

**Section 7: Handling and Storage**

Refer to section 8.

**Section 8: Exposure Controls/Personal Protection**

**Personal protective equipment:** Chemical safety gloves

**General Protective and Hygiene Measures:** Avoid contact with eyes, skin and clothing. Wash thoroughly with water after handling. Irritant. Keep tightly closed. Store in a cool dry place.

**Section 9: Physical and Chemical Properties**

**Form:** Liquid

**Color:** Colorless

**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**

**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.

**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**

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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

### **Material Safety Data Sheet**

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

**Name:** Positive Control Adenosylhomocysteine

**Catalog #:** a component of product Cat# 786-431 indicated above.

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

Contains Adenosylhomocysteine solution

#### **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

## **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

## **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:** 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**

**Form:** Liquid

**Color:** Clear

**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**

**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.

## **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** 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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## **Material Safety Data Sheet**

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

**Name:** SAMfluoro Enzyme Mix

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains proprietary Methyltransferase Enzyme mix

### **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

### **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

#### **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:** 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**

**Form:** Liquid

**Color:** Clear

**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**

**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.

#### **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**

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. 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*

*[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)*

### **Material Safety Data Sheet**

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

**Name:** S-Adenosylmethionine

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains S-Adenosylmethionine [CAS# 17176-17-9]

#### **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

#### **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

## **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:** 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**

**Form:** Liquid

**Color:** Clear

**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**

**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.

## **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**

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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## **Material Safety Data Sheet**

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

**Name:** SAM Methyltransferase Assay Buffer Additive

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains Manganese Chloride; MnCl<sub>2</sub>, [CAS# 13446-34-9]

### **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

### **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

### **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:** 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**

**Form:** Liquid

**Color:** Clear

**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**

**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.

### **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**

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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## **Material Safety Data Sheet**

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

**Name:** SAMfluoro Fluorometric Mix

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains proprietary non-hazardous fluorescent mix

### **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

### **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

### **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:** 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**

**Form:** Liquid

**Color:** Clear to slight orange

**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**

**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.

### **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**

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. 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*

*[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)*

## **Material Safety Data Sheet**

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

**Name:** Resorufin Standard

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains: Resorufin CAS# 635-78-9

### **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

### **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

**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:** 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**

**Form:** Liquid

**Color:** purple-red

**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**

**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.

**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** 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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

# **Material Safety Data Sheet**

## **Section 1: Chemical Identification**

**Name:** DMSO

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains: DMSO [CAS# 67-68-5]

## **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

## **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

## **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:** 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**

**Form:** Liquid

**Color:** Clear

**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**

**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.

**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** 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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

# **Material Safety Data Sheet**

## **Section 1: Chemical Identification**

**Name:** Resorufin Buffer

**Catalog #:** A component of product Cat# 786-431 indicated above.

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

Contains: Tris-HCl [CAS# 1185-53-1] <5%

## **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.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** None needed

**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:** Not known

## **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Chemical safety goggles, rubber boots and rubber gloves.

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

## **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:** 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**

**Form:** Liquid

**Color:** Clear

**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**

**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.

## **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** 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. 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*

[www.GBiosciences.com](http://www.GBiosciences.com); email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)