

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

Product: CyanoBlue Universal Dye

Catalog #: 786-101

[LU 07.15.09/JK-IA]

Components

Quantity

CyanoBlue Universal Dye

1.5ml/vial

Material Safety Data Sheet

Section 1: Chemical Identification

Name: CyanoBlue Universal Dye

Catalog #: a component of product Cat # 786-101 indicated above

Section 2: Composition/Information On Ingredients

One or more of the following items- Bromphenol Blue, Xylene cyanol, formide, ethidium bromide and Sodium dodecylsulfate.

Section 3: Hazards Identification

Label precautionary statements, Irritant, irritating to eyes, 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 contact, immediately flush eyes with copious amounts of water for at least 15 minutes. If swallowed, wash out mouth with water provided person is conscious. Call the physician. Wash contaminated clothing before reuse

Section 5: Fire Fighting Measures

Extinguishing media, 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.

Section 6: Accidental Release Measures

Sweep up, place in a bag and hold for waste disposal.

Section 7: Handling And Storage

Refer to section 8.

Section 8: Exposure Control And Personal Protection

Chemical safety goggles, Safety shower and eye bath. Mechanical exhaust required. Do not breathe dust & vapor. 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

Appearance and odor: dark color liquid

Section 10: Stability And Reactivity

STABLE AT 0-40°C TEMPERATURE, Avoid strong oxidizing agents. May produce carbon oxides, sulfur oxides and bromine.

Section 11: Toxicological Information

Known carcinogen, May be harmful by inhalation, ingestion, or skin absorption. Causes eye and skin irritation. To the best of our knowledge, the chemical, physical, and toxicological properties have not been thoroughly investigated.

Section 12: Ecological Information

Data not yet available

Section 13: Disposal Considerations

Dissolve or mix with plenty of water.

Observe all federal, state and local environmental regulations

Section 14: Transport Information

Contact g-biosciences for transportation information

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 complete 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.

G-Biosciences/Genotech St. Louis, MO, USA

Web: www.GBiosciences.com, email: technical@GBiosciences.com