

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

**Product:** Protein Standard – BSA

**Catalog #:** 786-006

[LU 4.15.09-IA/JK]

**Components**

**Quantity**

Protein Standard BSA (2mg/ml)

5ml

**Material Safety Data Sheet**

**Section 1: Chemical Identification**

**Name:** Protein Standard- BSA

**Catalog #:** a component of product Cat #786-006 indicated above

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

2mg/ml solution of Bovine Serum Albumin in dilute solution of azide, CAS# 9048-46-8

**Section 3: Hazards Identification**

In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

**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**

**Special Fire Fighting Procedures:**

Special fire fighting measures not needed.

**Section 6: Accidental Release Measures**

**Cleaning/Collecting Measures:**

Wash with water, 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:**

**General Protective and Hygiene Measures:**

Avoid contact with eyes, skin and clothing. Wash thoroughly after handling.

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. Avoid moisture, ignition sources, strong oxidizing agents and strong acids.

**Dangerous Reactions:** No dangerous reactions known. Stable.

**Dangerous Products of Decomposition:** Carbon oxides.

**Section 11: Toxicological Information**

**Primary Irritant Effects:** Harmful if inhaled or ingested.

**On the skin:** None

**On the eye:** None

**Sensitization:** No sensitizing effects known.

**Section 12: Ecological Information**

Data not yet available

**Section 13: Disposal Considerations**

**Recommendation:**

Dissolve or mix the material with water. Laboratory waste disposal sink may be used.

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*

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