

Material Safety Data Sheet

Section 1 Product and Company Identification

Product number: B011
Product name: D-Biocytin TMR
Unit Size: 5 mg
Manufacturer/Supplier: GeneCopoeia Inc
9620 Medical Center Drive, Suite 101, Rockville, MD 20850, USA
Tel: 1-301-762-0888, Fax: 1-301-762-3888
Web: www.abpbio.com

Section 2 Hazardous Ingredients/Identity Information

Chemical Name	CAS-No	EINECS-No	Weight %
D-Biocytin TMR	-	-	60-100

Section 3 Physical and Chemical Properties

Form	solid
Appearance	No information available
Odor	No information available
Boiling point	No data available
Melting point	No data available
Flash point	No data available
Autoignition temperature	No data available
Oxidizing properties	No information available
Water solubility	No information available

Section 4 Potential Health Hazard Effects

Eyes	May cause eye irritation with susceptible persons.
Skin	May cause skin irritation in susceptible persons.
Inhalation	May be harmful by inhalation.
Ingestion	May be harmful if swallowed.
Carcinogenic effects	not applicable.
Reproductive toxicity	not applicable.
Sensitization	not applicable.

Section 5 Stability and Reactivity Data

Stability	Stable under normal conditions.
Materials to avoid	No information available.
Hazardous decomposition	No information available.
Polymerization	Hazardous polymerization does not occur.

Section 6 Toxicological and Ecological Information

Chemical Name	LD50 (oral, rat/mouse)	LD50 (dermal, rat/rabbit)	LC50 (inhalation, rat/mouse)
D-Biocytin TMR	No data available	No data available	No data available

Ecotoxicity effects

No information available.

Mobility

No information available.

Biodegradation

Inherently biodegradable.

Bioaccumulation

Does not bioaccumulate.

Section 7 Handling and Storage

Handling

Always wear recommended Personal Protective Equipment. No special handling advice required.

Storage

Keep in a dry, cool and well-ventilated place.

Section 8 First Aid Measures

Skin contact

Rinse with plenty of water. If symptoms arise, call a physician.

Eye contact

Rinse immediately with plenty of water, also under the eyelids, for at least 15 minutes. If symptoms persist, call a physician.

Ingestion

Never give anything by mouth to an unconscious person. If symptoms persist, call a physician. Do not induce vomiting without medical advice.

Inhalation

Move to fresh air. If symptoms persist, call a physician. If not breathing, give artificial respiration.

Section 9 Disposal Considerations

Dispose of in accordance with local regulations.

Section 10 Other Information

For research use only. Not intended for human or animal diagnostic or therapeutic uses.

End of Material Safety Data Sheet