

Material Safety Data Sheet

Section 1 Product and Company Identification

Product number: B004
Product name: D-Biotin-X, succinimidyl ester
Unit Size: 100 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-Biotin-X, succinimidyl ester	72040-63-2	-	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-Biotin-X, succinimidyl ester	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