

## Material Safety Data Sheet

### Section 1 Product and Company Identification

---

**Product number:** B005  
**Product name:** D-Biotin-XX, succinimidyl ester  
**Unit Size:** 50 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](http://www.abpbio.com)

### Section 2 Hazardous Ingredients/Identity Information

---

Chemical Name	CAS-No	EINECS-No	Weight %
D-Biotin-XX, succinimidyl ester	89889-52-1	-	60-100

### Section 3 Physical and Chemical Properties

---

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

### Section 4 Potential Health Hazard Effects

---

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

### Section 5 Stability and Reactivity Data

---

<b>Stability</b>	Stable under normal conditions.
<b>Materials to avoid</b>	No information available.
<b>Hazardous decomposition</b>	No information available.
<b>Polymerization</b>	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-XX, succinimidyl ester	No data available	No data available	No data available

<b>Ecotoxicity effects</b>	No information available.
<b>Mobility</b>	No information available.
<b>Biodegradation</b>	Inherently biodegradable.
<b>Bioaccumulation</b>	Does not bioaccumulate.

## Section 7 Handling and Storage

---

<b>Handling</b>	Always wear recommended Personal Protective Equipment. No special handling advice required.
<b>Storage</b>	Keep in a dry, cool and well-ventilated place.

## Section 8 First Aid Measures

---

<b>Skin contact</b>	Rinse with plenty of water. If symptoms arise, call a physician.
<b>Eye contact</b>	Rinse immediately with plenty of water, also under the eyelids, for at least 15 minutes. If symptoms persist, call a physician.
<b>Ingestion</b>	Never give anything by mouth to an unconscious person. If symptoms persist, call a physician. Do not induce vomiting without medical advice.
<b>Inhalation</b>	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**