

Material Safety Data Sheet (MSDS)

Revision Number: 1.0

Last updated February 2004

Product and Company Identification

Product Name: DiD
[1,1'-Dioctadecyl-3,3,3',3'- tetramethylindodicarbocyanine iodide]
Catalog Number: 84903
Unit Size: 25 mg
Manufacturer/Supplier: AnaSpec, Inc.
2149 O'Toole Avenue
San Jose, CA 95131
Tel: 408-452-5055
Fax: 408-452-5059
<http://www.anaspec.com>
e-Mail: service@anaspec.com

Composition / Information on Ingredients

DiD
[1,1'-Dioctadecyl-3,3,3',3'- tetramethylindodicarbocyanine iodide]
Molecular Formula: C₆₁H₉₉IN₂
Molecular Weight: 987.3
CAS Number/Name: 127274-91-3

Hazards Identification

Emergency Overview: Avoid contact and inhalation. To our knowledge, the hazards of this material have not been thoroughly investigated. We recommend handling all chemicals with caution.

Potential Health Effects

<i>Inhalation:</i>	Not determined
<i>Ingestion:</i>	Not determined
<i>Skin:</i>	Not determined
<i>Eyes:</i>	Not determined
<i>Chronic Exposures:</i>	Not determined
<i>Target Organs:</i>	Not determined

Potentially harmful. Avoid prolonged or repeated exposure. Wash thoroughly after handling. If eye or skin contact occurs, immediately wash affected area with soap and copious amounts of water for at least 15 minutes. Assure adequate flushing by separating the eyelids with fingers. Call a physician to seek medical advice. If swallowed, wash out mouth with water provided person is conscious and call a physician to seek medical advice. If inhaled, move individual to fresh air and call a physician to seek medical advice.

Fire Fighting Measures

Extinguishing media:

Water spray, carbon dioxide, dry chemical powder or appropriate foam.

Special firefighting procedures:

Wear self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.

Unusual fire and explosions hazards:

Emits toxic fumes under fire conditions.

Accidental Release Measures

Use appropriate protective equipment and methods to clean up spilled substances promptly. Absorb spill onto an appropriate material. Collect and dispose of all waste in accordance with applicable laws.

Handling and Storage

Protect material from long-term exposure to light; may be exposed to light for short periods of time.

Exposure Controls / Personal Protection

Wear appropriate gloves, protective clothing and eyewear and follow safe laboratory practices.

ACGIH/OSHA Permissible Exposure Limit Data:

Not determined

Physical and Chemical Properties

Form:

Solid

Odor:

Not determined

Solubility in Water:

Low

Specific Gravity:

Not determined

pH:

Not determined

Boiling Point:

Not determined

Melting Point:

Not determined

Flash Point:

Not determined

Vapor Pressure:

Not determined

Stability and Reactivity

Thermal Decomposition:

No decomposition if used according to specifications.

Dangerous Products of Decomposition:

No dangerous decomposition products identified.

Dangerous Reactions:

No dangerous reactions identified.

Toxicological Information

RTECS Number:

None known

Toxicity:

We are not aware of any toxicity data for this product.

Health Hazards:

Not thoroughly investigated.

Potential Hazards:

The complete properties have not been investigated; however, similar compounds are known to be chemically reactive with proteins and other biochemicals and should be treated as potentially hazardous.

Carcinogenicity:

Not listed by NTP, IARC or OSHA.

Ecological Information

Do not allow product to reach ground water, water course, or sewage system.

Disposal Considerations

Consult local, state or national regulations for proper disposal.

Transport Information

<i>Hazard Class:</i>	Not determined
<i>Identification Number:</i>	Not determined
<i>Packing Group:</i>	Not classified
<i>Proper Shipping Name (Technical Name):</i>	Not determined

Regulations

<i>US Toxic Substances Control Act (TSCA):</i>	Not listed
<i>US Other:</i>	Not applicable
<i>EEC EINECS Number:</i>	Not identified
<i>Other Country Regulations:</i>	None identified
<i>EEC Risk Statements:</i>	Not determined

Other Information

This material is sold for research purposes only and is not required to appear on the TSCA inventory. It is not intended for food, drug, household, agricultural or cosmetic use. Its use must be supervised by a technically qualified individual experienced in handling potentially hazardous chemicals. The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. Users should make independent decisions regarding completeness of the information based on all sources available. HiLyte Biosciences shall not be held liable for any damage resulting from handling or from contact with the above product.