



Material Safety Data Sheet (MSDS)

Revision Number: 1.1		Last updated April 2010	
<u>Product and Company Identification</u>			
Product Name:		Fmoc - (N - β - Boc) - D - α , β - diaminopropionic acid	
Manufacturer/Supplier:		AnaSpec, Inc. 34801 Campus Drive, Fremont, CA 94555 Tel: 510-791-9560 Fax: 510-791-9573 http://www.anaspec.com	
<u>Composition / Information on Ingredients</u>			
Fmoc - (N - β - Boc) - D - α , β - diaminopropionic acid			
Fmoc - D - Dap(Boc) - OH			
Molecular Formula: C ₂₃ H ₂₆ N ₂ O ₆			
Molecular Weight: 426.5			
CAS Number: 198544 - 42 - 2			
<u>Hazards Identification</u>			
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			
Inhalation:		If inhaled, remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, give oxygen.	
Ingestion:		If swallowed, wash out mouth with water provided person is conscious. Call a physician.	
Skin:		In case of contact, immediately wash skin with soap and plenty of water.	
Eyes:		In case of contact, immediately flush eyes with copious amounts of water for at least 15 minutes.	
Chronic Exposures:		Not determined	
Target Organs:		Not determined	

AnaSpec Inc.

34801 Campus Drive, Fremont, CA 94555

- Toll-Free: 800-452-5530 • Tel: 510-791-9560 • Fax: 510-791-9572



<u>Fire Fighting Measures</u>	
<i>Extinguishing media:</i>	Water spray, carbon dioxide, dry chemical powder or appropriate foam.
<i>Special firefighting procedures:</i>	Wear self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.
<i>Unusual fire and explosions hazards:</i>	Emits toxic fumes under fire conditions.
<u>Accidental Release Measures</u>	
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. Avoid dust formation. Do not let product enter drains.	
<u>Handling and Storage</u>	
Store at 2-8 °C. Keep container tightly closed in a dry and well-ventilated place.	
<u>Exposure Controls / Personal Protection</u>	
Wear appropriate gloves, protective clothing and eyewear and follow safe laboratory practices.	
<u>ACGIH/OSHA Permissible Exposure Limit Data:</u>	
Not determined	
<u>Physical and Chemical Properties</u>	
<i>Form:</i>	Solid
<i>Color:</i>	White
<i>Solubility in Water:</i>	Not soluble
<i>Boiling Point:</i>	Not determined
<i>Melting Point:</i>	Not determined
<i>Flash Point:</i>	Not determined
<u>Stability and Reactivity</u>	
<i>Stability:</i>	Stable under recommended storage conditions
<i>Dangerous Products of Decomposition:</i>	Carbon monoxide, carbon dioxide and nitrogen oxides
<i>Materials to avoid:</i>	Strong oxidizing agents
<u>Toxicological Information</u>	
<i>Toxicity:</i>	We are not aware of any toxicity data for this product.
<i>Health Hazards:</i>	Skin Contact: May cause skin irritation. Skin absorption: May be harmful if absorbed through skin Eye Contact: May cause eye irritation. Inhalation: May be harmful if inhaled. Ingestion: May be harmful if swallowed.
<i>Carcinogenicity:</i>	Not listed by NTP, IARC or OSHA.

AnaSpec Inc.

34801 Campus Drive, Fremont, CA 94555

• Toll-Free: 800-452-5530 • Tel: 510-791-9560 • Fax: 510-791-9572

**Ecological Information**

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

Disposal Considerations

Consult local, state or national regulations for proper disposal.

Transport Information

This substance is considered to be non-hazardous for transport

Regulations

OSHA Hazard:	No known OSHA hazards
SARA 311/312 Hazard:	No SARA hazards
Other Country Regulations:	None identified
EEC Risk Statements:	Not determined

Other Information

It is not intended for food, drug, household, agricultural or cosmetic use. A technically qualified individual experienced in handling potentially hazardous chemicals must supervise its use. 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. Anaspec shall not be held liable for any damage resulting from handling or from contact with the above product.

AnaSpec Inc.

34801 Campus Drive, Fremont, CA 94555

• Toll-Free: 800-452-5530 • Tel: 510-791-9560 • Fax: 510-791-9572