

Material Safety Data Sheet (MSDS)

<i>Revision Number: 1.1</i>	<i>Last updated December 2010</i>
<u>Product and Company Identification</u>	
<i>Product Name:</i>	<i>ClearPoint™ Angiotensin II, human, 13C and 15N - labeled</i>
<i>Catalog Number:</i>	<i>AS-64805</i>
<i>Manufacturer/Supplier:</i>	AnaSpec, Inc. 34801 Campus Drive Fremont, CA 94555 http://www.anspec.com
<u>Composition / Information on Ingredients</u>	
<i>Sequence:</i> <i>H - Asp - Arg - Val - Tyr - *Ile - His - Pro - Phe - OH [Ile* = Ile(U¹³C6, ¹⁵N)]</i> <i>Molecular Weight: 1053.2</i> <i>CAS Number/Name: NA</i>	
<u>Hazards Identification</u>	
Avoid contact and inhalation. To our knowledge, the hazards of this material have not been thoroughly investigated. We recommend handling all chemicals with caution	
<i>First Aid Measures:</i>	
<i>Inhalation:</i>	If inhaled, remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, give oxygen.
<i>Ingestion:</i>	If swallowed, wash out mouth with water provided person is conscious. Call a physician.
<i>Skin:</i>	In case of contact, immediately wash skin with soap and plenty of water.
<i>Eyes:</i>	In case of contact, immediately flush eyes with copious amounts of water for at least 15 minutes.
<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.	

AnaSpec Inc.

34801 Campus Drive, Fremont, CA 94555

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

<u>Handling and Storage</u>	
Avoid contact with skin and eyes. Store at -20 °C.	
<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>	Lyophilized powder
<i>Odor:</i>	Not determined
<i>Boiling Point:</i>	Not determined
<i>Melting Point:</i>	Not determined
<i>Flash Point:</i>	Not determined
<u>Stability and Reactivity</u>	
<i>Materials to avoid:</i>	Strong acids, strong bases
<i>Dangerous Products of Decomposition:</i>	Carbon oxides, nitrogen oxides (NOx)
<i>Dangerous Reactions:</i>	No dangerous reactions identified.
<u>Toxicological Information</u>	
<i>Toxicity:</i>	We are not aware of any toxicity data for this product.
<i>Health Hazards:</i>	May cause skin irritation. May cause eye irritation. Material may be irritating to mucous membranes and upper respiratory tract. May be harmful if inhaled. May be harmful if swallowed.
<i>Target Organ:</i>	Not determined
<i>Carcinogenicity:</i>	Not listed by NTP, IARC or OSHA.
<u>Ecological Information</u>	
Do not allow product to reach ground water, watercourse, or sewage system.	
<u>Disposal Considerations</u>	
Consult local, state or national regulations for proper disposal.	
<u>Transport Information</u>	
This substance is considered to be non-hazardous for transport	
<u>Regulations</u>	
<i>OSHA hazard:</i>	Not determined
<i>SARA 311/312 hazard:</i>	Not determined
<i>Other Country Regulations:</i>	None identified
<i>EEC Risk Statements:</i>	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