

Product Information Sheet

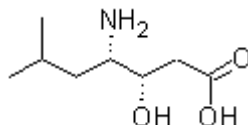
Ordering Information

Product Number: 5109
Product Name: Statine
Unit Size: 100 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 175.23

Chemical Structure:



Soluble in: DMF
Excitation Wavelength: NA
Emission Wavelength: NA

Application Notes

(3S,4S)-Statine is a novel amino acid of pepstatin, a low molecular weight inhibitor of acid proteases.