

Product Information Sheet

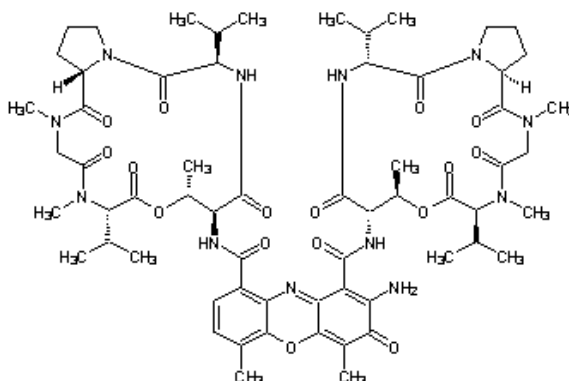
Ordering Information

Product Number: 17505
Product Name: Actinomycin D
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 1255.42

Chemical Structure:



Soluble in: DMSO

Application Notes

Actinomycin is a polypeptide antibiotic isolated from soil bacteria of the genus *Streptomyces*. It binds to DNA duplexes, thereby interfering with the action of enzymes engaged in replication and transcription. Actinomycin D is one of the old chemotherapy drugs which have been used in therapy for many years. Actinomycin D is also used as a selective antibiotic for resistant bacteria, generally at a concentration of about 1 µg/ml. This product is provided as a powder and should be made into a stock solution at 1 mg/ml in DMSO.