

Aclacinomycin A (Aclarubicin)

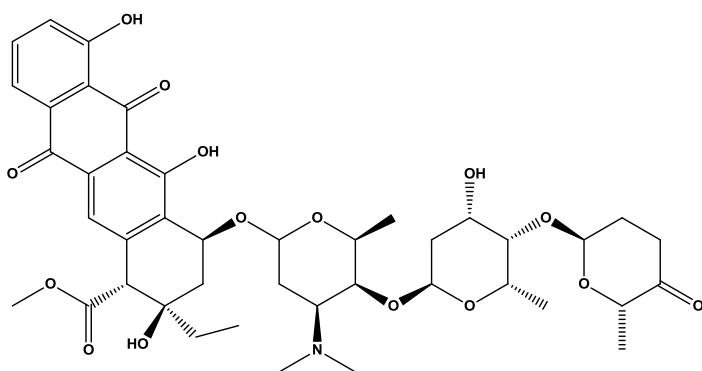
Cat. # SI9730

Background: Aclacinomycin A is specific inhibitor of the 20S proteasome chymotrypsin-like activity and calpain activity. It acts by forming a stable complex with DNA. It was originally identified as antitumor agent. Cell permeable.

Application: Inhibition of 20S proteasome activity.

Product Information

CAS No.	57576-44-0
Purity:	>95% (TLC); NMR (Conforms)
Molecular Weight:	811.9
Physical State:	Orange powder
Quantity:	10 mg
Solubility:	DMSO (30 mg/ml)
Storage:	Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to one month.



Formula: C₄₂H₅₃NO₁₅

References

1. Figueiroda-Pereira, ME *et al.* The antitumor drug aclacinomycin A, which inhibits the degradation of ubiquitinated proteins, shows selectivity for the chymotrypsin-like activity of the bovine pituitary 20 S proteasome. *J. Biol. Chem.* 1996, 271, 16455.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved