Aclacinomycin A

Originally identified as an anti-tumor drug, aclacinomycin A is the first described non-peptidic inhibitor showing discrete specificity for the chymotrypsin-like activity of the 20S proteasome. Aclacinomycin A has no effect on cathepsin B, it stimulates trypsin activity but inhibits the activity of both chymotrypsin and calpain.

Citations: 8

View Online »

Ordering Information

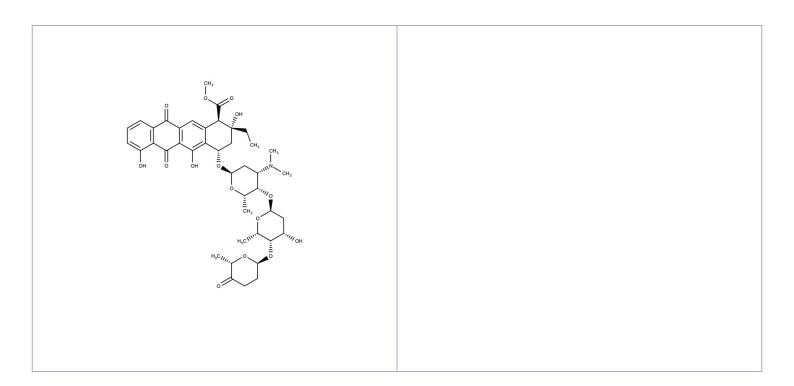
Order Online »

BML-AW8655-0005

5mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Aclarubicin

Appearance Yellow powder with orange cast.

CAS 57576-44-0

Couple Target Calpain, Chymotrypsin, Proteasome, Trypsin

Couple Type Activator, Inhibitor

Formula $C_{42}H_{53}NO_{15}$

MW 811.9

Purity ≥95% (TLC)

Soluble in DMSO (25mg/ml) or dimethyl formamide

(25mg/ml).

Last modified: May 29, 2024