

PRODUCT DATA SHEET



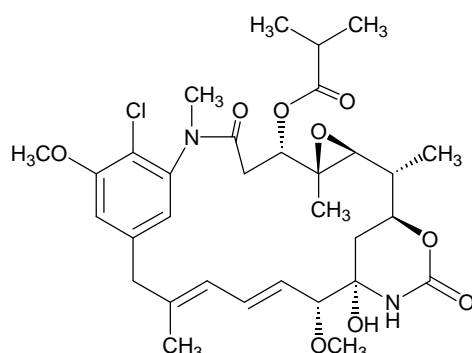
PRODUCT: Ansamitocin P-3

CAS NO: 66547-09-9

CATALOG NO.: T-121

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Lyophilized solid

MOLECULAR FORMULA: $C_{32}H_{43}ClN_2O_9$

MOLECULAR WEIGHT: 635.2

PURITY: 95%(TLC)

SOLUBILITY: Soluble in DMSO

STORAGE: Store, as supplied, at $-20^{\circ}C$ for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: A fungal metabolite from *Actinosynnema pretiosum* which binds to actin. Shares the same binding site as phalloidin and rhizoxin on porcine tubulin.¹ A maytansine analog which displays potent cytotoxicity against human solid tumor cell lines A-549 and HT-29.² Maytansine and analogs cause extensive disassembly of microtubules and totally prevent tubulin spiralization.³

- REFERENCES:
1. Y.Li *et al.* *Biochem.Biophys.Res.Commun.* 1992 **187** 722
 2. K.Suwanborirux *et al.* *Experientia* 1990 **46** 117
 3. K.Ootsu *et al.* *Cancer Res.* 1980 **40** 1707

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.