

# PRODUCT DATA SHEET



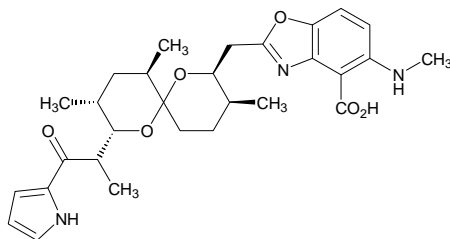
PRODUCT: Antibiotic A-23187  
Calcimycin

CAS NO: 52665-69-7

CATALOG NO.: CA-100

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Off-white solid (m.p. 181-182°C)

MOLECULAR FORMULA:  $C_{29}H_{37}N_3O_6$

MOLECULAR WEIGHT: 523.6

PURITY: 98% (TLC)

SOLUBILITY: Soluble in DMSO and ethanol

STORAGE: Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: Calcium ionophore. Causes rapid influx of calcium in various cells.

REFERENCES: 1. V.L.Lew and J.Garcia-Sancho *Cell.Calcium* 1985 **6** 15  
2. D.P.Negri and Y.Kishi *Tetrahedron.Lett.* 1987 **28** 1063  
3. T.A.Rao *et al. Life Sci.* 1993 **53** PL147

**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.**