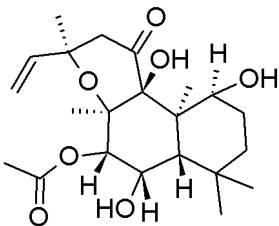


Coleonol, Colforsin, 7 β -Acetoxy-8,13-epoxy-1 α ,6 β ,9 α -trihydroxylabd-14-en-11-one

Instruction Manual

Catalog Number	PK-CA577-1531-5
Description	Activates adenylate cyclase leading to an increase in the intracellular concentration of cAMP.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	410.5 Da; C ₂₂ H ₃₄ O ₇
Chemical Structure	
Purity	≥98%
Appearance / Formulation / Solubility	Off-white powder; soluble in DMSO or 100% ethanol.
Storage & Stability	Store at -20°C. Protect from moisture.
Applications	see Description
References	NA
Caution	NA

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.