

PRODUCT INFORMATION



Kendomycin

Item No. 11037

CAS Registry No.: 183202-73-5
Formal Name: (13E)-6R,7S,8R,9R,10,11,12S,15,16S,17,18S,19S-dodecahydro-4,7,19-trihydroxy-2,6,8,12,14,16,18-heptamethyl-1,19:5R,9-diepoxybenzocyclooctadecen-3(5H)-one

Synonym: TAN 2162

MF: C₂₉H₄₂O₆

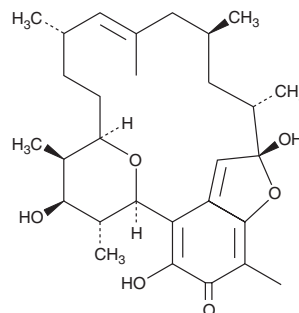
FW: 486.6

Purity: ≥90%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Kendomycin is a macrolide from *Streptomyces* which displays potent cytotoxicity against several carcinoma cell lines ($GI_{50} < 100$ nM).¹ It also acts as an endothelin receptor antagonist and has antiosteoporotic activities.¹ Certain effects are thought to be mediated through inhibition of proteasomal degradation of proteins.

Reference

1. Bahnck, K.B. and Rychnovsky, S.D. Formal synthesis of (-)-kendomycin featuring a Prins-cyclization to construct the macrocycle. *J. Am. Chem. Soc.* **130(39)**, 13177-13181 (2008).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/07/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM