

PRODUCT INFORMATION



PX-866-17OH

Item No. 21204

CAS Registry No.: 1012327-63-7
Formal Name: (1E,4S,4aR,5R,6aS,7S,9aR)-5-(acetyloxy)-1-[(di-2-propen-1-ylamino)methylene]-4a,5,6,6a,7,8,9,9a-octahydro-7,11-dihydroxy-4-(methoxymethyl)-4a,6a-dimethyl-cyclopenta[5,6]naphtho[1,2-c]pyran-2,10(1H,4H)-dione

MF: C₂₉H₃₇NO₈

FW: 527.6

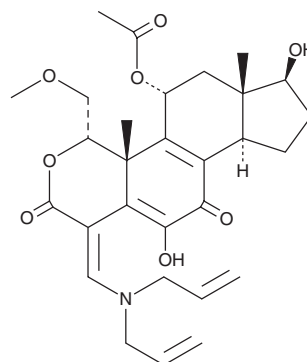
Purity: ≥98%

UV/Vis.: λ_{max}: 249, 316, 405 nm

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.

Laboratory Procedures

PX-866-17OH is supplied as a crystalline solid. A stock solution may be made by dissolving the PX-866-17OH in the solvent of choice, which should be purged with an inert gas. PX-866-17OH is soluble in the organic solvent DMSO.

Description

PX-866-17OH is an active metabolite of PX-866 (Item No. 13645).¹ It inhibits PI3Kα, PI3Kβ, PI3Kγ, and PI3Kδ (IC₅₀s = 14, 57, 131, and 148 nM, respectively).

Reference

1. Hausman, D., Kirkman, R., and Peterson, S. Combination therapy using PI3 kinase and braf inhibitors. *Oncothyreon Inc. WO 2014/031856 A1* (2014).

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, 12/16/2022

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