

PRODUCT INFORMATION

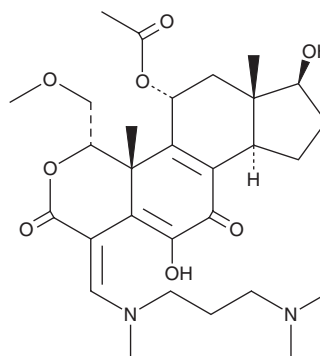


PX-13-17OH

Item No. 21203

CAS Registry No.: 884539-95-1
Formal Name: (1E,4S,4aR,5R,6aS,7S,9aR)-5-(acetyloxy)-1-[[[3-(dimethylamino)propyl]methylamino]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: C₂₉H₄₂N₂O₈
FW: 546.7
Purity: ≥98%
UV/Vis.: λ_{max}: 251, 319, 407 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-13-17OH is supplied as a crystalline solid. A stock solution may be made by dissolving the PX-13-17OH in the solvent of choice, which should be purged with an inert gas. PX-13-17OH is soluble in DMSO.

Description

PX-13-17OH is a PI3K inhibitor.¹ It selectively inhibits PI3K α , PI3K β , PI3K γ , and PI3K δ (IC₅₀s = 6.4, 13, 8, and 11 nM, respectively) over mTOR (IC₅₀ = 2.9 μ M). PX-13-17OH is greater than 420-fold selective for PI3K in a panel of 20 lipid and protein kinases. PX-13-17OH inhibits phosphorylation of Akt and S6 kinase (S6K) in PTEN-negative U87MG cells when used at concentrations ranging from 0.03 to 1 μ g/ml. It inhibits tumor growth in a U87MG mouse xenograft model when administered at doses ranging from 2.5 to 10 mg/kg.

References

1. Zask, A., Kaplan, J., Toral-Barza, L., *et al.* Synthesis and structure-activity relationships of ring-opened 17-hydroxywortmannins: Potent phosphoinositide 3-kinase inhibitors with improved properties and anticancer efficacy. *J. Med. Chem.* **51**(5), 1319-1323 (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, 12/05/2022

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