

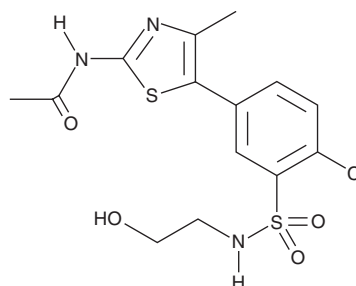
PRODUCT INFORMATION



PIK-93

Item No. 10009212

CAS Registry No.: 593960-11-3
Formal Name: N-[5-[4-chloro-3-[[[(2-hydroxyethyl)amino]sulfonyl]phenyl]-4-methyl-2-thiazolyl]-acetamide
MF: C₁₄H₁₆ClN₃O₄S₂
FW: 389.9
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 306 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

PIK-93 is supplied as a crystalline solid. A stock solution may be made by dissolving the PIK-93 in the solvent of choice. PIK-93 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of PIK-93 in ethanol is approximately 0.25 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

PIK-93 is an inhibitor of phosphatidylinositol 4-kinase IIIβ (PI4KIIIβ; IC₅₀ = 19 nM).¹ It is selective for PI4KIIIβ over PI4KIIIα and PI4KIIα (IC₅₀s = 1,100 and >100,000 nM, respectively). However, it also inhibits the PI3K catalytic subunits (IC₅₀s = 39, 590, 120, and 16 nM for p110α, p110β, p110δ, and p110γ, respectively). PIK-93 reduces enteroviral RNA replication when used at a concentration of 125 nM in HeLa cells.² It also inhibits the transport of ceramide from the endoplasmic reticulum to the Golgi apparatus.³

References

1. Knight, Z.A., Gonzalez, B., Feldman, M.E., *et al.* A pharmacological map of the PI3-K family defines a role for p110α in insulin signaling. *Cell* **125**(4), 733-747 (2006).
2. Hsu, N.Y., Illynska, O., Belov, G., *et al.* Viral reorganization of the secretory pathway generates distinct organelles for RNA replication. *Cell* **141**(5), 799-811 (2010).
3. Toth, B., Balla, A., Ma, H., *et al.* Phosphatidylinositol 4-kinase IIIβ regulates the transport of ceramide between the endoplasmic reticulum and Golgi. *J. Biol. Chem.* **281**(47), 36369-36377 (2006).

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, 11/09/2022

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