

PRODUCT INFORMATION



CAY10669

Item No. 10974

CAS Registry No.: 1243583-88-1

Formal Name: 2-hydroxy-6-[(1Z)-2-[4-(pentyloxy)phenyl]ethenyl]-benzoic acid

MF: C₂₀H₂₂O₄

FW: 326.4

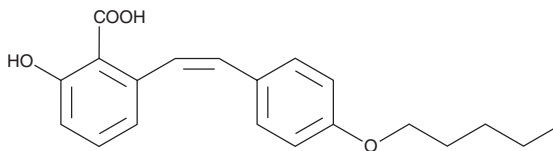
Purity: ≥98%

UV/Vis.: λ_{max}: 327 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

CAY10669 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10669 in the solvent of choice, which should be purged with an inert gas. CAY10669 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10669 in these solvents is approximately 30 mg/ml.

Description

CAY10669 is an inhibitor of the histone acetyltransferase PCAF (p300/CREB-binding protein-associated factor; IC₅₀ = 662 μM), displaying a 2-fold improvement in inhibitory potency over anacardic acid (Item No. 13144).¹ At 30-60 μM, CAY10669 can dose-dependently inhibit histone H4 acetylation in HepG2 cells *in vitro*.¹

Reference

1. Ghizzoni, M., Boltjes, A., Graaf, C., *et al.* Improved inhibition of the histone acetyltransferase PCAF by an anacardic acid derivative. *Bioorg. Med. Chem.* **18(16)**, 5826-5834 (2010).

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, 02/21/2024

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