

PRODUCT INFORMATION

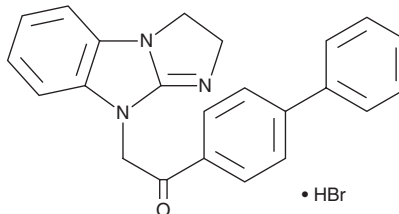


CCT031374 (hydrobromide)

Item No. 32866

CAS Registry No.: 1219184-91-4
Formal Name: 1-[1,1'-biphenyl]-4-yl-2-(2,3-dihydro-9H-imidazo[1,2-a]benzimidazol-9-yl)-ethanone, monohydrobromide

MF: C₂₃H₁₉N₃O • HBr
FW: 434.3
Purity: ≥98%
Supplied as: A 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

CCT031374 (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the CCT031374 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. CCT031374 (hydrobromide) is soluble in DMSO.

Description

CCT031374 is an inhibitor of β -catenin/transcription factor (TCF) complex signaling.¹ It inhibits TCF-dependent transcription induced by expression of Δ N- β -catenin, a constitutively active, stabilized form of β -catenin, but not expression of a constitutively active TCF-VP16 fusion protein in reporter assays at 30 μ M. CCT031374 (80 μ M) induces apoptosis in, and inhibits growth of, HCT116 colon cancer cells.

Reference

1. Ewan, K., Pajak, B., Stubbs, M., *et al.* A useful approach to identify novel small-molecule inhibitors of Wnt-dependent transcription. *Cancer Res.* **70(14)**, 5963-5973 (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, 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