

PRODUCT INFORMATION



ABC99

Item No. 25858

CAS Registry No.: 2331255-53-7
Formal Name: 7-(4-chlorobenzyl)-1,3-dioxohexahydroimidazo[1,5-a]pyrazin-2(3H)-yl 2,3-dihydro-4H-benzo[b][1,4]oxazine-4-carboxylate

MF: C₂₂H₂₁ClN₄O₅

FW: 456.9

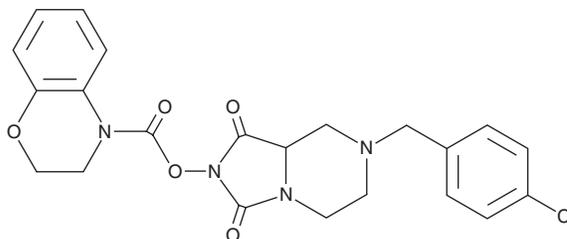
Purity: ≥95%

UV/Vis.: λ_{max}: 210, 282 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ABC99 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABC99 in the solvent of choice. ABC99 is soluble in organic solvents such as chloroform, acetonitrile, and methyl acetate, which should be purged with an inert gas. The solubility of ABC99 in these solvents is approximately 1 mg/ml.

Description

ABC99 is an N-hydroxyhydantoin (NHH) carbamate that inhibits the Wnt-deacylating enzyme NOTUM (IC₅₀ = 13 nM).¹ It preserves Wnt3A signaling in the presence of NOTUM (EC₅₀ = 89 nM in a cell-based reporter assay).

Reference

1. Suciu, R.M., Cognetta, A.B., III., Potter, Z.E., *et al.* Selective irreversible inhibitors of the Wnt-deacylating enzyme NOTUM developed by activity-based protein profiling. *ACS Med. Chem. Lett.* **9(6)**, 563-568 (2018).

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, 10/05/2020

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