

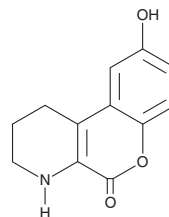
PRODUCT INFORMATION



CID-797718

Item No. 23428

CAS Registry No.: 370586-05-3
Formal Name: 1,2,3,4-tetrahydro-9-hydroxy-5H-[1]benzopyrano[3,4-b]pyridin-5-one
MF: C₁₂H₁₁NO₃
FW: 217.2
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 250, 347 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

CID-797718 is supplied as a crystalline solid. A stock solution may be made by dissolving the CID-797718 in the solvent of choice, which should be purged with an inert gas. CID-797718 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CID-797718 in these solvents is approximately 1, 20, and 30 mg/ml, respectively.

Description

CID-797718 is a protein kinase D1 (PKD1) inhibitor (IC₅₀ = 7 μM).¹ It also inhibits CDK activating kinase (CAK) and polo-like kinase 1 (PLK1) but has no activity at AKT (IC₅₀s = 8.4, 21.9, and >50 μM, respectively). CID-797718 inhibits PKD1 activity in LNCaP cells with an IC₅₀ value of 0.21 μM. CID-797718 is a byproduct in the synthesis of CID-755673 (Item No. 15924).²

References

1. Sharlow, E.R., Giridhar, K.V., LaValle, C.R., *et al.* Potent and selective disruption of protein kinase D functionality by a benzoxoloazepinolone. *J. Biol. Chem.* **283(48)**, 33516-33526 (2008).
2. Lavalle, C.R., Bravo-Altamirano, K., Gorodhar, K.V., *et al.* Novel protein kinase D inhibitors cause potent arrest in prostate cancer cell growth and motility. *BMC Chem Biol.* **10(5)**, (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, 04/13/2023

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