

PRODUCT INFORMATION



C-82

Item No. 42038

CAS Registry No.: 1422253-37-9
Formal Name: (6S,9S,9aS)-hexahydro-6-[(4-hydroxyphenyl)methyl]-2,9-dimethyl-4,7-dioxo-N-(phenylmethyl)-8-(8-quinolinylmethyl)-2H-pyrazino[2,1-c][1,2,4]triazine-1(6H)-carboxamide

MF: C₃₃H₃₄N₆O₄

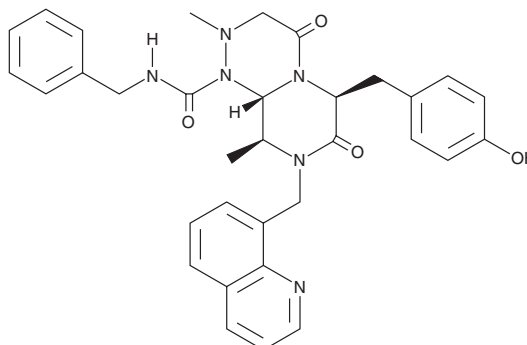
FW: 578.7

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

C-82 is supplied as a solid. A stock solution may be made by dissolving the C-82 in the solvent of choice, which should be purged with an inert gas. C-82 is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

C-82 is a β -catenin/CREB-binding protein (CBP) inhibitor and an active metabolite of PRI-724 (Item No. 41728).¹ It reduces the proliferation and protein levels of cyclin D1, α -smooth muscle actin (α -Sma), and Ki-67, a marker of cell proliferation, in primary mouse hepatic stellate cells (HSCs) when used at a concentration of 10 μ M. C-82 (10 μ M) decreases viability in continuous culture-activated primary mouse HSCs. It induces apoptosis when used in combination with nilotinib (Item Nos. 10010422 | 36628) in non-stem- and stem cell subpopulations of tyrosine kinase inhibitor-resistant, patient-derived, and blast crisis phase chronic myeloid leukemia (CML) cells with or without co-culture of primary human bone marrow mesenchymal stromal cells (MSCs).²

References

1. Osawa, Y., Oboki, K., Imamura, J., *et al.* Inhibition of cyclic adenosine monophosphate (cAMP)-response element-binding protein (CREB)-binding protein (CBP)/ β -catenin reduces liver fibrosis in mice. *EBioMedicine* **2**(11), 1751-1758 (2015).
2. Zhou, H., Mak, P.Y., Mu, H., *et al.* Combined inhibition of β -catenin and Bcr-Abl synergistically targets tyrosine kinase inhibitor-resistant blast crisis chronic myeloid leukemia blasts and progenitors *in vitro* and *in vivo*. *Leukemia* **31**(10), 2065-2074 (2017).

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, 11/18/2024

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