

PRODUCT INFORMATION

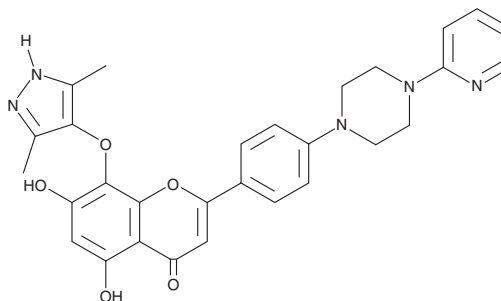


DC-U4106

Item No. 41723

CAS Registry No.: 2410534-62-0
Formal Name: 8-[(3,5-dimethyl-1H-pyrazol-4-yl)oxy]-5,7-dihydroxy-2-[4-[4-(2-pyridinyl)-1-piperazinyl]phenyl]-4H-1-benzopyran-4-one

MF: C₂₉H₂₇N₅O₅
FW: 525.6
Purity: ≥95%
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

DC-U4106 is supplied as a solid. A stock solution may be made by dissolving the DC-U4106 in the solvent of choice, which should be purged with an inert gas. DC-U4106 is slightly soluble (0.1-1 mg/ml) in the organic solvent DMSO.

Description

DC-U4106 is an inhibitor of ubiquitin-specific protease 8 (USP8; IC₅₀ = 1.2 μM).¹ It is selective for USP8 over USP2 and USP7 (IC₅₀s = 58.4 and >100 μM, respectively). DC-U4106 (1-5 μM) induces apoptosis in, and inhibits colony formation by, MCF-7 breast cancer cells. *In vivo*, DC-U4106 (20 mg/kg) reduces tumor volume in an MCF-7 mouse xenograft model.

Reference

1. Tian, Y., Liu, K., Liu, R., *et al.* Discovery of potent small-molecule USP8 inhibitors for the treatment of breast cancer through regulating ERα expression. *J. Med. Chem.* **65**(13), 8914-8932 (2022).

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, 10/22/2024

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