

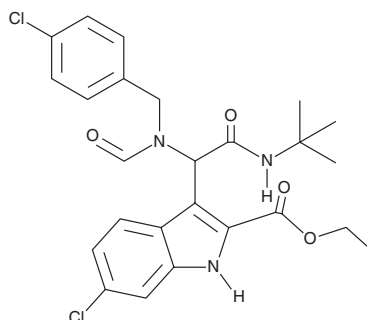
PRODUCT INFORMATION



YH 239-EE

Item No. 27678

CAS Registry No.: 1364488-67-4
Formal Name: 6-chloro-3-[1-[[[4-chlorophenyl)methyl]formylamino]-2-[(1,1-dimethylethyl)amino]-2-oxoethyl]-1H-indole-2-carboxylic acid, ethyl ester
MF: C₂₅H₂₇Cl₂N₃O₄
FW: 504.4
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 308 nm
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

YH 239-EE is supplied as a solid. A stock solution may be made by dissolving the YH 239-EE in the solvent of choice, which should be purged with an inert gas. YH 239-EE is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of YH 239-EE in these solvents is approximately 2, 5, and 10 mg/ml, respectively.

YH 239-EE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, YH 239-EE should first be dissolved in DMF and then diluted with the aqueous buffer of choice. YH 239-EE has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

YH 239-EE is an ethyl ester prodrug form of YH 239, a potent inhibitor of the p53-MDM2 protein-protein interaction.¹ YH 239-EE (20 μM) inhibits the growth of OCI-AML-3 cells and halts the cell cycle at the sub-G₁ phase and induces apoptosis in NB4 and MOLM-13 cells.

Reference

1. Huang, y., Wolf, S., Beck, B., *et al.* Discovery of highly potent p53-MDM2 antagonists and structural basis for anti-acute myeloid leukemia activities. *ACS Chem. Biol.* **9**(3), 802-811 (2014).

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, 12/02/2022

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