

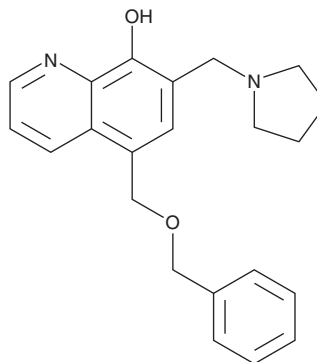
PRODUCT INFORMATION



UC-112

Item No. 45378

CAS Registry No.: 383392-66-3
Formal Name: 5-[(phenylmethoxy)methyl]-7-(1-pyrrolidinylmethyl)-8-quinolinol
MF: C₂₂H₂₄N₂O₂
FW: 348.4
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

UC-112 is supplied as a solid. A stock solution may be made by dissolving the UC-112 in the solvent of choice, which should be purged with an inert gas. UC-112 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

UC-112 is a Smac mimetic.¹ It reduces the level of survivin and induces apoptosis in A375 melanoma and PC3 prostate cancer cells and inhibits the growth of A375, M14 human melanoma, PC3 and DU145 prostate cancer cells (IC₅₀s = 1.6, 2.5, 0.7, and 3.4 μM, respectively). *In vivo*, UC-112 (20 and 40 mg/kg) reduces tumor growth and intratumoral levels of survivin in an A375 mouse xenograft model.

Reference

1. Wang, J. and Li, W. Discovery of novel second mitochondria-derived activator of caspase mimetics as selective inhibitor of apoptosis protein inhibitors. *J. Pharmacol. Exp. Ther.* **349**(3), 319-329 (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, 04/20/2026

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